



Ready-Load™ 100 bp DNA Ladder

Cat. No. 10380-012

Size: 100 applications

Conc. 0.1 µg/µl

Store at Room Temperature or +4°C

Description

Ready-Load™ 100 bp DNA Ladder is intended for use as a molecular weight standard in agarose or polyacrylamide gel electrophoresis. The ladder is supplied in gel loading buffer and can be applied directly to an agarose or polyacrylamide gel. These fragments can be visualized by ethidium bromide staining.

Ready-Load™ 100 bp DNA Ladder consists of 15 blunt-ended fragments between 100 and 1500 bp in multiples of 100 bp and an additional fragment at 2072 bp. The 600 bp band is approximately 2 to 3 times brighter than the other ladder bands to provide internal orientation. This ladder is not designed for quantitation.

Storage Buffer

10 mM Tris-HCl (pH 7.5)

10 mM EDTA

0.05% bromophenol blue

5% glycerol

Recommended Procedure

If product is stored at room temperature, briefly centrifuge prior to use. Apply 1 µl of ladder per mm lane width. **Do not** heat before loading.

Quality Control

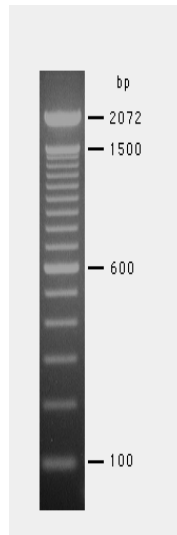
Agarose gel analysis shows that the bands between 100 to 1500 bp are distinguishable. The 600 bp band must be more intense than any other band except the band at 2072 bp.

Part no. 10380012.pps

Rev. Date: 08/14/03

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-Line® U.S.A. 800 955 6288



Note

During 2% agarose gel electrophoresis with tris-acetate (pH 7.6) as the running buffer, bromophenol blue migrates near the 100 bp fragment. The 100 bp band migrates behind the bromophenol blue marker on 6% polyacrylamide gels with tris-borate (pH 8.0) as the running buffer.

Part of the 600 bp band may migrate anomalously slowly in polyacrylamide gels (1,2). This band may appear as an extra band near or on top of the 700 bp band.

References

1. Hsieh, C., et al. (1991) *Mol. Gen. Genet.* 225, 25.
2. Stellwagen, N.C. (1983) *Biochemistry* 22, 6186.

100 bp DNA Ladder
0.5 µg/lane
2% agarose gel
stained with ethidium bromide