



Certificate of Analysis

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

Catalog #: 786-505

Quantity: 500ml

Lot#: 140610

Appearance: Clear, colorless solution

Tests:

	Min	Max	Actual	Units	Disposition
Conductivity		10.0	8.9	umhos	Passed
pH at 25 ⁰ C	REPORT		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: December 17, 2014

G-Biosciences / Genotech

9800 Page Avenue
St. Louis, MO 63132, USA
Fax: 314-991-1504

Acrylamide Bis Acrylamide 29:1 40% Sol., 786-505/C of A
Rev 05.04.12/IA

