

## 1 kb DNA Ladder

### 1 kb to 10 kb log scale

ready-to-use

Cat. No.	Volume	Amount
M-217	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

#### Concentration

105 ng/µl

#### Recommended Load

5 µl DNA ladder per lane

#### DNA Fragments

1, 2, 3, 4, **5**, 6, 7, 8, 9, 10 bp

#### Form

DNA fragment mix in 10 mM Tris-HCl (pH 7.5),  
10 mM EDTA, 10% (w/v) Glycerol, Bromophenol Blue  
and Xylene Cyanol FF

#### Description

1 kb DNA Ladder consist of 10 DNA fragments ranging from 1 kb to 10 kb and is supplied in ready-to-use format containing blue tracking dyes. 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.

#### 1 kb DNA ladder

