

RNAscope™ VS Universal AP Assay

For Ventana DISCOVERY ULTRA System

Document Number UM 323250

For Research Use Only. Not for use in diagnostic procedures.

Trademarks

RNAscope, BaseScope, PretreatPro, and miRNAscope are trademarks of Advanced Cell Diagnostics, Inc. Lunaphore, seqIF and COMET logo are trademarks of Lunaphore. All other trademarks belong to their respective owners.

Citing RNAscope in Publications

When describing a procedure for publication using this product, please refer to it as the RNAscope Assay and cite: Wang F, Flanagan J, Su N, Wang L-C, Bui S, Nielson A, Wu X, Vo H-T, Ma X-J and Luo Y. RNAscope: A Novel *In Situ* RNA Analysis Platform for Formalin-Fixed Paraffin-Embedded Tissues. J. Mol. Diagnostics, 2012, 14:22–29.

Disclaimers

Advanced Cell Diagnostics, Inc. reserves the right to change its products and services at any time to incorporate technological developments. This manual is subject to change without notice.

Although this manual has been prepared with every precaution to ensure accuracy, Advanced Cell Diagnostics, Inc. assumes no liability for any errors, omissions, or for any damages resulting from the use of this information.

Copyright

© 2024. Advanced Cell Diagnostics, Inc. All rights reserved.

Contents

Chapter 1. Product Information	6
About this guide	6
Product description	6
Background	6
Overview	6
Kit contents and storage	7
RNAscope VS Probes.....	7
RNAscope VS Control Slides.....	8
RNAscope VS Universal Reagents.....	8
Required materials from Roche Diagnostics.....	9
Roche materials for Sequential RNA-Protein Co-Detection.....	10
Instrument buffers	12
User-supplied materials.....	12
Chapter 2. Before You Begin.....	14
Important procedural guidelines.....	14
Chapter 3. Prepare Samples	15
Workflow	15
Materials required	16
Fix and embed samples.....	16
Prepare FFPE tissue sections	16
Chapter 4. Determine Pretreatment Conditions	18
Pretreat FFPE sections.....	18
Pretreatment workflow selection	18
Target retrieval and permeabilization.....	18
Tissue pretreatment recommendations for the protease-free workflow	19
Tissue pretreatment recommendations for the protease workflow	19
Chapter 5. Automated RNAscope VS Universal AP Assay	20
Workflow	20
Prepare materials	20
Materials required.....	20
Prepare the instrument.....	21
Dilute bulk reagents.....	21
Register new reagents	21
Prepare instrument reagents	21
Create an instrument protocol.....	22
Print the labels	25

Load the reagents	25
Complete the run.....	26
Wash and dry the slides	26
Mount the samples.....	26
Chapter 6: Protease-Free, Sequential RNA-Protein Co-Detection.....	28
Workflow	28
Prepare materials	28
Materials required.....	28
Prepare the instrument	29
Dilute instrument bulk reagents	29
Register new reagents.....	29
Prepare user-fillable reagents for RNA-Protein Co-Detection.....	29
Create an instrument protocol.....	30
Print the labels	36
Load the reagents	36
Complete the run.....	37
Wash and dry the slides	37
Mount the samples.....	38
Chapter 7. Evaluate the Results.....	39
Evaluate the samples	39
Troubleshooting.....	39
Tissue example	40
Appendix A. Semi-automated RNAscope VS Universal AP Assay	41
Workflow	41
Kit contents and storage	42
Prepare the materials.....	42
Materials required	42
Prepare the instrument.....	42
Dilute bulk reagents.....	43
Register new reagents.....	43
Prepare instrument reagents.....	43
Prepare deparaffinization reagents	43
Prepare 1X Target Retrieval	43
Create an instrument protocol.....	44
Print the labels	45
Manually pretreat the samples	45
Bake the slides	45
Deparaffinize FFPE sections	45
Pretreat the slides	46
Run the RNAscope VS Universal AP Assay	47
Load the reagents	47

Start the run	47
Complete the run.....	48
Wash the slides	48
Dry and mount the samples.....	48
Appendix B. Pretreatment Guidance for FFPE Samples	49
Tissue-specific pretreatment conditions	49
Appendix C. Safety.....	50
Chemical safety	50
Biological hazard safety	50
In the U.S.:	51
In the EU:.....	51
Documentation and Support	52
Obtaining SDSs	52
Obtaining support	52
Contact information	52
Limited product warranty	52

1

Chapter 1. Product Information



Before using this product, read and understand the information in **Appendix C. Safety** of this document.

IMPORTANT! We recommend reading the entire user manual before beginning any protocols.

About this guide

This user manual provides several versions of the RNAscope VS Universal AP Assay:

- Chapter 5. Automated RNAscope VS Universal AP Assay
- Chapter 6. Protease-Free, Sequential RNA-Protein Co-Detection
- Appendix A. Semi-automated RNAscope VS Universal AP Assay

Product description

Background

The RNAscope VS Universal Assays uses a novel and proprietary method of in situ hybridization (ISH) to visualize single RNA molecules per cell in formalin-fixed, paraffin-embedded (FFPE) tissue mounted on slides. The assays are based on Advanced Cell Diagnostic's patented signal amplification and background suppression technology and can detect RNA molecules in archival samples and partially degraded specimens. The RNAscope VS Universal Assay allows users to automate the highly sensitive RNAscope Assay using the Ventana DISCOVERY ULTRA System.

Overview

Figure 1 illustrates the RNAscope VS Universal Assay procedure, which can be completed on the instrument in ~11 hours. Starting with properly prepared samples, sections are first pretreated, and then RNA-specific probes are hybridized to target RNA. The signal is amplified using multiple steps, followed by hybridization to alkaline phosphatase (AP)-labeled probes and detection using a chromogenic substrate. A single RNA transcript appears as a distinct dot of chromogen precipitate visible using a common bright-field microscope.

Figure 1. Procedure overview



1. Tissue section	2. Hybridize to target RNA	3. Amplify	4. Image
Start with properly prepared sections and pretreat to allow access to target RNA.	Hybridize gene-specific probe pairs to the target mRNA.	Probes are hybridized to a cascade of signal amplification molecules, culminating in binding of AP-labeled probes. Add Fast Red substrate to detect target RNA.	Visualize target RNA using a standard bright field microscope.

Kit contents and storage

The RNAscope VS Universal Assay requires the RNAscope 2.5 VS Probes and the RNAscope VS Universal Reagents, available from Advanced Cell Diagnostics.

RNAscope VS Probes

The RNAscope 2.5 VS Probes consist of the user-specified Target Probe and the Positive and Negative Control Probes. Visit <https://acdbio.com/products> to find a gene-specific Target Probe or appropriate Control Probes. Each probe is sufficient for staining ~30 standard slides. The probes have a shelf life of two years from the manufacturing date when stored as indicated in the following table:

Target Probes					
<input checked="" type="checkbox"/>	Reagent	Cat. No.	Content	Quantity	Storage
	RNAscope 2.5 VS Target Probe ([species] – [gene])	Various	Probe targeting specific RNA	7 mL x 1 bottle	2–8°C
Control Probes					
<input checked="" type="checkbox"/>	Reagent	Cat. No.	Content	Quantity	Storage
	RNAscope 2.5 VS Positive Control Probe – [species] – PPIB)	Various	Probe targeting common housekeeping gene	7 mL x 1 bottle	2–8°C
	RNAscope 2.5 VS– Negative Control Probe – DapB	312039	Probe targeting bacterial gene dapB	7 mL x 1 bottle	2–8°C

RNAScope VS Control Slides

The RNAScope VS Control Slides (Cat. No. 310045 for Human control slide, HeLa; Cat. No. 310023 for Mouse control slide, 3T3) contain FFPE cell pellets sectioned and mounted on slides. The control slides can be used for assay control with the RNAScope 2.5 VS Positive Control Probe and RNAScope 2.5 VS Negative Control Probe. The slides have a shelf life listed on the label when stored at 2–8°C with desiccants.

RNAScope VS Universal Reagents

RNAScope VS Universal kits provide enough reagents to stain ~60 standard slides. You will receive four boxes when you order the RNAScope VS Universal AP Reagent Kit (Cat. No. 323250).

RNAScope VS Universal AP Reagents include:

- RNAScope VS Universal AP Detection Reagents (Cat. No. 323260)
 - RNAScope VS Universal AP Standard Reagents
 - RNAScope VS Pro Reagents
- RNAScope VS Universal Sample Prep Reagents v2 (Cat. No. 323740)
- RNAScope VS Accessory Kit (Cat. No. 320630)

The reagents are Ready-To-Use (RTU) and have shelf life listed on the label when stored as indicated in the following table:

RNAScope VS Universal AP Detection Reagents (Cat. No. 323260)				
RNAScope VS Universal AP Standard Reagents (Cat. No. 322040)				
<input checked="" type="checkbox"/>	Cat. No.	Reagent	Quantity	Storage
	323261	RNAScope VS Universal AP AMP 1	14 mL x 1 bottle	2–8°C
	323262	RNAScope VS Universal AP AMP 2	14 mL x 1 bottle	2–8°C
	323263	RNAScope VS Universal AP AMP 3	14 mL x 1 bottle	2–8°C
	323264	RNAScope VS Universal AP AMP 4	14 mL x 1 bottle	2–8°C
	323265	RNAScope VS Universal AP AMP 5	14 mL x 1 bottle	2–8°C
	323266	RNAScope VS Universal AP AMP 6	14 mL x 1 bottle	2–8°C
	323267	RNAScope VS Universal AP AMP 7	14 mL x 1 bottle	2–8°C
	323218	RNAScope VS Protease	14 mL x 1 bottle	2–8°C
RNAScope VS Pro Reagents (Cat. No. 322035)				
<input checked="" type="checkbox"/>	Cat. No.	Reagent	Quantity	Storage
	322031	RNAScope VS PretreatPro	18 mL x 1 bottle	2–8°C
	322032	RNAScope VS CoDetectPro	18 mL x 1 bottle	2–8°C

RNAscope VS Universal Sample Prep Reagents v2 (323740)				
<input checked="" type="checkbox"/>	Cat. No.	Reagent	Quantity	Storage
	323221	RNAscope VS Target Retrieval v2	14 mL x 2 bottles	Room Temp (15–30°C)
	323222	RNAscope VS Dewax	14 mL x 1 bottle	Room Temp (15–30°C)
RNAscope VS Accessory Kit (Cat. No. 320630)				
<input checked="" type="checkbox"/>	Cat. No.	Reagent	Quantity	Storage
	320631	RNAscope VS Hematoxylin — RTU	7 mL x 1 bottle	2–8°C
	320632	RNAscope VS Bluing Reagent — RTU	7 mL x 1 bottle	2–8°C

IMPORTANT! Ensure Dewax is in solution and at room temperature before use on the instrument. Warm the solution in your hand or place it at 37°C for 15 MIN prior to each use, regardless of previous storage conditions, as it may precipitate during shipment.

IMPORTANT! Do not interchange the reagent components of the reagent kits, even those having the same name.

Required materials from Roche Diagnostics

The RNAscope VS Universal Assay requires specific materials and equipment available only from Roche Diagnostics. Catalog numbers are valid in the U.S. only. For other regions, please check catalog or ordering numbers with your local lab supplier.

