

# PRODUCT DATA SHEET

## **cis-9,10-Methyleneoctadecanoic acid**

**Catalog No:** 1822

**Common Name:** Dihydrosterculic acid

**Source:** synthetic

**Solubility:** chloroform, methanol, ethanol,  
hexane

**CAS No:** 4675-61-0

**Molecular Formula:** C<sub>19</sub>H<sub>36</sub>O<sub>2</sub>

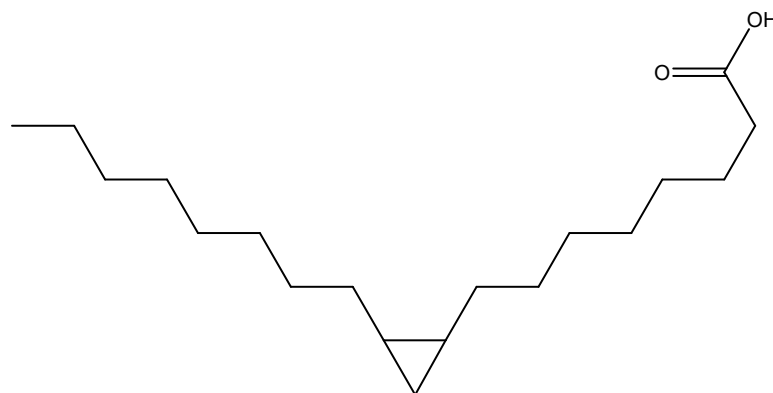
**Molecular Weight:** 297

**Storage:** -20°C

**Purity:** TLC > 98%, GC > 98%

**TLC System:** hexane/ethyl ether/acetic acid  
(80:20:1 by vol.)

**Appearance:** solid



### **Application Notes:**

This cyclopropanoid fatty acid is a major constituent of some seed oils and also occurs in some bacterial membranes but is not synthesized or used by humans.<sup>1</sup> Dihydrosterculic acid is a major constituent of the phospholipids of many trypanosomatid flagellates including some pathogenic species. 10-thiastearic acid has been found to be a potent inhibitor of dihydrosterculic acid synthesis and has been used as a therapeutic drug against these organisms.<sup>2</sup> The enzyme S-adenosylmethionine donates a methylene group to oleic acid in the sn-1 position of phosphatidylethanolamine to form dihydrosterculic acid. Dihydrosterculic acid is further desaturated to sterculic acid by cyclopropane desaturase.

### **Selected References:**

1. G. Knothe "NMR Characterization of Dihydrosterculic Acid and Its Methyl Ester" *Lipids*, Vol. 41(4) pp. 393-396, 2006
2. R. Pascal, Jr, S. Mannarelli, D. Ziering "10-Thiastearic acid inhibits both dihydrosterculic acid biosynthesis and growth of the protozoan *Crithidia fasciculata*" *The Journal of Biological Chemistry*, Vol. 261 pp. 12441-12443, 1986

This product is to be used for research only. It is not intended for drug or diagnostic use, human consumption or to be used in food or food additives. Matreya assumes no liability for any use of this product by the end user. We believe the information, offered in good faith, is accurate.