



Certificate of Analysis / Data Sheet

Product: Nourseothricin Sulfate

Catalog No: RC-188

Release Date: 1/24/2019

Size: 100mg

Retest Date: 1/17/2020

Lot No: 191701

Storage: 4°C

Specifications

CAS Number: 96736-11-7

Appearance: Beige Powder

Test	Specification	Result
Components D & F (HPLC analysis)	>85%	91.9%
Components C & E (HPLC analysis)	<15%	8.1%
Biological Activity (Assay)	≥100% standard NR. 220200	110% standard NR. 220200
Ash	= 2.0±1.0% of material	2.0% of material

This document has been produced electronically and is valid without signature

Date of Print: December 23, 2019

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP

