

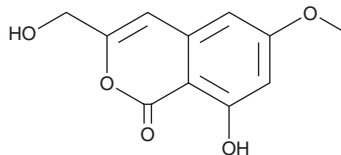
PRODUCT INFORMATION



Cytogenin

Item No. 40220

CAS Registry No.: 132971-59-6
Formal Name: 8-hydroxy-3-(hydroxymethyl)-6-methoxy-1H-2-benzopyran-1-one
MF: C₁₁H₁₀O₅
FW: 222.2
Purity: ≥98%
Supplied as: A powder
Storage: -20°C
Stability: ≥1 year
Item Origin: Bacterium/*Streptoverticillium eurocidicum*



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

Laboratory Procedures

Cytogenin is supplied as a powder. A stock solution may be made by dissolving the cytogenin in the solvent of choice, which should be purged with an inert gas. Cytogenin is soluble in acetone, acetonitrile, and DMSO.

Description

Cytogenin is a coumarin derivative originally found in *S. eurocidicum* that has diverse biological activities.^{1,2} It reduces tumor weight in an Ehrlich murine spontaneous adenocarcinoma model when administered at a dose of 25 mg/kg per day.¹ *In vivo*, cytogenin (30 and 100 mg/kg per day) reduces streptozotocin-induced decreases in body weight and increases in plasma glucose levels and macrophage pancreatic infiltration in a mouse model of diabetes.² It also decreases zymosan- and LPS-induced nitric oxide production, as well as LPS-induced increases in IL-6 levels, in isolated mouse macrophages after administration at a dose of 100 mg/kg per day.

References

1. Kumagai, H., Masuda, T., Ohsono, M., *et al.* Cytogenin, a novel antitumor substance. *J. Antibiot. (Tokyo)* **43(11)**, 1505-1507 (1990).
2. Kumagai, H., Wakazono, K., Agata, N., *et al.* Effect of cytogenin, a novel immunomodulator, on streptozotocin-induced diabetes in mice. *J. Antibiot. (Tokyo)* **58(3)**, 202-205 (2005).

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, 11/21/2023

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