Probe Dispensers (Cat. No. 960-761 to 960-780; for ordering code, contact local Roche representative)		
<input checked="" type="checkbox"/>	Component	Storage
	250 Test Probe #1–20 dispensers — fill dispensers with RNAscope 2.5 VS Probes. Use up to 20 probes at one time.	Room Temp (15–30°C)
mRNA Sample Prep Kit (Cat. No. 760-248; Ordering Code 08127166001)		
<input checked="" type="checkbox"/>	Component	Storage
	mRNA Target Retrieval dispenser — fill dispenser with RNAscope VS Universal Target Retrieval v2 (Optional: This dispenser is not needed for Protease-Free Workflow)	Room Temp (15–30°C)
	mRNA Dewax — fill dispenser with RNAscope VS Dewax	Room Temp (15–30°C)
	mRNA Protease dispenser — fill dispenser with RNAscope VS Protease (Optional: This dispenser is not needed for Protease-Free Workflow)	Room Temp (15–30°C)
mRNA RED Probe Amplification Kit (Cat. No. 760-236; Ordering Code 7095341001)		
<input checked="" type="checkbox"/>	Component	Storage
	mRNA AMP 1 dispenser — fill dispenser with VS Universal AP AMP 1	Room Temp (15–30°C)
	mRNA AMP 2 dispenser — fill dispenser with VS Universal AP AMP 2	Room Temp (15–30°C)
	mRNA AMP 3 dispenser — fill dispenser with VS Universal AP AMP 3	Room Temp (15–30°C)
	mRNA AMP 4 dispenser — fill dispenser with VS Universal AP AMP 4	Room Temp (15–30°C)
	mRNA AMP 5 dispenser — fill dispenser with VS Universal AP AMP 5	Room Temp (15–30°C)

	mRNA AMP 6 dispenser — fill dispenser with VS Universal AP AMP 6	Room Temp (15–30°C)		
	mRNA AMP 7 dispenser — fill dispenser with VS Universal AP AMP 7	Room Temp (15–30°C)		
mRNA RED Detection Kit (Cat. No. 760-234; Ordering Code 7099037001)				
<input checked="" type="checkbox"/>	Component	Storage		
	mRNA Inhibitor-prefilled	2–8°C		
	mRNA Activator dispenser-prefilled	2–8°C		
	mRNA Naphthol dispenser-prefilled	2–8°C		
	mRNA Fast Red dispenser-prefilled	2–8°C		
Additional Dispensers for RNAscope				
<input checked="" type="checkbox"/>	Component	Cat. No.	Ordering Code	Fill with:
	Pretreatment dispenser*	960-901 to 960-909	Contact local Roche representative	VS PretreatPro
	Counterstain 1 dispenser	771-741	05271720001	VS Hematoxylin
	Counterstain 2 dispenser	771-742	05271738001	VS Bluing Reagent

*Choose distinct barcode numbers for VS PretreatPro and VS CoDetectPro. Ensure that staining protocols reflect the barcode numbers assigned to each reagent.

Roche materials for Sequential RNA-Protein Co-Detection

Sequential RNA-Protein Co-Detection can be performed with either Roche RTU primary antibody or with your choice of primary antibody concentrate diluted for use on the DISCOVERY ULTRA. For a list of available Roche antibodies and ordering information, please contact your local Roche representative.

The following additional materials from Roche Tissue Diagnostics can be used for Sequential RNA-Protein Co-Detection on Ventana DISCOVERY ULTRA.

IHC assay-specific materials

Reagent Options for Ventana IHC AP Detection				
<input checked="" type="checkbox"/>	Component	Cat. No.	Ordering Code	Storage
	DISCOVERY UltraMap anti-Ms Alk Phos*	760-4312	05269687001	2–8°C
	DISCOVERY UltraMap anti-Rb Alk Phos*	760-4314	05269709001	2–8°C
	DISCOVERY Yellow Kit*	760-239	07698445001	2–8°C
	DISCOVERY Red Kit	760-228	07425333001	2–8°C
	DISCOVERY ChromoMap Red Kit*	760-160	05266653001	2–8°C

Reagent Options for Ventana IHC HRP Detection				
<input checked="" type="checkbox"/>	Product	Cat. No.	Ordering Code	Storage
	DISCOVERY UltraMap anti-Ms HRP*	760-4313	05269695001	2–8°C
	DISCOVERY UltraMap anti-Rb HRP*	760-4315	05269717001	2–8°C
	DISCOVERY UltraMap anti-Rat HRP*	760-4456	05891884001	2–8°C
	DISCOVERY UltraMap anti-Gt HRP*	760-4648	06607241001	2–8°C
	DISCOVERY Purple Kit*	760-229	07053983001	2–8°C
	DISCOVERY Green HRP Kit †	760-278	07053983001	2–8°C
	DISCOVERY Teal HRP Kit †	760-247	08254338001	2–8°C
	DISCOVERY ChromoMap DAB Kit †	760-159	05266645001	2–8°C
	DISCOVERY DCC Kit †	760-240	07988192001	2–8°C
	DISCOVERY FAM Kit †	760-243	07988150001	2–8°C
	DISCOVERY FITC Kit†	760-232	07259212001	2–8°C
	DISCOVERY Rhodamine Kit †	760-233	07259883001	2–8°C
	DISCOVERY Rhodamine 6G Kit †	760-244	07988168001	2–8°C
	DISCOVERY Red 610 Kit†	760-245	07988176001	2–8°C
	DISCOVERY Cy5 Kit †	760-238	07551215001	2–8°C

* Choose one secondary detection antibody depending on the primary antibody species and desired IHC chromogen.

† Choose one IHC chromogen per protein target.

Dispensers for Sequential RNA-Protein Co-Detection				
<input checked="" type="checkbox"/>	Component	Cat. No.	Ordering Code	Fill with:
	Antibody dispensers (<i>Optional*</i>)	770-001 to 770-099	Contact local Roche representative	User-sourced Primary Antibody Concentrate diluted in Co-Detection Antibody Diluent
	Pretreatment dispenser†	960-901 to 960-909	Contact local Roche representative	VS PretreatPro
	Pretreatment dispenser‡	960-901 to 960-909	Contact local Roche representative	VS CoDetectPro
	Counterstain 1 dispenser	771-741	05271720001	VS Hematoxylin
	Counterstain 2 dispenser	771-742	05271738001	VS Bluing Reagent
	Roche RTU Counterstain Reagents	Various	Contact local Roche representative	Pre-filled
	DISCOVERY Antibody Block	760-4204	05268869001	Pre-filled

*Sequential RNA-Protein Co-Detection can be performed with either Roche RTU primary antibody or your choice of primary antibody.

†Choose distinct barcode numbers for VS PretreatPro and VS CoDetectPro. Ensure that staining protocols reflect the barcode numbers assigned to each reagent.

Instrument buffers

<input checked="" type="checkbox"/>	Component	Cat. No./Ordering Code
	10X DISCOVERY Wash (RUO)	950-510 / 07311079001
	ULTRA LCS (Predilute)	650-210 / 05424534001
	SSC (10X)	950-110 / 5353947001
	Reaction Buffer (10X)	760-107 / 5266262001
	DISCOVERY CC1	950-500 / 6414575001

IMPORTANT! To run the VS Universal assay successfully, use DISCOVERY Wash (950-510). Do not use DISCOVERY EZ Prep. Place 2X SSC (950-110) in the SSC bulk container. You may fill the option bulk container with reaction buffer.

User-supplied materials

IMPORTANT! Do not substitute other materials for the SuperFrost® Plus Slides listed in the following table.

<input checked="" type="checkbox"/>	Description	Supplier	Cat. No.
	SuperFrost Plus Slides (required)	Fisher Scientific	12-550-15
	100% ethanol (EtOH)	American Master Tech Scientific/MLS	ALREAGAL
	Xylene (or Histo-Clear)	Fisher Scientific/MLS	X3P-1GAL
	10% neutral-buffered formalin (NBF)	MLS	—
	Paraffin wax	MLS	—
	1X PBS	MLS	—
	Microtome	MLS	—
	Drying oven, capable of holding temperature at 60 +/- 1°C	MLS	—
	EcoMount (if using Red detection)	Biocare or ACD	EM897L or 320409
	Tissue-Tek® Vertical 24 Slide Rack	American Master Tech Scientific/MLS	LWSRA24
	Tissue-Tek Staining Dishes	American Master Tech Scientific/MLS	LWT4457EA
	Tissue-Tek Clearing Agent Dishes, xylene resistant	American Master Tech Scientific/MLS	LWT4456EA
	Cover Glass 24 x 50 mm	Fisher Scientific/MLS	12-545F
	Distilled water	MLS	—
	Dawn detergent or similar detergent	MLS	—
	Fume hood	MLS	—

<input checked="" type="checkbox"/>	Description	Supplier	Cat. No.
	Optional: Glass beaker (1 or 2 L)	MLS	—
	Optional: Hot plate	Fisher Scientific/MLS	11-300-49SHP
	Primary antibody – RTU or concentrate	User	Various
	ProLong Gold Antifade Reagent	Life Technologies	P36930

* Major Laboratory Supplier in North America. For other regions, please check catalog numbers with your local lab supplier.

2

Chapter 2. Before You Begin

Prior to running the RNAscope VS Universal Assay on your samples for the first time, we recommend that you:

- Be familiar with the Ventana DISCOVERY ULTRA system. Refer to the *Ventana DISCOVERY ULTRA System User Manual*.
- Run the assay on FFPE RNAscope VS Control Slides (Cat. No. 310045 for Human control slide, HeLa; Catalog No. 310023 for Mouse control slide, 3T3) using the Positive and Negative Control Probes.

Important procedural guidelines

- Start with properly fixed and prepared sections. Refer to **Chapter 3. Prepare and Pretreat Samples, Recommended guidelines** and to our sample preparation and pretreatment user guides available at <https://acdbio.com/technical-support/user-manuals>.
- Regularly maintain and clean your automated staining instrument.
- Always run positive and negative control probes on your sample to assess sample RNA quality and background, respectively.
- Do not substitute required materials. Assay has been qualified with these materials only.
- Follow the protocol exactly for the best results. Do not substitute buffers or diluents.
- Store reagents accordingly right after use. To avoid any contamination or evaporation, do not leave reagents open for an extended period.
- Do not let your sections dry out during the procedure unless specified in the protocol.
- Use good laboratory practices and follow all necessary safety procedures. Refer to **Appendix C. Safety** for more information.

3

Chapter 3. Prepare Samples

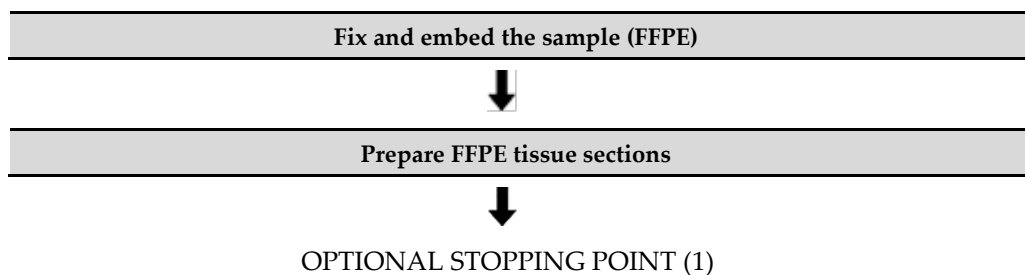
This chapter describes the formalin-fixed, paraffin-embedded (FFPE) sample preparation method and slide pre-processing procedure. If you have already embedded your FFPE samples, you can proceed directly to the **Prepare FFPE tissue sections** procedure.

For other sample types and preparation methods, contact support.acd@bio-techne.com for the latest protocols and guidelines.

IMPORTANT! We highly recommend following these guidelines. We cannot guarantee assay results with other preparation methods.

For samples treated differently from the following protocol, you may need to optimize pretreatment conditions. Contact support.acd@bio-techne.com if you need additional guidance.

Workflow



Materials required

-
- 10% Neutral Buffered Formalin (NBF)
 - 1X PBS
 - Paraffin wax
 - Tissue-Tek Clearing Agent Dishes
 - Tissue-Tek Staining Dishes
 - Tissue-Tek Vertical 24 Slide Rack
 - 100% alcohol (EtOH)
 - Histo-Clear or xylene
 - Microtome
 - Water bath
 - SuperFrost Plus slides
 - Drying oven
 - Fume hood
-

Fix and embed samples

1. Immediately following dissection, fix tissue in 10% NBF for **16–32 HRS at ROOM TEMPERATURE (RT)**. Fixation time will vary depending on tissue type and size.

 **CAUTION!** Handle biological specimens appropriately.

IMPORTANT! Fixation for <16 HRS or >32 HRS will impair the performance of the assay.

2. Wash sample with 1X PBS.
3. Dehydrate sample using a standard ethanol series, followed by xylene.

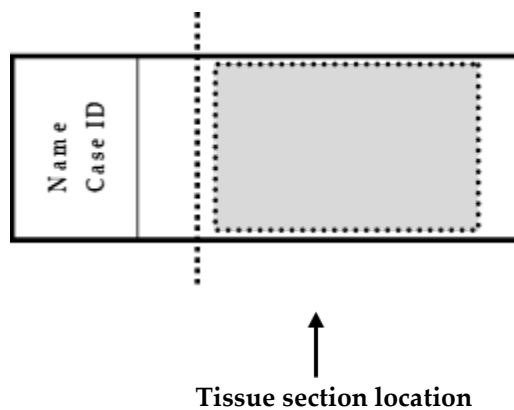
IMPORTANT! Use fresh reagents. Embed samples as quickly as possible to preserve RNA quality.

4. Embed sample in paraffin using standard procedures.

Note: Embedded samples can be stored at room temperature with desiccants. For best preservation of RNA quality over an extended period (>1 yr), storage at 2–8°C with desiccants is recommended.

