

# 12(S)- Hydroperoxyeicosa- 5Z,8Z,10E,14Z- tetraenoic acid

## (HPETE)

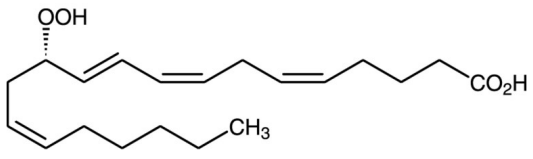
Primary metabolite of arachidonic acid via the 12-lipoxygenase pathway. Inhibits thromboxane synthase. Inhibits platelet aggregation *via* activation of guanylate cyclase (100% at 5 $\mu$ M). Stimulates leukotriene biosynthesis. **Thromboxane synthase inhibitor** Potent and selective inhibitor of Ca<sup>2+</sup>-calmodulin-dependent protein kinase II (IC<sub>50</sub>=0.7 $\mu$ M). Endogenous TRPV1 agonist (K<sub>i</sub>=0.35 $\mu$ M).

Citations: 6

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Stable for at least 6 months after receipt when stored at -80°C. Product is highly unstable in aqueous solution. We recommend using aqueous solutions within fifteen minutes. For experiments with cells or tissues, aliquots of ethanol solutions may be added
<b>Handling</b>	After opening, prepare aliquots and store at -80°C.
<b>Long Term Storage</b>	-80°C
<b>Shipping</b>	Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	12(S)-HPETE
<b>Appearance</b>	Liquid.
<b>CAS</b>	71774-10-2
<b>Couple Target</b>	CaM kinase, Guanylyl cyclase, Thromboxane synthase, TRP channel
<b>Couple Type</b>	Inhibitor, Ligand
<b>Formula</b>	$C_{20}H_{32}O_4$
<b>Formulation</b>	Solution in ethanol.
<b>MW</b>	336.5
<b>Purity</b>	≥98%
<b>Solubility</b>	Soluble in DMSO or dimethyl formamide; slightly soluble in PBS, pH 7.2 (0.8mg/ml) or 0.1M Na <sub>2</sub> CO <sub>3</sub> (2mg/ml).