

Single-Stranded DNA 7K Ladder Insert

STORE AT -25°C to -15°C UPON RECEIPT

P/N CLS157950: 1 vial of 40 µL of ssDNA 7K Ladder
For Research Use Only, not for use in diagnostic procedures

Properties: Colorless solution, volume per vial: 40 µL,
total concentration: 70-110 ng/µL.

Storage buffer components: 50 mM Potassium Acetate, 20mM Tris-acetate, 10mM Magnesium Acetate, 20mM EDTA, ~6% Glycerol, pH ~7.8.

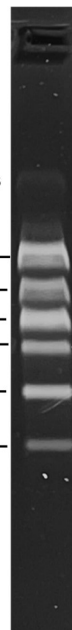
Storage: Store at -25°C to -15 °C. Avoid multiple freeze-thaw cycles. Aliquots of the product may be stored for up to one week at 2°C to 8°C. Do not store in a frost-free freezer.

Handling: Use DNase-free and RNase-free reagents, DNA Low binding tubes, and barrier pipette tips.

Thawing instructions: Thaw just to completion at up to 37°C, vortex a few seconds when fully thawed, and place on ice. To avoid multiple freeze-thaw cycles, make aliquots in DNase-free and RNase-free, DNA Low binding tubes with the required amount typical for daily use. The ssDNA 7K ladder has no obvious stability loss after 3 freeze-thaw cycles.

Shelf life upon receipt: With recommended storage, until expiration specified on the vial.

Mass (ng)	Kilobases
20	7.2
14.2	5.1
22.2	4
8.9	3.2
11.7	2.1
3.1	1.1



**0.8% 0.5 X TBE Agarose
~80 ng**

Figure 1: ssDNA 7K Ladder by Safe Stains (SYBR Gold) visualized on a 0.8% TBE agarose gel using Gel Imaging System with Blue Light Illumination. Mass values are for ~80 ng/lane. Suggested Load: 0.5 to 2 µl/lane.

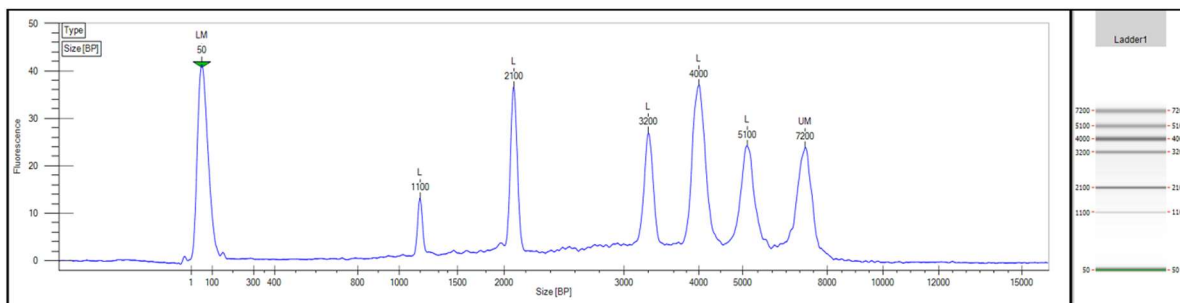


Figure 2: ssDNA 7K on LabChip® GXII Touch™ (P/N CLS138160). The sample (1µL ssDNA 7K Ladder + 39 µL nuclease-free H2O) was run by ssDNA 7K.asyx (P/N CLS157952) program using RNA Pico reagent (P/N CLS960012) or ssDNA 7K reagent (P/N CLS158169) and DNA 5K/RNA/CZE HT LabChip® (P/N 760435).