



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

Certificate of Analysis

Product: Sodium azide Solution, 1%

Catalog #: 786-750

Size: 500ml

Lot #: 201711

Re-test date: 11/2025

QC Tests:

Appearance - Clear Liquid	Passed	Yes / No
Specific gravity test	Passed	Yes / No
pH Test	Passed	Yes / No

Storage: Store at room temperature

For Research Use Only

This document has been produced electronically and is valid without signature

Date of Print: November 20, 2020

G-Biosciences / Genotech

9800 Page Avenue
St. Louis, MO 63132, USA
Ph: 314-991-6034, Fax: 314-991-1504

Sodium azide 1% Sol C of A/Com#3
Rev 08.25.16/MM



think proteins! think G-Biosciences!

www.GBiosciences.com