

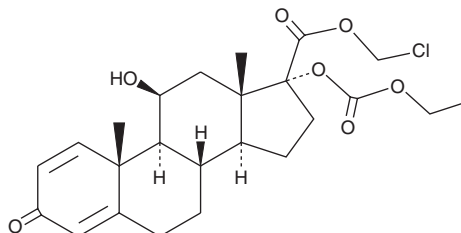
# PRODUCT INFORMATION



## Loteprednol etabonate

Item No. 23305

**CAS Registry No.:** 82034-46-6  
**Formal Name:** (11 $\beta$ ,17 $\alpha$ )-17-[(ethoxycarbonyl)oxy]-11-hydroxy-3-oxo-androsta-1,4-diene-17-carboxylic acid, chloromethyl ester  
**Synonym:** CDDD 5604  
**MF:** C<sub>24</sub>H<sub>31</sub>ClO<sub>7</sub>  
**FW:** 467.0  
**Purity:**  $\geq$ 98%  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 245 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

Loteprednol etabonate is supplied as a crystalline solid. A stock solution may be made by dissolving the loteprednol etabonate in the solvent of choice. Loteprednol etabonate is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of loteprednol etabonate in these solvents is approximately 30 mg/ml.

Loteprednol etabonate is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, loteprednol etabonate should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Loteprednol etabonate has a solubility of approximately 0.2 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Loteprednol etabonate is a glucocorticoid receptor agonist with an affinity 4.3 times that of dexamethasone (Item No. 11015).<sup>1</sup> It decreases wound healing time in mice and minimizes scarring of rabbit cornea wounds.<sup>2,3</sup> Formulations containing loteprednol etabonate are used in the treatment of inflammatory eye conditions.

### References

1. Druzgala, P., Hochhaus, G., and Bodor, N. Soft drugs--10. Blanching activity and receptor binding affinity of a new type of glucocorticoid: loteprednol etabonate. *J. Steroid Biochem. Mol. Biol.* **38(2)**, 149-154 (1991).
2. Bodor, N.S., Kiss-Buris, S.T., and Buris, L. Novel soft steroids: Effects on cell growth *in vitro* and on wound healing in the mouse. *Steroids* **56(8)**, 434-439 (1991).
3. Bodor, N. and Varga, M. Effect of a novel soft steroid on the wound healing of rabbit cornea. *Exp. Eye Res.* **50(2)**, 183-187 (1990).

#### 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, 10/24/2022

#### 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