

## miRNAscope<sup>™</sup> HD Reagent Kit – RED

Use the miRNAscope" HD Reagent Kit – RED (Cat.No.324500), along with the miRNAscope" Probe–New Target Description Probe, for highly sensitive and highly specific in situ detection of RNA in slide-mounted samples. Each kit contains sufficient reagents in a convenient, Ready-To-Use (RTU) format to perform 20 slide/tests. View the results of the assay with a standard brightfield microscope.

Contents and Storage The components of the miRNAscope<sup>™</sup> HD Reagent Kit – RED and quantities supplied are listed in the following table. The contents have a shelf life until the date of expiration indicated on the kit\_label. Follow the product storage conditions as specified on the label so that the reagents maintain their reliability.

Cat. No.	Sub-Kit	Component	Quantity	Storage
322381	RNAscope® H <sub>2</sub> O <sub>2</sub> and Protease Reagents	RNAscope <sup>®</sup> Hydrogen Peroxide	4 mL x 2 bottles	2–8°C
		RNAscope <sup>®</sup> Protease III*	4.5 mL x 2 bottles	2–8°C
		RNAscope <sup>®</sup> Protease Plus	4.5 mL x 2 bottle	2–8°C
		RNAscope <sup>®</sup> Protease IV	4.5 mL x 2 bottle	2–8°C
322000	RNAscope® Target Retrieval Reagents	RNAscope <sup>®</sup> Target Retrieval (10X)	70 mL x 4 bottles	15–30°C
324510	miRNAscope <sup>™</sup> HD Detection Reagents – RED	miRNAscope <sup>™</sup> HD AMP 1	3 mL x 1 bottle	2–8°C
		miRNAscope <sup>™</sup> HD AMP 2	4.5 mL x 1 bottle	2-8°C
		miRNAscope <sup>™</sup> HD AMP 3	3 mL x 1 bottle	2–8°C
		miRNAscope <sup>™</sup> HD AMP 4 - RED	4.5 mL x 1 bottle	2 <b>-</b> 8°C
		miRNAscope <sup>™</sup> HD AMP 5 - RED	4.5 mL x 1 bottle	2-8°C
		miRNAscope <sup>™</sup> HD AMP 6 - RED	3 mL x 1 bottle	2 <b>-</b> 8°C
		miRNAscope <sup>™</sup> Fast RED-A	3 mL x 1 bottle	2-8°C
		miRNAscope <sup>™</sup> Fast RED-B	50 µL x 1 bottle	2 <b>-</b> 8°C
310091	RNAscope® Wash Buffer Reagents	RNAscope® Wash Buffer (50X)	60 mL x 4 bottles	15–30°C

Pretreatment Reagents based on different tissue types. Do not interchange the reagent components of the Detection Reagents, even those having the same name. Refer to the latest version of the miRNAscope<sup>™</sup> HD (RED) Assay with Sample Preparation and Pretreatment User Manual (Doc. No. UM-324510) available online at https://acdbio.com/technical-support/user-manuals.

Handle biological specimens and other materials according to local laboratory guidelines and regulations. Avoid Precautions cross-contamination of samples and reagents to ensure quality results. Use equipment and reagents as specified by the manufacturer. For safety and biohazard guidelines, refer to the "Safety" section in the user manual. For every chemical, read the Safety Data Sheets (SDSs) and follow the handling instructions available at https://acdbio.com/documents/product-documents. Wear appropriate protective eyewear, clothing, and gloves.

Contact Information	Address:	7707 Gateway Blvd CA 94545 USA			
	Phone:	1-877-576-3636 option 1	Fax:	1-510-576-8801	
	Web:	www.acdbio.com	Email:	support.acd@bio-techne.com	

Copyright ©2020. Advanced Cell Diagnostics, Inc. All rights reserved.

Trademarks

miRNAscope and RNAscope are trademarks or registered trademarks of Advanced Cell Diagnostics, Inc. All other trademarks belong to their respective owners. Advanced Cell Diagnostics, Inc. reserves the right to change its products and services at any time to incorporate technological developments. This insert is subject to change without notice. Although this insert has been prepared with every precaution to ensure accuracy, Advanced Cell Diagnostics, Inc. assumes no liability for any errors, omissions, or for any damages resulting from the use of this information