

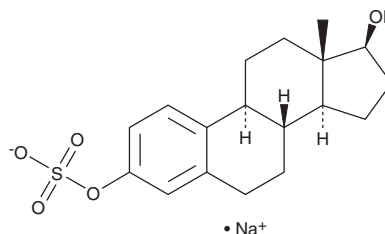
PRODUCT INFORMATION



17β-Estradiol 3-O-Sulfate (sodium salt)

Item No. 39169

CAS Registry No.: 4999-79-5
Formal Name: estra-1,3,5(10)-triene-3,17β-diol
3-(hydrogen sulfate), monosodium salt
Synonyms: Estradiol Sulfate, E2S
MF: C₁₈H₂₃O₅S • Na
FW: 374.4
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

17β-Estradiol 3-O-sulfate (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the 17β-estradiol 3-O-sulfate (sodium salt) in the solvent of choice, which should be purged with an inert gas. 17β-Estradiol 3-O-sulfate (sodium salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 17β-estradiol 3-O-sulfate (sodium salt) in these solvents is approximately 10 mg/ml.

Description

17β-Estradiol 3-O-sulfate is an inactive metabolite of the major premenopausal estrogen 17β-estradiol (Item No. 10006315).¹ 17β-Estradiol 3-O-sulfate is formed from 17β-estradiol via the sulfotransferase (SULT) isoforms SULT1A1 and SULT1A2.

Reference

1. Harris, R.M., Waring, R.H., Kirk, C.J., et al. Sulfation of "estrogenic" alkylphenols and 17β-estradiol by human platelet phenol sulfotransferases. *J. Biol. Chem.* **275**(1), 159-166 (2000).

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, 10/09/2023

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