



CE-LIF dsDNA 100 Base Pair Ladder

Catalog Number 148-2021

Contents	This size standard is a mixture of 10 double-stranded DNA fragments ranging in length from 100 bp to 1,000 bp in exact 100 bp increments.
Quantity	200 μ l supplied in TE (10 mM Tris-HCl, 1 mM EDTA, pH 8.0).
Concentration	1 μ g/ml
Storage	The size standard should be stored at 4 °C. The standard can be stored at -20 °C in aliquots for long term storage. Use only sterile pipet tips when removing aliquots. Introduction of nucleases will shorten shelf life.
Use	Under the conditions described in the instruction manual for CE dsDNA 1000 Fluorescent Detection and Analysis Kit, the standard will exhibit 10 major peaks, see Figure 1. Pipette 20–50 μ l of the CE-LIF dsDNA 100 bp Ladder into a 500 μ l sample vial.

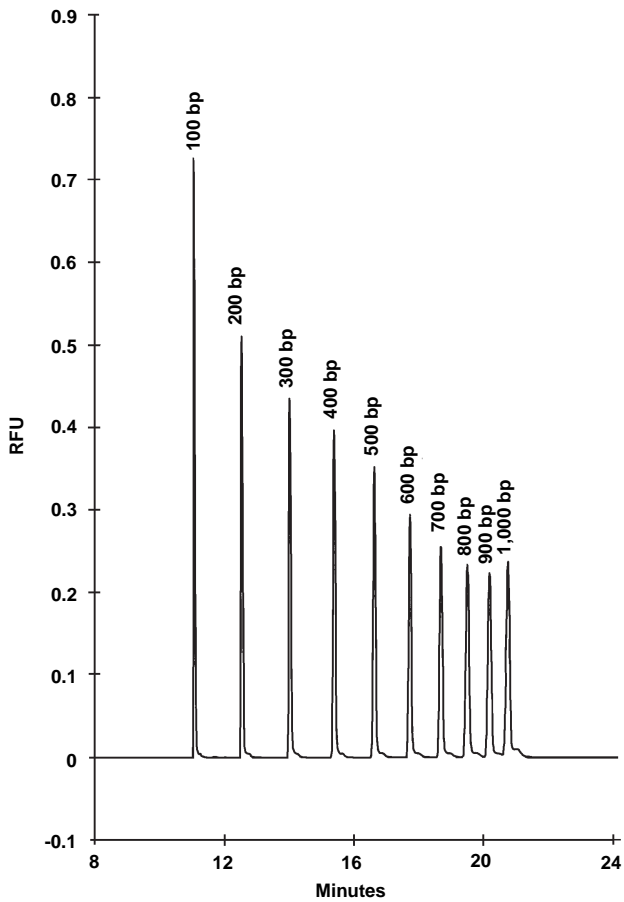


Fig. 1. Separation of 100 bp ladder.