



## Certificate of Analysis

**Product:** TBST [10X]

**Catalog #:** R042

**Size:** 1L

**Lot #:** 200212

**Re-test date:** 12/2023

### Tests:

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

**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

