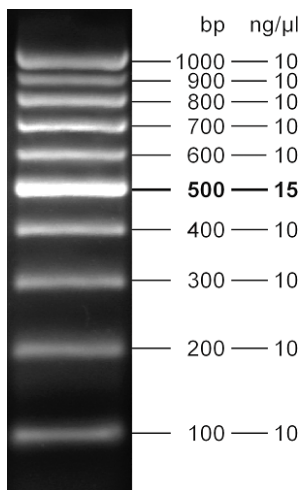




100 bp DNA Ladder

100 bp to 1 kb log scale
ready-to-use, orange / blue

Cat. No.	Amount
M-214S	500 µl (105 ng/µl)
M-214L	5 x 500 µl (105 ng/µl)



5 µl loaded onto 2.4 % 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: 105 ng/µl

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.

Recommended Load:

5 µl per lane

DNA Fragments:

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