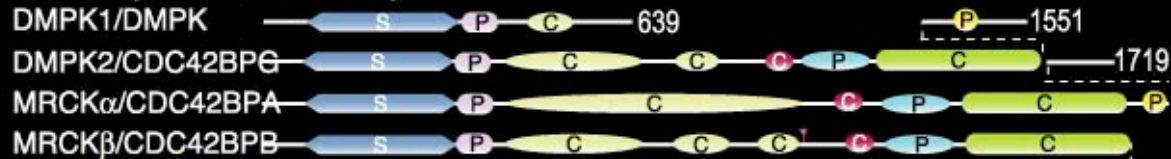


AGC Group - Human Serine Kinases - 1

AKT Family



DMPK Family, GEK Subfamily



DMPK Family, ROCK Subfamily



DMPK Family, Unique Subfamily



GRK Family, BARK Subfamily



GRK Family, GRK Subfamily



MAST Family

