

HT RNA Pico Reagent Kit

Product Number: CLS960012

Lot Number: CF05RK01

Expiration Date: 21 APR 2025

PARAMETER	SAMPLE	SPECIFICATION	RESULT
Marker Migration Time	Ladder Samples	Meets spec	Pass
Ladder Peak Migration Time	Ladder Samples	Meets spec	Pass
Ladder Concentration	Ladder Samples	Mean is within +/- 30% of nominal value for each ladder concentration	Pass
Ladder Purity	Ladder Samples	No extra peak taller than the last ladder for 10ng/ul, 20ng/ul ladder samples	Pass
Ladder Peak Ratio	Ladder Well	Average of ladder peak height ratio [200nt / 6000nt] is within 2.5 to 5.1	Pass
Quality Score	Rat Brain Sample	Min & Max values are within 10% of each other	Pass

Reagent	Vial	Quantity
RNA Dye Concentrate	Blue ●	1 vial, 0.5 mL
Chip Storage Buffer	White ○	5 vials, 1.8 mL each
RNA Pico Gel Matrix	Red ●	5 vials, 0.51 mL each
RNA Pico Ladder (packaged separately, Part Number CLS760652)	Yellow ●	2 vials, 0.04 mL
RNA Pico Marker	Green ●	1 vial, 0.8 mL
10X RNA Sample Buffer Concentrate	Purple ●	3 vials, 2 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.

CLS960012-R Rev01