



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:** Boric Acid  
**Size:** 500g  
**Catalog #:** RC-028  
**Lot#:** 131604  
**Appearance:** Fine white granular powder  
**Molecular Formula:** H<sub>3</sub>BO<sub>3</sub>  
**Mol Weight:** 58.44

Tests	Specifications	Results
Purity	99.5%	PASSED
Solubility	PASS	PASSED

**Storage:** Store at RT. Protect from moisture.

This document has been produced electronically and is valid without signature

**Date of Print:** April 17, 2013

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

9800 Page Avenue  
St. Louis, MO 63132, USA  
Ph: 314-991-6034, Fax: 314-991-1504

Boric acid, RC-028 C of A/Com#3/Rev 04.17.13/MM



think proteins! think G-Biosciences!

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