

Novo-TRX

Catalog #: 27500

Lot #: 120723

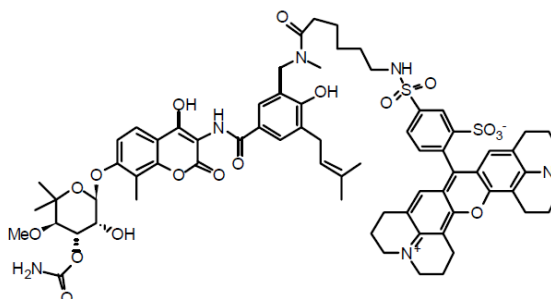
Size: 20 μ l

Structure:

CAS Registry #: 1476-53-5 (Novobiocin, sodium salt)

Purity: \geq 93.3%

Chemical Formula: C₇₀H₈₀N₆O₁₈S₂



Molecular Weight: 1357.5

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. Novo-TRX is Texas Red-labeled novobiocin, which can be used as a fluorescence probe in assays for screening DNA gyrase inhibitors.

Appearance: A red aqueous solution

Formulation: Supplied as a 10 μ M solution in anhydrous DMSO

Biological Activity: Novo-TRX has been shown to inhibit bacterial DNA gyrase activity with an IC₅₀ of 0.6 μ M.

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. Glaser BT, *et al.*, *J Biomol Screen.* 2011 Feb;**16(2)**:230-8.