

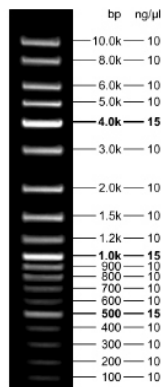


Long Range DNA Ladder

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

Cat. No.	Amount
M-205S	500 µl (205 ng/µl)
M-205L	5 x 500 µl (205 ng/µl)

Long Range DNA Ladder



1% Agarose

5 µl loaded onto 1 % 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, 8 % (w/v) Glycerol, Tartrazine and Xylene Cyanol FF

Color: green

Concentration: 205 ng/µl

Description:

Long Range DNA Ladder consists of 19 DNA fragments ranging from 100 bp to 10 kb. It is designed to show virtually uniform spacing over an extreme wide fragment range covering virtually the complete application spectrum of DNA ladders. 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.

Recommended Load:

5 µl per lane

DNA Fragments:

100, 200, 300, 400, **500**, 600, 700, 800, 900 bp, **1.0**, 1.2, 1.5, 2.0, 3.0, **4.0**, 5.0, 6.0, 8.0, 10.0 kb