



510PR-01

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

A Geno Technology, Inc. (USA) brand name

DNAmark™ 1kb Plus

(Cat. # 786-854)



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

INTRODUCTION

DNAMark™ 1kb Plus contains 15 DNA bands ranging from 100bp to 10Kb. The 500bp and 3,000bp 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 500bp and 3,000bp are 40ng. The 500bp and 3,000bp fragment are 100ng. This ladder is pre-mixed with loading dye and is ready to use.

ITEM(S) SUPPLIED (Cat. # 786-854)

Description	Size
DNAMark™ 1kb Plus	50µg

STORAGE CONDITION

Shipped on blue ice, upon arrival, store at -20°C.

CONCENTRATION

144ng/µl

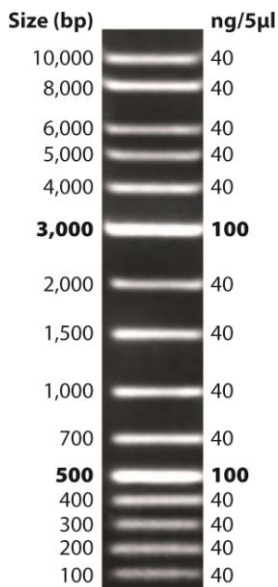
LOADING BUFFER COMPOSITION

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

PROTOCOL

Load 2-5µl DNAMark™ 1kb Plus Ladder directly to wells designated for markers.

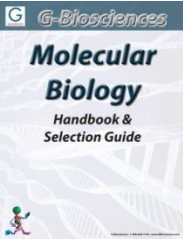
The volume loaded is dependent on the well size, gel thickness and concentration and level of dye used to visualize the bands.



DNAMark™ 1kb Plus Ladder resolved on a 1% 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: 4/30/2014 CMH



www.GBiosciences.com