



Certificate of Analysis / Data Sheet

Product: Potassium Phosphate (Dibasic)

Catalog No: RC-082

Release Date: 02/04/21

Size: 1kg

Retest Date: 02/04/26

Lot No: 210402

Storage: Room Temp.

Specifications

CAS Number: 7758-11-4

Appearance: White powder

Test	Specification	Result
Assay	≥ 98.0%	98.47%
pH (5% solution)	8.5 – 9.6	9.38
Insoluble matter	≤ 0.01%	0.004%
Moisture (LOD)	≤ 1.0%	0.069%
Chloride	≤ 0.003%	< 0.003%
Nitrogen compounds	≤ 0.001%	< 0.001%
Sulfate	≤ 0.005%	< 0.005%
Heavy metals	≤ 5ppm	< 5ppm
Iron	≤ 0.001%	< 0.001%
Sodium	≤ 0.05%	0.0477%

This document has been produced electronically and is valid without signature

Date of Print: February 23, 2021

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP

