

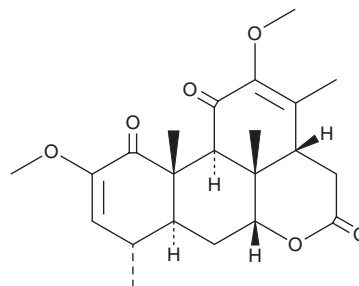
PRODUCT INFORMATION



Quassin

Item No. 34382

CAS Registry No.: 76-78-8
Formal Name: 2,12-dimethoxy-picrasa-2,12-diene-1,11,16-trione
Synonyms: NSC 36342, (+)-Quassin
MF: C₂₂H₂₈O₆
FW: 388.5
Purity: ≥98%
UV/Vis.: λ_{max}: 257 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Quassia amara*



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

Laboratory Procedures

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

Description

Quassin is a quassinoid that has been found in *Q. amara* and has diverse biological activities.¹⁻³ It is active against chloroquine-sensitive and -resistant strains of *P. falciparum* (IC₅₀ = 0.15 μM for both).¹ Quassin has antifeedant and insecticidal activity against the diamondback moth (*P. xylostella*) when applied to dietary cabbage leaf discs at 16, 31.9, or 63.7 μg/cm².² It increases viability of SH-SY5Y cells in a model of hydrogen peroxide-induced neuronal injury when used at concentrations of 25, 50, or 100 μM.³

References

1. Mishra, K., Chakraborty, D., Pal, A., et al. *Plasmodium falciparum*: in vitro interaction of quassin and neo-quassin with artesunate, a hemisuccinate derivative of artemisinin. *Exp. Parasitol.* **124(4)**, 421-427 (2010).
2. Daido, M., Fukamiya, N., Okano, M., et al. Antifeedant and insecticidal activity of quassinoids against diamondback moth (*Plutella xylostella*). *Biosci. Biotechnol. Biochem.* **57(2)**, 2440246 (1993).
3. He, C., Wang, Y., Yang, T., et al. Quassinoids with insecticidal activity against *Diaphorina citri* Kuwayama and neuroprotective activities from *Picrasma quassioides*. *J. Agric. Food. Chem.* **68(1)**, 117-127 (2020).

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