



253PR

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

A Geno Technology, Inc. (USA) brand name

# LongLife™ RNase

(Cat. # 786-040)



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

## INTRODUCTION

LongLife™ RNase is a ready to use endo-ribonuclease preparation, which efficiently hydrolyzes RNA contaminants, present in the sample preparation. The enzyme preparation is carefully optimized for use in molecular biology research, as it is free from DNase activity. The LongLife™ RNase is supplied in a buffer that ensures its full activity, while in storage. It is simple to use, just take an aliquot and add in your sample.

## ITEM(S) SUPPLIED (Cat. # 786-040)

Description	Size
LongLife™ RNase (5mg/ml; >60U/mg)	2 x 0.5ml

## STORAGE CONDITIONS

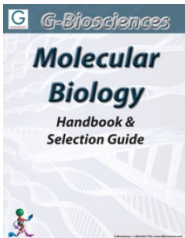
Shipped at ambient temperature. Store at -20°C upon arrival.

## INSTRUCTION FOR USE

1. LongLife™ RNase is supplied as **suspension** in a concentration of 5mg/ml (>60U/mg).
2. Thaw the tube and flick it briefly to mix the tube contents before taking an aliquot.
3. Add appropriate volume of LongLife™ RNase to the reaction mixture (e.g. 10µl/ml or as needed),
4. Vortex it briefly and incubate it at 25°C for 15-30 minutes or as needed.
5. Store the LongLife™ RNase tube at -20°C after its use.

## 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: 5/19/2015 CMH

*This page is intentionally left blank*



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