

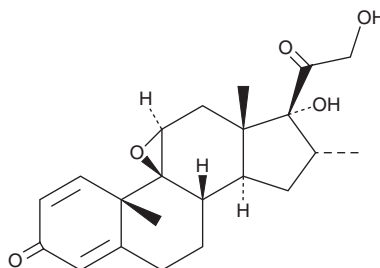
PRODUCT INFORMATION



Dexamethasone 9,11-epoxide

Item No. 36161

CAS Registry No.: 24916-90-3
Formal Name: 9 β ,11 β -epoxy-17,21-dihydroxy-16 α -methyl-pregna-1,4-diene-3,20-dione
MF: C₂₂H₂₈O₅
FW: 372.5
Purity: \geq 98%
UV/Vis.: λ_{max} : 249 nm
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

Dexamethasone 9,11-epoxide is supplied as a solid. A stock solution may be made by dissolving the dexamethasone 9,11-epoxide in the solvent of choice, which should be purged with an inert gas. Dexamethasone 9,11-epoxide is soluble in DMSO.

Description

Dexamethasone 9,11-epoxide is a potential impurity found in preparations of dexamethasone.¹ It is an autooxidative degradation product formed under alkaline conditions.

Reference

1. Li, M., Chen, B., Monteiro, S.G., *et al.* Mechanism of base-catalyzed autooxidation of corticosteroids containing 20-keto-21-hydroxyl side chain. *Tetrahedron Lett.* **50**(32), 4575-4581 (2009).

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, 11/04/2022

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