



G-Biosciences, St Louis, MO. USA ♦ 1-800-628-7730 ♦ 1-314- 991-6034 ♦ technical@GBiosciences.com

Certificate of Analysis

Product: TBE [10X] (0.9M Tris-borate, 20mM EDTA, pH8.3))

Catalog #: R025

Lot #: 210406

Re-test date: 06/2026

Size: 1L

Tests:

Appearance - Clear Liquid	Passed	Yes / No
Specific gravity test	Passed	Yes / No
pH test	Passed	Yes / No

Storage: Store at room temperature

Note: *If precipitate develops after a long-term storage, microwave the solution bottle for 3-5 minutes. Let it cool down then use. This wouldn't affect the product quality and functioning.*

This document has been produced electronically and is valid without signature

For Research Use Only

Date of Print: December 21, 2021

G-Biosciences / Genotech

9800 Page Avenue
St. Louis, MO 63132, USA
Ph: 314-991-6034, Fax: 314-991-1504

10X TBE Cat# R025 C of A/Com#3Rev 07/10/2014/IA



think proteins! think G-Biosciences!

www.GBiosciences.com