

# 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  $\text{Ca}^{2+}$  release: *trans*-Ned-19 ( $\text{IC}_{50} = 6\text{nM}$ ); *cis*-Ned-19 ( $\text{IC}_{50} = 800\text{nM}$ ). [ $^{32}\text{P}$ ] NAADP binding: *trans*-Ned-19 ( $\text{IC}_{50} = 0.4\text{nM}$ ); *cis*-Ned-19 ( $\text{IC}_{50} = 15\mu\text{M}$ ). Inhibits  $\text{Ca}^{2+}$  release from lysosomes. Noncompetitive antagonist. Fluorescent.

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

Citations: 28

[View Online »](#)

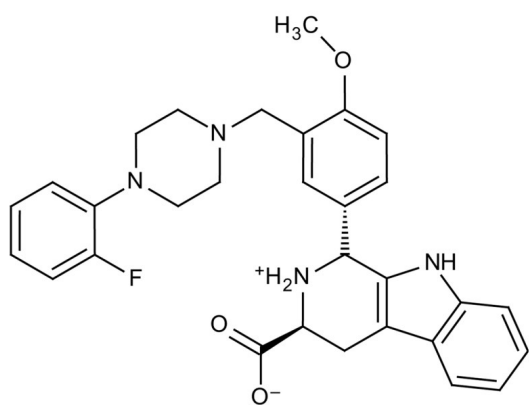
## Ordering Information

[Order Online »](#)

ALX-270-503-MC05	0.5mg
ALX-270-503-M001	1mg

## 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](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)

Last modified: May 29, 2024