



**G-Biosciences** ♦

♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

A Geno Technology, Inc. (USA) brand name

## Certificate of Analysis / Data Sheet

**Product:** Tris, pH 7.5 [1M]

**Catalog No:** 786-265

**Release Date:** 1/29/20

**Size:** 100ml

**Retest Date:** 1/29/23

**Lot No:** 202901

**Storage:** Room Temp.

### Specifications

**CAS Number:** N/A

**Appearance:** Clear, colorless liquid

Test	Specification	Result
pH	7.45-7.55	7.53

This document has been produced electronically and is valid without signature

**Date of Print:** February 5, 2020

**G-Biosciences/Geno Technology, Inc.**

Rev 10/25/18 BP



**think proteins! think G-Biosciences** [www.GBiosciences.com](http://www.GBiosciences.com)