



G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: SDS [20%]

Catalog No: 786-017

Release Date: 02/02/21

Size: 1 L

Retest Date: 02/02/25

Lot No: 210202

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: February 19, 2021

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP



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