

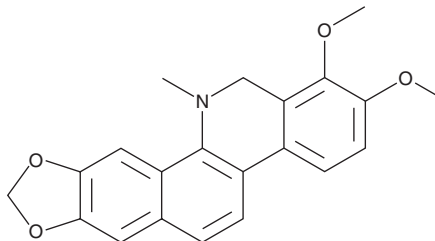
# PRODUCT INFORMATION



## Dihydrochelerythrine

Item No. 30118

**CAS Registry No.:** 6880-91-7  
**Formal Name:** 12,13-dihydro-1,2-dimethoxy-12-methyl-[1,3]benzodioxolo[5,6-c]phenanthridine  
**MF:** C<sub>21</sub>H<sub>19</sub>NO<sub>4</sub>  
**FW:** 349.4  
**Purity:** ≥95%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥2 years  
**Item Origin:** Plant/*Macleaya cordata* (Willd.) R. Br.



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

### Laboratory Procedures

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

### Description

Dihydrochelerythrine is a benzophenanthridine alkaloid that has been found in *Bocconia arborea* and has diverse biological activities.<sup>1,2,3</sup> It is active against *S. aureus*, *S. faecalis*, and *C. albicans* (MIC = 1.87 µg/ml for all).<sup>1</sup> Dihydrochelerythrine inhibits secretion of the hepatitis B virus (HBV) antigens HBsAg and HBeAg from HepG2 2.2.15 cells (IC<sub>50</sub>s = <0.05 µM for both).<sup>2</sup> It is cytotoxic to HCT8, BGC-823, and A2780 cancer cells (IC<sub>50</sub>s = 1.4, 0.4, and 3.5 µM, respectively).<sup>3</sup>

### References

1. Navarro, V. and Delgado, G. Two antimicrobial alkaloids from *Bocconia arborea*. *J. Ethnopharmacol.* **66(2)**, 223-226 (1999).
2. Wu, Y.R., Ma, Y.B., Zhao, Y.X., *et al.* Two new quaternary alkaloids and anti-hepatitis B virus active constituents from *Corydalis saxicol*. *Planta Med.* **73(8)**, 787-791 (2007).
3. Deng, A.J. and Qin, H.L. Cytotoxic dihydrobenzophenanthridine alkaloids from the roots of *Macleaya microcarp*. *Phytochemistry* **71(7)**, 816-822 (2010).

#### 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, 05/18/2020

#### 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