



G-Biosciences, St Louis, MO. USA ♦ 1-800-628-7730 ♦ 1-314- 991-6034 ♦ technical@GBiosciences.com

Certificate of Analysis

Product: *Longlife Zymolyase*

Catalog #: 786-036

Size: 2 x 0.5ml

Lot#: 180507

Test:

Zymolyase Enzyme activity – OD change/minute Passed Yes / ~~No~~
(Test Sample: Yeast – *S. cerevisiae*)

Storage: Store at –20°C.

Note: *Warm it to room temperature before opening the vial.*

This document has been produced electronically and is valid without signature

Date of Print: July 13, 2018

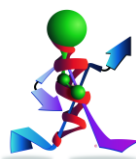
G-Biosciences / Genotech

9800 Page Avenue

St. Louis, MO 63132, USA

Phone: 314-991-6034, Fax: 314-991-1504

Longlife Zymolyase C of A/MN



think proteins! think G-Biosciences!

www.GBiosciences.com