



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

Certificate of Analysis

Product: DetergentOUT GBS-10-3000

Catalog #: 786-156

Size: Kit

Lot#: 200804B

Re-test date: 04/2025

Analysis

Passes

This kit performs as per corresponding and/or companion protocol

Yes/~~No~~

This document has been produced electronically and is valid without signature

Date of Print: January 20, 2022

For Research Use Only

G-Biosciences / Genotech

9800 Page Avenue

St. Louis, MO 63132, USA

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



think proteins! think G-Biosciences!

www.GBiosciences.com