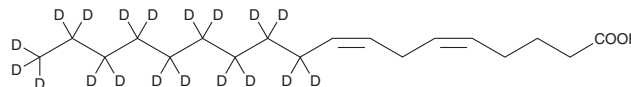


# PRODUCT INFORMATION



## Sebaleic Acid-d<sub>19</sub> Item No. 28228

**CAS Registry No.:** 2692624-10-3  
**Formal Name:** (5Z,8Z)-5,8-octadecadienoic-10,10,11,11,12,12,13,13,14,14,15,15,16,16,17,17,18,18,18-d<sub>19</sub> acid  
**Synonyms:** 5(Z),8(Z)-Octadecadienoic Acid-d<sub>19</sub>, C18:2-d<sub>19</sub>, *cis,cis*-5,8-Octadecadienoic Acid-d<sub>19</sub>, FA 18:2-d<sub>19</sub>  
**MF:** C<sub>18</sub>H<sub>13</sub>D<sub>19</sub>O<sub>2</sub>  
**FW:** 299.6  
**Chemical Purity:** ≥98% (Sebaleic Acid)  
**Deuterium Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>19</sub>); ≤1% d<sub>0</sub>  
**Supplied as:** A solution in ethanol, with 5 µg BHT  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Laboratory Procedures

Sebaleic acid-d<sub>19</sub> is intended for use as an internal standard for the quantification of sebaleic acid by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated *versus* unlabeled).

### Description

Sebaleic acid is a fatty acid that has been found in human sebum, hair, and nails.<sup>1</sup>

### Reference

1. Destailhats, F., Guitard, M., and Cruz-Hernandez, C. Identification of Δ<sup>6</sup>-monounsaturated fatty acids in human hair and nail samples by gas-chromatography-mass-spectrometry using ionic-liquid coated capillary column. *J. Chromatogr.* **8(52)**, 9384-9389 (2011).

**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, 03/25/2024

### 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