



G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: Tris, pH 8.5 [1M]

Catalog No: 786-269

Release Date: 12/29/20

Size: 100ml

Retest Date: 12/29/24

Lot No: 202912

Storage: Room Temp.

Specifications

CAS Number: N/A

Appearance: Clear, colorless liquid

Test	Specification	Result
pH	8.45 – 8.55	8.52

This document has been produced electronically and is valid without signature

Date of Print: January 27, 2021

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP



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