

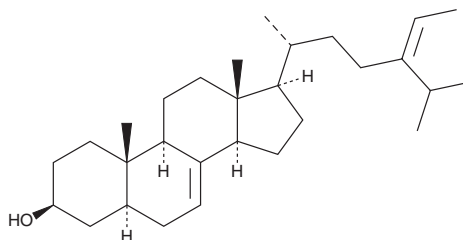
# PRODUCT INFORMATION



## $\Delta^7$ -Avenasterol

Item No. 30155

CAS Registry No.: 7212-91-1  
Formal Name: (3 $\beta$ ,5 $\alpha$ )-stigmasta-7,24(28)-dien-3-ol  
MF: C<sub>29</sub>H<sub>48</sub>O  
FW: 412.7  
Purity:  $\geq 95\%$  (mixture of isomers)  
Supplied as: A solid  
Storage: -20°C  
Stability:  $\geq 4$  years  
Item Origin: Synthetic



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

### Laboratory Procedures

$\Delta^7$ -Avenasterol is supplied as a solid. A stock solution may be made by dissolving the  $\Delta^7$ -avenasterol in the solvent of choice, which should be purged with an inert gas.  $\Delta^7$ -Avenasterol is slightly soluble in chloroform and methanol.

### Description

$\Delta^7$ -Avenasterol is a phytosterol that has been found in a variety of plant oils and grains, including olive and soybean oils and wheat, oats, and rye, and has antioxidant activities.<sup>1,2</sup> It reduces temperature-induced oxidation of safflower oil.<sup>3</sup>

### References

1. Yorulmaz, H.O. and Konuskan, D.B. Antioxidant activity, sterol and fatty acid compositions of turkish olive oils as an indicator of variety and ripening degree. *J. Food Sci. Technol.* **54**(12), 4067-4077 (2017).
2. Dutta, P.C. and Appelqvist, L.Å. Saturated sterols (stanols) in unhydrogenated and hydrogenated edible vegetable oils and in cereal lipids. *J. Sci. Fd. Agric.* **71**(3), 383-391 (1996).
3. Sims, R.J., Fioriti, J.A., and Kanuk, M.J. Sterol additives as polymerization inhibitors for frying oils. *J. Am. Oil Chem. Soc.* **49**, 298-301 (1972).

#### 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, 12/05/2022

#### 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