



255PR-01

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

A Geno Technology, Inc. (USA) brand name

Longlife™ Zymolyase®

(Cat. # 786-036, 786-914)



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

INTRODUCTION

A ready to use Zymolyase[®] preparation is provided, which can be used for spheroplast preparation, yeast plasmid isolation / lysis of yeast cells and further extraction of soluble protein from them. The size of the supplied Longlife[™] Zymolyase[®] is 1ml (2 x 0.5ml) in a concentration of 1500Units/ml.

APPLICATIONS

Spheroplast preparation / plasmid isolation, lysis and extraction of proteins from yeast cells

ITEM(S) SUPPLIED

Description	Cat. # 786-036	Cat. # 786-914
Longlife [™] Zymolyase [®]	2 x 0.5ml [1.5Units*/μl]	5ml [1.5Units*/μl]

*One unit will produce a change of 0.001 OD₈₀₀ per minute at 25°C, using a yeast suspension as substrate in 3ml reaction mixture.

STORAGE CONDITIONS

It is shipped at ambient temperature. Upon arrival, store Longlife[™] Zymolyase[®] at -20°C.

ADDITIONAL ITEMS REQUIRED

- β-mercaptoethanol

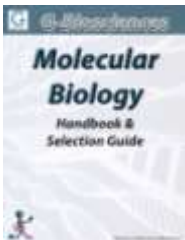
PROTOCOL

1. Suspend yeast cell pellet in an appropriate suspension buffer for efficient lysis. (G-Biosciences' Yeast Suspension Buffer is recommended). Add 1 μ l of β -mercaptoethanol per 100 μ l of yeast suspension buffer.
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 Zymolyase[®] to mix the solution. Add 5 μ l of Zymolyase[®] per 50 μ l of yeast cell pellet suspended in yeast 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 yeast cell suspension at 10,000g for 5 minutes and remove the supernatant.
6. Lysis: Suspend the yeast cell pellet obtained, in an appropriate lysis buffer (G-Bioscience's Yeast-PE LB[™] recommended) two to three times the volume added in step 1. 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.
7. Centrifuge at 1000g for 1h at 4°C and collect clear lysate. The lysate is now ready for protein purification and further analyses.

Zymolyase[®] is a registered trademark of Kirin Brewery Co. Ltd.

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.



www.GBiosciences.com