

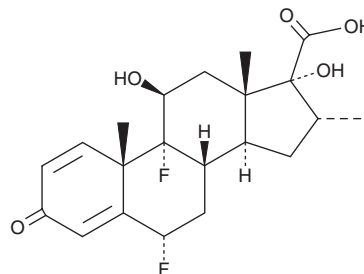
# PRODUCT INFORMATION



## Fluticasone 17 $\beta$ -Carboxylic Acid

Item No. 27855

**CAS Registry No.:** 28416-82-2  
**Formal Name:** (6 $\alpha$ ,11 $\beta$ ,16 $\alpha$ ,17 $\alpha$ )-6,9-difluoro-11,17-dihydroxy-16-methyl-3-oxo-androsta-1,4-diene-17-carboxylic acid  
**MF:** C<sub>21</sub>H<sub>26</sub>F<sub>2</sub>O<sub>5</sub>  
**FW:** 396.4  
**Purity:**  $\geq$ 98%  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 216, 242 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

Fluticasone 17 $\beta$ -carboxylic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the fluticasone 17 $\beta$ -carboxylic acid in the solvent of choice, which should be purged with an inert gas. Fluticasone 17 $\beta$ -carboxylic acid is soluble in the organic solvent DMSO at a concentration of approximately 5 mg/ml.

Fluticasone 17 $\beta$ -carboxylic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, fluticasone 17 $\beta$ -carboxylic acid should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Fluticasone 17 $\beta$ -carboxylic acid has a solubility of approximately 0.10 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Fluticasone 17 $\beta$ -carboxylic acid is an intermediate in the synthesis of the glucocorticosteroid fluticasone propionate (Item No. 20703).<sup>1</sup>

### Reference

1. Zhou, J., Jin, C., and Su, W. Improved synthesis of fluticasone propionate. *Org. Process Res. Dev.* **18**(8), 928-933 (2014).

#### 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, 10/06/2022

#### 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