



Certificate of Analysis

Product: Acrylamide/BisAcrylamide 37.5:1, 40% Solution

Catalog #: 786-502

Quantity: 500ml

Lot#: 110501

Appearance: Clear, colorless solution

Tests:

	Min	Max	Actual	Units	Disposition
Conductivity		10.0	3.9	umhos	Passed
pH at 25 ⁰ C	REPORT		7.00		Passed
Electrophoresis	PASS		PASS	P/F	Passed
Polymerization Time		20	<20	min	Passed
DNase		None	None	P/F	Passed
RNase		None	None	P/F	Passed
Protease		None	None	P/F	Passed

Storage Temp: Room Temp. Protect from Moisture

This document has been produced electronically and is valid without signature

Date of Print: January 5, 2011

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 40% Sol., 786-502 C of A
Rev 03.09.10/IA

