



G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: APG (p-Azidophenyl Glyoxal)

Catalog No: BC30

Release Date: 2/5/20

Size: 100mg

Retest Date: 2/5/21

Lot No: 200502

Storage: 4°C

Specifications

CAS Number: 1196151-49-1

Appearance: White fluffy powder

Test	Specification	Result
IR spectrum	Conforms to reference spectrum	PASS

This document has been produced electronically and is valid without signature

Date of Print: February 17, 2020

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP



think proteins! think G-Biosciences www.GBiosciences.com