

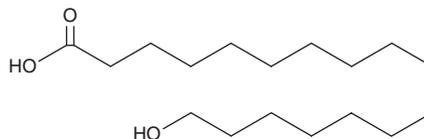
# PRODUCT INFORMATION



## 17-hydroxy Heptadecanoic Acid

Item No. 24650

**CAS Registry No.:** 13099-34-8  
**Formal Name:** 17-hydroxy-heptadecanoic acid  
**Synonyms:** FA 17:0;O, ω-hydroxy Heptadecanoic Acid  
**MF:** C<sub>17</sub>H<sub>34</sub>O<sub>3</sub>  
**FW:** 286.5  
**Purity:** ≥98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

17-hydroxy Heptadecanoic acid is supplied as a solid. A stock solution may be made by dissolving the 17-hydroxy heptadecanoic acid in the solvent of choice. 17-hydroxy Heptadecanoic acid is soluble in chloroform, ethanol (warmed), and ethyl ether.

### Description

17-hydroxy Heptadecanoic acid is a hydroxylated fatty acid that has been found in aeolian particles collected on the Frioul Islands and in the outer bark of *E. globulus*.<sup>1,2</sup> It has been used in the synthesis of various macrocyclic lactones.<sup>3</sup>

### References

1. Rontani, J.-F., Charriere, B., Vaultier, F., *et al.* Origin and degradation of lipids in aeolian particles from a coastal area of the north-western Mediterranean Sea. *Atmos. Environ.* **56**, 124-135 (2012).
2. Freire, C.S.R., Silvestre, A.J.D., Neto, C.P., *et al.* Lipophilic extractives of the inner and outer barks of *Eucalyptus globulus*. *Holzforschung* **56(4)**, 372-379 (2002).
3. Scott, L.T. and Naples, J.O. Macrolide synthesis by the polystyrene-mediated lactonization of ω-hydroxycarboxylic acids. *Synthesis* **11**, 738-739 (1976).

#### 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, 02/27/2024

#### 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