

Tosufloxacin . tosylate

Fluoroquinolone antibacterial agent

Fluoroquinolone antibacterial agent. Inhibits bacterial DNA gyrase (topoisomerase). Effective against Gram-positive and Gram-negative aerobic bacteria, anaerobic bacteria and *Chlamydia trachomatis*.

Citations: 8

[View Online »](#)

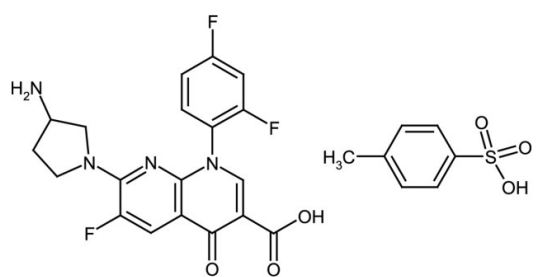
Ordering Information

[Order Online »](#)

ALX-380-303-M250	250mg
------------------	-------

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 7-(3-Amino-1-pyrrolidinyl)-1-(2,4-difluorophenyl)-6-fluoro-1,4-dihydro-4-oxo-1,8-naphthyridine-3-carboxylic acid . tosylate, A-64730

Appearance White to yellow crystalline solid.

CAS 115964-29-9

Couple Target DNA gyrase

Couple Type Inhibitor

Formula $C_{19}H_{15}F_3N_4O_3 \cdot C_7H_8O_3S$

MI 14: 9555

MW 404.4 . 172.2

Purity ≥98% (HPLC)

Solubility Slightly soluble 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-eu@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)