



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

Certificate of Analysis

Product: Immobilized EndotoxinOUT resin
(50% slurry in 20% ethanol)

Catalog #: 786-367

Size: 10 ml Resin

Lot#: 212709

Expiry: 09/2026

Test(s)	Specifications	Result
Appearance	White, free of flocculent material	Passed
Capacity	≥9995 endotoxin units (EU) removed by 1ml resin	Passed
Storage:	Store at 4°C. DO NOT FREEZE.	

This document has been produced electronically and is valid without signature

Date of Print: November 4, 2021

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