



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

## Certificate of Analysis

**Product:** Tris [1M] solution, pH 8.0

**Catalog #:** R002

**Lot #:** 191712

**Re-test date:** 12/2024

### Tests:

Appearance: Clear Liquid      Passed      Yes / ~~No~~

Specific gravity test      Passed      Yes / ~~No~~

pH Test      Passed      Yes / ~~No~~

**Storage:** Store at RT      *for Research Use Only*

This document has been produced electronically and is valid without signature

**Date of Print:** February 25, 2020

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

9800 Page Avenue

St. Louis, MO 63132, USA

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

Tris 1M, pH 8.0 C of A/01.31.07/Com#3

Rev 12.23.08-SA/MM/IA



think proteins! think G-Biosciences!

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