

Fluorescent 200 bp DNA Ladder

200 bp to 2 kb log scale

ready-to-use, blue / blue, green fluorescent

Cat. No.	Volume	Amount
M-235	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

200, 400, 600, 800, **1000**, 1200, 1400, 1600, 1800, 2000 bp

Form

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

Description

Fluorescent 200 bp DNA Ladder consist of 10 DNA fragments ranging from 200 bp to 2 kb. The prestained ladder is supplied in a ready-to-use format containing EvaGreen™ fluorescent DNA stain and blue tracking dyes. It allows sizing and concentration estimate of DNA fragments on agarose gels generated by PCR or restriction digest.

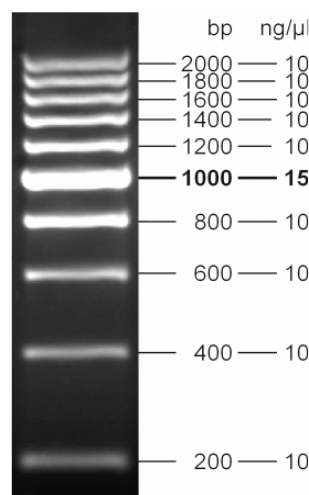
DNA log scale ladders are well suited for combination if small and large fragments should be analyzed on one gel.

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.

200 bp DNA Ladder



1.8%
 Agarose