



## Certificate of Analysis / Data Sheet

**Product:** Calcium chloride dihydrate

**Catalog No:** RC-030

**Release Date:** 6/8/2021

**Size:** 500g

**Retest Date:** 6/3/2026

**Lot No:** 210306

**Storage:** RT

### Specifications

**CAS Number:** 10035-04-8

**Appearance:** White colorless crystalline powder

**Grade:** ACS reagent

Test	Specification	Result
Solubility	326 g in 100 mL water at 60 °C	Pass
Heavy Metals	< 2 ppm	Pass
Anion	Sulfate (SO <sub>4</sub> <sup>2-</sup> ): < 0.005%	Pass
Cation	Ammonium (NH <sub>4</sub> ): < 0.001%	Pass
	Barium (Ba): < 0.001%	
	Iron (Fe): < 0.0005%	
	Magnesium (Mg): < 0.002%	
	Potassium (K): < 0.003%	
	Sodium (Na): < 0.007%	
	Strontium (Sr): < 0.02%	
pH	5.82	N/A
Oxidizable substances	0.0015%	N/A
Insoluble material	0.0012%	N/A
Assay by titration	99.98%	N/A

This document has been produced electronically and is valid without signature

**Date of Print:** January 28, 2022

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP



