

# Cyclin-Dependent Kinase Family Domains

## CDK Family, CDC2 Subfamily



## CDK Family, CDK10 Subfamily



## CDK Family, CDK4 Subfamily



## CDK Family, CDK5 Subfamily



## CDK Family, CDK7 Subfamily



## CDK Family, CDK8 Subfamily



## CDK Family, CDK9 Subfamily



## CDK Family, CRK7 Subfamily



## CDK Family, PITSLRE Subfamily



## CDK Family, TAIRE Subfamily

