



Certificate of Analysis / Data Sheet

Product: Potassium Phosphate (Monobasic)

Catalog No: RC-083

Release Date: 08/12/2021

Size: 500g

Retest Date: 08/12/2026

Lot No: 211208

Storage: Room Temp.

Specifications

CAS Number: 7778-77-0

Appearance: White crystalline powder

Test	Specification	Result
Heavy metals	< 0.001%	PASS
Chloride (Cl ⁻)	< 0.001%	PASS
Sulfate (SO ₄ ²⁻)	< 0.003%	PASS
Iron (Fe)	< 0.002%	PASS
Sodium (Na)	< 0.004%	PASS
pH	4.20-4.50	PASS
Loss on Drying	< 0.1%	PASS
Insoluble material	< 0.003%	PASS
Assay by titration	≥ 99.00%	PASS

This document has been produced electronically and is valid without signature

Date of Print: November 10, 2021



G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP