

# PRODUCT INFORMATION



## 1-Palmitoyl-2-10-OAHSA-3-Oleoyl-*sn*-glycerol

Item No. 29406

**Formal Name:** 16-(((R)-1-(oleoyloxy)-3-(stearoyloxy)propan-2-yl)oxy)-16-oxohexadecan-7-yl oleate

**Synonyms:** 16:0-10-OAHSA-18:1-TG, TG(16:0/10-OAHSA/18:1)

**MF:** C<sub>73</sub>H<sub>136</sub>O<sub>8</sub>

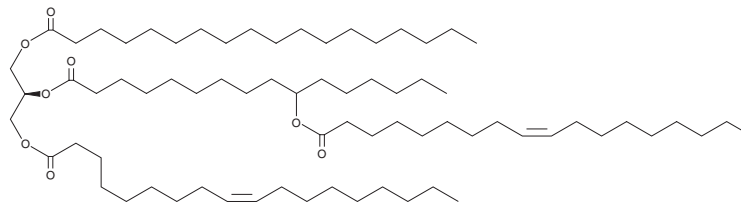
**FW:** 1,141.9

**Purity:** ≥95%

**Supplied as:** A solution in methyl acetate

**Storage:** -20°C

**Stability:** ≥2 years



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

### Laboratory Procedures

1-Palmitoyl-2-10-OAHSA-3-oleoyl-*sn*-glycerol is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of 1-palmitoyl-2-10-OAHSA-3-oleoyl-*sn*-glycerol in ethanol and DMF is approximately 20 mg/ml and approximately 15 mg/ml in DMSO.

1-Palmitoyl-2-10-OAHSA-3-oleoyl-*sn*-glycerol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 1-palmitoyl-2-10-OAHSA-3-oleoyl-*sn*-glycerol should be diluted with the aqueous buffer of choice. 1-Palmitoyl-2-10-OAHSA-3-oleoyl-*sn*-glycerol has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method.

### Description

1-Palmitoyl-2-10-OAHSA-3-oleoyl-*sn*-glycerol is a triacylglycerol that contains palmitic acid (Item No. 10006627) at the *sn*-1 position, the fatty acid ester of hydroxy fatty acid (FAHFA) 10-OAHSA (Item No. 19975) at the *sn*-2 position, and oleic acid (Item Nos. 90260 | 24659) at the *sn*-3 position. FAHFA-containing triacylglycerol levels are greater than 100-fold higher than non-esterified FAHFAs in mouse adipose tissues, indicating the triacylglycerols act as intracellular storage reservoirs for FAHFAs.<sup>1</sup>

### Reference

1. Tan, D., Ertunc, M.E., Konduri, S., *et al.* Discovery of FAHFA-containing triacylglycerols and their metabolic regulation. *J. Am. Chem. Soc.* **141**(22), 8798-8806 (2019).

#### 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, 11/10/2023

#### 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