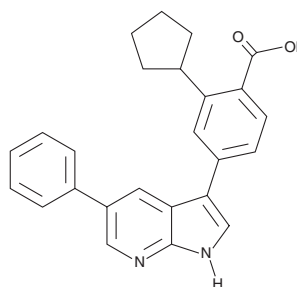


PRODUCT INFORMATION



GSK650394
Item No. 17001

CAS Registry No.: 890842-28-1
Formal Name: 2-cyclopentyl-4-(5-phenyl-1H-pyrrolo[2,3-b]pyridin-3-yl)-benzoic acid
MF: C₂₅H₂₂N₂O₂
FW: 382.5
Purity: ≥98%
UV/Vis.: λ_{max}: 241, 257, 295 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

GSK650394 is supplied as a crystalline solid. A stock solution may be made by dissolving the GSK650394 in the solvent of choice, which should be purged with an inert gas. GSK650394 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of GSK650394 in these solvents is approximately 30 mg/ml, respectively.

GSK650394 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, GSK650394 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. GSK650394 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

GSK650394 is an inhibitor of serum- and glucocorticoid-regulated kinase (IC₅₀s = 62 and 103 nM for SGK1 and SGK2, respectively).¹ It demonstrates greater than 30-fold selectivity for SGK1 over Akt and other related kinases.¹ GSK650394 can inhibit the androgen-stimulated growth of human prostate carcinoma LNCaP cells with an IC₅₀ value of ~ 1 μM.¹

Reference

1. Sherk, A.B., Frigo, D.E., Schnackenberg, C.G., *et al.* Development of a small-molecule serum- and glucocorticoid-regulated kinase-1 antagonist and its evaluation as a prostate cancer therapeutic. *Cancer Res.* **68**(18), 7475-7483 (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, 11/14/2022

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