

PRODUCT INFORMATION



PKC θ Inhibitor (hydrochloride)

Item No. 36165

CAS Registry No.: 2253640-49-0
Formal Name: N⁴-[[4-(aminomethyl)cyclohexyl]methyl]-5-nitro-N²-[[2-(trifluoromethoxy)phenyl]methyl]-2,4-pyrimidinediamine, monohydrochloride

Synonym: Protein Kinase C θ Inhibitor

MF: C₂₀H₂₅F₃N₆O₃ • HCl

FW: 490.9

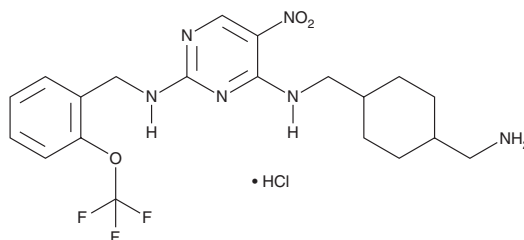
Purity: \geq 98%

UV/Vis.: λ_{max} : 268, 363 nm

Supplied as: A solid

Storage: -20°C

Stability: \geq 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

PKC θ inhibitor (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the PKC θ inhibitor (hydrochloride) in the solvent of choice, which should be purged with an inert gas. PKC θ inhibitor (hydrochloride) is soluble in the organic solvent DMSO at a concentration of approximately 1 mg/ml.

Description

PKC θ inhibitor is an inhibitor of PKC θ (IC₅₀ = 18 nM).¹

Reference

1. Cywin, C.L., Dahmann, G., Prokopowicz, A.S., III, *et al.* Discovery of potent and selective PKC- θ inhibitors. *Bioorg. Med. Chem. Lett.* **17(1)**, 225-230 (2007).

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