

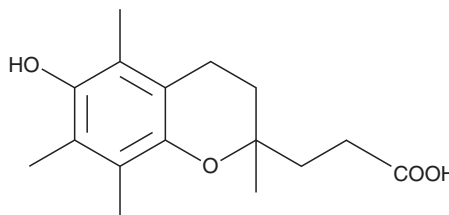
PRODUCT INFORMATION



α -CEHC

Item No. 10007705

CAS Registry No.: 4072-32-6
Formal Name: 3,4-dihydro-6-hydroxy-2,5,7,8-tetramethyl-2H-1-benzopyran-2-propanoic acid
MF: C₁₆H₂₂O₄
FW: 278.3
Purity: \geq 98%
UV/Vis.: λ_{max} : 206, 291 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

α -CEHC is supplied as a crystalline solid. A stock solution may be made by dissolving the α -CEHC in the solvent of choice, which should be purged with an inert gas. α -CEHC is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of α -CEHC in ethanol is approximately 10 mg/ml and approximately 20 mg/ml in DMSO and DMF.

α -CEHC is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, α -CEHC should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. α -CEHC has a solubility of approximately 0.25 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Molecules having vitamin E antioxidant activity include four tocopherols (α , β , δ , and γ) and four tocotrienols (α , β , δ , and γ).¹ α -Tocopherol is the major lipid soluble antioxidant *in vivo* and protects against lipid peroxidation.² α -CEHC is the major urinary metabolite of α -tocopherol following vitamin E supplementation.³ The concentration of α -CEHC in human serum is in the range of 5-10 pmol/ml but increases significantly up to 200 pmol/ml upon supplementation with RRR- α -tocopherol. About one-third of the α -CEHC circulating in the blood is present as a glucuronide conjugate.⁴ α -CEHC is excreted when a threshold concentration of 7-9 μ mol α -tocopherol/g total lipid in plasma is exceeded. Therefore, excretion of α -CEHC may be considered to be a marker of optimum vitamin E intake.⁵

References

1. Kamal-Eldin, A. and Appelqvist, L.-Å. The chemistry and antioxidant properties of tocopherols and tocotrienols. *Lipids* **31**, 671-701 (1996).
2. Burton, G.W., Joyce, A., and Ingold, K.U. Is vitamin E the only lipid-soluble, chain-breaking antioxidant in human blood plasma and erythrocyte membranes? *Arch. Biochem. Biophys.* **221(1)**, 281-290 (1983).
3. Traber, M.G., Elsner, A., and Brigelius-Flohé, R. Synthetic as compared with natural vitamin E is preferentially excreted as α -CEHC in human urine: Studies using deuterated α -tocopheryl acetates. *FEBS Lett.* **437**, 145-148 (1998).
4. Stahl, W.S., Graf, P., Brigelius-Flohé, R., et al. Quantification of the α - and γ -tocopherol metabolites 2,5,7,8-tetramethyl-2-(2'-carboxyethyl)-6-hydroxychroman and 2,7,8-trimethyl-2-(2'-carboxyethyl)-6-hydroxychroman in human serum. *Anal. Biochem.* **275**, 254-259 (1999).
5. Schultz, M., Leist, M., Petrzika, M., et al. Novel urinary metabolite of α -tocopherol, 2,5,7,8-tetramethyl-2(2'-carboxyethyl)-6-hydroxychroman, as an indicator of an adequate vitamin E supply. *Am. J. Clin. Nutr.* **62(6)**, 1527S-34S (1995).

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, 04/29/2024

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