



BIOSCIENCES®

G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: OneQuant™ sulfoSMCC

Catalog No: 786-072

Release Date: 8/7/20

Size: 5 x 5mg

Retest Date: 8/7/22

Lot No: 200708

Storage: -20°C

Specifications

CAS Number: 92921-24-9

Appearance: Off-white to tan powder

Test	Specification	Result
Infrared spectrum	Consistent with structure and reference spectrum	PASS
Solubility	10mm solution in water should be clear with no insoluble material	PASS
TLC	Show a positive amine reaction with one spot present	PASS
HPLC	≥97.5%	PASS

This document has been produced electronically and is valid without signature

Date of Print: August 18, 2020

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP



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