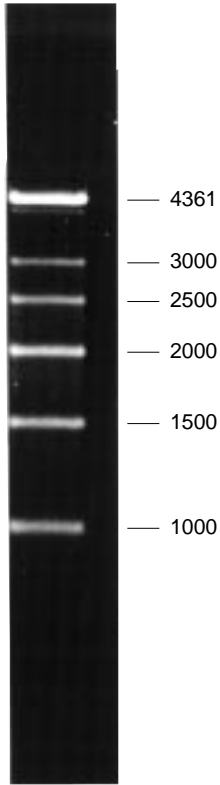


## **DNA Size Standards — 1-4.4 kb Ladder**

**Catalog Number 170-8210**

<b>Contents</b>	Proprietary blunt-ended double-stranded DNA of precise length and known sequence. Six fragments of 1000, 1500, 2000, 2500, 3000 and 4361 bp.
<b>Quantity</b>	250 $\mu$ l. Sufficient for 50 lanes when used at 5 $\mu$ l per lane.
<b>Concentration</b>	60 ng/ $\mu$ l
<b>Storage</b>	4 °C
<b>Shelf Life</b>	1 yr at 4 °C. <b>Use only sterile pipet tips when removing aliquots.</b> Introduction of nucleases will shorten shelf life.
<b>Use</b>	These standards are used to estimate molecular weights of DNA fragments separated on agarose gels. <b>Use only sterile solutions.</b>



**DNA Size Standards, 1-4.4 kb** separated on 1.5% Molecular Biology Certified Agarose with TAE buffer.