



## Certificate of Analysis

**Product:** Agarose I  
**Catalog #:** RC-122  
**Quantity:** 5g  
**Lot#:** 110411  
**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
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:** November 8, 2011

### **Genotech/GBiosciences**

92 Weldon Parkway, St. Louis, MO 63043, USA

Ph: 314-991-6034, Fax: 314-991-1504

Web: [www.GBiosciences.com](http://www.GBiosciences.com), Email: [technical@genotech.com](mailto:technical@genotech.com)

Agarose I, RC-122 C of A, Rev 12.16.08-SA/MM/IA



think proteins! think G Biosciences!