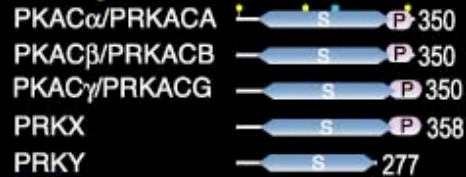


# AGC Group - Human Serine Kinases - 2

## NDR Family



## PKA Family



## PDK1 Family



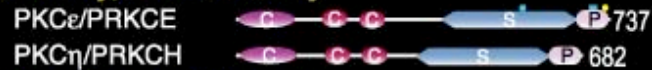
## PKC Family, Alpha Subfamily



## PKC Family, Delta Subfamily



## PKC Family, Eta Subfamily



## PKC Family, Iota Subfamily



## PKG Family



## PKN Family

