



G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: Tris-HCl [1M], pH 9.0

Catalog No: 786-476

Release Date: 12/17/19

Size: 100ml

Retest Date: 12/17/22

Lot No: 191712

Storage: Room temp.

Specifications

Appearance: Clear, colorless liquid

Test	Specification	Result
pH @ RT	8.90 – 9.10	9.00

This document has been produced electronically and is valid without signature

Date of Print: December 24, 2019

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP



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