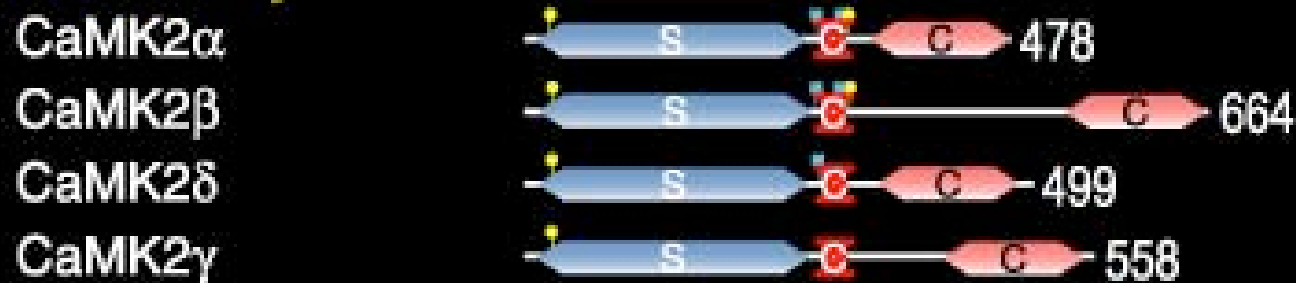


Ca²⁺/Calmodulin-Regulated Protein Kinases 1 + 2

CAMK1 Family



CAMK2 Family



Calmodulin-bind
CaMK2 assoc.
SerThrKinase

