

# PRODUCT INFORMATION



## GSK-F1

Item No. 41366

CAS Registry No.: 1402345-92-9

Formal Name: 5-[2-amino-3,4-dihydro-4-oxo-3-[2-(trifluoromethyl)phenyl]-6-quinazoliny]-N-(2,4-difluorophenyl)-2-methoxy-3-pyridinesulfonamide

MF:  $C_{27}H_{18}F_5N_5O_4S$

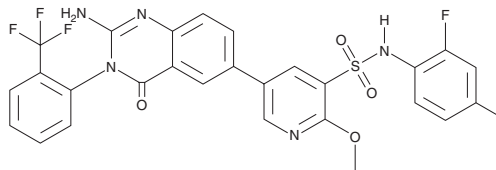
FW: 603.5

Purity:  $\geq 98\%$

Supplied as: A solid

Storage:  $-20^{\circ}\text{C}$

Stability:  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

GSK-F1 is supplied as a solid. A stock solution may be made by dissolving the GSK-F1 in the solvent of choice, which should be purged with an inert gas. GSK-F1 is soluble ( $\geq 10$  mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

### Description

GSK-F1 is a PI4KIII $\alpha$  inhibitor ( $IC_{50} = 5.01$  nM).<sup>1</sup> It is selective for PI4KIII $\alpha$  over PI4K $\beta$ , PI3K $\alpha$ , PI3K $\beta$ , PI3K $\delta$ , and PI4K $\gamma$  ( $IC_{50}$ s = 1,000, 2,512, 7,943, 2,512, and 2,512 nM, respectively). GSK-F1 inhibits the replication of hepatitis C virus (HCV) genotypes 1a, 1b, and 2a in a cell-based replicon assay ( $IC_{50}$ s = 12.6, 2.51, and 12.6 nM, respectively).

### Reference

- Leivers, A.L., Tallant, M., Shotwell, J.B., *et al.* Discovery of selective small molecule type III phosphatidylinositol 4-kinase  $\alpha$  (PI4KIII $\alpha$ ) inhibitors as anti hepatitis C (HCV) agents. *J. Med. Chem.* 57(5), 2091-2106 (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, 08/19/2024

#### 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