

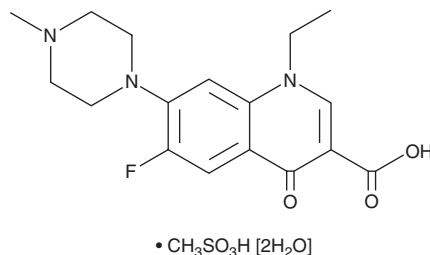
PRODUCT INFORMATION



Pefloxacin (mesylate hydrate)

Item No. 29103

CAS Registry No.: 149676-40-4
Formal Name: 1-ethyl-6-fluoro-1,4-dihydro-7-(4-methyl-1-piperazinyl)-4-oxo-3-quinolinecarboxylic acid, monomethanesulfonate, dihydrate
MF: $C_{17}H_{20}FN_3O_3 \cdot CH_3SO_3H [2H_2O]$
FW: 465.5
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 280, 318, 332 nm
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

Pefloxacin (mesylate hydrate) is supplied as a solid. A stock solution may be made by dissolving the pefloxacin (mesylate hydrate) in the solvent of choice, which should be purged with an inert gas. Pefloxacin (mesylate hydrate) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of pefloxacin (mesylate hydrate) in these solvents is approximately 5 and 3 mg/ml, respectively.

Pefloxacin (mesylate hydrate) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, pefloxacin (mesylate hydrate) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Pefloxacin (mesylate hydrate) has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Pefloxacin is a broad-spectrum and synthetic fluoroquinolone antibiotic and a methylated derivative of norfloxacin (Item No. 25975).^{1,2} It is active against Gram-negative and Gram-positive bacteria, including clinical isolates of *E. coli*, *K. pneumoniae*, *S. aureus*, and *G. vaginalis* (MIC_{90s} = 0.12, 0.5, 0.5, and 32 mg/L, respectively).^{1,2} Pefloxacin selectively inhibits *E. coli* topoisomerase I and DNA gyrase over calf thymus topoisomerase I (IC_{50s} = 45, 4.5, and >140 $\mu g/ml$, respectively) in cell-free assays.³ *In vivo*, pefloxacin is active against systemic *E. coli*, *K. pneumoniae*, and *P. aeruginosa* infections in mice with 50% of efficacy maximum (P_{50}) values of 6.5, 29.2, and 219 mg/kg, respectively.⁴

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. Jones, B.M., Geary, I., Lee, M.E., et al. Activity of pefloxacin and thirteen other antimicrobial agents *in vitro* against isolates from hospital and genitourinary infections. *J. Antimicrob. Chemother.* **17**(6), 736-746 (1986).
3. Tabary, X., Moreau, N., Dureuil, C., et al. Effect of DNA gyrase inhibitors pefloxacin, five other quinolones, novobiocin, and clorobiocin on *Escherichia coli* topoisomerase I. *Antimicrob. Agents Chemother.* **31**(12), 1925-1928 (1987).
4. Fantin, B., Leggett, J., Ebert, S., et al. Correlation between *in vitro* and *in vivo* activity of antimicrobial agents against gram-negative bacilli in a murine infection model. *Antimicrob. Agents Chemother.* **35**(7), 1413-1422 (1991).

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, 11/01/2022

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