

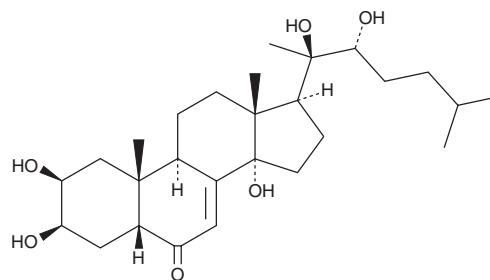
# PRODUCT INFORMATION



## Ponasterone A

Item No. 16386

**CAS Registry No.:** 13408-56-5  
**Formal Name:** (5 $\beta$ )-2 $\beta$ ,3 $\beta$ ,14,20,22R-pentahydroxy-cholest-7-en-6-one  
**Synonym:** 25-Deoxyecdysterone  
**MF:** C<sub>27</sub>H<sub>44</sub>O<sub>6</sub>  
**FW:** 464.6  
**Purity:**  $\geq$ 95%  
**Supplied as:** A powder  
**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

Ponasterone A is supplied as a powder. A stock solution may be made by dissolving the ponasterone A in the solvent of choice. Ponasterone A is soluble in organic solvents such as ethanol, methanol, and DMSO, which should be purged with an inert gas.

### Description

Ponasterone A is an analog of 20-hydroxy ecdysone (Item No. 16145), the insect steroid hormone that regulates the metamorphosis of *Drosophila*.<sup>1</sup> Ecdysteroids such as this compound have been employed as inducers of ecdysone-inducible mammalian expression systems.<sup>2,3</sup>

### References

1. Fallon, A.M. and Gerenday, A. Ecdysone and the cell cycle: Investigations in a mosquito cell line. *J. Insect Physiol.* **56(10)**, 1396-1401 (2010).
2. Saez, E., Nelson, M.C., Eshelman, B., *et al.* Identification of ligands and coligands for the ecdysone-regulated gene switch. *Proc. Natl. Acad. Sci. USA* **97(26)**, 14512-14517 (2000).
3. Karzenowski, D., Potter, D.W., and Padidam, M. Inducible control of transgene expression with ecdysone receptor: Gene switches with high sensitivity, robust expression, and reduced size. *Biotechniques* **39(2)**, 191-192 (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, 04/24/2018

#### 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