



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 7.0

Catalog #: R001

Size: 100ml

Lot #: 131211

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*

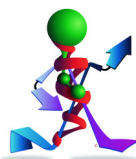
This document has been produced electronically and is valid without signature

Date of Print: November 14, 2013

G-Biosciences / Genotech

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

Tris 1M, pH 7.0 C of A/Com#3
Rev 04.26.10 /IA



think proteins! think G-Biosciences!

www.GBiosciences.com