

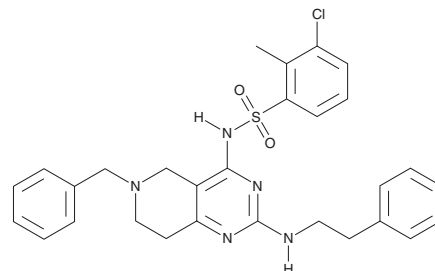
PRODUCT INFORMATION



CaMKII-IN-1

Item No. 37303

CAS Registry No.: 1208123-85-6
Formal Name: 3-chloro-2-methyl-N-[5,6,7,8-tetrahydro-2-[(2-phenylethyl)amino]-6-(phenylmethyl)pyrido[4,3-d]pyrimidin-4-yl]-benzenesulfonamide
Synonyms: Calcium/Calmodulin-dependent Protein Kinase II Inhibitor 1, CaMKII Inhibitor 1
MF: C₂₉H₃₀ClN₅O₂S
FW: 548.1
Purity: ≥95%
UV/Vis.: λ_{max}: 267 nm
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

CaMKII-IN-1 is supplied as a solid. A stock solution may be made by dissolving the CaMKII-IN-1 in the solvent of choice, which should be purged with an inert gas. CaMKII-IN-1 is soluble in DMSO.

Description

CaMKII-IN-1 is a calcium/calmodulin-dependent kinase II (CaMKII) inhibitor (IC₅₀ = 0.063 μM).¹ It is selective for CaMKII over CaMKIV, myosin light-chain kinase (MLCK), p38α, Akt1, and PKC (IC₅₀s = >60, 36, 11, 30, and 21 μM, respectively).

Reference

- Asano, S., Komiya, M., Koike, N., *et al.* 5,6,7,8-Tetrahydropyrido[4,3-d]pyrimidines as novel class of potent and highly selective CaMKII inhibitors. *Bioorg. Med. Chem. Lett.* **20(22)**, 6696-6698 (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, 04/18/2023

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