



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 7.5

Catalog #: 786-265

Lot #: 190504

Re-test date: 04-2024

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: December 17, 2019

G-Biosciences /Genotech

9800 Page Avenue, St. Louis, MO 63132, USA

Fax: 314-991-1504

Tris 1M , pH 7.5 C of A/Com#3
Rev 05.30.12/MM



think proteins! think G-Biosciences!

www.GBiosciences.com