

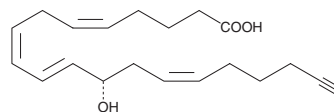
# PRODUCT INFORMATION



## 12(S)-HETE-19,20-alkyne

Item No. 9001937

<b>Formal Name:</b>	(S,5Z,8Z,10E,14Z)-12-hydroxyicosa-5,8,10,14-tetraen-19-ynoic acid
<b>Synonyms:</b>	Click Tag™ 12(S)-HETE-19,20-alkyne, 12(S)-Hydroxyeicosatetraenoic Acid-19,20-alkyne
<b>MF:</b>	C <sub>20</sub> H <sub>28</sub> O <sub>3</sub>
<b>FW:</b>	316.4
<b>Purity:</b>	≥97%
<b>UV/Vis.:</b>	λ <sub>max</sub> : 236 nm
<b>Supplied as:</b>	A solution in ethanol
<b>Storage:</b>	-20°C
<b>Stability:</b>	≥1 year



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

### Laboratory Procedures

12(S)-HETE-19,20-alkyne 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. 12(S)-HETE-19,20-alkyne is miscible in these solvents.

12(S)-HETE-19,20-alkyne is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 12(S)-HETE-19,20-alkyne should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. The solubility of 12(S)-HETE-19,20-alkyne in PBS (pH 7.2) is approximately 0.5 mg/ml. For greater aqueous solubility, 12(S)-HETE-19,20-alkyne can be directly dissolved in 0.1 M Na<sub>2</sub>CO<sub>3</sub> (solubility of 2 mg/ml) and then diluted with PBS (pH 7.2) to achieve the desired concentration or pH. We do not recommend storing the aqueous solution for more than one day.

### Description

12(S)-HETE (Item No. 34570) is the predominant lipoxygenase product of mammalian platelets.<sup>1</sup> It enhances tumor cell adhesion to endothelial cells, fibronectin, and the subendothelial matrix at 0.1 μM.<sup>2,3</sup> 12(S)-HETE-19,20-alkyne is a form of 12(S)-HETE with an ω-terminal alkyne. The terminal alkyne group can be used in click chemistry linking reactions, to tag 12(S)-HETE with fluorescent or biotinylated labels for analysis of its metabolism and biological activity.<sup>4,5</sup>

### References

1. Hamberg, M. and Samuelsson, B. Prostaglandin endoperoxides. Novel transformations of arachidonic acid in human platelets. *Proc. Natl. Acad. Sci. USA* **71(9)**, 3400-3404 (1974).
2. Grossi, I.M., Fitzgerald, L.A., Umbarger, L.A., *et al.* Bidirectional control of membrane expression and/or activation of the tumor cell IRGpIIb/IIIa receptor and tumor cell adhesion by lipoxygenase products of arachidonic acid and linoleic acid. *Cancer Res.* **49(4)**, 1029-1037 (1989).
3. Honn, K.V., Nelson, K.K., Renaud, C., *et al.* Fatty acid modulation of tumor cell adhesion to microvessel endothelium and experimental metastasis. *Prostaglandins* **44(5)**, 413-429 (1992).
4. Kolb, H.C. and Sharpless, K.B. The growing impact of click chemistry on drug discovery. *Drug Discov. Today* **8(24)**, 1128-1137 (2003).
5. Lutz, J.-F. and Zarafshani, Z. Efficient construction of therapeutics, bioconjugates, biomaterials and bioactive surfaces using azide-alkyne "click" chemistry. *Adv. Drug Deliv. Rev.* **60(9)**, 958-970 (2008).

#### 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, 08/24/2023

#### 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