

Clozapine

Dopamine receptor and serotonin receptor antagonist

Dopamine D₄ and D₂ receptor antagonist. High affinity for the cloned rat dopamine D4 receptor (K_i<20nM). Atypical neuroleptic agent. Antagonist at 5HT_{2A}, 5HT_{2C}, 5HT₃, 5HT₆ and 5HT₇ receptors.

Citations: 10

[View Online »](#)

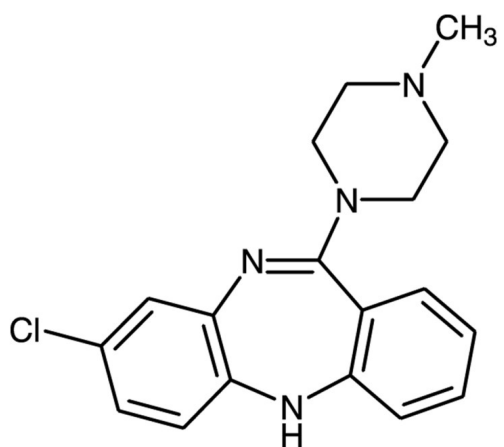
Ordering Information

[Order Online »](#)

BML-D107-0500	500mg
---------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 8-Chloro-11-(4-methyl-1-piperazinyl)-5H-dibenzo[b,e]-[1,4]diazepine

Appearance Yellow crystalline powder.

CAS 5786-21-0

Couple Target Dopamine receptor, Serotonin receptor

Couple Type Ligand

Formula $C_{18}H_{19}ClN_4$

Identity Determined by IR.

MI 14: 2423

MW 326.8

Purity ≥99% (HPLC)

RTECS HP1750000

Solubility Soluble in DMSO (30mg/ml) or dilute aqueous acid.

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)