

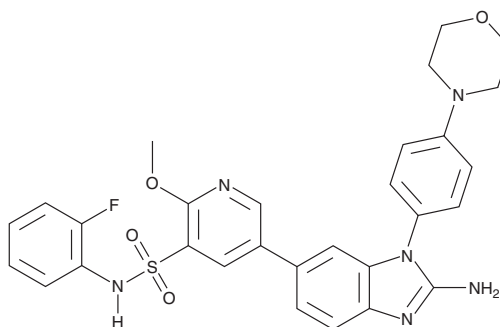
PRODUCT INFORMATION



GSK-A1

Item No. 34502

CAS Registry No.: 1416334-69-4
Formal Name: 5-[2-amino-1-[4-(4-morpholinyl)phenyl]-1H-benzimidazol-6-yl]-N-(2-fluorophenyl)-2-methoxy-3-pyridinesulfonamide
MF: C₂₉H₂₇FN₆O₄S
FW: 574.6
Purity: ≥98%
Supplied as: A 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

GSK-A1 is supplied as a solid. A stock solution may be made by dissolving the GSK-A1 in the solvent of choice, which should be purged with an inert gas. GSK-A1 is soluble in the organic solvent DMSO at a concentration of approximately 2 mg/ml.

Description

GSK-A1 is an inhibitor of phosphatidylinositol 4-kinase III α (PI4KIII α ; IC₅₀ = 3.16 nM).¹ It is selective for PI4KIII α over PI4KIII β , PI3K α , PI3K β , and PI3K δ (IC₅₀s = >50 nM) but also inhibits PI3K γ (IC₅₀ = 15.8 nM). GSK-A1 decreases the levels of phosphatidylinositol 4-phosphate (PtdIns-(4)-P₁) but not PtdIns-(4,5)-P₂ in HEK293 cells expressing the angiotensin II type 1 (AT₁) receptor (IC₅₀ = 3 nM). GSK-A1 inhibits hepatitis C virus (HCV) genotype 1a and 1b replication.

Reference

1. Bojjireddy, N., Botyanszki, J., Hammond, G., *et al.* Pharmacological and genetic targeting of the PI4KA enzyme reveals its important role in maintaining plasma membrane phosphatidylinositol 4-phosphate and phosphatidylinositol 4,5-bisphosphate levels. *J. Biol. Chem.* **289**(9), 6120-6132 (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, 12/06/2022

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