

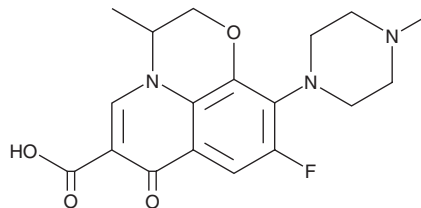
PRODUCT INFORMATION



Ofloxacin

Item No. 22891

CAS Registry No.: 82419-36-1
Formal Name: 9-fluoro-2,3-dihydro-3-methyl-10-(4-methyl-1-piperazinyl)-7-oxo-7H-pyrido[1,2,3-de]-1,4-benzoxazine-6-carboxylic acid
Synonyms: DL-8280, WPO405
MF: C₁₈H₂₀FN₃O₄
FW: 361.4
Purity: ≥95%
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.

Description

Ofloxacin is a broad-spectrum fluoroquinolone antibiotic that prevents supercoiling of bacterial chromosomes by DNA gyrase.^{1,2} It is active against Gram-positive and Gram-negative bacteria with MIC₉₀s ranging from 0.39 to 3.13 µg/ml for clinical isolates of *S. aureus*, *S. epidermidis*, *S. pyogenes*, and *S. faecalis* and ≤0.78 µg/ml for *N. gonorrhoeae* and various species of Enterobacteriaceae. Ofloxacin is active *in vivo*, with ED₅₀ values ranging from 0.7 to 75.1 mg/kg in infected mice. Formulations containing ofloxacin have been used to treat urinary tract infections, gonorrhea, prostatitis, and gastroenteritis.³

References

1. Sato, K., Matsuura, Y., Inoue, M., *et al.* *In vitro* and *in vivo* activity of DL-8280, a new oxazine derivative. *Antimicrob. Agents and Chemother.* **22(4)**, 548-553 (1982).
2. Smith, J.T. The mode of action of 4-quinolones and possible mechanisms of resistance. *J. Antimicrob. Chemother.* **18 (Suppl. D)**, 21-29 (1986).
3. Hooper, D.C., and Wolfson, J.S. The fluoroquinolones: pharmacology, clinical uses, and toxicities in humans. *Antimicrob. Agents Chemother.* **28(5)**, 716-721 (1985).

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, 02/24/2021

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