

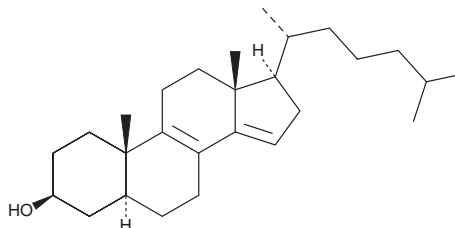
# PRODUCT INFORMATION



## 14-dehydro Zymostenol

Item No. 29534

CAS Registry No.: 19431-20-0  
Formal Name: (3 $\beta$ ,5 $\alpha$ )-cholesta-8,14-dien-3-ol  
Synonyms:  $\Delta^8$ ,14-Cholestandienol,  
Cholesta-8,14-dien-3 $\beta$ -ol  
MF: C<sub>27</sub>H<sub>44</sub>O  
FW: 384.6  
Purity:  $\geq 95\%$   
UV/Vis.:  $\lambda_{\text{max}}$ : 250 nm  
Supplied as: A crystalline solid  
Storage: -20°C  
Stability:  $\geq 2$  years



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

### Laboratory Procedures

14-dehydro Zymostenol is supplied as a crystalline solid. A stock solution may be made by dissolving the 14-dehydro zymostenol in the solvent of choice, which should be purged with an inert gas. 14-dehydro Zymostenol is soluble in ethanol and dimethyl formamide (DMF). The solubility of 14-dehydro zymostenol in these solvents is approximately 2 and 3 mg/ml, respectively.

### Description

14-dehydro Zymostenol is a precursor to cholesterol (Item No. 9003100).<sup>1</sup> It increases the percentage of myelin basic protein-positive (MBP<sup>+</sup>) oligodendrocytes formed from oligodendrocyte precursor cells when used at concentrations of 5.8 and 17  $\mu\text{M}$ .<sup>2</sup>

### References

1. Lutsky, B.N. and Schroepfer, G.J., Jr. Studies on the enzymic conversion of 5 $\alpha$ -cholesta-8,14-dien-3 $\beta$ -ol to cholesterol. *J. Biol. Chem.* **245**(23), 6449-6455 (1970).
2. Hubler, Z., Allimuthu, D., Bederman, I., *et al.* Accumulation of 8,9-unsaturated sterols drives oligodendrocyte formation and remyelination. *Nature* **560**(7718), 372-376 (2018).

#### 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, 05/07/2020

#### 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