

RNAscope[™] 2.5 LS and LSx Reagent Kits – **RED**

For use with BOND RX System, from Leica Biosystems

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

Trademarks

RNAscope is a trademark of Advanced Cell Diagnostics, Inc. BOND RX is a registered trademark of Leica Biosystems. 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

© 2025. Advanced Cell Diagnostics, Inc. All rights reserved.

Contents

Chapter 1. Product Information	6
About this guide	6
Product description	6
Background Overview	6 6
Kit contents and storage	7
RNAscope 2.5 LS Probes RNAscope 2.5 LS Reagent Kit – RED (Cat. No. 322150) RNAscope 2.5 LSx Reagent Kit – RED (Cat. No. 322750)	7 8 8
Additional reagents required for the protease-free workflow Required materials from Leica Biosystems for the BOND RX	9 9
Equipment User-supplied materials	10 10
Chapter 2. Before You Begin	12
Important procedural guidelines	12
Chapter 3. Prepare Samples	13
Prepare FFPE sections	13
Materials required Fix the sample Dehydrate, embed, and cut the sample	13 13 13
Prepare fixed-frozen sections	14
Materials required Fix samples	14 14
Freeze tissues Prepare sections Debydrate and dry the sections	14 15
Prepare fresh-frozen sections	16
Materials required Prepare fresh frozen tissue sections	16 16
Fix the sections Dehydrate the sections	16
Chapter 4. Determine Pretreatment Conditions	
Pretreat FFPE sections	18
Target retrieval	18
RNAscope 2.5 LS and LSx Reagent Kits – RED User Manual UM 322150/Rev D/ Effective Date (04/24/2025

Permeabilization	
Tissue pretreatment recommendations	20
Pretreat fixed-frozen sections	20
Target Retrieval	20
Permeabilization	20
Tissue pretreatment recommendations	
Pretreat fresh-frozen sections	21
Target Retrieval	
Tissue pretreatment recommendations	
issue prededukcia recontinentatuons	
Chapter 5. Staining Protocols and RNA Probes	
RNA probe(s)	22
Staining Protocols	22
Workflow	23
Edit *ACD 2.5 Red Rev B (first use only)	23
Create RNA probe(s)	24
Chapter 6. Preparing the reagents	
RNAscope 2.5 LS Reagent Kit reagents (Cat. No. 322150)	27
Materials required	
Prepare the instrument reagents	
RNAscope 2.5 LSx Reagent Kit-RED reagents (Cat. No. 322750)	29
Materials required	29
Prepare the instrument reagents	29
Chapter 7. Running the Assay	
Workflow	31
Prepare the instrument	
Set up a study	
Set up slides	
Start the run	
Complete the run	
Dry and mount the samples	
Chapter 8. Evaluate the Results	
Scoring guidelines	
Control example	
Troubleshooting	
-	

Appendix A. Pretreatment Guidance for FFPE Samples	40
Tissue-specific pretreatment conditions	40
Appendix B. Edit a Heat Treatment Protocol	42
Create a prestaining protocol	42
Appendix C. Edit an Enzyme Pretreatment Protocol	45
Appendix D. Slide Setup for Additional Tissue Types	47
Fixed-frozen tissues	47
Fresh-frozen tissues	47
Protease-free workflow	48
Protease workflow	48
Appendix E. Safety	49
Chemical safety	49
Biological hazard safety	49
In the U.S.:	49
In the EU:	50
Documentation and Support	51
Obtaining SDSs	51
Obtaining support	51
Contact information	51
Limited product warranty	51



Chapter 1. Product Information



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

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

About this guide

This user manual provides guidelines and protocols to use the RNAscope 2.5 LS and LSx Reagent Kits with the BOND RX Research Advanced Staining System, from Leica Biosystems. RNAscope 2.5 LS Assays are compatible with a variety of sample types.

Product description

Background

The RNAscope 2.5 LS Assays use a novel and proprietary method of *in situ* hybridization (ISH) to visualize single RNA molecules per cell in formalin-fixed, paraffin-embedded (FFPE), fixed-frozen, and fresh-frozen 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 2.5 LS Assay allows users to automate the highly sensitive RNAscope Assay using the BOND RX System.

Overview

Figure 1 illustrates the RNAscope 2.5 LS Assay procedure, which can be completed on the instrument in ~9–10 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 the hybridization of alkaline phosphatase (AP)-labeled probes and detection using the Fast Red chromogenic substrate. Each single RNA transcript appears as a distinct dot of chromogen precipitate visible using a common bright-field microscope.

Figure 1. Procedure overview

	ZZ Target RNA-Specific Oligo Prot 2 z Z Z Z Z Z Z 2 z Z Z Z Z Z Z 2 z Z Z Z Z Z Z Z Z Z Z Z Z Z Z Z Z Z Z	PreAMP AMP Labe	I Probe
1: Tissue section	2: Hybridize to target RNA	3: Amplify signal	4: Image
Start with properly prepared sections and load slides onto the instrument. Pretreat tissue to allow access to target RNA.	Hybridize gene-specific probe pairs to the target mRNA.	Apply 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 2.5 LS RED Assay requires the RNAscope 2.5 LS Probes and the RNAscope 2.5 LS Reagents available from Advanced Cell Diagnostics. The reagents are available in two different formats:

- RNAscope 2.5 LS Reagent Kit RED (Cat. No. 322150)
- RNAscope 2.5 LSx Reagent Kit RED (Cat. No. 322750) and RNAscope 2.5 LS Pro Reagents (Cat. No. 322020)

RNAscope 2.5 LS Probes

The RNAscope 2.5 LS Probes consist of the user-specified Target Probe and the Positive and Negative Control Probes. Visit **https://www.bio-techne.com/reagents/rnascope-ish-technology** to find a gene-specific Target Probe or control probes or order a custom probe.

RNAscope 2.5 LS singleplex assays can only be used with C1 probes. These probes either have a C1 designation in their name or no designation at all. A probe with a C2, C3 or C4 designation in its name is not compatible with this chemistry.

Each probe is sufficient for staining ~30 standard slides. The probes have a shelf life of two years from date of bulk manufacturing when stored as indicated in the following table:

Target Probes						
V	Reagent	Cat. No.	Content	Quantity	Storage	
	RNAscope 2.5 LS Target Probe – [species] – [gene]	Various	Probe targeting specific RNA	16 mL x 1 bottle	2–8°C	
Control Probes						
Ø	Reagent	Cat. No.	Content	Quantity	Storage	
V	Reagent RNAscope 2.5 LS — Positive Control Probe – [<i>species</i>] – PPIB	Cat. No. Various	Content Probe targeting common housekeeping gene	Quantity 16 mL x 1 bottle	Storage 2–8°C	

RNAscope 2.5 LS Reagent Kit – RED (Cat. No. 322150)

The RNAscope 2.5 LS Reagent Kit – RED (Cat. No. 322150) contains 2 boxes (322030 and 322020) all the reagents needed to run the RNAscope 2.5 LS Assay on the BOND RX System, except for the RNA-specific probes. The kits provide enough reagents to stain ~60 standard slides.

The reagents are shipped in Ready-To-Use (RTU) bottles and must be transferred to BOND Open Containers by the user for use on the BOND RX instrument. Use the reagents in conjunction with the BOND Polymer Refine RED Detection Kit (DS9390). The RNAscope reagents are stored as indicated in the following table:

RNAscope 2.5 LS Reagent Kit – RED (Cat. No. 322150)				
Ø	Reagent	Quantity	Storage	
	RNAscope 2.5 LS Hydrogen Peroxide	21 mL x 1 bottle	2–8°C	
	RNAscope 2.5 LS Protease III	21 mL x 1 bottle	2–8°C	
	RNAscope 2.5 LS PretreatPro™	28 mL x 1 bottle	2–8°C	
	RNAscope 2.5 LS AMP 1	21 mL x 1 bottle	2–8°C	
	RNAscope 2.5 LS AMP 2	21 mL x 1 bottle	2–8°C	
	RNAscope 2.5 LS AMP 3	21 mL x 1 bottle	2–8°C	
	RNAscope 2.5 LS AMP 4 - RED	21 mL x 1 bottle	2–8°C	
	RNAscope 2.5 LS AMP 5 – RED	21 mL x 1 bottle	2–8°C	
	RNAscope 2.5 LS AMP Pro TM	21 mL x 1 bottle	2–8°C	
	RNAscope 2.5 LS AMP 6 – RED	21 mL x 1 bottle	2–8°C	
	RNAscope 2.5 LS Rinse	29 mL x 2 bottles	2–8°C	
	RNAscope 2.5 LS Bluing Reagent*	21 ml x 1 bottles	2–8°C	

* Bluing is optional.

IMPORTANT! Use only RNAscope 2.5 LS Probes. Do not substitute the reagent components of the RNAscope 2.5 LS Reagent Kit with those of any other RNAscope Reagent Kits.

