

## HT DNA NGS 3K Reagent Kit

Product Number: CLS960013  
 Lot Number: CK24RK01  
 Expiration Date: 24 JUL 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	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