



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

Certificate of Analysis

Product: Boric Acid
Size: 1Kg
Catalog #: RC-029
Lot#: 122709
Appearance: Fine white granular powder
Molecular Formula: H_3BO_3
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: September 27, 2012

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 09.19.12/MM



think proteins! think G-Biosciences!

www.GBiosciences.com