



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 sulfate

**Catalog #:** RC-019

**Size:** 1kg

**Lot#:** 160103

**Mol Formula:** H<sub>8</sub>N<sub>2</sub>O<sub>4</sub>S

**MW:** 132.14

### Tests:

	<b>Min</b>	<b>Max</b>	<b>Actual</b>	<b>Units</b>
Purity (Dry Basis)	99.5		99.9	%
pH (5%, Water) at 25C	5.0	6.0	5.12	
Insolubles		0.005%	<0.005	%
Phosphate		0.0005	<0.0002	%
Chloride		0.0005	<0.0003	%
Residue on Ignition		0.005%	<0.005	%
Heavy Metals (as Pb)		0.0002%	<0.0002	%
Iron		0.0005%	0.0001	%
Selenium		0.0020	0.0005	%
Magnesium Test	PASS		PASS	P/F
Arsenic		0.0001	<0.00002	%

This document has been produced electronically and is valid without signature

**Storage Temp:** Room Temp. Protect from Moisture

**Date of Print:** November 11, 2016

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

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

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

Amm sulfate RC-019 C of A, Rev 07.16.12/MM



think proteins! think G-Biosciences!

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