

Ciprofloxacin

Fluoroquinolone antibacterial agent

Fluoroquinolone antibacterial agent. Inhibits bacterial DNA gyrase (topoisomerase). Inhibits cell division and causes double-strand breaks in the bacterial chromosome. Broad-spectrum antibiotic effective against Gram-positive, Gram-negative and anaerobic bacteria. Effective against *Mycobacterium tuberculosis*. Potentially anxiogenic.

Citations: 10

[View Online »](#)

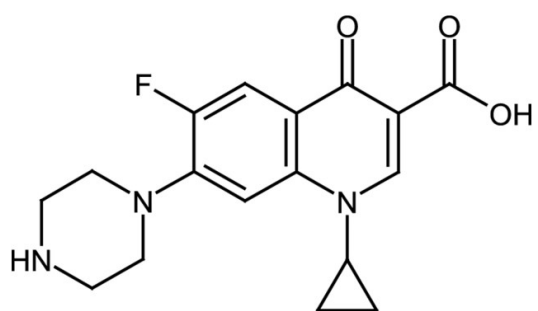
Ordering Information

[Order Online »](#)

ALX-380-287-G005	5g
ALX-380-287-G025	25g

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 1-Cyclopropyl-6-fluoro-1,4-dihydro-4-oxo-7-(1-piperazinyl)-3-quinolinecarboxylic acid

Appearance White to light yellow crystalline powder

CAS 85721-33-1

Couple Target DNA gyrase

Couple Type Inhibitor

Formula $C_{17}H_{18}FN_3O_3$

MI **14:** 2314

MW 331.3

Purity ≥98% (HPLC)

RTECS VB1993800

Solubility Soluble in dilute hydrochloric acid; insoluble in water.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)