RNAscope 2.5 LSx Reagent Kit – RED (Cat. No. 322750)

RNAscope 2.5 LS and LSx Reagent Kits - RED User Manual

The RNAscope 2.5 LSx Reagent Kit – RED (Cat. No. 322750) contains all the reagents needed to run the RNAscope 2.5 LS Assay on the BOND RX System, except for the RNA-specific probes and RNAscope 2.5 LS Pro Reagents (Cat. No. 322020). The kit includes two reagent trays consisting of containers pre-filled with the RNAscope reagents. The kits provide enough reagents to stain ~60 standard slides. The LSx RED Kit replaces the BOND Polymer Refine RED kit. To use the existing protease-free assay workflow, the RNAscope 2.5 LS Pro Reagents (Cat. No. 322020) kit must be transferred to BOND Open containers by the user.

RNAscope 2.5 LSx Reagent Kit – RED (Cat. No. 322750)				
V	Reagent Name	Abbreviated Name	Volume (mL)	Storage
	RNAscope 2.5 LSx H2O2	Tray 2, position 1	9	2–8°C
	RNAscope 2.5 LSx Protease	Tray 2, position 8	12	2–8°C
	RNAscope 2.5 LSx AMP 1 RED	Tray 2, position 2	18	2–8°C
	RNAscope 2.5 LSx AMP 2 RED	Tray 2, position 3	18	2–8°C
	RNAscope 2.5 LSx AMP 3 RED	Tray 2, position 4	18	2–8°C
	RNAscope 2.5 LSx AMP 4 RED	Tray 2, position 5	18	2–8°C
	RNAscope 2.5 LSx AMP 5 RED	Tray 2, position 6	18	2–8°C
	RNAscope 2.5 LSx AMP 6 RED	Tray 2, position 7	18	2–8°C
	RNAscope 2.5 LSx Rinse	Tray 1, position 1-2	27 x 2	2–8°C
	RNAscope 2.5 LSx Bluing Reagent*	Tray 1, position 6	9	2–8°C
	RNAscope 2.5 LSx Hematoxylin	Tray 1, position 3	9	2–8°C
	RNAscope 2.5 LSx Red Part A	Tray 1, position 4	4.5	2–8°C
	RNAscope 2.5 LSx Red Part B	Tray 1, position 5	1	2–8°C
	RNAscope 2.5 LSx Red Part C	Tray 1, position 6	1	2–8°C
	RNAscope 2.5 LSx Red Part D	Tray 1, position 7	30.5	2–8°C
* D1				

Store the RNAscope 2.5 LSx Reagent Kit – RED at 2–8°C.

* Bluing is optional.

IMPORTANT! Use only RNAscope 2.5 LS Probes. Do not substitute the reagent components of the RNAscope 2.5 LSx Red Reagent Kit with those of any other RNAscope Reagent Kits.

Additional reagents required for the protease-free workflow

RNAscope 2.5 LS Pro Reagents Kit (Cat. No. 322020)			
V	Reagent	Quantity	Storage
	RNAscope 2.5 LS PretreatPro	28 mL x 1 bottle	2–8°C
	RNAscope 2.5 LS AMP Pro	21 mL x 1 bottle	2–8°C

Required materials from Leica Biosystems for the BOND RX.

The RNAscope 2.5 LS Assay requires specific materials and equipment available only from Leica Biosystems.

\checkmark	Component	Cat. No.	Storage
	BOND Open Containers 30 mL	OP09700	Room temp (20–25°C)
	BOND Universal Covertiles (pack of 160)	S21.4611	Room temp (20–25°C)
	BOND Epitope Retrieval Solution 1-1L (RTU)	AR9961	2–8°C
	BOND Epitope Retrieval Solution 2-1L (RTU)	AR9640	2–8°C
	BOND Dewax Solution – 1L (RTU)	AR9222	2–26°C
	BOND Wash Solution 10X Concentrate – 1L	AR9590	2–8°C
	BOND Polymer Refine Red Detection*	DS9390	2–8°C
	BOND Aspirating Probe Cleaning System	CS9100	2–8°C
	BOND Mixing Stations	S21.1971	Room temp (20–25°C)

* Do not substitute with any other chromogen kit.

Equipment

Ø	Component	Cat. No.
	BOND RX System – automated slide stainer	_

User-supplied materials

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

V	Description	Supplier	Cat. No.
	SuperFrost Plus Slides (required)	Fisher Scientific	12-550-15
	95% Ethanol (EtOH)	American Master Tech Scientific/MLS*	ALREA95
	Xylene	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 (optional)	MLS	_
	Water bath or incubator, capable of holding temperature at 40 +/- 1°C	MLS	_
	BioCare EcoMount or Vector Labs Vectamount	ACD (or directly from manufacturer)	320409 or 321584
	Tissue-Tek® Vertical 24 Slide Rack	American Master Tech Scientific/MLS	LWS2124
	Tissue-Tek Staining Dish (4 required)	American Master Tech Scientific/MLS	LWS20WH
	Tissue-Tek Clearing Agent Dish, xylene resistant (2 required)	American Master Tech Scientific/MLS	LWS20GR
	Cover Glass 24 x 50 mm	Fisher Scientific/MLS	12-545-F
	Distilled water	MLS	_

RNAscope 2.5 LS and LSx Reagent Kits - RED User Manual

ß	Description	Supplier	Cat. No.
	Fume hood	MLS	_

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



Chapter 2. Before You Begin

Prior to running the RNAscope 2.5 LS Assay on your samples for the first time, we recommend that you:

- Become familiar with BOND RX Research Advanced Staining System. Refer to the *BOND RX 7 User Manual*.
- Run the assay on RNAscope Control Slides (Cat. No. 310045 for Human HeLa Cell Pellet, and Cat. No. 310023 for Mouse 3T3 Cell Pellet) using the RNAscope 2.5 LS Positive and Negative Control Probes.

Important procedural guidelines

- Start with properly fixed and prepared sections. Refer to Chapter 3. Prepare Samples for preparation of FFPE, fixed-frozen or fresh-frozen slides. For preparation of other sample types, contact support.acd@bio-techne.com.
- Regularly maintain and clean your automated staining instrument.
- Always run positive and negative control probes on your sample to assess sample RNA quality and optimal permeabilization.
- Do not substitute required materials. Assay has been validated with these materials only.
- Follow the protocol exactly for the best results.
- Use good laboratory practices and follow all necessary safety procedures. Refer to **Appendix E. Safety** for more information.



Chapter 3. Prepare Samples

IMPORTANT! The following protocols describe formalin-fixed, paraffin-embedded (FFPE), fixed frozen and fresh frozen sample preparation. We highly recommend following these guidelines. We cannot guarantee assay results with other preparation methods.

Prepare FFPE sections

Materials required

- 10% neutral buffered formalin (NBF)
- 1X PBS
- Paraffin wax
- 95% Ethanol (EtOH)
- Xylene
- Microtome
- Water bath
- SuperFrost Plus slides

Fix the sample

- 1. Immediately following dissection cut the tissue into blocks of 3–4 mm in thickness.
- 2. Place the tissue blocks into fixative within **1 HR** of biopsy.
- 3. Fix the 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 RNAscope 2.5 LS Assay.

Dehydrate, embed, and cut the sample

IMPORTANT! Use fresh reagents.

- 1. Wash sample with 1X PBS.
- 2. Dehydrate sample using a standard ethanol series, followed by xylene.
- 3. Embed sample in paraffin using standard procedures.

Note: Embedded samples may be stored at room temperature with desiccation. To better preserve RNA quality over a long period (>1 yr), storing at 2–8°C with desiccation is recommended.

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



Tissue section location

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

6. Air dry slides **OVERNIGHT** at **RT**.

OPTIONAL STOPPING POINT. Use sectioned tissue within three months. Store sections with desiccants at room temperature.

Prepare fixed-frozen sections

Materials required

- 1X PBS
- 10% Neutral Buffered Formalin (NBF) or 4% Paraformaldehyde (PFA)
- 100% alcohol (EtOH)
- Tissue-Tek Vertical 24 Slide Rack
- Tissue-Tek Staining Dishes
- Drying oven

Fix samples

- 1. If needed, perfuse the tissue with freshly prepared 4% paraformaldehyde (PFA) in 1X PBS, or go directly to Step 2.
- 2. Dissect the tissue and fix in freshly prepared 4% PFA for 24 HRS at 4°C.

Freeze tissues

- 1. Immerse the tissue in 10% sucrose in 1X PBS at 4°C until the tissue sinks to the bottom of the container (approximately **18 HRS** for brain tissue).
- 2. Repeat this step with 20% sucrose in 1X PBS, followed by 30% sucrose in 1X PBS, each time allowing the tissue to sink to the bottom of the container.

UM 322150/Rev D/ Effective Date 04/24/2025

3. Freeze the tissue in Optimal Cutting Temperature (OCT) embedding media with dry ice or liquid nitrogen, and store it in an airtight container at **-80°C**.

Prepare sections

Before tissue sectioning, equilibrate the tissue blocks at -20°C for at least 1 HR in a cryostat.

 Section the blocks by cutting 7–15 μm thick sections. 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.

2. Air dry the slides for 60 –120 MIN at –20°C.

OPTIONAL STOPPING POINT (1). Use sectioned tissue within **3 MONTHS**. Store sections with desiccants at **-80°C**.

- 3. Wash the slides with 200 mL 1X PBS in a Tissue-Tek slide rack for **5 MIN** while moving the rack up and down to remove OCT.
- 4. Bake the slides for 60 MIN at 60°C.
- 5. Immediately post-fix the slides by immersing them in prechilled 10% NBF or 4% PFA in 1X PBS for **60 MIN** at **RT**.

