

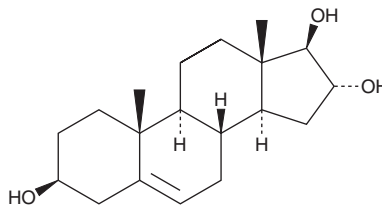
PRODUCT INFORMATION



5-Androstenetriol

Item No. 35273

CAS Registry No.: 4150-30-5
Formal Name: androst-5-ene-3 β ,16 α ,17 β -triol
Synonym: Δ^5 -AT
MF: C₁₉H₃₀O₃
FW: 306.4
Purity: $\geq 98\%$
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

5-Androstenetriol is supplied as a solid. A stock solution may be made by dissolving the 5-androstenetriol in the solvent of choice, which should be purged with an inert gas. 5-Androstenetriol is soluble in ethanol and DMSO.

Description

5-Androstenetriol is a steroid and an active metabolite of dehydroepiandrosterone (DHEA; Item No. 15728).¹ *In vivo*, 5-androstenetriol (0.75 mg/25 g body weight) reduces radiation-induced lethality and induces hematopoietic recovery from surviving stem cells in mice. It increases survival in a rat model of surgical trauma-induced sepsis when administered at a dose of 40 mg/kg.² Urinary levels of 5-androstenetriol are increased in women with androgenic alopecia.³

References

1. Loria, R.M., Conrad, D.H., Huff, T., *et al.* Androstenetriol and androstenediol. Protection against lethal radiation and restoration of immunity after radiation injury. *Ann. N. Y. Acad. Sci.* **917**(1), 860-867 (2000).
2. Marcu, A.C., Kielar, N.D., Paccione, K.E., *et al.* Androstenetriol improves survival in a rodent model of traumatic shock. *Resuscitation* **71**(3), 379-386 (2006).
3. Juricksay, S. and Telegdy, E. Urinary steroids in women with androgenic alopecia. *Clin. Biochem.* **33**(2), 97-101 (2000).

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, 10/12/2022

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