

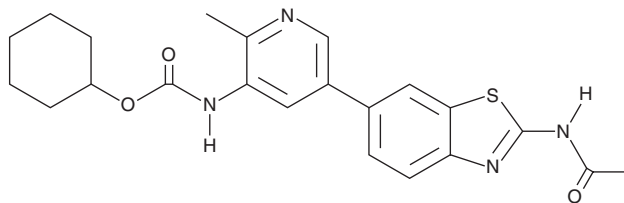
PRODUCT INFORMATION



PK68

Item No. 36653

CAS Registry No.: 2173556-69-7
Formal Name: N-[5-[2-(acetylamino)-6-benzothiazolyl]-2-methyl-3-pyridinyl]-carbamic acid, cyclohexyl ester
MF: C₂₂H₂₄N₄O₃S
FW: 424.5
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

PK68 is supplied as a solid. A stock solution may be made by dissolving the PK68 in the solvent of choice, which should be purged with an inert gas. PK68 is soluble in the organic solvent DMSO. PK68 is slightly soluble in methanol and acetonitrile.

Description

PK68 is an inhibitor of receptor-interacting serine/threonine kinase 1 (RIPK1; IC₅₀ = 90 nM).¹ It is selective for RIPK1 over RIPK3 and a panel of 369 kinases at 1,000 nM. PK68 inhibits necroptosis induced by TNF-α, a Smac mimetic, and z-VAD in HT-29 colon cancer cells (EC₅₀ = 23 nM). It reduces the number of lung metastases in a B16/F10 murine melanoma model when administered at a dose of 5 mg/kg. PK68 (1 mg/kg) increases survival and prevents decreases in body temperature and increases in serum II-1β levels in a mouse model of Tnf-α-induced systemic inflammatory response syndrome (SIRS).

Reference

1. Hou, J., Ju, J., Zhang, Z., *et al.* Discovery of potent necroptosis inhibitors targeting RIPK1 kinase activity for the treatment of inflammatory disorder and cancer metastasis. *Cell Death Dis.* **10(7)**, 493 (2019).

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, 11/14/2023

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