

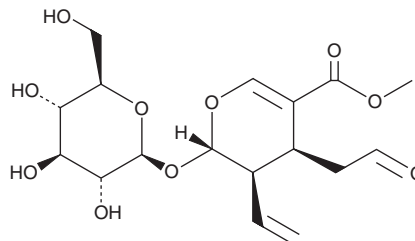
PRODUCT INFORMATION



Secologanin

Item No. 20647

CAS Registry No.: 19351-63-4
Formal Name: (2S,3R,4S)-3-ethenyl-2-(β-D-glucopyranosyloxy)-3,4-dihydro-4-(2-oxoethyl)-2H-pyran-5-carboxylic acid, methyl ester
Synonym: NSC 640525
MF: C₁₇H₂₄O₁₀
FW: 388.4
Purity: ≥90%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Secologanin is supplied as a solid. Secologanin is slightly soluble in aqueous solutions. We do not recommend storing the aqueous solution for more than one day.

Description

Secologanin is a monoterpene that has been found in *V. rosea* and is an intermediate in the synthesis of monoterpene indole alkaloids from geraniol.^{1,2} It is a metabolite of loganin (Item No. 19997) formed by the cytochrome P450 (CYP) isoform CYP72A1, also known as secologanin synthase.³

References

1. Battersby, A.R., Burnett, A.R., and Parsons, P.G. Alkaloid biosynthesis. XIV. Secologanin: Its conversion into ipecoside and its role as biological precursor of the indole alkaloids *J. Chem. Soc. C Organic* **8**, 1187-1192 (1969).
2. Miettinen, K., Dong, L., Navrot, N., *et al.* The seco-iridoid pathway from *Catharanthus roseus*. *Nat. Commun.* **5:3606**, (2014).
3. Irmeler, S., Schröder, G., St-Pierre, B., *et al.* Indole alkaloid biosynthesis in *Catharanthus roseus*: New enzyme activities and identification of cytochrome P450 CYP72A1 as secologanin synthase. *Plant J.* **24(6)**, 797-804 (2000).

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, 12/21/2022

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