



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

A Geno Technology, Inc. (USA) brand name

CERTIFICATE OF ANALYSIS

Product Name: [10X] femto™ TBST

Cat #: 786-161

Lot #: 201712

Re-test date: 12/2024

Size: 250ml

Test Procedure:

Each batch of [10X] femto™ TBST is checked for correct molar specification of the buffer components and the pH.

Specification for 10X femto™ TBST to pass:

For [10X] femto™ TBST to pass, it must have correct molar specification of the buffer components and the pH.

Molar Specification	Passed	Yes / No
pH Specification	Passed	Yes / No

Storage: Store at 4 degrees C.

This document has been produced electronically and is valid without a signature

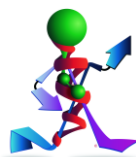
Date of print: December 29, 2020

G-Biosciences / Genotech

St. Louis. MO. USA

www.GBiosciences.com

10X femto TBST 786-161/C of A, Rev 01.27.09-SA/MM/IA



think proteins! think G-Biosciences!

www.GBiosciences.com