

## cfDNA Reagent Kit

Product Number: CLS157242  
 Lot Number: CE21RK01  
 Expiration Date: 06 FEB 2025

PARAMETER	SAMPLE	SPECIFICATION	RESULT
Sizing Accuracy (1)	200 bp DNA fragment	180-220 nt	Pass
Sizing Accuracy (2)	Gene Ruler	Mean of each peak $\pm 10\%$ to expected	Pass
Concentration Accuracy	200 bp DNA fragment	$\pm 20$ to Qubit measurement	Pass
Low Marker Migration Time	All wells	16.5-18.5 s	Pass
Upper Marker Migration Time	All wells	28-36 s	Pass

Reagent	Vial	Quantity
cfDNA Dye Concentrate	Blue ●	1 vial, 0.09 mL
Chip Storage Buffer	White ○	9 vials, 1.8 mL each
cfDNA Gel Matrix	Red ●	5 vials, 1.1 mL each
cfDNA Ladder	Yellow ●	1 vial, 0.26 mL
cfDNA Marker	Green ●	1 tube, 4.5 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.

CLS157242-R Rev01