

N^G-Hydroxy-L-arginine . monoacetate

Intermediate in NO biosynthesis

Intermediate in the biosynthesis of nitric oxide from L-arginine by nitric oxide synthases (NOS). Inhibitor of arginases.

Citations: 15

[View Online »](#)

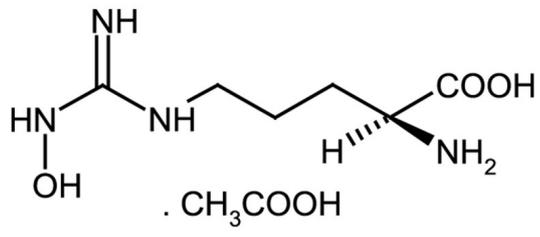
Ordering Information

[Order Online »](#)

| | |
|------------------|------|
| ALX-106-004-M025 | 25mg |
|------------------|------|

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

| | |
|-------------------|--|
| Use/Stability | As indicated on product label or CoA when stored as recommended. |
| Long Term Storage | +4°C |
| Shipping | Ambient Temperature |

Regulatory Status

RUO - Research Use Only

Product Details

| | |
|------------------|----------------------------------|
| Alternative Name | L-HOArg . AcOH |
| Appearance | White to off-white solid. |
| CAS | 53598-01-9 |
| Couple Target | Arginase, NOS |
| Couple Type | Inhibitor, Substrate |
| Formula | $C_6H_{14}N_4O_3 \cdot CH_3COOH$ |
| MW | 190.2 . 60.1 |
| Purity | ≥98% (HPLC) |
| Solubility | Soluble in water (50mg/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-
be@enzolifesciences.com](mailto:info-be@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)