

Occurrence of Selected Binding Domains in Protein Kinases

Domain	SH2	FHA	Polo Box	WD40	SH3	C1	C2	PH	PY
Domain Specificity	Phos-Tyr	Phos-Ser	Phos-Ser	Phos-Ser	Poly-Pro	Phospho-lipid	Phospho-lipid	Phospho-lipid	Phospho-lipid
AGC	0	0	0	0	0	15	8	12	2
ATYPICAL	0	0	0	0	0	0	1	1	0
CAMK	0	1	0	0	4	3	0	6	0
CK1	0	0	0	0	0	0	0	0	0
CMGC	0	0	0	0	0	0	0	0	0
OTHER	0	0	3	1	0	0	0	0	1
RGC	0	0	0	0	0	0	0	0	0
STE	0	0	0	0	0	0	0	0	0
TK	28	0	0	0	21	0	0	4	0
TKL	0	0	0	1	4	5	0	0	0
Separate proteins	28	1	3	2	29	23	9	23	3
Total number	30	1	6	13	29	32	9	24	3
Average length (aa)	86	64	68	45	59	52	114	121	34
Shortest length (aa)	76	64	64	32	53	46	104	90	31
Longest length (aa)	106	64	72	61	68	57	137	130	39

