

PRODUCT INFORMATION



PKC δ Substrate (trifluoroacetate salt)

Item No. 37513

Formal Name:	L-alanyl-L-arginyl-L-arginyl-L-lysyl-L-arginyl-L-lysylglycyl-L-seryl-L-phenylalanyl-L-phenylalanyl-L-tyrosylglycylglycine	
Synonym:	Protein Kinase C δ Substrate	H—Ala—Arg—Arg—Lys—Arg—Lys—Gly—Ser—Phe—
Peptide Sequence:	ARRKRKGSFFYGG-OH	Tyr—Gly—Gly—OH
MF:	C ₆₉ H ₁₀₈ N ₂₄ O ₁₆ • XCF ₃ COOH	
FW:	1,529.8	• XCF ₃ COOH
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

PKC δ Substrate (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the PKC δ substrate (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

PKC δ substrate is a peptide substrate for PKC δ .¹ It has been used to monitor PKC δ activity in cell-free and cell-based assays.

Reference

1. Miralem, T., Lerner-Marmarosh, N., Gibbs, P.E.M., *et al.* The human biliverdin reductase-based peptide fragments and biliverdin regulate protein kinase C δ activity: The peptides are inhibitors or substrate for the protein kinase C. *J. Biol. Chem.* **287(29)**, 24698-24712 (2012).

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, 10/11/2023

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