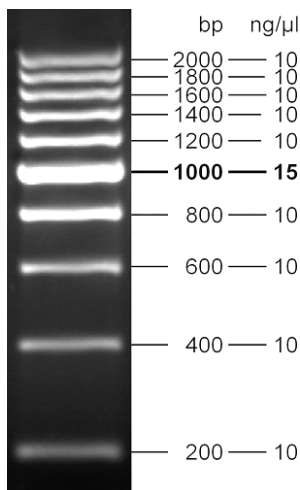




200 bp DNA Ladder

200 bp to 2 kb log scale
ready-to-use, blue / blue

Cat. No.	Amount
M-215S	500 µl (105 ng/µl)
M-215L	5 x 500 µl (105 ng/µl)



5 µl loaded onto 1.8 % 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: 105 ng/µl

Description:

200 bp DNA Ladder consist of 10 DNA fragments ranging from 200 bp to 2 kb and is supplied in ready-to-use format containing 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:

200, 400, 600, 800, **1000**, 1200, 1400, 1600, 1800, 2000 bp