



**G-Biosciences** ♦

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

A Geno Technology, Inc. (USA) brand name

## Certificate of Analysis / Data Sheet

**Product:** SDS [20%]

**Catalog No:** 786-017

**Release Date:** 10/27/20

**Size:** 1 L

**Retest Date:** 10/27/24

**Lot No:** 202710

**Storage:** Room Temp.

### Specifications

**CAS Number:** N/A

**Appearance:** Clear, colorless liquid. Some white precipitant may form.

Test	Specification	Result
Solubility	Soluble at 20%	PASS
Concentration specs.	Meets specs.	PASS

This document has been produced electronically and is valid without signature

**Date of Print:** November 6, 2020

**G-Biosciences/Geno Technology, Inc.**

Rev 10/25/18 BP



**think proteins! think G-Biosciences** [www.GBiosciences.com](http://www.GBiosciences.com)