

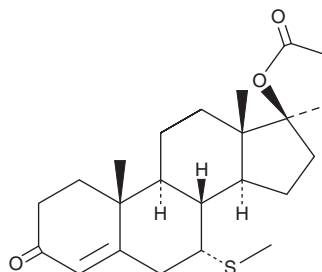
PRODUCT INFORMATION



7 α -Thiomethylspironolactone

Item No. 34526

CAS Registry No.: 38753-77-4
Formal Name: (7 α ,17 α)-17-hydroxy-7-(methylthio)-3-oxo-pregn-4-ene-21-carboxylic acid, γ -lactone
Synonym: SC 26519
MF: C₂₃H₃₂O₃S
FW: 388.6
Purity: \geq 95%
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 -Thiomethylspironolactone is supplied as a solid. A stock solution may be made by dissolving the 7 α -thiomethylspironolactone in the solvent of choice, which should be purged with an inert gas. 7 α -Thiomethylspironolactone is slightly soluble in chloroform and ethyl acetate.

Description

7 α -Thiomethylspironolactone is a major metabolite of the synthetic steroid spironolactone.^{1,2}

References

1. Jankowski, A., Skorek-Jankowska, A., and Lamparczyk, H. Simultaneous determination of spironolactone and its metabolites in human plasma. *J. Pharm. Biomed. Anal.* **14(8-10)**, 1359-1365 (1996).
2. Gardiner, P., Schrode, K., Quinlan, D., et al. Spironolactone metabolism: Steady-state serum levels of the sulfur-containing metabolites. *J. Clin. Pharmacol.* **29(4)**, 342-347 (1989).

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, 11/22/2022

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