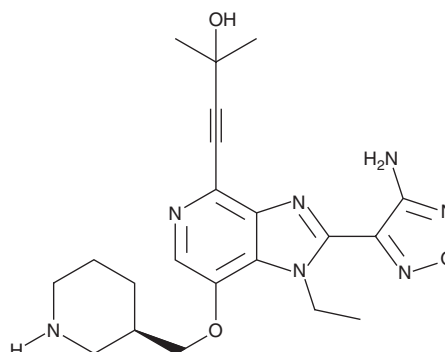


PRODUCT INFORMATION



GSK690693
Item No. 16891

CAS Registry No.: 937174-76-0
Formal Name: 4-[2-(4-amino-1,2,5-oxadiazol-3-yl)-1-ethyl-7-[(3S)-3-piperidinylmethoxy]-1H-imidazo[4,5-c]pyridin-4-yl]-2-methyl-3-Butyn-2-ol
MF: C₂₁H₂₇N₇O₃
FW: 425.5
Purity: ≥95%
UV/Vis.: λ_{max}: 215, 250, 315 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

GSK690693 is supplied as a crystalline solid. A stock solution may be made by dissolving the GSK690693 in the solvent of choice, which should be purged with an inert gas. GSK690693 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of GSK690693 in these solvents is approximately 25 mg/ml.

GSK690693 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, GSK690693 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. GSK690693 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

GSK690693 is an ATP-competitive, pan-Akt kinase inhibitor with IC₅₀ values of 2, 13, and 9 nM for Akt1, 2, and 3, respectively.¹ It also inhibits other members of the AGC kinase family, including AMPK, DAPK3, PAKs, PKA, PKC, PKG, and PrkX (IC₅₀ values range from 2-81 nM).¹ By preventing the phosphorylation of downstream Akt substrates, GSK690693 has been shown to inhibit proliferation and to induce apoptosis in various human tumor cells *in vitro* and in xenografts in immunocompromised mice.¹

Reference

1. Rhodes, N., Heerding, D.A., Duckett, D.R., *et al.* Characterization of an Akt kinase inhibitor with potent pharmacodynamic and antitumor activity. *Cancer Res.* **68**(7), 2366-2374 (2008).

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, 11/10/2022

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