

PRODUCT INFORMATION

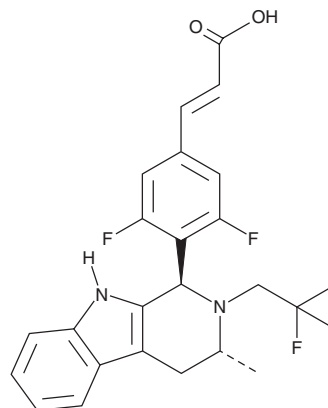


AZD 9496

Item No. 21805

CAS Registry No.: 1639042-08-2
Formal Name: (2E)-3-[3,5-difluoro-4-[(1R,3R)-2-(2-fluoro-2-methylpropyl)-2,3,4,9-tetrahydro-3-methyl-1H-pyrido[3,4-b]indol-1-yl]phenyl]-2-propenoic acid

MF: C₂₅H₂₅F₃N₂O₂
FW: 442.5
Purity: ≥98%
UV/Vis.: λ_{max}: 225, 276 nm
Supplied as: A crystalline 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

AZD 9496 is supplied as a crystalline solid. A stock solution may be made by dissolving the AZD 9496 in the solvent of choice, which should be purged with an inert gas. AZD 9496 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of AZD 9496 in these solvents is approximately 30 mg/ml.

AZD 9496 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AZD 9496 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. AZD 9496 has a solubility of approximately 0.14 mg/ml in a 1:6 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

AZD 9496 is a potent and selective estrogen receptor downregulator (SERD) with an IC₅₀ value of 0.138 nM for estrogen receptor α (ERα) downregulation.¹ It is selective for ERα compared to other nuclear hormone receptors with IC₅₀ values of 0.0008, 0.54, 9.2, and 30 μM for ERα, progesterone receptor (PR), glucocorticoid receptor (GR), and androgen receptor (AR), respectively. AZD 9496 decreases ERα activity (IC₅₀ = 0.282 nM), measured *via* quantification of downstream PR activity, and reduces proliferation of MCF-7 breast cancer cells (IC₅₀ = 0.0398 nM). It also inhibits MCF-7 xenograft growth in mice in a dose-dependent manner.

Reference

1. De Savi, C., Bradbury, R.H., Rabow, A.A., *et al.* Optimization of a novel binding motif to (E)-3-(3,5-difluoro-4-((1R,3R)-2-(2-fluoro-2-methylpropyl)-3-methyl-2,3,4,9-tetrahydro-1H-pyrido[3,4-b]indol-1-yl)phenyl) acrylic acid (AZD9496), a potent and orally bioavailable selective estrogen receptor downregulator and antagonist. *J. Med. Chem.* **58**(20), 8128-8140 (2015).

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, 11/14/2022

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