



G-Biosciences, St Louis, MO. USA ♦ 1-800-628-7730 ♦ 1- 314- 991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

## Certificate of Analysis

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

**Catalog #:** 786-503

**Quantity:** 40g

**Lot#:** 130807

**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:** July 8, 2013

### **Genotech/GBiosciences**

9800 Page Avenue  
St. Louis, MO 63132, USA  
Ph: 314-991-6034, Fax: 314-991-1504  
[www.GBiosciences.com](http://www.GBiosciences.com)

Acrylamide Bis Acrylamide 29:1, 786-507 C of A  
Rev 07.08.2013/JK



think proteins! think G-Biosciences!

[www.GBiosciences.com](http://www.GBiosciences.com)