

Mid Range DNA Ladder 100 to 3000 bp linear scale ready-to-use

Cat. No.	Volume	Amount
M-203	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, orange G and xylene cyanol FF

Concentration

130 ng/µl

Description

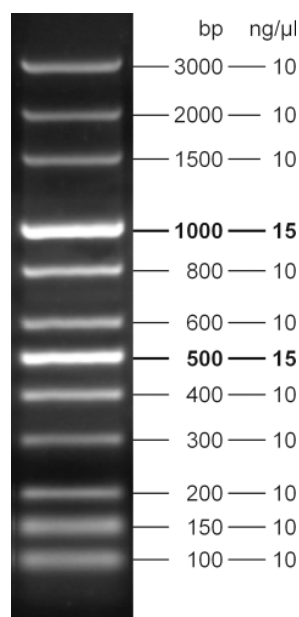
Mid Range DNA Ladder consists of 12 DNA fragments ranging from 100 bp to 3000 bp. It is designed to show virtually uniform spacing over a wide fragment range. The ladder allows sizing and concentration estimate of DNA fragments on agarose gels generated by PCR or restriction digest. The fragment mix is supplied in ready-to-use format containing orange and blue tracking dyes. It yields a ladder of the following double-stranded DNA fragments:

100, 150, 200, 300, 400, **500**, 600, 800, **1000**, 1500, 2000, 3000 bp

Recommended Load

5 µl DNA ladder per lane

Mid Range DNA Ladder



1.8%
Agarose