



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

Certificate of Analysis

Product: Sodium Phosphate, monobasic (anhydrous)

Catalog #: RC-098

Size: 1kg

Lot#: 192609

Mol Formula: NaH₂PO₄

MW: 119.98

Tests:

	<u>Min</u>	<u>Max</u>	<u>Result</u>
Purity	99.5%		Passed
Moisture		0.3%	Passed
pH (1%, water) @25C	4.4	4.7	Passed
Insolubles		0.03%	Passed
Heavy Metals (as Pb)		0.001%	Passed

Storage Temp: Room Temp. Protect from Moisture

This document has been produced electronically and is valid without signature

Date of Print: October 15, 2019

G-Biosciences / Genotech

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

Sodium Phosphate, monobasic C of A/Com#3
Rev 04.28.15/IA



think proteins! think G-Biosciences!

www.GBiosciences.com