



G-Biosciences, St Louis, MO. USA ♦ 1-800-628-7730 ♦ 1-314- 991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

## Certificate of Analysis

**Product:** APG; p-Azidophenylglyoxal monohydrate

**MW:** 193.16

**Catalog #:** BC30

**Quantity:** 100mg

**Lot#:** 123010

**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:** October 30, 2012

### **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](http://www.GBiosciences.com)