



G-Biosciences, St Louis, MO. USA ♦ 1-800-628-7730 ♦ 1-314- 991-6034 ♦ technical@GBiosciences.com

Certificate of Analysis

Product: N-succinimidyl-3-(4-hydroxyphenyl) propionate
(*Bolton-Hunter Reagent*)

Catalog #: BC84

Quantity: 1.0 gm

Lot#: 122302

Appearance: White powder

MW: 263.25

Tests:

Infrared Spectrum – Consistent with structure and reference spectrum Passed Yes / ~~No~~

This document has been produced electronically and is valid without signature

Date of Print: February 23, 2012

G-Biosciences/Genotech

92 Weldon Parkway
St. Louis, MO 63043, USA
Ph: 314-991-6034, Fax: 314-991-1504

Bolton-Hunter Reagent C of A/Com#3
Rev 01.12.11/IA



think proteins! think G-Biosciences!

www.GBiosciences.com