

PRODUCT INFORMATION



Δ^{14} -Triamcinolone acetonide

Item No. 9000087

CAS Registry No.: 1260149-96-9
Formal Name: (6aS,6bR,7S,8aS,8bS,11aR,12bS)-6b-fluoro-7-hydroxy-8b-(2-hydroxyacetyl)-6a,8a,10,10-tetramethyl-1,2,6a,6b,7,8,8a,8b,11a,12b-decahydro-4H-naphtho[2',1':4,5]indeno[1,2-d][1,3]dioxol-4-one

Synonyms: 14,15-dehydro Triamcinolone acetonide, Triamcinolone acetonide Impurity B

MF: C₂₄H₂₉FO₆

FW: 432.5

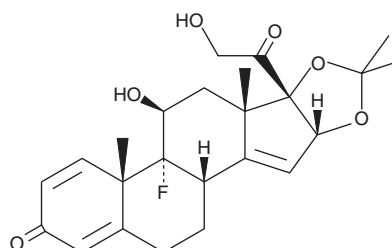
Purity: ≥96%

UV/Vis.: λ_{max}: 238 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

Δ^{14} -Triamcinolone acetonide is a potential impurity found in commercial preparations of triamcinolone acetonide.¹ Triamcinolone acetonide is a synthetic corticosteroid.² It decreases cytokine levels, the firing rate of sensory neurons, and mechanical hypersensitivity in a rat spinal nerve ligation model when used at a dose of 1.5 mg/kg prior to and following surgery for three days. Triamcinolone acetonide also decreases outflow facility in a mouse model of steroid-induced glaucoma when 20 μ l of a 40 mg/ml suspension is administered subconjunctivally.³ Formulations containing triamcinolone acetonide are used in the treatment of diabetic macular edema.

References

1. Cavina, G., Alimenti, R., Gallinella, B., *et al.* The identification of related substances in triamcinolone acetonide by means of high-performance liquid chromatography with diode array detector and mass spectrometry. *J. Pharm. Biomed. Anal.* **10(9)**, 685-692 (1992).
2. Li, H., Xie, W., Strong, J.A., *et al.* Systemic antiinflammatory corticosteroid reduces mechanical pain behavior, sympathetic sprouting, and elevation of proinflammatory cytokines in a rat model of neuropathic pain. *Anesthesiology* **107(3)**, 469-477 (2007).
3. Kumar, S., Shah, S., Deutsch, E.R., *et al.* Triamcinolone acetonide decreases outflow facility in C57BL/6 mouse eyes. *Invest. Ophthalmol. Vis. Sci.* **54(2)**, 1280-1287 (2013).

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, 05/31/2018

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