

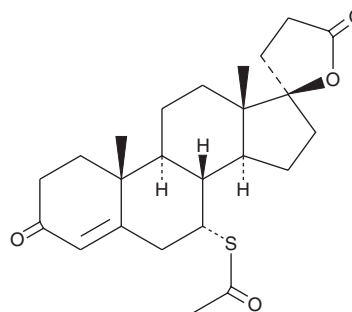
PRODUCT INFORMATION



Spironolactone

Item No. 9000324

CAS Registry No.: 52-01-7
Formal Name: (7 α ,17 α)-7-(acetylthio)-17-hydroxy-3-oxo-pregn-4-ene-21-carboxylic acid- γ -lactone
Synonyms: NSC 150399, SC-9420
MF: C₂₄H₃₂O₄S
FW: 416.6
Purity: \geq 98%
UV/Vis.: λ_{\max} : 236 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

Spironolactone (Item No. 9000324) is an analytical reference standard categorized as a steroid.^{1,2} It is a diuretic that has been abused as a performance-enhancing drug and masking agent in sports doping.³ This product is intended for research and forensic applications

References

1. Chun, T.Y., Bloem, L.J., and Pratt, J.H. Aldosterone inhibits inducible nitric oxide synthase in neonatal rat cardiomyocytes. *Endocrinology* **144**(5), 1712-1717 (2003).
2. 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).
3. 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, 07/12/2023

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