

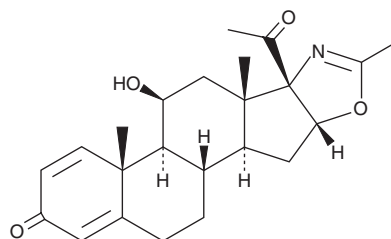
PRODUCT INFORMATION



21-deacetoxy Deflazacort

Item No. 43019

CAS Registry No.: 13649-88-2
Formal Name: 11 β -hydroxy-2'-methyl-5'H-pregna-1,4-dieno[17,16 β -d]oxazole-3,20-dione
MF: C₂₃H₂₉NO₄
FW: 383.5
Purity: $\geq 95\%$
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

21-deacetoxy Deflazacort is supplied as a solid. A stock solution may be made by dissolving the 21-deacetoxy deflazacort in the solvent of choice, which should be purged with an inert gas. 21-deacetoxy Deflazacort is slightly soluble (0.1-1 mg/ml) in ethanol and sparingly soluble (1-10 mg/ml) in DMSO.

Description

21-deacetoxy Deflazacort is a glucocorticoid.¹ It induces hepatic gluconeogenesis in adrenalectomized rats when administered at a dose of 0.2 mg/kg.

Reference

1. Nathansohn, G., Pasqualucci, C.R., Radaelli, P., *et al.* Steroids possessing nitrogen atoms. V. (1) Some pharmacological and chemico-physical aspects of a new class of active corticoids: [17 α ,16 α -d]-oxazolino-steroids. *Steroids* **13**(3), 365-381 (1969).

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, 03/31/2025

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