

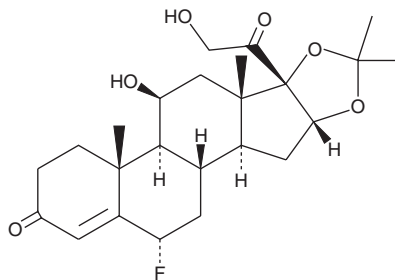
PRODUCT INFORMATION



Flurandrenolide

Item No. 20933

CAS Registry No.: 1524-88-5
Formal Name: (6 α ,11 β ,16 α)-6-fluoro-11,21-dihydroxy-16,17-[(1-methylethylidene)bis(oxy)]-pregn-4-ene-3,20-dione
Synonyms: Fludroxycortide, Flurandrenolone, L-33,379
MF: C₂₄H₃₃FO₆
FW: 436.5
Purity: \geq 98%
UV/Vis.: λ_{max} : 235 nm
Supplied as: A crystalline 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

Flurandrenolide is supplied as a crystalline solid. A stock solution may be made by dissolving the flurandrenolide in the solvent of choice, which should be purged with an inert gas. Flurandrenolide is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of flurandrenolide in ethanol and DMF is approximately 15 mg/ml and approximately 1 mg/ml in DMSO.

Description

Flurandrenolide is a topical corticosteroid with anti-inflammatory actions.¹ It has also been shown to activate EGFR with an EC₅₀ value of 23 nM.²

References

1. Boris, A., and Hurley, J.F. Assessment of topical anti-inflammatory activity in rats with cantharidin-induced inflammation. *J. Invest. Dermatol.* **68(3)**, 161-164 (1977).
2. Antczak, C., Mahida, J.P., Bhinder, B., *et al.* A high-content biosensor-based screen identifies cell-permeable activators and inhibitors of EGFR function: Implications in drug discovery. *Journal of Biomolecular Screening* **17(7)**, 885-899 (2012).

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, 10/10/2022

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