

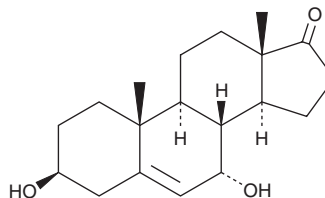
PRODUCT INFORMATION



7 α -hydroxy Dehydroepiandrosterone

Item No. 34268

CAS Registry No.: 53-00-9
Formal Name: 3 β ,7 α -dihydroxy-androst-5-en-17-one
Synonym: 7 α -hydroxy DHEA
MF: C₁₉H₂₈O₃
FW: 304.4
Purity: \geq 98%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

7 α -hydroxy Dehydroepiandrosterone (7 α -hydroxy DHEA) is supplied as a solid. A stock solution may be made by dissolving the 7 α -hydroxy DHEA in the solvent of choice, which should be purged with an inert gas. 7 α -hydroxy DHEA is soluble in the organic solvent ethanol.

Description

7 α -hydroxy DHEA is an active metabolite of the endogenous steroid hormone dehydroepiandrosterone (Item No. 15728).¹ 7 α -hydroxy DHEA is formed from dehydroepiandrosterone via 7-hydroxylation by cytochrome P450s (CYPs).² It increases antigen-induced production of anti-tetanus toxoid and anti-pertussis antigen IgGs in co-cultures of human tonsil-derived B and T cells when used at a concentration of 1 μ M.³ 7 α -hydroxy DHEA (6.25 mg/kg) increases serum concentrations of anti-lysozyme IgG in mice.¹

References

1. Morfin, R. and Courchay, G. Pregnenolone and dehydroepiandrosterone as precursors of native 7-hydroxylated metabolites which increase the immune response in mice. *J. Steroid Biochem. Mol. Biol.* **50(1-2)**, 91-100 (1994).
2. Hampl, R., Morfin, R., and Starka, L. 7-Hydroxylated derivatives of dehydroepiandrosterone: What are they good for? *Endocr. Regul.* **31(4)**, 211-218 (1997).
3. Lafaye, P., Chmielewski, V., Nato, F., et al. The 7 α -hydroxysteroids produced in human tonsils enhance the immune response to tetanus toxoid and *Bordetella pertussis* antigens. *Biochim. Biophys. Acta* **1472(1-2)**, 222-231 (1999).

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, 01/03/2023

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