

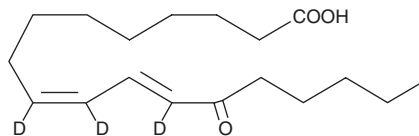
# PRODUCT INFORMATION



## 13-OxoODE-d<sub>3</sub>

Item No. 338620

**Formal Name:** 13-oxo-9Z,11E-9,10,12,-d<sub>3</sub>-octadecadienoic acid  
**Synonym:** 13-KODE-d<sub>3</sub>  
**MF:** C<sub>18</sub>H<sub>27</sub>D<sub>3</sub>O<sub>3</sub>  
**FW:** 297.5  
**Chemical Purity:** ≥98% 13-OxoODE  
**Deuterium Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>3</sub>); ≤1% d<sub>0</sub>  
**UV/Vis.:** λ<sub>max</sub>: 275 nm ε: 22,000  
**Supplied as:** A solution in acetonitrile  
**Storage:** -80°C  
**Stability:** ≥1 year



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

### Laboratory Procedures

13-OxoODE-d<sub>3</sub> is intended for use as an internal standard for the quantification of 13-oxoODE (Item No. 38620) 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).

13-OxoODE-d<sub>3</sub> is supplied as a solution in acetonitrile. To change the solvent, simply evaporate the acetonitrile 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. The solubility of 13-oxoODE-d<sub>3</sub> in these solvents is approximately 50 mg/ml.

### Description

13-OxoODE is produced from 13-HODE by a NAD<sup>+</sup>-dependent dehydrogenase present in rat colonic mucosa.<sup>1</sup> 13-OxoODE has been shown to stimulate cell proliferation when instilled intra-rectally in rats.<sup>2</sup> 13-OxoODE has also been detected in preparations of rabbit reticulocyte plasma and mitochondrial membranes, mostly still esterified to phospholipids. Production of 13-oxoODE is putatively linked to the maturation of reticulocytes to erythrocytes through the activity of 15-lipoxygenase.<sup>3,4</sup>

### References

1. Earles, S.M., Bronstein, J.C., Winner, D.L., *et al.* Metabolism of oxidized linoleic acid: Characterization of 13-hydroxyoctadecadienoic acid dehydrogenase activity from rat colonic tissue. *Biochim. Biophys. Acta* **1081**, 174-180 (1991).
2. Bull, A.W., and Bronstein, J.C. Production of unsaturated carbonyl compounds during metabolism of hydroperoxy fatty acids by colonic homogenates. *Carcinogenesis* **11**, 1699-1704 (1990).
3. Kühn, H., Belkner, J., Wiesner, R., *et al.* Occurrence of 9- and 13-keto-octadecadienoic acid in biological membranes oxygenated by the reticulocyte lipoxygenase. *Arch. Biochem. Biophys.* **279**, 218-224 (1990).
4. Kühn, H., Belkner, J., and Wiesner, R. Subcellular distribution of lipoxygenase products in rabbit reticulocyte membranes. *Eur. J. Biochem.* **191**, 221-227 (1990).

#### 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, 11/08/2017

#### 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