



G-Biosciences, St Louis, MO. USA ♦ 1-800-628-7730 ♦ 1-314- 991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

## Certificate of Analysis

**Product:** Citric Acid Trisodium Salt Dihydrate

**Catalog #:** RC-040

**Quantity:** 500g

**Lot#:** 142503

**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	Colorless
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:** March 26, 2014

### **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](http://www.GBiosciences.com)