



## Certificate of Analysis

**Product:** Acrylamide/BisAcrylamide 37.5:1 Premixed Powder

**Catalog #:** 786-503

**Quantity:** 40g

**Lot#:** 112302

**Appearance:** White free-flowing powder

### Tests:

	<b>Min</b>	<b>Max</b>	<b>Actual</b>	<b>Units</b>	<b>Disposition</b>
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:** February 24, 2011

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

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

Acrylamide Bis Acrylamide 29:1, 786-507 C of A  
Rev 02.24.11/IA

