

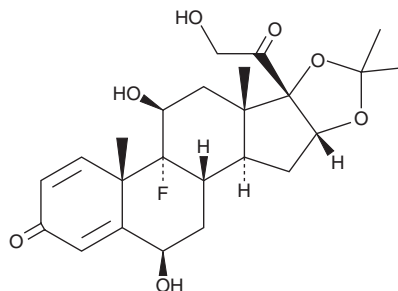
PRODUCT INFORMATION



6 β -hydroxy Triamcinolone acetonide

Item No. 40122

CAS Registry No.: 3869-32-7
Formal Name: 9-fluoro-6 β ,11 β ,21-trihydroxy-16 α ,17-
[[1-methylethylidene]bis(oxy)]-pregna-
1,4-diene-3,20-dione
Synonym: 6 β -hydroxy TrA
MF: C₂₄H₃₁FO₇
FW: 450.5
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

6 β -hydroxy Triamcinolone acetonide is supplied as a solid. A stock solution may be made by dissolving the 6 β -hydroxy triamcinolone acetonide in the solvent of choice, which should be purged with an inert gas. 6 β -hydroxy Triamcinolone acetonide is slightly soluble in the organic solvent DMSO.

Description

6 β -hydroxy Triamcinolone acetonide is a metabolite of the synthetic corticosteroid triamcinolone acetonide (Item No. 18026).¹ It is formed from triamcinolone acetonide by the cytochrome P450 (CYP) isoforms CYP3A4, CYP3A5, and CYP3A7. 6 β -hydroxy Triamcinolone acetonide has been found in wastewater effluent.²

References

1. Moore, C.D., Roberts, J.K., Orton, C.R., *et al.* Metabolic pathways of inhaled glucocorticoids by the CYP3A enzymes. *Drug Metab. Dispos.* **41(2)**, 379-389 (2013).
2. Weizel, A., Schlüsener, M.P., Dierkes, G., *et al.* Occurrence of glucocorticoids, mineralocorticoids, and progestogens in various treated wastewater, rivers, and streams. *Environ. Sci. Technol.* **52(9)**, 5296-5307 (2018).

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, 12/12/2023

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