



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 #: 190403

Size: 100ml

Tests:

Appearance - Clear Liquid	Passed	Yes / No
Specific gravity test	Passed	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: September 12, 2019

Expiry: 3-04-2023

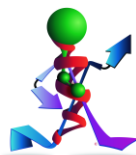
G-Biosciences / Genotech

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