



G-Biosciences, St Louis, MO. USA ♦ 1-800-628-7730 ♦ 1-314- 991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

### Certificate of Analysis

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

**Catalog #:** R004

**Size:** 100ml

**Lot #:** 190801

**Re-test date:** 01/2024

**Appearance:** Clear Liquid

**Tests:**

Magnesium chloride 1M solution	Passed	Yes / <del>No</del>
--------------------------------	--------	---------------------

Specific gravity test	Passed	Yes / <del>No</del>
-----------------------	--------	---------------------

**Storage:** Store it at RT.

This document has been produced electronically and is valid without signature
---

**Date of Print:** November 15, 2019

**G-Biosciences / Genotech**

9800 Page Avenue

St. Louis, MO 63132, USA

Ph: 314-991-6034, Fax: 314-991-1504



think proteins! think G-Biosciences!

[www.GBiosciences.com](http://www.GBiosciences.com)