

PRODUCT INFORMATION



Estrogen Receptor Antagonist 2

Item No. 43587

CAS Registry No.: 2222844-89-3

Formal Name: N-[1-(3-fluoropropyl)-3-azetidiny]-6-
[(6S,8R)-6,7,8,9-tetrahydro-8-methyl-7-
(2,2,2-trifluoroethyl)-3H-pyrazolo[4,3-f]
isoquinolin-6-yl]-3-pyridinamine

Synonym: AZD9833

MF: $C_{24}H_{28}F_4N_6$

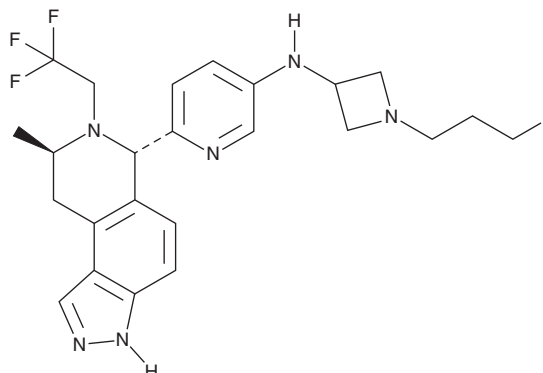
FW: 476.5

Purity: $\geq 95\%$

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

Estrogen receptor antagonist 2 is supplied as a solid. A stock solution may be made by dissolving the estrogen receptor antagonist 2 in the solvent of choice, which should be purged with an inert gas. Estrogen receptor antagonist 2 is soluble (≥ 10 mg/ml) in ethanol and DMSO.

Description

Estrogen receptor antagonist 2 is a degrader of estrogen receptor α (ER α).¹ It induces degradation of ER α in MCF-7 breast cancer cells ($IC_{50} = 0.16$ nM). Estrogen receptor antagonist 2 inhibits estradiol-induced growth in MCF-7 cells ($IC_{50} = 5.01$ nM). *In vivo*, estrogen receptor antagonist 2 (0.5-50 mg/kg per day) decreases tumor volume in an MCF-7 mouse xenograft model.

Reference

1. Scott, J.S., Moss, T.A., Balazs, A., *et al.* Discovery of AZD9833, a potent and orally bioavailable selective estrogen receptor degrader and antagonist. *J. Med. Chem.* **63**(23), 14530-14559 (2020).

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, 10/29/2025

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