Note: If you are experiencing issues with sample detachment, extending the post-fixation and baking times may be helpful.

Dehydrate and dry the sections

Reagents may be prepared ahead of time. Ensure all containers remain covered.

- 1. Prepare 200 mL 50% ethanol, 200 mL 70% ethanol, and 2X 200 mL 100% ethanol in Tissue Tek Staining Dishes.
- 2. Remove the slides from the 10% NBF or 4% PFA, and immerse them in 50% EtOH for **5 MIN** at **RT**.
- 3. Place the slides in 70% ethanol for 5 MIN at RT.
- 4. Place the slides in 100% ethanol for **5 MIN** at **RT**.
- 5. Place slides in fresh 100% ethanol for **5 MIN** at **RT**.
- 6. Remove slides from ethanol, and let them dry for 5 MIN at RT.

Prepare fresh-frozen sections

Materials required

- 1X PBS
- 10% Neutral Buffered Formalin (NBF) or 4% Paraformaldehyde (PFA)
- 100% alcohol (EtOH)
- Tissue-Tek Vertical 24 Slide Rack
- Tissue-Tek Staining Dishes

Prepare fresh frozen tissue sections

Remove tissue and cut to fit into cryomolds.

CAUTION! Handle biological specimens appropriately.

- 1. Freeze the specimen within 5 MIN of tissue harvest.
- 2. Embed the frozen tissue in cryo-embedding medium (OCT):
 - a. Add two drops of OCT into a cryomold.
 - b. Place the frozen tissue on the OCT in the correct orientation for cutting.
 - c. Add more OCT to fill the cryomold. Do not allow any air bubbles to form.
 - d. Hold the block with forceps on the surface of the liquid nitrogen or isopentane cooled by dry ice or liquid nitrogen or place the cryomold on dry ice.
- 3. Store the frozen block in an air-tight container at **-80°C** prior to sectioning.

Embedded tissue may be stored for up to three months.

OPTIONAL STOPPING POINT (1). Section tissue within 3 MONTHS.

- 4. Section the block:
 - a. Equilibrate block to -20°C in a cryostat ~1 HR.
 - b. Cut 10-20 µm thick sections and mount onto SUPERFROST® PLUS SLIDES.
 - c. Dry the sections at 60 –120 MIN at –20°C to retain tissue adherence.
- 5. Store the sections in slide boxes wrapped air-tight with aluminum foil or zip-lock bags at **-80°C** until use.

Sections may be stored for up to three months.

IMPORTANT! Do not fix the slides prior to this step.

OPTIONAL STOPPING POINT (2). Use sectioned tissue within 3 MONTHS.

Fix the sections

- 1. Remove fresh-frozen tissue slides from 80°C and place in a Tissue Tek Slide Rack.
- 2. Immediately immerse the slides in 200 mL of 10% NBF or freshly prepared 4% PFA.
- 3. Incubate the slides for at least **90 MIN** at **ROOM TEMPERATURE (RT)**.



Note: Formalin that has been stored for more than six months, exposed to air for more than a week, or used repeatedly may result in suboptimal tissue fixation. 4% PFA must be freshly prepared for each experiment.

Dehydrate the sections

Reagents may be prepared ahead of time. Ensure all containers remain covered.

- 1. Prepare 200 mL 50% ethanol, 200 mL 70% ethanol, and 2X 200 mL 100% ethanol in Tissue Tek Staining Dishes.
- 2. Place the slides in 50% ethanol for 5 MIN at RT.
- 3. Place the slides in 70% ethanol for **5 MIN** at **RT**.
- 4. Place the slides in 100% ethanol for **5 MIN** at **RT**.
- 5. Place slides in fresh 100% ethanol for **5 MIN** at **RT**.

Note: If needed, slides can be stored in 100% EtOH at -20°C for up to 1 week. Prolonged storage may degrade sample RNA.

6. Remove slides from ethanol, and let them dry for 5 MIN at RT.



Chapter 4. Determine Pretreatment Conditions

The following protocols describe formalin-fixed, paraffin-embedded (FFPE), fixed frozen and fresh-frozen 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.

Pretreat FFPE sections

Target retrieval

FFPE samples must be de-crosslinked with a target retrieval step. The RNAscope 2.5 LS Assays uses the BOND RX's ER2 solution exclusively for this step.

Permeabilization

Two options are available:

- Protease-based permeabilization is recommended for experiments that stain only RNA. This option uses LS Protease III or RNAScope 2.5 LSx Protease depending on the kit you are using.
- Protease-free permeabilization uses the LS PretreatPro reagent which is free of protease enzyme. This allows co-detection of RNA and proteins that were previously incompatible with protease on the same tissue section using immunohistochemistry (IHC).



To determine the correct permeabilization option, please refer to the following flowchart:

** For optimizing ISH signal strength for low RNA expressors in the protease free workflow.

Tissue pretreatment recommendations

Use these conditions as a starting point when tissues are prepared as described in **Chapter 3**. Depending on your tissue type, vary the amount of time for the ER2 and Protease or LS PretreatPro until positive RNA control signal is maximized with minimal/no negative RNA control signal (see **Appendix A** for a list of tissues).

Reagent	Mild Protease Workflow	Standard Protease Workflow	Protease-Free Workflow
BOND ER2*	15 MIN at 88°C	15 MIN at 95°C	15 MIN at 95°C
LS PretreatPro†			30 MIN at 95°C
LS Protease III OR RNAscope 2.5 LSx Protease‡	15 MIN at 40°C	15 MIN at 40°C	

* Sample types, such as certain xenografts and cell pellets, might require shorter incubation time. For these tissue types, reduce the BOND ER2 incubation time. Protease/ LS PretreatPro incubation times can also be adjusted but is rarely needed. Contact ACD Support at **support.acd@bio-techne.com** if you need guidance on your tissue type.

+ For ACD Control Cell pellets, we recommend starting with ER2 at 92°C for 15mins when using the LS PretreatPro workflow.

‡ Depending on which assay kit format you are using; user-filled OR pre-filled.

Pretreat fixed-frozen sections

Target Retrieval

Fixed-frozen samples must be gently de-crosslinked with a target retrieval step. The RNAscope 2.5 LS Assay uses the BOND RX's ER2 solution exclusively for this step.

Permeabilization

Only protease has been tested for use with fixed-frozen sections. Check with ACD Support for any updates.

Tissue pretreatment recommendations

Use these conditions as a starting point when tissues are prepared as described in **Chapter 3**. Depending on your tissue type, vary the amount of time for the ER2 and/or Protease until positive control RNA signal is maximized with minimal or no negative RNA control signal (see **Appendix E** for details).

Reagent	Standard
BOND ER2*	5 MIN at 95°C
LS Protease III or RNAscope 2.5 LSx Protease [†]	15 MIN at 40°C

* Sample types, such as certain xenografts and cell pellets, might require shorter incubation time. For these tissue types, reduce the BOND ER2 incubation time. Protease time can also be adjusted but is rarely needed. Contact ACD Support at **support.acd@bio-techne.com** if you need guidance on your tissue type.

+ You might need to create this enzyme treatment protocol. Please refer to **Appendix B** for further instructions.

Pretreat fresh-frozen sections

Target Retrieval

The retrieval requirements for fresh-frozen sections are determined by the permeabilization approach that is followed. Target retrieval is not needed when using protease. However, a short target retrieval is required to optimize the signal when using PretreatPro.

Permeabilization

ACD recommends starting with the PretreatPro workflow as all reagents are included in the kit purchases. If protease treatment is needed, then LS Protease IV needs to be purchased separately for optimal assay performance.

Tissue pretreatment recommendations

Use these conditions as a starting point when tissues are prepared as described in Chapter 3.

Protease-Free Workflow

Reagent	Standard
BOND ER2	5 MIN at 95°C*
LS PretreatPro	30 MIN at 40°C

* You might need to create this heat pretreatment protocol. Please refer to **Appendix B** for further instructions.

Protease Workflow

Reagent	Standard
LS Protease IV (ACD Part Number 322140)	30 MIN at ambient temperature*

*You might need to create this enzyme treatment protocol. Please refer to **Appendix C** for further instructions.



Chapter 5. Staining Protocols and RNA Probes

Use the instructions in this chapter to set up the RNAscope 2.5 LS Assay using software version 6.0 and above.

IMPORTANT! BXD42 or higher is required to run the following setup on software version 6.0 and above. Please contact your Leica FAS to upgrade to BXD42 before starting the assay.

IMPORTANT! We strongly recommend you run the Control Slides (Cat. No. 310045 or Cat. No. 310023) using the RNAscope 2.5 LS positive and negative control probes along with your samples in every run.

RNA probe(s)

This workflow uses the standard BOND RX software setup. Probes are manually registered in the software as 'Probe RNA' (not as 'Ancillary'), and a pre-defined staining protocol will be selected as the default protocol for this probe. Follow the steps in this chapter to enable the workflow. Your ACD Field Application Specialist (FAS) can help you implement this procedure.

