

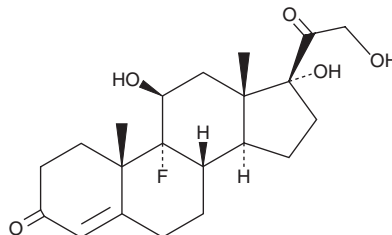
PRODUCT INFORMATION



Fludrocortisone

Item No. 29391

CAS Registry No.: 127-31-1
Formal Name: (11 β)-9-fluoro-11,17,21-trihydroxy-pregn-4-ene-3,20-dione
Synonyms: Astonin-H, 9 α -Fludrocortisone, Fluohydrocortisone, Fluorohydrocortisone, NSC 11318
MF: C₂₁H₂₉FO₅
FW: 380.5
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

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

Description

Fludrocortisone is a synthetic mineralocorticoid.¹ Dietary administration of fludrocortisone (10 μ g/kg per day) reduces tympanic membrane thickness and fluid area in a mouse model of *S. pneumoniae*-induced middle ear inflammation. Formulations containing fludrocortisone have been used in the treatment of Addison's disease.

Reference

1. MacArthur, C.J., DeGagne, J.M., Kempton, J.B., *et al.* Steroid control of acute middle ear inflammation in a mouse model. *Arch. Otolaryngol. Head Neck Surg.* **135(5)**, 453-457 (2009).

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, 12/15/2022

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