

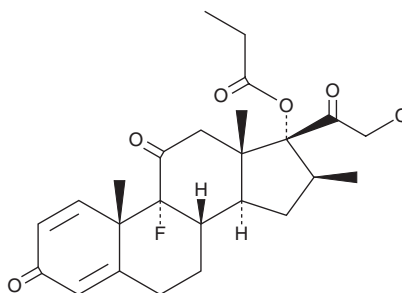
# PRODUCT INFORMATION



## Clobetasone Propionate

Item No. 35747

**CAS Registry No.:** 25122-56-9  
**Formal Name:** 21-chloro-9-fluoro-16 $\beta$ -methyl-17-(1-oxopropoxy)-pregna-1,4-diene-3,11,20-trione  
**Synonym:** Clobetasone 17-Propionate  
**MF:** C<sub>25</sub>H<sub>30</sub>ClFO<sub>5</sub>  
**FW:** 465.0  
**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

Clobetasone propionate is supplied as a solid. A stock solution may be made by dissolving the clobetasone propionate in the solvent of choice, which should be purged with an inert gas. Clobetasone propionate is slightly soluble in acetone, chloroform, and DMSO.

### Description

Clobetasone propionate is a corticosteroid.<sup>1</sup> Formulations containing clobetasol propionate have been used in the treatment of eczema, psoriasis, and lupus.

### Reference

1. Brandi, N., Starace, M., Alessandrini, A., *et al.* Tinea versicolor of the neck as side effect of topical steroids for alopecia areata. *J. Dermatolog. Treat.* **30(8)**, 757-759 (2019).

#### 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, 12/13/2022

#### 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