



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

Certificate of Analysis

Product: Acrylamide/BisAcrylamide 37.5:1 Premixed Powder

Catalog #: 786-503

Quantity: 40g

Lot#: 090911

Appearance: White free-flowing powder

Tests:

	Min	Max	Actual	Units	Disposition
Conductivity		10.0	<10.0	umhos	Passed
Solubility	PASS		PASSED	P/F	Passed
Electrophoresis	PASS		PASSED	P/F	Passed

This document has been produced electronically and is valid without signature

Storage Temp: Room Temp. Protect from Moisture

Date of Print: November 29, 2010

G-Biosciences / Genotech

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

Acrylamide Bis Acrylamide 29:1, 786-507 CofA Rev 12.16.08-SA/MM/IA



think proteins! think G-Biosciences!

www.GBiosciences.com