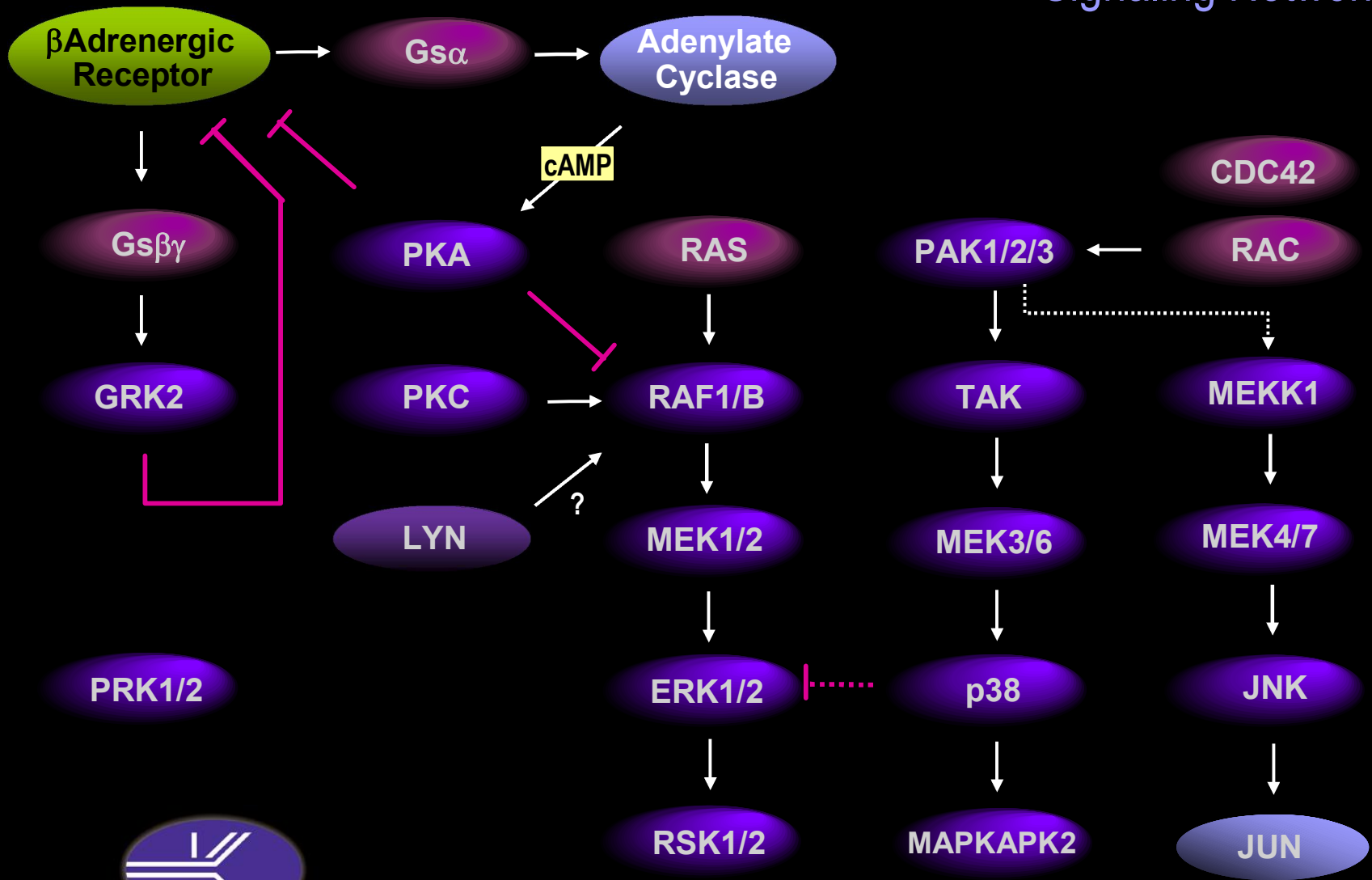


Signaling Networks



KINEXUS