



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

Certificate of Analysis

Product: Trichloroacetic acid (TC) Solution, 10% w/v

Catalog #: 786-886

Size: 500ml

Lot#: 210303

Re-test date: 03/2025

Tests:

Appearance – Clear colorless liquid	Passed	Yes / No
Specific gravity	Passed	Yes / No
pH @ 25°C	Passed	Yes / No

Storage: Store at room temperature.

This document has been produced electronically and is valid without signature

Date of Print: January 14, 2022

G-Biosciences/ Genotech

9800 Page Avenue
St. Louis, MO 63132, USA
Ph: 314-991-6034, Fax: 314-991-1504

Trichloroacetic acid (TC) Solution, 10% w/v C of A/Com#3/Rev 04.29.16/IA



think proteins! think G-Biosciences!

www.GBiosciences.com