



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

A Geno Technology, Inc. (USA) brand name

CERTIFICATE OF ANALYSIS

Product Name: Potassium phosphate Monobasic [1M]

Cat #: 786-488

Lot #: 201111

Size: 1L

Re-test date: 11/2025

Test Procedure:

Each batch of Potassium phosphate Monobasic [1M] solution is checked for correct molar specification of the buffer and its pH.

Specification for Potassium phosphate Monobasic [1M] to pass:

For Potassium phosphate Monobasic [1M] to pass, it must have correct molar specification of the buffer and the pH.

Molar Specification	Passed	Yes / No
pH Specification	Passed	Yes / No

Storage: Store at room temperature.

This document has been produced electronically and is valid without a signature

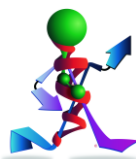
Date of print: November 20, 2020

G-Biosciences / Genotech

St. Louis, MO, USA

www.GBiosciences.com

Potassium phosphate Monobasic [1M] C of A
Rev 12.21.10/IA



think proteins! think G-Biosciences!

www.GBiosciences.com