

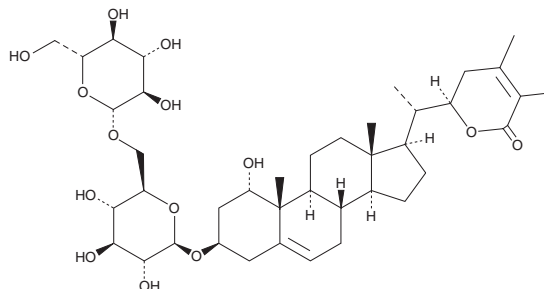
PRODUCT INFORMATION



Withanoside-V

Item No. 11806

CAS Registry No.: 256520-90-8
Formal Name: (1 α ,3 β ,22R)-3-[(6-O- β -D-glucopyranosyl- β -D-glucopyranosyl)oxy]-ergosta-5,24-dien-26-oic acid, δ -lactone
MF: C₄₀H₆₂O₁₄
FW: 766.9
Purity: $\geq$ 98%
Supplied as: A solid
Storage: -20°C
Stability: $\geq$ 4 years
Item Origin: Plant/*Withania somnifera*



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

Laboratory Procedures

Withanoside-V is supplied as a solid. A stock solution may be made by dissolving the withanoside-V in the solvent of choice, which should be purged with an inert gas. Withanoside-V is soluble ($\geq$ 10 mg/ml) in DMSO.

Description

Withanoside-V is a withanolide glycoside that has been found in *W. somnifera* and has diverse biological activities.¹⁻³ It inhibits iron-induced lipid peroxidation by 82.5% in a cell-free assay.¹ Withanoside-V (100 μ M) also inhibits the aggregation of amyloid- β (1-42) (A β 42) in a cell-free assay.² It decreases angiotensin-converting enzyme 2 (ACE2) levels in MCF-7 cells when used at a concentration of 5 μ M.³

References

1. Jayaprakasam, B., Stasburg, G.A., and Nair, M.G. Potent lipid peroxidation inhibitors from *Withania somnifera* fruits. *Tetrahedron* **60**(13), 3109-3121 (2004).
2. Dubey, S., Kallubai, M., and Subramanyam, R. Improving the inhibition of β -amyloid aggregation by withanolide and withanoside derivatives. *Int. J. Biol. Macromol.* **173**, 56-65 (2021).
3. Kalra, R.S., Kumar, V., Dhanjal, J.K., et al. COVID19-inhibitory activity of withanolides involves targeting of the host cell surface receptor ACE2: Insights from computational and biochemical assays. *J. Biomol. Struct. Dyn.* **40**(17), 7885-7895 (2021).

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, 05/06/2025

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