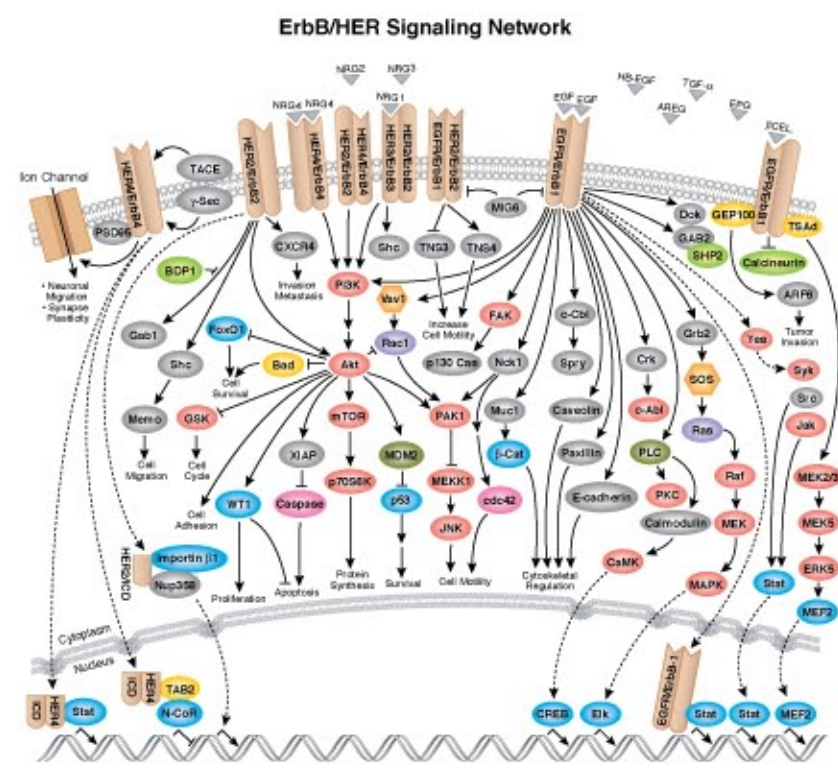


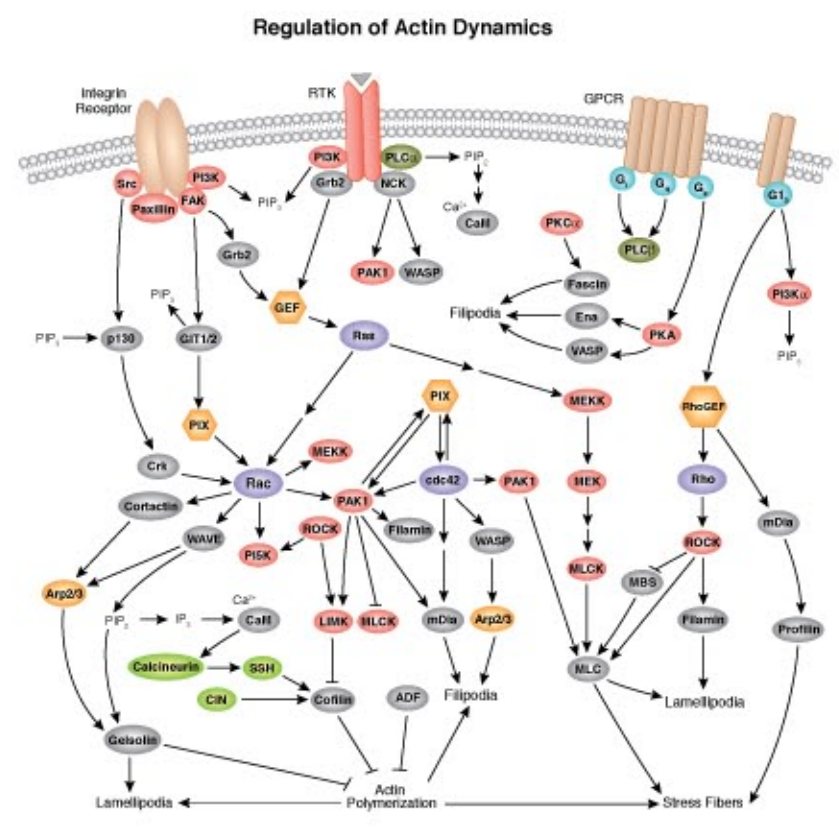


Signal Transduction Pathways

