



DNA Size Standards—5 Kb Ladder

Catalog Number 170-3624

- Use:** The 5 kb ladder DNA size standard is used to estimate molecular weights of DNA samples separated on agarose gels. This standard allows very accurate measurements of sizes from 5 kb to about 100 kb because of the predictable spacing between bands. Always use sterile solutions, pipet tips, and tubes. Approximately 8-10 μ l of the sample per lane is sufficient to see all the bands by ethidium bromide staining.
- Contents:** The 5 kb ladder DNA size standard consists of ligated pBR328 DNA partially digested with EcoRI. The standards are supplied in 28 mM Tris, pH 8.0, 50 mM EDTA, 28 mM NaCl, 0.05% SDS.
- Quantity:** 20 μ g
- Concentration:** 0.1 μ g/ μ l
- Storage:** 4 °C for short term storage. -20 °C for long term storage.
- Shelf life:** One year at -20 °C. Six months at 4 °C. Use only sterile pipet tips when removing aliquots. Introduction of nucleases can shorten the shelf life.
- Sizes:** A ladder starting from 4.9 kilobases¹ and extending to approximately 98 kilobases.
- Reference:** 1. Soberon, X., *et al.*, *Gene*, **9**, 287-305 (1980).

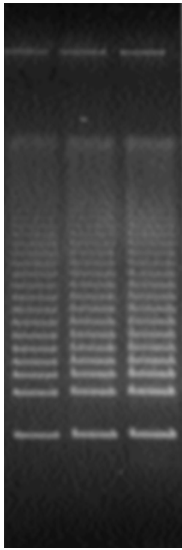


Fig. 1.5kb Ladder. Electrophoresis was done on the CHEF-DR® II system at 6 V/cm, 1% Molecular Biology Grade Agarose, 0.5x TAE buffer, and temperature at 14 °C. Switch time was a linear ramp from 1 second to 6 seconds with a run time of 11 hours.