

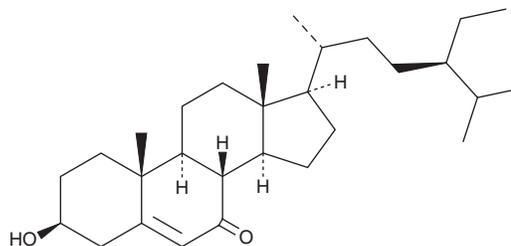
# PRODUCT INFORMATION



## 7-Ketositosterol

Item No. 37189

CAS Registry No.:	2034-74-4
Formal Name:	3 $\beta$ -hydroxy-stigmast-5-en-7-one
Synonyms:	7-oxo- $\beta$ -Sitosterol, 7-Oxositosterol
MF:	C <sub>29</sub> H <sub>48</sub> O <sub>2</sub>
FW:	428.7
Purity:	≥98%
UV/Vis.:	$\lambda_{\text{max}}$ : 238 nm
Supplied as:	A solid
Storage:	-20°C
Stability:	≥4 years
Item Origin:	Synthetic



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

### Laboratory Procedures

7-Ketositosterol is supplied as a solid. A stock solution may be made by dissolving the 7-ketositosterol in the solvent of choice, which should be purged with an inert gas. 7-Ketositosterol is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of 7-ketositosterol in these solvents is approximately 3 and 12 mg/ml, respectively.

### Description

7-Ketositosterol is a phytosterol and phytosterol oxidation product (POP).<sup>1,2</sup> It is a thermal oxidation product of sitosterol (24 $\alpha$ -ethyl cholesterol; Item No. 11756).<sup>2</sup> 7-Ketositosterol inhibits the proliferation of HL-60 acute myeloid leukemia (AML) cells (IC<sub>50</sub> = 45.6  $\mu$ M).<sup>3</sup>

### References

1. Yu, J.-D., Chen, M.-X., Wu, H., *et al.* Phytochemical and chemotaxonomic study on *Pholidota pallida* lindl. *Biochem. Syst. Ecol.* **97**, 104290 (2021).
2. McDonald, K., Langenbahn, H.J., Miller, J.D., *et al.* Phytosterol oxidation products from coffee silverskin. *J. Food Sci.* **87(2)**, 728-737 (2022).
3. Oyama, M., Tokiwano, T., Ota, H., *et al.*  $\beta$ -sitosterol oxides from the dried stem of *salix gilgiana* inhibit the proliferation of HL-60 leukemic cells. *Curr. Bioact. Compd.* **16(3)**, 329-334 (2020).

#### 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, 11/10/2022

#### 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