

Levofloxacin Hydrochloride Monohydrate

**Fluoroquinolone, S-(-) isomer of ofloxacin;
topoisomerase IV and bacterial DNA gyrase**

Inhibitor. It is the optical S-(-) isomer of the fluoroquinolone antibacterial, ofloxacin. It has twice the potency of ofloxacin.

Ordering Information

[Order Online »](#)

LKT-L1786-G005	5g
----------------	----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name S-(-)-Ofloxacin, Levofloxacin

Couple Type Inhibitor

Formula $C_{18}H_{20}FN_3O_4 \cdot HCl \cdot H_2O$

MW 415.84

Purity $\geq 90\%$

Solubility Water(25mg/mL)



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com