

Human Protein Phosphatome

Group	Number in group	Number with TM domains	Number with two catalytic domains
DUAL SPECIFICITY	34	0	0
RECEPTOR-TYROSINE	21	21	11
NON-RECEPTOR TYROSINE	28	4	2
SERINE/THREONINE	41	0	0
IPP-RELATED	14	0	0
Total	138	26	13

