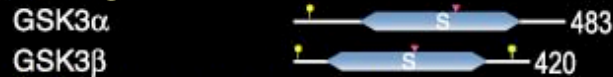
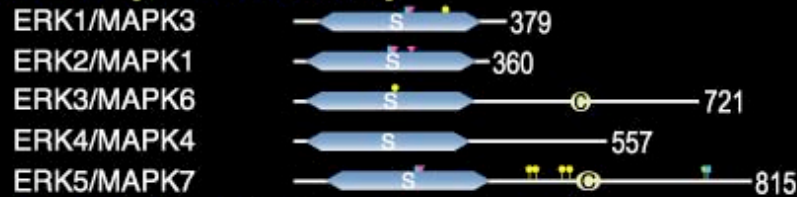


MAP Kinase Family Domains

GSK Family



MAPK Family, ERK Subfamily



MAPK Family, ERK7 Subfamily



MAPK Family, JNK Subfamily



MAPK Family, nmo Subfamily



MAPK Family, p38 Subfamily



RCK Family



SRPK Family

