



G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: HOOK™-PFP-Biotin

Catalog No: BG-10

Release Date: 5/18/2018

Size: 50mg

Retest Date: 5/18/2019

Lot No: 181505

Storage: 4°C

Specifications

CAS Number: NA

Appearance: White powder

Test	Specification	Result
Melting Point	195°C	Conforms
Identity Test	Passes (¹ H NMR)	Conforms
T.L.C.	One spot Support: Silica Gel 60; E. Merck #5715-7 Developing Solvent: [20:10:1, toluene, dioxane, acetic acid]; R _f = 0.58	
Assay	100% (300 MHz ¹ H NMR; DMSO- <i>d</i> ₆ ; Int Std) Microanalysis: C 46.83%; H 3.65%; N 6.81% Theory: C 46.83%; H 3.68%; N 6.82%	

This document has been produced electronically and is valid without signature

Date of Print: January 9, 2019

G-Biosciences/Geno Technology, Inc.
9800 Page Avenue, St. Louis, MO 63132, USA
Ph: 314-991-6034, Fax: 314-991-1504

Rev 05/18/18 BP



think proteins! think G-Biosciences www.GBiosciences.com