Prepare FFPE tissue sections

1. Trim paraffin blocks as needed and cut embedded tissue into 5 +/- 1 µm sections using a microtome.
2. Place paraffin ribbon in a **40–45 °C** water bath and mount the sections on **SUPERFROST PLUS SLIDES**. Place tissue as shown for optimal staining:



IMPORTANT! Do not mount more than one section per slide. Place sections in the center of the slide.

3. Air dry the slides **OVERNIGHT** at **RT**. Do NOT bake slides unless they will be used for RNAscope within 1 week.

OPTIONAL STOPPING POINT (1). You can store sections with desiccants at room temperature. For best preservation of RNA quality, storage at 2–8°C with desiccants is recommended. Use sectioned tissue within 3 months.

4

Chapter 4. Determine Pretreatment Conditions

The following protocols describe sample pretreatment for formalin-fixed, paraffin-embedded (FFPE) samples. For other sample types and preparation methods, contact support.acd@biotechne.com for the latest protocols and guidelines.

IMPORTANT! We highly recommend following these guidelines. We cannot guarantee assay results with other preparation methods.

Pretreat FFPE sections

Pretreatment workflow selection

FFPE samples must be de-crosslinked with a target retrieval step, and then permeabilized for best RNA access. The RNAscope VS Universal RNAscope AP Assay is compatible with both protease-based and protease-free pretreatment approaches. We recommend using a protease-free pretreatment if you plan to combine RNA with protein detection in future downstream applications.

Target retrieval and permeabilization

Two options are available:

- Protease-free permeabilization is recommended when there is interest in combining RNA with Protein detection downstream. This approach uses an IHC-style cell conditioning for Target Retrieval, using Roche CC1 bulk reagent, combined with permeabilization using ACD's VS PretreatPro reagent. This allows co-detection of RNA and proteins that were previously incompatible with protease on the same tissue section using the same cell conditioning as immunohistochemistry (IHC). We recommend the protease-free pretreatment to be considered as the default pretreatment.
- Protease-based pretreatment is recommended for experiments that stain only RNA for established protocols when maintenance of performance compared to historical ISH experiments is critical. This option uses ACD Target Retrieval v2 and VS Protease.

Tissue pretreatment recommendations for the protease-free workflow

Start with the following conditions when tissues are prepared as described in **Chapter 3**. Depending on the tissue type, you may need to adjust the time and temperature of CC1 to maximize the positive RNA control signal while minimizing or eliminating the negative RNA control signal.

Recommended pretreatment selections:

Reagent	Mild	Standard	Extended
CC1 Cell Conditioning – Protease-Free	32 MIN @95C	64 MIN @95C	92 MIN @95C
VS PretreatPro	32 MIN @40C	32 MIN @40C	32 MIN @40C

*Sample types, such as certain xenografts and cell pellets, might require shorter incubation time. For these tissue types, the CC1 time and or temperature can be reduced. VS PretreatPro incubation time is adjustable but is rarely needed. Refer to the table below for recommendations for optimization and to **Appendix B** for recommended pretreatment strength by tissue type. If you have a tissue type not listed, contact ACD Support at support.acd@bio-technne.com.

Optimization guidance (if needed):

Optimization	
Troubleshooting:	Recommended Optimization
Tissue detachment	Reduce CC1 Target Retrieval Time
Low ISH signal	Extend CC1 Target Retrieval Time
High dot background	Reduce CC1 Target Retrieval Time
Low staining patches	Reduce CC1 Target Retrieval Temperature

Tissue pretreatment recommendations for the protease workflow

Start with the following conditions when tissues are prepared as described in **Chapter 3**. Depending on the tissue type, you may need to adjust the time and temperature of Target Retrieval v2 (Cell Conditioning) to maximize the positive RNA control signal while minimizing or eliminating the negative RNA control signal.

Recommended pretreatment selections:

Reagent	Mild	Standard	Extended
ACD VS Target Retrieval v2	16 MIN @97C	24 MIN @97C	48 MIN @97C
RNAScope VS Protease	16 MIN @37C	16 MIN @37C	16 MIN @37C

*Sample types, such as certain xenografts and cell pellets, might require shorter incubation time. For these tissue types, reduce the target retrieval incubation time. RNAScope VS Protease incubation time can also be adjusted but is rarely needed. Refer to **Appendix B** for recommended pretreatment strength by tissue type. If you have a tissue type not listed, contact ACD Support at support.acd@bio-technne.com.

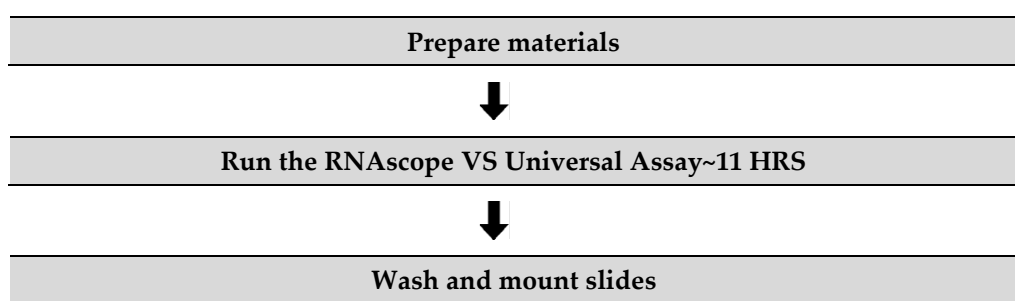
5

Chapter 5. Automated RNAscope VS Universal AP Assay

IMPORTANT! We strongly recommend you run the RNAscope VS Control Slides (Cat. No. 310045 or 310023) using the RNAscope 2.5 VS Positive and Negative Control Probes along with your samples in every run.

Appendix A. Semi-automated RNAscope VS Universal AP Assay describes an offline boiling procedure for use with Cat. No. 322000 (RNAscope Target Retrieval Reagents).

Workflow



Prepare materials

Materials required

Materials Provided by Advanced Cell Diagnostics	Materials Provided by Roche Tissue Diagnostics	Other Materials and Equipment
<ul style="list-style-type: none"> • RNAscope 2.5 VS Target Probe • RNAscope 2.5 VS Positive Control Probe • RNAscope 2.5 VS Negative Control Probe • RNAscope VS Dewax • RNAscope VS Protease† • VS PretreatPro* • RNAscope VS Target Retrieval† • RNAscope VS Universal AP AMP 1 • RNAscope VS Universal AP AMP 2 • RNAscope VS Universal AP AMP 3 • RNAscope VS Universal AP AMP 4 • RNAscope VS Universal AP AMP 5 	<ul style="list-style-type: none"> • DISCOVERY ULTRA — automated slide stainer • DISCOVERY Wash Buffer 10X • ULTRA LCS (Predilute) • SSC Buffer 10X • DISCOVERY CC1 • Reaction Buffer 10X • Probe dispensers • mRNA RED Probe Amplification Kit • mRNA Universal Sample Prep Kit • mRNA RED Detection Kit • User fillable dispensers 	<ul style="list-style-type: none"> • Distilled water • Dawn detergent or similar detergent • Fume hood • Xylene • Tissue-Tek Staining Dish • Tissue-Tek Clearing Agent Dish, xylene-resistant • Tissue-Tek Vertical 24 Slide Rack • EcoMount • Cover Glass, 24 mm x 50 mm

Materials Provided by Advanced Cell Diagnostics	Materials Provided by Roche Tissue Diagnostics	Other Materials and Equipment
<ul style="list-style-type: none"> • RNAscope VS Universal AP AMP 6 • RNAscope VS Universal AP AMP 7 • RNAscope VS Hematoxylin • RNAscope VS Bluing Reagent 		

* This dispenser is only used for the protease-free workflow

† These dispensers are used for the protease workflow

Prepare the instrument

Most tissue types can be fully automated using the DISCOVERY ULTRA Kits. Manual pretreatment may give a better result in some cases. Use the semi-automated procedure in **Appendix A** for tissues that do not have a satisfactory result when using the fully automated procedure.

If the instrument has not been used for ≥ 1 week, follow the guidelines for instrument maintenance in the *Ventana DISCOVERY ULTRA System User Manual*.

Dilute bulk reagents

Please prepare the bulk fluids per manufacturer instructions.

Register new reagents

Reagent dispensers come with appropriate barcode labels and registration buttons for dispensing RNAscope VS Universal Reagents. Refer to the *Ventana DISCOVERY ULTRA System User Manual* for details. To register reagents:

1. Log all ACD reagents and probes into the software as “log user-fillable reagents” or “log user-fillable probes”.
2. Use the wand or barcode scanner that comes with the instrument to register new reagent kits.

Prepare instrument reagents

After selecting the pretreatment workflow, refer to the dispenser table in **Chapter 1** to determine the proper dispenser for each reagent.

1. Transfer VS Protease and both bottles of VS Target Retrieval v2 **OR** VS PretreatPro to the corresponding labeled dispenser.
2. For RNAscope VS Universal AP reagents AMP 1 to AMP 7, transfer the entire volume of each component into the correspondingly labeled dispenser.
3. Transfer the VS Dewax, VS Hematoxylin, and VS Bluing Reagent to the correspondingly labeled dispenser.
4. Use probe dispensers for 2.5 VS Target Probe, 2.5 VS Positive Control Probe, 2.5 VS Negative Control Probe.

IMPORTANT! Avoid cross contamination between reagents. Dewax must be warmed to room temperature and be completely in solution before use.

- Follow the dispenser product insert instructions to properly prime and handle the dispensers.
- Store the tightly-capped mRNA Dewax and Target Retrieval v2 dispensers at **15–30°C**. Store all other capped dispensers at **4°C** when not in use.
- Check solution levels: DISCOVERY Wash, 2X SSC, Reaction Buffer, ULTRA LCS (Predilute), and CC1. Refill if they are less than half full. To run a full tray of 30 slides, completely fill the 2X SSC and Reaction Buffer containers.

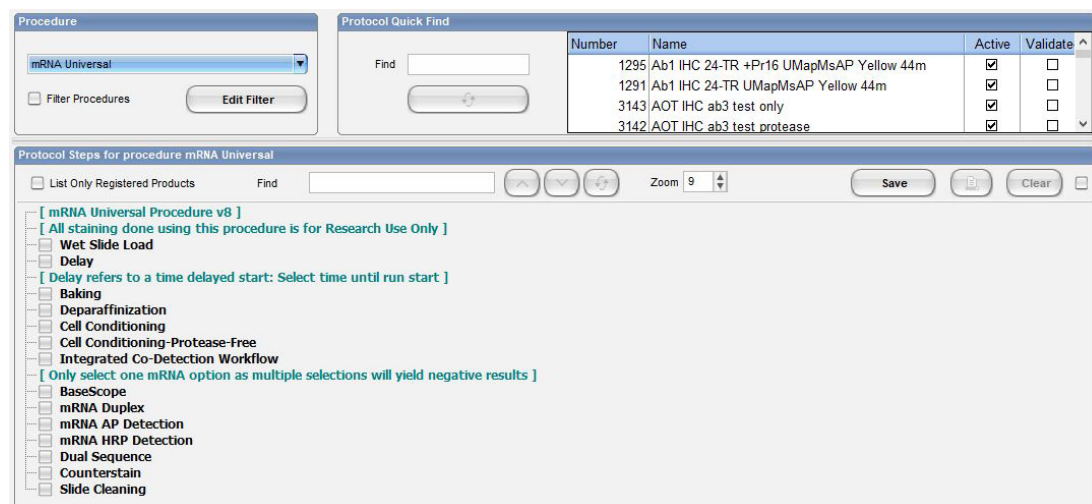
IMPORTANT! Do not use expired reagents.

- Empty the waste bottle if needed.

