



Certificate of Analysis / Data Sheet

Product: SB3-10

Catalog No: DG054

Size: 1g

Lot No: 191212

Release Date: 12/12/19

Retest Date: 12/12/23

Storage: Room Temp.

Specifications

CAS Number: 15163-36-7

Appearance: White powder

Test	Specification	Result
Solubility (color)	Colorless	Colorless
Solubility (Turbidity)	Clear	Clear
<i>100mg/ml in water</i>		
Conductivity (μ ohms/cm)	≤ 20	1
<i>0.1M aqueous solution @ 24°C</i>		
Infrared Spectrum	Conforms to structure	PASS
Purity (TLC)	$\geq 98\%$	100%
Carbon	57.4 – 59.8%	58.5%
Nitrogen	4.3 – 4.7%	4.6%

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

