

200 bp DNA Ladder 200 bp to 2 kb log scale ready-to-use

Cat. No.	Volume	Amount
M-215	500 μ l	100 lanes

For *in vitro* use only
Quality guaranteed for 12 months
Store at -20°C , short term storage (up to 4 weeks) at 4°C possible

Form

Liquid. DNA fragment mix in 10 mM Tris-HCl (pH 7.5), 10 mM EDTA, 10% (w/v) glycerol, bromophenol blue and xylene cyanol FF

Concentration

105 ng/ μ l

Description

200 bp DNA Ladder consist of 10 DNA fragments ranging from 200 bp to 2 kb and is supplied in ready-to-use format containing blue tracking dyes. It allows sizing and concentration estimate of DNA fragments on agarose gels generated by PCR or restriction digest. The fragment mix yields a ladder of the following double-stranded DNA fragments:

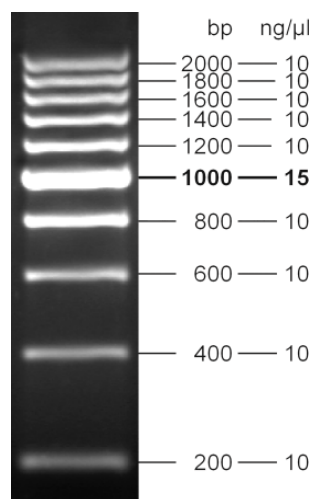
200, 400, 600, 800, **1000**, 1200, 1400, 1600, 1800, 2000 bp

DNA log scale ladders are well suited for combination if small and large fragments should be analyzed on one gel.

Recommended Load

5 μ l DNA ladder per lane

200 bp DNA Ladder



1.8%
Agarose