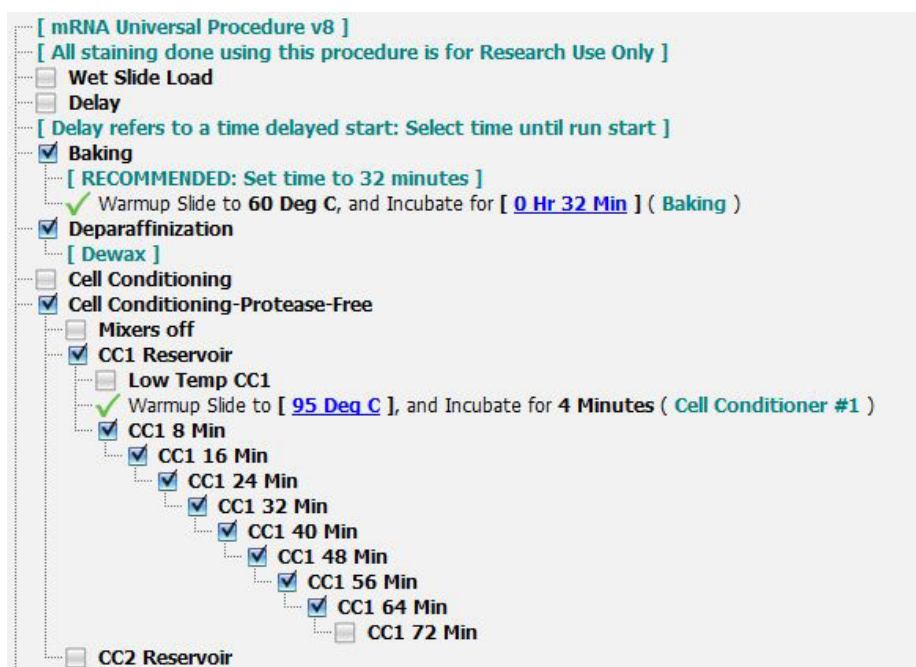
Create an instrument protocol

- Open the VS software and click on the **Protocol** button.
- Click on **Create/Edit Protocols**, go to the Procedure drop down menu and select **mRNA Universal**.

Main protocol steps appear as shown:



- Do one of the following, depending on your selected workflow:
 - For target retrieval using the **Protease-Free Workflow**, select the appropriate pretreatment conditions according to the following screenshot. After selecting the main steps, drop-down menus become available. We recommend selecting **Cell-Conditioning-Protease-Free** with CC1 for the time and temperature specified in **Chapter 4** according to tissue-specific recommendations found in **Appendix B**.



Note: The duration of tissue cell conditioning corresponds to the longest selected time. It is not the sum of all selected times.

- For target retrieval using the **Protease Workflow**, select the appropriate pretreatment conditions according to the following screenshot. After selecting the main steps, drop-down menus become available. We recommend selecting **Cell Conditioning** for the time and temperature specified in **Chapter 4** according to tissue-specific recommendations found in **Appendix B**.

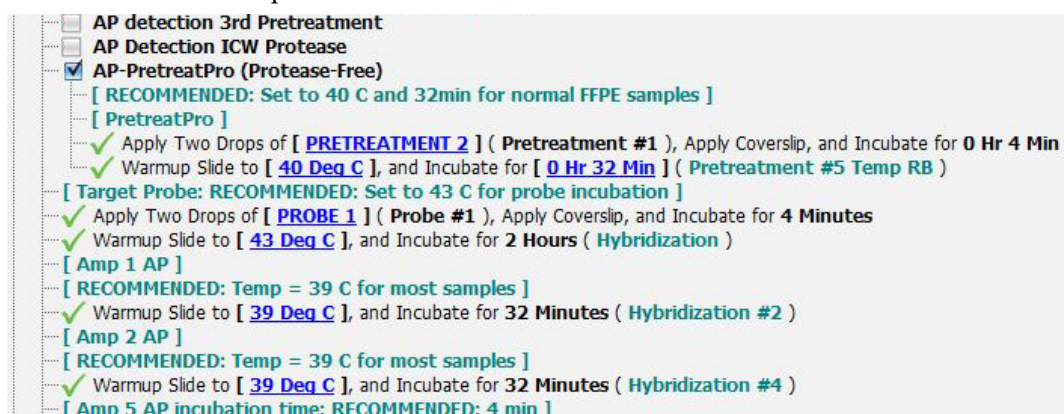


Note: The duration of tissue cell conditioning corresponds to the longest selected time. It is not the sum of all selected times.

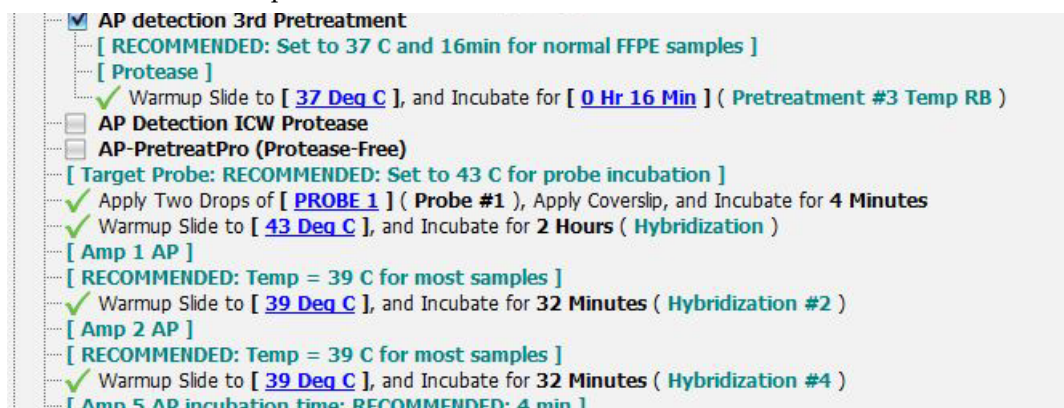
- For ISH staining with RNAscope VS Universal AP, select **mRNA AP Detection**.

Note: Select only one mRNA option per protocol. Choosing multiple mRNA options gives negative results.

5. Do one of the following, depending on your selected workflow:
 - If using the **Protease-Free Workflow**, select **AP-PretreatPro** to apply and incubate the VS PretreatPro reagent. We recommend incubating this reagent for **32 MIN** at **40°C**.
 - If using the **Protease Workflow**, select **AP detection 3rd Pretreatment** to apply and incubate the VS Protease reagent. We recommend incubating this reagent for **16 MIN** at **37°C**.
6. Select the remaining ISH staining conditions according to the following screenshots and table, depending on your selected workflow.
 - ISH Selections for protease-free workflow:



- ISH Selections for protease workflow:



Standard Times/Temperatures for mRNA AP Detection	
Cell Conditioning*	24 MIN @97°C
Cell-Conditioning-Protease-Free (CC1)*	64 MIN @95°C
VS Protease temperature and time*	37°C, 16 MIN
AP-PretreatPro (protease-free) temperature and time*	40°C, 32 MIN
Suggested probe temperature	Single Probes - 43°C HPV Pooled Probes - 50°C
Suggested RNAscope AP AMP 1 and AMP 2 temperature	39°C†
RNAscope AP AMP 5 incubation time	Instrument titrated ‡

*Choose either Cell Conditioning/VS Protease or Cell-Conditioning–Protease-Free/VS PretreatPro depending on the workflow. Do not select both within the same protocol. Refer to **Chapter 4** for pretreatment recommendations.

† If very high nuclear background is observed, you can raise temperatures to 40 or 41°C. However, some signal sensitivity may be lost. Please contact support or your FAS before making these changes.

‡ VS Universal AP Amp 5 incubation time is determined by instrument calibration. Use the instrument setting previously optimized for the mRNA Universal software. For assistance, consult your local ACD FAS for more information.

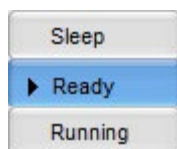
7. Select your preferred counterstain and post-counterstain settings. For optimal visualization of multiplex chromogenic staining, a light counterstain is recommended.
8. At the top of the Protocol Steps window, click **Save As**, then select a unique protocol number from the drop-down menu and choose a protocol name.
9. Click **Active**, specify any relevant comments in the available field, and then click **Save**.
10. Click **Close** to go back to the main screen.
11. Assign a probe number from the list to each probe of interest. For each probe selected, assign a protocol.

Print the labels

1. Select the **Print Label** icon from the upper right corner of the home screen.
2. Select your preferred template or create a new template. To create a new template, refer to the *Ventana DISCOVERY ULTRA System User Manual* for details.
3. Click **Protocol**.
4. Select the protocols you created in the section above. Click the **Add** button.
5. When the protocols for all slides have been assigned, click **Close/Print**.
6. Fill in the template for each slide. Click **Print** when completed.

Load the reagents

1. Remove the nozzle caps from the filled dispensers and place each cap on the post located on the back of the dispenser.
2. Prime the user-fillable dispensers. For guidance, refer to the instructions provided by Roche Tissue Diagnostics.
3. If needed, remove any air bubbles at the nozzle tip by pushing down on the nozzle until the liquid reaches the tip of the nozzle or forms a small meniscus at the tip of the nozzle.
4. Remove the yellow locking ring from the dispensers in all the prefilled dispensers. Refer to the instructions provided by Roche Tissue Diagnostics.
5. Load the dispensers onto the reagent racks.
6. Load the reagent racks onto the reagent carousel.
7. Select the **Ready** button.

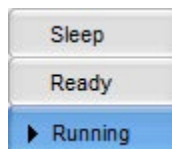


8. Open the slide drawers.

- Load each slide onto a heater pad with the label facing upward and inward. Ensure that the slides sit securely on the pads.

IMPORTANT! Prior to loading the slides, ensure heater pads are completely dry. Wipe off any liquid using laboratory tissue paper.

- Close the slide drawers.
- Select the **Running** button.



- The assay duration varies based on assay selections, approximately **10–13 HRS**.

IMPORTANT! Before leaving the instrument unattended, ensure all reagents and slides are successfully registered and the instrument is running.

Complete the run

- After the run is complete, remove the Dewax reagent and Target Retrieval v2, place nozzle caps on the dispensers, and store at room temperature.
- For the remaining reagents, place nozzle caps back on the dispensers and place racks onto magnet locking tray and Store at **4°C**.

Wash and dry the slides

- Prepare 200 mL of diluted detergent by adding 1–2 drops detergent to 200 mL distilled water in a container with a cap.
- Mix well by inverting the container 4–5 times.
- Add diluted detergent to a Tissue-Tek Staining Dish.

Note: Store diluted detergent at **RT**.

- Submerge a Tissue-Tek Slide Rack into the Tissue-Tek Staining Dish containing 200 mL diluted detergent.
- Open the instrument slide drawers and unload slides.
- Decant solution on the slides into the slide drawer, then *immediately* load slides into the Tissue-Tek Slide Rack submerged in detergent.
- Rinse oil off the slides by moving the slide rack up and down in the dish 10 times.
- Replace the detergent with distilled water and rinse slides by moving the slide rack up and down a minimum of **10** times.
- Repeat Step 8 three to five times.
- Transfer the slides into a Tissue-Tek Staining Dish containing **200 mL** distilled water.
- Place slides in a drying oven at **60°C** for at least **30 MIN**.

Mount the samples

- In a fume hood, fill two clearing agent dishes with ~200 mL fresh xylene.

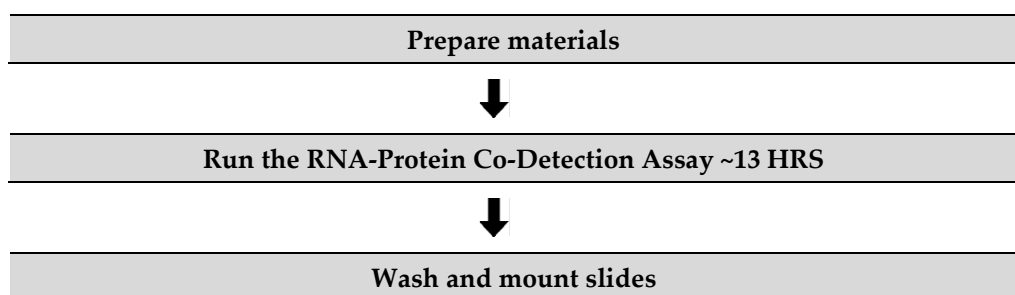
2. Once slides are dry, move the Tissue-Tek Slide rack into the staining dish containing xylene for **1 MIN** with occasional agitation.
3. Move the Tissue-Tek Slide rack into the staining dish containing xylene for **1 MIN** with occasional agitation.
4. Lay each slide flat with the sections facing up in the fume hood then add 1–2 drops of EcoMount or other chromogen-compatible xylene-based mounting medium. Carefully place a 24 mm x 50 mm coverslip over the section and avoid trapping air bubbles.
5. Air dry slides for at least **15 MIN** before evaluation.
6. Proceed to **Chapter 7. Evaluate the Results.**

6

Chapter 6: Protease-Free, Sequential RNA-Protein Co-Detection

In the protease-free workflow for sequential RNA-Protein Co-Detection, tissue samples undergo pretreatment using Roche IHC-style cell conditioning with CC1, followed by further digestion using ACD's VS PretreatPro. This is followed by ISH detection, and then IHC detection and counterstaining. We recommend first establishing successful detection of RNA and IHC separately in your tissue of interest before proceeding to combined detection.

