



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

Certificate of Analysis

Product: IgG elution buffer

Catalog #: 786-545

Size: 100 ml

Lot #: 202011

Re-test date: 11/2026

QC/QA Tests:

Appearance: Clear Liquid Passed Yes / ~~No~~

Specific gravity Test Passed Yes / ~~No~~

pH Test [2.5-3.5] Passed Yes / ~~No~~

Storage: Store at RT.

This document has been produced electronically and is valid without signature

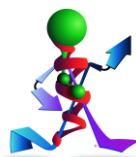
For Research Use Only

Date of Print: December 2, 2020

Genotech / G-Biosciences

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

Rev 04.19.13/IA



think proteins! think G-Biosciences!

www.GBiosciences.com