

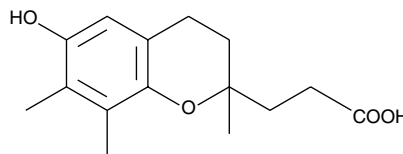
Product Information



γ -CEHC

Item No. 89630

CAS Registry No.: 178167-75-4
Formal Name: 3,4-dihydro-6-hydroxy-2,7,8-trimethyl-2H-1-Benzopyran-2-propanoic acid
Synonyms: γ -Tocopherol Metabolite, GTM, 2,7,8-trimethyl-2-(β -Carboxy-Ethyl)-6-Hydroxychroman
MF: C₁₅H₂₀O₄
FW: 264.3
Purity: $\geq 98\%$
Stability: ≥ 1 year at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max} : 296 nm



Laboratory Procedures

For long term storage, we suggest that γ -CEHC be stored as supplied at -20°C . It should be stable for at least one year. γ -CEHC is supplied as a crystalline solid. A stock solution may be made by dissolving the γ -CEHC in an organic solvent purged with an inert gas. γ -CEHC is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of γ -CEHC in these solvents is approximately 25 mg/ml.

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 γ -CEHC can be prepared by directly dissolving the crystalline compound in aqueous buffers. The solubility of γ -CEHC in PBS (pH 7.2) is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

The major tocopherol obtained from natural dietary sources is γ -tocopherol, whereas α -tocopherol is the form of Vitamin E typically obtained from synthetic supplements. γ -CEHC is a β -oxidized metabolite of dietary γ -tocopherol that functions as a natriuretic hormone.¹ It was initially purified and characterized from the urine of uremic patients, but it has since been found in urine from both control patients and those with congestive heart failure.² γ -CEHC is also anti-inflammatory, reducing 8-isoprostane and inflammatory eicosanoid synthesis in rat models.³

References

1. Christen, S., Woodall, A.A., Shigenaga, M.K., *et al.* γ -Tocopherol traps mutagenic electrophiles such as NO_x and complements α -tocopherol: Physiological implications. *Proc. Natl. Acad. Sci. USA* **94**, 3217-3222 (1997).
2. Wechter, W.J., Kantoci, D., Murray, J.E.D., *et al.* A new endogenous natriuretic factor: LLU- α . *Proc. Natl. Acad. Sci. USA* **93**, 6002-6007 (1996).
3. Jiang, Q., Elson-Schwab, I., Courtemanche, C., *et al.* γ -Tocopherol and its major metabolite, in contrast to α -tocopherol, inhibit cyclooxygenase activity in macrophages and epithelial cells. *Proc. Natl. Acad. Sci. USA* **97**, 11494-11499 (2000).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/89630

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY; NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's exclusive remedy and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 10/22/2013