

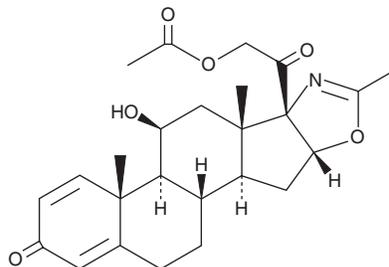
PRODUCT INFORMATION



Deflazacort

Item No. 20386

CAS Registry No.: 14484-47-0
Formal Name: 21-(acetyloxy)-11 β -hydroxy-2'-methyl-5'H-pregna-1,4-dieno[17,16 β -d]oxazole-3,20-dione
Synonyms: DL-458IT, L-5,458, MDL 458
MF: C₂₅H₃₁NO₆
FW: 441.5
Purity: \geq 98%
UV/Vis.: λ_{max} : 242 nm
Supplied as: A crystalline 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

Deflazacort is supplied as a crystalline solid. A stock solution may be made by dissolving the deflazacort in the solvent of choice, which should be purged with an inert gas. Deflazacort is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of deflazacort in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Deflazacort is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, deflazacort should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Deflazacort has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Deflazacort is a prodrug for the active glucocorticoid 21-desacetyl deflazacort (Item No. 20316). Deflazacort and related glucocorticoids have anti-inflammatory and immunosuppressant effects, as well as benefits in certain muscular dystrophies.¹⁻³

References

1. Canonica, G.W., Ciprandi, G., Passalacqua, G., *et al.* Molecular events in allergic inflammation: Experimental models and possible modulation. *Allergy* **52(Suppl. 34)**, 25-30 (1997).
2. Czaja, A.J. Promising pharmacological, molecular and cellular treatments of autoimmune hepatitis. *Curr. Pharm. Des.* **17(29)**, 3120-3140 (2011).
3. Matthews, E., Brassington, R., Kuntzer, T., *et al.* Corticosteroids for the treatment of Duchenne muscular dystrophy. *Cochrane Database Syst. Rev.* **5**, CD003725 (2016).

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