

Norfluoxetine . HCl

Fluoxetine Metabolite

Fluoxetine metabolite that induces long QT syndrome via slow delayed rectifier potassium current block. It is a time-dependent inhibitor of CYP2C19 and CYP3A4. The active N-demethylated metabolite is a potent and selective inhibitor of serotonin uptake in vivo. Also inhibits coxsackievirus replication acting via inhibition of viral RNA synthesis. State-dependent TREK-2 blocker.

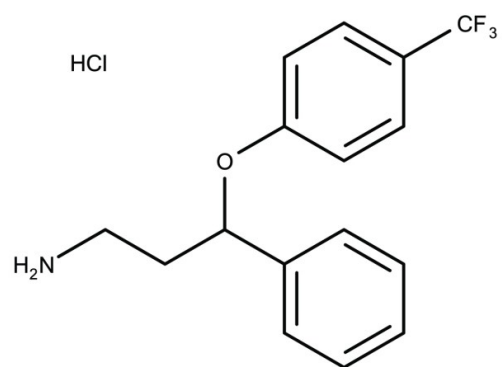
Ordering Information

[Order Online »](#)

ENZ-CHM269-0025	25mg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored unopened at -20°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Desmethylfluoxetine
Appearance	Beige waxy solid.
CAS	57226-68-3
Couple Target	Cytochrome P450
Couple Type	Inhibitor
Formula	$C_{16}H_{16}F_3NO \cdot HCl$
Identity	Identity determined by NMR.
MW	331.8
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (25 mg/ml).

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)