

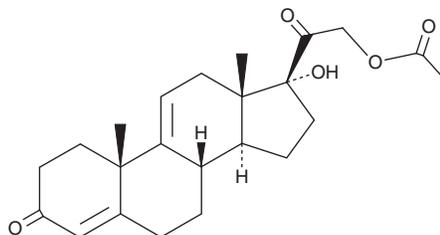
# PRODUCT INFORMATION



## Anecortave Acetate

Item No. 36441

**CAS Registry No.:** 7753-60-8  
**Formal Name:** 21-(acetyloxy)-17-hydroxy-pregna-4,9(11)-diene-3,20-dione  
**Synonyms:** AL 3789, NSC 15475, NSC 24345  
**MF:** C<sub>23</sub>H<sub>30</sub>O<sub>5</sub>  
**FW:** 386.5  
**Purity:** ≥98%  
**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

Anecortave acetate is supplied as a solid. A stock solution may be made by dissolving the anecortave acetate in the solvent of choice, which should be purged with an inert gas. Anecortave acetate is soluble in the organic solvent chloroform at a concentration of approximately 10 mg/ml.

### Description

Anecortave acetate is a sterol lipid.<sup>1</sup> It inhibits VEGF-induced increases in vascular generation, number, and length in isolated rat retinas when used at a concentration of 1 μM. Sub-tenon administration of anecortave acetate (7.5% w/v) prevents and reduces increases in intraocular pressure induced by prednisolone (Item No. 20866) in sheep.<sup>2</sup> Subconjunctival administration of anecortave acetate (15% w/v), in combination with 2-deoxyglucose, decreases tumor growth in a Tg(TagRb)1Plm mouse model of retinoblastoma.<sup>3</sup>

### References

1. Unoki, N., Murakami, T., Ogino, K., *et al.* Time-lapse imaging of retinal angiogenesis reveals decreased development and progression of neovascular sprouting by anecortave desacetate. *Invest. Ophthalmol. Vis. Sci.* **51(5)**, 2347-2355 (2010).
2. Candia, O.A., Gerometta, R., Millar, J.C., *et al.* Suppression of corticosteroid-induced ocular hypertension in sheep by anecortave. *Arch. Ophthalmol.* **128(3)**, 338-343 (2010).
3. Houston, S.K., Pina, Y., Murray, T.G., *et al.* Novel retinoblastoma treatment avoids chemotherapy: The effect of optimally timed combination therapy with angiogenic and glycolytic inhibitors on LH<sub>BETA</sub> T<sub>AG</sub> retinoblastoma tumors. *Clin. Ophthalmol.* **5**, 129-137 (2011).

#### 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, 05/10/2023

#### 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