

PRODUCT INFORMATION

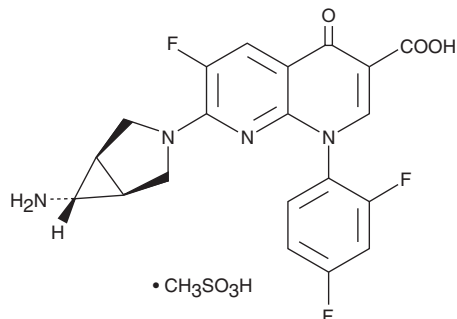


Trovafloxacin (mesylate)

Item No. 9000303

CAS Registry No.: 147059-75-4
Formal Name: 7-[(1 α ,5 α ,6 α)-6-amino-3-azabicyclo[3.1.0]hex-3-yl]-1-(2,4-difluorophenyl)-6-fluoro-1,4-dihydro-4-oxo-1,8-naphthyridine-3-carboxylic acid, monomethanesulfonate

Synonym: CP 99,219
MF: C₂₀H₁₅F₃N₄O₃ • CH₄SO₃
FW: 512.5
Purity: ≥98%
UV/Vis.: λ_{\max} : 269, 342 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Trovafloxacin (mesylate) is supplied as a crystalline solid. A stock solution may be made by dissolving the trovafloxacin (mesylate) in the solvent of choice, which should be purged with an inert gas. Trovafloxacin (mesylate) is soluble in the organic solvent DMSO at a concentration of approximately 25 mg/ml. It is also soluble in water at a concentration of approximately 20 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Trovafloxacin is a broad spectrum antibiotic that inhibits the uncoiling of supercoiled DNA in bacteria by blocking the activity of topoisomerase IV (IC₅₀ = 3.02 μ g/ml) and DNA gyrase (IC₅₀ = 7.13 μ g/ml).¹

Reference

1. Takei, M., Fukuda, H., Kishii, R., *et al.* Target preference of 15 quinolones against *Staphylococcus aureus*, based on antibacterial activities and target inhibition. *Antimicrob. Agents Chemother.* **45(12)**, 3544-3547 (2001).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 04/10/2024

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM