



~
-

Certificate of Analysis

Product: Magnesium chloride [1M]

Catalog #: R004

Size: 100ml

Lot #: 201009

Re-test date: 09/2025

Appearance: Clear Liquid

Tests:

Magnesium chloride 1M solution Passed Yes / ~~No~~

pH test 6.45-7.1 Actual(6.83) Passed Yes / ~~No~~

Storage: Store it at RT.

This document has been produced electronically and is valid without signature

Date of Print: January 25, 2022

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