

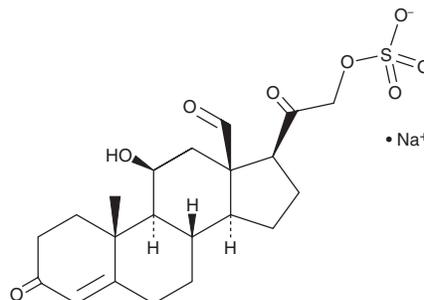
PRODUCT INFORMATION



Aldosterone 21-sulfate (sodium salt)

Item No. 9003450

Formal Name: (11 β)-11-hydroxy-3,20-dioxo-21-(sulfooxy)-pregn-4-en-18-al, monosodium salt
Synonym: Aldosterone 21-sulphate
MF: C₂₁H₂₇O₈S • Na
FW: 462.5
Purity: \geq 98%
UV/Vis.: λ_{max} : 240 mn
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 2 years



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

Laboratory Procedures

Aldosterone 21-sulfate (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the aldosterone 21-sulfate (sodium salt) in the solvent of choice, which should be purged with an inert gas. Aldosterone 21-sulfate (sodium salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of aldosterone 21-sulfate (sodium salt) in these solvents is approximately 1 mg/ml in ethanol and 15 mg/ml in DMSO and DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of aldosterone 21-sulfate (sodium salt) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of aldosterone 21-sulfate (sodium salt) in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Aldosterone 21-sulfate is a metabolite of the steroid hormone aldosterone (Item No. 15273).¹ It reduces the sodium-to-potassium ratio in rat urine when administered at doses ranging from 10 to 100 μ g, indicating mineralocorticoid activity that is less than 1% that of aldosterone.²

References

1. Grose, J.H., Nowaczynski, W., Kuchel, O., *et al.* Isolation of aldosterone urinary metabolites, glucuronides and sulfate. *J. Steroid Biochem.* **4(6)**, 551-566 (1973).
2. Jowett, T.P. and Slater, J.D.H. Aldosterone 21-sulphate in man. *J. Steroid. Biochem.* **18(4)**, 471-479 (2019).

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, 09/03/2024

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