Staining Protocols

Staining protocols have been pre-defined in the software (*) for the four different RNAscope RED assays, depending on the reagent kit and enzyme retrieval used. A summary of each staining protocol follows:

- *ACD 2.5 Red Rev B: This protocol has been pre-defined for use with:
 - RNAscope 2.5 LS Reagent Kit -RED (Cat. No. 322150) user-filled
 - BOND Polymer Refine Red Detection (DS9390)
 - Protease Enzyme Retrieval
- *ACD 2.5 Red Rev B_Protease-Free: This protocol has been pre-defined for use with:
 - RNAscope 2.5 LS Reagent Kit-RED (Cat. No. 322150) user-filled
 - BOND Polymer Refine Red Detection (DS9390)
 - PretreatPro (protease-free) Enzyme Retrieval
- *RNAscope 2.5 LSx RED ISH: This protocol has been pre-defined for use with:
 - RNAscope 2.5 LSx RED ISH Kit (Cat. No. 322750)
 - Protease Enzyme Retrieval
- *RS2.5 LSx RED_Protease-Free: This protocol has been pre-defined for use with:
 - RNAscope 2.5 LSx RED ISH Kit (Cat. No. 322750)
 - PretreatPro (protease-free) Enzyme Retrieval



Slide setup and supporting protocols for HIER, EIER and Hybridization are described in more detail in **Chapter 6**.

Workflow



Edit *ACD 2.5 Red Rev B (first use only)

This edit removes the first 14 reagent and wash steps that are no longer needed in the staining protocol when using software version 6.0 or greater.

- 1. In the Protocol setup screen, select **Staining** from the Protocol group menu.
- 2. Highlight the ***ACD 2.5 Red Rev B** protocol. Select **Copy**.
- 3. Change the protocol name for your first probe to **ACD 2.5 Red Rev B Protease** in the Name text box, **25RRBPt** in the Abbreviated name text box, and **ACD 2.5 Red Rev B protocol Protease** in the Description text box.
- 4. Highlight the default probe reagent lines (Lines 1-3: *ACD 2.5 P1) and the following 11 wash steps and select **Delete step** at the bottom right of the window. Steps must be deleted one at a time. Once this is complete, *ACD Amp1 should be the first step of the protocol.
- 5. Select **Preferred** in the bottom right corner of the window.
- 6. Select Save.
- 7. Click on **Next** to proceed. Acknowledge any pop-up warnings that may appear on the screen by clicking **OK**.

The following screenshot shows the protocol after editing. Moving forward, use this new protocol, which means you can deselect the ***ACD 2.5 Red Rev B** protocol as the preferred protocol.

New protocol properties

Name:	ACD 2.5 Red Rev B Protease					
Abbreviated name:	2.5RRBPt					
Description:	ACD 2.5 Red Rev B protocol	Protease				
Staining method:	< Single < Preliminar	ry 📝 Final			✓ Prefer	red
BOND RX ^m	BOND RX			Import protocol	Protocol type: ISH detect	tion
Preferred detection s	Bond Polymer Ref	fine Red Detection				
Step N° Was	h Reagent	Supplier	Ambient Temperature	Inc. (min)	Dispense type	
1	*ACD Amp 1	Advanced Cell Diagnostics	42	1:00	150 µL	Î
2	*ACD Amp 1	Advanced Cell Diagnostics	42	30:00	150 µL	1
3 🗸	*Bond Wash Solution	Leica Microsystems	4	0:00	150 µL	
4 🖌	*Bond Wash Solution	Leica Microsystems	4	0:00	150 µL	
5 🖌	*Bond Wash Solution	Leica Microsystems	~	0:00	150 μL	
6 🗸	*Bond Wash Solution	Leica Microsystems	*	3:00	150 µL	
7 🖌	*Bond Wash Solution	Leica Microsystems	4	3:00	150 µL	
8 🗸	*Bond Wash Solution	Leica Microsystems	~	0:00	150 µL	
۵ ،	*D	Laina Minanuntama		0.00	450	*
Show wash ste	ps			Insert wash	Insert reagent Delete s	tep

Create RNA probe(s)

For each probe not already in the database, create it with the default workflow protocols that are appropriate for your needs. To ensure efficiency, select the protocols you anticipate using most frequently as your defaults. These defaults can be modified during slide setup if any slides require different protocols.

See the following workflow examples:

Name:	ACD Probe 1 LS	
Abbreviated name:	ACDP1 LS	
Туре:	Probe RNA	*
Supplier:	ACD	
Staining method:	Single/Sequentia	al multiplex
Single P	reliminary Fin	al
Default staining protoc	:ol:	ACD 2.5 Red Rev B Protease
Default HIER protocol:		*ACD HIER 15 min with ER2 (95)
Default enzyme protoc	col:	*ACD 15 min Protease
Default denaturation p	rotocol:	*****
Default hybridization p	n protocol: *RNAscope 2.5 LSx Hybridization	
Compatible bulks:		
BWash		
Preferred	Hazardous	
	Caus	Canaal

User-filled RNAscope 2.5 LS reagents using protease pretreatment

User-filled RNAscope 2.5 LS reagents using protease-free pretreatment

		Add reagent		
Name:	ACD Probe 1 LS			
Abbreviated name:	ACDP1 LS			
Туре:	Probe RNA	A 👻		
Supplier:	ACD			
Staining method:	Single/Sequential multiplex			
Single	Preliminary	Final		
Default staining proto	col:	*ACD 2.5 Red_Protease-Free		
Default HIER protocol:		*ACD HIER 15min with ER2 (95)		
Default enzyme proto	col:	*ACD PretreatPro		
Default denaturation	protocol:	*		
Default hybridization	*RNAscope 2.5 LSx Hybridization			
Compatible bulks:				
*BWash				
✓ Preferred ✓	Hazardous			
	s	ave Cancel		

RNAscope 2.5 LSx Reagents kit using protease pretreatment

		Add reagent	
Name:	ACD Probe 1 LSx		
Abbreviated name:	ACDP1LSx		
Туре:	Probe RNA	↓	
Supplier:	ACD		
Staining method:	Single/Sequential multiplex -		
Single P	reliminary	Final	
Default staining protoc	ol:	*RNAscope 2.5 LSx RED ISH	•
Default HIER protocol:		*RNAscope 2.5 LSx Target Retrieval (95)	
Default enzyme protoc	ol:	*RNAscope 2.5 LSx Enzyme	
Default denaturation p	*		•
Default hybridization p	*RNAscope 2.5 LSx Hybridization		•
Compatible bulks:			
*BWash			
Preferred	Hazardous		
	S	ave Cancel	

RNAscope 2.5 LSx Reagents using protease-free pretreatment

	1	Add reagent	
Name:	ACD Probe 1 LSx		
Abbreviated name:	ACDP1LSx		
Туре:	Probe RNA	-	
Supplier:	ACD		
Staining method:	Single/Sequential multiplex		
Single P	reliminary	Final	
Default staining protoc	:ol:	*RS 2.5 LSx Red_Protease-Free	
Default HIER protocol:		RNAscope 2.5 LSx Target Retrieval (95)	
Default enzyme protoc	col:	*RS 2.5 LSx_PreTPro	
Default denaturation p	rotocol:	*	
Default hybridization p	rotocol:	*RNAscope 2.5 LSx Hybridization	
Compatible bulks:			
*BWash			
Preferred	Hazardous		
	S	ave	
	Sa	ave Cancel	



Chapter 6. Preparing the reagents

RNAscope 2.5 LS Reagent Kit reagents (Cat. No. 322150)

Materials required

	Materials provided by Advanced Cell Diagnostics	Materials provided by Leica Biosystems	Materials provided by user
· · · · · · ·	Materials provided by Advanced Cell DiagnosticsRNAscope 2.5 LS Target ProbeRNAscope 2.5 LS Positive Control ProbeRNAscope 2.5 LS Negative Control ProbeRNAscope 2.5 LS Hydrogen PeroxideRNAscope 2.5 LS Protease IIIRNAscope 2.5 LS Protease IIIRNAscope 2.5 LS AMP 1RNAscope 2.5 LS AMP 2RNAscope 2.5 LS AMP 3RNAscope 2.5 LS AMP 4 - REDRNAscope 2.5 LS AMP 5 - REDRNAscope 2.5 LS AMP 6 - RED	Materials provided by Leica Biosystems BOND RX System • Stainer Bulk Reagents • BOND Wash Solution, 10X • BOND Dewax Solution • BOND Epitope Retrieval Solution 1 • BOND Epitope Retrieval Solution 2 Reagents • BOND Polymer Refine Red Detection	 Materials provided by user Distilled water Drying oven Fume hood Tissue-Tek Staining Dish Tissue-Tek Clearing Agent Dish, xylene-resistant (2) Tissue-Tek Vertical 24 Slide Rack Drying oven EcoMount or VectaMount Cover glass, 24 mm x 50 mm
• •	RNAscope 2.5 LS AMP Pro RNAscope 2.5 LS Rinse RNAscope 2.5 Bluing Reagent		

Prepare the instrument reagents

1. Label the BOND Open Containers as shown in the following table.

Note: The table lists the required container names. If unsure, check your protocols to confirm the reagent names and label your containers accordingly.

