

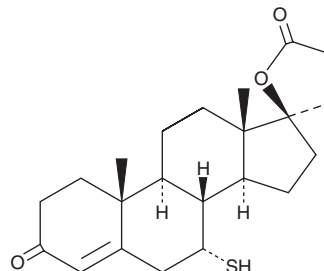
PRODUCT INFORMATION



7 α -Thiospironolactone

Item No. 34868

CAS Registry No.: 38753-76-3
Formal Name: 17 α -hydroxy-7 α -mercapto-3-oxo-pregn-4-ene-21-carboxylic acid, γ -lactone
Synonyms: Deacetylspironolactone, Mercaptospironolactone, SC-24813, 7 α -Thio-SL
MF: C₂₂H₃₀O₃S
FW: 374.5
Purity: $\geq 90\%$
UV/Vis.: λ_{max} : 237 nm
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

7 α -Thiospironolactone (7 α -Thio-SL) is supplied as a solid. A stock solution may be made by dissolving the 7 α -thio-SL in the solvent of choice, which should be purged with an inert gas. 7 α -Thio-SL is soluble in the organic solvent ethanol.

Description

7 α -Thio-SL is an active metabolite of the synthetic steroid spironolactone.¹ *In vivo*, 7 α -thio-SL (100 mg/kg) decreases plasma testosterone and estradiol levels in male dogs.² It reduces mercuric chloride-induced nephrocalcinosis and mortality in rats.³

References

1. Los, L.E., Coddington, A.B., Ramjit, H.G., *et al.* Identification of spironolactone metabolites in plasma and target organs of guinea pigs. *Drug Metab. Dispos.* **21**(6), 1086-1090 (1993).
2. Menard, R.H., Loriaux, D.L., Bartter, F.C., *et al.* The effect of the administration of spironolactone on the concentration of plasma testosterone, estradiol and cortisol in male dogs. *Steroids* **31**(6), 771-782 (1978).
3. Szabo, S. and Gagnon, M.J. The protective effect of sulfur-containing steroids against nephrocalcinosis induced by mercuric chloride in rats. *Gen. Pharmacol.* **14**(2), 269-272 (1983).

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, 11/04/2022

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