

# NG, NG-Dimethyl-L-arginine . dihydrochloride

## NOS inhibitor

Endogenous nitric oxide synthase (NOS) inhibitor.

Citations: 1

[View Online »](#)

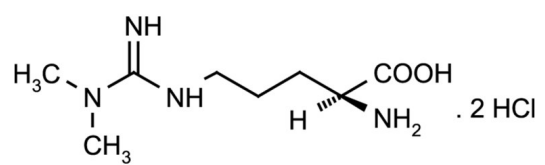
## Ordering Information

[Order Online »](#)

|                  |      |
|------------------|------|
| ALX-106-006-M005 | 5mg  |
| ALX-106-006-M025 | 25mg |

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

|                          |  |
|--------------------------|--|
| <b>Use/Stability</b>     | As indicated on product label or CoA when stored as recommended. Stock solutions are stable for up to 4 months when stored at -20°C. |
| <b>Handling</b>          | Hygroscopic.   |
| <b>Long Term Storage</b> | -20°C  |
| <b>Shipping</b>          | Ambient Temperature  |

## Regulatory Status

RUO - Research Use Only

## Product Details

|                         |  |
|-------------------------|--|
| <b>Alternative Name</b> | H-Arg(di-Me)-OH . 2HCl (asymmetrical), ADMA . 2HCl |
| <b>Appearance</b>       | White to off-white crystalline solid.              |
| <b>CAS</b>              | 220805-22-1  |
| <b>Couple Target</b>    | NOS  |
| <b>Couple Type</b>      | Inhibitor  |
| <b>Formula</b>          | $C_8H_{18}N_4O_2 \cdot 2 HCl \cdot H_2O$           |
| <b>MW</b>               | 202.2 . 73.0 . 18.0                                |
| <b>Purity</b>           | ≥97% (TLC)   |
| <b>Solubility</b>       | 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-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)