

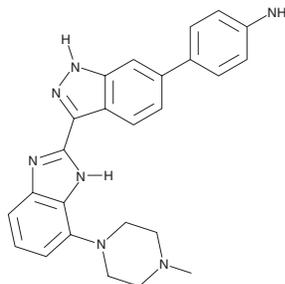
PRODUCT INFORMATION



PKC-9

Item No. 25556

CAS Registry No.: 1132609-87-0
Formal Name: 4-[3-[7-(4-methyl-1-piperazinyl)-1H-benzimidazol-2-yl]-1H-indazol-6-yl]-benzenamine
MF: C₂₅H₂₅N₇
FW: 423.5
Purity: ≥95%
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

PKC-9 is supplied as a solid. A stock solution may be made by dissolving the PKC-9 in the solvent of choice. PKC-9 is soluble in the organic solvent DMSO, which should be purged with an inert gas.

Description

PKC-9 is an inhibitor of PKC ζ that is 8,000-, 20,671-, 711-, 918-, 1,895-, 12,424-, 10-, and 1,218-fold selective for PKC ζ over PKC α , PKC β II, PKC γ , PKC δ , PKC ϵ , PKC η , PKC ι , and PKC θ , respectively.¹ It also inhibits 27 additional kinases in a panel of 37 kinases when used at a concentration of 10 μ M.

Reference

1. Trujillo, J.I., Kiefer, J.R., Huang, W., *et al.* 2-(6-Phenyl-1H-indazol-3-yl)-1H-benzo[d]imidazoles: Design and synthesis of a potent and isoform selective PKC- ζ inhibitor. *Bioorg. Med. Chem. Lett.* **19(3)**, 908-911 (2009).

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, 12/19/2022

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