

PRODUCT INFORMATION



6 β -hydroxy Budesonide

Item No. 34979

Formal Name: (2R,6aR,6bS,7S,8aS,8bS,11aR,12aS,12bS)-2,7-dihydroxy-8b-(2-hydroxyacetyl)-6a,8a-dimethyl-10-propyl-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: C₂₅H₃₄O₇

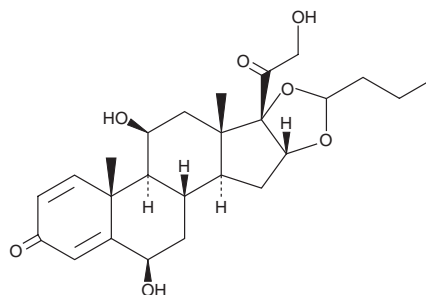
FW: 446.5

Purity: \geq 95%

Supplied as: A 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

6 β -hydroxy Budesonide is supplied as a solid. A stock solution may be made by dissolving the 6 β -hydroxy budesonide in the solvent of choice, which should be purged with an inert gas. 6 β -hydroxy Budesonide is soluble in organic solvents such as methanol, DMSO, acetonitrile, and dimethyl formamide.

Description

6 β -hydroxy Budesonide is a metabolite of the glucocorticoid budesonide (Item No. 15407).¹

Reference

1. Matabosch, X., Pozo, O.J., Pérez-Mañá, C., *et al.* Identification of budesonide metabolites in human urine after oral administration. *Anal. Bioanal. Chem.* **404**(2), 325-340 (2012).

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, 12/02/2022

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