

PRODUCT INFORMATION



(+)- α -Tocopherol Acetate

Item No. 31795

CAS Registry No.: 58-95-7
Formal Name: (2R)-3,4-dihydro-2,5,7,8-tetramethyl-2-[(4R,8R)-4,8,12-trimethyltridecyl]-2H-1-benzopyran-6-ol, 6-acetate
Synonyms: D- α -Tocopherol Acetate, D-Vitamin E Acetate, RRR- α -Tocopherol Acetate

MF: C₃₁H₅₂O₃

FW: 472.7

Purity: \geq 95%

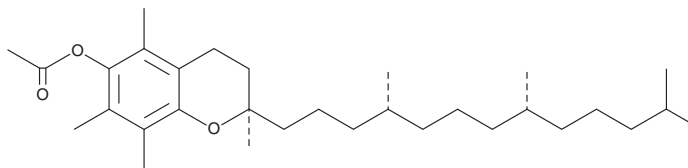
Supplied as: A semi-solid

Storage: 22°C

Stability: \geq 4 years

Special conditions: Protect from light and oxygen

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



Laboratory Procedures

(+)- α -Tocopherol acetate is supplied as a semi-solid. A stock solution may be made by dissolving the (+)- α -tocopherol acetate in the solvent of choice, which should be purged with an inert gas. (+)- α -Tocopherol acetate is sparingly soluble (1-10 mg/ml) in ethanol.

Description

(+)- α -Tocopherol acetate is an acetate form of (+)- α -tocopherol, a natural form of vitamin E.¹ It reduces liver necrosis and increases survival in male rats in a model of hepatotoxicity induced by carbon tetrachloride (CCl₄) when administered at a dose of 450 mg/kg three times at 72, 48, and 24 hours before CCl₄ administration.² (+)- α -Tocopherol acetate (20 mg/kg in the diet) decreases abdominal fat percentage and breast muscle malondialdehyde (MDA) levels and increases breast muscle total antioxidant capacity, total superoxide dismutase (SOD) activity, and glutathione peroxidase (GPX) activity in broiler chickens.³

References

1. Jensen, S.K. and Lauridsen, C. α -tocopherol stereoisomers. *Vitamins and Hormones*. Litwack, G., editor, Elsevier (2007).
2. Cawthorne, M.A., Bunyan, J., Sennitt, M.V., et al. Vitamin E and hepatotoxic agents. 3. Vitamin E, synthetic antioxidants and carbon tetrachloride toxicity in the rat. *Br. J. Nutr.* **24(2)**, 357-384 (1970).
3. Hu, Z.P., Wang, T., Ahmad, H., et al. Effects of different formulations of α -tocopherol acetate (vitamin E) on growth performance, meat quality and antioxidant capacity in broiler chickens. *Br. Poult. Sci.* **56(6)**, 687-695 (2015).

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, 07/10/2025

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