

Protein Kinase Dimerization

Group	Examples of Dimerized Human Protein Kinases
AGC	DMPK1, GRPK6, MSK1+2, PKA, PKB, PKG, ROCK2, RSK1-4,
ATYPICAL	ATM, ATR, DNAPK
CAMK	CHK2, DAPK1-3, PHK γ , smMLCK
CK1	N.D. reported
CMGC	ERK1, ERK2, ERK3, JNK, p38
OTHER	CK2, GCN2, HRI, IKK $\alpha+\beta$, IRAK, IRE1+2, NEK2, PEK, PKR, Plk4, SgK069
RGC	ANPa+b, CYGD+F, HSER
STE	ASK1, MEKK2+4, MST1+2, PAK1, SLK, TAO1-3,
TK	All 58 receptor-tyrosine kinases, JAK1-3, Tyk, Abl, FAK, Bmx, Btk
TKL	All 12 receptor-serine kinases, DLK, Ksr1, LZK, MLK3, Raf1, RafA, RafB
Total	Over 145 reported human protein kinases

- Most protein kinases can dimerize or oligomerize
- N.D.able exceptions appear to be the Src, PKC and CDK families