



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:** 1Kg  
**Catalog #:** RC-029  
**Lot#:** 101603  
**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:** November 30, 2010

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

92 Weldon Parkway, St. Louis, MO 63043, USA  
Ph: 314-991-6034, Fax: 314-991-1504

Boric acid, RC-028 C of A/Com#3/Rev 02.16.10/IA



**think proteins! think G-Biosciences!**

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