

Novobiocin, sodium salt

Catalog #: 27501

Lot #: 120910

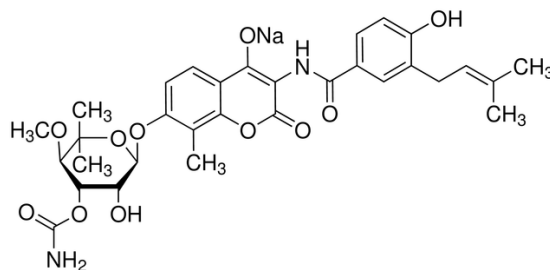
Size: 250 µl

Structure:

CAS Registry #: 1476-53-5

Purity: ≥93%

Chemical Formula: C₃₁H₃₅N₂NaO₁₁



Molecular Weight: 634.61

Description: Novobiocin is an aminocoumarin antibiotic that is produced by *Streptomyces niveus*. Along with other aminocoumarin antibiotics, novobiocin is a very potent inhibitor of bacterial DNA gyrase. It works by targeting the GyrB subunit of the enzyme involved in energy transduction.

Appearance: A beige powder

Formulation: Supplied as a 400 µM solution in anhydrous DMSO

Biological Activity: Novobiocin has been shown to inhibit bacterial DNA gyrase supercoiling with an IC₅₀ of 0.08 µM.

Storage/Stability: Store at or below -20 °C.

Quality Control: The purity was determined by HPLC analysis.

Reference(s):

1. Lamour V, *et al.*, *J Biol Chem.* 2002 May 24; **277**(21):18947-53.
2. Alt, S., *et al.*, *J Antimicrob Chemother.* 2011 Jun 11; doi:10.1093/jac/dkr247.