Workflow



Prepare materials

Materials required

Materials Provided by Advanced Cell Diagnostics	Materials Provided by Roche Tissue Diagnostics	Other Materials and Equipment
<ul style="list-style-type: none"> • RNAscope 2.5 VS Target Probe • RNAscope 2.5 VS Positive Control Probe • RNAscope 2.5 VS Negative Control Probe • RNAscope VS Dewax • VS PretreatPro* • RNAscope VS Target Retrieval† • RNAscope VS Universal AP AMP 1 • RNAscope VS Universal AP AMP 2 • RNAscope VS Universal AP AMP 3 • RNAscope VS Universal AP AMP 4 • RNAscope VS Universal AP AMP 5 • RNAscope VS Universal AP AMP 6 • RNAscope VS Universal AP AMP 7 	<ul style="list-style-type: none"> • DISCOVERY ULTRA — automated slide stainer • DISCOVERY Wash Buffer 10X • ULTRA LCS (Predilute) • SSC Buffer 10X • DISCOVERY CC1 • Reaction Buffer 10X • Probe dispensers • mRNA RED Probe Amplification Kit • mRNA Universal Sample Prep Kit • mRNA RED Detection Kit • User fillable dispensers • IHC detection options (see Chapter 1) 	<ul style="list-style-type: none"> • Distilled water • Dawn detergent or similar detergent • Fume hood • Xylene • Tissue-Tek Staining Dish • Tissue-Tek Clearing Agent Dish, xylene-resistant • Tissue-Tek Vertical 24 Slide Rack • EcoMount • Cover Glass, 24 mm x 50 mm • Primary antibody • ProLong Gold Antifade Reagent

Materials Provided by Advanced Cell Diagnostics	Materials Provided by Roche Tissue Diagnostics	Other Materials and Equipment
<ul style="list-style-type: none"> • RNAscope VS Hematoxylin (optional) • RNAscope VS Bluing Reagent (optional) 		

* This dispenser is only used for the protease-free workflow

† These dispensers are used for the protease workflow

Prepare the instrument

If the instrument has not been used for > 1 week, follow guidelines for instrument maintenance from Roche Tissue Diagnostics. Before use, empty the waste carboys if needed.

Dilute instrument bulk reagents

1. Prepare the instrument bulk fluids according to the manufacturer's instructions.
2. Fill bulk solution containers for 1X DISCOVERY Wash, ULTRA LCS (predilute), and CC1 (predilute) to be at least half full. Completely fill bulk solution containers for 2X SSC and 1X Reaction Buffer.

IMPORTANT! Do not use expired reagents.

Register new reagents

Reagent dispensers come with appropriate barcode labels and registration buttons for dispensing RNAscope VS Universal Reagents. Refer to the *Ventana DISCOVERY ULTRA System User Manual* for details. To register reagents:

- Log all ACD reagents and probes into the software as **log user-fillable reagents** and **log user-fillable probes**, respectively.
- Use the reagent registration wand that comes with the instrument to register new reagent kits from Roche Tissue Diagnostics

Prepare user-fillable reagents for RNA-Protein Co-Detection

Refer to **Chapter 1** to determine the proper dispenser for each reagent.

IMPORTANT! Avoid cross contamination between reagents. Dewax must be warmed to room temperature and be completely in solution before use.

1. Transfer the entire volume of each AMP component of the Detection Kit to the corresponding labeled dispenser from the appropriate mRNA Amplification kit.
2. Fill the mRNA Sample Prep Kit by transferring the VS Dewax reagent from the RNAscope VS Sample Prep Reagent Kit v2 to the mRNA Dewax Dispenser.

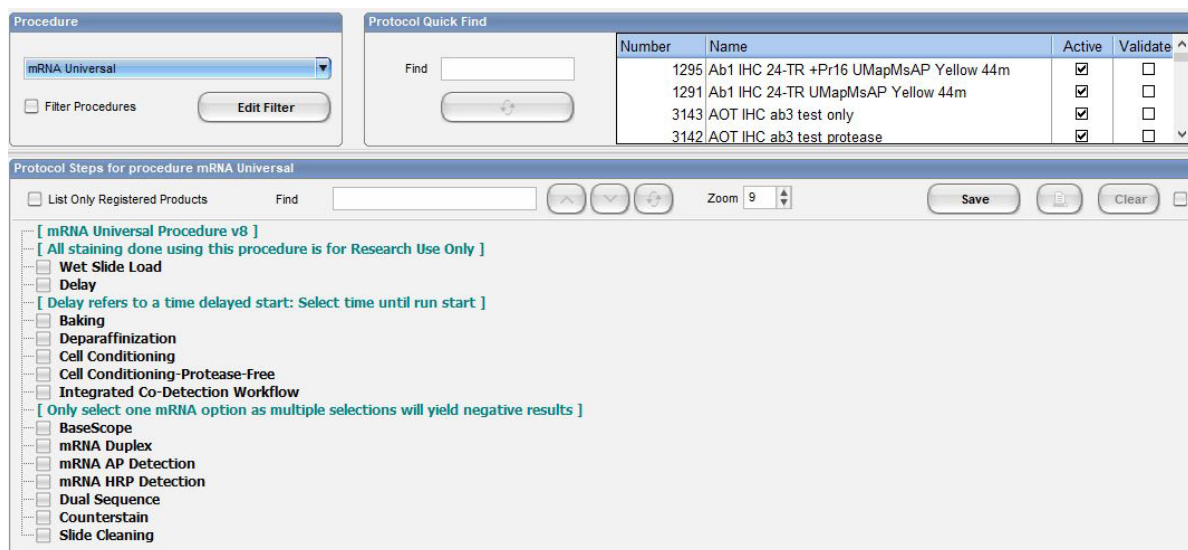
Note: mRNA Target Retrieval and mRNA Protease dispensers are not needed for the protease-free workflow.

3. Fill the following dispensers:

- a. Transfer the RNAscope 2.5 VS Target Probe and control probes to the corresponding probe dispensers.
- b. Transfer the VS PretreatPro and VS CoDetectPro from the RNAscope VS Pro Reagents to the open pretreatment dispensers. Ensure the barcode number aligns with the protocol used.
- c. Transfer the VS Hematoxylin and VS Bluing to the Counterstain 1 and Counterstain 2 dispensers.
4. Follow the dispenser product insert instructions to properly prime and handle the dispensers.
5. Store tightly capped dispensers at 4°C when not in use.
6. Store mRNA Dewax and Target Retrieval v2 dispenser at room temperature when not in use.

Create an instrument protocol

1. Open the NexES software and click the **Protocol** button.
2. Click **Create/Edit Protocols**, then select **mRNA Universal** from the Procedure drop-down menu.
3. Confirm that “mRNA Universal v8” is displayed. The main protocol selections appear as shown:



4. Select the appropriate pretreatment conditions according to the following screenshot. After selecting the main steps, drop-down menus become available. We recommend selecting **Cell-Conditioning-Protease-Free** with CC1 for the time and temperature specified in **Chapter 4** according to tissue-specific recommendations found in **Appendix B**.

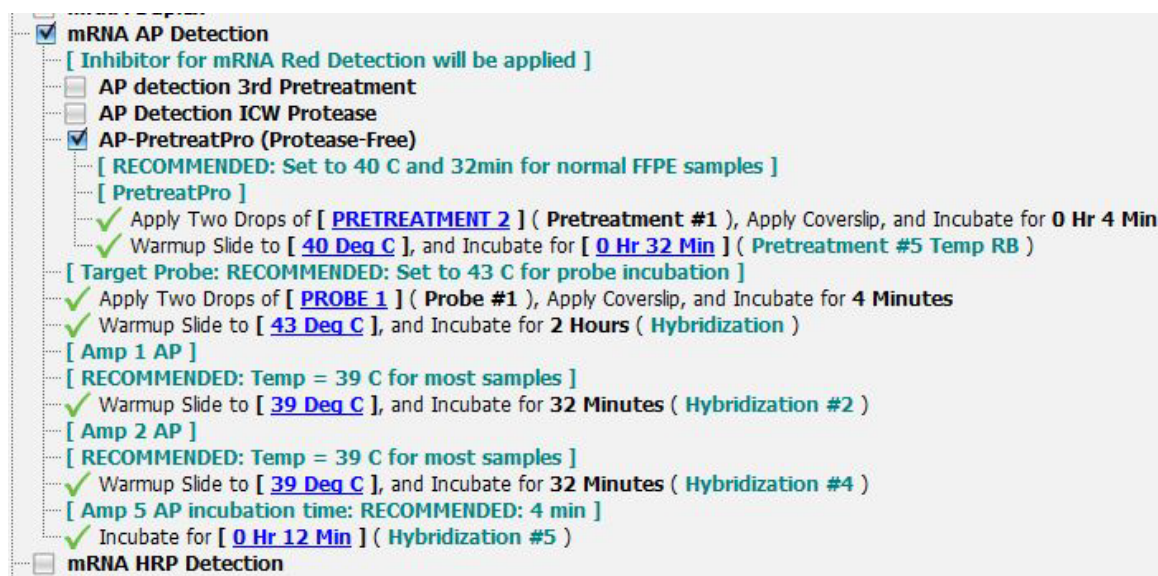


Note: The duration of tissue cell conditioning corresponds to the longest selected time. It is not the sum of all selected times.

- To enable ISH staining with RNAscope VS Universal AP, select **mRNA AP Detection**.

Note: Select only one mRNA option per protocol. Choosing multiple mRNA options gives negative results.

- Select **AP-PretreatPro** to apply and incubate the VS PretreatPro reagent. We recommend incubating this reagent for **32 MIN** at **40°C**.
- Select the remaining ISH staining conditions according to the following screenshot and table:



Standard Times/Temperatures for mRNA AP Detection	
Cell-Conditioning–Protease-Free (CC1) *	64 MIN @95°C
VS Protease temperature and time*	37°C, 16 MIN
AP-PretreatPro (protease-free) temperature and time*	40°C, 32 MIN

Standard Times/Temperatures for mRNA AP Detection	
Suggested probe temperature	Single Probes - 43°C HPV Pooled Probes - 50°C
Suggested RNAscope AP AMP 1 and AMP 2 temperature	39°C†
RNAscope AP AMP 5 incubation time	Instrument titrated ‡

* Refer to **Chapter 4** for the CC1 time and temperature recommended for your tissue type.

† If very high nuclear background is observed, you can raise temperatures to 40 or 41°C. However, some signal sensitivity may be lost. Please contact support or your FAS before making these changes.

‡ VS Universal AP Amp 5 incubation time is determined by instrument calibration. Use the instrument setting previously optimized for the mRNA Universal software. For assistance, consult your local ACD FAS for more information.

8. Select **Dual Sequence** to enable IHC following ISH detection.
9. Apply VS CoDetectPro from the RNAscope VS Pro Reagents by selecting **DS Pretreatment**, **DS 2nd Pretreatment** and **Use DW for DS 2nd Pretreatment**. Select a pretreatment barcode to correspond to the VS CoDetectPro reagent. We recommend incubating at **40°C** for **32 MIN** prior to protein detection to quench RNAscope enzymes prior to IHC detection.

IMPORTANT! Make sure that VS CoDetectPro uses a unique Pretreatment number, different from VS PretreatPro, to ensure the protocol runs correctly.

☒ **Dual Sequence**

- ☐ **Antibody Denaturation**
- ☒ **DS Pretreatment**
 - ☐ **DS 1st Pretreatment**
 - ☐ **Protease-ISH**
 - ☐ **DS Cell Conditioning**
 - ☒ **DS 2nd Pretreatment**
 - ☐ **Use RB for DS 2nd Pretreatment**
 - ☒ **Use DW for DS 2nd Pretreatment**
 - ☐ **DS Disable heat PT2-EZ**
- [RECOMMENDED: If using temperatures above 42°C, set incubation for less than 1 hour]
- ✓ Warmup Slide to [40 Deg C], and Incubate for 4 Minutes (DS Pretreatment #2 Temp DW)
- ✓ Apply One Drop of [PRETREATMENT 1] (DS Pretreatment #2), and Incubate for [0 Hr 32 Min]
- ☐ **DS Enzyme**
 - ☐ **DS Option**
 - ☐ **DS A/B Block**
 - ☐ **DS Inhibitor**

10. To apply primary antibody, select **DS Antibody** and your desired incubation time and temperature, using previously optimized IHC-only conditions. See the following for an example.

