

PRODUCT INFORMATION



2,3-Oxidosqualene

Item No. 36187

CAS Registry No.: 7200-26-2

Formal Name: 2,2-dimethyl-3-[(3E,7E,11E,15E)-3,7,12,16,20-pentamethyl-3,7,11,15,19-heneicosapentaen-1-yl]-oxirane

Synonyms: (3R,S)-Oxidosqualene, Squalene 2,3-oxide

MF: $C_{30}H_{50}O$

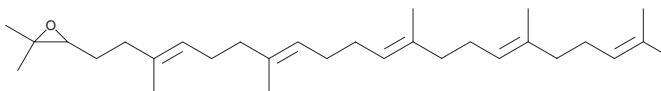
FW: 426.7

Purity: $\geq 90\%$

Supplied as: A neat oil

Storage: $-80^{\circ}C$

Stability: ≥ 2 years



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

Laboratory Procedures

2,3-Oxidosqualene is supplied as a neat oil. A stock solution may be made by dissolving the 2,3-oxidosqualene in the solvent of choice, which should be purged with an inert gas. 2,3-Oxidosqualene is slightly soluble in chloroform.

Description

2,3-Oxidosqualene is an intermediate in the biosynthesis of sterols such as lanosterol (Item No. 19521), cycloartenol, and cholesterol (Item No. 9003100).^{1,2} It is also an intermediate in the biosynthesis of the triterpenoid β -amyrin (Item No. 20949).³

References

1. Nes, W.D. Biosynthesis of cholesterol and other sterols. *Chem. Rev.* **111**(10), 6423-6451 (2011).
2. Willett, J.D., Sharpless, K.B., Lord, K.E., *et al.* Squalene-2,3-oxide, an intermediate in the enzymatic conversion of squalene to lanosterol and cholesterol. *J. Biol. Chem.* **242**(18), 4182-4191 (1967).
3. Kajikawa, M., Yamato, K.T., Fukuzawa, H., *et al.* Cloning and characterization of a cDNA encoding beta-amyrin synthase from petroleum plant *Euphorbia tirucalli* L. *Phytochemistry* **66**(15), 1759-1766 (2005).

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, 01/27/2022

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