

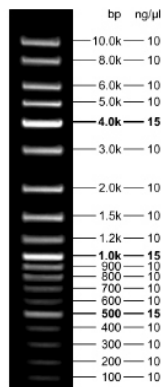


Fluorescent Long Range DNA Ladder

100 bp to 10 kb linear scale
ready-to-use, orange / blue, green fluorescent

Cat. No.	Amount
M-225S	500 µl (205 ng/µl)
M-225L	5 x 500 µl (205 ng/µl)

Long Range DNA Ladder



1% Agarose

5 µl loaded onto 1 % agarose

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Additional Storage Conditions: store dark

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, 10 % (w/v) Glycerol, SYBR® Orange G and Xylene Cyanol FF

Color: green

Concentration: 205 ng/µl

Description:

Fluorescent Long Range DNA Ladder consists of 19 DNA fragments ranging from 100 bp to 10 kb. It is designed to show virtually uniform spacing over an extreme wide fragment range covering virtually the complete application spectrum of DNA ladders. The ladder allows sizing and concentration estimate of DNA fragments on agarose gels generated by PCR or restriction digest. The prestained fragment mix is supplied in ready-to-use format containing SYBR® Green fluorescent DNA stain and orange / blue tracking dyes.

Fluorescent DNA Stain:

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

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 (#PCR-274 - #PCR-276).

The ladders provide highest convenience during routine handling and avoid commonly used gel stains like Ethidium Bromide.

Recommended Load:

5 µl per lane

DNA Fragments:

100, 200, 300, 400, **500**, 600, 700, 800, 900 bp,
1.0, 1.2, 1.5, 2.0, 3.0, **4.0**, 5.0, 6.0, 8.0, 10.0 kb