

# PRODUCT INFORMATION



## (S)-Canadine

Item No. 33805

**CAS Registry No.:** 5096-57-1  
**Formal Name:** 5,8,13,13aS-tetrahydro-9,10-dimethoxy-6H-benzo[g]-1,3-benzodioxolo[5,6-a]quinolizine

**Synonyms:** (-)-Canadine,  
(S)-Tetrahydroberberine

**MF:** C<sub>20</sub>H<sub>21</sub>NO<sub>4</sub>

**FW:** 339.4

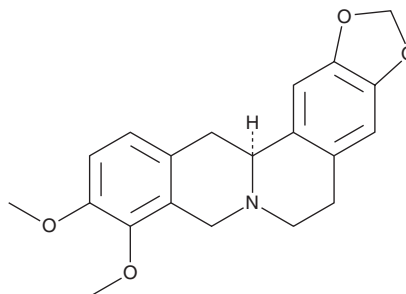
**Purity:** ≥98%

**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

(S)-Canadine is supplied as a solid. A stock solution may be made by dissolving the (S)-canadine in the solvent of choice, which should be purged with an inert gas. (S)-Canadine is slightly soluble in chloroform and is very slightly soluble in methanol (warmed).

### Description

(S)-Canadine is an alkaloid and intermediate in the biosynthesis of berberine (Item No. 10006427) that has been found in *H. canadensis*.<sup>1-3</sup>

### References

1. Malhotra, S., Taneja, S.C., and Dhar, K.L. Minor alkaloid from *Coscinium fenestratum*. *Phytochem.* **28(7)**, 1998-1999 (1989).
2. Galanie, S., and Smolke, C.D. Optimization of yeast-based production of medicinal protoberberine alkaloids. *Microb. Cell Fact.* **14**, 144 (2015).
3. Leyte-Lugo, M., Britton, E.R., Foil, D.H., *et al.* Secondary metabolites from the leaves of the medicinal plant goldenseal (*Hydrastis canadensis*). *Phytochem. Lett.* **20**, 54-60 (2017).

#### 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, 08/25/2023

#### 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