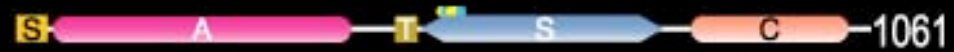


RGC Group - Human Receptor Serine Kinases

RGC Family

ANP α /NPR1



ANP β /NPR2



CYGD/GUCY2D



CYGF/GUCY2F



HSER/GUCY2C

