



Product Data Sheet

Product Name:	Protein Kinase C ϵ Peptide Substrate	
Catalog Number:	AS-27183 (1 mg)	Lot Number: See label on vial
Sequence:	H-Glu-Arg-Met-Arg-Pro-Arg-Lys-Arg-Gln-Gly-Ser-Val-Arg-Arg-Arg-Val-OH (3-letter code) ERM RPRKRQGSVRRRV (1-letter code)	
Molecular Weight:	2067.5	
Peptide Purity:	>95%	
Appearance:	Lyophilized white powder	

Peptide Reconstitution: Protein Kinase C ϵ peptide is freely soluble in water.

Storage: Protein Kinase C ϵ peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at -20°C or lower. Reconstituted peptide can be aliquoted and stored at -20°C or lower.

Description: The peptide sequence, ERM RPRKRQGSVRRRV (cat# AS-27183), corresponds to pseudosubstrate region of the ϵ -isotype of protein kinase C (PKC ϵ) with an alanine to serine substitution. PKC ϵ exhibits an apparent specificity for the native peptide ($K_m = 68 \mu\text{M}$). Ref: Schaap, D. and P.J. Parker, J. Biol. Chem. 265, 7301 (1990); Schaap, D. et al. FEBS Lett. 243, 351 (1989).

Related Products:

For Research Use Only