



**G-Biosciences** ♦

♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

A Geno Technology, Inc. (USA) brand name

## Certificate of Analysis / Data Sheet

**Product:** Magnesium Chloride [1M]

**Catalog No:** R004

**Release Date:** 11/8/19

**Size:** 100mL

**Retest Date:** 11/8/22

**Lot No:** 190811

**Storage:** Room temp.

### Specifications

**Appearance:** Clear, colorless solution

| Test    | Specification | Result |
|---------|---------------|--------|
| pH test | 6.45-7.10     | 6.81   |

This document has been produced electronically and is valid without signature

**Date of Print:** November 18, 2019

**G-Biosciences/Geno Technology, Inc.**

Rev 10/25/18 BP



**think proteins! think G-Biosciences** [www.GBiosciences.com](http://www.GBiosciences.com)