

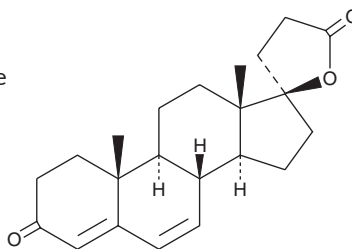
PRODUCT INFORMATION



Canrenone

Item No. 21307

CAS Registry No.: 976-71-6
Formal Name: 17 α -hydroxy-3-oxo-pregna-4,6-diene-21-carboxylic acid γ -lactone
Synonyms: Aldadiene, NSC 261713, SC-9376, SC-14266
MF: C₂₂H₂₈O₃
FW: 340.5
Purity: \geq 98%
UV/Vis.: λ_{max} : 210, 281 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 5 years



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

Description

Canrenone (Item No. 21307) is an analytical reference standard categorized as a steroid.¹ It is an active metabolite of spironolactone (Item No. 9000324) that acts as a diuretic. Diuretics, including canrenone, have been abused as performance-enhancing drugs and masking agents in sports doping.² This product is intended for research and forensic applications.

References

1. Larik, F.A., Saaed, A., Shahzad, D., *et al.* Synthetic approaches towards the multi target drug spironolactone and its potent analogues/derivatives. *Steroids* **118**, 76-92 (2017).
2. Cadwallader, A.B., de la Torre, X., Tieri, A., *et al.* The abuse of diuretics as performance-enhancing drugs and masking agents in sport doping: Pharmacology, toxicology and analysis. *Br. J. Pharmacol.* **161**(1), 1-16 (2010).

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, 06/28/2023

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