



006PR-03

**G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)**

A Geno Technology, Inc. (USA) brand name

# DNAmark™ 500bp Ladder

(Cat. # 786-462)



**think proteins! think G-Biosciences [www.GBiosciences.com](http://www.GBiosciences.com)**

## INTRODUCTION

The DNAmark™ 500bp ladder is a ready-to-use preparation that consists of 10 DNA fragments ranging from 500bp to 5.0kb. The 2,000bp 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 2,000bp are 40ng. The 2,000bp fragment is 100ng. This ladder is pre-mixed with loading dye and is ready to use.

## ITEM(S) SUPPLIED

Cat. #	Description	Size
786-462	DNAmark™ 500bp ladder	50µg

## STORAGE CONDITION

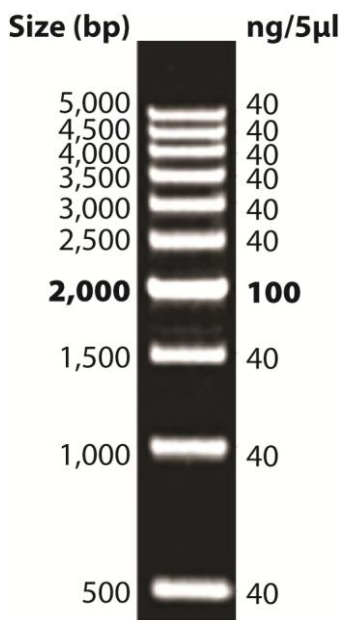
Shipped at ambient temperature. Upon arrival, store at -20°C.

## CONCENTRATION

92ng/µl

## LOADING BUFFER COMPOSITION

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



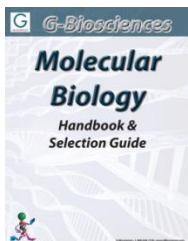
## PROTOCOL

Load 2-5µl DNAmark™ 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.

*DNAmark™ 500bp 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](http://www.GBiosciences.com) or contact us.

Last saved: 4/30/2014 SM



**[www.GBiosciences.com](http://www.GBiosciences.com)**