

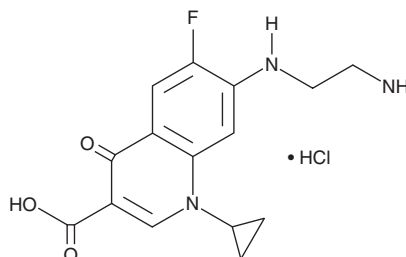
PRODUCT INFORMATION



Desethylene Ciprofloxacin (hydrochloride)

Item No. 22123

CAS Registry No.: 528851-31-2
Formal Name: 7-[(2-aminoethyl)amino]-1-cyclopropyl-6-fluoro-1,4-dihydro-4-oxo-3-quinolinecarboxylic acid, monohydrochloride
MF: C₁₅H₁₆FN₃O₃ • HCl
FW: 341.8
Purity: ≥98%
UV/Vis.: λ_{max}: 278 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

Desethylene ciprofloxacin is a major metabolite of ciprofloxacin (Item No. 14286).¹ It is also a degradation product of ciprofloxacin that can be created through advanced oxygenation processes as a potential way to remove ciprofloxacin from wastewater.² It is also created during degradation of ciprofloxacin by chlorination.³

References

1. Bergan, T., Thorsteinsson, S.B., Rohwedder, R., *et al.* Elimination of ciprofloxacin and three major metabolites and consequences of reduced renal function. *Chemotherapy* **35(6)**, 393-405 (1989).
2. De Witte, B., Dewulf, J., Demeestere, K., *et al.* Ozonation and advanced oxidation by the peroxide process of ciprofloxacin in water. *J. Hazard. Mater.* **161(2-3)**, 701-708 (2009).
3. Wang, H., Hu, C., Liu, L., *et al.* Interaction of ciprofloxacin chlorination products with bacteria in drinking water distribution systems. *J. Hazard. Mater.* **339**, 174-181 (2017).

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