

## Manual assay: RNAscope® Multiplex Fluorescent Assay v2

Doc. No. CHK 47-007



## Reagents/Materials required from ACD

<ul> <li>RNAscope® Target Probes in Channel C1, Channel C2, and Channel C3 (C4 Channel probes are required to perform the 4-plex assay).</li> <li>Species-specific RNAscope® 3-plex Positive Control Probes- POLR2A (Channel C1), PPIB (Channel C2), UBC (Channel C3: Human Cat.No.320861; Mouse Cat.No.320881; Rat CatNo.320891). Probes target common housekeeping genes to help qualify samples and interpret results.</li> <li>RNAscope® 3-plex Negative Control Probe (Cat. No. 320871). The probe targets the bacterial dapB gene to control for background noise, and to help interpret assay results.</li> <li>RNAscope® Multiplex Fluorescent Reagent Kit v2 (Cat. No. 323100). The kit provides enough reagents to stain ~20 tissue sections, each with an area of approximately 20 mm x 20 mm (0.75in x 0.75in). The reagent kit contains pretreatment kit, detection kit, TSA buffer, and wash buffer.</li> <li>HybEZ™ Hybridization System with ACD EZ-Batch Slide System I (Cat. Nos. 321461, 110VAC and 321462, 220VAC) or II (Cat. Nos. 321711, 110VAC and 321721, 220VAC). The system maintains optimum humidity and temperature at 40°C during RNAscope® hybridization.</li> <li>RNAscope® Control Slides (Cat. No. 310045 for Human Cell Pellet; Cat. No. 310023 for Mouse 3T3 Cell Pellet).</li> <li>ImmEdge™ Hydrophobic Barrier Pen (Cat. No. 310018) is the only pen that will maintain a hydrophobic barrier throughout the RNAscope® procedure. Do not use any other pen as the barrier may not stay intact throughout the workflow).</li> </ul>	
<ul> <li>C2), UBC (Channel C3: Human Cat.No.320861; Mouse Cat.No.320881; Rat CatNo.320891). Probes target common housekeeping genes to help qualify samples and interpret results.</li> <li>□ RNAscope® 3-plex Negative Control Probe (Cat. No. 320871). The probe targets the bacterial dapB gene to control for background noise, and to help interpret assay results.</li> <li>□ RNAscope® Multiplex Fluorescent Reagent Kit v2 (Cat. No. 323100). The kit provides enough reagents to stain ~20 tissue sections, each with an area of approximately 20 mm x 20 mm (0.75in x 0.75in). The reagent kit contains pretreatment kit, detection kit, TSA buffer, and wash buffer.</li> <li>□ HybEZ™ Hybridization System with ACD EZ-Batch Slide System I (Cat. Nos. 321461, 110VAC and 321462, 220VAC) or II (Cat. Nos. 321711, 110VAC and 321721, 220VAC). The system maintains optimum humidity and temperature at 40°C during RNAscope® hybridization.</li> <li>□ RNAscope® Control Slides (Cat. No. 310045 for Human Cell Pellet; Cat. No. 310023 for Mouse 3T3 Cell Pellet).</li> <li>□ ImmEdge™ Hydrophobic Barrier Pen (Cat. No. 310018) is the only pen that will maintain a hydrophobic barrier throughout the RNAscope® procedure. Do not use any other pen as the</li> </ul>	· · · · · · · · · · · · · · · · · · ·
<ul> <li>dapB gene to control for background noise, and to help interpret assay results.</li> <li>RNAscope® Multiplex Fluorescent Reagent Kit v2 (Cat. No. 323100). The kit provides enough reagents to stain ~20 tissue sections, each with an area of approximately 20 mm x 20 mm (0.75in x 0.75in). The reagent kit contains pretreatment kit, detection kit, TSA buffer, and wash buffer.</li> <li>HybEZ™ Hybridization System with ACD EZ-Batch Slide System I (Cat. Nos. 321461, 110VAC and 321462, 220VAC) or II (Cat. Nos. 321711, 110VAC and 321721, 220VAC). The system maintains optimum humidity and temperature at 40°C during RNAscope® hybridization.</li> <li>RNAscope® Control Slides (Cat. No. 310045 for Human Cell Pellet; Cat. No. 310023 for Mouse 3T3 Cell Pellet).</li> <li>ImmEdge™ Hydrophobic Barrier Pen (Cat. No. 310018) is the only pen that will maintain a hydrophobic barrier throughout the RNAscope® procedure. Do not use any other pen as the</li> </ul>	C2), UBC (Channel C3: Human Cat.No.320861; Mouse Cat.No.320881; Rat CatNo.320891).
reagents to stain ~20 tissue sections, each with an area of approximately 20 mm x 20 mm (0.75in x 0.75in). The reagent kit contains pretreatment kit, detection kit, TSA buffer, and wash buffer.  □ HybEZ™ Hybridization System with ACD EZ-Batch Slide System I (Cat. Nos. 321461, 110VAC and 321462, 220VAC) or II (Cat. Nos. 321711, 110VAC and 321721, 220VAC). The system maintains optimum humidity and temperature at 40°C during RNAscope® hybridization.  □ RNAscope® Control Slides (Cat. No. 310045 for Human Cell Pellet; Cat. No. 310023 for Mouse 3T3 Cell Pellet).  □ ImmEdge™ Hydrophobic Barrier Pen (Cat. No. 310018) is the only pen that will maintain a hydrophobic barrier throughout the RNAscope® procedure. Do not use any other pen as the	
<ul> <li>and 321462, 220VAC) or II (Cat. Nos. 321711, 110VAC and 321721, 220VAC). The system maintains optimum humidity and temperature at 40°C during RNAscope® hybridization.</li> <li>□ RNAscope® Control Slides (Cat. No. 310045 for Human Cell Pellet; Cat. No. 310023 for Mouse 3T3 Cell Pellet).</li> <li>□ ImmEdge™ Hydrophobic Barrier Pen (Cat. No. 310018) is the only pen that will maintain a hydrophobic barrier throughout the RNAscope® procedure. Do not use any other pen as the</li> </ul>	reagents to stain ~20 tissue sections, each with an area of approximately 20 mm x 20 mm (0.75in x 0.75in). The reagent kit contains pretreatment kit, detection kit, TSA buffer, and wash
3T3 Cell Pellet).  □ ImmEdge <sup>™</sup> Hydrophobic Barrier Pen (Cat. No. 310018) is the only pen that will maintain a hydrophobic barrier throughout the RNAscope® procedure. Do not use any other pen as the	and 321462, 220VAC) or II (Cat. Nos. 321711, 110VAC and 321721, 220VAC). The system
hydrophobic barrier throughout the RNAscope® procedure. Do not use any other pen as the	·
	hydrophobic barrier throughout the RNAscope® procedure. Do not use any other pen as the



