



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:** Ammonium acetate [5M]

**Catalog #:** R012

**Lot #:** 120805

### Tests:

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

**Storage:** Store at room temperature

*For Research Use Only*

This document has been produced electronically and is valid without signature

**Date of Print:** May 9, 2012

### **G-Biosciences / Genotech/**

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

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

Ammonium acetate 5M Cat# R012 C of A, Rev 05.09.12/MM



think proteins! think G-Biosciences!

[www.GBiosciences.com](http://www.GBiosciences.com)