

PRODUCT INFORMATION

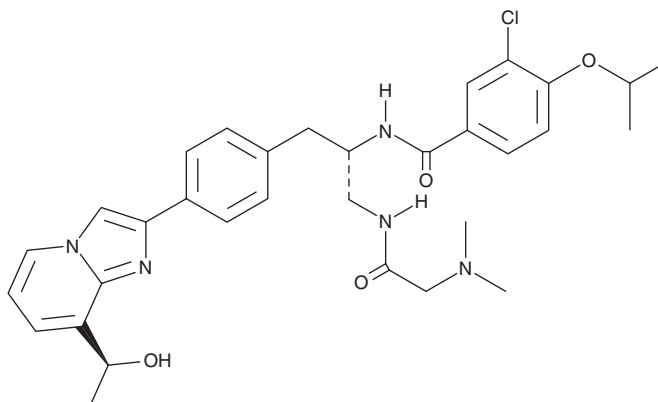


GSK923295

Item No. 18389

CAS Registry No.: 1088965-37-0
Formal Name: 3-chloro-N-[(1S)-2-[[2-(dimethylamino)acetyl]amino]-1-[[4-[8-[(1S)-1-hydroxyethyl]imidazo[1,2-a]pyridin-2-yl]phenyl]methyl]ethyl]-4-(1-methylethoxy)-benzamide

Synonym: GSK923295A
MF: C₃₂H₃₈ClN₅O₄
FW: 592.1
Purity: ≥98%
UV/Vis.: λ_{max}: 209, 251, 317 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

GSK923295 is supplied as a crystalline solid. A stock solution may be made by dissolving the GSK923295 in the solvent of choice. GSK923295 is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of GSK923295 in these solvents is approximately 30 mg/ml.

GSK923295 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, GSK923295 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. GSK923295 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

GSK923295 is a potent inhibitor of centromere-associated protein E (CENP-E; K_i = 3.2 nM), a kinesin motor protein involved in mitotic checkpoint signaling.^{1,2} GSK923295 causes mitotic cell cycle delay, leading to apoptosis in various cancer cell lines.¹ It induces apoptosis in cancer cells in mice bearing xenografts of COLO 205 cells.^{1,2} GSK923295 can be used with AZ3146, an inhibitor of the spindle checkpoint kinase Mps1, to generate aneuploidy in cells.³

References

1. Wood, K. W., Lad, L., Luo, L., *et al.* Antitumor activity of an allosteric inhibitor of centromere-associated protein-E. *PNAS* **107**(13), 5839-5844 (2010).
2. Qian, X., McDonald, A., Zhou, H.-J., *et al.* Discovery of the first potent and selective inhibitor of centromere-associated protein E: GSK923295. *ACS Med. Chem. Lett.* **1**, 30-34 (2010).
3. Bennett, A., Bechi, B., Tighe, A., *et al.* Cenp-E inhibitor GSK923295: Novel synthetic route and use as a tool to generate aneuploidy. *Oncotarget* **6**(25), 20921-20932 (2015).

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, 11/21/2022

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