

Human Protein Kinase Families

Group	Number in group	Number with TM domains	Number with two catalytic domains	Number pseudo genes	Known Pho-Ser sites	Known Pho-Thr sites	Known Pho-Tyr sites
AGC	63	0	0	6	44	38	17
ATYPICAL	38	2	0	1	15	2	6
CAMK	74	0	8	39	24	31	3
CK1	12	1	0	5	0	0	0
CMGC	61	0	0	12	41	10	18
OTHER	82	4	1	21	17	11	6
RGC	5	5	0	3	0	0	0
STE	47	0	0	6	29	19	5
TK	90	55	4	5	38	5	215
TKL	43	12	0	6	39	9	5
Total	515	79	13	104	247	125	275

