

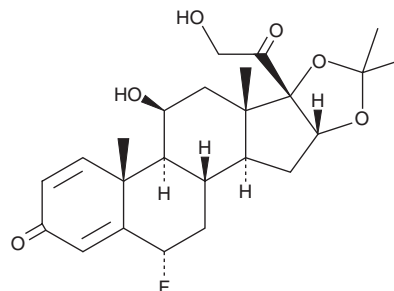
PRODUCT INFORMATION



Flunisolide

Item No. 23842

CAS Registry No.: 3385-03-3
Formal Name: 6 α -fluoro-11 β ,21-dihydroxy-16 α ,17-[(1-methylethylidene)bis(oxy)]-pregna-1,4-diene-3,20-dione
Synonym: RS 3999
MF: C₂₄H₃₁FO₆
FW: 434.5
Purity: \geq 95%
UV/Vis.: λ_{max} : 238 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

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

Flunisolide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, flunisolide should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Flunisolide has a solubility of approximately 0.2 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Flunisolide is a fluorinated corticosteroid receptor agonist that induces reversible pyloric hyalinization in CD-1 mice following oral administration at a dose of 200 μ g/kg per day.¹ Formulations containing flunisolide have been used for the treatment of asthma.²

References

- McKevitt, T.P., Giffen, P., Woodfine, J.A., *et al.* Hyalinization of the pyloric stomach in CD-1 mice following oral (dietary) administration of the corticosteroid agonists mometasone furoate, budesonide, and flunisolide. *Toxicol. Pathol.* **39**(6), 958-968 (2011).
- O'Callaghan, C., White, J.A., and Kantar, A. Nebulization of corticosteroids to asthmatic children: Large variation in dose inhaled. *Respirology* **19**(2), 276-279 (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, 11/14/2022

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