

# PRODUCT INFORMATION



## all-*cis*-1,2,3-Docosahexaenoyl Glycerol

Item No. 10009667

CAS Registry No.: 124596-98-1

Formal Name: (4Z,4'Z,4''Z,7Z,7'Z,7''Z,10Z,10'Z,10''Z,13Z,13'Z,13''Z,16Z,16'Z,16''Z,19Z,19'Z,19''Z)-4,7,10,13,16,19-docosahexaenoic acid, 1,1',1''-(1,2,3-propanetriyl) ester

Synonyms: all-*cis*-1,2,3-Docosahexaenoyl-*rac*-glycerol,  
all-*cis*-Glycerol Tridocosahexaenoate,  
all-*cis*-Tridocosahexaenoic Acid,  
all-*cis*-Tridocosahexaenoylglycerol,  
Glycerol Tridocosahexaenoyl

MF: C<sub>69</sub>H<sub>98</sub>O<sub>6</sub>

FW: 1,023.5

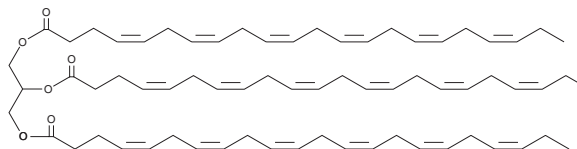
Purity: ≥98%

UV/Vis.: λ<sub>max</sub>: 205 nm

Supplied as: A solution in ethanol

Storage: -20°C

Stability: ≥2 years



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

### Laboratory Procedures

all-*cis*-1,2,3-Docosahexaenoyl glycerol is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as dimethyl formamide purged with an inert gas can be used. The solubility of all-*cis*-1,2,3-docosahexaenoyl glycerol in dimethyl formamide is approximately 10 mg/ml.

all-*cis*-1,2,3-Docosahexaenoyl glycerol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of all-*cis*-1,2,3-docosahexaenoyl glycerol should be diluted with the aqueous buffer of choice. all-*cis*-1,2,3-Docosahexaenoyl 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

all-*cis*-1,2,3-Docosahexaenoyl glycerol is a polyunsaturated triacylglycerol that contains all-*cis* docosahexaenoic acid (DHA; Item No. 90310) at the *sn*-1, *sn*-2, and *sn*-3 positions.

#### 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, 06/04/2020

#### 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