



G-Biosciences, St Louis, MO. USA ♦ 1-800-628-7730 ♦ 1-314- 991-6034 ♦ technical@GBiosciences.com

Certificate of Analysis

Product: Citric Acid Trisodium Salt Dihydrate

Catalog #: RC-040

Quantity: 500g

Lot#: 121707

Mol Formula: $C_6H_5Na_3O_7 \cdot 2H_2O$

MW: 294.1

Tests:

	<u>Specification</u>	<u>Result</u>
Appearance (Color)	White	White
Appearance (Form)	Powder	Powder
Solubility (Color)	Colorless	Cloroless
Sodium (Na)	21.1-24.6%	22.5%
Heavy Metals	< = 5ppm	< = 5ppm
Purity (Titration)	≥ 98.000%	99.953%

Storage Temp: Room Temp. Protect from Moisture

Date of Print: July 17, 2012

G-Biosciences / Genotech

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

Citric Acid Trisodium Salt Dihydrate CofA/Com#3
Rev 07.12.12/MM



think proteins! think G-Biosciences!

www.GBiosciences.com