

FFA120 Ladder Insert

STORE AT -25°C to -15°C UPON RECEIPT

P/N CLS158127: 1 vial of 120 µL of FFA120 Ladder

For Research Use Only, not for use in diagnostic procedures

Properties: Colorless solution, volume per vial: 120 µL.

Storage buffer components: < 20mM Tris, < 10mM other salts, pH ~8.0.

Storage: Store at -25°C to -15 °C. Avoid multiple freeze-thaw cycles. 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 the FFA120 Ladder on ice or at 4 °C. Vortex the thawed FFA120 ladder for 10-15 seconds to mix and then spin down for a few seconds before each use. To avoid multiple freeze-thaw cycles, make 65 µL aliquots in DNase-free and RNase-free, DNA Low binding tubes for daily use. Store the aliquots at -20 °C until the expiration date. The FFA120 ladder has no obvious stability loss after 6 freeze-thaw cycles.

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

Ladder gel image on LabChip® GXII Touch™ (P/N CLS138160) using Small RNA Reagent (P/N CLS153530) on the chip of DNA 5K/RNA/CZE HT LabChip® (P/N 760435) using FFA120.asyx program.

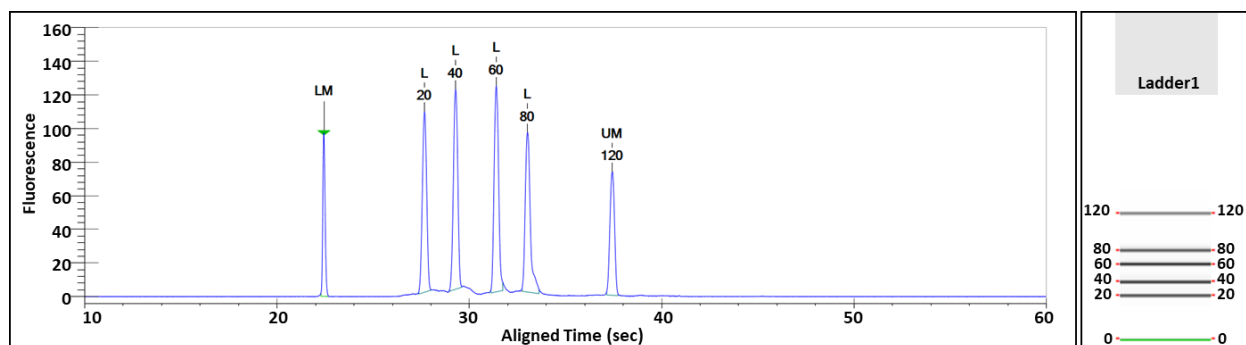


Figure: Example of ladder performance on LabChip® GXII Touch™ with 1:10 dilution by DMSO.