



## Certificate of Analysis

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

**Catalog #:** R004

**Size:** 100ml

**Lot #:** 191401

**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
---

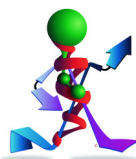
**Date of Print:** February 4, 2019

### **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)