

trans-Ned-19

Cell permeable, selective and potent inhibitor of nicotinic acid adenine dinucleotide phosphate (NAADP) receptor in mammalian cells and sea urchin eggs. Blocks NAADP signaling. Inhibits Ca^{2+} release: *trans*-Ned-19 ($\text{IC}_{50} = 6\text{nM}$); *cis*-Ned-19 ($\text{IC}_{50} = 800\text{nM}$). [^{32}P] NAADP binding: *trans*-Ned-19 ($\text{IC}_{50} = 0.4\text{nM}$); *cis*-Ned-19 ($\text{IC}_{50} = 15\mu\text{M}$). Inhibits Ca^{2+} release from lysosomes. Noncompetitive antagonist. Fluorescent.

NAADP (nicotinic acid adenine dinucleotide phosphate) is an emerging Ca^{2+} releasing second messenger in several mammalian tissues including pancreatic, brain and cardiac tissue.

Citations: 28

[View Online »](#)

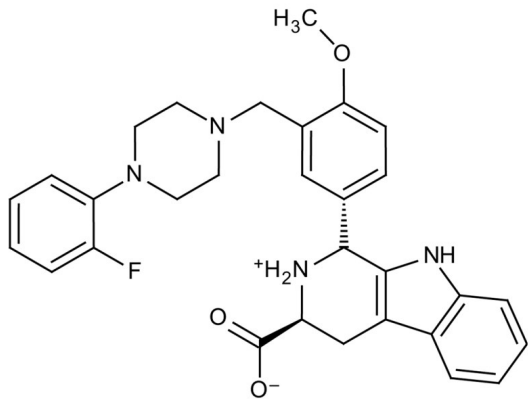
Ordering Information

[Order Online »](#)

ALX-270-503-MC05	0.5mg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	1-(3-((4-(2-Fluorophenyl)piperazin-1-yl)methyl)-4-methoxyphenyl)-2,3,4,9-tetrahydro-1H-pyrido[3,4-b]indole-3-carboxylic acid
Appearance	White to off-white powder.
Couple Target	NAADP receptor
Couple Type	Inhibitor
Formula	$C_{30}H_{31}FN_4O_3$
Identity	Determined by 1H-NMR and MS.
MW	514.6
Purity	≥97% (HPLC)
Solubility	Soluble in DMSO or 100% ethanol.
Source	Synthetic
Technical Info / Product Notes	Fluorescence of Ned-19 is compatible with visualization of the NAADP receptor in intact cells via confocal microscopy using an ultraviolet Argon ion laser (excitation 351 and 365 nm).



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-fr@enzolifesciences.com

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

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

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

Last modified: May 29, 2024