

# PRODUCT INFORMATION



## Allocryptopine

Item No. 34751

**CAS Registry No.:** 485-91-6  
**Formal Name:** 5,7,8,15-tetrahydro-3,4-dimethoxy-6-methyl-[1,3]benzodioxolo[5,6-e][2]benzazecin-14(6H)-one

**Synonyms:**  $\alpha$ -Allocryptopine,  $\alpha$ -Fagarine,  $\beta$ -Homochelidonine

**MF:** C<sub>21</sub>H<sub>23</sub>NO<sub>5</sub>

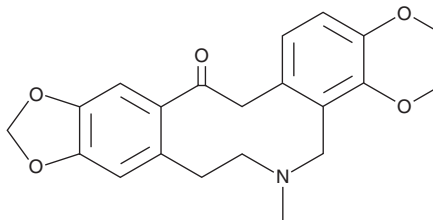
**FW:** 369.4

**Purity:**  $\geq$ 95%

**Supplied as:** A solid

**Storage:** -20°C

**Stability:**  $\geq$ 4 years



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

### Laboratory Procedures

Allocryptopine is supplied as a solid. A stock solution may be made by dissolving the allocryptopine in the solvent of choice, which should be purged with an inert gas. Allocryptopine is soluble in ethanol.

### Description

Allocryptopine is an isoquinoline alkaloid that has been found in *M. cordata* and has antimicrobial activity.<sup>1-3</sup> It is active against *S. aureus*, *P. aeruginosa*, and *E. coli* (MICs = 250, 125, and 125  $\mu$ g/ml, respectively), the protozoa *P. falciparum* and *E. histolytica* (IC<sub>50</sub>s = 5.08 and 135  $\mu$ M, respectively), as well as a panel of seven phytopathogenic fungi, including *F. graminearum* and *F. oxysporum*, with IC<sub>50</sub> values ranging from 15.2 to 42.4  $\mu$ g/ml.

### References

1. Kosina, P., Gregorova, J., Gruz, J., *et al.* Phytochemical and antimicrobial characterization of *Macleaya cordata* herb. *Fitoterapia* **81(8)**, 1006-1012 (2010).
2. Wright, C.W., Marshall, S.J., Russell, P.F., *et al.* In vitro antiplasmodial, antiamebic, and cytotoxic activities of some monomeric isoquinoline alkaloids. *J. Nat. Prod.* **63(12)**, 1638-1640 (2000).
3. Liu, H., Wang, J., Zhao, J., *et al.* Isoquinoline alkaloids from *Macleaya cordata* active against plant microbial pathogens. *Nat. Prod. Commun.* **4(11)**, 1557-1560 (2009).

#### 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, 11/04/2022

#### 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