

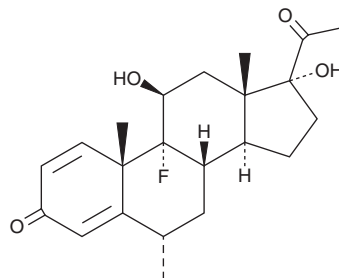
# PRODUCT INFORMATION



## Fluorometholone

Item No. 9000157

**CAS Registry No.:** 426-13-1  
**Formal Name:** 9-fluoro-11 $\beta$ ,17-dihydroxy-6 $\alpha$ -methyl-pregna-1,4-diene-3,20-dione  
**Synonyms:** NSC 33001, U-8614  
**MF:** C<sub>22</sub>H<sub>29</sub>FO<sub>4</sub>  
**FW:** 376.5  
**Purity:**  $\geq$ 98%  
**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

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

### Description

Fluorometholone is a synthetic glucocorticoid that binds to glucocorticoid receptors ( $IC_{50}$  = 1.5 nM;  $K_d$  = 2.8 nM in a radioligand binding assay).<sup>1,2</sup> Fluorometholone (0.01%) decreases survival of cultured corneal epithelial cells and inhibits their migration in a scratch assay following 12 hours of treatment.<sup>3</sup> Formulations containing fluorometholone have been used for the treatment of eye inflammation.

### References

1. Southren, A.L., Dominguez, M.O., Gordon, G.G., *et al.* Nuclear translocation of the cytoplasmic glucocorticoid receptor in the iris-ciliary body and adjacent corneoscleral tissue of the rabbit following topical administration of various glucocorticoids. A rapid screening method for glucocorticoid activity. *Invest. Ophthalmol. Vis. Sci.* **24**(2), 147-152 (1983).
2. Lin, M.T. Demonstration of specific glucocorticoid binding sites in bovine cornea. *Exp. Eye Res.* **38**(4), 333-339 (1984).
3. Lee, L.S., Kim, Y.H., and Park, Y.M. The toxicity of nonsteroidal anti-inflammatory eye drops against human corneal epithelial cells in vitro. *J. Korean Med. Sci.* **30**(12), 1856-1864 (2015).

#### 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, 11/09/2022

#### 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](http://WWW.CAYMANCHEM.COM)