

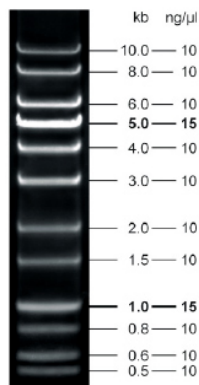


High Range DNA Ladder

500 bp to 10 kb linear scale
ready-to-use, blue / blue

Cat. No.	Amount
M-204S	500 µl (130 ng/µl)
M-204L	5 x 500 µl (130 ng/µl)

High Range DNA Ladder



0.6% Agarose

5 µl loaded onto 0.6 % agarose

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Short term storage (up to 1 month) at 4 °C possible.

Shelf Life: 12 months

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

Color: blue

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.

Recommended Load:

5 µl per lane

DNA Fragments:

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