

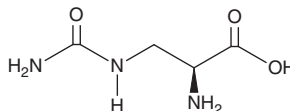
PRODUCT INFORMATION



L-Albizziin

Item No. 36665

CAS Registry No.: 1483-07-4
Formal Name: 3-[(aminocarbonyl)amino]-L-alanine
Synonyms: NSC 132089,
L- α -amino- β -Ureidopropionic Acid,
L- β -Ureidoalanine
MF: $C_4H_9N_3O_3$
FW: 147.1
Purity: $\geq 95\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

L-Albizziin is supplied as a solid. Aqueous solutions of L-albizziin can be prepared by directly dissolving the solid in aqueous buffers. The solubility of L-albizziin in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

L-Albizziin is a non-proteinogenic amino acid and metabolite of 5-aminouracil.¹ It is formed from 5-aminouracil via a 5-amino-5,6-dihydrouracil intermediate by dihydropyrimidinase.¹ L-Albizziin inhibits 4-methyleneglutaminase ($IC_{50} = 20$ mM for the peanut enzyme).²

References

1. Brown, E.G. and Turan, Y. Pyrimidine metabolism and secondary product formation; biogenesis of albizziine, 4-hydroxyhomoarginine and 2,3-diaminopropanoic acid. *Phytochem.* **40(3)**, 763-771 (1995).
2. Ibrahim, S.A., Lea, P.J., and Fowden, S.L. Purification and properties of 4-methyleneglutaminase from the leaves of peanut (*Arachis hypogaea*). *Phytochem.* **23(8)**, 1545-1549 (1984).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 08/29/2023

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM