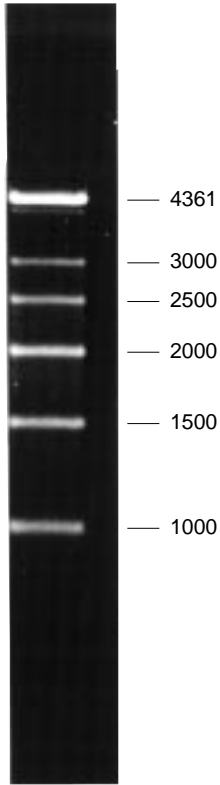


DNA Size Standards — 1-4.4 kb Ladder

Catalog Number 170-8210

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



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