



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: 50X TAE

Size: 1 gallon

Cat #: R024

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 / No
pH Specification	Passed	Yes / No

Storage: Store at room temp.

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

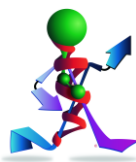
Date of print: November 6, 2020

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

