

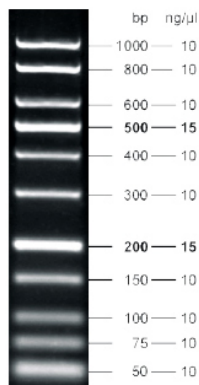


## Low Range DNA Ladder

50 bp to 1 kb linear scale  
ready-to-use, orange / blue

| Cat. No. | Amount                 |
|----------|------------------------|
| M-202S   | 500 µl (120 ng/µl)     |
| M-202L   | 5 x 500 µl (120 ng/µl) |

Low Range DNA Ladder



3.0 % Agarose

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, 8 % (w/v) Glycerol, Tartrazine and Xylene Cyanol FF

**Color:** green

**Concentration:** 120 ng/µl

**Description:**

Low Range DNA Ladder consists of 11 DNA fragments ranging from 50 bp to 1 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:**

50, 75, 100, 150, 200, 300, 400, 500, 600, 800, 1000 bp