



## Certificate of Analysis

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

**Catalog #:** 786-502

**Quantity:** 500ml

**Lot#:** 132203

**Appearance:** Clear, colorless solution

### Tests:

	<b>Min</b>	<b>Max</b>	<b>Actual</b>	<b>Units</b>	<b>Disposition</b>
Conductivity		10.0	3.9	umhos	Passed
pH at 25 <sup>0</sup> 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:** March 27, 2013

### **G-Biosciences / Genotech**

9800 Page Avenue  
St. Louis, MO 63132, USA  
Ph: 314-991-6034, Fax: 314-991-1504

Acrylamide Bis Acrylamide 37.5:1 40% Sol., 786-502 C of A  
Rev 03.27.13/MM

