

## Fluorescent 20 bp DNA Ladder

### 20 bp to 200 bp log scale

ready-to-use, orange, green fluorescent

Cat. No.	Volume	Amount
M-232	500 µl	100 lanes

For *in vitro* use only  
Quality guaranteed for 12 months  
Store at -20°C in the dark  
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

20, 40, 60, 80, **100**, 120, 140, 160, 180, 200 bp

#### Form

DNA fragment mix in 10 mM Tris-HCl (pH 7.5),  
10 mM EDTA, 10% (w/v) Glycerol, EvaGreen™,  
Orange G

#### Description

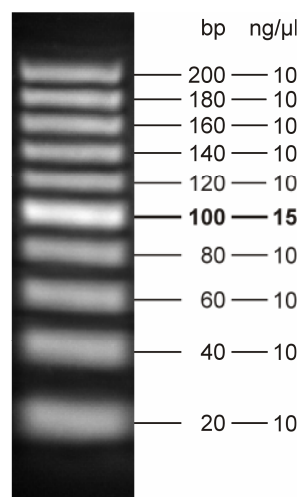
Fluorescent 20 bp DNA Ladder consists of 10 DNA fragments ranging from 20 bp to 200 bp. The prestained ladder is supplied in a ready-to-use format containing EvaGreen™ fluorescent DNA stain and an orange tracking dye. It allows sizing and concentration estimate of DNA fragments on agarose gels generated 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.

#### Fluorescent DNA Stain

DNA Ladders with Fluorescent Stain contain EvaGreen™ a fluorescent DNA intercalator dye specially developed for DNA analysis applications. High quantum yield and excellent stability makes EvaGreen™ the ideal fluorophore for DNA staining applications and a superior replacement for the widely used dyes Ethidium Bromide or SYBR® Green.

DNA Ladders with Fluorescent Stain are optimized for direct loading into unstained agarose gels and are recommended for use in combination with Gel Loading Buffer with DNA Stain (Cat.-No. PCR-255). The ladders provide highest convenience during routine handling and avoid commonly used gel staining procedures with Ethidium Bromide or SYBR® Green.

#### 20 bp DNA Ladder



3.6%  
Agarose