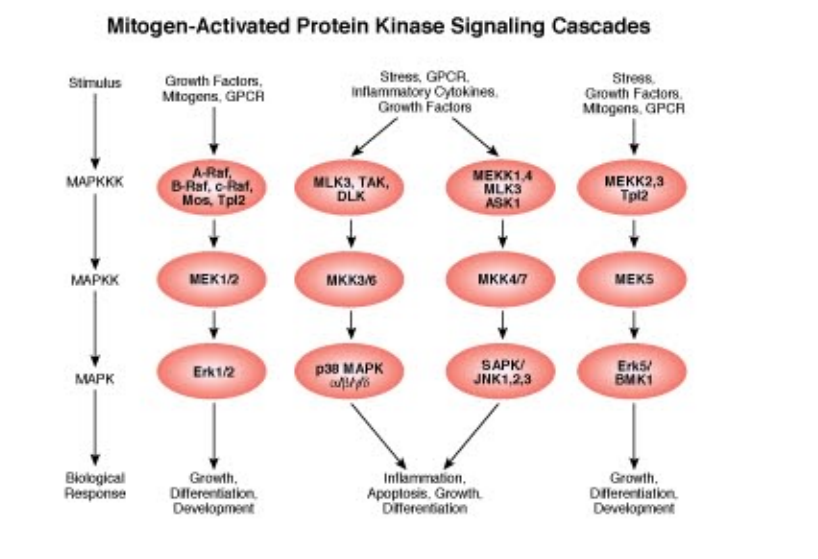
RECEPTOR TYROSINE KINASE



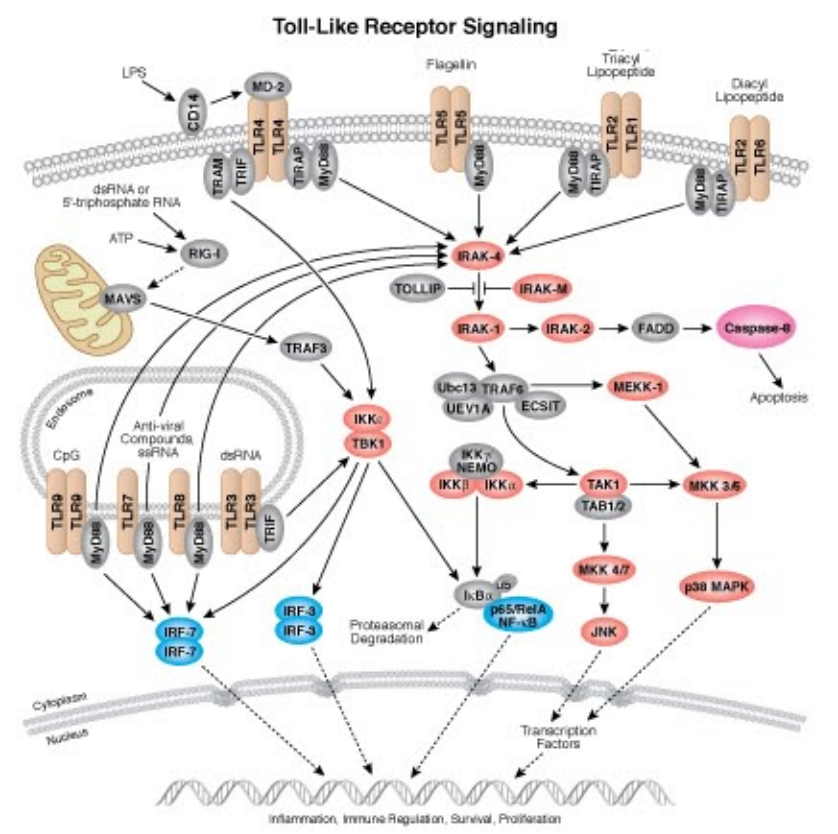
