



Certificate of Analysis

Product: TBST [10X]

Catalog #: R043

Size: 1 gallon

Lot #: 200212

Re-test date: 12/2023

Tests:

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

Storage: Store at room temperature

For Research Use Only

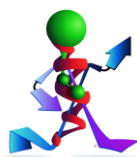
This document has been produced electronically and is valid without signature

Date of Print: December 21, 2020

G-Biosciences / Genotech

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

10X TBS Cat# R029 C of A/Com#3
Rev 09.21.12/MM



think proteins! think G-Biosciences!

www.GBiosciences.com