



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:** Calcium Chloride, Dihydrate

**Catalog #:** RC-030

**Quantity:** 500g

**Lot#:** 121503

**Molecular Formula:** CaCl<sub>2</sub>.2H<sub>2</sub>O

**Molecular Weight:** 147.0

**Grade:** Biotechnology Grade

**Appearance:** Fine white powder

**Tests:**

Assay	(Max 105.0%)	Passed	Yes / <del>No</del>
Heavy Metals	(Max 0.0005%)	Passed	Yes / <del>No</del>
pH (5% water) @ 25°C	(Max 8.5)	Passed	Yes / <del>No</del>

**Storage:** Store at room temperature. Protect from moisture.

This document has been produced electronically and is valid without signature

**Date of Print:** March 15, 2012

**G-Biosciences/ Genotech**

92 Weldon Parkway, St. Louis, MO 63043, USA  
Ph: 314-991-6034, Fax: 314-991-1504

Calcium Chloride, Dihydrate RC-030 CofA/Com#3  
Rev 12.9.10/IA



think proteins! think G-Biosciences!

[www.GBiosciences.com](http://www.GBiosciences.com)