



Certificate of Analysis

Product: Calcium Chloride, Dihydrate

Catalog #: RC-031

Quantity: 1kg

Lot#: 160709

Molecular Formula: CaCl₂.2H₂O

Molecular Weight: 147.0

Grade: Biotechnology Grade

Appearance: Fine white powder

Tests:

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

Storage: Store at room temperature. Protect from moisture.

This document has been produced electronically and is valid without signature

Date of Print: March 27, 2017

G-Biosciences/ Genotech

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

Calcium Chloride, Dihydrate RC-031 CofA/Com#3
Rev 09.12.12/MM

