

L-Thiocitrulline . dihydrochloride

NOS inhibitor

Inhibitor of neuronal (nNOS; NOS I; $K_i=60\text{nM}$) and inducible (iNOS; NOS II) nitric oxide synthase.

Citations: 2

[View Online »](#)

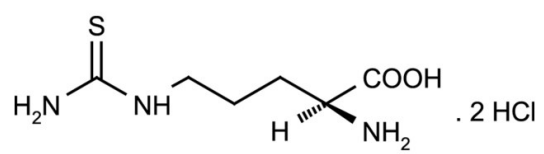
Ordering Information

[Order Online »](#)

ALX-106-013-M010	10mg
ALX-106-013-M050	50mg

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

Appearance White solid.

CAS 212051-53-1

Couple Target NOS

Couple Type Inhibitor

Formula $C_6H_{13}N_3O_2S \cdot 2HCl$

MW 191.3 . 72.9

Purity ≥97% (TLC)

Solubility Soluble in water or DMSO.

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)