



A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: Novobiocin Sodium Salt

Catalog No: RC-736

Release Date: 10/22/2018

Size: 5 g

Retest Date: 06/22/2020

Lot No: 182210

Storage: -20°C

Specifications

CAS Number: 1476-53-5

Appearance: Off-white powder

Test	Specification	Result
Specific Optical Rotation [alpha]D	-58° to -50° (5% in MeOH/HCl, 25°C)	Complies
Activity	≥850 µg/mg	940 µg/mg
Sulfated Ash	≤ 15.0%	11.2%
Water	≤5.0%	0.4%
Assay (from N)	≥95.0%	>95.0%

This document has been produced electronically and is valid without signature

Date of Print: January 9, 2019

G-Biosciences/Geno Technology, Inc.
9800 Page Avenue, St. Louis, MO 63132, USA
Ph: 314-991-6034, Fax: 314-991-1504

Rev 02/13/17 CCH

