

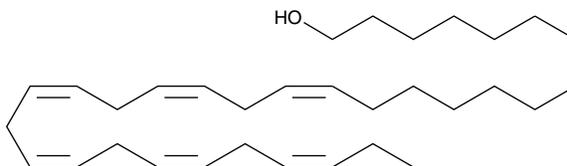
# Product Information



## CAY10631

Item No. 10562

**Formal Name:** 14Z,17Z,20Z,23Z,26Z,29Z-dotriaconta-hexaen-1-ol  
**MF:** C<sub>32</sub>H<sub>54</sub>O  
**FW:** 454.8  
**Purity:** ≥95%  
**Stability:** ≥1 year at -20°C  
**Supplied as:** A solution in ethanol



### Laboratory Procedures

For long term storage, we suggest that CAY10631 be stored as supplied at -20°C. It should be stable for at least one year.

CAY10631 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 DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of CAY10631 in these solvents is approximately 10 mg/ml.

If aqueous stock solutions are required for biological experiments, they can best be prepared by diluting the organic solvent into aqueous buffers or isotonic saline. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Very long chain fatty acids derived from docosahexaenoic acid by the action of fatty acid elongases are important in retinal development and macular diseases.<sup>1-3</sup> These fatty acids are commonly incorporated into the *sn*-1 position of membrane phospholipids, predominantly phosphatidylcholine, and are involved in binding the photoreceptor rhodopsin.<sup>4</sup> CAY10631 is an alcohol related to a C32:6 very long chain fatty acid found in the retina. Its biological actions are unknown.

### References

1. Agbaga, M.-P., Mandal, M.N.A., and Anderson, R.E. Retinal very long-chain PUFAs: New insights from studies on ELOVL4 protein. *J. Lipid Res.* **51**, 1624-1642 (2010).
2. McMahon, A., Jackson, S.N., Woods, A.S., *et al.* A Stargardt disease-3 mutation in the mouse Elov4 gene causes retinal deficiency of C32-C36 acyl phosphatidylcholines. *FEBS Lett.* **581**, 5459-5463 (2007).
3. Bedell, M., Harkewicz, R., Wang, X., *et al.* Focus on molecules: ELOVL4. *Exp. Eye Res.* **90**, 476-477 (2010).
4. Aveldaño, M.I. Phospholipid species containing long and very long polyenoic fatty acids remain with rhodopsin after hexane extraction of photoreceptor membranes. *Biochemistry* **27**, 1229-1239 (1988).

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/10562](http://www.caymanchem.com/catalog/10562)

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY. NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

#### SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery.**

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog.**

Copyright Cayman Chemical Company, 07/23/2014

### Cayman Chemical

#### Mailing address

1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

#### Phone

(800) 364-9897  
(734) 971-3335

#### Fax

(734) 971-3640

#### E-Mail

[custserv@caymanchem.com](mailto:custserv@caymanchem.com)

#### Web

[www.caymanchem.com](http://www.caymanchem.com)