

500 bp DNA Ladder 500 bp to 5 kb log scale ready-to-use

Cat. No.	Volume	Amount
M-216	500 µl	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/µl

Description

500 bp DNA Ladder consists of 10 DNA fragments ranging from 500 bp to 5 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:

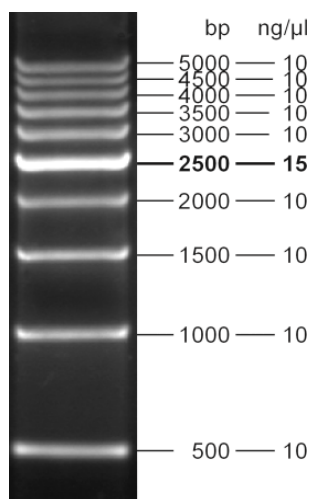
500, 1000, 1500, 2000, **2500**, 3000, 3500, 4000, 4500, 5000 bp

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

Recommended Load

5 µl DNA ladder per lane

500 bp DNA Ladder



1.2%
Agarose