



508PR-01

G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

DNAMark™ 100bp Ladder

(Cat. # 786-855)



think proteins! think G-Biosciences www.GBiosciences.com

INTRODUCTION

DNAmark™ 100bp Ladder consists of 11 fragments ranging from 100-1500 base pairs (bp). The 500bp fragment is present at increased intensity to allow easy identification. All fragments are precisely quantified and mixed during the production. For 5µl loading, all fragments except 500bp are 40ng. The 500bp fragment is 100ng. This ladder is pre-mixed with loading dye and is ready to use.

ITEMS SUPPLIED (Cat. # 786-855)

Description	Size
DNAmark™ 100bp Ladder	50µg

STORAGE CONDITION

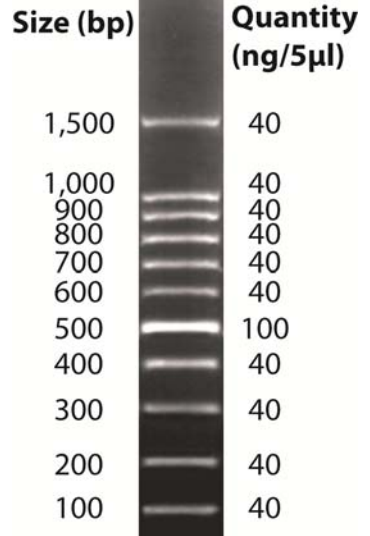
Upon arrival, store at -20°C.

CONCENTRATION

100ng/µl

LOADING BUFFER COMPOSITION

10mM Tris-HCL (pH7.6)
0.03% Bromophenol Blue
0.03% Xylene Cyanol
60% Glycerol
60mM EDTA



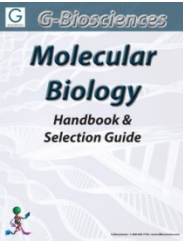
DNAmark™ 100bp Ladder resolved on a 1.7% Agarose-TBE gel and stained with LabSafe™ Nucleic Acid Stain

PROTOCOL

Load 2-5µl DNAmark™ 100bp Ladder directly to wells designated for markers. The volume loaded is dependent on the well size, gel thickness and concentration and level of dye used to visualize the bands.

RELATED PRODUCTS

Download our Molecular Biology Handbook.



<http://info.gbiosciences.com/complete-molecular-biology-handbook>

For other related products, visit our website at www.GBiosciences.com or contact us.

Last saved: 6/12/2013 CMH



www.GBiosciences.com