



G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: TBE Buffer [10X] (890mM Tris, 890mM Boric acid, 20mM EDTA)

Catalog No: R025

Release Date: 10/19/20

Size: 1L

Retest Date: 10/19/22

Lot No: 201910

Storage: Room Temp.

Specifications

CAS Number: N/A

Appearance: Clear, colorless liquid. If precipitation occurs in solution, microwave for 3-5 min.

Test	Specification	Result
pH	8.20 – 8.40	8.33

This document has been produced electronically and is valid without signature

Date of Print: October 28, 2020

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP



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