

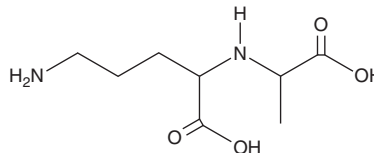
PRODUCT INFORMATION



Octopinic Acid

Item No. 33093

CAS Registry No.: 20197-09-5
Formal Name: N²-[(1R)-1-carboxyethyl]-L-ornithine
Synonym: (+)-Octopinic Acid
MF: C₈H₁₆N₂O₄
FW: 204.2
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



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

Laboratory Procedures

Octopinic acid is supplied as a solid. A stock solution may be made by dissolving the octopinic acid in the solvent of choice, which should be purged with an inert gas. Octopinic acid is soluble in methanol and DMSO.

Description

Octopinic acid is an opine that has been found in crown gall tumors induced by the plant pathogenic bacterium *A. tumefaciens*.^{1,2}

References

1. Scott, I.M. Opine content of unorganised and teratomatous tobacco crown gall tissues. *Plant Sci. Lett.* **16(2-3)**, 239-248 (1979).
2. Menage, A. and Morel, G. Presence of a new amino acid in crown gall tissue. *C. R. Acad. Sci.* **261(9)**, 2001-2002 (1965).

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, 11/11/2022

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