



G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: Nourseothricin Sulfate

Catalog No: RC-190

Release Date: 4/13/20

Size: 500mg

Retest Date: 3/9/22

Lot No: 201304

Storage: 4°C

Specifications

CAS Number: 96736-11-7

Appearance: White Powder

Test	Specification	Result
Carbon (C)	33.0-34.0%	33.2%
Hydrogen (H)	6.0-6.5%	6.36%
Nitrogen (N)	15.0-16.0%	15.1%
Purity (HPLC)	≥ 85%	> 85%
Sulfur (S)	7.0-8.0%	7.2%

This document has been produced electronically and is valid without signature

Date of Print: April 15, 2020

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP



think proteins! think G-Biosciences www.GBiosciences.com