

# miRNAscope<sup>™</sup> HD Reagent Kit (RED)

### Introduction

This quick guide is intended for advanced users who are familiar with the procedures in the *miRNAscope™ HD (RED) Assay User Manual* (Document No. UM 324510). Refer to the user manual for safety guidelines. For every chemical, read the Safety Data Sheet (SDS) and follow handling instructions. Wear appropriate protective eyewear, clothing, and gloves. For the latest services and support information, go to: https://acdbio.com/technical-support/support-overview.

Part 1: Prepare and Pretreat Samples

	Workflow Steps			
PREPARE SLIDES ~1.5 HOURS	Bake Slides  1. Bake slides in a dry oven for 1 HR at 60°C.			
Bake Slides	OPTIONAL STOPPING POINT (1). Use sectioned tissue within 1 week. Store sections with desiccants at room temperature (RT).			
Deparaffinize FFPE Sections	Deparaffinize FFPE Sections  1. In a fume hood:  • Fill two Tissue-Tek® Clearing Agent dishes with ~200 mL fresh xylene.  • Fill two Tissue-Tek® Staining dishes with ~200 mL fresh 100% alcohol.  2. Place slides in a Tissue-Tek® Slide Rack in xylene 2 x 5 MIN.  3. Incubate slides in 100% alcohol 2 x 2 MIN.  4. Remove slides from rack. Dry slides for 5 MIN at 60°C or until completely dry.			
POST-FIX SAMPLES 2 HOURS or OVERNIGHT	<ol> <li>In a fume hood, do one of the following:         <ul> <li>Incubate slides in fresh 10% NBF OVERNIGHT (~16 HRS) at RT or</li> <li>Incubate slides in fresh 12% formaldehyde for 2 HRS at RT</li> </ul> </li> <li>Remove slides from fix and wash them for 2 MIN in distilled water.</li> <li>Dry slides for 5 MIN at 60°C or until completely dry.</li> </ol>			
PRETREAT SAMPLES ~1-2 HOURS	Prepare Oven and Reagents (30 MIN at 40°C)  4. Set HybEZ <sup>™</sup> oven to 40°C and warm HybEZ <sup>™</sup> Humidity Control Tray containing wet humidifying paper for 30 MIN before use.			
Prepare Oven and Reagents ↓	5. Prepare 200 mL fresh 1X Target Retrieval Reagent by adding 180 mL distilled water to 20 mL of 10X Target Retrieval Reagents. Mix well.			
Apply Hydrogen Peroxide ↓	6. Turn on the steamer, and set the heating time to <b>95 MIN</b> . Allow the temperature to rise to at least <b>99°C</b> .			
Apply Target Retrieval  ↓ Create Barrier  ↓ Apply Protease III	<ol> <li>Apply Hydrogen Peroxide (10 MIN at RT)</li> <li>Add ~5-8 drops of hydrogen peroxide to each section for 10 MIN at RT.</li> <li>Remove the hydrogen peroxide, and place the slides in a slide rack.</li> <li>Wash the slides by moving the rack up and down 3-5 times in a staining dish filled with distilled water. Repeat with fresh water.</li> <li>Apply Target Retrieval (15-30 MIN at boiling point)</li> </ol>			
	<ol> <li>With a pair of forceps slowly submerge the slide rack into boiling 1X Target retrieval solution for 15 MIN. Refer to Appendix A of the user manual (Doc. No. UM 324510) to determine the appropriate conditions for your tissue type.</li> <li>Wash the slides by moving the rack up and down 3–5 times in a staining dish filled with distilled water. Repeat with fresh water.</li> </ol>			



3. Wash slides in 100% alcohol, then air dry.

#### Create Barrier

Once the slides are dry, draw 2-4 times around the tissue using the Immedge<sup>™</sup> hydrophobic barrier pen. Dry completely ≥ 5 MIN at RT.

OPTIONAL STOPPING POINT (2). Dry slides overnight for use the following day or proceed directly to the next section.

### Apply Protease III (15-30 MIN at 40°C)

- Place the slides in the ACD EZ-Batch<sup>™</sup> Slide Holder, and then add ~4 drops of protease to cover each section.
- 2. Place the slide holder into the pre-warmed humidity control tray. Seal the tray.
  - For HeLa cell pellets, incubate slides for 15 MIN at 40°C in the HybEZ™ oven.
  - For cell pellets, incubate slides for 15 MIN at RT.
  - For tissues, incubate slides for 30 MIN at 40°C in the HybEZ<sup>™</sup> oven.
- 3. Place the slide holder with the slides into the ACD EZ-Batch™ Wash Tray filled with fresh distilled water. Wash the slides with slight agitation.

