



**G-Biosciences, St Louis, MO. USA ♦ 1-800-628-7730 ♦ 1-314- 991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)**

A Geno Technology, Inc. (USA) brand name

### **CERTIFICATE OF ANALYSIS**

**Product Name:** 50X TAE

**Size:** 500ml

**Cat #:** 786-060

**Lot #:** 202910

**Re-test date:** 10/2023

**Test Procedure:**

Each batch of 50X TAE is checked for correct molar specification of the buffer components and the pH.

**Specification for 50X TAE Buffer to pass:**

It must have correct molar specification of the buffer components and the pH.

Molar Specification	Passed	Yes / <del>No</del>
pH Specification	Passed	Yes / <del>No</del>

**Storage:** Store at room temp.

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

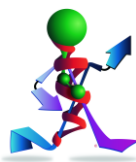
**Date of print:** November 6, 2020

**G-Biosciences / Genotech**

9800 Page Avenue, St. Louis, MO 63132, USA

Ph: 314-991-6034, Fax: 314-991-1504

50X TAE 786-060 C of A  
Rev 04/26/2011/CMH



think proteins! think G-Biosciences!

[www.GBiosciences.com](http://www.GBiosciences.com)

