



CE-LIF dsDNA Mid Range Size Standard

Catalog Number 148-2020

Contents	This double-stranded DNA size standard is a proprietary mixture of blunt-ended dsDNA of precise length and known sequence. There are six fragments of 1,000, 1,500, 2,000, 2,500, 3,000, and 4,361 bp.
Quantity	200 μ l supplied in TE (10 mM Tris-HCl, 1 mM EDTA, pH 8.0)
Concentration	1 ng/fragment/ μ l
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 4000 Fluorescent Detection and Analysis Kit, the standard will exhibit 6 major peaks, see Figure 1. Pipette 20–50 μ l of the CE-LIF dsDNA Mid Range Size Standard into a 500 μ l sample vial.

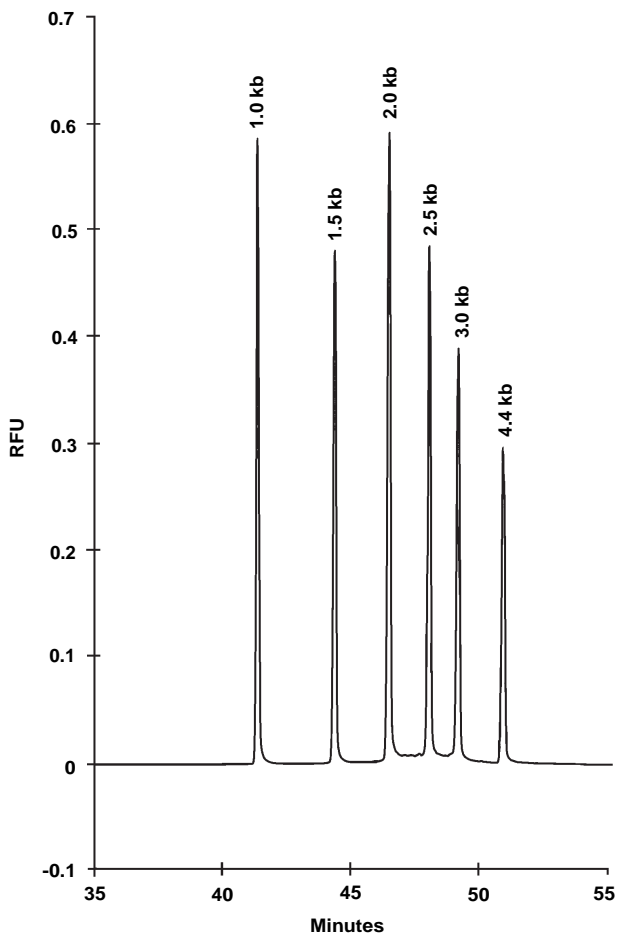


Fig. 1. Separation of the mid range standard.