



Certificate of Analysis / Data Sheet

Product: Penicillin G Sodium Salt, USP grade***Catalog No: RC-741****Release Date: 08/05/21****Size: 50g****Retest Date: 08/05/23****Lot No: 210508****Storage: -20°C****Specifications****CAS Number: 69-57-8****Appearance: White crystalline powder**

Test	Specification	Result
Assay	98.0 – 102.0 %	99.5%
Crystallinity	PASS	PASS
Identification	PASS	PASS
Loss on drying	≤ 1.0%	0.1%
pH (6% solution @ 25°C)	5.5-7.5	7.2
Potency	1440 – 1680 U/mg	1585 U/mg
Residual solvent (ethyl acetate)	≤ 0.2%	0.003%
Residual solvent (N-butyl alcohol)	≤ 0.2%	0.01%

*Conforms to USP32 specification

This document has been produced electronically and is valid without signature

Date of Print: December 29, 2021**G-Biosciences/Geno Technology, Inc.**

Rev 10/25/18 BP

