



Certificate of Analysis / Data Sheet

Product: BSOCOES

Catalog No: BC01

Release Date: 10/21/20

Size: 100mg

Retest Date: 10/21/22

Lot No: 202110

Storage: 4°C

Specifications

CAS Number: 57683-72-4

Chemical Formula: C₁₄ H₁₆ N₂ O₁₈ S₃

M.W.: 436.35

Appearance: White powder

Test	Specification	Result
Infrared	Conforms to structure and reference spectrum	PASS
Solubility	Clear Solution	PASS

This document has been produced electronically and is valid without signature

Date of Print: November 20, 2020

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP

