



## Certificate of Analysis / Data Sheet

**Product:** Penicillin G Sodium Salt

**Catalog No:** RC-742

**Release Date:** 12/12/19

**Size:** 100g

**Retest Date:** 10/30/22

**Lot No:** 191212

**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%     |

This document has been produced electronically and is valid without signature

**Date of Print:** December 24, 2019

**G-Biosciences/Geno Technology, Inc.**

Rev 10/25/18 BP

