



G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: Sodium Azide Solution [1%]

Catalog No: 786-750

Release Date: 02/04/21

Size: 500ml

Retest Date: 02/04/26

Lot No: 210402

Storage: Room Temp.

Specifications

CAS Number: N/A

Appearance: Clear, colorless liquid

| Test | Specification | Result |
|------|---------------|--------|
| pH | 9.30 to 9.90 | 9.83 |

This document has been produced electronically and is valid without signature

Date of Print: February 11, 2021

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP



think proteins! think G-Biosciences www.GBiosciences.com