

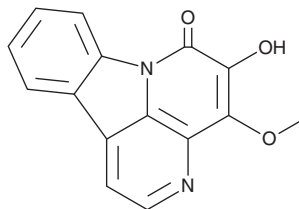
PRODUCT INFORMATION



4-methoxy-5-Hydroxycanthin-6-one

Item No. 40234

CAS Registry No.: 18110-86-6
Formal Name: 5-hydroxy-4-methoxy-6H-indolo[3,2,1-de][1,5] naphthyridin-6-one
Synonym: Nigakinone
MF: C₁₅H₁₀N₂O₃
FW: 266.3
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

4-methoxy-5-Hydroxycanthin-6-one is supplied as a solid. A stock solution may be made by dissolving the 4-methoxy-5-hydroxycanthin-6-one in the solvent of choice, which should be purged with an inert gas. 4-methoxy-5-Hydroxycanthin-6-one is soluble in methanol.

Description

4-methoxy-5-Hydroxycanthin-6-one is an alkaloid that has been found in *P. quassioides* and has anti-inflammatory activity.^{1,2} It reduces LPS-induced TNF- α release from, and nitric oxide (NO) production in, RAW 264.7 macrophages when used at concentrations of 10, 30, and 100 μ M.¹ *In vivo*, 4-methoxy-5-hydroxycanthin-6-one (9 and 27 mg/kg) reduces paw swelling in a rat model of arthritis induced by complete Freund's adjuvant (CFA). It also reduces colonic inflammatory cytokine levels and maintains mucosal barrier function in wild-type, but not *Fxr*^{-/-}, mice in a model of colitis induced by the sulfated polysaccharide dextran sulfate sodium (DSS; Item No. 23250) when administered at doses ranging from 25 to 100 mg/kg.²

References

1. Fan, H., Qi, D., Yang, M., *et al.* *In vitro* and *in vivo* anti-inflammatory effects of 4-methoxy-5-hydroxycanthin-6-one, a natural alkaloid from *Picrasma quassioides*. *Phytomedicine* **20**(3-4), 319-323 (2013).
2. Liu, F., Yao, Y., Wang, Q., *et al.* Nigakinone alleviates DSS-induced experimental colitis *via* regulating bile acid profile and FXR/NLRP3 signaling pathways. *Phytother. Res.* **37**(1), 15-34 (2023).

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, 02/13/2024

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