RNAscope Reagents	BOND RX Container Name
RNAscope 2.5 LS Hydrogen Peroxide	*Open 0 Haz
RNAscope 2.5 LS Protease III (for protease retrieval only)	*ACD Enzyme
RNAscope LS PretreatPro (protease-free retrieval only)	*RNAscope LS PreteatPro
RNAscope 2.5 LS AMP 1	*ACD Amp 1
RNAscope 2.5 LS AMP 2	*ACD Amp 2
RNAscope 2.5 LS AMP 3	*ACD Amp 3
RNAscope 2.5 LS AMP 4	*ACD Amp 4

RNAscope 2.5 LS and LSx Reagent Kits - RED User Manual

RNAscope Reagents	BOND RX Container Name
RNAscope 2.5 LS AMP 5 – RED (for protease assay only)	*ACD Amp 5 Red
RNAscope 2.5 LS AMP Pro (For protease-free assay only)	*RNAscope 2.5 LS Amp Pro
RNAscope 2.5 LS AMP 6 – RED	*ACD Amp 6 Red
RNAscope 2.5 LS Rinse	*LS Rinse
RNAscope 2.5 LS Bluing Reagent ⁺	*ACD Blue
RNAscope 2.5 LS Target Probe	Name of your choice

⁺Bluing is optional

Note: Leica BOND Polymer Refine Red Detection (DS9390) is a pre-filled BOND Detection System. The protocols use the RED Chromogen and Hematoxylin from this kit.

2. Carefully transfer all the RNAscope LS reagents into the labeled 30 mL BOND Open containers.

3. Determine the volume of probe reagent needed for the run.

Note: If using other open containers to deliver your probe, you must account for the dead volume required in each container. Suggested volumes can be found in the following table.

Open Container	Suggested Dead-Volume
30 mL	2.5 mL
7 mL	1 mL
6 mL	600 μL

- 4. Using the barcode scanner, scan the front barcode on the labeled 30 mL BOND Open container. A window will appear.
- 5. From the drop-down menu, select the corresponding name of the reagent as shown in the table above under **BOND RX Container Name**:
- 6. Enter the RNAscope 2.5 LS Reagent Kit lot number and the expiration date in their respective fields. Select **OK**.
- 7. Scan the two side barcodes on the tray for BOND Polymer Refine Red Detection. When the window opens, select **OK** to register the kit.

IMPORTANT! Do not introduce bubbles into the solutions by shaking the containers. To mix reagents, gently invert the containers several times. If bubbles are present, leave the containers out at room temperature until the bubbles dissipate.

RNAscope 2.5 LSx Reagent Kit-RED reagents (Cat. No. 322750)

Materials required

	Materials provided by Advanced Cell Diagnostics	Materials provided by Leica Biosystems		Materials provided by user
•	RNAscope 2.5 LS Target Probe	BOND RX System	•	Distilled water
•	RNAscope 2.5 LS Positive Control Probe	• Stainer	•	Drying oven
•	RNAscope 2.5 LS Negative Control Probe	Bulk Reagents	•	Fume hood
•	RNAscope 2.5 LS PretreatPro	BOND Wash Solution, 10X	•	Tissue-Tek Staining Dish
•	RNAscope 2.5 LS AMP Pro	BOND Dewax SolutionBOND Epitope Retrieval Solution 1	•	Tissue-Tek Clearing Agent Dish. xvlene-resistant (2)
•	KNAScope 2.5 LSX keagent Kit - KED	BOND Epitope Retrieval Solution 2	•	Tissue-Tek Vertical 24 Slide Rack
			•	Drying oven
			•	EcoMount or VectaMount
			•	Cover glass, 24 mm x 50 mm

Prepare the instrument reagents

1. Label the BOND Open Containers for your RNA Probes (including controls).

Note: The table lists the required container names. If unsure, check your protocols to confirm the reagent names and label your containers accordingly.

RNAscope Reagents	BOND RX Container Name
RNAscope 2.5 LS Target Probe	Name of your choice
RNAscope LS PretreatPro (for protease-free enzyme retrieval only)	*RNAscope LS PretreatPro
RNAscope 2.5 LS AMP Pro (for protease-free assay only)	*RNAscope 2.5 LS AMP Pro

2. Determine the volume of probe reagent needed for the run.

Note: If using other open containers to deliver your probe, you must account for the dead volume required in each container. Suggested volumes can be found in the following table.

Open Container	Suggested Dead-Volume
30 mL	2.5 mL
7 mL	1 mL
6 mL	600 μL

3. Using the barcode scanner, scan the front barcode on the labeled 30 mL BOND Open container. A window will appear.

- 4. From the drop-down menu, select the corresponding name of the RNA probe created in Chapter 5. Enter in a Lot# and expiration date in their respective fields. Select **OK**.
- 5. If running the protease-free assay, label BOND Open Containers for PretreatPro and AMP Pro. Scan the front barcode on these containers and select the corresponding name from the drop-down list. Select **OK**.
- 6. Transfer the RNA Probes, (and if running the protease-free assay) the LS PretreatPro and LS 2.5 AMP Pro reagents into the corresponding BOND Open Containers.
- 7. Scan the two side barcodes on Tray #1 for the RNAscope 2.5 LSx Reagent Kit-RED kit. When the window opens, select **OK** to register the tray.
- 8. Next scan the front barcode for the first reagent in Tray #2 (ie, AMP reagents). Select **OK** and then repeat for all containers in Tray #2.



Chapter 7. Running the Assay

Workflow



Prepare the instrument

1. Fill the large containers located in the bottom of the instrument with the BOND RX bulk reagents. Dilute Bond Wash Solution 10X Concentrate 1:10, with DI Water to make the 1X Bond Wash Solution (working solution).

Note: Insufficient bulk reagent volumes may lead to run failure.

IMPORTANT! Do not introduce bubbles into the solutions by shaking the containers. To mix reagents, gently invert the containers several times. If bubbles are present, leave the containers out at room temperature until the bubbles dissipate.

- 2. Use clean, dry covertiles for every run. Clean used covertiles with water, bleach, and ethanol. Air dry before reuse. See Leica Biosystems documentation for details.
- 3. Before starting a run, empty bulk waste containers. Discard waste according to all local, state/provincial, and/or national regulations.

Set up a study

1. To build a study, select the **Slide setup** icon at the top of the screen.



2. Select Add study and enter a name in the Study ID field (keep the Dispense volume at 150 μ l as shown).

Slide setup	Protocal setup	Reagent setup		Search	Help	Log out +			Leica
	Slide se	etup						Add study Edit study Delate	ludy Copy study
		Study	ID			Study name	Researcher name	Slides Slides	12 Add panel
					Sh Sh Re Sh Dis	ndy ID: rdy name: rdy comments: searcher: rdy N ⁴ : pense volume: paration protocol:	Add study *		

- 3. For FFPE tissues, select ***Bake and Dewax** as the Preparation protocol. (For alternative tissue preps, see **Appendix D**).
- 4. Select OK.

Set up slides

The slide setup differs based on the assay (protease or protease-free) and the reagents (user-filled RNAscope 2.5 LS reagents or RNAscope 2.5 LSx Reagents – RED kit).

The following screen shots indicate the protocols to select for the slide setup for each of the assay and reagent variables. Ensure you select the correct protocols.

- 1. Select **Add slide** to assign a protocol to each slide.
- 2. Enter the tissue type and probe name in the Comments field.
- 3. Choose the Staining, HIER, Enzyme and Hybridization protocols depending on the reagent kit and enzyme method you are using. Ensure you change Probe Application and Probe Removal protocols from *Default to the *RNAscope 2.5 LSx choice.

See the following four options.

CD Testing		Sinde Comments						
lesearcher:	Tissue type: Dis	pense volume:						
	V Test tissue	100 μL						
ilide ID:	Negative tissue	🧭 150 μL						
tudy N°:	Positive tissue							
tudy comments:	Staining mode:							
ate created:	Single	Routine						
/28/2025 1:41:34 PM	Single							
	Process: IHC	s 📀 ISH						
	Marker: ACD P	robe 1 LS (ACD)						
	Protocols							
	Staining:	ACD 2.5 Red Rev B Protease						
	Preparation:	*Bake and Dewax						
	HIER:	*ACD HIER 15 min with ER2 (95)						
	Enzyme:	*ACD 15 min Protease						
larkor IIPI	Probe Application:	*RNAscope 2.5 LSx Probe Application						
Auto	Denaturation:	*						
staction System UPI:	Hybridization:	*RNAscope 2.5 LSx Hybridization						
Auto	Probe Removal:	*RNAscope 2.5 LSx Probe Removal						

User-filled RNAscope 2.5 LS reagents using protease pretreatment

User-filled RNAscope 2.5 LS reagents using protease-free pretreatment

D Testing	Tiesuo tupo:	Dieponeo	so volumo:
esearcher: 	Test tissue		100 µL
lide ID:	Negative tissue	1	150 µL
tudy N°: 33	Positive tissue		
udy comments:	Staining mode:		
ate created:	Single	▼ R	Routine
/28/2025 1:41:34 PM	Single		
	Process:	нс	📀 ISH
	Marker:	ACD Probe 1	1 LS (ACD)
	Protocols		
	Staining:		*ACD 2.5 Red_Protease-Free
	Preparation:		*Bake and Dewax
	HIER:		*ACD HIER 15min with ER2 (95)
	Enzyme:		*ACD PretreatPro
larker UPI:	Probe Application:		*RNAscope 2.5 LSx Probe Application
Auto	Denaturation:		*****
etection System UPI:	Hybridization:		*RNAscope 2.5 LSx Hybridization
Auto and a system of it.	Probe Removal:		*RNAscope 2.5 LSx Probe Removal

