

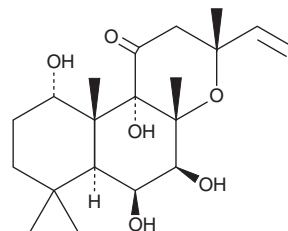
PRODUCT INFORMATION



Deacetylforskolin

Item No. 11701

CAS Registry No.: 64657-20-1
Formal Name: (3R,4aR,5S,6S,6aS,10S,10aR,10bS)-3-ethenyldodecahydro-5,6,10,10b-tetrahydroxy-3,4a,7,7,10a-pentamethyl-1H-naphtho[2,1-b]pyran-1-one
Synonyms: 7-Deacetylcoleonol, Forskolin D
MF: C₂₀H₃₂O₆
FW: 368.5
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Coleus forskohlii*



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

Laboratory Procedures

Deacetylforskolin is supplied as a solid. A stock solution may be made by dissolving the deacetylforskolin in the solvent of choice, which should be purged with an inert gas. Deacetylforskolin is slightly soluble in chloroform and methanol.

Description

Deacetylforskolin is a diterpene and a derivative of forskolin (Item No. 11018) that has been found in *C. forskohlii* and has diverse biological activities.¹⁻⁴ It activates rat adipocyte adenylyl cyclase (IC₅₀ = 20 μM) and inhibits glucose transport in rat adipocyte plasma membranes.³ Deacetylforskolin (30-1,000 μg/kg) reduces blood pressure in spontaneously hypertensive rats.⁴ It also attenuates hypercapnia-induced impairments in the passive avoidance response in mice.²

References

1. Gabetta, B., Zini, G., and Danieli, B. Minor Diterpenoids of *Coleus forskohlii*. *Phytochemistry* **28(3)**, 859-862 (1989).
2. McCulloch, A.J., Thomson, T.A., and Deacon, R. Hypoxic amnesia and its reversal with forskolin. *Biochem. Soc. Trans.* **17(1)**, 212-213 (1988).
3. Joost, H.G., Habberfield, A.D., Simpson, I.A., et al. Activation of adenylate cyclase and inhibition of glucose transport in rat adipocytes by forskolin analogues: structural determinants for distinct sites of action. *Mol. Pharmacol.* **33(4)**, 449-453 (1988).
4. Bhat, S.V., Dohadwalla, A.N., Bajwa, B.S., et al. The antihypertensive and positive inotropic diterpene forskolin: Effects of structural modifications on its activity. *J. Med. Chem.* **26(4)**, 486-492 (1983).

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, 11/07/2022

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