

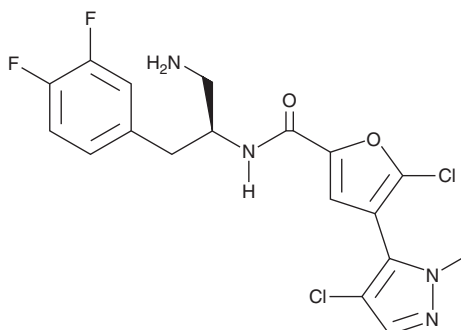
PRODUCT INFORMATION



Uprosertib

Item No. 19904

CAS Registry No.: 1047634-65-0
Formal Name: N-[(1S)-2-amino-1-[(3,4-difluorophenyl)methyl]ethyl]-5-chloro-4-(4-chloro-1-methyl-1H-pyrazol-5-yl)-2-furancarboxamide
Synonyms: GSK795, GSK2141795
MF: C₁₈H₁₆Cl₂F₂N₄O₂
FW: 429.2
Purity: ≥98%
UV/Vis.: λ_{max}: 239 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

Uprosertib is supplied as a crystalline solid. A stock solution may be made by dissolving the uprosertib in the solvent of choice, which should be purged with an inert gas. Uprosertib is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of uprosertib in these solvents is approximately 1 mg/ml.

Description

Uprosertib is a selective, orally bioavailable inhibitor of Akt (IC₅₀s = 180, 328, and 38 nM for Akt1, Akt2, and Akt3, respectively).¹ Uprosertib preferentially inhibits the proliferation of human cancer cell lines with Akt pathway activation via PI3K/PTEN mutation or loss.² It can cause cell cycle arrest and consequent tumor growth inhibition in mice bearing either BT474 breast tumor or SKOV3 ovarian tumor xenografts.²

References

1. Pachl, F., Plattner, P., Ruprecht, B., *et al.* Characterization of a chemical affinity probe targeting Akt kinases *J. Proteome Res.* **12**(8), 3792-3800 (2013).
2. Dumble, M., Crouthamel, M.-C., Zhang, S.-Y., *et al.* Discovery of novel AKT inhibitors with enhanced anti-tumor effects in combination with the MEK inhibitor. *PLoS One* **9**(6), (2014).

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, 03/09/2023

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