RNAscope 2.5 LS and LSx Reagent Kits - RED User Manual

Tinnun funna Dinnan						
Tissue type: Dispen	se volume:					
V Test tissue	100 µL					
Negative tissue	150 µL					
Positive tissue						
Staining mode:						
Single	Routine					
Single						
Process: IHC	✓ ISH					
Marker: ACD Probe	⇒ 1 LSx (ACD) 🗸					
Protocols						
Staining:	*RNAscope 2.5 LSx RED ISH					
Preparation:	*Bake and Dewax					
HIER:	*RNAscope 2.5 LSx Target Retrieval (95)					
Enzyme:	*RNAscope 2.5 LSx Enzyme					
Probe Application:	*RNAscope 2.5 LSx Probe Application					
Denaturation:	*					
Hybridization:	*RNAscope 2.5 LSx Hybridization 🗸					
	Tissue type: Dispension Image: Image: Image: Image:					

RNAscope 2.5 LSx Reagents - RED kit using protease pretreatment

RNAscope 2.5 LSx Reagents - RED kit using protease-free pretreatment

CD Testing	Tissue type: D)ispense volume:
esearcher:	Test tissue	00 μL
lide ID:	Negative tissue	
tudy N°: 33	Positive tissue	
udy comments:	Staining mode:	
ate created:	Single	Routine
/28/2025 1:41:34 PM	Single	
	Process:	нс 🥑 ізн
	Marker: ACD	Probe 1 LSx (ACD)
	Protocols	
	Staining:	*RS 2.5 LSx Red_Protease-Free
	Preparation:	*Bake and Dewax
	HIER:	RNAscope 2.5 LSx Target Retrieval (95)
	Enzyme:	*RS 2.5 LSx_PreTPro
Marker UPI:	Probe Application:	*RNAscope 2.5 LSx Probe Application
Auto	Denaturation:	*
	Hybridization:	*RNAscope 2.5 LSx Hybridization
etection System IIPI-	Dub Durit	*RNAscope 2.5 LSx Probe Removal

RNAscope 2.5 LS and LSx Reagent Kits - RED User Manual

UM 322150/Rev D/ Effective Date 04/24/2025



Note: HIER protocol time and temperature varies depending on tissue type.

- 4. Select Add slide.
- 5. After adding all the slides to the study, select Close to return to the Slide setup screen.
- 6. Select **Print labels** to print barcodes to attach to the slides.



Start the run

1. Attach the barcode labels to the slides and add the slides to the slide tray with the label sides facing up.

Note: Add a covertile on top of each slide. The rectangular-shaped neck of the covertile should fit into the groove of the slide tray. Verify placement and seating of covertile.

- 2. Place the tray in the BOND RX and press the button to load the tray onto the machine.
- 3. Once the slides have been scanned, select the **PLAY** (triangular) button on the screen located under the start tray to start the run. Alternatively, right-click on scanned label images and select **Delayed Start** to start the run at a future time.

IMPORTANT! Before leaving the instrument unattended, ensure that the instrument is running successfully. In the event of a problem, please contact **support.acd@bio-techne.com** or your Field Application Scientist.

Complete the run

- 1. After the run is complete, press the button on the front of the instrument to unload the slides.
- 2. Place the slides onto the Tissue-Tek Slide Rack and move the rack into a staining dish containing distilled water.
- 3. Flush the slides with flowing tap water for a few minutes.

UM 322150/Rev D/ Effective Date 04/24/2025



4. (Optional) Perform offline bluing of choice or use the RNAscope 2.5 LS Bluing Reagent included in the kit by pipetting a small amount onto each slide for **1–2 MIN** followed by rinsing.

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 or VectaMount on the slide before the xylene dries.

IMPORTANT! Use EcoMount (Biocare Medical) or VectaMount (VectaLabs) mounting media.

- 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 5 MIN.
- 7. Proceed to Chapter 8. Evaluate the Results.



Chapter 8. Evaluate the Results

Examine tissue sections under a standard bright field microscope at 20-40X magnification:

- Assess tissue and cell morphology.
- Assess positive control signal strength. Positive control signal should be visible as punctate dots within the cell at 20–40X magnification.
- Assess negative control background. One dot to every 10 cells displaying background staining per 20X microscope field is acceptable.
- Evaluate target probe signal using the scoring guidelines in the next section.

Scoring guidelines

The RNAscope 2.5 LS Assay enables a semi-quantitative scoring guideline utilizing the estimated number of punctate dots present within each cell boundary.

An example of how to develop such a guideline for semi-quantitative assessment of RNAscope staining intensity is presented below for a gene with expression level varying between 1 to > 10 copies per cell.

Note: If your gene expression level is higher or lower than this range, you may need to scale the criteria accordingly.

Categorize staining into five grades: 0, 1+, 2+, 3+, and 4+ according to the following table:

Staining Score	Microscope Objective Scoring*
0	No staining or less than 1 dot for every 10 cells (40X magnification)
1	1–3 dots/cell (visible at 20–40X magnification)
2	4–9 dots/cell. No or very few dot clusters (visible at 20–40X magnification)
3	10–15 dots/cell and/or < 10% positive cells have dot in clusters (visible at 20X magnification)
4	>15 dots/cell and/or >10% positive cells have dot in clusters (visible at 20X magnification)

* Discount cells with artificially high nuclear background staining.

biotechne[®] / A@D

Control example

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



Figure 2. RNAscope 2.5 Assay detection of PPIB mRNA in HeLa cells

Troubleshooting

If you obtain less than satisfactory results, troubleshoot your assay by following these simple guidelines:

- For the protease assay, if you observe the presence of background staining, increase the Epitope Retrieval 2 (ER2) in increments of five minutes and/or increase the protease time in increments of ten minutes (see **Appendix B** and **C** for instructions on editing protocols).
- For the protease-free assay, if you observe lower signal intensity, increase the ER2 temperature from 88°C to 92°C for cell pellets or from 95°C to 100°C for tissues to obtain better signal intensity.
- Use the previous process for over-fixed tissues.
- The RNAscope 2.5 LS RED and LS BROWN assays use Leica Biosystems' BOND Polymer Refine Detection and BOND Polymer Refine Red Detection kits, respectively. Do not use any other chromogen kits.
- Do not shake the contents in the dispensers as this will form bubbles and may lead to weak or no staining. If bubbles are present, leave the containers out at room temperature until the bubbles dissipate.
- Do not alter the staining protocol in any way except for the hematoxylin incubation time. The parameters in the staining protocol have been optimized to run the RNAscope assay on the instrument.

For troubleshooting information, please contact technical support at support.acd@bio-techne.com.



Appendix A. Pretreatment Guidance for FFPE Samples

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

- Any new or previously untested FFPE tissue types
- For specific guidance on other sample preparations contact ACD Support at **support.acd@bio-techne.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-techne.com**.

Species	Tissue	ae Pathology Pretreatment Species		Tissue Type	Pathology	Pretreatment	
	Туре		Condition				Condition
Mouse/	Intestine	Normal	Standard	Human	Neck	Cancer	Standard
Rat	Intestine	Tumor	Standard		Liver	Cancer	Standard
	Embryo	Normal	Standard		Liver	Normal	Standard
	Brain	Normal	Standard		Heart	Normal	Standard
	Spleen	Normal	Standard		GI tract	Normal	Standard
	Eye/Retina	Normal	Extended		Kidney	Normal	Standard
	Liver	Normal	Standard		Skin	Normal	Standard
	Kidney	Normal	Standard		Lymphoma	Cancer	Standard
Human	Breast	Tumor	Standard		Thymus	Normal	Mild/Standard
	Colon	Tumor	Standard		Melanoma	Tumor	Standard
	Colon	Normal	Standard		Nevus	Benign	Standard
	Lung	Tumor	Standard		Placenta	Normal	Standard
	Lung	Normal	Standard		Skin (TMA*)	Normal	Standard
	Prostate	Tumor	Standard		Breast (TMA*)	Normal	Standard
	Prostate	Normal	Standard		Melanoma	Normal	Standard
	Lymph	Tumor	Standard		(TMA*)		
	node				Nevus (TMA)	Benign	Standard
	Lymph	Normal	Mild		Stomach (TMA)	Normal	Standard
	node				Stomach (TMA)	Tumor	Standard
	Tonsil	Normal	Mild/Standard		Cell pellets,	—	Mild
	Pancreas	Normal	Standard		fixed with 10%		
	Cervical	Cancer	Standard		NBF		
	Cervical	Normal	Standard		HeLa or 3T3	—	Mild
	Cervical	Abnormal	Standard		cells, fixed with		
	dysplasia				10%		
	Brain	Tumor	Standard		Formaldehyde		
	Brain	Normal	Standard		/PBS/ACD		
	Cancer	Standard	Head		Control		A (*1.1
					Xenograft	—	Milld
					tissue		

