

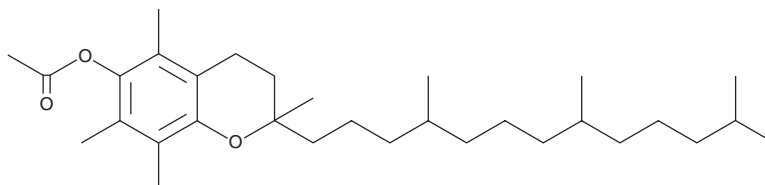
PRODUCT INFORMATION



(±)-α-Tocopherol Acetate

Item No. 28399

CAS Registry No.: 7695-91-2
Formal Name: 3,4-dihydro-2,5,7,8-tetramethyl-2-(4,8,12-trimethyltridecyl)-2H-1-benzopyran-6-ol, 6-acetate
Synonyms: all-*rac*-α-Tocopherol Acetate, DL-α-Tocopherol Acetate, DL-Vitamin E acetate
MF: C₃₁H₅₂O₃
FW: 472.7
Purity: ≥95%
Supplied as: A liquid
Storage: -20°C
Stability: ≥2 years
Item Origin: Synthetic



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 liquid. 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 soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide.

(±)-α-Tocopherol acetate is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, (±)-α-tocopherol acetate should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. (±)-α-Tocopherol acetate has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

(±)-α-Tocopherol acetate is an acetate form of (±)-α-tocopherol (Item No. 25985), a biologically active form of vitamin E. It prevents reductions in the percentage of embryos developing into blastocysts and the levels of glutathione (GSH), as well as prevents increases in oxidative stress, in zygotes from mice exposed to heat stress prior to mating when administered to dams prior to heat stress at a dose of 1 g/kg.¹ (±)-α-Tocopherol acetate (>1,050 IU/kg in the diet) prevents body weight loss and prevents decreases in pulmonary dynamic compliance in mice exposed to sidestream cigarette smoke but not in sham exposure animals.² However, it also increases the levels of IL-1β, IL-4, and IFN-γ in the bronchoalveolar lavage fluid (BALF) of sidestream cigarette smoke-exposed, but not sham exposure, mice.

References

1. Sakamoto, N., Ozawa, M., Yokotani-Tomita, K., *et al.* DL-α-Tocopherol acetate mitigates maternal hyperthermia-induced pre-implantation embryonic death accompanied by a reduction of physiological oxidative stress in mice. *Reproduction* **135**(4), 489-496 (2008).
2. Wang, S., Sun, N.N., Zhang, J., *et al.* Immunomodulatory effects of high-dose α-tocopherol acetate on mice subjected to sidestream cigarette smoke. *Toxicology* **175**(1-3), 235-245 (2002).

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, 11/03/2022

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