

# Propenyl-L-NIO . hydrochloride

## iNOS inhibitor

Potent, selective inhibitor of iNOS (NOS II). The  $K_i$  values for inhibition of iNOS (NOS II), nNOS (NOS I), and eNOS (NOS III) by ENIPO are 17, 10.3 and 58.2  $\mu$ M, respectively, as determined using initial rate binding kinetics. However, ENIPO is a time-dependent inhibitor that, with longer incubations, demonstrates reversible tight binding inhibition that is selective for iNOS (NOS II) over nNOS (NOS I) or eNOS (NOS III). This selectivity results from iNOS (NOS II) exhibiting a 4-fold faster binding and 10-fold slower dissociation with ENIPO compared to nNOS (NOS I).  $K_i$  values for ENIPO binding to iNOS (NOS II) and nNOS (NOS I) are 0.56 and 6  $\mu$ M, respectively. Accurate determination of the  $K_i$  value of ENIPO for eNOS (NOS III) has not been determined but is expected to be 50-fold less than iNOS (NOS II).

Citations: 1

[View Online »](#)

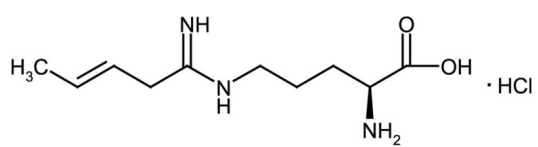
## Ordering Information

[Order Online »](#)

ALX-270-497-M025	25mg
------------------	------

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 2 years after receipt when stored at -20°C. We do not recommend storing aqueous solutions for more than one day.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	N5-[(3E)-1-Imino-3-pentenyl]-L-ornithine . HCl, ENIPO . HCl
Appearance	White to off-white crystalline powder.
Couple Target	NOS
Couple Type	Inhibitor
Formula	$C_{10}H_{19}N_3O_2 \cdot HCl$
Identity	Determined by IR, MS and <sup>1</sup> H-NMR.
MW	213.3 . 36.5
Purity	≥97% (TLC)
Solubility	Soluble in PBS, pH 7.2 (10mg/ml), 100% ethanol (10mg/ml), DMSO (2mg/ml) or dimethyl formamide (2mg/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)