



245PR

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

A Geno Technology, Inc. (USA) brand name

LongLife™ DNase

(Cat. # 786-041)



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

INTRODUCTION

LongLife™ DNase is a ready to use DNase preparation provided with a co-factor bivalent metal ion. LongLife™ DNase is an endonuclease, which splits phosphodiester linkages, yielding 5'- phosphate terminated polynucleotides with a free hydroxyl group on position 3'. LongLife™ DNase is supplied in 0.5ml volume (10 Units/μl) in a buffer that ensures its full activity and is RNase free. Simple to use, take an aliquot and add in your reaction sample.

ITEM(S) SUPPLIED (Cat. # 786-041)

Description	Size
LongLife™ DNase (10 Units/μl)	1 x 0.5ml

STORAGE CONDITIONS

It is shipped on blue ice. Upon arrival, store at -20°C.

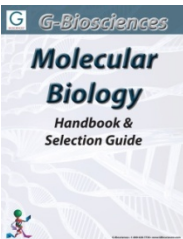
INSTRUCTION FOR USE

NOTE: LongLife™ DNase is supplied as **suspension** 10 Units/μl.

1. Flick the tube briefly before taking out an aliquot.
2. Add appropriate volume of LongLife™ DNase in the reaction mix (e.g. 10μl/ml)
3. Vortex briefly and incubate at 25°C for 30-45 minutes or as needed.
4. Store the unused LongLife™ DNase at -20°C.

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: 11/4/2013 CMH

This page is intentionally left blank



www.GBiosciences.com