

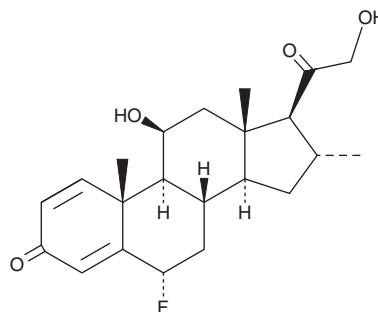
PRODUCT INFORMATION



Fluocortolone

Item No. 35246

CAS Registry No.: 152-97-6
Formal Name: 6 α -fluoro-11 β ,21-dihydroxy-16 α -methyl-pregna-1,4-diene-3,20-dione
Synonym: SH 742
MF: C₂₂H₂₉FO₄
FW: 376.5
Purity: $\geq 95\%$
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

Fluocortolone is supplied as a solid. A stock solution may be made by dissolving the fluocortolone in the solvent of choice, which should be purged with an inert gas. Fluocortolone is soluble in organic solvents such as chloroform and methanol. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

Fluocortolone is a glucocorticoid.^{1,2} It inhibits progesterone 3 α -hydroxysteroid dehydrogenase in rat cortical homogenates (IC₅₀ = 40 μ M).¹ Fluocortolone (3, 6, and 12 mg/kg) decreases adjuvant-induced paw edema in rats.²

References

1. Stürenburg, H.J., Fries, U., and Kunze, K. Glucocorticoids and anabolic/androgenic steroids inhibit the synthesis of GABAergic steroids in rat cortex. *Neuropsychobiology* **35**(3), 143-146 (1997).
2. Bhargava, A.S. Effect of anti-inflammatory agents on adjuvant-induced edema modified for primary test. *Pharmacol. Res. Commun.* **3**(1), 83-91 (1971)

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, 11/03/2022

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