

PRODUCT INFORMATION

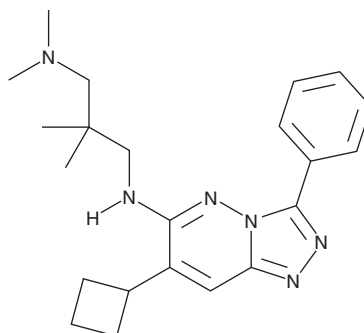


Akt-I-1

Item No. 41016

CAS Registry No.: 473382-39-7
Formal Name: N³-(7-cyclobutyl-3-phenyl-1,2,4-triazolo[4,3-b]pyridazin-6-yl)-N¹,N¹,2,2-tetramethyl-1,3-propanediamine

Synonym: Akt-IN-1
MF: C₂₂H₃₀N₆
FW: 378.5
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

Akt-I-1 is supplied as a solid. A stock solution may be made by dissolving the Akt-I-1 in the solvent of choice, which should be purged with an inert gas. Akt-I-1 is sparingly soluble (1-10 mg/ml) in ethanol.

Description

Akt-I-1 is an inhibitor of Akt1 (IC₅₀ = 4.6 μM).¹ It is selective for Akt1 over Akt2, Akt3, and serum/glucocorticoid-regulated kinase (SGK; IC₅₀s = >250 μM for all).

Reference

1. Barnett, S.F., Defeo-Jones, D., Fu, S., *et al.* Identification and characterization of pleckstrin-homology-domain-dependent and isoenzyme-specific Akt inhibitors. *Biochem. J.* **385**(Pt. 2), 399-408 (2005).

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, 06/26/2024

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