



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

Certificate of Analysis

Product: RBC Lysis Buffer

Catalog No.: 786-649

SIZE: 100ml

LOT NO.: 190309

Re-test date: 09/2024

SPECIFICATIONS:

Appearance: Liquid Clear liquid

Specific Gravity: Passed

pH test Passed

STORAGE: Ambient

SHIPPING: Material shall be packaged and shipped in a manner that will insure protection of the material from the adverse effects of moisture, dirt and other forms of contamination or the adverse effects of temperature.

This document has been produced electronically and is valid without signature

Date of Print: August 7, 2020

G-Biosciences / Genotech

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

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

Last saved: 8/7/2020 IA



think proteins! think G-Biosciences!

www.GBiosciences.com