



Certificate of Analysis

Product: Acrylamide

Cat #: RC-002

Size: 500g

Lot#: 130708

Mol Formula: C₃H₅NO

MW: 71.08

Tests:

	Min	Max	Actual	Disposition
Purity (%)	98.0%		99.96	Passed
Melting Point (Degree C)	84	86	85.2	Passed
pH (10%, 0.1M NaCl) @ 25C	5.5	6.5	6.05	Passed
Insoluble		<0.005%	<0.005	Passed
Free Acrylic acid (%)		<0.0001	<0.0001	Passed
Conductivity (40%, Water, umhos)		5	2.2	Passed
Heavy Metals (as Pb)		0.005%		Passed
Iron		<0.0001%	0.0001	Passed
Magnesium		<0.0001%	<0.0001	Passed
RNase		None	None	Passed
DNase		None	None	Passed
Protease		None	None	Passed

Storage Temp: Store at Room Temp. Protect from Moisture

This document has been produced electronically and is valid without signature

Date of Print: February 2, 2016

G-Biosciences / Genotech

9800 Page Avenue
St. Louis, MO 63132, USA
Ph: 314-991-6034, Fax: 314-991-1504

Acrylamide RC-002 C of A
Rev 04.17.13/MM

