

Clozapine N-oxide

Major metabolite of clozapine

Clozapine N-oxide (CNO) is a major metabolite of clozapine (Prod. No. BML-D107). CNO activates designer receptors exclusively activated by designer drugs (DREADDs) that are derived from human muscarinic acetylcholine receptors (M₁₋₅D receptors).

Clozapine is an antipsychotic medication used in the treatment of schizophrenia, and is also sometimes used off-label for the treatment of bipolar disorder and borderline personality disorder. It is on the World Health Organization's List of Essential Medicines, a list of the most important medication needed in a basic health system.

Citations: 159

[View Online »](#)

Ordering Information

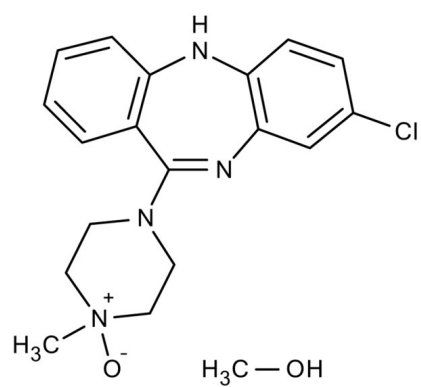
[Order Online »](#)

BML-NS105-0005	5mg
BML-NS105-0025	25mg

Manuals, SDS & CofA

[View Online »](#)

- High purity
- Active *in vitro* and *in vivo*
- Possesses little-to-no activity at 5-HT receptors



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 CNO

Appearance Yellow solid.

CAS 34233-69-7

Formula $C_{18}H_{19}ClN_4O$

Formulation Contains 1 equivalent of methanol.

MW 342.8

Purity $\geq 99\%$ (HPLC)

Solubility Soluble in water (5mg/ml).

Source Synthetic.

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)