



**G-Biosciences** ♦

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

A Geno Technology, Inc. (USA) brand name

## Certificate of Analysis / Data Sheet

**Product:** OneQuant™ sulfoSMCC

**Catalog No:** 786-082

**Release Date:** 08/03/2021

**Size:** 8 x 5mg

**Retest Date:** 08/03/2023

**Lot No:** 210308

**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

This document has been produced electronically and is valid without signature

**Date of Print:** November 4, 2021

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

Rev 10/25/18 BP



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