

PNU-120596

Nicotinic acetylcholine receptor activator

Type II positive allosteric modulator (PAM) of $\alpha 7$ neuronal nicotinic acetylcholine receptor (nAChR) (EC_{50} =216nM). Shows no detectable effect on $\alpha 4\beta 2$, $\alpha 3\beta 4$ and $\alpha 9\alpha 10$ receptors.

Citations: 3

[View Online »](#)

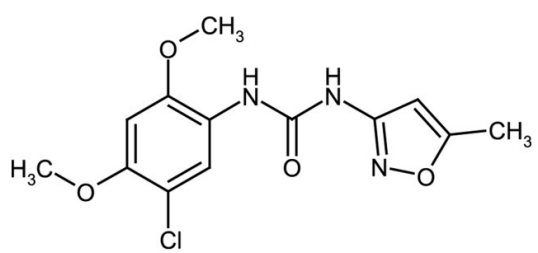
Ordering Information

[Order Online »](#)

ALX-550-526-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name NSC 216666, N-(5-Chloro-2,4-dimethoxyphenyl)-N'-(5-methyl-3-isoxazolyl)urea

Appearance White to off-white solid.

CAS 501925-31-1

Couple Target Acetylcholine receptor

Couple Type Activator

Formula $C_{13}H_{14}ClN_3O_4$

Identity Determined by MS and NMR.

MW 311.7

Purity ≥98% (HPLC)

Solubility Soluble in DMSO (50mg/ml); slightly soluble in methanol (1mg/ml) or acetonitrile (0.5mg/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)