

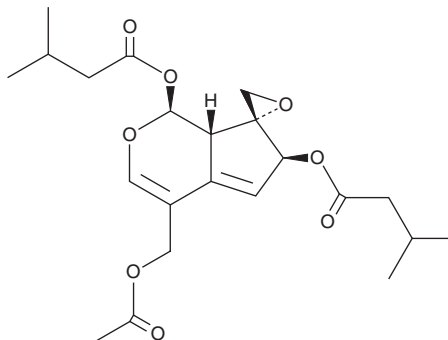
PRODUCT INFORMATION



Valtrate

Item No. 36522

CAS Registry No.: 18296-44-1
Formal Name: butanoic acid, 3-methyl-1'-
[(1S,2'R,6S,7aS)-4-[(acetyloxy)methyl]-
6,7a-dihydrospiro[cyclopenta[c]pyran-
7(1H),2'-oxirane]-1,6-diyl]ester
Synonym: Halazuchrome B
MF: C₂₂H₃₀O₈
FW: 422.5
Purity: ≥98%
UV/Vis.: λ_{max}: 256 nm
Supplied as: A 10 mg/ml solution in chloroform
Storage: -20°C
Stability: ≥2 years
Item Origin: Plant/*Valeriana officinalis* L.



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

Laboratory Procedures

Valtrate is supplied as a solution in chloroform. To change the solvent, simply evaporate the chloroform under a gentle stream of nitrogen and immediately add the solvent of choice. Valtrate is soluble in DMSO. Valtrate is slightly soluble in methanol.

Description

Valtrate is a valepotriate that has been found in *Valeriana* and has diverse biological activities.¹⁻³ It inhibits nuclear export of influenza virus ribonucleoprotein complex (vRNP) in HeLa cells when used at a concentration of 3.6 μM and inhibits influenza virus infection in MDCK cells (IC₅₀ = 0.19 μM).¹ Valtrate is cytotoxic to GLC-4 lung and COLO 320 colorectal cancer cells (IC₅₀s = 1.4 and 3 μM, respectively).² It increases the number of entries into, and percentage of time spent in, the open arms of the elevated plus maze, indicating anxiolytic-like activity, and decreases serum corticosterone levels in rats when administered at a dose of 10 mg/kg.³

References

1. Watanabe, K., Takatsuki, H., Sonoda, M., *et al.* Anti-influenza viral effects of novel nuclear export inhibitors from *Valerianae Radix* and *Alpinia galanga*. *Drug Discov. Ther.* **5**(1), 26-31 (2016).
2. Bos, R., Hendriks, H., Scheffer, J.J., *et al.* Cytotoxic potential of valerian constituents and valerian tinctures. *Phytomedicine* **5**(3), 219-225 (1998).
3. Shi, S.-N., Shi, J.-L., Liu, Y., *et al.* The anxiolytic effects of valtrate in rats involves changes of corticosterone levels. *Evid. Based Complement. Alternat. Med.* **325948** (2014).

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, 06/25/2026

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