

100 bp DNA Ladder

100 bp to 1 kb log scale

ready-to-use

Cat. No.	Volume	Amount
M-214	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

Concentration

105 ng/µl

Recommended Load

5 µl DNA ladder per lane

DNA Fragments

100, 200, 300, 400, **500**, 600, 700, 800, 900, 1000 bp

Form

DNA fragment mix in 10 mM Tris-HCl (pH 7.5), 10 mM EDTA, 10% (w/v) Glycerol, Orange G and Xylene Cyanol FF

Description

100 bp DNA Ladder consists of 10 DNA fragments ranging from 100 bp to 1 kb and is supplied in ready-to-use format containing orange and blue tracking dyes. It allows sizing and concentration estimate of DNA fragments on agarose gels generated by PCR or restriction digest. DNA log scale ladders are well suited for combination if small and large fragments should be analyzed on one gel.

100 bp DNA Ladder

