

PRODUCT INFORMATION



N-desmethyl Levofloxacin (hydrochloride)

Item No. 27178

Formal Name: (3S)-9-fluoro-2,3-dihydro-3-methyl-7-oxo-10-(1-piperazinyl)-7H-pyrido[1,2,3-de]-1,4-benzoxazine-6-carboxylic acid, hydrochloride

Synonym: Desmethyl Levofloxacin

MF: C₁₇H₁₈FN₃O₄ • HCl

FW: 347.3

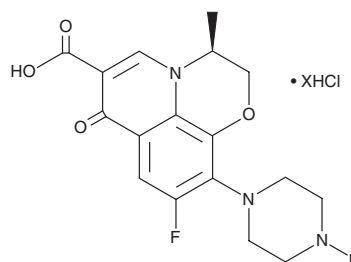
Purity: ≥98%

UV/Vis.: λ_{max}: 223, 294, 326 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

N-desmethyl Levofloxacin (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the N-desmethyl levofloxacin (hydrochloride) in the solvent of choice, which should be purged with an inert gas. N-desmethyl Levofloxacin (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of N-desmethyl levofloxacin (hydrochloride) in these solvents is approximately 1, 25, and 30 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 N-desmethyl levofloxacin (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of N-desmethyl levofloxacin (hydrochloride) in PBS, pH 7.2, is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

N-desmethyl Levofloxacin is an active metabolite of the fluoroquinolone antibiotic levofloxacin (Item No. 20382).¹ It is active against *S. aureus*, *S. epidermidis*, *B. subtilis*, *E. coli*, *P. aeruginosa*, and *K. pneumoniae* (MICs = 4, 1, 1, 0.012, >4, and 0.25 µg/ml, respectively).

Reference

1. Mohammadhosseini, N., Alipanahi, Z., Alipour, E., *et al.* Synthesis and antibacterial activity of novel levofloxacin derivatives containing a substituted thienylethyl moiety. *Daru*. **20(1)**, (2012).

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, 12/21/2023

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