



G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: Tris, pH8.0 [1M]

Catalog No: 786-815

Release Date: 11/15/2021

Size: 500ml

Retest Date: 11/4/2023

Lot No: 210411

Storage: Room temp

Specifications

Test	Specification	Result
Appearance	Clear Liquid	Passed
Sterility Test Passed	Yes / No	Yes
pH Test Passed	Yes / No	Yes

This document has been produced electronically and is valid without signature

Date of Print: January 28, 2022

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP



think proteins! think G-Biosciences www.GBiosciences.com