

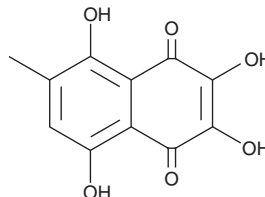
PRODUCT INFORMATION



Methylspinazarin

Item No. 29898

CAS Registry No.: 41768-12-1
Formal Name: 2,3,5,8-tetrahydroxy-6-methyl-1,4-naphthalenedione
Synonym: Methylspinazaline
MF: C₁₁H₈O₆
FW: 236.2
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Bacterium/*Streptomyces* sp.



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

Laboratory Procedures

Methylspinazarin is supplied as a solid. A stock solution may be made by dissolving the methylspinazarin in the solvent of choice, which should be purged with an inert gas. Methylspinazarin is soluble in ethanol, methanol, DMSO, and dimethyl formamide.

Description

Methylspinazarin is a naphthoquinone bacterial metabolite that has been found in *Streptomyces* and is an inhibitor of catechol O-methyltransferase (COMT; IC₅₀ = 0.8 µg/ml).¹ It is selective for COMT over tyrosine hydroxylase, DOPA decarboxylase, and dopamine-β-hydroxylase at 100 µg/ml. Methylspinazarin decreases blood pressure in spontaneously hypertensive rats when administered at a dose of 50 mg/kg.

Reference

1. Chimura, H., Sawa, T., Takita, T., *et al.* Methylspinazarin and dihydromethylspinazarin, catechol-O-methyl transferase inhibitors produced by *Streptomyces*. *J. Antibiot.* **26(2)**, 112-114 (1973).

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, 10/21/2022

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