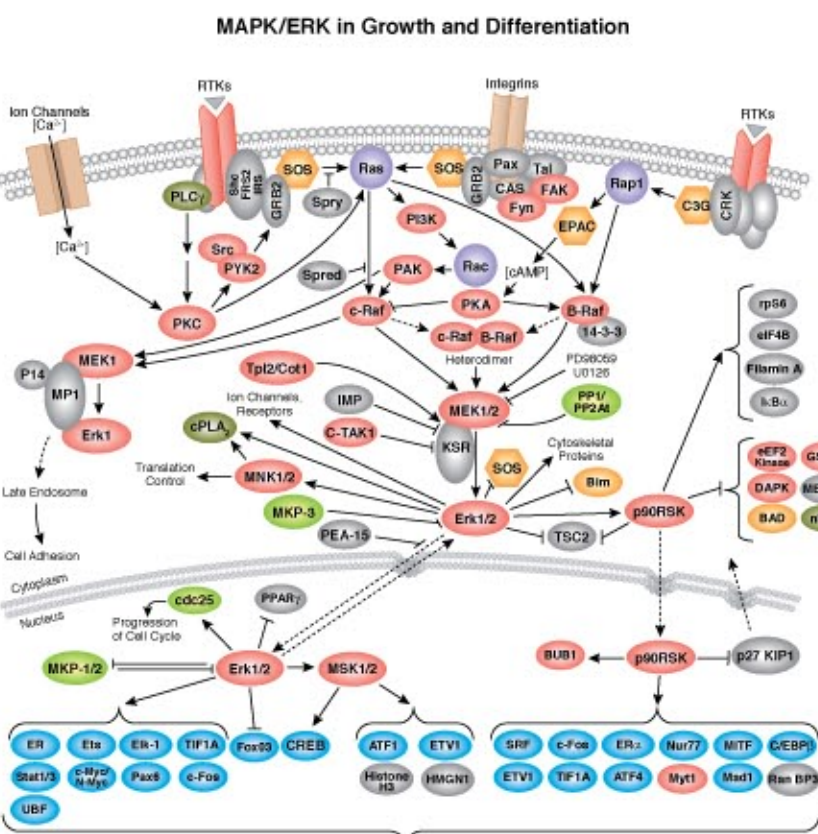
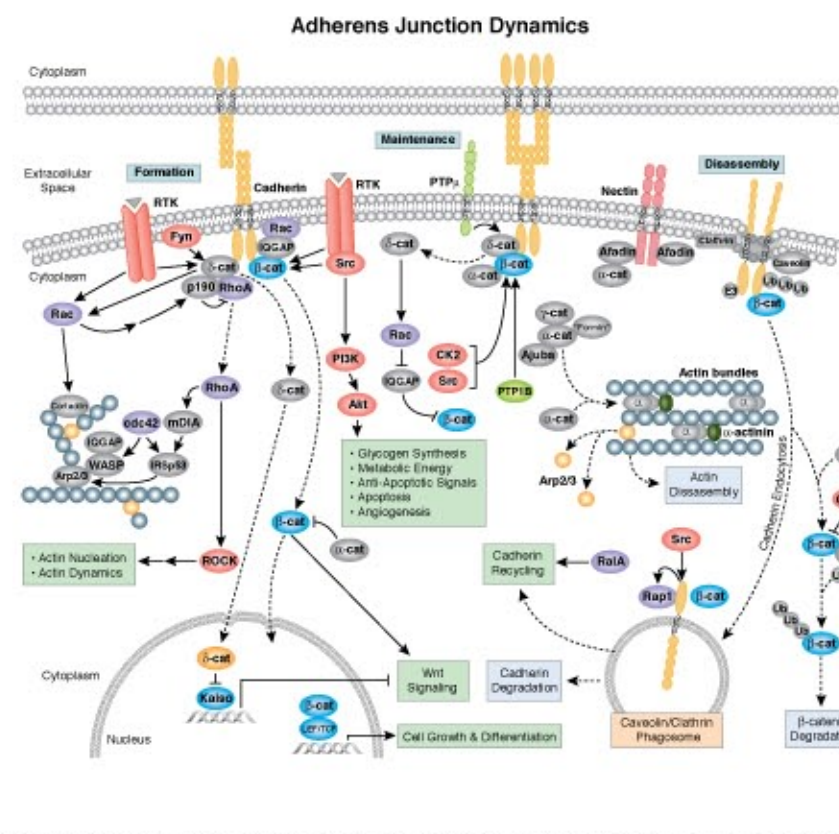
