



247PR

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

A Geno Technology, Inc. (USA) brand name

LongLife™ Lysozyme

(Cat. # 786-037)



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

INTRODUCTION

A ready to use lysozyme preparation is provided, which can be used for spheroplast preparation, lysis of bacterial cells and further extraction of protein and DNA from bacterial cells. Simply take an aliquot and add to your reaction. The total volume of the supplied LongLife™ Lysozyme is 1ml (2 x 0.5ml) in a concentration of 1,500 Units/μl.

APPLICATIONS:

Spheroplast preparation, lysis and extraction of protein and DNA from bacterial cells.

ITEM(S) SUPPLIED (Cat. # 786-037)

Description	Size
LongLife™ Lysozyme (1,500 units/μl)	1ml

STORAGE CONDITIONS

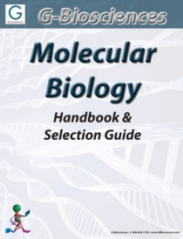
It is shipped at ambient temperature. Store at -20°C, upon arrival.

PROTOCOL

1. Suspend bacterial cell pellet in an appropriate suspension buffer for efficient lysis. (*G-Biosciences' Bacterial Suspension Buffer is recommended*).
2. Vortex for approximately 1 minute until the cell suspension becomes homogenous. Incubate the suspension for 5 minutes in cold and vortex again to suspend the cells.
3. Flick the vial containing lysozyme to mix the solution. Add 10μl of Lysozyme per 100μl of bacterial cell pellet suspended in suspension buffer. Gently mix the content.
4. Incubate the cell suspension at 37°C for 30-60 minutes.
5. At the end of incubation, centrifuge the bacterial cell suspension at 10,000g for 5 minutes and remove the supernatant.
6. Lysis: Suspend the spheroplast obtained in an appropriate lysis buffer (*G-Biosciences' Bacterial-PE LB™ recommended if extraction of protein is required*) two to three times the volume added in step 1.
7. Pipette the suspension up and down a few times then vortex and incubate on ice for 30 minutes. The lysis may be further facilitated by incubating the cells at 37°C for 1-3 minutes.

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: 8/7/2012 CMH



www.GBiosciences.com