

## Low Range DNA Ladder

### 50 bp to 1 kb linear scale

### ready-to-use

Cat. No.	Volume	Amount
M-202	500 µl	100 lanes

For *in vitro* use only  
 Quality guaranteed for 12 months  
 Store at -20°C, short term storage (up to 4 weeks) at 4°C possible

#### Form

Liquid. DNA fragment mix in 10 mM Tris-HCl (pH 7.5), 10 mM EDTA, 10% (w/v) glycerol, orange G and xylene cyanol FF

#### Concentration

120 ng/µl

#### Description

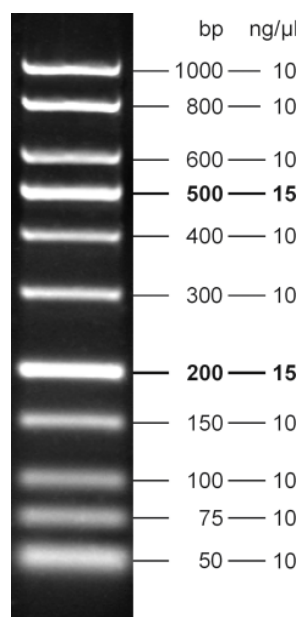
Low Range DNA Ladder consists of 11 DNA fragments ranging from 50 bp to 1 kb. It is designed to show virtually uniform spacing over a wide fragment range. The ladder allows sizing and concentration estimate of DNA fragments on agarose gels generated by PCR or restriction digest. The fragment mix is supplied in ready-to-use format containing orange and blue tracking dyes. It yields a ladder of the following double-stranded DNA fragments:

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

#### Recommended Load

5 µl DNA ladder per lane

#### Low Range DNA Ladder



3.0%  
Agarose