

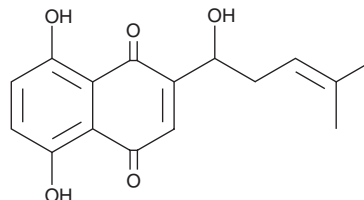
PRODUCT INFORMATION



Shikalkin

Item No. 34355

CAS Registry No.: 54952-43-1
Formal Name: 5,8-dihydroxy-2-(1-hydroxy-4-methyl-3-penten-1-yl)-1,4-naphthalenedione
Synonyms: (±)-Alkannin, (±)-Shikonin, NSC 291845
MF: C₁₆H₁₆O₅
FW: 288.3
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Arnebia euchroma*



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

Laboratory Procedures

Shikalkin is supplied as a solid. A stock solution may be made by dissolving the shikalkin in the solvent of choice, which should be purged with an inert gas. Shikalkin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of shikalkin in these solvents is approximately 1, 5, and 10 mg/ml, respectively.

Description

Shikalkin is a naphthoquinone and racemic mixture of alkannin (Item No. 25058) and shikonin (Item No. 14751) that has been found in *A. hispidissima*.¹

Reference

1. Singh, B., Sharma, M.K., Meghwal, P.R., *et al.* Anti-inflammatory activity of shikonin derivatives from *Arnebia hispidissima*. *Phytomedicine* **10**(5), 375-380 (2003).

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, 11/22/2022

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