

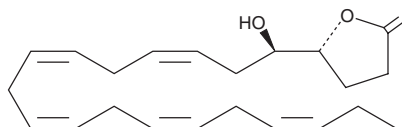
PRODUCT INFORMATION



(±)4(5)-DiHDPA lactone

Item No. 18344

CAS Registry No.: 845673-68-9
Formal Name: dihydro-5-[(3Z,6Z,9Z,12Z,15Z)-1-hydroxy-3,6,9,12,15-octadecapentaen-1-yl]-2(3H)-furanone
MF: C₂₂H₃₂O₃
FW: 344.5
Purity: ≥95%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



NOTE: Relative stereochemistry shown in chemical structure

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

Laboratory Procedures

(±)4(5)-DiHDPA lactone 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 ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of (±)4(5)-DiHDPA lactone in ethanol and DMF is approximately 30 mg/ml and approximately 25 mg/ml in DMSO.

(±)4(5)-DiHDPA lactone is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of (±)4(5)-DiHDPA lactone should be diluted with the aqueous buffer of choice. (±)4(5)-DiHDPA lactone has a solubility of approximately 0.3 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

(±)5(6)-DiHET lactone (Item No. 51212) is a 1,5 cyclic ester derived from (±)5(6)-DiHET (Item No. 51211), which, in turn, is a potential derivative of epoxidation of arachidonic acid (Item No. 90010) at the α-5 double bond. (±)4(5)-DiHDPA lactone is a derivative of docosahexaenoic acid (DHA; Item No. 90310) that is analogous to (+)5(6)-DiHET lactone. It is the 1,4 cyclic ester derived from (±)4(5)-DiHDPA, which is produced by epoxidation of DHA at the α-4 double bond. Its biological activity is unknown.

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, 04/16/2024

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