

High Range DNA Ladder

500 bp to 10 kb linear scale

ready-to-use

Cat. No.	Volume	Amount
M-204	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, 10 mM Tris-HCl (pH 7.5), 10 mM EDTA, 10% (w/v) glycerol, bromophenol blue and xylene cyanol FF

Concentration

130 ng/µl

Description

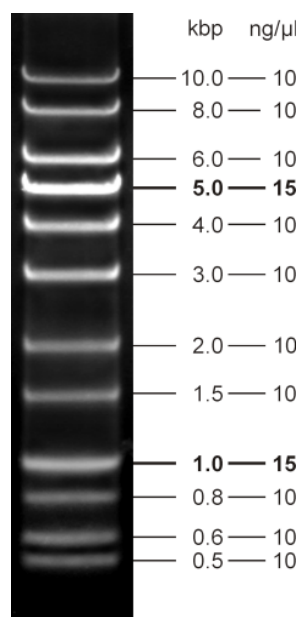
High Range DNA Ladder consists of 12 DNA fragments ranging from 500 bp to 10 kb. It is designed to show virtually uniform spacing over a wide fragment range. The ladder allows sizing and concentration estimate of DNA fragments on agarose gels generated by PCR or restriction digest. The fragment mix is supplied in ready-to-use format containing blue tracking dyes. It yields a ladder of the following double-stranded DNA fragments:

0.5, 0.6, 0.8, **1**, 1.5, 2, 3, 4, **5**, 6, 8, 10 kb

Recommended Load

5 µl DNA ladder per lane

High Range DNA Ladder



0.6%
Agarose