



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

Certificate of Analysis

Product: Boric Acid
Size: 500g
Catalog #: RC-028
Lot#: 101603
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: 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