

HT RNA Reagent Kit

Product Number: CLS960010

Lot Number: CC01RK01

Expiration Date: 28 JAN 2025

PARAMETER	SAMPLE	SPECIFICATION	RESULT
Marker Migration	Ladder Well	Meets spec	Pass
Ladder Peak Migration	Ladder Well	Meets spec	Pass
Peak Height to Marker	Ladder Sample	2 to 5	Pass
Ladder Concentration	Ladder Sample	432 to 528	Pass
Concentration Reproducibility	Ladder Sample	Tracking only	Pass
Ladder Peak Migration	Ladder Sample	Within 10% of migration time of ladder well	Pass
Ladder Purity	Ladder Sample	No extra peaks in ladder	Pass
28S/18S Ratio	RB Sample	Meets spec	Pass
Peak height Linearity	RB Sample	No Value deviates more than 20% from line	Pass
Marker Purity	Blank Sample	No extra peaks from marker	Pass
False Peaks	Blank Sample	No more than 1	Pass
Quality Score	RB Sample	No value less than 6.5	Pass

Reagent	Vial	Quantity
RNA Dye Concentrate	Blue ●	1 vial of 0.5 mL
Chip Storage Buffer	White ○	5 vials, 1.8 mL each
RNA Gel Matrix	Red ●	5 vials, 0.51 mL each
RNA Ladder (packed separately)	Yellow ●	2 vial, 0.03 mL
RNA Marker	Green ●	1 vial, 0.8 mL
10X RNA Sample Buffer Concentrate	Purple ●	6 vials, 2.0 mL each
Consumable Items	PN 701672	1

The information provided in this document is valid for the specified lot number and date of analysis. This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.

CLS960010-R Rev01