

Sparfloxacin

Fluoroquinolone antibacterial agent

Fluoroquinolone with broad spectrum antibacterial activity against Gram-positive and Gram-negative bacteria. Inhibits bacterial DNA gyrase (topoisomerase). Effective against *Mycobacterium tuberculosis*. Shows photosensitizing effects.

Citations: 7

[View Online »](#)

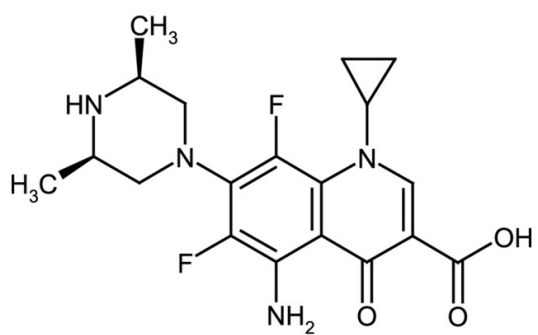
Ordering Information

[Order Online »](#)

ALX-380-302-G005	5g
------------------	----

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 rel-5-Amino-1-cyclopropyl-7-[(3R,5S)-3,5-dimethyl-1-piperazinyl]-6,8-difluoro-1,4-dihydro-4-oxo-3-quinolinecarboxylic acid, AT-4140

Appearance White to yellow crystalline solid.

CAS 110871-86-8

Couple Target DNA gyrase

Couple Type Inhibitor

Formula $C_{19}H_{22}F_2N_4O_3$

Identity Determined by IR.

MI 14: 8735

MW 392.4

Purity ≥98% (Assay)

RTECS VB1986500

Solubility Soluble in 100% ethanol or 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-
be@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)