

AGC Group - Human Serine Kinases - 3

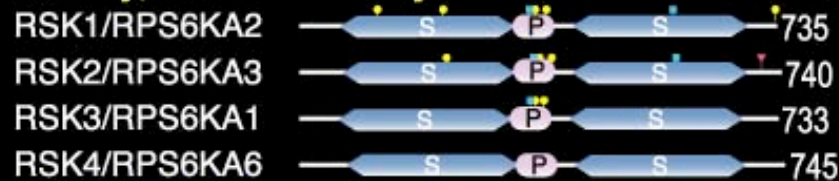
RSK Family, MSK Subfamily



RSK Family, p70 Subfamily



RSK Family, RSK Subfamily



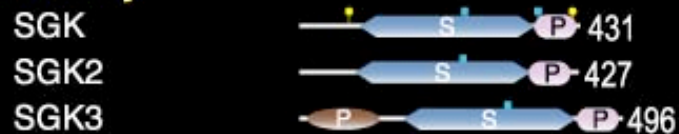
RSK Family, Unique Subfamily



RSKL Family



SGK Family



YANK Family

