



BIOSCIENCES®

G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: Magnesium Chloride Hexahydrate**Catalog No:** RC-069**Release Date:** 11/19/19**Size:** 500g**Retest Date:** 12/15/21**Lot No:** 191911**Storage:** RT

Specifications

CAS Number: 7791-18-6**Appearance:** White, odorless deliquescent crystals

Test	Specification	Result
Assay	99.0% - 101.0%	99.34%
Identification	PASS	PASS
Solubility (10% in water)	Clear & colorless	PASS
pH (5% in water)	4.5 – 7.0	5.57
Moisture	51.0% - 55.0%	52.8%
Insoluble Matter	≤ 0.005%	0.003%
Acidity/Alkalinity	PASS	PASS
Ammonium	≤ 0.002%	< 0.002%
Nitrate	≤ 0.001%	< 0.001%
Phosphate	≤ 0.0005%	< 0.0005%
Sulfate	≤ 0.002%	< 0.002%
Barium	PASS	PASS
Bromide	≤ 500ppm	< 500ppm
Calcium	≤ 0.01%	0.007%
Manganese	≤ 5ppm	3ppm
Potassium	≤ 0.005%	0.0015%
Sodium	≤ 0.005%	0.0017%



Strontium	$\leq 0.005\%$	0.0009%
Lead	$\leq 5\text{ppm}$	$< 5\text{ppm}$
Iron	$\leq 5\text{ppm}$	

This document has been produced electronically and is valid without signature

Date of Print: December 2, 2019

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP