

Nadifloxacin

Fluoroquinolone antibacterial agent

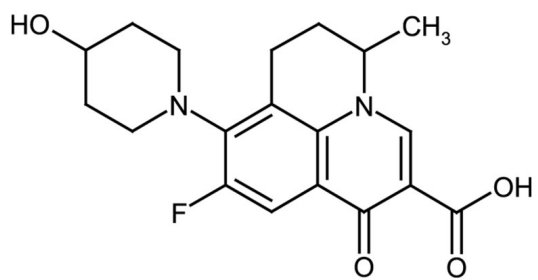
Fluoroquinolone with broad spectrum activity against Gram-negative and Gram-positive bacteria. Inhibits the generation of O_2^- and OH radicals by neutrophils. Inhibits the production of inflammatory cytokines.

Citations: 8

[View Online »](#)

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	9-Fluoro-6,7-dihydro-8-(4-hydroxypiperidino)-5-methyl-1-oxo-1H,5H-benzol[ij]quinolizine-2-carboxylic acid, Jinofloxacin, OPC-7251
Appearance	White to yellow crystalline solid.
CAS	124858-35-1
Couple Target	DNA gyrase
Couple Type	Inhibitor
Formula	$C_{19}H_{21}FN_2O_4$
Identity	Determined by IR.
MI	14: 6345
MW	360.4
Purity	≥98% (Assay)
Solubility	Soluble in DMSO.

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-
be@enzolifesciences.com](mailto:info-be@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)