



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

Certificate of Analysis

Product: Tris [1M] solution, pH 8.0

Catalog #: R002

Lot #: 201603

Re-test date: 03/2025

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*

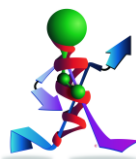
This document has been produced electronically and is valid without signature

Date of Print: July 30, 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