

PRODUCT INFORMATION



Androst-5-enediol Sulfate (sodium salt)

Item No. 43450

Formal Name: androst-5-ene-3 β ,17 β -diol, hydrogen sulfate, sodium salt

Synonym: 5-Androstenediol Sulfate

MF: C₁₉H₂₉O₅S • Na

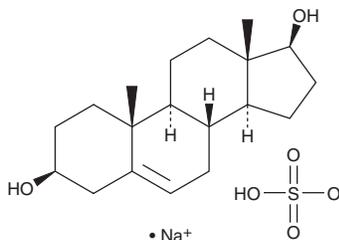
FW: 392.5

Purity: \geq 95% (mixture of isomers)

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

Androst-5-enediol sulfate (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the androst-5-enediol sulfate (sodium salt) in the solvent of choice, which should be purged with an inert gas. Androst-5-enediol sulfate (sodium salt) is soluble (\geq 10 mg/ml) in DMSO. Androst-5-enediol sulfate (sodium salt) is slightly soluble (0.1-1 mg/ml) in ethanol.

Description

Androst-5-enediol sulfate is a metabolite of 5-androstenediol.¹

Reference

1. Raps, M., Kleider, C., and Lehmann, L. Quantitative analysis of 34 sex (pro)hormones, conjugates and bioactive oxidation products thereof in human plasma by GC- and LC-MS/MS and systematic investigation of overestimations of analyte concentrations not accounted for by method validation. *Steroids* **208:109441**, (2024).

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, 08/18/2025

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