

Bupropion . HCl

Inhibitor of dopamine and norepinephrine transporters

Inhibits dopamine and norepinephrine transporters ($K_i=2.8$ and $1.4\mu\text{M}$ respectively). Inhibits the serotonin transporter at higher concentration ($K_i=45\mu\text{M}$). In clinical use as an antidepressant and for smoking cessation. Also acts as a nicotinic receptor antagonist.

Citations: 5

[View Online »](#)

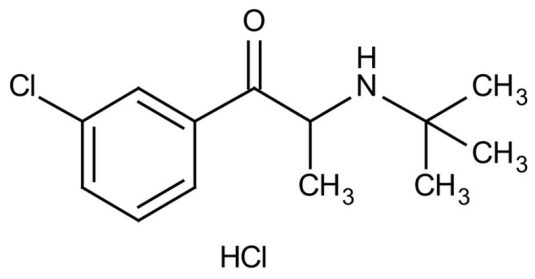
Ordering Information

[Order Online »](#)

BML-NS111-0050	50mg
BML-NS111-0250	250mg

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

Appearance	White solid.
CAS	31677-93-7
Couple Target	Dopamine transporter, Norepinephrine transporter
Couple Type	Inhibitor
Formula	$C_{13}H_{18}ClNO \cdot HCl$
MW	276.2
Purity	≥98% (HPLC)
Solubility	Soluble in water (>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-
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)