

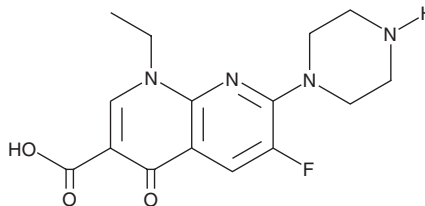
PRODUCT INFORMATION



Enoxacin

Item No. 16956

CAS Registry No.: 74011-58-8
Formal Name: 1-ethyl-6-fluoro-1,4-dihydro-4-oxo-7-(1-piperazinyl)-1,8-naphthyridine-3-carboxylic acid
Synonym: NSC 629661
MF: C₁₅H₁₇FN₄O₃
FW: 320.3
Purity: ≥98%
UV/Vis.: λ_{max}: 220, 270, 340 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

Enoxacin is supplied as a crystalline solid. A stock solution may be made by dissolving the enoxacin in the solvent of choice, which should be purged with an inert gas. Enoxacin is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of enoxacin in these solvents is approximately 0.1 and 15 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of enoxacin can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of enoxacin in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Enoxacin is a fluoroquinolone antibiotic.¹⁻⁴ It is active against clinical isolates of a variety of Gram-positive and Gram-negative bacteria, including *S. aureus*, *E. coli*, *K. pneumoniae*, *P. aeruginosa*, and *S. marcescens* (MIC_{50s} = 1, 0.12, 0.25, 0.5, and 1 mg/L, respectively).¹ Enoxacin inhibits *S. aureus* DNA gyrase supercoiling activity and topoisomerase IV DNA decatenation (IC_{50s} = 126 and 26.5 μg/ml, respectively).² It increases survival in mouse models of systemic *S. aureus*, *E. coli*, *K. pneumoniae*, *P. aeruginosa*, and *S. marcescens* infection with ED₅₀ values of 15.1, 2.2, 4.1, 120.3, and 7.6 mg/kg, respectively.³ Enoxacin (4 and 8 mg/kg per day) also reduces tumor growth in a 143B human osteosarcoma mouse xenograft model.⁴ Formulations containing enoxacin have previously been used in the treatment of urinary tract infections and gonorrhea.

References

1. Clarke, A.M., Zemcov, S.J., and Campbell, M.E. *In-vitro* activity of pefloxacin compared to enoxacin, norfloxacin, gentamicin and new β-lactams. *J. Antimicrob. Chemother.* **15(1)**, 39-44 (1985).
2. 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).
3. Ozaki, M., Matsuda, M., Tomii, Y., *et al.* In vivo evaluation of NM441, a new thiazeto-quinoline derivative. *Antimicrob. Agents Chemother.* **35(12)**, 2496-2499 (1991).
4. Luo, X., Liu, X., Tao, Q., *et al.* Enoxacin inhibits proliferation and invasion of human osteosarcoma cells and reduces bone tumour volume in a murine xenograft model. *Oncol. Lett.* **20(2)**, 1400-1408 (2020).

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, 11/30/2022

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