Sparfloxacin

Fluoroquinolone antibacterial agent

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

Citations: 7

View Online »

Ordering Information

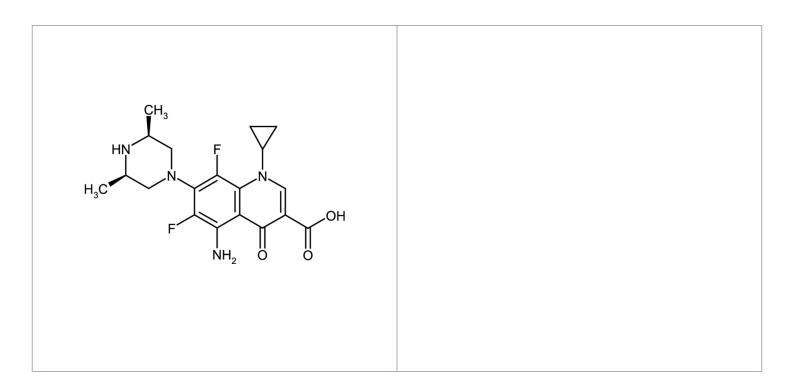
Order Online »

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

Soluble in 100% ethanol or water.