

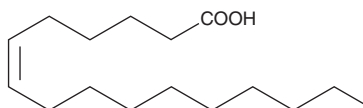
# PRODUCT INFORMATION



## *cis*-6-Hexadecenoic Acid

Item No. 9001845

**CAS Registry No.:** 17004-51-2  
**Formal Name:** 6Z-hexadecenoic acid  
**Synonyms:** C16:1(6Z), C16:1Δ6, C16:1ω10, FA 16:1, Sapienic Acid  
**MF:** C<sub>16</sub>H<sub>30</sub>O<sub>2</sub>  
**FW:** 254.4  
**Purity:** ≥99%  
**Supplied as:** A solution in ethanol  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Laboratory Procedures

*cis*-6-Hexadecenoic acid 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 methanol, chloroform, and ether purged with an inert gas can be used.

### Description

*cis*-6-Hexadecenoic acid is a monounsaturated fatty acid and is one of the primary fatty acids in human skin.<sup>1</sup> *cis*-6-Hexadecenoic acid levels are increased in isolated sebum from the face and back of patients with acne.<sup>2</sup> In contrast, levels are decreased in the non-lesional skin and isolated sebum of atopic dermatitis patients, which correlates with an increase in *S. aureus* in the sebum.<sup>1</sup> It is active against *S. aureus in vitro* when used at a concentration of 5 μg/ml at pH 5.5.<sup>3</sup> *cis*-6-Hexadecenoic acid disrupts membrane integrity, the proton motive force, increases membrane fluidity, and inhibits the electron transport chain in *S. aureus*.

### References

1. Takigawa, H., Nakagawa, H., Kuzukawa, M., *et al.* Deficient production of hexadecenoic acid in the skin is associated in part with the vulnerability of atopic dermatitis patients to colonization by *Staphylococcus aureus*. *Dermatology* **211(3)**, 240-248 (2005).
2. Li, W.-H., Zhang, Q., Flach, C.R., *et al.* In vitro modeling of unsaturated free fatty acid-mediated tissue impairments seen in acne lesions. *Arch. Dermatol. Res.* **309(7)**, 529-540 (2017).
3. Cartron, M.L., England, S.R., Chiriac, A.I., *et al.* Bactericidal activity of the human skin fatty acid *cis*-6-hexadecanoic acid on *Staphylococcus aureus*. *Antimicrob. Agents Chemother.* **58(7)**, 3599-3609 (2014).

#### 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, 02/15/2024

#### 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