



BIOSCIENCES®

G-Biosciences ♦

♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

A Geno Technology, Inc. (USA) brand name

## Certificate of Analysis / Data Sheet

**Product:** Kanamycin Sulfate, USP Grade

**Catalog No:** RC-183

**Release Date:** 10/27/20

**Size:** 10g

**Retest Date:** 2/13/23

**Lot No:** 202710

**Storage:** 4°C

### Specifications

**CAS Number:** 25389-94-0

**Appearance:** White powder

Test	Specification	Result
Chromatographic Purity	PASS/FAIL	PASS
Identification	PASS/FAIL	PASS
Loss on Drying	≤ 1.5%	0.6%
pH (1% solution)	6.5-8.5	8.1
Potency	≥ 750µg/mg	798µg/mg
Residue on Ignition	≤ 1.0%	0.2%
Specific Rotation	+116° to +123°	+121°

This document has been produced electronically and is valid without signature

**Date of Print:** November 6, 2020

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

Rev 10/25/18 BP



**think proteins! think G-Biosciences [www.GBiosciences.com](http://www.GBiosciences.com)**