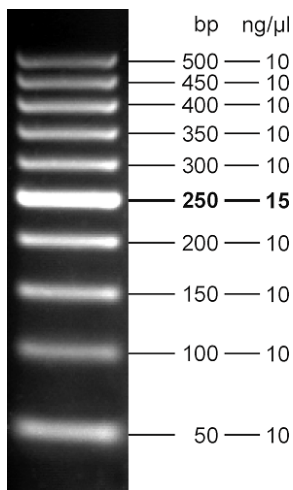




50 bp DNA Ladder

50 bp to 500 bp log scale
ready-to-use, orange

Cat. No.	Amount
M-213S	500 µl (105 ng/µl)
M-213L	5 x 500 µl (105 ng/µl)



5 µl loaded onto 3.0 % agarose

For *in vitro* use only!

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 and Orange G

Color: orange

Concentration: 105 ng/µl

Description:

50 bp DNA Ladder consists of 10 DNA fragments ranging from 50 bp to 500 bp and is supplied in ready-to-use format containing an orange tracking dye. 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:

50, 100, 150, 200, **250**, 300, 350, 400, 450, 500 bp