## Additional Reagents/Materials required from ACD for the RNAscope 4-plex assay

RNAscope® 4-Plex Ancillary Kit for Multiplex Fluorescent Kit v2 (Cat. No. 323120). The kit is
designed to work with the RNAscope® Multiplex Fluorescent Reagent Kit v2 and C1, C2, C3, and
C4 Channel Target Probes for 4-plex fluorescent in situ hybridization. It contains C4 HRP and HRP
blocker.

□ Species specific RNAscope® 4-plex Positive Control Probes for RNAscope Multiplex Fluorescent Assay- Polr2a (Channel C1), PPIB (Channel C2), UBC (Channel C3), and HPRT1 (Channel C4). UBC has the highest relative expression level; PPIB is considered to be a moderate-high expressor, while HPRT1 and POLR2A are moderate-to low expressor targets (Human Cat.No.321801; Mouse Cat.No.321818; Rat CatNo.321821).



□ RNAscope® 4-plex Multiplex Negative Control Probe (Cat. No. 321831). The probe targets the bacterial *dapB* gene to control for background noise, and help interpret assay results.



## Reagents/Materials provided by Researcher

□ Supertrost® Plus slides (Fisher Scientitic Part No.12-550-15) are required. Other slide types r result in tissue detachment.	may
□ ProLong® Gold Antifade Reagent (Life Technologies, Part No. P36930).	
□ Fresh reagents (e.g. ethanol and xylene).	
□ 20X SSC (final concentration needed is 5X SSC).	
□ Fresh 10% NBF (neutral-buffered formalin).	
□ Hotplate, drying oven, water bath, thermometer, and microscope.	
□ Paraffin wax, microtome, and fume hood.	
☐ Purchased from Akoya Biosciences. For 3-plex assays, we recommend the following dyes at	a

Opal 520 Reagent Pack (Part No. FP1487001KT)

dilution range of 1:750 – 1:3000:

- Opal 570 Reagent Pack (Part No. FP1488001KT)
- Opal 690 Reagent Pack (Part No. FP1497001KT)
- □ Purchased from Akoya Biosciences. For 4-plex assays, we recommend the following dyes at a dilution range of 1:750 1:3000:
  - Opal 520 Reagent Pack (Part No. FP1487001KT)
  - Opal 570 Reagent Pack (Part No. FP1488001KT)
  - Opal 620 Reagent Pack (Part No. FP1495001KT)
  - Opal 690 Reagent Pack (Part No. FP1497001KT)

**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® and HybEZ™ are trademarks of Advanced Cell Diagnostics, Inc. All other trademarks belong to their respective owners.