

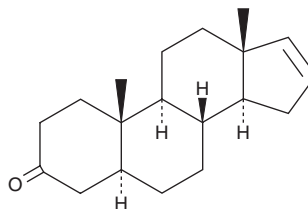
PRODUCT INFORMATION



5 α -Androst-16-en-3-one

Item No. 18233

CAS Registry No.: 18339-16-7
Formal Name: (5 α)-androst-16-en-3-one
Synonym: Androstenone
MF: C₁₉H₂₈O
FW: 272.4
Purity: $\geq 98\%$
Supplied as: A crystalline 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

5 α -Androst-16-en-3-one is supplied as a crystalline solid. A stock solution may be made by dissolving the 5 α -androst-16-en-3-one in the solvent of choice. 5 α -Androst-16-en-3-one is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of 5 α -androst-16-en-3-one in these solvents is approximately 10, 15, and 25 mg/ml, respectively.

5 α -Androst-16-en-3-one is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 5 α -androst-16-en-3-one should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. 5 α -Androst-16-en-3-one has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

5 α -Androst-16-en-3-one is a mammalian pheromone released as a volatile chemical cue in boar saliva to facilitate social and sexual interactions.¹ It has been used to prime sexual behavior of sows in estrus for mating or artificial insemination. 5 α -Androst-16-en-3-one is also found in human sweat and urine and has been used to study receptor-mediated odorant detection and the genetic basis for anosmias.²

References

1. Dehnard, M., Rohrmann, H., and Kauffold, J. Measurement of 16-androstenes (5 α -androst-16-en-3-one, 5 α -androst-16-en-3 α -ol, 5 α -androst-16-en-3 β -ol) in saliva of German landrace and Göttingen minipig boars, Chapter 30, in Chemical Signals in Vertebrates 12. East, M.L. and Dehnard, M., editors, 12, Springer-Verlag, New York, 381-390 (2013).
2. Wang, L., Chen, L., and Jacob, T. Evidence for peripheral plasticity in human odour response. *J. Physiol.* **554.1**, 236-244 (2003).

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, 10/31/2022

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