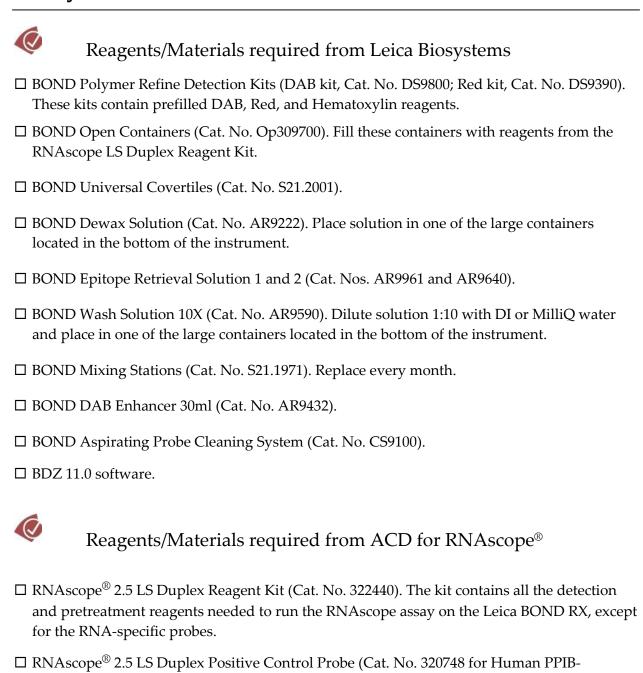


Checklist for the RNAscope® Automated Assay on Leica Biosystems' BOND RX System RNAscope®2.5 LS Duplex

Assay Doc. No. CHK 47-001



C1/Polr2A-C2; Cat No. 320768 for Mouse PPIB-C1/POLR2A-C2). Probes target common

□ RNAscope[®] 2.5 LS Duplex Negative Control Probe (Cat. No. 312758). The probe targets the bacterial *dapB* gene in two channels to control for background noise, and to help interpret

housekeeping genes to help qualify your samples and interpret your results.

assay results (dapB C1/dapB-C2).



□ RNAscope [®] 2.5 LS Target Probes target specific RNAs of interest in Channel C1 and Channel C2.
□ RNAscope [®] Control Slides (Cat. No. 310045 for Human Cell Pellet; Cat. No. 310023 for Mouse 3T3 Cell Pellet).
$\hfill\square$ Software protocols created for each probe (workaround solution).
Reagents/Materials provided by Researcher
☐ Superfrost® Plus slides (Fisher Scientific): Superfrost Plus slides are required. Other slide types may result in tissue detachment.
☐ Fresh reagents (e.g. ethanol and xylene).
☐ Fresh 10% NBF (neutral-buffered formalin).
☐ Tissue-Tek® slide rack and staining dishes.
☐ EcoMount or PERTEX®.
☐ Hotplate, drying oven, water bath, thermometer, and microscope.
☐ Paraffin wax, microtome, and fume hood.
Note: Always refer to the latest version of the user manuals for the complete set of materials required to run the RNAscope assay, available at http://www.acdbio.com/technical-support/user-manuals.

© 2020. Advanced Cell Diagnostics. All rights reserved. RNAscope $^{\circledcirc}$ is a trademark of Advanced Cell Diagnostics, Inc. Leica BOND RX is a registered trademark of Leica Biosystems. All other trademarks belong to their respective