

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 μM , 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 μM , 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

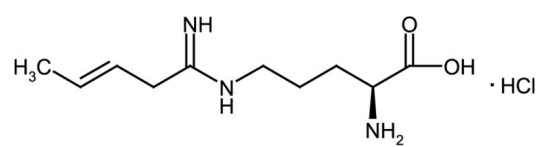
[Order Online »](#)

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 1H-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

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