

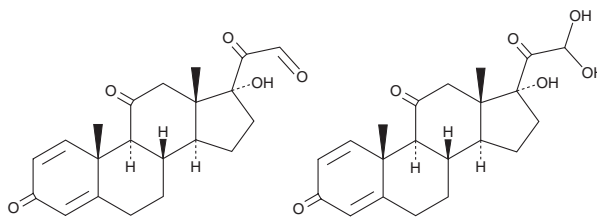
PRODUCT INFORMATION



Prednisone 21-aldehyde/22-hydroxy Prednisone

Item No. 22570

Formal Name: 17-hydroxy-3,11,20-trioxo-pregna-1,4-dien-21-al; 17,21,21-trihydroxy-pregna-1,4-diene-3,11,20-trione
MF: $C_{21}H_{24}O_5 \bullet C_{21}H_{26}O_6$
FW: 730.9
Purity: $\geq 85\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

Prednisone 21-aldehyde/22-hydroxy Prednisone is supplied as a solid. A stock solution may be made by dissolving the prednisone 21-aldehyde/22-hydroxy prednisone in the solvent of choice, which should be purged with an inert gas. Prednisone 21-aldehyde/22-hydroxy Prednisone is slightly soluble in chloroform, DMSO, and methanol.

Description

Prednisone 21-aldehyde is a derivative of the synthetic corticosteroid prednisone (Item No. 20677) that has antiviral activity against influenza virus A-PR8 in embryonated chicken eggs.¹ 22-hydroxy Prednisone is an unstable adduct of prednisone.² It has been observed following incubation of prednisone with nornicotine. This product is a mixture of prednisone 21-aldehyde and 22-hydroxy prednisone.

Reference

1. Cavallini, G., Massarani, E., and Nardi, D. Antiviral Compounds. IX. Steroid Derivatives. *J. Med. Chem.* **7(5)**, 673-675 (1964).
2. Dickerson, T.J., and Janda, K.D. A previously undescribed chemical link between smoking and metabolic disease. *Proc. Natl. Acad. Sci. U.S.A.* **99(23)**, 15084-15088 (2002).

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, 10/21/2022

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