

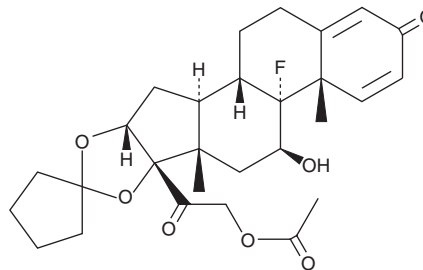
# PRODUCT INFORMATION



## Amcinonide

Item No. 23903

**CAS Registry No.:** 51022-69-6  
**Formal Name:** (11 $\beta$ ,16 $\alpha$ )-21-(acetyloxy)-16,17-[cyclopentylidenebis(oxy)]-9-fluoro-11-hydroxy-pregna-1,4-diene-3,20-dione  
**Synonym:** CL 34,699  
**MF:** C<sub>28</sub>H<sub>35</sub>FO<sub>7</sub>  
**FW:** 502.6  
**Purity:**  $\geq$ 98%  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 236 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

Amcinonide is supplied as a crystalline solid. A stock solution may be made by dissolving the amcinonide in the solvent of choice, which should be purged with an inert gas. Amcinonide is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of amcinonide in these solvents is approximately 10, 25, and 20 mg/ml, respectively.

Amcinonide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, amcinonide should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Amcinonide has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Amcinonide is a synthetic corticosteroid.<sup>1</sup> Topical administration of amcinonide reduces inflammation in rat models of carrageenan-induced plantar edema and air pouch granulomas (ID<sub>50</sub>s = 1.85 and 4.2 mg/kg, respectively).

### Reference

1. Wepierre, J. and Bogaievsky, Y. Comparative pharmacological activity of triamcinolone acetone and of a new topical steroid: Amcinonide, after percutaneous or oral administration (author's transl). *J. Pharmacol.* **11**(4), 449-462 (1980).

#### 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, 12/06/2022

#### 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