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 »

Ordering Information

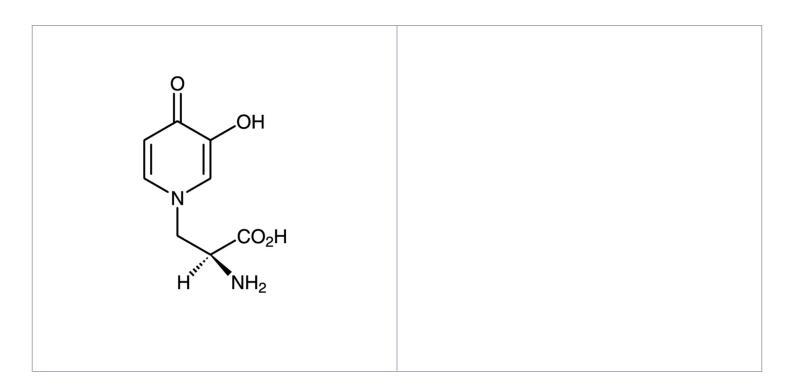
Order Online »

BML-CC102-0500

500mg

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 ≥98%

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

Last modified: May 29, 2024