



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

## Certificate of Analysis

*For Research Use Only*

**Product:** Zymogram Sample Buffer

**Catalog #:** 786-483

**Lot #:** 183003

**Re-test date:** 03/2023

### Tests:

Specific gravity test	Passed	Yes / <del>No</del>
pH test	Passed	Yes / <del>No</del>

**Storage:** Store at room temperature

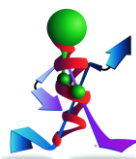
This document has been produced electronically and is valid without signature

**Date of Print:** June 1, 2012

### **G-Biosciences / Genotech**

9800 Page Ave  
St. Louis, MO 63132, USA  
Ph: 314-991-6034, Fax: 314-991-1504

Zymogram Sample Buffer Cat# 786-483 CofA/Com#3, Rev 12.24.08-SA/MM/IA



think proteins! think G-Biosciences!

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