RNAscope 2.5 LS and LSx Reagent Kits - RED User Manual

Species	Tissue Type	Pathology	Pretreatment	Species	Tissue	Pathology	Pretreatment
Cyno	Spleen	Normal	Mild	Dog	Spleen	Normal	Mild
monkey	Lymph Node	Normal	Mild	Dog	Lymph	Mild	Mild
	Tonsil	Normal	Mild		Node	wind	Wind
	Thymus	Normal	Mild		Tonsil	N.A.	N.A.
	Retina	Normal	Mild		Thymus	Mild	Mild
	Prostate	Normal	Standard/Mild		Retina	Mild	Mild
	Gland				Prostate	Mild	Mild
	Epididymis	Normal	Mild/Standard		Gland		
	Testis	Normal	Mild/Standard		Epididy	Mild	Mild
	Ovary	Normal	Mild/Standard		mis		
	Duodenum	Normal	Mild/Standard		Testis	Mild/Stand	Mild/Standard
	Jejunum	Normal	Mild/Standard			ard	
	Colon	Normal	Standard		Ovary	Mild/Stand	Mild/Standard
	Adrenal	Normal	Mild/Standard			ard	
	Gland				Duoden	Normal	Mild
Rat	Spleen	Normal	Mild		um		
	Lymph Node	Normal	Mild		Jeiunum	Normal	Mild
	Tonsil	Normal	N.A.		Jejunum	ivorinai	Wind
	Thymus	Normal	Mild		Colon	Normal	Mild
	Retina	Normal	Mild		Adrenal	Normal	Standard/Mild
	Prostate	Normal	Standard/Mild		Gland		
	Gland						
	Epididymis	Normal	Standard				
	Testis	Normal	Standard				
	Ovary	Normal	Standard				
	Duodenum	Normal	Standard/Mild				
	Jejunum	Normal	Standard				
	Colon	Normal	Standard				
	Adrenal	Normal	N.A				
	Giana	<u> </u>	l				



Appendix B. Edit a Heat Treatment Protocol

The following example shows how to edit the Epitope Retrieval procedure from within the software.

Create a prestaining protocol

1. Open the Leica BOND software and click on the Protocol setup icon as shown.



2. Select **Prestaining** under the Protocol group menu and **Heat pretreatment** under the Protocol type menu to access the heat pretreatment protocols.

Slide setup	Protocol setup Reagent setup Silde history	Search Help	Log out ◆Ĵ		•	<u>feica</u>
3219600	Protocol setup			Copy Open	Delete	Report
	Protocol name +	Protocol type	Description	Modified by	Mod. date	Pref.
	*AccuCyte CTC HIER ER2 8 mins	Heat pretreatment	AccuCyte CTC HIER ER2 8 mins	Leica	9/1/2020	× 1
	*ACD HIER 15 min with ER2 (95)	Heat pretreatment	ACD RNAscope heat pretreatment	Leica	9/1/2020	~
	*E10 and H2 20 HIER	Heat pretreatment	10 min enzyme & 20 min heat retrieval with ER2 variant 1	Leica	9/1/2020	
	*E10 and H2 20 HIER2	Heat pretreatment	10 min enzyme & 20 min heat retrieval with ER2 variant 2	Leica	9/1/2020	
	"HIER 10 min with ER1	Heat pretreatment	10 min Heat Retrieval using ER1	Leica	9/1/2020	~
	"HIER 10 min with ER2	Heat pretreatment	10 min Heat Retrieval using ER2	Leica	9/1/2020	~
	*HIER 20 min with ER1	Heat pretreatment	20 min Heat Retrieval using ER1	Leica	9/1/2020	~
	*HIER 20 min with ER1 (95)	Heat pretreatment	20 min Heat Retrieval using ER1 at 95	Leica	9/1/2020	
	*HIER 20 min with ER2	Heat pretreatment	20 min Heat Retrieval using ER2	Leica	9/1/2020	~
	*HIER 25 min with ER1 (97)	Heat pretreatment	25 min Heat Retrieval using ER1	Leica	9/1/2020	~
	'HIER 30 min with ER1	Heat pretreatment	30 min Heat Retrieval using ER1	Leica	9/1/2020	~
	*HIER 30 min with ER2	Heat pretreatment	30 min Heat Retrieval using ER2	Leica	9/1/2020	~
	*HIER 40 min with ER1	Heat pretreatment	40 min Heat Retrieval using ER1	Leica	9/1/2020	~
	*HIER 40 min with ER1 (98)	Heat pretreatment	40 min Heat Retrieval using ER1 at 98	Leica	9/1/2020	
	*HIER 40 min with ER2	Heat pretreatment	40 min Heat Retrieval using ER2	Leica	9/1/2020	~
	*HIER 5 min with ER1	Heat pretreatment	5 min Heat Retrieval using ER1	Leica	9/1/2020	~
	*RNAscope 2.5 LSx Target Retrieval (88)	Heat pretreatment	RNAscope 2.5 LSx heat retrieval 88C	Leica	9/1/2020	~
	*RNAscope 2.5 LSx Target Retrieval (95)	Heat pretreatment	RNAscope 2.5 LSx heat pretreatment 95C	Leica	9/1/2020	~
	*ViewRNA HIER 10 min, ER1 (95)	Heat pretreatment	10 min heat pretreatment, ER1 @ 95C for Affymetrix ViewRNA eZ Assay	Leica	9/1/2020	~
	*ViewRNA HIER 10 min, ER2 (90)	Heat pretreatment	10 min heat pretreatment, ER2 @ 90C for Affymetrix ViewRNA eZ Assay	Leica	9/1/2020	~
	ACD HIER 10 min with ER2 (88)	Heat pretreatment	ACD RNAscope heat pretreatment 10 min at 88	bondpoweruser	10/18/2021	~
	Protocol group:	Protocol type:	Strining status: Protocol origin:	Preferred status:		-
	Prestaining	Heat pretreatment	All	All		•

3. Highlight a protocol, for example, *ACD HIER 15 min with ER2 (95). Select Copy.

Note: ER2 = Epitope Retrieval 2.

- 4. Rename the protocol as ACD HIER 15 min with ER2 (88).
- 5. Rename the Abbreviated name as ER2-88.
- 6. Rename the Description to ACD RNAscope heat pretreatment 88.
- 7. Highlight the third *BOND ER Solution 2 step (see above) and change temperature to 88°C.

		New proto	ocol proper	ties		
Name:	ACD HIER 15 min with ER2 (8	38)				
Abbreviated name:	ACD88					
Description:	ACD RNAscope heat pretreat	ment 88				
						Preferred
BOND RX					Import protocol	Protocol type: Heat pretreatment
Step N° Was	sh Reagent	Supplier	Ambient	Temperature	Inc. (min)	Dispense type
1	*Bond ER Solution 2	Leica Microsystems	~		0:00	150 µL
2	*Bond ER Solution 2	Leica Microsystems	~		0:00	150 µL
3	*Bond ER Solution 2	Leica Microsystems		88	5:00	Intermediate
4	*Bond ER Solution 2	Leica Microsystems	~		0:00	150 µL
Show wash ste	eps					
		Savo	Canco			
		Save	Cance			

RNAscope 2.5 LS and LSx Reagent Kits - RED User Manual

UM 322150/Rev D/ Effective Date 04/24/2025

- 8. For RNAscope, ER 2 temperature varies between **95°C** and **88°C** depending on the tissue type used. Please see **Appendix A** for a list of tissues.
- 9. Select **Save** to create a protocol for ER2 pretreatment at **88°C**.
- 10. If needed, repeat Steps 1–8 to create new heating protocols for different incubation times (for example, ACD 25minER2).



Appendix C. Edit an Enzyme Pretreatment Protocol

The following example shows how to edit the protease procedure from within the software.

- 1. Select **Enzyme Pretreatment** under the Protocol type menu (bottom left).
- 2. Highlight the *ACD 15min Protease protocol. Select Copy.

