
Checklist for Manual BaseScope[™] Detection Reagent Kit v2 – RED

Doc. No. CHK 47-015



Reagents/Materials required from ACD

- BaseScope[™] Positive Control Probes targeting common housekeeping genes to help qualify RNA quality of the samples and interpret assay results.
- BaseScope[™] Negative Control Probe targeting the bacterial gene *dapB* to control for background noise and to help interpret assay results.
- BaseScope[™] Reagent Kit v2 RED Cat No. (323900) consisting of a pretreatment kit, a detection kit and wash buffer.

ALTERNATIVE:

- BaseScope[™] Introductory Pack for Reagent Kit v2 Red. Designed for first-time users of the RNAscope assay (Cat. No. 323970 Human or Cat. No. 323971 Mouse), and providing a complete set of reagents including controls probes, slides, hydrophobic barrier pen, and the BaseScope[™] Reagent Kit v2 – RED.
- BaseScope[™] Target Probes C1 Channel probe.
- HybEZ[™] Hybridization System with ACD EZ-Batch Slide System (321710 or, 321720) is required for RNAscope hybridization to maintain optimum humidity and temperature at 40°C.
- HybEZ[™] Hybridization System with ACD EZ-Batch Slide System I or II (Cat. Nos. 321461 (110VAC) and 321462 (220VAC); Cat. Nos. 321711 (110VAC) and 321721 (220VAC)). The system maintains optimum humidity and temperature at 40°C during RNAscope hybridization.
- RNAscope[®] Control Slides (Human Cell Pellet, Cat. No. 310045; Mouse 3T3, Cell Pellet Cat. No. 310023).
- ImmEdge Hydrophobic Barrier Pen (Cat. No. 310018) is the only pen that will maintain a hydrophobic barrier throughout the RNAscope[®] procedure.
- Vectamount* Mounting media (Cat. No. 321584) or Ecomount (Cat. No. 320409) can be used with the BaseScope[™] HD Red assay.



Reagents/Materials provided by Researcher

- Superfrost[™] Plus slides (Fisher Scientific): Superfrost Plus slides are required. Other slide types may result in tissue detachment.
- Coverlips (Fisher Scientific; Cat. No. 12-545F)
- Gill's Hematoxylin - I (American Master Tech Scientific; Cat. No. HXGHE1LT)
- Fresh reagents (e.g. ethanol and xylene).
- Fresh 10% NBF (neutral-buffered formalin).
- Ammonium Hydroxide, 28–30%.
- Hotplate, drying oven, water bath, thermometer, microscope
- Paraffin wax, microtome, fume hood.

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