

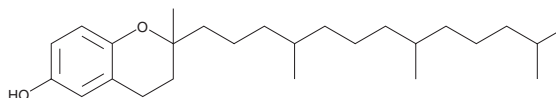
# PRODUCT INFORMATION



## (±)-Tocol

Item No. 27195

**CAS Registry No.:** 119-98-2  
**Formal Name:** 3,4-dihydro-2-methyl-2-(4,8,12-trimethyltridecyl)-2H-1-benzopyran-6-ol  
**Synonyms:** DL-Tocol, *rac*-Tocol  
**MF:** C<sub>26</sub>H<sub>44</sub>O<sub>2</sub>  
**FW:** 388.6  
**Purity:** ≥95%  
**Supplied as:** A solution in hexane  
**Storage:** -20°C  
**Stability:** ≥2 years  
**Concentration:** 50 mg/ml



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

### Laboratory Procedures

(±)-Tocol is supplied as a solution in hexane. To change the solvent, simply evaporate the hexane under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as methanol and ethanol purged with an inert gas can be used.

### Description

(±)-Tocol is a synthetic vitamin E derivative.<sup>1</sup> Unlike (±)- $\alpha$ -tocopherol (Item No. 25985), (±)-tocol does not suppress retinol-induced erythrocyte hemolysis or increase microviscosity of rat liver phosphatidylcholine (PC) liposomes.

### Reference

1. Urano, S., Inomori, Y., Sugawara, T., *et al.* Vitamin E: Inhibition of retinol-induced hemolysis and membrane-stabilizing behavior. *J. Biol. Chem.* **267**(26), 18365-18370 (1992).

#### 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, 11/15/2023

#### 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](http://WWW.CAYMANCHEM.COM)