
RNAscope™ VS Duplex Assay Automated Assay on the Ventana™ DISCOVERY™ ULTRA System

Doc. No. CHK 47-011



Reagents/Materials required from Ventana/Roche Diagnostics

- mRNA Duplex Amplification Kit, mRNA Sample Prep Kit, mRNA Detection Kits, and mRNA Link Kits contain prefilled mRNA Detection Reagents and the user-fillable empty dispensers for all ACD reagents except probes.
 - mRNA Sample Prep Kit (Cat. No. 760-248; Ordering Code 08127166001)
 - mRNA Duplex Amp Kit (Cat. No. 760-249; Ordering Code 08127174001)
 - mRNA RED Detection Kit (Cat. No. 760-234; Ordering Code 7099037001)
 - mRNA DAB Detection Kit (Cat. No. 760-224; Ordering Code 06614353001) OR mRNA TEAL Detection Kit (Cat. No. 760-256; Ordering Code 08352941001)
 - mRNA Link (Cat. No. 760-6014; Ordering Code 08127115001)

- DISCOVERY™ Wash Buffer (Cat. No. 950-510; Ordering Code 07311079001).
- ULTRA LCS (Predilute) (Cat No. 650-210; Ordering Code 5424534001).
- Reaction Buffer (10X) (Cat No. 950-300; Ordering Code 05353955001).
- DISCOVERY CC1 (Cat No. 950-500; Ordering Code 6414575001).
- 10X SSC Solution (Cat. No. 950-110; Ordering Code 05353947001). Required to run mRNA procedures on the DISCOVERY™ ULTRA.
- User-fillable probe dispensers (Cat. No. 960-761 to 960-785) for negative control probe, positive control probe, and target probe. Additional dispensers can be ordered for more target probes. See the Ventana™ website for part numbers.
- mRNA Universal Protocol (which can be installed by Roche or ACD FAS)
- Counterstaining reagents.
 - Hematoxylin II (Cat. No. 760-2208; Ordering Code 05277965001)
 - Bluing (Cat. No. 760-2037; Ordering Code 05266769001)

Notes:

- If ACD's or one's own hematoxylin and bluing reagents is used, Counterstain 1 (Cat. No. 771-741) and Counterstain 2 (Cat. No. 771-742) user-fillable dispensers from Ventana™ will need to be ordered.
- The Ventana™ hematoxylin and bluing reagent can be used with the RNAscope® VS Universal Assay.



Reagents/Materials required from ACD for RNAscope

- RNAscope™ VS Duplex Reagent Kit (Cat. No. 323300). Contains the following:
 - RNAscope™ VS Duplex Detection Reagents (Cat. No. 323310)
 - RNAscope™ VS Universal Sample Prep Reagents v2 (Cat. No. 323740)
 - RNAscope™ VS Accessory Kit (Cat. No. 320630)
- RNAscope™ 2.5 VS Duplex Positive Control Probe: PPIB-C1 and Polr2A-C2 (Cat No. 320749 for Human; Cat No. 320769 for Mouse).
- RNAscope™ 2.5 VS Duplex Negative Control Probe: DapB-C1 and DapB-C2 (Cat. No 320759). Targets the bacterial *dapB* gene to control for background noise, and to help interpret assay results.
- RNAscope™ 2.5 VS Target Probes targeting specific RNAs of interest.
- RNAscope™ Control Slides (Cat. No. 310045 for Human Cell Pellet; Cat. No. 310023 for Mouse 3T3 Cell Pellet).



Reagents/Materials provided by Researcher

- Superfrost™ Plus slides (Fisher Scientific): Superfrost Plus slides are required. Other slide types may result in tissue detachment.
- EcoMount is required for use with the RNAscope® VS Duplex Assay.
- Fresh reagents (e.g. ethanol and xylene)
- Dawn detergent
- Fresh 10% NBF (neutral-buffered formalin)
- Tissue-Tek® slide rack and staining dishes
- 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 <https://acdbio.com/technical-support/user-manuals>.