BOND - (bondpo	weruser) - Internet E	xplorer							-	ø×
Slide setup	Protocol setup	Reagent setup	Slide history	Search	Help	Log out				<u>feica</u>
3219600	Protoco	ol setup						Copy Open	Delete	Report
		Protocol	name 🔺	1	Protocol type	Description		Modified by	Mod. date	Pref.
	*ACD 15 r	min Protease		Er	nzyme pretreatment	ACD RNAscope enzyme pretreatment		Leica	9/1/2020	× 1
	*BaseSco	pe LSx Enzyme		Er	nzyme pretreatment	15 min Enzyme Pretreatment using BaseScope LSx Protease		Leica	5/5/2021	
	*E10 and	H2 20		Er	nzyme pretreatment	10 min enzyme & 20 min heat retrieval with ER2		Leica	9/1/2020	
	*Enzyme	1 (20)		Er	nzyme pretreatment	Enzyme 1 pretreatment (20 min) for ViewRNA eZ-LAssay		Leica	9/1/2020	
	*Enzyme	1 for 10 min		Er	nzyme pretreatment	10 min Enzyme Pretreatment using Enzyme 1		Leica	9/1/2020	~
	*Enzyme	1 for 15 min		Er	nzyme pretreatment	15 min Enzyme Pretreatment using Enzyme 1		Leica	9/1/2020	~
	*Enzyme	1 for 5 min		Er	nzyme pretreatment	5 min Enzyme Pretreatment using Enzyme 1		Leica	9/1/2020	~
	*Enzyme	2 (20)		Er	nzyme pretreatment	Enzyme 2 pretreatment (20 min) for ViewRNA eZ-LAssay		Leica	9/1/2020	
	*Enzyme	2 for 10 min		Er	nzyme pretreatment	10 min Enzyme Pretreatment using Enzyme 2		Leica	9/1/2020	~
	*Enzyme	2 for 15 min		Er	nzyme pretreatment	15 min Enzyme Pretreatment using Enzyme 2		Leica	9/1/2020	~
	*Enzyme	3 (20)		Er	nzyme pretreatment	Enzyme 3 pretreatment (20 min) for ViewRNA eZ-L Assay		Leica	9/1/2020	
	*Enzyme	3 for 10 min		Er	nzyme pretreatment	10 min Enzyme Pretreatment using Enzyme 3		Leica	9/1/2020	~
	*Enzyme	3 for 15 min		Er	nzyme pretreatment	15 min Enzyme Pretreatment using Enzyme 3		Leica	9/1/2020	~
	*Enzyme	5 for 25 min		Er	nzyme pretreatment	25 min Enzyme Pretreatment using Enzyme 5		Leica	9/1/2020	~
	*Protease	20 min and fix		Er	nzyme pretreatment	Protease 20 min and fix		Leica	9/1/2020	~
	"RNAscop	e 2.5 LSx Enzyme		Er	nzyme pretreatment	15 min Enzyme Pretreatment using RNAscope 2.5 LSx Proteas	50 50	Leica	9/1/2020	~
	*ViewRNA	Enzyme 1 (20)		Er	nzyme pretreatment	20 min enzyme pretreatment for Affymetrix ViewRNA eZ Assay		Leica	9/1/2020	~
	"ViewRNA	A Enzyme 2 (20)		Er	nzyme pretreatment	20 min enzyme pretreatment for Affymetrix ViewRNA eZ Assay		Leica	9/1/2020	~
	*ViewRNA	Enzyme 3 (20)		Er	nzyme pretreatment	20 min enzyme pretreatment for Affymetrix ViewRNA eZ Assay		Leica	9/1/2020	~
	2hr 54C p	retreat		Er	nzyme pretreatment	ACD RNAscope enzyme pretreatment 2hr 54C		bondpoweruser	2/21/2023	~
	30min 400	C pretreat		Er	nzyme pretreatment	ACD RNAscope enzyme pretreatment 30min 40c		bondpoweruser	7/31/2023	× .
	Protocol grou	p:		Protocol type:		Staining status:	Protocol origin:	Preferred status:		
	Prestaining		•	Enzyme pretre	eatment	•	All	All		•

- 3. Rename the protocol to ACD 25min Protease.
- 4. Rename the Abbreviated name to **25minPro**.
- 5. Rename the Description to ACD RNAscope 25min enzyme.
- 6. Highlight the second ***ACD Enzyme** step. Keep the temperature at **40°C** and set the enzyme incubation time to desired time (for example, 25min).

	ACD 25 min Protease						
Abbreviated name:	25mPro						
Description:	ACD RNAscope 25min enzy	yme pretreatment					
							 Prefe
BOND RX				Import p	rotocol	Protocol type: Enzyr	ne pretreatr
Step N° Wash	Reagent	Supplier	Ambient	Temperature	Inc. (min)	Dispense type	
2	*ACD Enzyme	Advanced Cell Diagnostics		40	0:00	150 µL	
3	*ACD Enzyme	Advanced Cell Diagnostics		40	25:00	50 µL	
7	*Open 0 Haz	User	~		10:00	150 µL	
Show wash steps	s						

- 7. Select Save.
- 8. If needed, repeat Steps 1–7 to create a new protease protocol for different sample types (for example, ACD 10min Protease or ACD 15min Protease at ambient temperature).
- 9. A similar workflow can be used to edit the *ACD PretreatPro and *RS 2.5 LSx PreTreatPro protocols if needed.
- * Indicates reagent is hard coded in software by Leica Biosystems.

×



Appendix D. Slide Setup for Additional Tissue Types

Alternatively prepared samples can be stained on the BOND RX using the following slide setup parameters.

Note: The following guidelines are not for use with the older mock probe workflow. If you need to use the older workflow and need assistance, please contact **support.acd@bio-techne.com** or your ACD FAS.

Fixed-frozen tissues

As described in **Chapter 4**, fixed-frozen tissues have only been optimized using the Protease Enzyme method. Start with the following conditions:

- ER2 retrieval of 5 minutes at 95°C
- 15 minutes of ACD's Protease treatment at 40°C
- 1. Confirm that you have a heat pretreatment available for the 5-minute treatment. See **Appendix B** to create this protocol if needed.
- 2. In Slide setup, select the following:
 - a. Staining: Choose the appropriate staining protocol for the chemistry and workflow you are using; LS 2.5 or LSx assay kits
 - b. Preparation: Select *----.
 - c. HIER: Choose ACD HIER 5 min with ER2 (95) or an edited RNAscope 2.5 LSx Target Retrieval protocol.
 - d. Enzyme: Select the appropriate protocol for the chemistry and workflow you are using; ***ACD 15min Protease** or **RNAscope 2.5 LSx Enzyme**.
 - e. Probe Application: Select ***RNAscope 2.5 LSx Probe Application**.
 - f. Denaturation: Choose *....
 - g. Hybridization: Choose the appropriate protocol for the chemistry and workflow you are using; **RNAscope 2.5 LSx Hybridization**.
 - h. Probe Removal: Select ***RNAscope 2.5 LSx Probe Removal**.

Note: When the run is complete, the BOND RX rinses the slides every 10 minutes which can impact the counterstain. Set up the instrument as late in the day as possible. Rinsing does not affect the RNAscope signal and counterstaining can be repeated offline in the morning if needed.

Start your run immediately after setting it up. Do not use a delayed start. This causes poor protease spreadability and negatively impacts results.

Fresh-frozen tissues

As described in Chapter 4, there are 2 options for processing Fresh-frozen tissues on the BOND RX: a protease-free option and a protease-based option that requires RNAscope LS Protease IV



Cat. No. 322140 which must be ordered separately. See below for both slide set ups. We recommend starting the protease-free workflow.

Protease-free workflow

With this method, a short retrieval step is applied followed by treatment with LS PretreatPro.

- 1. In Slide setup, please skip the Bake and Dewax steps
 - a. Staining: Select the appropriate protocol depending on the kit format you have: LS 2.5 or LSx assay kits.
 - b. Preparation: Choose *----.
 - c. HIER: Choose ACD HIER 5 min with ER2 (95) or edited RNAscope 2.5 LSx Target Retrieval (for 5mins) protocol depending on the kit format you have (See Appendix B to create this protocol if needed)
 - d. Enzyme: *ACD PretreatPro OR *RS 2.5 LSx_PreTPro depending on the kit format you have
 - e. Probe Application: Select *RNAscope 2.5 LSx Probe Application.
 - f. Denaturation: Select *....
 - g. Hybridization: Select *RNAscope 2.5 LSx Hybridization.
 - h. Probe Removal: Select *RNAscope 2.5 LSx Probe Removal.

Note: Start your run immediately after setting it up. Do not use a delayed start. This causes poor protease spreadability and negatively impacts results.

Protease workflow

With this method tissues do NOT need a target retrieval. Instead, permeabilize the tissue at ambient temperature with a stronger protease; RNAscope LS Protease IV Cat. No. 322140

- 1. In Slide setup, please skip the following steps: 1) Bake or Bake and Dewax 2) Heat retrieval. Choose the following instead:
 - a. Staining: Select the appropriate protocol depending on the kit format you have: LS 2.5 or LSx assay kits.
 - b. Preparation: Choose *----.
 - c. HIER: Choose *----.
 - d. Enzyme: Create and select a protocol with the following conditions **ACD 30min RT** with LS Protease IV (see Appendix C. Edit an Enzyme Pretreatment Protocol
 - e. Probe Application: Select *RNAscope 2.5 LSx Probe Application.
 - f. Denaturation: Select *....
 - g. Hybridization: Choose *RNAscope 2.5 LSx Hybridization.
 - h. Probe Removal: Select *RNAscope 2.5 LSx Probe Removal.

Note: Start your run immediately after setting it up. Do not use a delayed start. Otherwise, your protease will not distribute equally on the slide which can result in poor permeabilization.



Appendix E. 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), found at: www.access.gpo.gov/nara/cfr/waisidx_01/%2029cfr1910a_01.html
- 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, found at:

http://www.who.int/csr/resources/publications/biosafety/WHO_CDS_CSR_LYO_2004_11/en/

• Information about the Registration, Evaluation, Authorisation and Restriction of Chemicals (REACH) can be found at: http://echa.europa.eu/regulations/reach

Documentation and Support

Obtaining SDSs

Safety Data Sheets (SDSs) are available **in the documents download section of individual product pages at www.bio-techne.com**. For the SDSs of chemicals not distributed by Advanced Cell Diagnostics, contact the chemical manufacturer.

Obtaining support

For the latest support information, go to: https://www.bio-techne.com/resources or Contact us at https://www.bio-techne.com/support/contact-us.

Contact information

Advanced Cell Diagnostics, Inc.

7707 Gateway Blvd

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: order.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 at **https://www.bio-techne.com/terms-and-conditions**. If you have any questions, please contact Advanced Cell Diagnostics at **https://www.bio-techne.com/support/contact-us**

