



Certificate of Analysis

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

Catalog #: BC29

Quantity: 100mg

Lot#: 111711

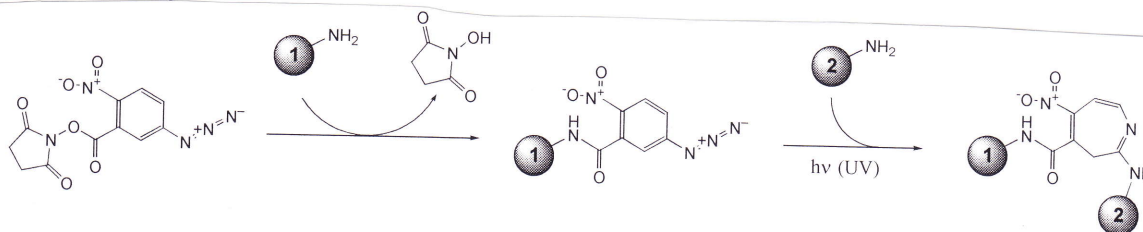
MW: 305.2

Appearance: Bright yellow powder

Tests:

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

Reaction Pathway for ANB-NOS:



This document has been produced electronically and is valid without signature

Date of Print: August 28, 2012

G-Biosciences/Geno Technology Inc

9800 Page Avenue

St. Louis, MO 63132, USA

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

ANB-NOS CofA/Com/Temp/Rev 07.01.12/IA



think proteins! think G-Biosciences!

www.GBiosciences.com