



249PR

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

A Geno Technology, Inc. (USA) brand name

LongLife™ Nuclease

Ready To Use, 100X Concentration

(Cat. # 786-039)



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

INTRODUCTION

LongLife™ Nuclease is a cocktail preparation containing DNase and RNase as well as co-factors necessary for optimal activity. The LongLife™ Nuclease is ready-to-use solution that has a long shelf life (> 1 year). Simply add directly to your sample for high efficiency removal of RNA and DNA. LongLife™ Nuclease can be used in conjunction with protease inhibitor cocktails that do not contain EDTA, a nuclease inhibitor, e.g. FOCUS ProteaseArrest™.

A 1X LongLife™ Nuclease solution contains 80 units DNase and 1.2 Units RNase from bovine pancreas and 0.5mM cofactors.

ITEM(S) SUPPLIED (Cat. # 786-039)

Description	Size
LongLife™ Nuclease [100X]	0.5ml

STORAGE CONDITIONS

Shipped at ambient temperature. Store at -20°C upon arrival.

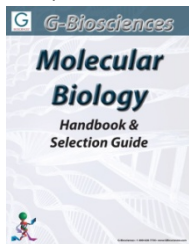
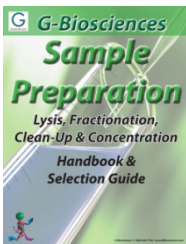
INSTRUCTIONS FOR USE

1. LongLife™ Nuclease is supplied as a 100X suspension, vortex before use.
2. Use 10µl LongLife™ Nuclease for every 1ml reaction.
3. Briefly, vortex the reaction and incubate at room temperature for 30-45 minutes.

NOTE: Protease inhibitors may be added in the same reaction mix.

RELATED PRODUCTS

Download our Sample Preparation or Molecular Handbooks.



<http://info.gbiosciences.com/complete-protein-sample-preparation-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: 2/22/2013 CMH

This page is intentionally left blank



www.GBiosciences.com