



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

Certificate of Analysis

For Research Use Only

Product: TBE Urea Sample Buffer

Catalog #: 786-474

Lot #: 212306

Tests:

Appearance – Blue Liquid	Passed	Yes / No
Specific gravity test	Passed	Yes / No
pH test	Passed	Yes / No

Storage: Store at room temperature

This document has been produced electronically and is valid without signature

Date of Print: October 12, 2021

G-Biosciences / Genotech

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

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

TBE-Urea Sample Buffer Cat# 786-474 C of A/Com#3/Rev 10.12.12/MM



think proteins! think G-Biosciences!

www.GBiosciences.com