

Canavanine . sulfate

iNOS inhibitor

The non-protein amino acid L-canavanine is an analog of L-arginine. Selective inducible nitric oxide synthase (iNOS; NOS II) inhibitor. Induces apoptotic cell death and shows antiproliferative and immunotoxic effects.

Citations: 15

[View Online »](#)

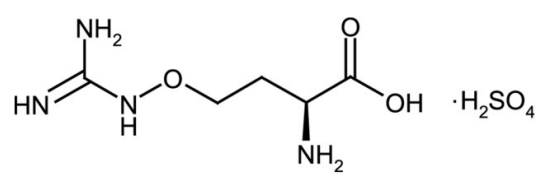
Ordering Information

[Order Online »](#)

ALX-350-002-M100	100mg
ALX-350-002-M500	500mg

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- α -Amino- γ -(guanidinoxy)-n-butyric acid . sulfate

Appearance White to-off white powder.

CAS 2219-31-0

Couple Target NOS

Couple Type Inhibitor

Formula $C_5H_{12}N_4O_3 \cdot H_2SO_4$

Identity Identity determined by IR.

MI **14:** 1737

MW 176.2 . 98.1

Purity $\geq 99\%$ (TLC)

Solubility Soluble in water (100mg/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)