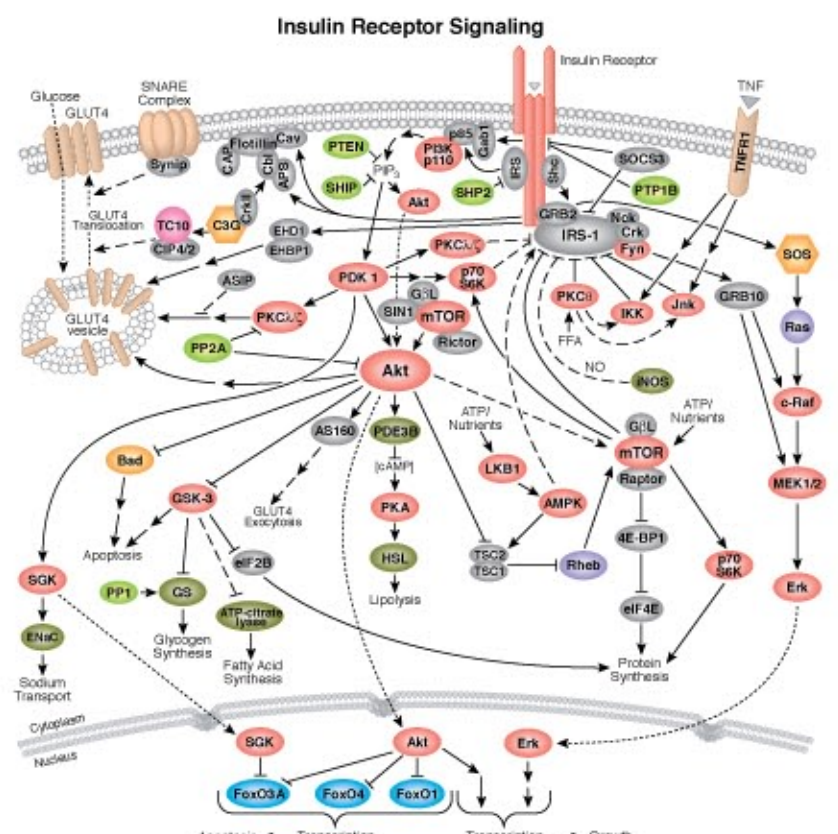
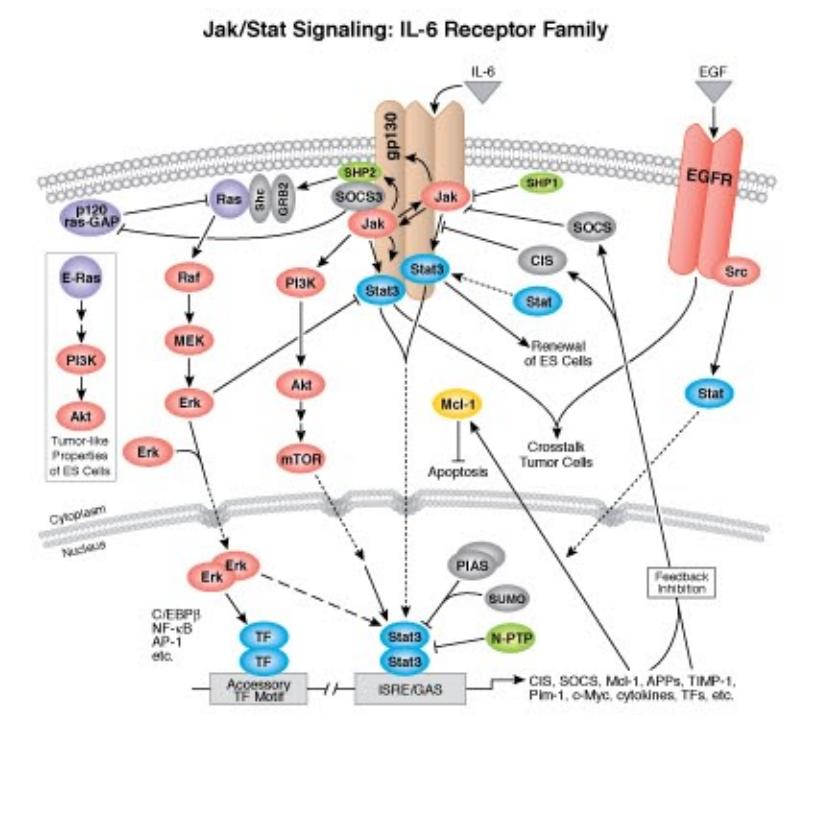
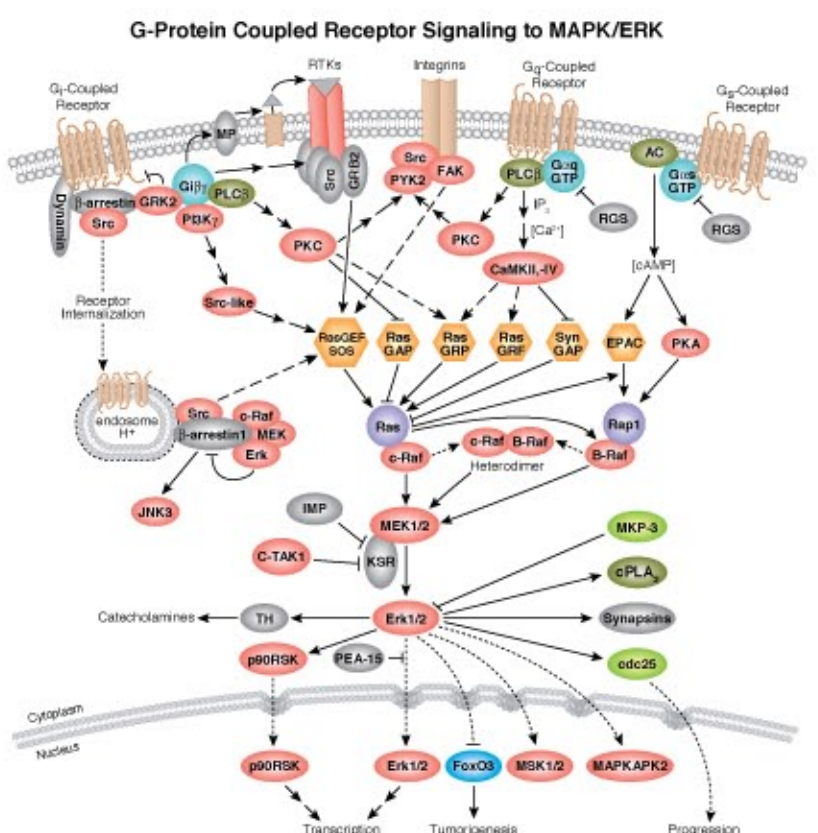
