

## HT Small RNA Reagent Kit

Product Number: CLS153530  
 Lot Number: CL20RK01  
 Expiration Date: 20 AUG 2025

PARAMETER	SAMPLE	SPECIFICATION	RESULT
Lower Marker Migration Time	Ladder Samples	25.5 – 30 s	Pass
Ladder Peak Migration Time	Ladder Samples	41 – 47 s	Pass
Resolution	RNA sample mix	Achieve greater than FWHM separation at: 5 nt difference for RNA ( 20-100 nt) 10% difference in RNA MW for RNA (100-150 nt)	Pass
Ladder Purity	Ladder Samples	No extra peak taller than the 30% of 40 nt ladder peak height	Pass
Limit of Detection	LOD test sample (on-plate 50 pg/μL)	Peak height > 1.2 rfu	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 CLS153594)	Yellow ●	2 vials, 0.05 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

Additional Small RNA Ladder can be ordered separately using Part Number CLS153594.

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.

CLS153530-R Rev01