☒ **Dual Sequence**
☐ Antibody Denaturation
☐ DS Pretreatment
☐ DS Option
☐ DS A/B Block
☐ DS Inhibitor
 [DISC Inhibitor will not be applied]
☐ DS ISH
☒ **DS Antibody**
☐ DS Antibody Manual Application
☐ DS Antibody Blocking
☐ Disable heat for DS Antibody
☐ DS High Temp Ab incubation
☒ Warmup Slide to [37 Deg C] from Very Low Temperatures (DS Primary Antibody)
☐ DS Extended Ab incubation
☒ Apply One Drop of [ANTIBODY 1] (DS Antibody), and Incubate for [32 Minutes]
☐ Disable heat for 2nd Fixative

11. Select the settings for secondary detection:

Note: To prevent non-specific cross-detection of RNAscope, we recommend using Roche's Antibody Block reagent immediately before applying the secondary antibody. Skipping the Antibody Block step could cause a hue shift in the RNA dots, making it difficult to accurately interpret RNA-protein colocalization.

- Do the following for **HRP-based secondary detection**:
 - a. Choose a Roche Multimer HRP reagent corresponding to the primary host species.
 - b. To enable secondary incubation, select **DS Multimer HRP**.
 - c. To apply Antibody Block, select **DS Multimer HRP Blocker**, then select Antibody Block reagent.
 - d. Select your Multimer HRP reagent of choice and desired incubation time based on previously optimized IHC-only conditions.

☐ Disable heat for 2nd Detection
☒ **DS Multimer HRP**
☒ **DS Multimer HRP Blocker**
 [Select Multimer blocker and Multimer species]
 [Note: Recommended Multimer HRP Reagent incubation time for ICW is 32min]
☒ Apply One Drop of [Antibody Block] (DS Multimer HRP Blocking), No Coverslip and Incubate for 32 Minutes
☒ Apply One Drop of [Umap anti-Ms HRP] (DS Multimer HRP), and Incubate for [32 Minutes]
☐ DS Multimer AP
☐ DS Proximity Detection
☐ DS Enzyme conjugate
☐ DS DISCOVERY Amplification
☐ DS DAB

- Do the following for **AP-based secondary detection**:
 - a. Choose a Roche Multimer AP reagent corresponding to the primary host species.
 - b. To enable secondary incubation, select **DS Multimer AP**.
 - c. To apply Antibody Block, select **DS Multimer AP Blocker**, then select Antibody Block reagent.

- d. Select your Multimer AP reagent of choice and desired incubation time based on previously optimized IHC-only conditions.

☐ Disable heat for 2nd Detection
☐ DS Multimer HRP
☒ DS Multimer AP

☒ DS Multimer AP Blocker

[Select Multimer blocker and Multimer species]
[Note: Recommended Multimer AP Reagent incubation time for ICW is 32min]
☒ Apply One Drop of [Antibody Block] (DS Multimer AP Blocking #1), No Coverslip and Incubate for 32 Minutes
☒ Apply One Drop of [UMap anti-Ms AP] (DS Multimer AP), and Incubate for [32 Minutes]
[Select Multimer blocker and Multimer species]

☐ DS Proximity Detection
☐ DS Enzyme conjugate
☐ DS DISCOVERY Amplification
☐ DS DAB

8. Select chromogens for IHC detection using the following recommendations:

☐ DS Enzyme conjugate
☐ DS DISCOVERY Amplification
☒ DS DAB
☐ DS Silver
☒ DS DISCOVERY Purple
☐ DS DISCOVERY Yellow HRP
☒ DS DISCOVERY Teal HRP
☒ DS DISCOVERY Green HRP
☐ DS DISCOVERY Red HRP
☐ DS DISCOVERY Blue HRP
☒ DS Red
☒ DS DISCOVERY Red
☒ DS DISCOVERY Yellow
☐ DS Blue
☒ DS Rhodamine
☒ DS FITC
☒ DS Cy5
☒ DS DCC
☒ DS FAM
☒ DS Red 610
☒ DS Rhodamine 6G
☒ DS Open Detection Kit

HRP Chromogens

AP Chromogens

DISCOVERY HRP Fluorescent Kits

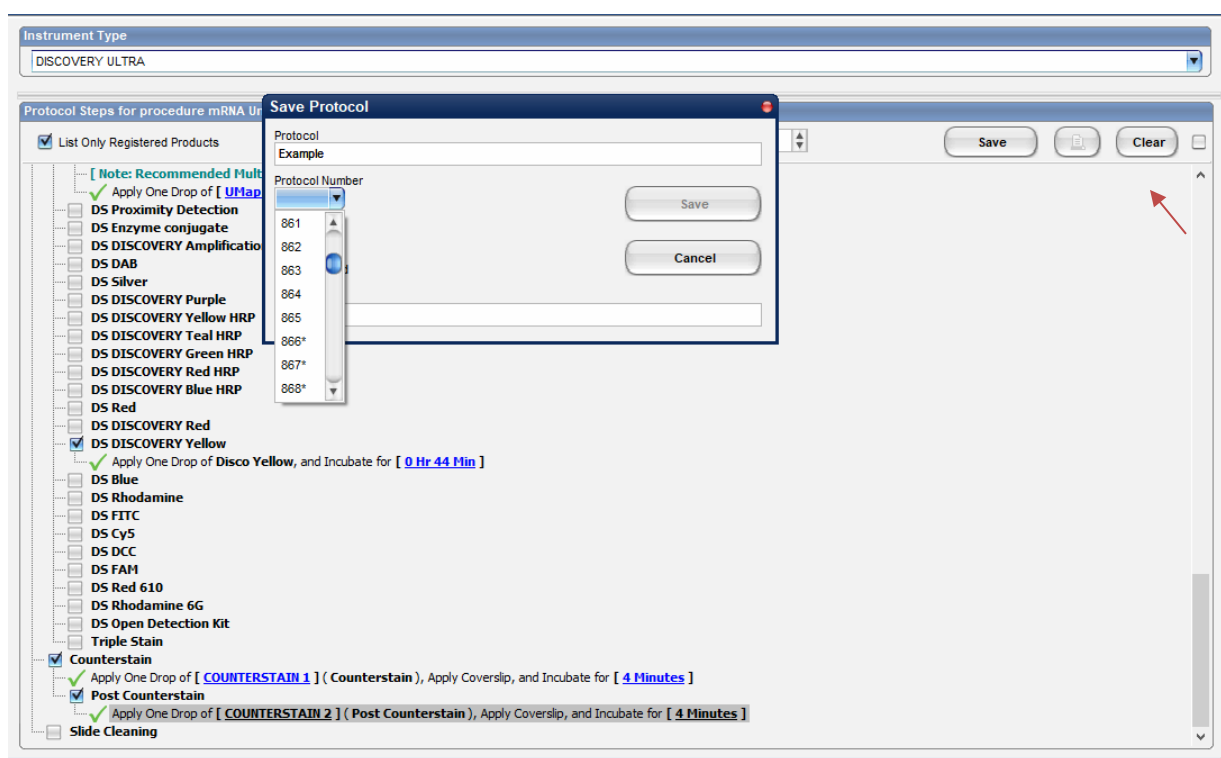
IHC Chromogen Selection and Settings		
IHC Enzyme	IHC Chromogen	Recommended Chromogen Settings*
AP	DISCOVERY Yellow	44 MIN – 2 HRS
	DISCOVERY Red	12 MIN†
	DS Red / ChromoMap Red‡	Default
HRP (chromogenic)	DS DAB / ChromoMap DAB	Default
	DISCOVERY Purple	40 MIN
	DISCOVERY Teal HRP	DISCOVERY Teal H ₂ O ₂ – 16–32 MIN DISCOVERY Teal Act – 16 MIN
	DISCOVERY Green HRP	DISCOVERY Green H ₂ O ₂ – 16–32 MIN DISCOVERY Green Act – 16 MIN
HRP (fluorescent)	DISCOVERY DCC Kit	32 MIN
	DISCOVERY FAM Kit	20 MIN
	DISCOVERY FITC Kit	20 MIN
	DISCOVERY Rhodamine Kit	32 MIN
	DISCOVERY Rhodamine 6G Kit	32 MIN
	DISCOVERY Red 610 Kit	32 MIN
	DISCOVERY Cy5 Kit	40 MIN

*We recommend using the same IHC chromogen settings for sequential ISH-IHC as optimized for IHC alone.

†Extending DISCOVERY Red incubation is not recommended, as it can result in a dot-like background.

‡For stronger AP-based Red IHC detection, ChromoMap Red is recommended (select **DS Red**).

- Select your preferred counterstain and post-counterstain settings. A light counterstain is recommended for best visualization of multiplex chromogenic staining.
- At the top of the Protocol Steps window, click **Save As**, then select a unique protocol number from the drop-down menu and choose a protocol name.
- Click **Active**, specify any relevant comments in the available field, and then click **Save**.



12. Create a new protocol for each probe/antibody/chromogen combination. Click **Save**.

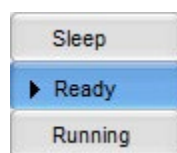
Print the labels

1. Select the **Print Label** icon from the upper right corner of the home screen.
2. Select your preferred template or create a new template. To create a new template, refer to the *Ventana DISCOVERY ULTRA System User Manual* for details.
3. Click **Protocol**.
4. Select the protocols you created in the section above. Click the **Add** button.
5. When the protocols for all slides have been assigned, click **Close/Print**.
6. Fill in the template for each slide. Click **Print** when complete.

Load the reagents

Remove the nozzle caps from the filled dispensers and place each cap on the post located on the back of the dispenser.

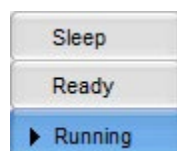
1. Prime the user-fillable dispensers. For guidance, refer to the instructions provided by Roche Tissue Diagnostics.
2. If needed, remove any air bubbles at the nozzle tip by pushing down on the nozzle until the liquid reaches the tip of the nozzle or forms a small meniscus at the tip of the nozzle.
3. Remove the yellow locking ring from the dispensers in all the prefilled dispensers. Refer to the instructions provided by Roche Tissue Diagnostics.
4. Load the dispensers onto the reagent racks.
5. Load the reagent racks onto the reagent carousel.
6. Select the **Ready** button.



7. Open the slide drawers.
8. Load each slide onto a heater pad with the label facing upward and inward. Ensure that the slides sit securely on the pads.

IMPORTANT! Prior to loading the slides, ensure heater pads are completely dry. Wipe off any liquid using laboratory tissue paper.

9. Close the slide drawers.
10. Select the **Running** button.



11. The assay duration varies based on assay selections, approximately **15 – 20 HRS**.

IMPORTANT! Before leaving the instrument unattended, ensure all reagents and slides are successfully registered and the instrument is running.

Complete the run

1. After the run is complete, remove the Dewax reagent and Target Retrieval v2, place nozzle caps on the dispensers, and store at room temperature.
2. For the remaining reagents, place nozzle caps back on the dispensers and place racks onto magnet locking tray and Store at **4°C**.

Wash and dry the slides

1. Prepare 200 mL of diluted detergent by adding 1–2 drops detergent to 200 mL distilled water in a container with a cap.
2. Mix well by inverting the container 4–5 times.
3. Add diluted detergent to a Tissue-Tek Staining Dish.

Note: Store diluted detergent at **RT**.

4. Submerge a Tissue-Tek Slide Rack into the Tissue-Tek Staining Dish containing 200 mL diluted detergent.
5. Open the instrument slide drawers and unload slides.
6. Decant solution on the slides into the slide drawer, then *immediately* load slides into the Tissue-Tek Slide Rack submerged in detergent.
7. Rinse oil off the slides by moving the slide rack up and down in the dish 10 times.
8. Replace the detergent with distilled water and rinse slides by moving the slide rack up and down a minimum of **10** times.
9. Repeat Step 8 three to five times.
10. Transfer the slides into a Tissue-Tek Staining Dish containing **200 mL** distilled water.
11. Place slides in a drying oven at **60°C** for at least **30 MIN**.

Mount the samples

1. In a fume hood, fill two clearing agent dishes with ~200 mL fresh xylene.
2. Once slides are dry, move the Tissue-Tek Slide rack into the staining dish containing xylene for **1 MIN** with occasional agitation.
3. Move the Tissue-Tek Slide rack into the staining dish containing xylene for **1 MIN** with occasional agitation.
4. Lay each slide flat with the sections facing up in the fume hood then add 1–2 drops of EcoMount or other chromogen-compatible xylene-based mounting medium. Carefully place a 24 mm x 50 mm coverslip over the section and avoid trapping air bubbles.
5. Air dry slides for at least **15 MIN** before evaluation.

IMPORTANT! mRNA Teal, mRNA Green, DISCOVERY Teal HRP and DISCOVERY Green HRP chromogens are light sensitive and can fade over time. For best results, protect stored slides from the light and image within one week of staining.

