

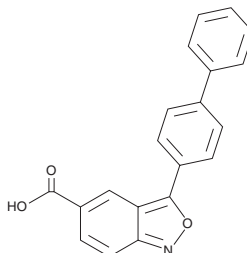
PRODUCT INFORMATION



UNC7467

Item No. 41597

CAS Registry No.: 2922283-43-8
Formal Name: 3-[1,1'-biphenyl]-4-yl-2,1-benzisoxazole-5-carboxylic acid
MF: $C_{20}H_{13}NO_3$
FW: 315.3
Purity: $\geq 98\%$
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

UNC7467 is supplied as a solid. A stock solution may be made by dissolving the UNC7467 in the solvent of choice, which should be purged with an inert gas. UNC7467 is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

UNC7467 is an inhibitor of inositol hexakisphosphate kinase 1 (IP6K1) and IP6K2 (IC_{50} s = 8.9 and 4.9 nM, respectively).¹ It is selective for IP6K1 and IP6K2 over IP6K3 and inositol-triphosphate 3-kinase A (ITPKA; IC_{50} s = 1,320 and 832 nM, respectively), as well as a panel of 30 other kinases at 1 μM , but does inhibit inositol phosphate multi-kinase (IPMK; IC_{50} = 24.4 nM). *In vivo*, UNC7467 (5 mg/kg per day) reduces body weight, fat mass, epididymal- and inguinal-adipose tissue deposit weights, and liver weight, as well as blood glucose levels in a pyruvate tolerance test after four weeks, in a mouse model of diet-induced obesity.

Reference

1. Zhou, Y., Mukherjee, S., Huang, D., *et al.* Development of novel IP6K inhibitors for the treatment of obesity and obesity-induced metabolic dysfunctions. *J. Med. Chem.* **65**(9), 6869-6887 (2022).

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, 08/26/2024

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