



G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: Sodium Azide, 1%

Catalog No: 786-750

Size: 500ml

Lot No: 191411

Release Date: 11/14/19

Retest Date: 11/14/24

Storage: Room temp.

Specifications

Appearance: Clear, colorless liquid

Test	Specification	Result
pH	9.3-9.7	9.34

This document has been produced electronically and is valid without signature

Date of Print: December 2, 2019

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP



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