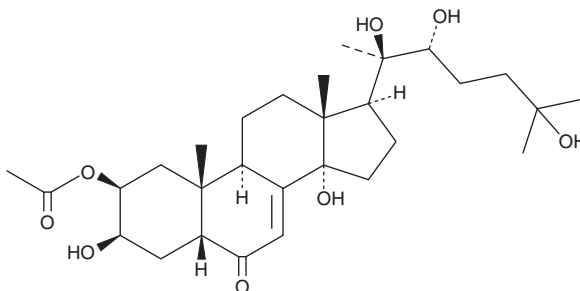


PRODUCT INFORMATION

2-O-Acetyl-20-hydroxyecdysone

Item No. 28063

CAS Registry No.: 19536-25-5
Formal Name: (2 β ,3 β ,5 β ,22R)-2-(acetyloxy)-3,14,20,22,25-pentahydroxycholest-7-en-6-one
Synonym: β -Ecdysone 2-acetate
MF: C₂₉H₄₆O₈
FW: 522.7
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 244 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Plant/*Cyanotis arachnoides*



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

Laboratory Procedures

2-O-Acetyl-20-hydroxyecdysone is supplied as a solid. A stock solution may be made by dissolving the 2-O-acetyl-20-hydroxyecdysone in the solvent of choice, which should be purged with an inert gas. 2-O-Acetyl-20-hydroxyecdysone is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 2-O-acetyl-20-hydroxyecdysone in DMSO and DMF is approximately 30 mg/ml and approximately 20 mg/ml in ethanol.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of 2-O-acetyl-20-hydroxyecdysone can be prepared by directly dissolving the solid in aqueous buffers. The solubility of 2-O-acetyl-20-hydroxyecdysone in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

2-O-Acetyl-20-hydroxyecdysone is an ecdysone steroid hormone that has been found in plants and animals.¹ It completely inhibits the growth of the bacteria *S. aureus*, *E. coli*, and *P. vulgaris*, as well as the fungi *C. albicans*, *A. alternata*, and *F. solani* when used at a concentration of 10 $\mu\text{g/ml}$. 2-O-Acetyl-20-hydroxyecdysone (50 μM) increases aggregation of amyloid- β (1-42) (A β 42), increases fibril formation by 60%, and decreases oligomer levels in cell-free assays.² It reduces A β 42-induced cytotoxicity of SH-SY5Y cells when used at concentrations of 5 and 10 μM .

References

- Shirshova, T.I., Politova, N.K., Burtseva, S.A., *et al.* Antimicrobial activity of natural ecdysteroids from *Serratula coronata* L. and their acyl derivatives. *Pharm. Chem. J.* **40**(5), 268-271 (2006).
- Yang, S.-G., Zhang, X., Sun, X.-S., *et al.* Diverse ecdysterones show different effects on amyloid- β 42 aggregation but all uniformly inhibit amyloid- β 42-induced cytotoxicity. *J. Alzheimers. Dis.* **22**(1), 107-117 (2010).

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, 12/22/2022

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