



Certificate of Analysis / Data Sheet

Product: Streptomycin Sulfate**Catalog No: RC-196****Release Date: 9/17/20****Size: 100g****Retest Date: 11/3/20****Lot No: 201709****Storage: -20°C****Specifications****CAS Number: 3810-74-0****Appearance: Fine White Powder**

| Test | Specification | Result |
|------------------------|--------------------------|----------------------|
| Appearance of Solution | Clear | Clear |
| Abnormal Toxicity | Meets Requirements | PASS |
| Bacterial Endotoxins | ≤ 0.25 IU/mg | < 0.25 IU/mg |
| Colorimetric | $\geq 90.0\%$ | 99.5% |
| Identification | Complies | PASS |
| Loss on Drying (EP) | $\leq 7.0\%$ | 4.9% |
| Loss on Drying (USP) | $\leq 5.0\%$ | 4.9% |
| Methanol (GC) | $\leq 0.3\%$ | $< 0.3\%$ |
| pH | 4.5-7.0 | 6.3 |
| Potency (Dried basis) | $\geq 720\mu\text{g/mg}$ | 738 $\mu\text{g/mg}$ |
| Residual Solvents | None present | Conforms |
| Streptomycin B | $\leq 3.0\%$ | $< 3.0\%$ |
| Sulfate | 18.0-21.5% | 19.8% |
| Sulfated Ash | $\leq 1.0\%$ | 0.5% |
| BSE/TSE | BSE/TSE Free | PASS |



This document has been produced electronically and is valid without signature

Date of Print: October 9, 2020

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP