



Certificate of Analysis

Product: Agarose I
Catalog #: RC-005
Quantity: 25g
Lot#: 131906
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/cm2	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: June 24, 2013

Genotech/GBiosciences
 9800 Page Avenue
 St. Louis, MO 63132, USA
 Ph: 314-991-6034, Fax: 314-991-1504
www.GBiosciences.com

Agarose I, RC-005 C of A/Rev 08.27.12/MM



think proteins! think G Biosciences!