

# RNAscope® 2.5 LSx Reagent Kit (BROWN) Quick Guide

### For FFPE Tissues

#### Introduction

This quick guide is intended for advanced users who are familiar with the procedures in the RNAscope® 2.5 LSx Reagent Kit-BROWN User Manual (Document No. 322700-USM). 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: Reagents and Software

Workflow Steps	
MATERIALS FROM ACD	<ul> <li>The RNAscope® 2.5 LSx Reagent Kit – BROWN (Cat. No. 322700) consisting of two trays of reagents in labeled Leica BOND compatible containers</li> <li>The first tray has two tray barcodes and a label on the handle, and contains the LS Rinse, DAB chromogen, and counterstaining reagents</li> <li>The second tray contains the pretreatment and signal amplification reagents</li> <li>RNAscope® 2.5 LS control probes and custom probes</li> </ul>
MATERIALS FROM LEICA	<ul> <li>Leica Biosystems' BOND RX System — automated slide stainer</li> <li>BOND Open Containers 30 mL</li> <li>BOND Universal Covertiles 100 pack</li> <li>BOND bulk buffers, including ER buffers, wash buffer, and dewax buffer</li> </ul>
VERIFY THE SOFTWARE VERSION	Update the BDZ/BXD software to database version 15 before running the RNAscope® 2.5 LSx Brown Assay.      Note: Contact your ACD FAS for help with the 5.2 software database upgrade. To upgrade 4.0 software, contact a Leica field service scientist.      Note: The LSx assay does not need a workaround protocol or mock probe. Each probe is used as a marker.

Part 2: Set up RNAscope® 2.5 LSx Assay

Workflow Steps	
REGISTER THE LSx REAGENTS ~5-10 MIN	1. Register the reagents on the first tray by scanning the two tray barcodes sequentially with the Barcode Scanner. Select <b>OK</b> .
	2. Register each component separately on the second tray by scanning the barcode on each reagent container with the Barcode Scanner. Select <b>OK</b> .
REGISTER PROBES ~10-20 MIN	Register each probe as a marker for the ISH procedure:  1. Go to the <b>Reagent Setup</b> screen, and select <b>Add</b> to enter the probe name, type, and other information.

	Workflow Steps	
	<ol> <li>For software version 4.0, select Probe in the Type drop down menu and RNA for Probe Type. For software version 5.2, select Probe RNA in the Type drop down menu.</li> <li>Select the default protocols you want to use for the probe.</li> <li>Select Hazardous, then Save.</li> </ol>	
SET UP A STUDY ~10-30 MIN	<ol> <li>Go to the Slide setup screen, select Add Study, and enter information.</li> <li>For each slide and probe used in the assay:         <ul> <li>Go to the Add slide screen, select ISH and the probe of interest under the Marker drop down menu.</li> <li>In the Protocols tab, choose the following LSx protocols. Do not use use protocols from other LS assays.</li></ul></li></ol>	
START THE ASSAY ~10-30 MIN	<ol> <li>Prepare the instrument, including filling the bulk reagents and emptying the waste container.</li> <li>Load the two registered 2.5 LSx reagent trays onto the instrument.</li> <li>Empty each target probe into a new 30mL Open container. Register each probe container by scanning the barcode in the front and selecting the corresponding name of the probe. Load probes onto the instrument.</li> <li>Go to the Slide setup screen.</li> <li>Attach the barcodes to the slides, add the covertiles, and place the slides on the slide tray with the label sides facing up.</li> <li>Place the tray in the Leica BOND RX, and press the button to load the tray onto the machine.</li> <li>After scanning the slides, select the PLAY (triangular) button on the screen located under the start tray. You may also right-click on the scanned label images, and select Delayed Start to start the run at a future time.</li> </ol>	

## **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: PLEASE REFER TO THE RNASCOPE® 2.5 LS REAGENT KIT- BROWN USER MANUAL FOR LIMITED USE LABEL LICENSE OR DISCLAIMER INFORMATION. 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.

© 2018 Advanced Cell Diagnostics. All rights reserved. RNAscope® is a registered trademark of Advanced Cell Diagnostics, Inc. Leica BOND RX is a registered trademark of Leica Biosystems. All other trademarks belong to their respective owners.

#### Headquarter:

