

Fluoxetine hydrochloride

FIASMA, SERT inhibitor, 5-HT and $\sigma 1$ antagonist.

A specific serotonin uptake inhibitor. It was found to exert inhibitory or stimulating effects on T-lymphocyte proliferation depending on the concentration.

Ordering Information

[Order Online »](#)

LKT-F4780-G005	5g
----------------	----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

CAS 59333-67-4

Couple Type Inhibitor

Formula $C_{17}H_{18}F_3NO \cdot HCl$

MI 13: 4211

MW 345.79

Purity $\geq 98\%$

Solubility Soluble in methanol, ethanol, acetone, acetonitrile and chloroform. Water to 10 mM, DMSO to 100 mM. DMF to 16 mg/mL. Ethanol to 69mg/mL



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)