

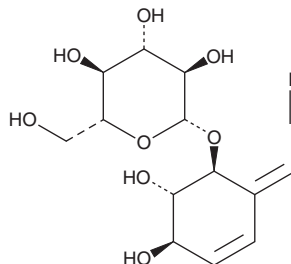
PRODUCT INFORMATION



Lithospermoside

Item No. 34771

CAS Registry No.: 63492-69-3
Formal Name: 2Z-[(4R,5S,6S)-6-(β-D-glucopyranosyloxy)-4,5-dihydroxy-2-cyclohexen-1-ylidene]-acetonitrile
Synonyms: Griffonin, NSC 290807
MF: C₁₄H₁₉NO₈
FW: 329.3
Purity: ≥98%
UV/Vis.: λ_{max}: 260 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Semiaquilegia adoxoides*



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

Laboratory Procedures

Lithospermoside is supplied as a solid. Aqueous solutions of lithospermoside can be prepared by directly dissolving the solid in aqueous buffers. The solubility of lithospermoside in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Lithospermoside is a nitrile glucoside that has been found in *Cowania mexicana* and has antitumor activity.¹ It inhibits activation of Epstein-Barr virus early antigen (EBV-EA) induced by phorbol 12-myristate 13-acetate (TPA; Item No. 10008014), a marker of tumor formation, in Raji cells.

Reference

1. Ito, H., Miyake, M., Nishatani, E., *et al.* Anti-tumor promoting activity of polyphenols from *Cowania mexicana* and *Coleogyne ramosissima*. *Cancer Lett.* **143(1)**, 5-13 (1999).

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