

ssDNA 7K Assay Reagent Kit

Product Number: CLS158169
 Lot Number: BM05RK01
 Expiration Date: 28NOV2024

PARAMETER	SAMPLE	SPECIFICATION	RESULT
Lower marker migration time	ssDNA 7K ladder sample	Mean is within [31, 37] seconds	Pass
7200nt ladder migration time	ssDNA 7K ladder sample	Mean is within [73, 92] seconds	Pass
Peak quality of 1.1-4kb	ssDNA 7K ladder sample	Mean of FWHM (sec) within 1.50 sec	Pass
LM peak height	ssDNA 7K ladder sample	Mean of LM peak height \geq 15 RFU	Pass
Detection sensitivity of 1.1kb ladder	ssDNA 7K ladder sample	Mean of 1.1kb peak height is \geq 5 RFU for all ladder samples with 1:40 dilution in nuclease-free water	Pass
Ladder quality	ssDNA 7K ladder	All 6 peaks detected	Pass
Ladder purity	ssDNA 7K ladder sample	No extra peak (with peak height repeatedly more than 5 RFU), and only six visible peaks	Pass

Reagent	Vial	Quantity
ssDNA 7K Ladder	● Yellow	1 vial, 0.04 mL
ssDNA Chip Storage Buffer	○ White	5 vials, 1.8mL
ssDNA Gel Matrix	● Red	5 vials, 1.1mL
ssDNA Marker	● Green	1 vial, 0.8 mL
10X Sample Buffer	● Purple	3 vials, 2.0mL
ssDNA Dye Concentrate	● Blue	1 vial, 0.5 mL
Consumable Items	PN 701672	1

Additional ssDNA 7K Ladder can be ordered separately using Part Number CLS157950.

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.

CLS158169-R Rev01