

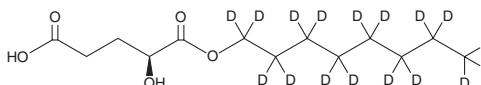
# PRODUCT INFORMATION



## (2S)-Octyl- $\alpha$ -hydroxyglutarate-d<sub>17</sub>

Item No. 19695

**CAS Registry No.:** 2748489-76-9  
**Formal Name:** 2S-hydroxy-pentanedioic acid,  
1-octyl-d<sub>17</sub> ester  
**Synonym:** (2S)-Octyl-2-HG-d<sub>17</sub>  
**MF:** C<sub>13</sub>H<sub>7</sub>D<sub>17</sub>O<sub>5</sub>  
**FW:** 277.4  
**Chemical Purity:** ≥95% ((2S)-Octyl- $\alpha$ -hydroxyglutarate)  
**Deuterium**  
**Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>17</sub>); ≤1% d<sub>0</sub>  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Laboratory Procedures

(2S)-Octyl- $\alpha$ -hydroxyglutarate-d<sub>17</sub> contains 17 deuterium atoms located on the octyl group. It is intended for use as an internal standard for the quantification of (2S)-octyl- $\alpha$ -hydroxyglutarate (Item No. 16367) 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).

(2S)-Octyl- $\alpha$ -hydroxyglutarate-d<sub>17</sub> is supplied as a crystalline solid. A stock solution may be made by dissolving the (2S)-octyl- $\alpha$ -hydroxyglutarate-d<sub>17</sub> in the solvent of choice. (2S)-Octyl- $\alpha$ -hydroxyglutarate-d<sub>17</sub> is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of (2S)-octyl- $\alpha$ -hydroxyglutarate-d<sub>17</sub> in ethanol is approximately 20 mg/ml and approximately 10 mg/ml in DMSO and DMF.

### Description

(2S)-Octyl- $\alpha$ -hydroxyglutarate is a cell-permeable derivative of the L-isomer of  $\alpha$ -hydroxyglutarate (Item No. 16374).

#### 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, 08/17/2022

#### 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