



G-Biosciences, St Louis, MO. USA ♦ 1-800-628-7730 ♦ 1-314- 991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

## Certificate of Analysis

**Product:** Sodium Phosphate Monobasic, Anhydrous

**Catalog #:** RC-097

**Quantity:** 500g

**Lot#:** 192609

**Mol Formula:** NaH<sub>2</sub>PO<sub>4</sub>

**MW:** 119.98

**Tests:**

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

**Storage Temp:** Room Temp. Protect from Moisture

This document has been produced electronically and is valid without signature

**Date of Print:** October 9, 2019

**G-Biosciences / Genotech**

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

Sodium Phosphate, monobasic RC-097CofA/Com#3  
Rev 05.24.12/MM



think proteins! think G-Biosciences!

[www.GBiosciences.com](http://www.GBiosciences.com)