PRODUCT INFORMATION



Ascorbyl Palmitate

Item No. 29604

CAS Registry No.:	137-66-6	
Formal Name:	L-ascorbic acid, 6-hexadecanoate	ОН
Synonyms:	Ascorbylpalmitic Acid, NSC 402451	
MF:	$C_{22}H_{38}O_7$	
FW:	414.5	
Purity:	≥95%	
UV/Vis.:	λ _{max} : 246 nm	Í Č Č Č Ŭ Ĭ Ĥ
Supplied as:	A solid	ОН
Storage:	-20°C	\checkmark \checkmark \land
Stability:	≥4 years	
Information represent	the product expecifications. Patch expecific an	adutical results are provided on each cortificate of analysis

Laboratory Procedures

Ascorbyl palmitate is supplied as a solid. A stock solution may be made by dissolving the ascorbyl palmitate in the solvent of choice, which should be purged with an inert gas. Ascorbyl palmitate is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of ascorbyl palmitate in ethanol is approximately 10 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Ascorbyl palmitate is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ascorbyl palmitate should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Ascorbyl palmitate has a solubility of approximately 0.5 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

Ascorbyl palmitate is a lipophilic derivative of ascorbic acid (Item No. 14656) with antioxidant and antiproliferative activities.¹⁻³ It scavenges hydroxyl radicals in cell-free assays.² Ascorbyl palmitate (0.01%) reduces the rate of autoxidation of soybean, safflower, sunflower, peanut, and corn oil.³ It inhibits increases in epidermal ornithine decarboxylase activity and DNA synthesis induced by phorbol 12-myristate 13-acetate (TPA; Item No. 10008014) in mice in a concentration-dependent manner when applied topically.¹ Ascorbyl palmitate (0.8 and 4 μ mol per 200 μ l of acetone, applied topically) reduces the number of tumors per mouse and the percentage of mice with tumors in a mouse skin two-stage model of tumor formation initiated and promoted by 7,12-dimethylbenz[a]anthracene (DMBA) and TPA, respectively. Formulations containing ascorbyl palmitate have been used as antioxidants and preservatives in foods, pharmaceuticals, and cosmetics.

References

- 1. Smart, R.C. and Crawford, C.L. Effect of ascorbic acid and its synthetic lipophilic derivative ascorbyl palmitate on phorbol ester-induced skin-tumor promotion in mice. Am. J. Clin. Nutr. 54(6 Suppl.), 1266S-1273S (1991).
- 2. Perricone, N., Nagy, K., Horváth, F., et al. The hydroxyl free radical reactions of ascorbyl palmitate as measured in various in vitro models. Biochem. Biophys. Res. Commun. 262(3), 661-665 (1999).
- 3. Cort, W.M. Antioxidant activity of tocopherols, ascorbyl palmitate, and ascorbic acid and their mode of action. J. Am. Oil Chem. Soc. 51(7), 321-325 (1974).

WARNING THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFFTY 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, 01/24/2025

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