



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 #: 786-504

Quantity: 200g

Lot#: 110604

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

Storage Temp: Room Temp. Protect from Moisture

This document has been produced electronically and is valid without signature

Date of Print: October 12th, 2010

G-Biosciences / Genotech

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

Acrylamide Bis Acrylamide 37.5:1, 786-504 C of A
Rev 07.21.10/IA



think proteins! think G-Biosciences!

www.GBiosciences.com