

PRODUCT INFORMATION

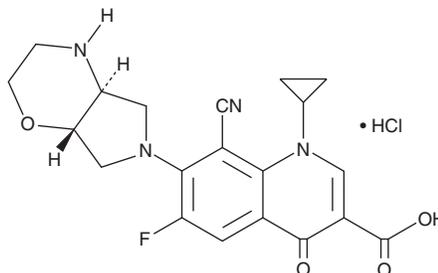


Finafloxacin (hydrochloride)

Item No. 22410

CAS Registry No.: 209342-41-6
Formal Name: 8-cyano-1-cyclopropyl-6-fluoro-7-[[4a*S*,7a*S*]-hexahydropyrrolo[3,4-*b*]-1,4-oxazin-6(2*H*)-yl]-1,4-dihydro-4-oxo-3-quinolinecarboxylic acid, monohydrochloride

Synonym: BAY35-337
MF: C₂₀H₁₉FN₄O₄ • HCl
Purity: ≥98%
FW: 434.9
UV/Vis.: λ_{max}: 245, 286, 371 nm
Supplied as: A 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

Finafloxacin (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the finafloxacin (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Finafloxacin (hydrochloride) is soluble in DMSO.

Description

Finafloxacin is a fluoroquinolone antibiotic.¹ It is active against *S. pneumoniae* and *P. aeruginosa*, as well as methicillin-resistant *S. aureus* (MRSA) that is sensitive or resistant to fluoroquinolones (MIC₉₀s = 2, 2, 0.125, and 4 mg/L, respectively). Finafloxacin (23.1 mg/kg) improves survival in *F. tularensis*- or *Y. pestis*-infected mice in models of tularemia or plague, respectively.² Formulations containing finafloxacin have been used in the treatment of acute otitis externa.

References

1. Kocsis, B., Domokos, J., and Szabo, D. Chemical structure and pharmacokinetics of novel quinolone agents represented by avarofloxacin, delafloxacin, finafloxacin, zabofloxacin and nemonoxacin. *Ann. Clin. Microbiol. Antimicrob.* **15**(1), 34 (2016).
2. Barnes, K.B., Richards, M.I., Laws, T.R., et al. Finafloxacin is an effective treatment for inhalational tularemia and plague in mouse models of infection. *Antimicrob. Agents Chemother.* **65**(6), e02294-20 (2021).

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, 10/19/2022

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