



**G-Biosciences** ♦

♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

A Geno Technology, Inc. (USA) brand name

## Certificate of Analysis / Data Sheet

**Product:** TE Buffer [100X]

**Catalog No:** R003

**Release Date:** 8/7/20

**Size:** 100mL

**Retest Date:** 8/7/32

**Lot No:** 200708

**Storage:** Room Temp.

### Specifications

**CAS Number:** N/A

**Appearance:** Clear, colorless liquid

Test	Specification	Result
DNase Activity	None Detected	PASS
RNase Activity	None Detected	PASS
pH (1X solution)	7.96 – 8.04	8.04

This document has been produced electronically and is valid without signature

**Date of Print:** September 2, 2020

**G-Biosciences/Geno Technology, Inc.**

Rev 10/25/18 BP



**think proteins! think G-Biosciences** [www.GBiosciences.com](http://www.GBiosciences.com)