



Certificate of Analysis

Product: Agarose I
Catalog #: RC-122
Quantity: 5g
Lot#: 090911
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
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 29, 2010

Genotech/GBiosciences

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

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

Web: www.GBiosciences.com, Email: technical@genotech.com

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



think proteins! think G Biosciences!