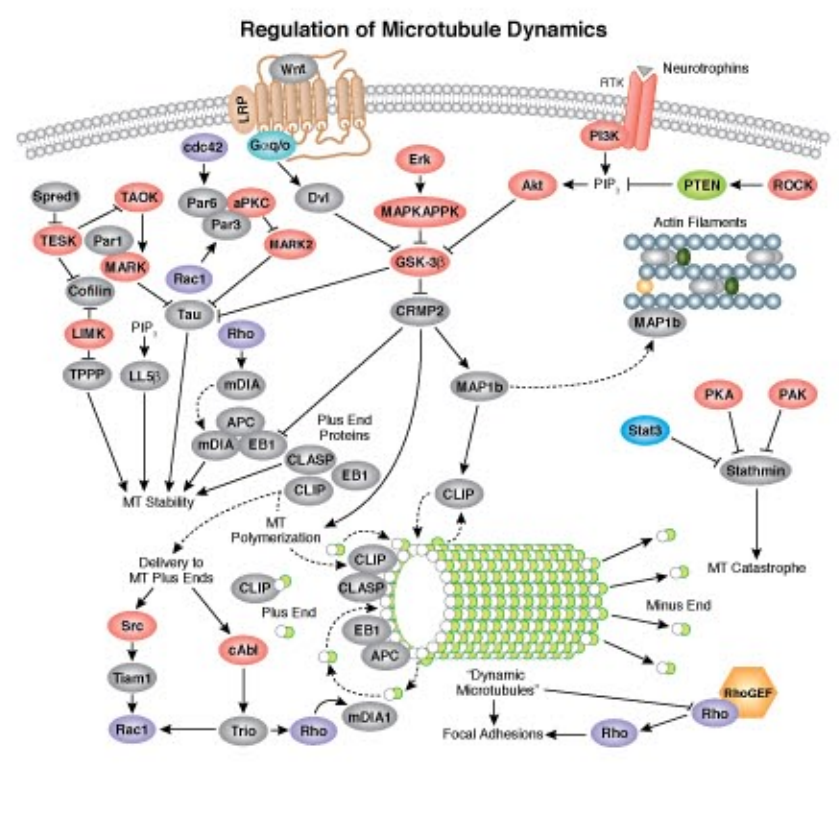
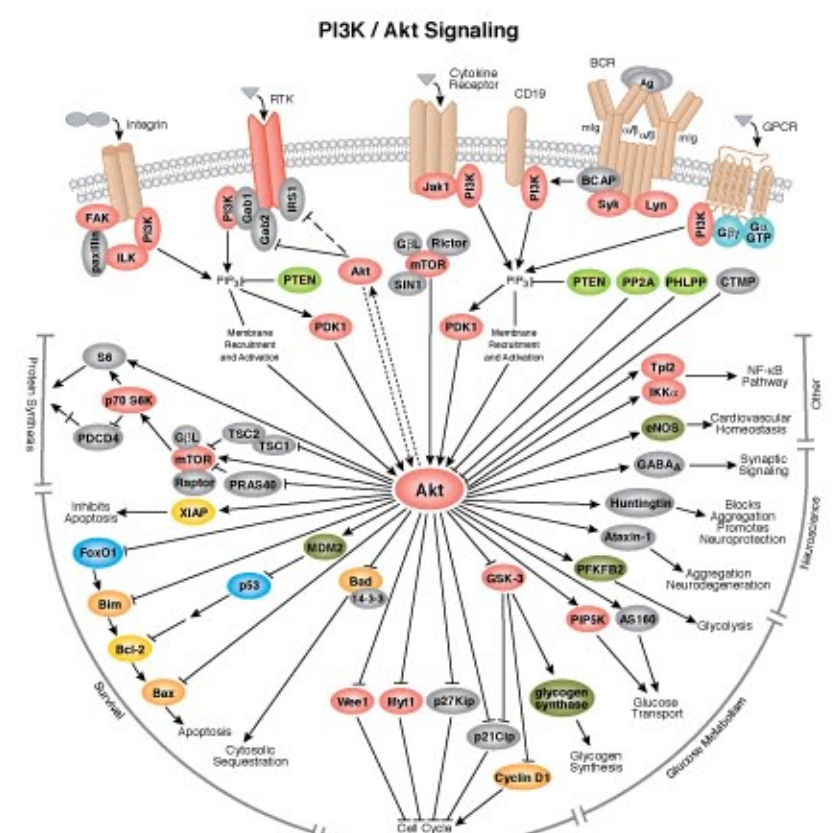
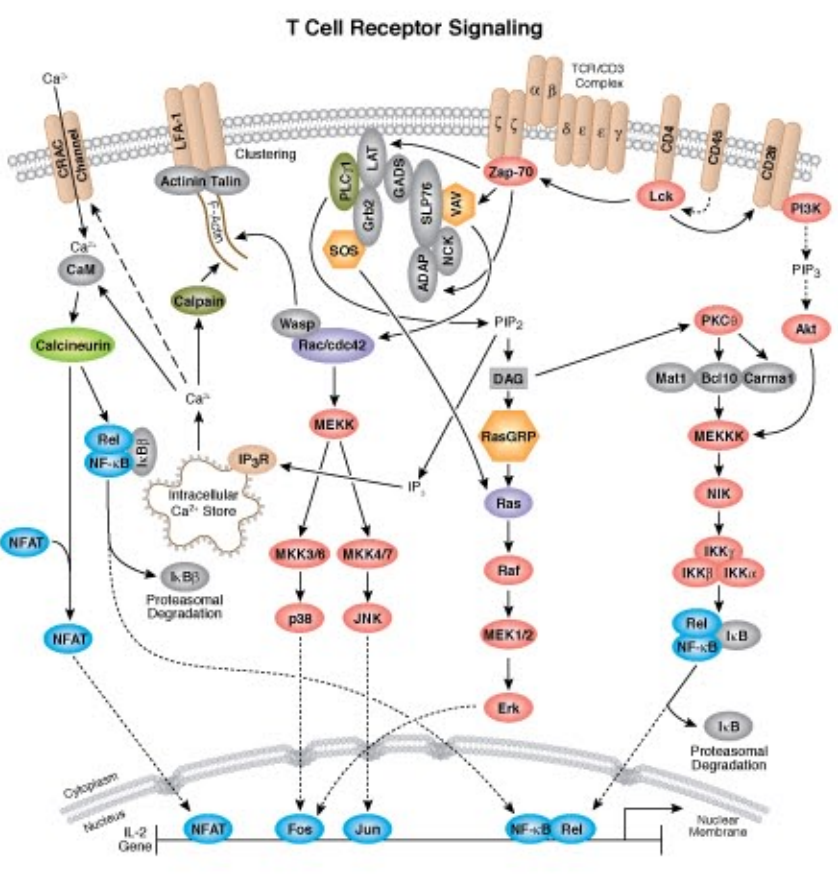
