

ssDNA 7K Ladder

Product Number: CLS157950
 Lot Number: 123263
 Expiration Date: 22 SEP 2024

PARAMETER	SAMPLE	SPECIFICATION	RESULT
Ladder Concentration	Ladder Samples	Measured concentration by Nanodrop $\geq 80\text{ng}/\mu\text{L}$	Pass
Ladder Peak Size Call	Ladder Samples	1 kb-5.1kb peaks are called within 10%, 7.2 bp peak is called within 15% of expected size on GX	Pass
Ladder Height	Ladder Samples	All 6 peaks are $\geq 80\%$ of peak height compared to reference ladder or all peaks in test ladder are ≥ 10 RFU	Pass
Ladder Purity	Ladder Samples	No extra peak taller ≥ 5 RFU and six visible peaks	Pass

Reagent	Vial	Quantity
ssDNA 7K Ladder	● Yellow	1 vial, 0.04 mL

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.

CLS157950-R Rev01