

PRODUCT INFORMATION



PKCθ Pseudosubstrate Inhibitor

Item No. 17483

Formal Name: H-leucyl-histidyl-glutamyl-arginyl-glycyl-alanyl-
isoleucyl-lysyl-glutamyl-alanyl-arginyl-arginyl-
tryptophyl-arginyl-lysyl-leucine

Synonym: Protein Kinase C θ Pseudosubstrate Inhibitor

MF: C₉₅H₁₆₅N₃₉O₂₀

FW: 2,173.6

Purity: ≥95%

Supplied as: A lyophilized powder

Storage: -20°C

Stability: ≥4 years

H—Leu—His—Gln—Arg—Arg—Gly—Ala—Ile—Lys—
Gln—Ala—Arg—Arg—Trp—Arg—Lys—Leu—OH

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

Description

PKCθ pseudosubstrate inhibitor is a synthetic peptide that corresponds to amino acid residues of the pseudosubstrate domain of protein kinase Cθ (PKCθ). It selectively, reversibly, and substrate-competitively inhibits PKCθ activity and, thus, is used to delineate the signaling functions of PKCθ, which are thought to be involved in several autoimmune disorders.¹

Reference

1. Hage-Sleiman, R. Hamze, A.B. Reslan, L., *et al.* The novel PKCθ from benchtop to clinic. *J. Immunol. Res.* 348798 (2015).

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, 10/25/2022

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