

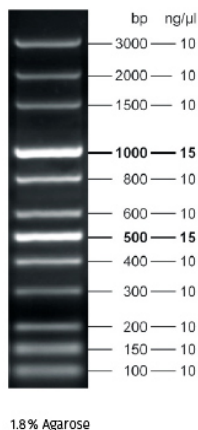


## Mid Range DNA Ladder

100 bp to 3 kb linear scale  
ready-to-use, orange / blue

| Cat. No. | Amount                 |
|----------|------------------------|
| M-203S   | 500 µl (130 ng/µl)     |
| M-203L   | 5 x 500 µl (130 ng/µl) |

Mid Range DNA Ladder



5 µl loaded onto 1.8 % 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, 8 % (w/v) Glycerol, Tartrazine and Xylene Cyanol FF

**Color:** green

**Concentration:** 130 ng/µl

### Description:

Mid Range DNA Ladder consists of 12 DNA fragments ranging from 100 bp to 3 kb. It is designed to show virtually uniform spacing over a wide fragment range. 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, 150, 200, 300, 400, **500**, 600, 800, **1000**, 1500, 2000, 3000 bp