

Phospholipase A₂ Catalytic Domain Alignment

		●●		●●●●	●	●●●●●●●●●●●●●●●●●●●●				
PLA2-IIA	1	MKTL	LL	LAVIMIFGLLQAHG--	NLVNFHRMIK-LTTG	KEAALS	YGFY	GCHCGV	GGRG	SPKDATDR
PLA2-IIIE	1	MKSP	HVLVFLCLL-VALVTG--	NLVQFGVMIE-KMTG-	KSALQY	NDYGCY	CGIGGS	HWPVDQTDW		
PLA2-G1B	1	MKLL	VLA	LLTVAAADSGISPR	AVWQFRKMIKCVIP	GSDFLE	YNNYGCY	CGLGGS	SGTPVDEL	DK
PLA2-IID	1	MELAL	LCGLVVMAGVIPIQG--	GILN	LNMVK-QVTG	KMPILSY	WPYGCH	CGLGGR	GQPKDATDW	
PLA2-V	1	MKGL	LPLAWFLACSVPAVQG--	GLLDL	KSMIE-KVTG	KNALTNY	GFYGCY	CGWGGR	GT	PKDGTDW

		●●●●●●●●		●●●●●●●●	●●●●●●●●	●●●●●●●●			
PLA2-IIA	63	CCVTHDCCYKRL	E	KRG-----	CGTKFLSYKFSNS	SGSRITCA-KQD	SCRSQ	LC	CDKAA
PLA2-IIIE	61	CCHAHDCCYGRLE	E	KL-----	CEPKLEKYLFSV	SERGI	FCA-GRTT	CQRLT	CECDKRA
PLA2-G1B	66	CCQTHDNCYDQAK	KL	LDSCFLLDNPY	HTYSYSCSGSAIT	CSSKNKE	CEAFI	CNCDR	NA
PLA2-IID	63	CCQTHDCCYDHL	K	TQG-----	CSIYKDYRYNFS	QGNIH	CSDKGS	WCEQQL	CACDKEV
PLA2-V	63	CCWAHDHCYGRLE	E	EKG-----	CNIRTQSYKYR	FAWGV	VTCE-PGPF	CHVNL	CACDRKL

		●●●●	●●	
PLA2-IIA	115	ATCFARNKTT	YNKKYQ	YYSNKHCRGS 140
PLA2-IIIE	113	ALCFRRNLGTY	NRKYAHYP	NKLCTGP 138
PLA2-G1B	125	AICFSK--	APYNKAHKNL	DTKKYCQS 148
PLA2-IID	116	AFCLKRNLDTY	QKRLRFYWR	PHCRGQ 141
PLA2-V	115	VYCLKRNLRSYN	PQYQYFPN	ILCS-- 138

Human Phospholipase
A2 - IIA
(residues 21-144, white)
PDB: 1BBC

