

Paroxetine . HCl

Highly potent and selective serotonin uptake inhibitor

Paroxetine is a highly potent and selective serotonin (5-HT) uptake inhibitor. It acts by binding to the serotonin transporter (SERT) with high affinity ($K_i = 05 \text{ nM}$). *In vitro* paroxetine competitively inhibits the uptake of 5-HT by rat hypothalamic and cortical synaptosomes with only very weak effects on noradrenaline and dopamine uptake. Paroxetine is in clinical use for the treatment of depression, anxiety and other disorders.

Citations: 4

[View Online »](#)

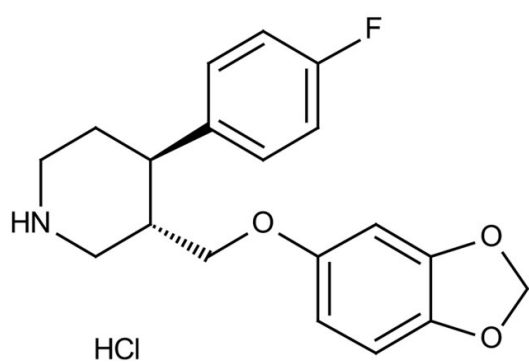
Ordering Information

[Order Online »](#)

BML-NS710-0100	100mg
----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Paxil®, Paroxetine . HCl . hemihydrate

Appearance White to off-white crystalline solid.

CAS 110429-35-1; 78246-49-8; 61869-08-7 (free base)

Couple Target Serotonin transporter

Couple Type Inhibitor

Formula $C_{19}H_{20}FNO_3 \cdot HCl \cdot 1/2 H_2O$

MW 374.8

Purity ≥98% (UHPLC)

Solubility Soluble in DMSO (25mg/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-
fr@enzolifesciences.com](mailto:info-fr@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)