



A Geno Technology, Inc. (USA) brand name

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

CERTIFICATE OF ANALYSIS

Product Name: 50X TAE

Size: 1L

Cat #: R023

Lot #: 212607

Re-test date: 07/2026

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: December 21, 2021

G-Biosciences / Genotech

St. Louis. MO. USA

www.GBiosciences.com

50X TAE C R023 of A/Common3
Rev 01.05.10/IA



think proteins! think G-Biosciences!

www.GBiosciences.com