



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

Certificate of Analysis

Product: ANB-NOS; N-5-azido-2-nitrobenzoyloxysuccinimide

MW: 305.2

Catalog #: BC29

Quantity: 100mg

Lot#: 030902

Appearance: Bright yellow powder

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: April 19, 2006

G-Biosciences

92 Weldon Parkway, St. Louis, MO 63043, USA

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

ANB-NOS CofA/Com/Temp, Rev 12.18.08-SA/MM/IA



think proteins! think G-Biosciences!

www.GBiosciences.com