

PI-Specific Phospholipase C Catalytic Domain Alignment

PLCδ1	297	MGQPLSHYLVSSSHNTYLLEDQLAGPSSTEAYIRALCKGCRCLELDCWDG--PNQEPIIYHGTYFTSKIF				
PLCδ4	292	MTQPLNHYFICSSHNTYLVGDQLCGQSSVEGYIRALKRGCRCEVDVWDG--PSGEPVYVHGHTLTSRIL				
PLCL2	302	MKQPLSHYFINSSHNTYLIEDQFRGPSDITGYIRALKMGCRSVELDVWDG--PDNEPVIYTGHTMTSQIV				
PLCγ1	322	MNNPLSHYWISSSHNTYLTGDQFSSESSLEAYARCLRMGCRCELDLDCWDG--PDGMPVIYHGHTLTTKIK				
PLCβ1	318	MSQPLSHYFINSSHNTYLTAGQLAGNSSVEMYRQVLLSGCRCELDLDCWKGRTAEEEEPVITHGFTMTTEIS				
PLCβ4	315	MDHPLAHYFISSSHNTYLTGRQFSGKSSVEMYRQVLLAGCRCELDLDCWDGKGEDQEPITTHGKAMCTDIL				
PLCδ1	366	FCDVLRAIRDYAFKASPYPVILSLENHC-TLEQQRVMARHLHAILGPMLLNRPLDGVT----NSLP				
PLCδ4	360	FKDVVATVAQYAFQTSYYPVILSLETHC-SWEQQQTMARHLTEILGEQLLSTTLDGVLV---TQLP				
PLCL2	370	FRSVIDIINKYAFFASEYPVILCLENHC-SIKQQKVMVQHMKKLLGDKLYTTPNVEE----SYLP				
PLCγ1	390	FSDVLHTIKEHAFVASEYPVILSIEDHC-SIAQQRNMAQYFKKVLGDTLLTKPVEISA----DGLP				
PLCβ1	388	FKEVIEAIAECAFKTSPFPILLSFENHVDSPKQQAKMAEYCRLIFGDALLMEPLEKYPLESGVPLP				
PLCβ4	385	FKDVIQAIKETAFVTSEYPVILSLENHC-SKYQQYKvSKYCEDLFGDLLLKQALESHPLEPGRALP				
PLCδ1	427	SPEQLKKGILLKGGKLL	442	PLCδ1	492	LSDMVIYCKSVHFGGFSSPGTPGQA
PLCδ4	422	SPEELRRKILVKGGKLL	437	PLCδ4	493	LSSLVIYLKSVSFRSFTHS-KEHYH
PLCL2	431	SPDVLKKGILIKAKKLL	446	PLCL2	492	LSELVSICKSVQFKEFQMS-FQVQK
PLCγ1	451	SPNQLKRKILIKHKKLL	466	PLCγ1	953	LSELvVYCRPVPFDEEKIG-TERAC
PLCβ1	454	SPMDLMYKILVKNKKK	469	PLCβ1	540	MSNLVNYIQPVKFESFEIS-KKRNK
PLCβ4	450	SPNDLKRKILIKNKRL	465	PLCβ4	565	LSTMINYAQPVKvQGFHVA-EERNI
PLCδ1	517	FYEMASFSENRALRLLQES-GNGFVRHNVGHL	SRIYYPAGWRTDSSSNYSPVEMWN			
PLCδ4	517	FYEISSFSETKAKRLIKEA-GNEFVQHNTWQL	SRVYPSGLRTDSSSNYNPQELWN			
PLCL2	516	YWEVCSFNEVLASKYANEN-PGDFVNYNKRFL	ARVFPSPMRIDSSNMNPQDFWK			
PLCγ1	977	YRDMSSFPEPKAEKYVNKAKGKKFLQYNRLQ	LRIYYPKGRDSSSNYDPLPMWI			
PLCβ1	564	SFEMSSFVETKGLEQLTKS-PVEFVEYNKMQL	SRIYYPKGRVDSNYPQLFWN			
PLCβ4	589	HYNMSSFNESVGLGYLKTH-AIEFVNYNKRQ	MSRIYYPKGRVDSNYPQIFWN			
PLCδ1	570	GGCQIVALNFQTPGPEMDVYQDRFQDNGACGYV	LKPAFLRD	610		
PLCδ4	570	AGCQMVAMNMQTAGLEMDICDGHFRQNGGCGYV	LKPDFLRD	610		
PLCL2	569	CGCQIVAMNFQTPGLMMDLNIWFQNGNCGYV	LRPAIMRE	609		
PLCγ1	1031	CGSQLVALNFQTPDKPMDMNQALFMTGRHCGYV	LQPSTMRD	1071		
PLCβ1	617	AGCQMVVALNFQTMDLAMQINMGMYEYNGKSGYR	LKPEFMRR	657		
PLCβ4	643	AGCQMVSLNYQTPDLAMQLNQGKFEYNGSCGYL	LKPDFMRR	682		

Rat
Phosphatidylinositide-Specific Phospholipase C-delta1
 (residues 133-756, white)
 Inositol-1,4,5-trisphosphate
 (black)
PDB: 1DJX

