

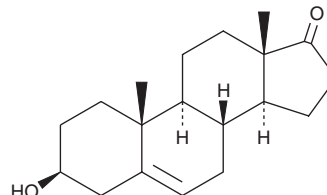
PRODUCT INFORMATION



Dehydroepiandrosterone

Item No. 15728

CAS Registry No.: 53-43-0
Formal Name: 5-androsten-3 β -ol-17-one
Synonyms: Androstenedione, *trans*-Dehydroandrosterone, DHEA, Diandron, NSC 9896, Prasterone, Psicosterone
MF: C₁₉H₂₈O₂
FW: 288.4
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

Dehydroepiandrosterone (DHEA) is supplied as a crystalline solid. A stock solution may be made by dissolving the DHEA in the solvent of choice, which should be purged with an inert gas. DHEA is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of DHEA in these solvents is approximately 10, 15, and 25 mg/ml, respectively.

DHEA is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, DHEA should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. DHEA 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

DHEA is an endogenous steroid hormone that is secreted primarily by the adrenal gland and is the most abundant sex steroid.¹⁻³ Metabolites of DHEA include androstenedione (Item No. 9001311), which subsequently may be metabolized to testosterone (Item Nos. ISO60154 | 15645) or estrone (Item No. 10006485), which is an estradiol (Item Nos. ISO60155 | 10006315) precursor.¹⁻³ In addition to serving as an intermediate in the biosynthesis of sex steroids, DHEA directly modulates a number of cellular and nuclear receptors.¹⁻³

References

1. Kim, C. and Halter, J.B. Endogenous sex hormones, metabolic syndrome, and diabetes in men and women. *Curr. Cardiol. Rep.* **16**(4), (2014).
2. Mannic, T., Viguie, J., and Rossier, M.F. *In vivo* and *in vitro* evidences of dehydroepiandrosterone protective role on the cardiovascular system. *Int. J. Endocrinol. Metab.* **13**(2), (2015).
3. Samaras, N., Papadopoulou, M.-A., Samaras, D., *et al.* Off-label use of hormones as an antiaging strategy: A review. *Clin. Interv. Aging* **9**, 1175-1186 (2014).

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, 02/06/2024

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