

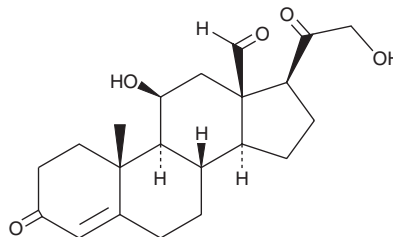
PRODUCT INFORMATION



Aldosterone

Item No. 15273

CAS Registry No.: 52-39-1
Formal Name: 11 β ,21-dihydroxy-3,20-dioxo-pregn-4-en-18-al
Synonym: 18-Aldocorticosterone, NSC 73856, Reichstein X
MF: C₂₁H₂₈O₅
FW: 360.4
Purity: \geq 98%
UV/Vis.: λ_{max} : 240 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

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

Aldosterone is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, aldosterone should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Aldosterone has a solubility of approximately 0.2 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Aldosterone is a steroid hormone secreted by the adrenal cortex and is the principle mineralocorticoid controlling sodium and potassium balance.^{1,2} Its primary role is to promote unidirectional salt reabsorption in epithelial tissues. Aldosterone is synthesized from cholesterol in the zona glomerulosa of the adrenal cortex and its secretion is regulated *via* the renin-angiotensin system.³

References

1. Agarwal, M.K. and Mirshahi, M. General overview of mineralocorticoid hormone action. *Pharmacol. Toxicol.* **84**, 273-326 (1999).
2. Rogerson, F.M. and Fuller, P.J. Mineralocorticoid action. *Steroids.* **65**, 61-73 (2000).
3. Lumbers, E.R. Angiotensin and aldosterone. *Regul. Pept.* **80**, 91-100 (1999).

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, 10/25/2022

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