

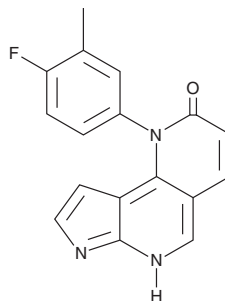
# PRODUCT INFORMATION



## STK16-IN-1

Item No. 27370

CAS Registry No.: 1223001-53-3  
Formal Name: 1-(4-fluoro-3-methylphenyl)-1,7-dihydro-2H-pyrrolo[2,3-h]-1,6-naphthyridin-2-one  
MF:  $C_{17}H_{12}FN_3O$   
FW: 293.3  
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

STK16-IN-1 is supplied as a solid. A stock solution may be made by dissolving the STK16-IN-1 in the solvent of choice, which should be purged with an inert gas. STK16-IN-1 is soluble in the organic solvent DMSO.

### Description

STK16-IN-1 is an inhibitor of the serine/threonine kinase STK16 ( $IC_{50} = 295 \text{ nM}$ ).<sup>1</sup> It is selective for STK16 over mTOR, PI3K $\delta$ , and PI3K $\gamma$  in an enzyme assay ( $IC_{50}$ s = 5,560, 856, and 867 nM, respectively) and over 440 additional kinases in a KinomeScan profiling assay. It decreases the growth of MCF-7 cells ( $GI_{50} = \sim 10 \text{ }\mu\text{M}$ ) and increases the number of binucleated cells. STK16-IN-1 (2.5-5  $\mu\text{M}$ ) potentiates the antiproliferative effects of colchicine (Item No. 9000760), paclitaxel (Item No. 10461), doxorubicin (Item No. 15007), and cisplatin (Item No. 13119) in MCF-7 cells.

### Reference

1. Liu, F., Wang, J., Yang, X., *et al.* Discovery of a highly selective STK16 kinase inhibitor. *ACS Chem. Biol.* **11**(6), 1537-1543 (2016).

#### 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, 11/07/2022

#### 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