

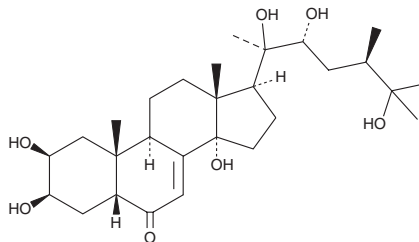
PRODUCT INFORMATION



Makisterone A

Item No. 11739

CAS Registry No.: 20137-14-8
Formal Name: (2 β ,3 β ,5 β ,22R,24R)-2,3,14,20,22,25-hexahydroxy-ergost-7-en-6-one
MF: C₂₈H₄₆O₇
FW: 494.7
Purity: $\geq 95\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years



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

Laboratory Procedures

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

Description

Makisterone A is an ecdysteroid that drives molting in insects.^{1,2} It binds the heterodimeric ecdysone receptor with nanomolar affinity.³

References

1. Cottam, D. M. and Milner, M. J. The effects of several ecdysteroids and ecdysteroid agonists on two *Drosophila* imaginal disc cell lines. *Cell. Mol. Life Sci.* **53**(7), 600-603 (1997).
2. Lavrynenko, O., Rodenfels, J., Carvalho, M., et al. The ecdysteroidome of *Drosophila*: Influence of diet and development. *Development* **142**(21), 3758-3768 (2015).
3. Tohidi-Esfahani, D., Graham, L. D., Hannan, G. N., et al. An ecdysone receptor from the pentatomomorph, *Nezara viridula*, shows similar affinities for moulting hormones makisterone A and 20-hydroxyecdysone. *Insect. Biochem. Mol. Biol.* **41**(2), 77-89 (2011).

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, 11/14/2022

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