



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

Certificate of Analysis

Product: Acrylamide/BisAcrylamide 29:1, 40% Solution

Catalog #: 786-505

Quantity: 500ml

Lot#: 100608

Appearance: Clear, colorless solution

Tests:

	Min	Max	Actual	Units	Disposition
Conductivity		10.0	8.9	umhos	Passed
pH at 25 ⁰ C	REPOERT		6.98		Passed
Electrophoresis	PASS		PASS	P/F	Passed

Storage Temp: Store cold. Keep dry. Warm it to room temperature before opening.

This document has been produced electronically and is valid without signature

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

crylamide Bis Acrylamide 29:1 40% Sol., 786-505/C of A
Rev 08.06.10/IA



think proteins! think G-Biosciences!

www.GBiosciences.com