

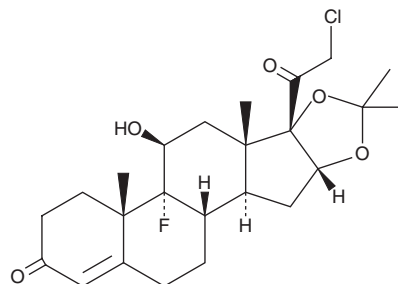
# PRODUCT INFORMATION



## Halcinonide

Item No. 23780

**CAS Registry No.:** 3093-35-4  
**Formal Name:** 21-chloro-9-fluoro-11 $\beta$ -hydroxy-16 $\alpha$ ,17-[(1-methylethylidene)bis(oxy)]-pregn-4-ene-3,20-dione  
**Synonym:** SQ 18,566  
**MF:** C<sub>24</sub>H<sub>32</sub>ClFO<sub>5</sub>  
**FW:** 455.0  
**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

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

### Description

Halcinonide is a synthetic glucocorticoid with anti-inflammatory properties.<sup>1,2</sup> In a rabbit model of edema, topical application of halcinonide (0.1% w/w) decreases edema by 55% in a reversed passive Arthus skin test (RPA).<sup>1</sup> Halcinonide (0.1% w/w) inhibits croton oil-induced ear edema by 70.1% and homologous passive cutaneous anaphylaxis (PCA) by 88.0% in rats dosed three hours prior to induction of inflammation.<sup>2</sup> Application of halcinonide (0.1% w/w) to the abraded surface of cecum in rats inhibits leukocyte adhesion (43.3% incidence *versus* 100% for controls) and lowers the total number of intraperitoneal leukocytes from 270 million to 121 million 3 days post surgery.<sup>3</sup> It also acts as an agonist at the smoothened (Smo) receptor with an EC<sub>50</sub> value of 1,100 nM.<sup>4</sup> Formulations containing halcinonide have been used to treat inflammatory and corticosteroid-responsive dermatoses.

### References

1. Goldlust, M.B., Palmer, D.M., and Augustine, M.A. Evaluation of topical anti-inflammatory steroid formulations in an Arthus model of inflammation. *J. Invest. Dermatol.* **66**(3), 157-160 (1976).
2. Iizuka, Y., Endo, Y., and Misawa, Y. Two simple methods for the evaluation of topically active anti-inflammatory steroidal ointments. *Agents Actions* **11**(3), 254-259 (1981).
3. Zhang, Y.-D., Yao, W., Wu, C.-X., *et al.* Topical application of halcinonide cream reduces the severity and incidence of intraperitoneal adhesions in a rat model. *Am. J. Surg.* **184**(1), 74-77 (2002).
4. Wang, J., Lu, J., Bond, M.C., *et al.* Identification of select glucocorticoids as smoothened agonists: Potential utility for regenerative medicine. *Proc. Nat. Acad. Sci. USA* **107**(20), 9323-9328 (2010).

#### 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, 11/01/2022

#### 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)