

## HT DNA 5K Reagent Kit

Product Number: CLS760675  
 Lot Number: CA18RK01  
 Expiration Date: 17OCT2024

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 +/- 30% of nominal	Pass
Marker Migration Time	All Wells	Meets spec	Pass
Marker Height Ratio (LM/UM)	All Wells	Ratio must be within 2 to 5	Pass

Reagents	Label	Quantity
DNA Dye Concentrate	Blue ●	1 vial, 0.08 mL
Chip Storage Buffer	White ○	7 vials, 1.8 mL each
DNA Gel Matrix	Red ●	5 vials, 1.1 mL each
10X DNA Ladder	Yellow ●	1 vial, 0.265 mL
DNA Marker	Green ●	1 vial, 1.2 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.

CLS760675-R Rev01