



## Certificate of Analysis / Data Sheet

**Product:** HOOK-Sulfo-NHS-Biotin

**Catalog No:** BS-06

**Release Date:** 2/5/20

**Size:** 25mg

**Retest Date:** 2/5/21

**Lot No:** 200502

**Storage:** 4°C

### Specifications

**CAS Number:** 119616-38-5

**Appearance:** White powder

Test	Specification	Result
IR Spectrum	Conforms to reference	PASS
Solubility (10mM)	No insolubles	PASS
TLC	Show positive amine reaction with one spot present	PASS
Elemental	C – 37.92; N – 9.48; H – 4.10	C – 38.32; N – 10.02; H – 4.55

This document has been produced electronically and is valid without signature

**Date of Print:** February 17, 2020

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

Rev 10/25/18 BP

