



Certificate of Analysis

Product: Calcium Chloride, Dihydrate

Catalog #: RC-030

Quantity: 500g

Lot#: 141009

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: October 10, 2014

G-Biosciences/ Genotech

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

Calcium Chloride, Dihydrate RC-030 CofA/Com#3
Rev 07.25.12/MM

