

# 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 (M<sub>1-5</sub>D 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

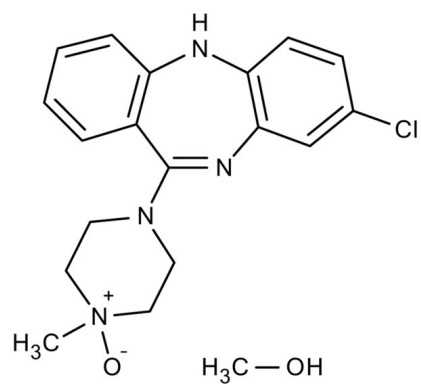
[Order Online »](#)

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-  
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)