7

Chapter 7. Evaluate the Results

Evaluate the samples

Examine the experimental results.

- Assess tissue and cell morphology.
Assess the negative control background first.
- Assess positive control signal strength. Positive control signal should be visible as punctate dots within the cell. Highly expressed targets might appear as clusters.
- Determine which pretreatment condition yielded the highest positive control signal and lowest negative control signal.
- If none of the conditions are satisfactory, contact technical support at **support.acd@bio-technne.com**.

Staining Score	Scoring Definition
0	No staining or <1 dot/10 cells
1	1 – 3 dots/cell
2	4 – 9 dots/ cell or very few dot clusters
3	10 – 15 dots/ cell and <10% dots are in clusters
4	>15 dots/ cell and >10% dots are in clusters

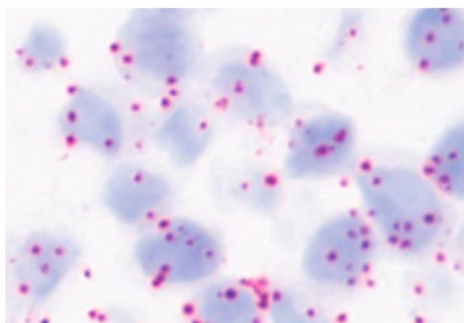
Troubleshooting

For troubleshooting information, please contact technical support at **support@acdbio.com**.

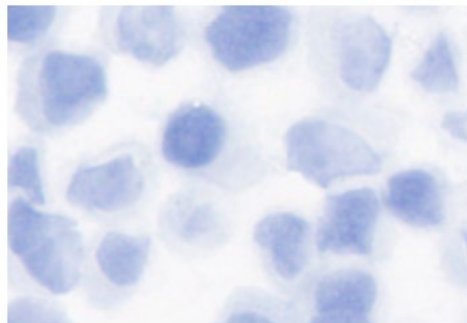
Tissue example

If the assay is successful, the staining should look like the following images:

Figure 2. RNAscope VS Universal AP Assay results in HeLa cells



Hs-TBP (Positive Control)



DapB (Negative Control)



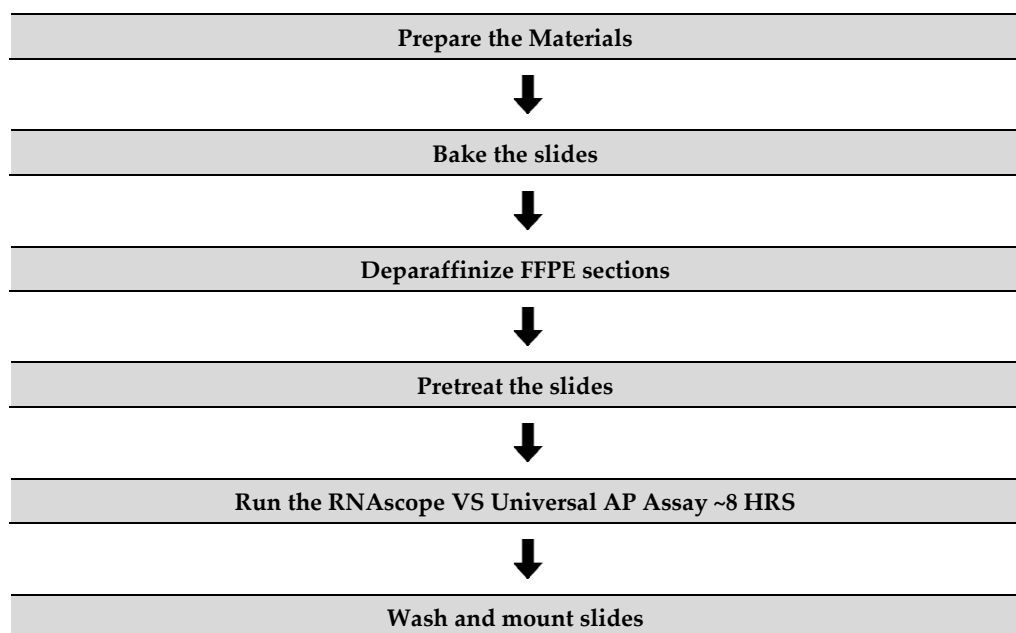
Appendix A. Semi-automated RNAscope VS Universal AP Assay

Most sample types can be fully automated on the Discovery ULTRA. However, manual pretreatment may yield better results in some cases. For tissues that do not achieve satisfactory results with the fully automated procedure and conditions in **Chapter 4**, use the semi-automated procedure.

The following protocols describe formalin-fixed, paraffin-embedded (FFPE) sample pretreatment. For other sample types and preparation methods, contact support.acd@bio-techne.com for the latest protocols and guidelines.

IMPORTANT! We highly recommend following these guidelines. We cannot guarantee assay results with other preparation methods.

Workflow



Kit contents and storage

For Offline Boiling: RNAscope Target Retrieval Reagents				
<input checked="" type="checkbox"/>	Cat. No.	Reagent	Quantity	Storage
	322000	RNAscope Target Retrieval Reagents	70 mL x 4 bottles	Room Temp (15–30°C)

IMPORTANT! Do not substitute the reagent components of the RNAscope VS Universal Reagent Kit with those of other RNAscope Reagent Kits, even those having the same name. The Target Retrieval v2 solution in the RNAscope VS Universal AP Kit CANNOT be used for offline boiling. Please use the RNAscope Target Retrieval Reagents (Cat. No. 322000) to boil samples off the instrument.

Prepare the materials

Materials can be prepared ahead of time or while baking the slides, unless otherwise stated.

Materials required

Materials provided by Advanced Cell Diagnostics	Materials Provided by Roche Tissue Diagnostics	Other Materials and Equipment
<ul style="list-style-type: none"> • RNAscope 2.5 VS Target Probe • RNAscope 2.5 VS Positive Control Probe • RNAscope 2.5 VS Negative Control Probe • RNAscope Target Retrieval Reagents (Offline) • RNAscope VS Protease • RNAscope VS Universal AP AMP 1 • RNAscope VS Universal AP AMP 2 • RNAscope VS Universal AP AMP 3 • RNAscope VS Universal AP AMP 4 • RNAscope VS Universal AP AMP 5 • RNAscope VS Universal AP AMP 6 • RNAscope VS Universal AP AMP 7 • RNAscope VS Hematoxylin • RNAscope VS Bluing Reagent 	<ul style="list-style-type: none"> • DISCOVERY ULTRA — automated slide stainer • DISCOVERY Wash 10x • ULTRA LCS (Predilute) • SSC Buffer 10X • Reaction Buffer 10X • Probe dispensers • mRNA RED Probe Amplification Kit • mRNA RED Detection Kit • User fillable dispensers • mRNA Universal Sample Prep Kit • CCI Buffer 	<ul style="list-style-type: none"> • Distilled water • Glass beaker (1 or 2 L) • Drying oven • Hot plate • Dawn detergent or similar detergent • Fume hood • Xylene • 100% ethanol (EtOH) • Tissue-Tek Staining Dishes • Tissue-Tek Clearing Agent Dishes, xylene-resistant • Tissue-Tek Vertical 24 Slide Rack • EcoMount • Cover Glass, 24 mm x 50 mm

Prepare the instrument

If the instrument has not been used for ≥ 1 week, follow the guidelines for instrument maintenance in the *Ventana DISCOVERY ULTRA System User Manual*.

Dilute bulk reagents

Please prepare the bulk fluids per manufacturer instructions.

Register new reagents

Reagent dispensers come with appropriate barcode labels and registration buttons for dispensing RNAscope VS Reagents. Refer to the *Ventana DISCOVERY ULTRA System User Manual* for details.

To register reagents:

1. Log all ACD reagents and probes into the software as “log user-fillable reagents” or “log user-fillable probes”.
2. Use the wand that comes with the instrument to register *new* reagent kits.

Prepare instrument reagents

Refer to **Chapter 1** to determine the proper dispenser for each reagent.

1. For RNAscope VS Universal AP AMP 1 to AMP 7, transfer the entire volume of each component into the correspondingly labeled dispenser.
2. Transfer the VS Protease, VS Hematoxylin, and VS Bluing Reagent to the correspondingly labeled dispenser.
3. Use Roche Tissue Diagnostics probe dispensers for 2.5 VS Target Probe, 2.5 VS Positive Control Probe, 2.5 VS Negative Control Probe. Follow the dispenser product insert instructions to properly prime and handle the dispensers.
4. Store capped dispensers at 4°C when not in use.
5. Check solution levels: DISCOVERY Wash, 2X SSC, Reaction Buffer, ULTRA LCS (Predilute), and CC1. Refill if they are less than half full. To run a full tray of 30 slides, completely fill the 2X SSC and Reaction Buffer containers.

IMPORTANT! Do not use expired reagents.

6. Empty the waste carboy if needed.

Prepare deparaffinization reagents

- In a fume hood, fill two clearing agent dishes with ~200 mL fresh xylene.
- In a fume hood, fill two staining dishes with ~200 mL fresh 100% EtOH.

Note: Ensure all containers remain covered when not in use.

Prepare 1X Target Retrieval

1X Target Retrieval is used in manual cell conditioning (CC).

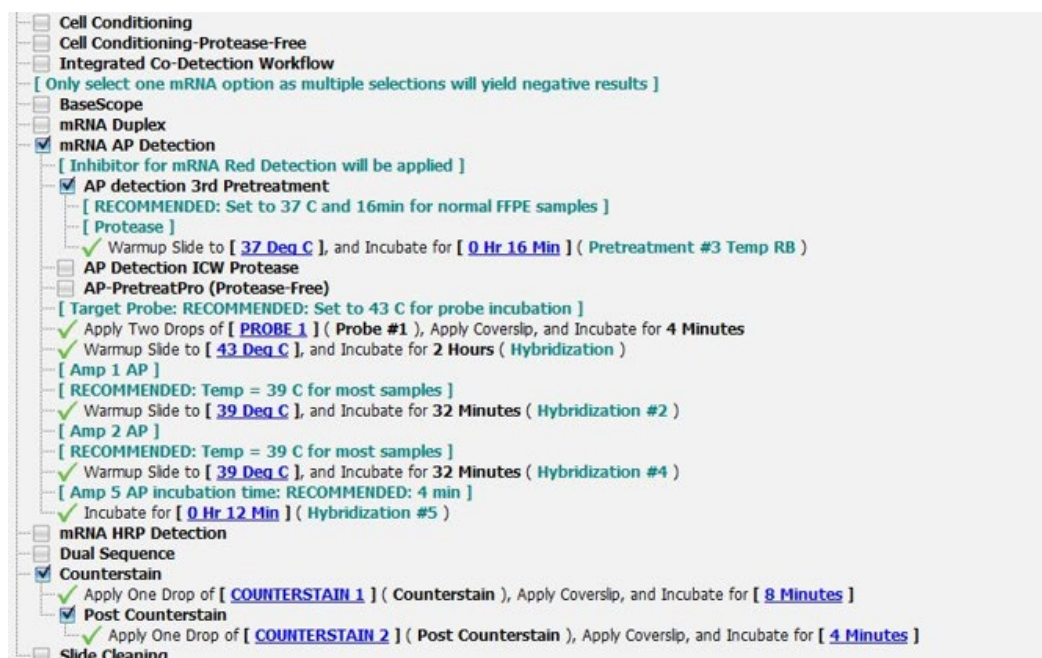
1. Prepare 700 mL of fresh 1X Target Retrieval **by** adding 630 mL distilled water to 1 bottle (70 mL) 10X Target Retrieval solution in the beaker.
2. Mix well and cover the beaker with foil.

IMPORTANT! Do not use RNAscope VS Target Retrieval v2 for offline boiling.

Create an instrument protocol

1. Open the VS software and click on the **Protocol** button.
2. Click on **Create/Edit Protocols**, go to the Procedure drop down menu and select **mRNA Universal**.

Main protocol steps appear as shown:



IMPORTANT! Do not select Baking, Deparaffinization, or Cell Conditioning.

3. After selecting the main protocol steps, drop down menus become available. Select the appropriate protocol steps by clicking on the associated check boxes as shown above.
4. Select the appropriate assay conditions from the drop-down menus according to the following table:

Standard Temperatures/Times	
VS Protease*	37°C
Standard probe temperatures	Single Probes 43°C
	HPV Pooled Probes - 50°C
Standard AMP 1 and AMP 2 temperatures*	39°C†
RNAScope AP AMP 5 incubation time	Instrument titrated‡

* Refer to **Chapter 4** for the CC1 time and temperature recommended for your tissue type

† If very high nuclear background is observed, you can raise temperatures to 40 or 41°C. However, some signal sensitivity may be lost. Please contact support or your FAS before making these changes.

