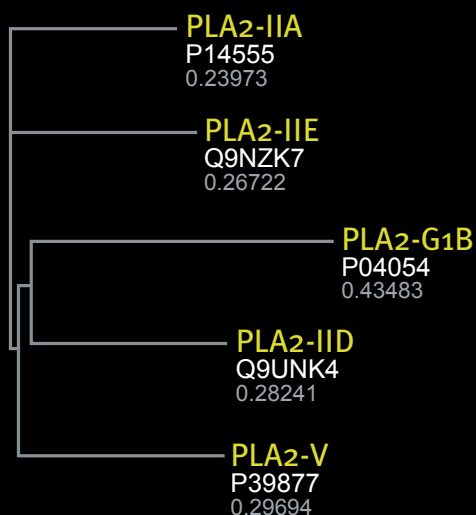




www.kinexus.ca
1-866-kinexus TOLL FREE PHONE

Human Phospholipase A2 - IIA
(Residues 21-144)
PDB: 1BBC

Phospholipase A2 Catalytic Domains



PLA2-IIA	1	MKTLALLAVIMIFGLLQAHG--NLVNFHRMIK-LTTGKEAALSYGFYGCCHCGVGGRGSPKDADR
PLA2-IIIE	1	MKSPHVLVFLCLL-VALVTG--NLVQFGVMIE-KMTG-KSALQYNDYGCYCGIGGSHWPVDQTDW
PLA2-G1B	1	MKLLVLAVLLTVAAADSGISPRVWQFRKMIKCVIPGSDPFLEYNNYGCYCGLGSGTVPDELDK
PLA2-IIID	1	MELALLCGLVVMAGVIPIQG--GILNLNKMVK-QVTGKMPILSWPYGCHCGLGGRGQPKDADR
PLA2-V	1	MKGLLPLAWFLACSVPAVQG--GLLDLKSME-KVTGKNALTNYGFYGCYCGWGGRGTPKDGTDW
PLA2-IIA	63	CCVTHDCCYKRLEKRG-----CGTKFLSYKFSNSGSRITCA-KQDSCRSQLECDKAA
PLA2-IIIE	61	CCHADCCYGRLEKLG-----CEPKLEKYLFSVSESGIFCA-GRITTCQRLTCECDKRA
PLA2-G1B	66	CCQTHDNCYDQAKKLDSCFLLDNPYTHYYSYCSGSAITCSSLNKECEAFICNCDRNA
PLA2-IIID	63	CCQTHDCCYDHLKTQG-----CSIYKDYRYNFSQGNITHCSDKGSWCEQQLCACDKEV
PLA2-V	63	CCWAHDHCYGRLEEKG-----CNIRTQSYKYRFAWGVVTC-PGPFCHVNLACDRKL
PLA2-IIA	115	ATCFARNKTTYNKKYQYYSNKHCRGS 140
PLA2-IIIE	113	ALCFRRNLGTYNRYAHYPNKLCTGP 138
PLA2-G1B	125	AICFSK--APYNKAHKNLDTKKYCQS 148
PLA2-IIID	116	AFCLKRNLDTYQKRLRFYWRPHCRGQ 141
PLA2-V	115	VYCLKRNLRSYNPQYQYFPNILCS-- 138