

PRODUCT INFORMATION

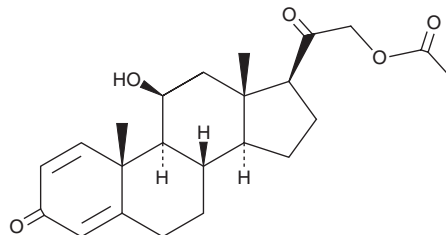


1-Dehydrocorticosterone 21-acetate

Item No. 36298

CAS Registry No.: 58652-04-3
Formal Name: 21-(acetyloxy)-11 β -hydroxy-pregna-1,4-diene-3,20-dione
Synonyms: 1-DHC 21-acetate,
11 β ,21-Dihydroxypregna-1,4-diene-3,20-dione Acetate

MF: C₂₃H₃₀O₅
FW: 386.5
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 244 nm
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

1-Dehydrocorticosterone 21-acetate is supplied as a solid. A stock solution may be made by dissolving the 1-dehydrocorticosterone 21-acetate in the solvent of choice, which should be purged with an inert gas. 1-Dehydrocorticosterone 21-acetate is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 1-dehydrocorticosterone 21-acetate in these solvents is approximately 2, 12, and 16 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of 1-dehydrocorticosterone 21-acetate can be prepared by directly dissolving the solid in aqueous buffers. The solubility of 1-dehydrocorticosterone 21-acetate in PBS (pH 7.2) is approximately 0.30 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

1-Dehydrocorticosterone 21-acetate is an acetylated derivative of corticosterone (Item No. 16063) that is dehydrogenated at the one position. It has been used as a precursor in the synthesis of 21-hydroxypregna-1,4-diene-3,11,20-trione, which is a metabolite of 1 α -hydroxycorticosterone in the skate.¹

Reference

1. Truscott, B., Kane, K.M., and Idler, D.R. 21-Hydroxypregna-1,4-diene-3,11,20-trione: A biliary metabolite of a cartilaginous fish, *Raja* sp. *Steroids* **31**(4), 573-582 (1978).

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, 12/19/2022

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