‡ VS Universal AP Amp 5 incubation time is determined by instrument calibration. Use the instrument setting previously optimized for the mRNA Universal software. For assistance, consult your local ACD FAS for more information.

5. Click **Save As**, then select a protocol number from the drop-down menu and choose a protocol name for each probe.
6. Click **Save**.
7. Click **Close** to go back to the main screen.
8. Assign a probe number from the list to each probe of interest. For each probe selected, assign a protocol.

Print the labels

1. Select the **Print Label** icon from the bottom of the home page screen.
2. Select your preferred template or create a new template. To create a new template, refer to the *Ventana DISCOVERY ULTRA System User Manual* for details.
3. Select the protocol you created for the RNAscope VS Universal Assay.
4. Click on **Protocol** to add and print the label.

Manually pretreat the samples

Bake the slides

1. Bake slides in a dry oven for **30-60 MIN** at **60°C**.

OPTIONAL STOPPING POINT Use immediately or store at **RT** with desiccants for ≤1 week. Prolonged storage may degrade sample RNA.

Deparaffinize FFPE sections

Note: If you have not done so already, create a protocol for your instrument and print slide labels during this procedure.

1. Place slides in a Tissue-Tek Slide Rack and submerge in the first xylene-containing clearing agent dish in the fume hood.
2. Incubate the slides in xylene for **5 MIN** at **RT**. Agitate the slides by occasionally lifting the slide rack up and down in the clearing agent dish.
3. Remove the slide rack from the first xylene-containing dish and *immediately* place in the second xylene-containing clearing agent dish in the fume hood.
4. Repeat Step 2.
5. Remove the slide rack from the second xylene-containing dish and immediately place in the staining dish containing 100% EtOH.
6. Incubate the slides in 100% EtOH for 1 MIN at RT with agitation.
7. Repeat Step 6 with fresh 100% EtOH.
8. Remove the slides from the rack, and place on absorbent paper with the section face-up. Air dry for **5 MIN** at **RT**.
9. While slides are drying, place printed labels on the slides.

IMPORTANT! Labels must be in place prior to the next section.

10. Insert the slides into a Tissue-Tek Slide Rack and proceed to condition the slides.

Pretreat the slides

Begin heating 1X Target Retrieval Buffer while FFPE slides are baking at 60°C or during the previous section.

IMPORTANT! Do not boil 1X Target Retrieval Buffer for more than 30 MIN before use.

1. Heat 1X Target Retrieval Buffer to 98–104°C:
 - a. Place the beaker containing 1X Target Retrieval Buffer on the hot plate. Cover the beaker with foil and turn the hot plate on high for 10–15 MIN.
 - b. Once 1X Target Retrieval Buffer reaches a slow boil (98–104°C), turn the hot plate to a lower setting to maintain the correct temperature. Check the temperature with a thermometer.
2. With a pair of forceps *very slowly* submerge the slide rack containing the slides into the boiling 1X Target Retrieval Buffer solution. Cover the beaker with foil and boil the slides for the time indicated below, according to target retrieval strength recommended in the following tables:

Extra Mild	Mild	Standard	Extended
7 MIN	10 MIN	15 MIN	30 MIN

Tissue-specific pretreatment recommendations:

Tissue Type	Offline Target Retrieval Time
Brain and spinal cord	Standard
Breast cancer	Standard
Cell lines	Mild
Colon	Standard
GI tract	Standard
Head and neck cancer	Standard
Heart	Standard
Kidney	Standard
Liver	Extended
Lung	Standard
Lymphoma	Mild
Placenta	Standard
Prostate	Standard
Skin	Standard
Stomach	Standard
Thymus	Mild
Tonsil	Mild
Xenograft derived from cell lines	Extra Mild
Xenograft derived from primary tumor	Standard
ACD control cell pellet	Mild

3. Immediately transfer the hot slide rack from the 1X Target Retrieval Buffer to a staining dish containing distilled water. Do not let the slides cool in Target Retrieval.
4. Wash slides 3–5 times by moving the Tissue-Tek Slide Rack up and down in the distilled water.
5. Repeat Step 4 with fresh distilled water.
6. Proceed directly to **Load the reagents** in the next section.

Run the RNAscope VS Universal AP Assay

Load the reagents

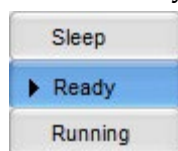
1. Remove the nozzle caps from the filled dispensers and place each cap on the post located on the back of the dispenser.
2. If needed, remove any air bubbles at the nozzle tips by pushing down on the nozzle until the liquid reaches the tip of the nozzle or forms a small meniscus at the tip of the nozzle.

IMPORTANT! Do not dispense any drops as this could compromise your drop inventory.

3. Load dispensers onto the reagent racks.
4. Remove the yellow locking ring from the dispensers in the prefilled **mRNA RED Detection Kit**. Refer to the instructions provided by Roche Diagnostics.
5. Load the reagent racks onto the reagent carousel.

Start the run

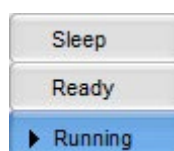
1. Select the **Ready** button.



2. Eject slide drawers.
3. Load each slide onto a heater pad with the label facing upward and inward. Ensure that the slides sit securely on the pads.

IMPORTANT! Prior to loading the slides, ensure the heater pads are completely dry. Wipe off any solution using laboratory tissue paper.

4. Close slide drawers.
5. Select the **Running** button. Automated assay will finish in **~8 HRS**.



IMPORTANT! Before leaving the instrument unattended, ensure all reagents and slides are successfully registered and the instrument is running.

Complete the run

After the run is complete, remove reagents, place nozzle caps back on the dispensers and place racks onto magnet locking tray.

IMPORTANT! Store reagent racks at 4°C until next use.

Wash the slides

1. Prepare 200 mL of diluted detergent by adding 1–2 drops detergent to 200 mL distilled water in a container with a cap.
2. Mix well by inverting the container 4–5 times.
3. Add diluted detergent to a Tissue-Tek Staining Dish.

Note: Store diluted detergent at RT.

4. Submerge a Tissue-Tek Slide Rack into the Tissue-Tek Staining Dish containing 200 mL diluted detergent.
5. Open the instrument slide drawers and unload slides.
6. Decant solution on the slides into the slide drawer, then *immediately* load slides into the Tissue-Tek Slide Rack submerged in detergent.
7. Rinse oil off the slides by moving the slide rack up and down in the dish 10 times.
8. Replace the detergent with distilled water and rinse slides by moving the slide rack up and down a minimum of 10 times.
9. Repeat Step 8 three to five times.

Dry and mount the samples

1. Remove the slide rack from the staining dish and dry slides in a 60°C dry oven for 30 MIN.

IMPORTANT! The RED substrate is alcohol sensitive. Do not dehydrate the slides in alcohol.

2. Cool the slides for 5 MIN at RT.
3. Briefly dip one slide into fresh pure xylene and immediately place 1–2 drops of EcoMount on the slide before the xylene dries.

IMPORTANT! Use the EcoMount mounting medium only.

4. Carefully place 24 mm x 50 mm coverslip over the tissue section. Avoid trapping air bubbles.
5. Repeat steps 2 and 3 for each slide.
6. Air dry slides for at least 5 MIN.
7. Proceed **Chapter 7. Evaluate the Results.**

B

Appendix B. Pretreatment Guidance for FFPE Samples

Follow the recommended pretreatment conditions based on your tissue type for:

- Any new or previously untested FFPE tissue types
- Samples prepared differently than the sample preparation protocol found in **Chapter 3**.
- For specific guidance on other sample preparations contact ACD Support at support.acd@bio-technne.com

Tissue-specific pretreatment conditions

Refer to the following table for tissue specific FFPE pretreatment conditions. For information about species or tissue type not listed here, contact support at support.acd@bio-technne.com.

Tissue Type	Protease-Free Workflow Recommended CC1 Target Retrieval (online)	Protease Workflow Recommended ACD Target Retrieval (online)
Cell Pellet	Mild	Mild
Brain	Standard	Mild
Spleen	Standard	Standard
Tonsil	Standard	Standard
Breast	Mild*	Standard
Prostate	Mild*	Standard
Intestine	Standard	Standard
Colon	Standard	Standard
Stomach	Standard	Standard
Kidney	Standard	Standard
Bladder	Standard	Standard
Heart	Standard	Standard
Lung	Standard	Standard
Skin	Standard	Standard
Head and Neck	Standard	Standard
Ovary	Standard	Standard
Cervical	Standard	Standard
Pancreas	Standard	Standard
Liver	Standard	Standard to Extended

*Mild pretreatment to reduce tissue detachment




Appendix C. Safety

Chemical safety

 **WARNING! GENERAL CHEMICAL HANDLING.** To minimize hazards, ensure laboratory personnel read and practice the general safety guidelines for chemical usage, storage, and waste provided below, and consult the relevant SDS for specific precautions and instructions:

- Read and understand the Safety Data Sheets (SDSs) before you store, handle, or work with any chemicals or hazardous materials. To obtain SDSs, see <http://www.acdbio.com/technical-support/user-manuals>
- Minimize contact with chemicals. Wear appropriate personal protective equipment when handling chemicals (for example, safety glasses, gloves, or protective clothing).
- Minimize the inhalation of chemicals. Do not leave chemical containers open. Use only with adequate ventilation (for example, fume hood).
- Characterize (by analysis if necessary) the waste generated by the particular applications, reagents, and substrates used in your laboratory.
- Ensure that the waste is stored, transferred, transported, and disposed of according to all local, state/provincial, and/or national regulations.
- **IMPORTANT!** Radioactive or biohazardous materials may require special handling, and disposal limitations may apply.

Biological hazard safety

 **WARNING! BIOHAZARD.** Biological samples such as tissues, body fluids, infectious agents, and blood of humans and other animals have the potential to transmit infectious diseases. Follow all applicable local, state/provincial, and/or national regulations. Wear appropriate protective equipment, which includes but is not limited to protective eyewear, face shield, clothing/lab coat, and gloves. All work should be conducted in properly equipped facilities using the appropriate safety equipment (for example, physical containment devices). Individuals should be trained according to applicable regulatory and company/institution requirements before working with potentially infectious materials. Read and follow the applicable guidelines and/or regulatory requirements in the following:

In the U.S.:

- U.S. Department of Health and Human Services guidelines published in Biosafety in Microbiological and Biomedical Laboratories found at www.cdc.gov/biosafety
 - Occupational Safety and Health Standards, Bloodborne Pathogens (29 CFR§1910.1030)
 - Your company's/institution's Biosafety Program protocols for working with/handling potentially infectious materials
 - Additional information about biohazard guidelines is available at www.cdc.gov/

In the EU:

- Check local guidelines and legislation on biohazard and biosafety precaution and refer to the best practices published in the World Health Organization (WHO) Laboratory Biosafety Manual, third edition
- Registration, Evaluation, Authorization and Restriction of Chemicals (REACH)

Documentation and Support

Obtaining SDSs

Safety Data Sheets (SDSs) are available at: <https://acdbio.com/documents/product-documents>. For the SDSs of chemicals not distributed by Advanced Cell Diagnostics, contact the chemical manufacturer.

Obtaining support

For the latest services and support information, go to: <https://acdbio.com/technical-support/support-overview>.

At the website, you can:

- Access telephone and fax numbers to contact Technical Support and Sales facilities.
- Search through frequently asked questions (FAQs).
- Submit a question directly to Technical Support.
- Search for user documents, SDSs, application notes, citations, training videos, and other product support documents.
- Find out information about customer training events.

Contact information

Advanced Cell Diagnostics, Inc.
 7707 Gateway Blvd Suite 200
 Newark, CA 94560
 Toll Free: 1-877-576-3636
 Direct: 1-510-576-8800
 Fax: 1-510-576-8801
 Information: info.acd@bio-techne.com
 Orders: orders.acd@bio-techne.com
 Support Email : support.acd@bio-techne.com

Limited product warranty

Advanced Cell Diagnostics, Inc. and/or its affiliate(s) warrant their products as set forth in the ACD General Terms and Conditions of Sale found on the ACD website. If you have any questions, please contact Advanced Cell Diagnostics at: <https://acdbio.com/about/contact>.

Headquarters

7707 Gateway Blvd Suite 200, Newark, CA 94560
For support, email support.acd@bio-technne.com
www.acdbio.com

Phone 1-510-576-8800

Toll Free 1-877-576-3636

