

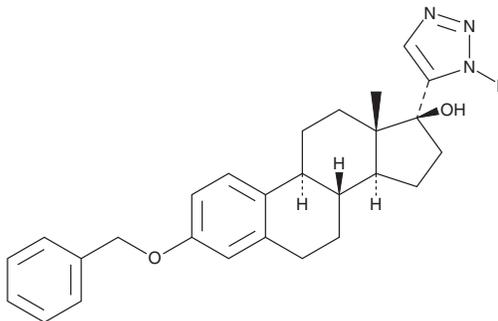
PRODUCT INFORMATION



Fz25

Item No. 37752

CAS Registry No.: 2762498-70-2
Formal Name: (17 β)-3-(phenylmethoxy)-17-(1H-1,2,3-triazol-5-yl)-estra-1,3,5(10)-trien-17-ol
MF: C₂₇H₃₁N₃O₂
FW: 429.6
Purity: \geq 98%
UV/Vis.: λ_{max} : 280 nm
Supplied as: A solid
Storage: -20°C
Stability: \geq 2 years



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

Laboratory Procedures

Fz25 is supplied as a solid. A stock solution may be made by dissolving the Fz25 in the solvent of choice, which should be purged with an inert gas. Fz25 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of Fz25 in DMSO and DMF is approximately 25 mg/ml and approximately 5 mg/ml in ethanol.

Description

Fz25 is a triazole-modified sterol and an anticancer agent.¹ It inhibits the proliferation of A2780 ovarian and HT-29 colorectal cancer cells (IC₅₀s = 15.29 and 15.98 μ M, respectively). Fz25 decreases the phosphorylation of EGFR and STAT3 in A2780 cells when used at a concentration of 30.58 μ M and phosphorylation of ERK, STAT3, and mammalian target of rapamycin (mTOR) in HT-29 cells at 31.96 μ M. It reduces the efflux activity of P-glycoprotein (P-gp), also known as multidrug resistance protein 1 (MDR1), in a reporter assay using HEK293 cells when used at a concentration of 10 μ M.

Reference

1. Ostlund, T., Alotaibi, F., Kyeremateng, J., et al. Triazole-estradiol analogs: A potential cancer therapeutic targeting ovarian and colorectal cancer. *Steroids* **177**, 108950 (2022).

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, 04/20/2023

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