

# PRODUCT INFORMATION



## Desmethyl Ofloxacin (hydrochloride)

Item No. 33807

**CAS Registry No.:** 402930-70-5  
**Formal Name:** 9-fluoro-2,3-dihydro-3-methyl-7-oxo-10-(1-piperazinyl)-7H-pyrido[1,2,3-de]-1,4-benzoxazine-6-carboxylic acid, monohydrochloride

**MF:** C<sub>17</sub>H<sub>18</sub>FN<sub>3</sub>O<sub>4</sub> • HCl

**FW:** 383.8

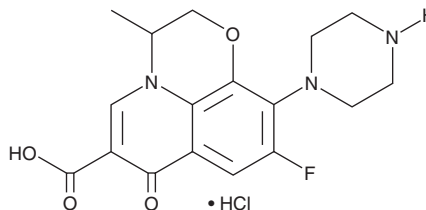
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 294 nm

**Supplied as:** A 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

Desmethyl ofloxacin (hydrochloride) is supplied as a solid. Aqueous solutions of desmethyl ofloxacin (hydrochloride) can be prepared by directly dissolving the solid in aqueous buffers. Desmethyl ofloxacin (hydrochloride) is slightly soluble in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

### Description

Desmethyl ofloxacin is a metabolite of the fluoroquinolone antibiotic ofloxacin (Item No. 22891).<sup>1</sup>

### Reference

1. Sudo, K., Okazaki, O., Tsumura, M., *et al.* Isolation and identification of metabolites of ofloxacin in rats, dogs and monkeys. *Xenobiotica* **16(8)**, 725-732 (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, 10/19/2022

#### 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