

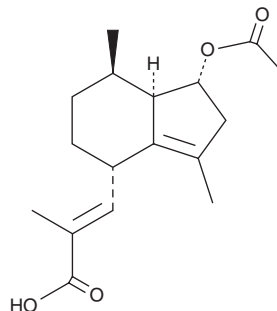
PRODUCT INFORMATION



Acetoxyvalerenic Acid

Item No. 31516

CAS Registry No.: 84638-55-1
Formal Name: (2E)-3-[(1R,4S,7R,7aR)-1-(acetyloxy)-2,4,5,6,7,7a-hexahydro-3,7-dimethyl-1H-inden-4-yl]-2-methyl-2-propenoic acid
Synonym: Acetylvalerenolic Acid
MF: C₁₇H₂₄O₄
FW: 292.4
Purity: ≥90%
UV/Vis.: λ_{max}: 217 nm
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years
Item Origin: Synthetic



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

Laboratory Procedures

Acetoxyvalerenic acid is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as methanol and chloroform purged with an inert gas can be used. Acetoxyvalerenic acid is slightly soluble in methanol and sparingly soluble in chloroform.

Description

Acetoxyvalerenic acid is a sesquiterpene acid that has been found in *Valeriana* roots.¹ It is cytotoxic to GLC-4 and COLO 320 cells (IC₅₀s = 111 and 124 μM, respectively).

Reference

1. Bos, R., Hendriks, H., Scheffer, J.J., *et al.* Cytotoxic potential of valerian constituents and valerian tinctures. *Phytomedicine* **5(3)**, 219-225 (1998).

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, 11/08/2022

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