

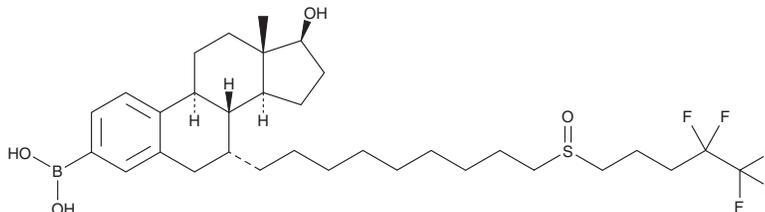
PRODUCT INFORMATION



ZB716

Item No. 42238

CAS Registry No.: 1853279-29-4
Formal Name: B-[(7 α ,17 β)-17-hydroxy-7-[9-[(4,4,5,5,5-pentafluoropentyl)sulfinyl]nonyl]estra-1,3,5(10)-trien-3-yl]-boronic acid
Synonym: Fulvestrant-3 Boronic Acid
MF: C₃₂H₄₈BF₅O₄S
FW: 634.6
Purity: \geq 95%
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

ZB716 is supplied as a solid. A stock solution may be made by dissolving the ZB716 in the solvent of choice, which should be purged with an inert gas. ZB716 is soluble in the organic solvent DMSO.

Description

ZB716 is a degrader of estrogen receptor α (ER α ; IC₅₀ = 4.1 nM).¹ It inhibits estradiol-induced ER α transcriptional activity in a reporter assay using T47D cells when used at concentrations of 1 or 10 nM. ZB716 decreases the proliferation of MCF-7 and T47D breast cancer cells and tamoxifen-resistant MCF-7 and T47D cells (IC₅₀s = 0.0032, 0.0062, 0.069, and 0.037 nM, respectively).

Reference

1. Liu, J., Zheng, S., Akerstrom, V.L., et al. Fulvestrant-3 boronic acid (ZB716): An orally bioavailable selective estrogen receptor downregulator (SERD). *J. Med. Chem.* **59**(17), 8134-8140 (2016).

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, 10/15/2024

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