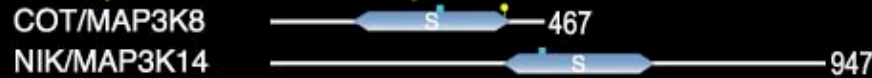


# Ste20 Group - Human Protein-Serine Kinases - 1

## STE Family, Unique Subfamily



## STE11 Family



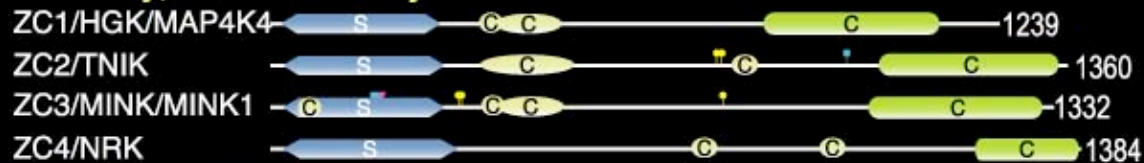
## STE20 Family, FRAY Subfamily



## STE20 Family, KHS Subfamily



## STE20 Family, MSN Subfamily



## STE20 Family, MST Subfamily

