





Protein Kinase C Isoforms Domain Structures

		Mr	Activators
cPKC		α	Ca^{2+}
		β	DAG
		β II	PS
		γ	lysoPC FFA
nPKC		δ	PS
		θ	DAG FFA
nPKC		η	PS
		ϵ	FFA
aPKC		ζ	PS
		λ	
		ι	