ADHESION AND MOBILITY



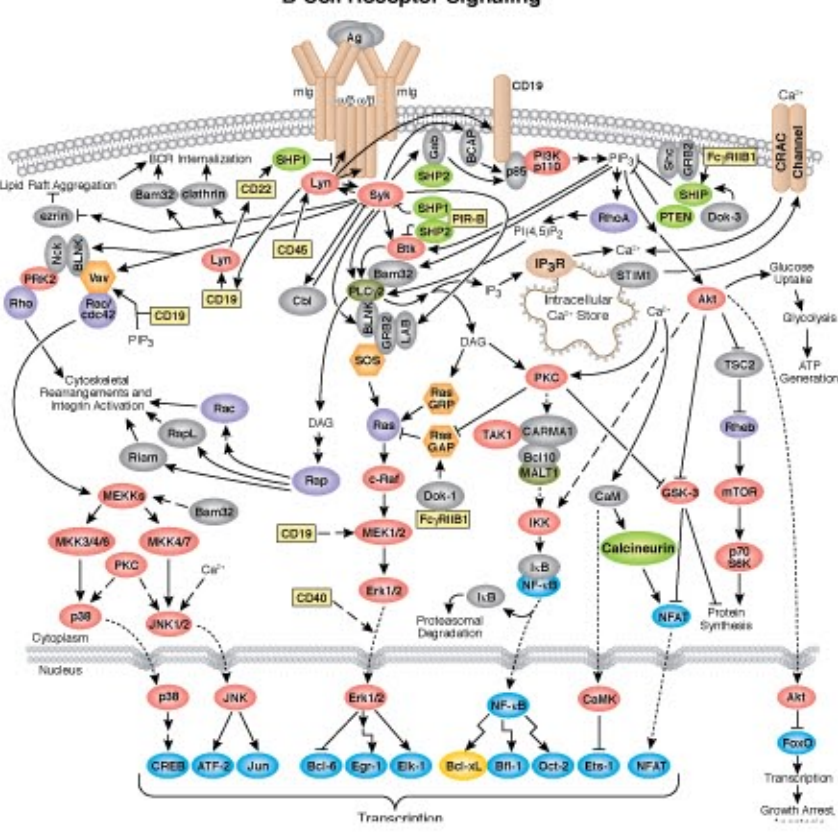
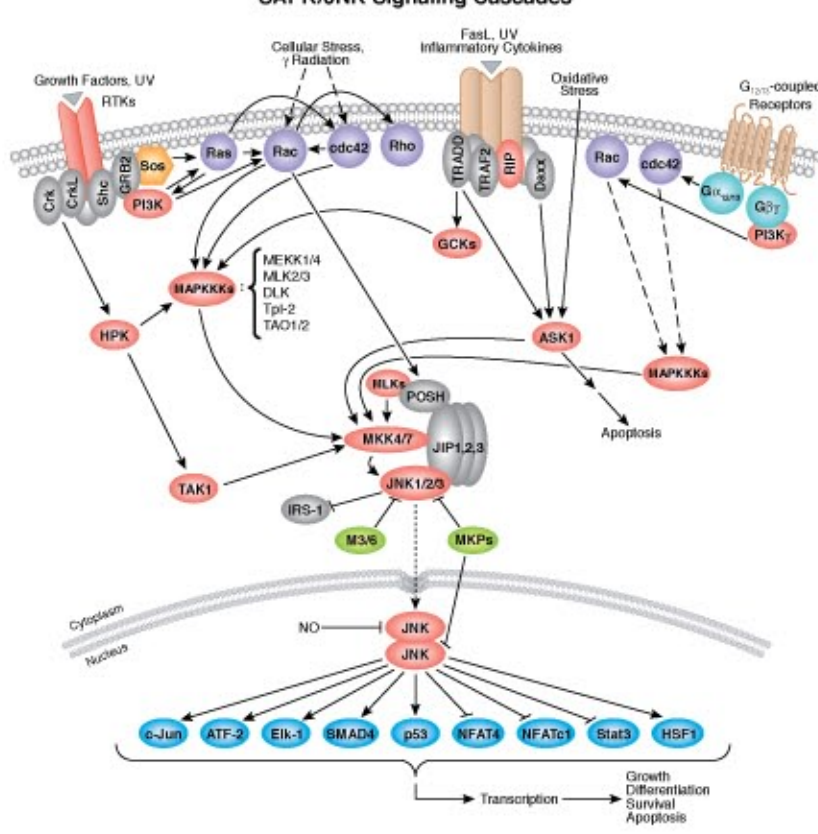
