

L-Mimosine

Cell cycle blocker

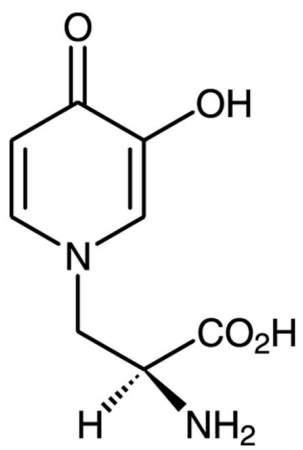
A rare plant amino acid that reversibly blocks the cell cycle at G1-phase. Inhibits the hydroxylation of deoxyhypusine and thus inhibits eukaryote initiation factor 5-A (eIF5-A). Inhibitor of mammalian tyrosinase and dopamine β -hydroxylase. Also chelates iron *via* the kojic acid metal-binding domain. Arrests DNA synthesis by inhibiting deoxyribonucleotide metabolism.

Citations: 5

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance Pink to brown powder.

CAS 500-44-7

Couple Target Dopamine beta-hydroxylase, eIF, Tyrosinase

Couple Type Inhibitor

Formula $C_8H_{10}N_2O_4$

Identity Determined by 1H-NMR.

MW 198.2

Purity $\geq 98\%$

Solubility Soluble in 10% sodium bicarbonate (15mg/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)