

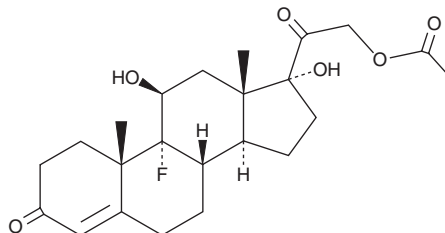
PRODUCT INFORMATION



Fludrocortisone Acetate

Item No. 24032

CAS Registry No.: 514-36-3
Formal Name: 21-(acetyloxy)-9-fluoro-11 β ,17-dihydroxy-pregn-4-ene-3,20-dione
Synonyms: 9 α -fluoro Hydrocortisone Acetate, NSC 15186, U 4845
MF: C₂₃H₃₁FO₆
FW: 422.5
Purity: \geq 98%
UV/Vis.: λ_{max} : 239 nm
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

Fludrocortisone acetate is supplied as a solid. A stock solution may be made by dissolving the fludrocortisone acetate in the solvent of choice, which should be purged with an inert gas. Fludrocortisone acetate is slightly soluble in chloroform and methanol.

Description

Fludrocortisone acetate is the acetate form of the synthetic corticosteroid fludrocortisone, which is a mineralocorticoid receptor agonist.¹ Fludrocortisone acetate (0.1-5 μ g/animal) promotes sodium retention in rats in a dose-dependent manner with 37% of sodium excreted compared with control when used at a dose of 5 μ g/animal.² Formulations containing fludrocortisone acetate have been used in the treatment of Addison's disease.

References

1. Agarwal, M.K., Coupry, F. and Philippe, M. Physiological activity and receptor binding of 9 α fluorohydrocortisone. *Biochem. Biophys. Res. Commun.* **78(2)**, 747-753 (1977).
2. Stafford, R.O., Barnes, L.E., Bowman, B.J., *et al.* Glucocorticoid and mineralocorticoids activities of Delta1-fluorohydrocortisone. *Proc. Soc. Exp. Biol. Med.* **89(3)**, 371-374 (1955).

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, 11/14/2022

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