

Expand your capabilities with RNAscope™ ISH Assays on the BOND RX from Leica Biosystems

RNAscope is the gold-standard for RNA *in situ* hybridization, trusted by top academic institutions, pharmaceutical companies and service providers around the world. Backed by over 8,000 peer reviewed publications, RNAscope delivers industry leading single-molecule sensitivity with unrivaled target specificity in morphological context.

Getting Started is Easy

Extensive Probe Catalog

Select probes to your genes of interest from an extensive online catalog with over 45,000 RNA probes to over 400 species.

Fast and Flexible Custom Probe Design

If we don't have the probe you need, we can easily design it for you to virtually ANY gene in ANY species for use in ANY tissue.

Expert Technical Support

Experienced field applications specialists in RNAscope ISH assay development on the BOND RX from Leica Biosystems will help you get up and running smoothly with each new application.

Visualize more Targets

Difficult Antigens

- No antibody available.
- Poor antibody performance.
- Secreted proteins including cytokines, chemokines, and growth factors.

RNA Biomarkers

- Any expressed mRNA, non-coding RNA, ASO, siRNA, miRNA, splice variant.
- RNA and protein co-detection.
- Viruses and microbial agents.

Engineered Cells and Vectors

- CAR-T and TCR-T cells.
- Gene therapy vectors (AAV, Lentivirus).
- Transgenes and reporters.

“RNAscope assay on BOND RX works beautifully. The support is everything you could possibly want. I highly recommend this combination!”

-Scientist II, Bristol-Myers Squibb Company

Simple RNAscope Assay Workflow

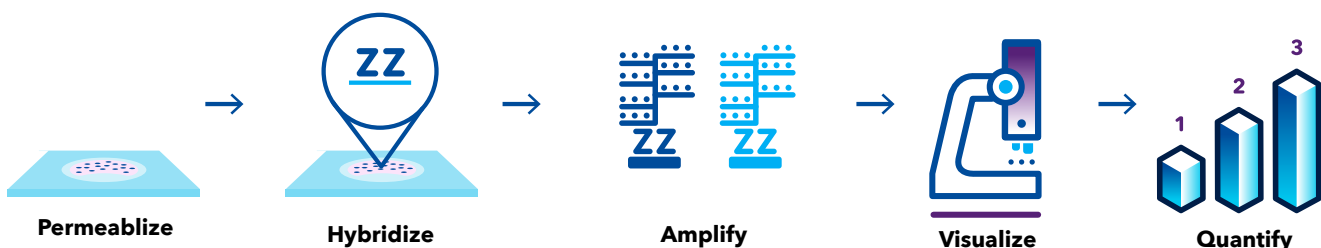
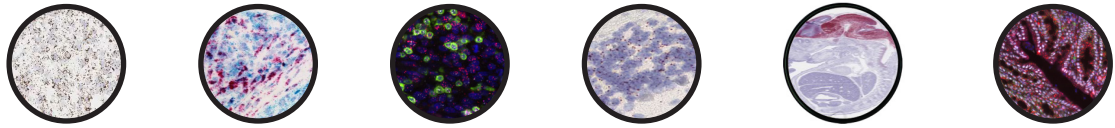


Fig 1. Get robust and reliable same-day results in 5 easy steps. Gold-standard RNAscope signal amplification and background suppression on the BOND RX research staining platform from Leica Biosystems delivers excellent specificity and sensitivity with scalability to over 100 slides per week.

Comprehensive Automated Assay Portfolio



	RNAscope™ (Single-Plex)	RNAscope™ (DuPlex)	RNAscope™ (Multi-Plex V2)	BaseScope™ (Single-Plex)	miRNAscope™ (Single-Plex)	RNAscope™ Plus smRNA-RNA
Visualization	Chromogenic	Chromogenic	Fluorescence	Chromogenic	Chromogenic	Fluorescence
'Plexing' Capability	Single Plex	Duplex	4-Plex	Single Plex	Single Plex	4-Plex: One small RNA and up to three mRNAs
Chromogen/Fluorophore Employed	DAB - HRP or Fast Red - AP	Green - HRP Fast Red - AP	ACD's Vivid™ 520, 570 and 650 (and Akoya's Opal™ 690, for 4-Plex assays)	Fast Red - AP	Fast Red - AP	ACD's Vivid™ 520, 570 and 650 (and Akoya's Opal™ 690, for 4-Plex assays)
Target size	>300 nt			50-300 nt	17-50 nt	

Target Type



Fig 2. Optimized RNAscope assays designed for use on the BOND RX research staining platform from Leica Biosystems. Easily select from a menu of chromogenic and fluorescent assays to visualize your desired targets.

Order Information

Catalog	Product Name	Catalog	Product Name
323600	BaseScope™ LS Reagent Kit	322700	RNAscope™ 2.5 LSx Reagent Kit-BROWN
324600	miRNAscope LS Reagent Kit - RED	322750	RNAscope™ 2.5 LSx Reagent Kit-RED
322440	RNAscope™ 2.5 LS Duplex Reagent Kit-(RED/BROWN)	322800	RNAscope™ LS Multiplex Fluorescent Reagent Kit
322550	RNAscope™ 2.5 LS Green Accessory Pack	323275	RNAscope™ LS Multiplex Fluorescent Reagent Kit with TSA Vivid Dyes
322100	RNAscope™ 2.5 LS Reagent Kit-BROWN	322830	RNAscope™ LS 4-Plex Ancillary Kit for LS Multiplex Fluorescent
322150	RNAscope™ 2.5 LS Reagent Kit-RED	322786	RNAscope™ Plus smRNA-RNA LS Reagents Kit

For Research Use Only. Not for use in diagnostic procedures



For more information Visit Us here!

Bio-Techne® | R&D Systems™ Novus Biologicals™ Tocris Bioscience™ ProteinSimple™ ACD™ ExosomeDx™ Asuragen®

For research use purposes only. Not for use in diagnostic procedures. Trademarks and registered trademarks are the property of their respective owners.

4757728105_SBD_FL_Leica-Sales_TA MK 51-183 Rev A

