



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

Certificate of Analysis

Product: Ammonium sulfate

Catalog #: RC-018

Size: 500g

Lot#: 132506

Mol Formula: H₈N₂O₄S

MW: 132.14

Tests:

	Min	Max	Actual	Units
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: June 26, 2013

G-Biosciences / Genotech

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

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

Amm sulfate RC-018 C of A, Rev 08.22.12/MM



think proteins! think G-Biosciences!

www.GBiosciences.com