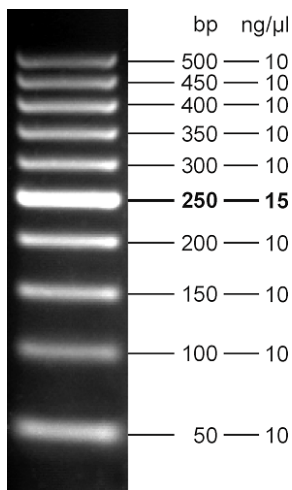




## 50 bp DNA Ladder

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 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 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