



Certificate of Analysis / Data Sheet

Product: DTSSP (3,3'-Dithiobis(sulfosuccinimidyl propionate))
Catalog No: BC-08 **Release Date:** 12/12/19
Size: 100mg **Retest Date:** 12/12/21
Lot No: 191212 **Storage:** 4°C

Specifications

CAS Number: 81069-02-5

Appearance: Off-white to light tan solid

Test	Specification	Result
Infrared Spectrum	Conforms to structure	PASS
Solubility	Clear 10mM solution	PASS
TLC	Show a positive amine reaction with one spot present	PASS
HPLC	≥ 94.0%	PASS

This document has been produced electronically and is valid without signature

Date of Print: December 24, 2019

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP

