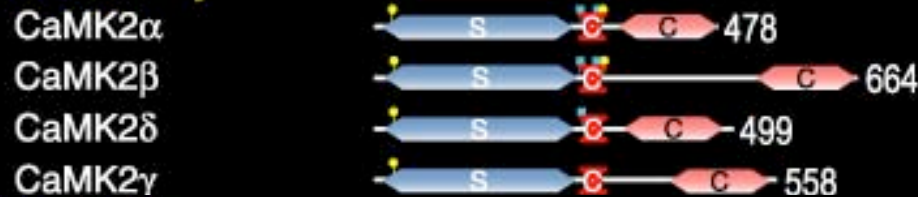


# Ca<sup>2+</sup>/Calmodulin-Regulated Protein Kinases Domains

## CAMK1 Family



## CAMK2 Family



## DAPK Family



## MLCK Family



## PHK Family

