

## HT DNA 12K Reagent Kit

Product Number: 760569  
 Lot Number: CD23RK01  
 Expiration Date: 22 JAN 2025

| PARAMETER                   | SAMPLE      | SPECIFICATION  | RESULT |
|-----------------------------|-------------|--|--------|
| Sizing                      | PhiX Sample | Mean of each of peaks 3-12 must be within +/- 10% of nominal | Pass   |
| Resolution                  | PhiX Sample | 10% resolution $\geq 0.83$                                   | Pass   |
| Concentration               | Gene Ruler  | Mean of each of peaks 2-10 must be within +/- 40% of nominal | Pass   |
| Marker Migration Time       | All Wells   | Meets spec   | Pass   |
| Marker Height Ratio (LM/UM) | All Wells   | Ratio must be within 2 to 6                                  | Pass   |

| Reagent             | Vial      | Quantity             |
|---------------------|-----------|----------------------|
| DNA Dye Concentrate | Blue ●    | 1 vial of 0.08 mL    |
| Chip Storage Buffer | White ○   | 5 vials, 1.8 mL each |
| DNA Gel Matrix      | Red ●     | 3 vials, 1.6 mL each |
| 10X DNA Ladder      | Yellow ●  | 1 vial, 0.15 mL      |
| DNA Marker          | Green ●   | 1 vial, 1.5 mL       |
| Consumable Items    | PN 701553 | 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.

760569-R Rev01