

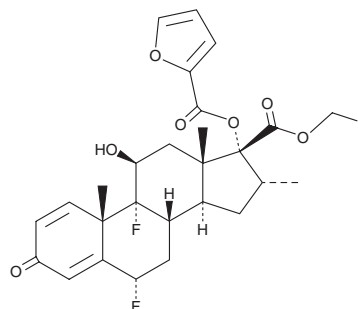
PRODUCT INFORMATION



Fluticasone O-fluoro methyl ester

Item No. 45756

CAS Registry No.: 572880-17-2
Formal Name: 6 α ,9-difluoro-17-[(2-furanylcarbonyl)oxy]-11 α -hydroxy-16 α -methyl-3-oxo-androsta-1,4-diene-17 α -carboxylic acid, fluoromethyl ester
Synonym: VSG158
MF: C₂₇H₂₉F₃O₇
FW: 522.5
Purity: $\geq 95\%$
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

Fluticasone O-fluoro methyl ester is supplied as a solid. A stock solution may be made by dissolving the fluticasone O-fluoro methyl ester in the solvent of choice, which should be purged with an inert gas. Fluticasone O-fluoro methyl ester is soluble (≥ 10 mg/ml) in DMSO.

Description

Fluticasone O-fluoro methyl ester is a synthetic glucocorticoid.¹ It binds to the glucocorticoid receptor ($K_i = 1.7$ nM) and inhibits transcription mediated by activator protein-1 (AP-1) in AD-293 cells. Intranasal administration of fluticasone O-fluoro methyl ester (0.125 mg/kg) represses airway hyperresponsiveness in a mouse model of ovalbumin-induced acute asthma. It reduces lung levels of inflammatory cytokines and bronchoalveolar lavage fluid (BALF) eosinophil infiltration in a mouse model of steroid-susceptible inflammation and reduces BALF neutrophil and CD4⁺ T cell infiltration in a mouse model of steroid-resistant inflammation.

Reference

1. He, Y., Shi, J., Nguyen, Q.T., *et al.* Development of highly potent glucocorticoids for steroid-resistant severe asthma. *Proc. Natl. Acad. Sci. USA* **116**(14), 6932-6937 (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, 07/13/2026

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