



## CE-LIF dsDNA Mid Range Size Standard

**Catalog Number 148-2020**

<b>Contents</b>	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.
<b>Quantity</b>	200 $\mu$ l supplied in TE (10 mM Tris-HCl, 1 mM EDTA, pH 8.0)
<b>Concentration</b>	1 ng/fragment/ $\mu$ l
<b>Storage</b>	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.
<b>Use</b>	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 $\mu$ l of the CE-LIF dsDNA Mid Range Size Standard into a 500 $\mu$ l sample vial.

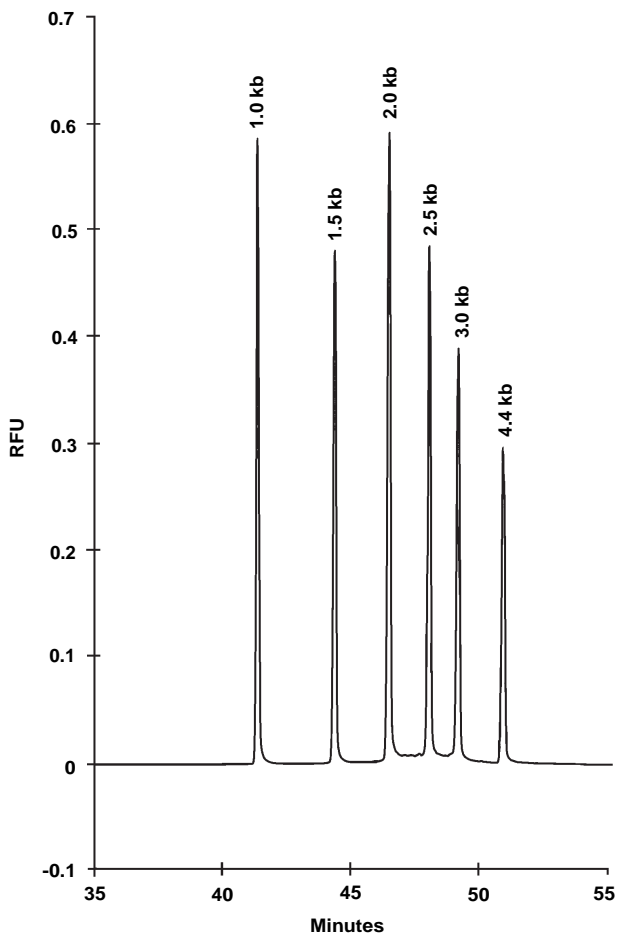


Fig. 1. Separation of the mid range standard.