



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

CERTIFICATE OF ANALYSIS

Product: LabSafe Gel Blue

Catalog #: 786-35G

Size: 1 Gallon

Lot Number: 212101

Re-test date: 01/2026

TESTS:

QC GEL STAINING-

SDS-PAGE

Gradient 15 well Gel 4-20%, SDS-Tris-Glycine

Protein BSA

BSA Loaded/Well:

1µg, 0.5µg, 0.25µg, 0.125µg, 0.062µg, 0.031µg, 0.015µg(15ng), 0.0078µg(8ng), 0.0039µg(4ng)

BAND VISIBILITY TEST:

1µg - 0.031µg bands visible within 2-3 minutes OK Yes / ~~No~~

1µg - 0.015µg bands visible within 5 minutes OK Yes / ~~No~~

8ng - 4ng bands visible within 30 minutes OK Yes / ~~No~~

Band visibility after washing 1µg - 8ng OK Yes / ~~No~~

Storage: Store at Room Temp

This document has been produced electronically and is valid without a signature

Date of print: February 11, 2021

9800 Page Avenue

St. Louis, MO 63132, USA

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

LabSafe Gel Blue C of A/Rev 06.20.12/MM/Common3



think proteins! think G-Biosciences!

www.GBiosciences.com