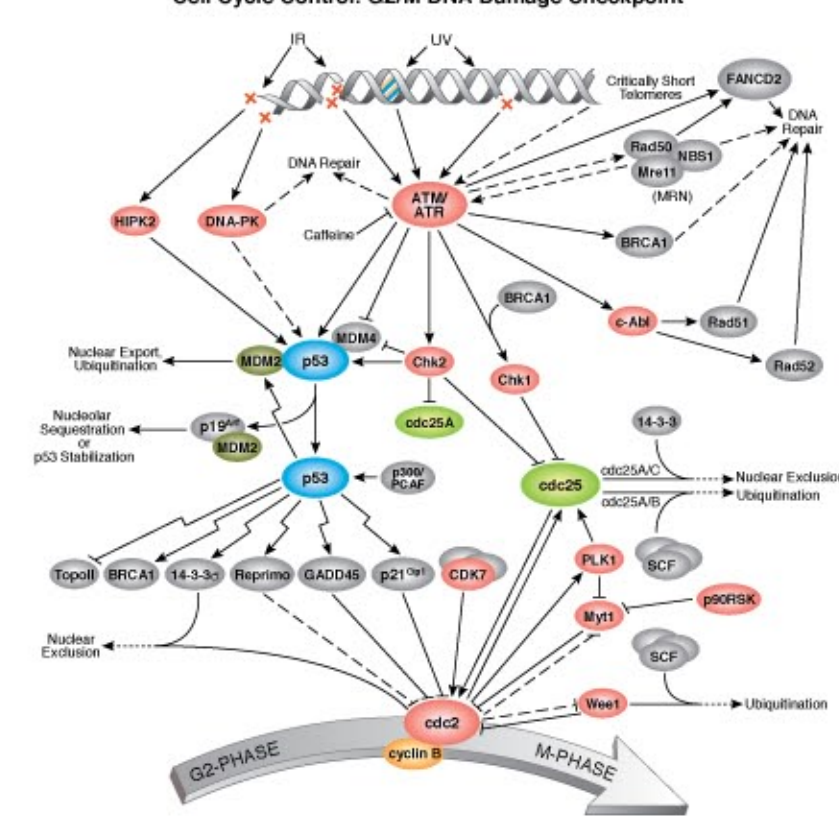
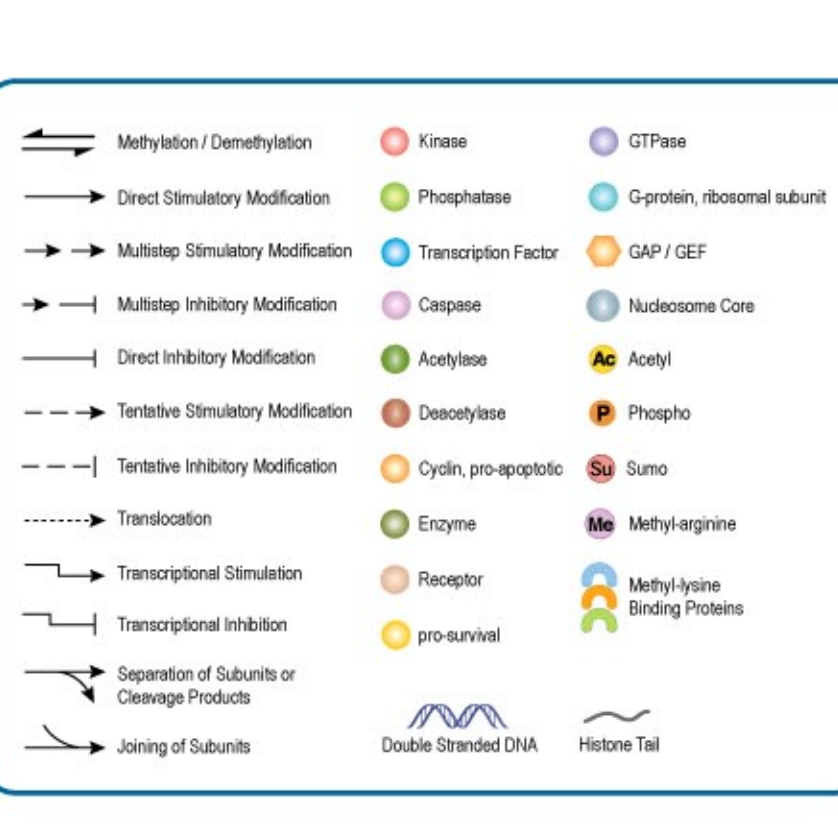
