



G-Biosciences, St Louis, MO. USA ♦ 1-800-628-7730 ♦ 1-314- 991-6034 ♦ technical@GBiosciences.com

Certificate of Analysis

Product: Potassium chloride; KCl [1M]

Catalog #: R005

Lot #: 212810

Size: 100ml

Re-test date: 10-2026

Tests:

Appearance - Clear Liquid
Specific gravity test

Passed
Passed

Yes / ~~No~~
Yes / ~~No~~

Storage: Store at room temperature

This document has been produced electronically and is valid without signature

For Research Use Only

Date of Print: May 16, 2022

G-Biosciences / Genotech

9800 Page Avenue
St. Louis, MO 63132, USA
Ph: 314-991-6034, Fax: 314-991-1504

KCl 1M C of A/Com#3, Rev 06.12.12/MM



think proteins! think G-Biosciences!

www.GBiosciences.com