

PRODUCT INFORMATION

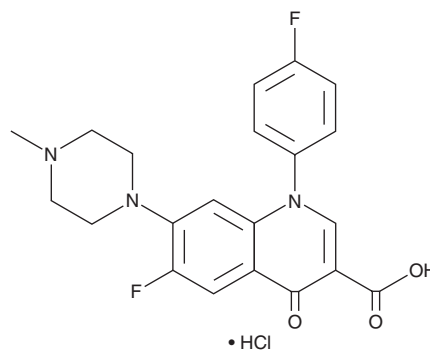


Difloxacin (hydrochloride)

Item No. 29000

CAS Registry No.: 91296-86-5
Formal Name: 6-fluoro-1-(4-fluorophenyl)-1,4-dihydro-7-(4-methyl-1-piperazinyl)-4-oxo-3-quinolinecarboxylic acid, monohydrochloride

Synonym: A-56619
MF: C₂₁H₁₉F₂N₃O₃ • HCl
FW: 435.9
Purity: ≥98%
UV/Vis.: λ_{max}: 281, 319, 332 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.

Description

Difloxacin is a fluoroquinolone antibiotic.¹ It is active against isolates of anaerobic bacteria, including *B. fragilis*, as well as *Fusobacterium* and *Actinomyces* species with MIC values ranging from ≤0.125 to 8 µg/ml. It eliminates *E. coli* and *B. fragilis* infection in a rat intra-abdominal abscess model when administered at a dose of 40 mg/kg three times per day.²

References

1. Bansal, M.B. and Thadepalli, H. Activity of difloxacin (A-56619) and A-56620 against clinical anaerobic bacteria in vitro. *Antimicrob. Agents Chemother.* **31(4)**, 619-621 (1987).
2. Thadepalli, H., Gollapudi, S.V., and Chuah, S.K. Therapeutic evaluation of difloxacin (A-56619) and A-56620 for experimentally induced *Bacteroides fragilis*-associated intra-abdominal abscess. *Antimicrob. Agents Chemother.* **30(4)**, 574-576 (1986).

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, 11/15/2022

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