

N-Desmethyl Clozapine

Serotonin receptor ligand

Major metabolite of clozapine. Potent serotonin 5-HT_{1c} receptor antagonist.

Citations: 5

[View Online »](#)

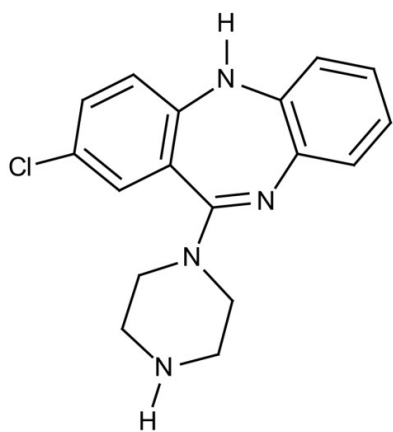
Ordering Information

[Order Online »](#)

BML-NS106-0025	25mg
----------------	------

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 Norclozapine

Appearance Yellow solid.

CAS 6104-71-8

Couple Target Serotonin receptor

Couple Type Ligand

Formula $C_{17}H_{17}ClN_4$

Identity Determined by ¹H-NMR.

MW 312.8

Purity ≥99% (HPLC)

Solubility Soluble in ethanol (25 mg/ml) or DMSO (>44 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-
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)