

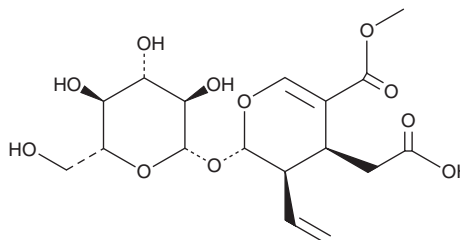
PRODUCT INFORMATION



Secoxyloganin

Item No. 30267

CAS Registry No.: 58822-47-2
Formal Name: (2S,3R,4S)-3-ethenyl-2-(β-D-glucopyranosyloxy)-3,4-dihydro-5-(methoxycarbonyl)-2H-pyran-4-acetic acid
MF: C₁₇H₂₄O₁₁
FW: 404.4
Purity: ≥95%
UV/Vis.: λ_{max}: 232 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years
Item Origin: Plant/*Lonicerae flos*



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

Laboratory Procedures

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

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of secoxyloganin can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of secoxyloganin in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Secoxyloganin is a secoiridoid glycoside that has been found in *L. japonica* and has diverse biological activities.¹⁻³ It is active against *E. coli* and *S. aureus* in a disc assay when used at a concentration of 2 mg/disc.¹ Secoxyloganin is cytotoxic to human dermal fibroblasts (IC₅₀ = 78.1 μM).² *In vivo*, secoxyloganin (10 mg/ml) prevents hen egg white lysozyme-induced decreases in tail vein blood flow, a marker of allergic inflammation, in hen egg white lysozyme-sensitized mice.³

References

1. Xiong, J., Li, S., Wang, W., *et al.* Screening and identification of the antibacterial bioactive compounds from *Lonicera japonica* Thunb. leaves. *Food Chem.* **138(1)**, 327-333 (2013).
2. Zebiri, I., Haddad, M., Duca, L., *et al.* Biological activities of triterpenoids from *Poraqueiba sericea* stems. *Nat. Prod. Res.* **31(11)**, 1333-1338 (2017).
3. Oku, H., Ogawa, Y., Iwaoka, E., *et al.* Allergy-preventive effects of chlorogenic acid and iridoid derivatives from flower buds of *Lonicera japonica*. *Biol. Pharm. Bull.* **34(8)**, 1330-1333 (2011).

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, 05/20/2020

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