



538PR-01

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

A Geno Technology, Inc. (USA) brand name

DNAmark™ 50bp Plus Ladder

(Cat. # 786-853)



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

INTRODUCTION

DNAmark™ 50bp Plus Ladder consists of 13 fragments ranging from 50-1000 base pairs (bp). This molecular weight marker is specially designed for identifying small DNA fragments and allowing for quantitative estimation of the DNA fragments in the gel at the same time. The 250bp and 500bp fragment are present at increased intensity to allow easy identification. All fragments are precisely quantified and mixed during the production. For 5µl loading, all fragments, except 250bp and 500bp, are 40ng. The 250bp and 500bp fragment are 100ng. This ladder is pre-mixed with loading dye and is ready to use.

ITEMS SUPPLIED (Cat. # 786-853)

Description	Size
DNAmark™ 50bp Plus Ladder	50µg

STORAGE CONDITION

Upon arrival, store at -20°C.

CONCENTRATION

128ng/µl

LOADING BUFFER COMPOSITION

10mM Tris-HCl

1mM EDTA (pH 8.0)

0.02% Bromophenol blue

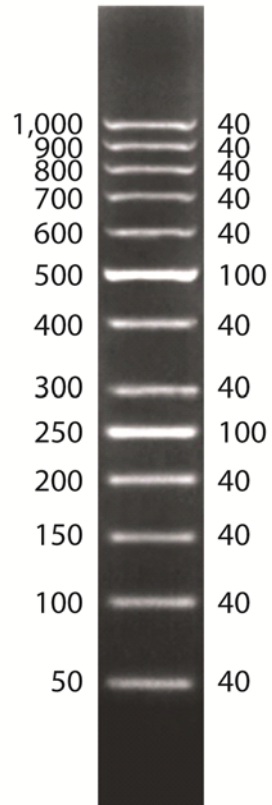
0.02% Xylene cyanol

5% Glycerol

PROTOCOL

Load 2-5µl DNAmark™ 50bp Plus 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

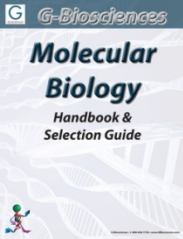
Size (bp) Quantity (ng)



5µl DNAmark™ 50bp Plus Ladder resolved on a 3% Agarose-TAE gel and stained with LabSafe™ Nucleic Acid Stain

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: 6/11/2013 CMH



www.GBiosciences.com