

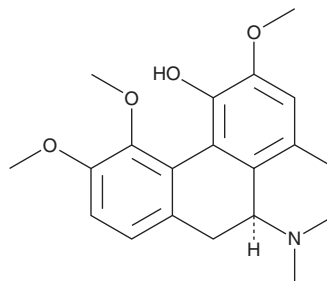
PRODUCT INFORMATION



Corydine

Item No. 35028

CAS Registry No.: 476-69-7
Formal Name: 5,6,6aS,7-tetrahydro-2,10,11-trimethoxy-6-methyl-4H-dibenzo[de,g]quinolin-1-ol
Synonym: (+)-Corydine
MF: $C_{20}H_{23}NO_4$
FW: 341.4
Purity: $\geq 95\%$
UV/Vis.: λ_{\max} : 224 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Plant/*Corydalis edulis* Maxim.



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

Laboratory Procedures

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

Description

Corydine is an aporphine alkaloid that has been found in *Corydalis* and has diverse biological activities.¹⁻³ It inhibits the proliferation of murine L1210 lymphocytic leukemia, B16/F10 melanoma, and P388 leukemia cells when used at concentrations ranging from 10 to 100 $\mu\text{g}/\text{ml}$.¹ Corydine binds to μ -opioid receptors ($K_i = 2.82 \mu\text{M}$) and induces [^{35}S]GTP γ S binding in CHO cells expressing human μ -opioid receptors ($\text{EC}_{50} = 0.51 \mu\text{M}$).² It also is active against *L. donovani* promastigotes ($\text{IC}_{50} = 26.7 \mu\text{M}$).³

References

1. Kondo, Y., Imai, Y., Hojo, H., *et al.* Suppression of tumor cell growth and mitogen response by aporphine alkaloids, dicentrine, glaucine, corydine, and apomorphine. *J. Pharmacobiodyn.* **13**(7), 426-431 (1990).
2. Kaserer, T., Steinacher, T., Kainhofer, R., *et al.* Identification and characterization of plant-derived alkaloids, corydine and corydaline, as novel mu opioid receptor agonists. *Sci. Rep.* **10**(1), 13804 (2020).
3. del Rayo Camacho, M., Kirby, G.C., Warhurst, D.C., *et al.* Oxoaporphine alkaloids and quinones from *Stephania dinklagei* and evaluation of their antiprotozoal activities. *Planta Med.* **66**(5), 478-480 (2000).

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, 12/08/2022

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