# **PRODUCT** INFORMATION



5(S),15(S)-DiHETE

Item No. 35280

CAS Registry No.:	82200-87-1	
Formal Name:	5S,15S-dihydroxy-6E,8Z,10Z,13E	-
	eicosatetraenoic acid	OH V
MF:	C <sub>20</sub> H <sub>32</sub> O <sub>4</sub>	СООН
FW:	336.5	
Purity:	≥95%	
UV/Vis.:	λ <sub>max</sub> : 243 nm	
Supplied as:	A solution in ethanol	
Storage:	-20°C	ÓН
Stability:	≥2 years	
Special Conditions: Oxygen and light sensitive		

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

# Laboratory Procedures

5(S),15(S)-DiHETE 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 5(S),15(S)-DiHETE in these solvents is approximately 50 mg/ml.

5(S),15(S)-DiHETE is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 5(S),15(S)-DiHETE should be diluted with the aqueous buffer of choice. The solubility of 5(S),15(S)-DiHETE in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

# Description

5(S),15(S)-DiHETE is synthesized by 15-lipoxygenase from 5(S)-HETE,<sup>1</sup> 5(S),15(S)-DiHETE potentiates the degranulation of human polymorphonuclear leukocytes in response to platelet activating factor, but not f-Met-Leu-Phe, calcium ionophore A23187, or Leukotriene B4.2 5(S),15(S)-DiHETE is chemotactic for eosinophils with an  $ED_{50}$  value of 0.3  $\mu$ M.<sup>3</sup>

# References

- 1. Green, F.A. Transformations of 5-HETE by activated keratinocyte 15-lipoxygenase and the activation mechanism. Lipids 25, 618-623 (1990).
- 2. O'Flaherty, J.T. and Thomas, M.J. Effect of 15-lipoxygenase-deprived arachidonate metabolites on human neutrophil degranulation. Prostaglandins Leukot. Med. 17, 199-212 (1985).
- 3. Morita, E., Schroder, J.M., and Christophers, E. Identification of a novel and highly potent eosinophil chemotactic lipid in human eosinophils treated with arachidonic acid. J. Immunol. 144, 1893-1900 (1990).

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

uyer 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, 01/16/2024

# 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