



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

Certificate of Analysis

Product: Ammonium acetate [5M]

Catalog #: R012

Lot #: 092910

Tests:

Appearance - Clear Liquid	Passed	Yes / No
Specific gravity test	Passed	Yes / No
pH test	Passed	Yes / No

Storage: Store at room temperature

For Research Use Only

This document has been produced electronically and is valid without signature

Date of Print: November 29, 2010

G-Biosciences / Genotech/

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

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

Ammonium acetate 5M Cat# R012 C of A, Rev 12.18.08-SA/MM/IA



think proteins! think G-Biosciences!

www.GBiosciences.com