

HT RNA Pico Reagent Kit

Product Number: CLS960012
 Lot Number: CK28RK01
 Expiration Date: 28 OCT 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