

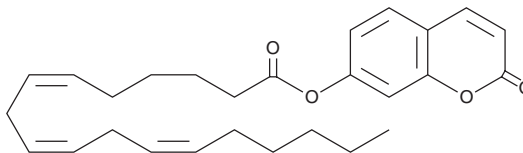
PRODUCT INFORMATION



7-hydroxycoumarinyl- γ -Linolenate

Item No. 10556

CAS Registry No.: 161180-12-7
Formal Name: 6Z,9Z,12Z-octadecatrienoic acid, 2-oxo-2H-1-benzopyran-7-yl ester
Synonym: Umbelliferol- γ -Linolenate
MF: C₂₇H₃₄O₄
FW: 422.6
Purity: \geq 98%
UV/Vis.: λ_{max} : 282, 311 nm
Supplied as: A solution in ethanol
Storage: -20°C
Stability: \geq 2 years



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

Laboratory Procedures

7-hydroxycoumarinyl- γ -Linolenate is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of 7-hydroxycoumarinyl- γ -linolenate in these solvents is approximately 25 and 50 mg/ml, respectively.

7-hydroxycoumarinyl- γ -Linolenate is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 7-hydroxycoumarinyl- γ -linolenate should be diluted with the aqueous buffer of choice. 7-hydroxycoumarinyl- γ -Linolenate has a solubility of approximately 50 μ g/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

7-hydroxycoumarinyl- γ -Linolenate is a γ -linolenic acid ester of 7-hydroxycoumarin (umbelliferone) that behaves as a substrate for cPLA₂. Hydrolysis of 7-hydroxycoumarinyl- γ -linolenate by phospholipase results in the release of the fluorescent compound 7-hydroxycoumarin which can be monitored spectrophotometrically (excitation at 335 nm, emission at 450 nm).¹

Reference

1. Huang, Z., Laliberté, F., Tremblay, N.M., *et al.* A continuous fluorescence-based assay for the human high-molecular-weight cytosolic phospholipase A₂. *Anal. Biochem.* **222**, 110-115 (1994).

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, 04/05/2024

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