



Certificate of Analysis

Product: ABH; p-Azidobenzoyl Hydrazide

Catalog #: BC28

Quantity: 100mg

Lot#: 05082

MW: 177.2

Appearance: Tan to off-white

Tests:

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

This document has been produced electronically and is valid without signature

Date of Print: November 1, 2012

G-Biosciences

9800 Page Avenue

St. Louis, MO 63132, USA

Ph: 314-991-6034, Fax: 314-991-1504

ANB-NOS C of A/Rev 11.01.12/IA/Com#3

