

HT DNA NGS 3K Reagent Kit

Product Number: CLS960013
 Lot Number: BL17RK01
 Expiration Date: 16AUG2024

| PARAMETER | SAMPLE | SPECIFICATION | RESULT |
|---------------------------------------|-------------------------------------|--|--------|
| Sizing | PhiX Sample | Mean of each of peaks 3-12 must be within +/- 10% of nominal | Pass |
| Resolution | PhiX Sample | Value > 1.5 for 10% size difference | Pass |
| Concentration | Gene Ruler | Mean of each of peaks 2-9 must be within +/- 30% of nominal | Pass |
| Upper Marker Migration Time | All Samples | Mean across all samples within 50-62 seconds | Pass |
| Marker/Sample Peak Area Ratio (UM/9A) | All Wells in row A (All PhiX wells) | Ratio must be >1 | Pass |

| Reagents | Label | Quantity |
|----------------------------|-----------|----------------------|
| DNA NGS 3K Dye Concentrate | Blue ● | 1 vial, 0.09 mL |
| Chip Storage Buffer | White ○ | 9 vials, 1.8 mL each |
| DNA HiSens Gel Matrix | Red ● | 5 vials, 1.1 mL each |
| 10X DNA NGS 3K Ladder | Yellow ● | 1 vial, 0.26 mL |
| DNA NGS 3K Marker | Green ● | 1 tube, 45 mL |
| 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.

CLS960013-R Rev01