

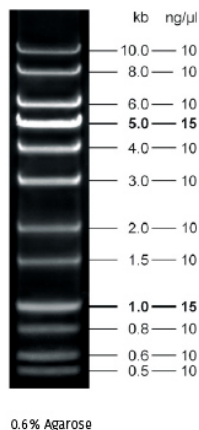


Fluorescent High Range DNA Ladder

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

Cat. No.	Amount
M-224S	500 µl (130 ng/µl)
M-224L	5 x 500 µl (130 ng/µl)

High Range DNA Ladder



0.6% Agarose

5 µl loaded onto 0.6 % 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, green-fluorescent DNA stain, Bromophenol Blue and Xylene Cyanol FF

Color: blue

Concentration: 130 ng/µl

Description:

Fluorescent High Range DNA Ladder consists of 12 DNA fragments ranging from 500 bp to 10 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 prestained fragment mix is supplied in ready-to-use format containing a green-fluorescent DNA stain and blue tracking dyes.

Fluorescent DNA Stain:

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

0.5, 0.6, 0.8, 1, 1.5, 2, 3, 4, 5, 6, 8, 10 kb