

Clozapine N-oxide

Inert clozapine analog, activates specialized GPCRs.

Ordering Information

[Order Online »](#)

LKT-C4759-M005	5mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name CNO

CAS 34233-69-7

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

MW 342.82

Purity $\geq 98\%$

Solubility Water to 10 mM, DMSO to 10 mg/mL, acetone, ether, methanol.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com