

PRODUCT INFORMATION

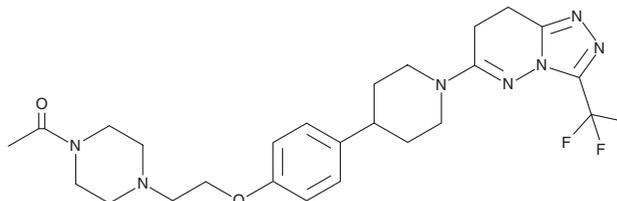


AZD 3514

Item No. 33222

CAS Registry No.: 1240299-33-5
Formal Name: 1-[4-[2-[4-[1-[7,8-dihydro-3-(trifluoromethyl)-1,2,4-triazolo[4,3-b]pyridazin-6-yl]-4-piperidinyl]phenoxy]ethyl]-1-piperazinyl]-ethanone

MF: C₂₅H₃₂F₃N₇O₂
FW: 519.6
Purity: ≥98%
UV/Vis.: λ_{max}: 222, 272 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 3514 is supplied as a crystalline solid. A stock solution may be made by dissolving the AZD 3514 in the solvent of choice, which should be purged with an inert gas. AZD 3514 is soluble in the organic solvent DMSO.

Description

AZD 3514 is a selective androgen receptor downregulator (SARD).¹ It inhibits AR nuclear translocation in LNCaP human prostate cancer cells with an IC₅₀ value of 1.78 μM. AZD 3514 (10 μM) inhibits dihydrotestosterone-induced proliferation of LNCaP and LAPC4 prostate cancer cells.² It reduces testosterone propionate-induced increases in seminal vesicle weight in castrated rats. AZD 3514 (50 mg/kg) decreases tumor growth of Dunning R3327H prostate tumor implants in rats.

References

1. Bradbury, R.H., Acton, D.G., Broadbent, N.L., *et al.* Discovery of AZD3514, a small-molecule androgen receptor downregulator for treatment of advanced prostate cancer. *Bioorg. Med. Chem. Lett.* **23(7)**, 1945-1948 (2013).
2. Loddick, S.A., Ross, S.J., Thomason, A.G., *et al.* AZD3514: A small molecule that modulates androgen receptor signaling and function *in vitro* and *in vivo*. *Mol. Cancer Ther.* **12(9)**, 1715-1727 (2013).

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, 10/17/2022

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