

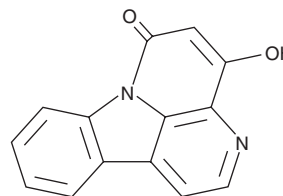
PRODUCT INFORMATION



4-Hydroxycanthin-6-one

Item No. 31646

CAS Registry No.: 106941-27-9
Formal Name: 4-hydroxy-6H-indolo[3,2,1-de][1,5]naphthyridin-6-one
Synonym: 4-Hydroxycanthinone
MF: C₁₄H₈N₂O₂
FW: 236.2
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Pentaceras australis*



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

Laboratory Procedures

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

Description

4-Hydroxycanthin-6-one is an alkaloid that has been found in *A. altissima*.¹

Reference

1. Kim, H.M., Lee, J.S., Sezirahiga, J., *et al.* A new canthinone-type alkaloid isolated from *Ailanthus altissima* swingle. *Molecules* **21(5)**, 642 (2016).

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, 12/05/2023

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