



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

Certificate of Analysis

Product: APG; p-Azidophenylglyoxal monohydrate

MW: 193.16

Catalog #: BC30

Quantity: 100mg

Lot#: 143005

Appearance: White fluffy 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: May 30, 2014

G-Biosciences / Genotech

9800 Page Avenue, St. Louis, MO 63132, USA

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

APG CofA, Rev 06.11.12/MM



think proteins! think G-Biosciences!

www.GBiosciences.com