# Part 2: Run the Assay

Workflow Steps					
PREPARE THE MATERIALS ~10-30 MIN	<ol> <li>Warm 50X Wash Buffer for 10–20 MIN before preparing 1X Wash Buffer solution.</li> <li>Prepare 5 L of 1X Wash Buffer by adding 4.90 L distilled water and 100 mL of 50X Wash Buffer to a large carboy. Mix well.</li> <li>Prepare 50% Gill's Hematoxylin-I and 0.02% Ammonium Hydroxide.</li> <li>Equilibrate reagents:         <ul> <li>Place Amps 1–6 at RT.</li> <li>Equilibrate probes for 15 MIN at 40°C.</li> </ul> </li> </ol>				
RUN THE ASSAY ~8 HOURS	Hybridize Probe (2 HRS at 40°C)  1. Remove excess liquid from slides, place in the ACD EZ-Batch™ Slide Holder, and add ~4 drops of probe to each section.				
Hybridize Probe ↓ Hybridize Amp 1 ↓ Hybridize Amp 2	<ol> <li>Insert sealed tray containing the slide holder into the HybEZ<sup>™</sup> Oven for 2 HRS at 40°C.</li> <li>Wash slides in 1X Wash Buffer for 2 MIN at RT. Repeat with fresh 1X Wash Buffer.     Hybridize Amp 1 (30 MIN at 40°C)</li> <li>Remove excess liquid from slides, place in the ACD EZ-Batch<sup>™</sup> Slide Holder, and add ~4 drops Amp 1 to each section.</li> </ol>				
↓ Hybridize Amp 3	<ol> <li>Insert sealed tray containing the slide holder into the HybEZ<sup>™</sup> Oven for 30 MIN at 40°C. Remove slide rack.</li> </ol>				
↓ Hybridize Amp 4 ↓	3. Wash slides in 1X Wash Buffer for <b>2 MIN</b> at <b>RT</b> . Repeat with fresh 1X Wash Buffer. Hybridize Amp 2 (15 MIN at 40°C)				
Hybridize Amp 5 ↓	<ol> <li>Remove excess liquid from slides, place in the ACD EZ-Batch™ Slide Holder, and add ~4 drops Amp 2 to each section.</li> </ol>				
Hybridize Amp 6 ↓	<ol> <li>Insert sealed tray containing the slide holder into the HybEZ<sup>™</sup> Oven for 15 MIN at 40°C. Remove slide rack.</li> </ol>				
Detect the Red Signal  3. Wash slides in 1X Wash Buffer for 2 MIN at RT. Repeat with fresh but Hybridize Amp 3 (30 MIN at 40°C)					
Counterstain the Slides	<ol> <li>Remove excess liquid from slides, place in the ACD EZ-Batch™ Slide Holder, and add ~4 drops Amp 3 to each section.</li> </ol>				



### Workflow Steps

## Evaluate the Results

- Insert sealed tray containing the slide holder into the HybEZ<sup>™</sup> Oven for 30 MIN at 40°C. Remove slide rack.
- 3. Wash slides in 1X Wash Buffer for **2 MIN** at **RT**. Repeat with fresh buffer.

#### Hybridize Amp 4 (15 MIN at 40°C)

- Remove excess liquid from slides, place in the ACD EZ-Batch<sup>™</sup> Slide Holder, and add ~4 drops Amp 4 to each section.
- Insert sealed tray containing the slide holder into the HybEZ<sup>™</sup> Oven for 15 MIN at 40°C.
- 3. Wash slides in 1X Wash Buffer for 2 MIN at RT. Repeat with fresh buffer.

**IMPORTANT!** Do not insert tray into HybEZ<sup>TM</sup> Oven for the rest of the procedure.

#### Hybridize Amp 5 (30 MIN at RT)

- 1. Remove excess liquid from slides, place in the ACD EZ-Batch™ Slide Holder, and add ~4 drops Amp 5 to each section.
- 2. Incubate the sealed tray containing the slide holder for 30 MIN at RT.
- 3. Wash slides in 1X Wash Buffer for **2 MIN** at **RT**. Repeat with fresh buffer.

#### Hybridize Amp 6 (15 MIN at RT)

- Remove excess liquid from slides, place in the ACD EZ-Batch<sup>™</sup> Slide Holder, and add ~4 drops Amp 6 to each section.
- 2. Incubate the sealed tray containing the slide holder for for 15 MIN at RT.
- 3. Wash slides in 1X Wash Buffer for **2 MIN** at **RT**. Repeat with fresh buffer.

#### Detect the Red Signal (10 MIN at RT)

- 1. Briefly spin Red-B and mix 1 volume of Red -B to 60 volumes of Red -A (use within **3–5 MIN**).
- 2. Remove excess liquid from slides, place in the ACD EZ-Batch™ Slide Holder, and pipette ~120 µL of Red solution onto each tissue section.
- Incubate sealed tray containing the slide holder for 10 MIN at RT.
- 4. Place slides in a Tissue-Tek® Slide Rack, and wash the slides by moving the rack up and down 3–5 times in a staining dish filled with distilled water.

#### Counterstain the Slides

- 1. Place slides in 50% Hematoxylin for 2 MIN at RT.
- Wash the slides by moving the rack up and down 3-5 times in distilled water.Repeat with fresh distilled water.
- Place slides in 0.02% Ammonium Hydroxide for 10 SEC.
- 4. Wash the slides by moving the rack up and down 3-5 times in distilled water.

#### Mount the Slides

- 1. Dry slides in a 60°C dry oven for 15-30 MIN.
- 2. Dip the slides into fresh pure xylene until the hydrophobic barrier disappears and immediately place 1–2 drops of EcoMount on the slides before the xylene dries. Place a coverslip over the section.
- Before viewing under the microscope, dry slides for at least 30 MIN or until the xylene completely evaporates.

#### Evaluate the Results

 Examine tissue sections under a standard bright field microscope at 20–40X magnification.



# **Troubleshooting**

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

For Research Use Only. Not for Diagnostic Use.

NOTICE TO PURCHASER: 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.

© 2020 Advanced Cell Diagnostics. All rights reserved. RNAscope<sup>™</sup>, miRNAscope<sup>™</sup> and HybEZ<sup>™</sup> are trademarks of Advanced Cell Diagnostics, Inc. All other trademarks belong to their respective owners.

