

Genomic DNA Reagent Kit

Product Number: CLS760685
 Lot Number: CB06RK01
 Expiration Date: 05 NOV 2024

PARAMETER	SAMPLE	SPECIFICATION	RESULT
Sizing Reproducibility	100 bp-7k bp Fragments	Within +/- 20%	Pass
Sizing Accuracy	100 bp-7k bp Fragments	Within +/- 20%	Pass
Concentration Reproducibility	500 bp-7k bp Fragments 10k bp-20/40k bp Fragments	Within +/- 20% Within +/- 30%	Pass
Concentration Accuracy	500 bp-7k bp Fragments 10k bp-20/40k bp Fragments	Within +/- 30% Within +/- 50%	Pass
Lower Marker Migration Time	Lower Marker from all Wells	Meets spec	Pass
Last Peak Migration Time	20k/40k Fragment	Meets spec	Pass
MigTime Reproducibility	10k bp - 20k/40k bp Fragments	< 5%	Pass

Reagents	Label	Quantity
DNA Dye Concentrate	Blue ●	1 vial, 0.09 mL
Chip Storage Buffer	White ○	9 vials, 1.8 mL each
Genomic DNA Gel Matrix	Red ●	5 vials, 1.1 mL each
10X Genomic DNA Ladder	Yellow ●	1 vial, 0.26 mL
Genomic DNA Marker	Green ●	1 vial, 1.5 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.

CLS760685-R Rev01