

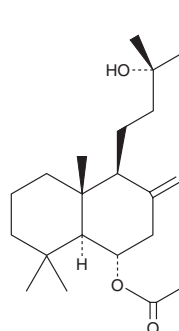
PRODUCT INFORMATION



Larixyl Acetate

Item No. 35309

CAS Registry No.: 4608-49-5
Formal Name: (α S,1S,4S,4aS,8aR)-4-(acetyloxy)- α -ethenyldecahydro- α ,5,5,8a-tetramethyl-2-methylene-1-naphthalenepropanol
Synonym: Larixol Acetate
MF: C₂₂H₃₆O₃
FW: 348.5
Purity: \geq 98%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Synthetic



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

Laboratory Procedures

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

Description

Larixyl acetate is a diterpenoid that has been found in *Larix* and has transient receptor potential canonical 6 (TRPC6) inhibitory and analgesic activities.^{1,2} It inhibits the activity of TRPC6 (IC₅₀ = 0.58 μ M) and is selective for TRPC6 over TRPC3 (IC₅₀ = 6.83 μ M), as well as over nine additional TRP channels at 10 μ M.¹ Intrathecal administration of larixyl acetate reduces mechanical and cold allodynia in a rat model of neuropathic pain induced by spared nerve injury (SNI).

References

1. Urban, N., Wang, L., Kwiek, S., *et al.* Identification and validation of larixyl acetate as a potent TRPC6 inhibitor. *Mol. Pharmacol.* **89**(1), 197-213 (2016).
2. Wang, J., Zhao, M., Jia, P., *et al.* The analgesic action of larixyl acetate, a potent TRPC6 inhibitor, in rat neuropathic pain model induced by spared nerve injury. *J. Neuroinflammation* **17**(1), 118 (2020).

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, 11/30/2022

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