

PRODUCT INFORMATION



[Ser¹⁵⁹]-PKCε (149-164) (trifluoroacetate salt)

Item No. 37508

Formal Name:	L-α-glutamyl-L-arginyl-L-methionyl-L-arginyl-L-prolyl-L-arginyl-L-lysyl-L-arginyl-L-glutamylglycyl-L-seryl-L-valyl-L-arginyl-L-arginyl-L-arginyl-L-valine, trifluoroacetate salt	
Synonyms:	[Ser ¹⁵⁹]-PKCε Pseudosubstrate Peptide (149-164), [Ser ¹⁵⁹]-Protein Kinase Cε Inhibitor Peptide (149-164)	H—Glu—Arg—Met—Arg—Pro—Arg—Lys—Arg—Gln—Gly— Ser—Val—Arg—Arg—Arg—Val—OH
Peptide Sequence:	ERMRRPKRQGSVRRRV-OH	
MF:	C ₈₃ H ₁₅₅ N ₃₉ O ₂₁ S • XCF ₃ COOH	• XCF ₃ COOH
FW:	2,067.4	
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

[Ser¹⁵⁹]-PKCε (149-164) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the [Ser¹⁵⁹]-PKCε (149-164) (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

[Ser¹⁵⁹]-PKCε (149-164) is a peptide substrate for PKCγ, PKCδ, PKCε, PKCη, and PKCξ, serine/threonine kinases involved in various cell processes such as proliferation, differentiation, and survival.¹⁻³ It has been used as a target substrate for the quantification of PKCγ, PKCδ, PKCε, PKCη, or PKCξ activity in cell-free or cell-based assays.

References

1. Nakanishi, H., Brewer, K.A., and Exton, J.H. Activation of the ζ isozyme of protein kinase C by phosphatidylinositol 3,4,5-trisphosphate. *J. Biol. Chem.* **268**(1), 13-16 (1993).
2. Luo, Y., Smith, R.A., Guan, R., *et al.* Pseudosubstrate peptides inhibit Akt and induce cell growth inhibition. *Biochemistry* **43**(5), 1254-1263 (2004).
3. Lee, A.W. The role of atypical protein kinase C in CSF-1-dependent Erk activation and proliferation in myeloid progenitors and macrophages. *PLoS One* **6**(10), e25580 (2011).

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, 10/30/2023

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