



Certificate of Analysis

Product: Agarose I
Catalog #: RC-122
Quantity: 5g
Lot#: 131107
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 |
| 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 15, 2013

Genotech/GBiosciences

9800 Page Avenue, St. Louis, MO 63132, 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 04.25.13/MM



think proteins! think G Biosciences!