



Certificate of Analysis

Product: Agarose I; Electrophoresis Matrix

Cat #: RC-149

Quantity: 500g

Lot#: 111207

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 ²	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: July 13, 2011

Genotech/GBiosciences

92 Weldon Parkway
St. Louis, MO 63043, USA
Ph: 314-991-6034, Fax: 314-991-1504
Web: www.GBiosciences.com
Email: technical@GBiosciences.com

Agarose I, 129A-C C of A, Rev 07.13.11/IA



think proteins! think G Biosciences!