

cfDNA Reagent Kit

Product Number: CLS157242
 Lot Number: DB06RK01
 Expiration Date: 06-Nov-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 | ± 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