



G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: ND SB 201 (Non-detergent Sulfofetaine)

Catalog No: DG080

Release Date: 6/18/21

Size: 25g

Retest Date: 5/8/24

Lot No: 211606

Storage: RT

Specifications

CAS Number: 15471-17-7

Appearance: Off-white crystalline powder

Test	Specification	Result
Assay	>99%	99.2%

This document has been produced electronically and is valid without signature

Date of Print: January 28, 2022

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP



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