

# Rufloxacin . hydrochloride

Fluoroquinolone antibacterial agent

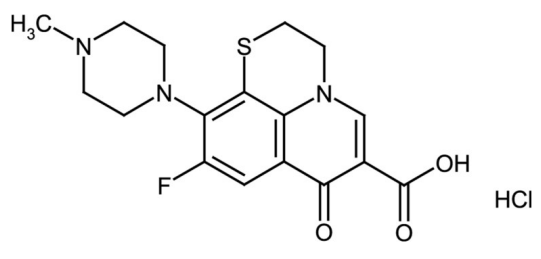
Fluoroquinolone antibacterial agent. Structurally similar to ofloxacin. Specific inhibitor of bacterial DNA gyrase (topoisomerase). Inhibits B-cell differentiation. Shows photosensitizing effects.

Citations: 10

[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** 6-Fluoro-2,3-dihydro-10-(4-methyl-1-piperazinyl)-7-oxo-7H-pyrido[1,2,3-de]-1,4-benzothiazine-6-carboxylic acid . hydrochloride, MF-934 . hydrochloride

**Appearance** Off-white to light green crystalline solid.

**CAS** 106017-08-7

**Couple Target** DNA gyrase

**Couple Type** Inhibitor

**Formula**  $C_{17}H_{18}FN_3O_3S \cdot HCl$

**Identity** Identity determined by IR.

**MI** 14: 8294

**MW** 363.4 . 36.5

**Purity** ≥98% (Assay)

**Solubility** Sparingly soluble in water; very slightly soluble in methanol; almost insoluble in chloroform or acetic acid.

**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)