



G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: Potassium Chloride [1M]

Catalog No: R005

Release Date: 03/24/2022

Size: 100 mL

Retest Date: 03/24/2024

Lot No: 222403

Storage: Room Temperature

Specifications

Appearance: Clear Liquid

Test	Specification	Result
pH	5.0-6.3	Pass

This document has been produced electronically and is valid without signature

Date of Print: July 27, 2023

G-Biosciences/Geno Technology, Inc.

KCl 1M C of A/Com#3, Rev 11.28.22/LO



think proteins! think G-Biosciences www.GBiosciences.com