

BaseScope[™] LS Automated Assay on Leica Biosystems' BOND RX System Doc. No. CHK 47-013



Reagents/Materials required from Leica Biosystems

- □ BOND Polymer Red Refine Detection Kit (Cat. No. DS9390)
 This kit contains prefilled Red chromogen and Hematoxylin reagents
 □ BOND 30ml Open Containers (Cat. No. Op309700) -minimum 20 required
 Fill these containers with reagents from the BaseScope™ LS Reagent Kit
 □ BOND Universal Covertiles (Cat. No. S21.2001)
 □ BOND Dewax Solution (Cat. No. AR9222)
 □ BOND Epitope Retrieval Solution 2 (Cat. No. AR9640)
 □ BOND Epitope Retrieval Solution 1 (Cat. No. AR9961)
 □ BOND Wash Solution 10X (Cat. No. AR9590)
 Dilute the solution 1:10 and place in Bond Wash bulk container located in the bottom of the instrument.
 □ BOND Aspirating Probe Cleaning System (Cat. No. CS9100).
 □ BOND Mixing Stations. (Cat. No. S21.1971). Replace every month.

Reagents/Materials required from ACD for BaseScope

☐ BaseScope[™] LS Reagent Kit (Cat. No. 323600)

☐ BDZ/BXD Software version 15.

- Contains detection and pretreatment reagents that are needed to run the assay on the Leica BOND RX, except for probes.
- ☐ BaseScope[™] LS Target Probes targeting specific RNAs of interest
- ☐ BaseScope[™] LS Positive Control Probe targeting common housekeeping genes.
 - For 1ZZ target probes, choose 1ZZ control probes. For 2ZZ and 3ZZ target probes, choose 3ZZ control probes.



- Human: BaseScope™ LS Probe BA-Hs-PPIB-1ZZ (Cat No. 710178); PPIB-3ZZ (Cat. No. 701038)
- Mouse: BaseScope™ LS Probe BA-Mm-Ppib-1 ZZ (Cat No. 712358); Ppib-3ZZ (Cat. No. (701078).
- □ BaseScope[™] LS Negative Control Probe. Targets the bacterial *dapB* gene to control for background noise, and to help interpret assay results.
 - BaseScope[™] LS Negative Control Probe DapB-1ZZ (Cat No. 701028)
 - BaseScope[™] LS Negative Control Probe DapB-3ZZ (Cat No. 701018)
- ☐ Probe Packs, available for Human and Mouse:

Cat# 323631 BaseScope™ LS Control Probe Pack-Human

- Hs-PPIB-1ZZ (Cat No. 710178)
- Hs-PPIB-3ZZ (Cat. No. 701038)
- DapB-1ZZ (Cat No. 701028)
- DapB-3ZZ (Cat No. 701018)

Cat# 323632 BaseScope™ LS Control Probe Pack-Mouse

- Mm- Ppib-1ZZ (Cat No. 712358)
- Mm-Ppib-3ZZ (Cat. No. 701078)
- DapB-1ZZ (Cat No. 701028)
- DapB-3ZZ (Cat No. 701018)
- ☐ RNAscope® Control Slides. Choose the species needed:
 - Human Cell Pellet (Cat. No. 310045)
 - Mouse 3T3 Cell Pellet (Cat. No. 310023)

Ø

Reagents/Materials provided by Researcher

\square Superfrost TM 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 (required)
☐ Paraffin wax, microtome, fume hood, drying oven, and microscope.

Note: Always refer to the latest version of the user manuals for the complete set of materials required to run the BaseScope™ LS Assay available at https://acdbio.com/technical-support/user-manuals.

© 2021 Advanced Cell Diagnostics. All rights reserved. RNAscope and BaseScope are trademarks and registered trademarks of Advanced Cell Diagnostics, Inc. Leica BOND RX is a registered trademark of Leica Biosystems. All other trademarks belong to their respective owners.