PRODUCT INFORMATION



(±)17-HETE

Item No. 10010636

CAS Registry No.:	128914-47-6	
Formal Name:	(±)17-hydroxy-5Z,8Z,11Z,14Z-	
	eicosatetraenoic acid	СООН
Synonym:	(±)17-Hydroxyeicosatetraenoic Acid	
MF:	$C_{20}H_{32}O_{3}$	
FW:	320.5	
Purity:	≥98%	
Supplied as:	A solution in ethanol	ОП
Storage:	-20°C	
Stability:	≥2 years	
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.		

Laboratory Procedures

(±)17-HETE 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 purged with an inert gas can be used. (±)17-HETE is miscible in these solvents. (±)17-HETE is soluble in 0.1 m Na₂CO₃ at a concentration of approximately 2 mg/ml.

(±)17-HETE is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of (±)17-HETE should be diluted with the aqueous buffer of choice. The solubility of (\pm) 17-HETE in PBS (pH 7.2) is approximately 0.8 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Electrolyte and fluid transport in the kidney are regulated in part by arachidonic acid and its metabolites. (±)17-HETE is the racemic version of a cytochrome P450 (CYP450) metabolite of arachidonic acid that has stereospecific effects on sodium transport in the kidney. At a concentration of 2 μ M the (S)-enantiomer of 17-HETE inhibits proximal tubule ATPase activity by as much as 70%, whereas the (R)-isomer is inactive.¹

Reference

1. Carroll, M.A., Balazy, M., Margiotta, P., et al. Cytochrome P-450-dependent HETEs: Profile of biological activity and stimulation by vasoactive peptides. Am. J. Physiol. 271(4 Pt 2), R863-R869 (1996).

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

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, 04/22/2024

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