



**G-Biosciences** ♦

♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

A Geno Technology, Inc. (USA) brand name

## Certificate of Analysis / Data Sheet

**Product:** Sodium Azide, 1%

**Catalog No:** 786-299

**Size:** 100ml

**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](http://www.GBiosciences.com)