

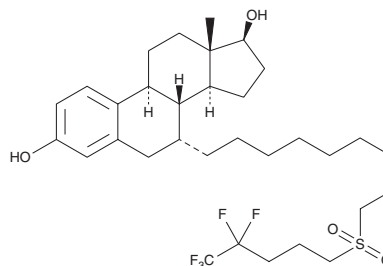
PRODUCT INFORMATION



Fulvestrant 9-sulfone

Item No. 35070

CAS Registry No.: 98008-06-1
Formal Name: 7 α -[9-[(4,4,5,5,5-pentafluoropentyl)sulfonyl]nonyl]-estra-1,3,5(10)-triene-3,17 β -diol
MF: C₃₂H₄₇F₅O₄S
FW: 622.8
Purity: \geq 98%
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

Fulvestrant 9-sulfone is supplied as a solid. A stock solution may be made by dissolving the fulvestrant 9-sulfone in the solvent of choice, which should be purged with an inert gas. Fulvestrant 9-sulfone is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of fulvestrant 9-sulfone in DMF is approximately 30 mg/ml and approximately 10 mg/ml in ethanol and DMSO.

Description

Fulvestrant 9-sulfone is an estrogen receptor antagonist and a derivative of the selective estrogen receptor degrader (SERD) fulvestrant (Item No. 10011269).¹ It binds to estrogen receptor α (ER α) and ER β and inhibits estrogen-induced gene expression in a reporter assay in Ishikawa endometrial cells expressing an estrogen response element (ERE) reporter when used at concentrations ranging from 0.1 to 100 μ M.

Reference

1. Nuttall, M.E., Pendrak, I., Emery, J.G., *et al.* Antagonism of oestrogen action in human breast and endometrial cells in vitro: Potential novel antitumour agents. *Cancer Chemother. Pharmacol.* **47(5)**, 437-443 (2001).

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, 12/09/2022

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