

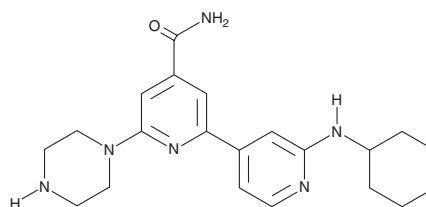
PRODUCT INFORMATION



BPKDi

Item No. 33707

CAS Registry No.: 1201673-28-0
Formal Name: 2'-(cyclohexylamino)-6-(1-piperazinyl)-[2,4'-bipyridine]-4-carboxamide
MF: C₂₁H₂₈N₆O
FW: 380.5
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

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

Description

BPKDi is an inhibitor of protein kinase D (PKD; IC₅₀s = 1, 9, and 1 nM for PKD1, PKD2, and PKD3, respectively).¹ It is selective for PKD over PKCδ and PKCε, as well as over a panel of additional calcium/calmodulin-dependent protein kinase (CaMK) superfamily members, at 1 μM. It inhibits phenylephrine-induced phosphorylation and nuclear export of histone deacetylase 4 (HDAC4) and HDAC5 in, as well reduces phenylephrine-induced hypertrophy of, isolated neonatal rat ventricular myocytes when used at a concentration of 1 μM.

Reference

1. Monovich, L., Vega, R.B., Meredith, E., *et al.* A novel kinase inhibitor establishes a predominant role for protein kinase D as a cardiac class IIa histone deacetylase kinase. *FEBS Lett.* **584(3)**, 631-637 (2010).

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, 12/05/2022

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