



## Certificate of Analysis

**Product:** Agarose I; Electrophoresis Matrix

**Cat #:** RC-149

**Quantity:** 500g

**Lot#:** 132404

**Grade:** Biotechnology Grade

**Appearance:** Fine White powder

<u>Assay:</u>	<u>Min</u>	<u>Max</u>	<u>Actual</u>	<u>Units</u>	<u>Disposition</u>
Gel Strength (1.5%)	1200		1280	g/cm <sup>2</sup>	Passed
Gelling Range (1.5%)	36	29	39	Degrees C	Passed
Melting Range (1.5%)	87	89	88.8	Degree C	Passed
EEO (-Mr)		0.10	0.06		Passed
Sulfate		0.15	0.06	%	Passed
RNase		NONE	NONE	P/F	Passed
DNase		NONE	NONE	P/F	Passed
Protease		NONE	NONE	P/F	Passed
Moisture		<10	7.1	%	Passed

**Storage:** Store at room temperature. Protect from moisture.

This document has been produced electronically and is valid without signature

**Date of Print:** April 25, 2013

### **Genotech/GBiosciences**

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

Agarose I, 129A-C C of A, Rev 08.27.12/MM



think proteins! think G Biosciences!