

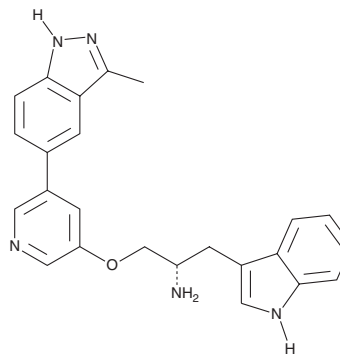
PRODUCT INFORMATION



A-443654

Item No. 16499

CAS Registry No.: 552325-16-3
Formal Name: αS-[[[5-(3-methyl-1H-indazol-5-yl)-3-pyridinyl]oxy)methyl]-1H-indole-3-ethanamine
MF: C₂₄H₂₃N₅O
FW: 397.5
Purity: ≥98%
UV/Vis.: λ_{max}: 222, 291 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

A-443654 is supplied as a crystalline solid. A stock solution may be made by dissolving the A-443654 in the solvent of choice, which should be purged with an inert gas. A-443654 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of A-443654 in these solvents is approximately 10, 15, and 20 mg/ml, respectively.

A-443654 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, A-443654 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. A-443654 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Three related forms of the kinase Akt (1, 2, 3, also known as protein kinase B isoforms PKBα, β, γ) modulate cell proliferation, metabolism, and survival as well as angiogenesis. A-443654 is an inhibitor of Akt (K_i = 160 pmol/L for all three isoforms) that interferes with mitotic progression and bipolar spindle formation.^{1,2} It induces G₂/M accumulation, defects in centrosome separation, and formation of either monopolar arrays or disorganized spindles.² A-443654 has been reported to slow the progression of Akt-dependent tumors in *in vivo* mouse models.¹ In response to an A-443654-induced decrease in phosphorylation of Akt targets, a concomitant increase in Thr³⁰⁸ and Ser⁴⁷³ phosphorylation of Akt has been observed in human cancer cell lines.³ A-443654 has been used to examine the mechanism of this rapid feedback reaction.

References

1. Luo, Y., Shoemaker, A.R., Liu, X., *et al.* Potent and selective inhibitors of Akt kinases slow the progress of tumors *in vivo*. *Mol. Cancer Ther.* **4(6)**, 977-968 (2005).
2. Liu, X., Shi, Y., Woods, K.W., *et al.* Akt inhibitor a-443654 interferes with mitotic progression by regulating aurora a kinase expression. *Neoplasia* **10(8)**, 828-837 (2008).
3. Han, E.K., Levenson, J.D., McGonigal, T., *et al.* Akt inhibitor A-443654 induces rapid Akt Ser-473 phosphorylation independent of mTORC1 inhibition. *Oncogene* **26(38)**, 5655-5661 (2007).

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