



High Molecular Weight DNA Markers

Cat. No. 15618-010

Conc.: 40 µg/ml

Size: 50 µg

Store at 4°C.

Description:

High Molecular Weight (MW) DNA Markers are prepared by digesting phage λ DNA with different restriction enzymes. They are suitable for sizing linear double-stranded DNA from 8.3 to 48.5 kb. These fragments can be visualized by ethidium bromide staining.

Storage Buffer:

10 mM Tris-HCl (pH 7.4)

50 mM NaCl

1 mM EDTA

Recommended Procedure:

Apply approximately 40 ng of markers per mm lane width. Heat to 65°C for no longer than 10 min. immediately before loading. Electrophoresis for 18 hours at 25 volts at room temperature results in optimal separation of high molecular weight DNA.

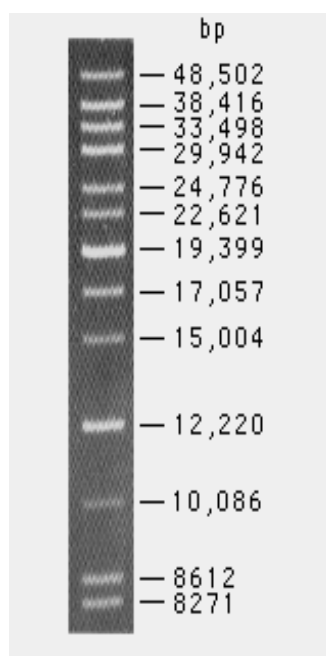
Note:

The migration and resolution of the DNA fragments will vary under different electrophoretic conditions. For the electrophoresis conditions outlined, DNA amounts of < 300 ng per well are recommended. Larger amounts of DNA will result in poor resolution. Because of this low DNA amount, it is necessary to take photographs with short wave (254 nm) transilluminators. Longer wave (302 or 366 nm) transilluminators show only faint bands. The top four DNA bands of the High Molecular Weight DNA Markers do not separate with shorter electrophoresis times.

Doc. Rev: 05/09/03

This product is distributed for laboratory research only. CAUTION: Not for diagnostic use. The safety and efficacy of this product in diagnostic or other clinical uses has not been established.

For technical questions about this product, call the Invitrogen Tech-LineSM U.S.A. 800 955 6288



Quality Control:

Upon electrophoresis of 100-300 ng of marker fragments under standard conditions and after ethidium bromide staining, all DNA bands are sharp and clearly separated.

Reference:

Uher, L. (1986) *FOCUS*® 8:1, 10.

High Molecular Weight DNA Markers
0.2 µg/lane
0.4% agarose gel
stained with ethidium bromide

Cat. No. 15618-010