



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

Certificate of Analysis

Product: SDS-PAGE Sample Loading Buffer [6X]

Catalog #: 786-701

Size: 25 ml

Lot #: 212312

Re-test date: 12/2025

Tests:

| | | |
|--------------------------------|--------|---------------------|
| Appearance – Blue color liquid | Passed | Yes / No |
| BSA Standard Protein Gel | Passed | Yes / No |
| pH Test | Passed | Yes/ No |

Storage: Store at room temp.

This document has been produced electronically and is valid without signature

Date of Print: January 13, 2022

G-Biosciences / Genotech

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

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

SDS-PAGE Sample Loading Buffer 6X, 786-701 C of A/Com#3
Rev 08.22.14/CM



think proteins! think G-Biosciences!

www.GBiosciences.com