

PRODUCT INFORMATION



Triamcinolone acetonide-d₇ Item No. 25236

Formal Name: (6aS,6bR,7S,8aS,8bS,11aR,12aS,12bS)-6b-fluoro-7-hydroxy-8b-(2-hydroxyacetyl-2-d)-6a,8a-dimethyl-10,10-bis(methyl-d₃)-1,2,6a,6b,7,8,8a,8b,11a,12,12a,12b-dodecahydro-4H-naphtho[2',1':4,5]indeno[1,2-d][1,3]dioxol-4-one

MF: C₂₄H₂₄D₇FO₆

FW: 441.5

Chemical Purity: ≥95% (Triamcinolone acetonide)

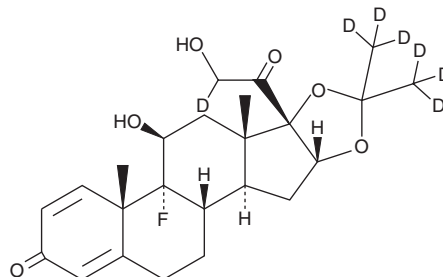
Deuterium

Incorporation: ≥99% deuterated forms (d₁-d₇); ≤1% d₀

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

Triamcinolone acetonide-d₇ is intended for use as an internal standard for the quantification of triamcinolone acetonide (Item No. 18026) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Triamcinolone acetonide-d₇ is supplied as a solid. A stock solution may be made by dissolving the triamcinolone acetonide-d₇ in the solvent of choice, which should be purged with an inert gas. Triamcinolone acetonide-d₇ is slightly soluble in chloroform and methanol.

Description

Triamcinolone acetonide is a synthetic corticosteroid.¹ 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 µl of a 40 mg/ml suspension is administered subconjunctivally.² Formulations containing triamcinolone acetonide are used in the treatment of diabetic macular edema.

References

1. 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).
2. 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, 12/06/2022

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