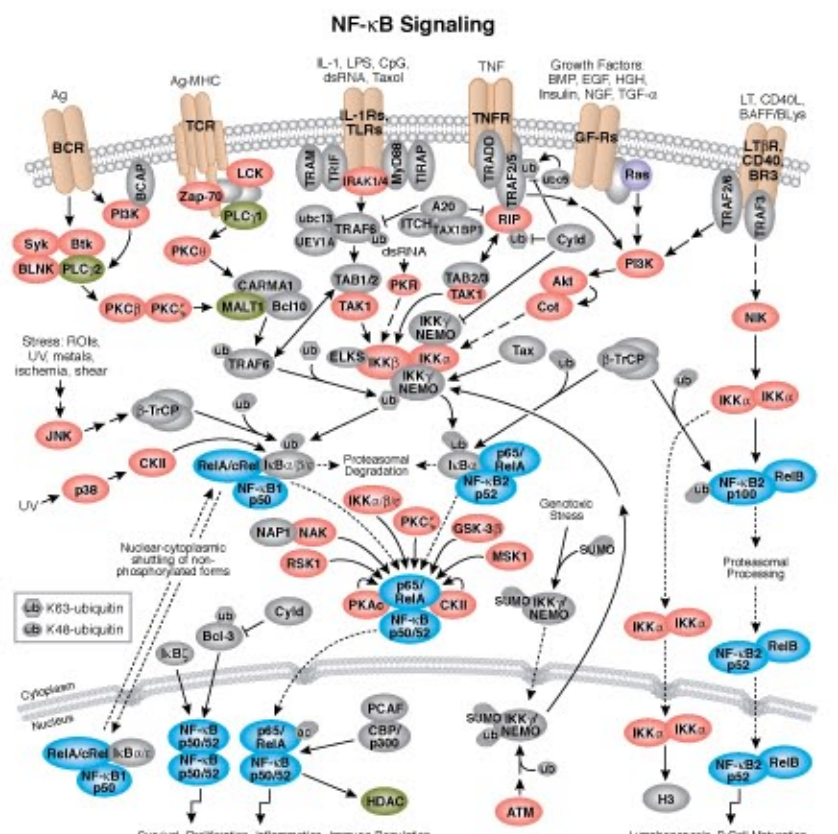
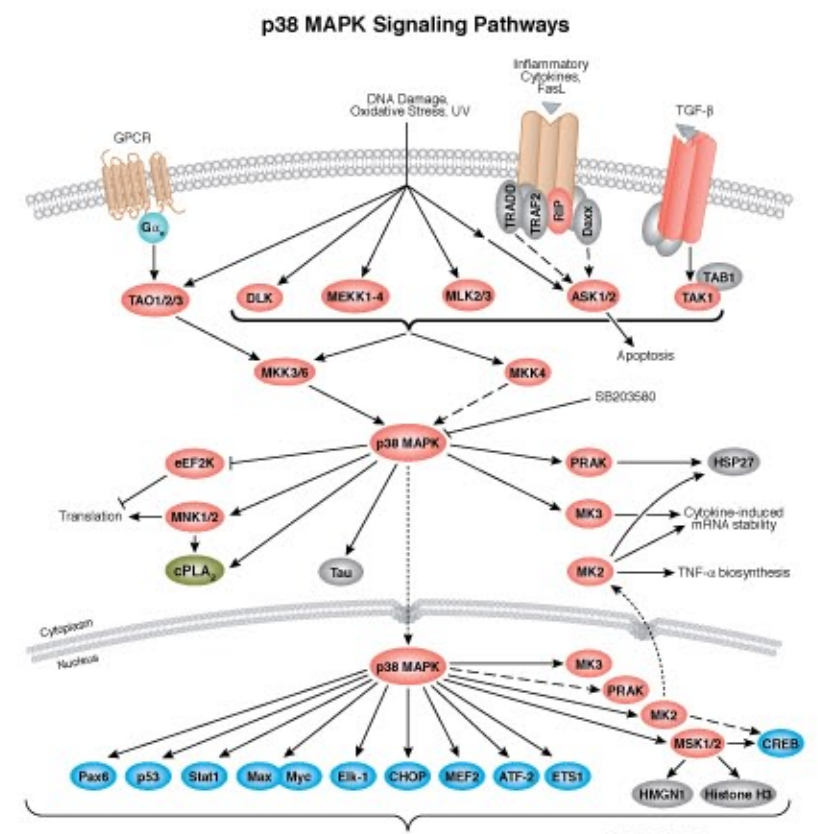
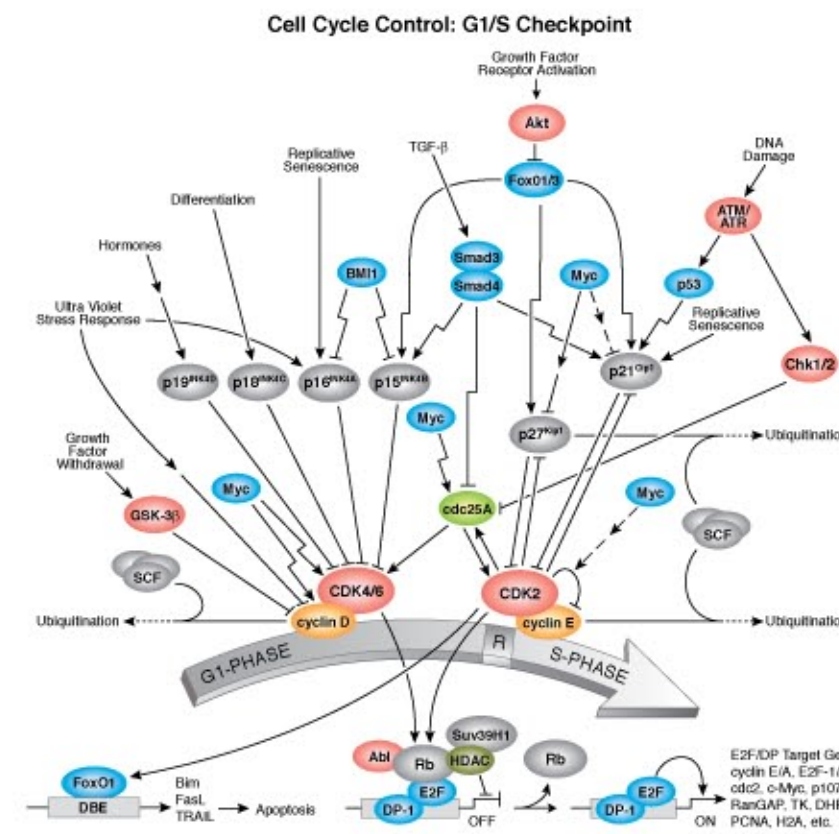
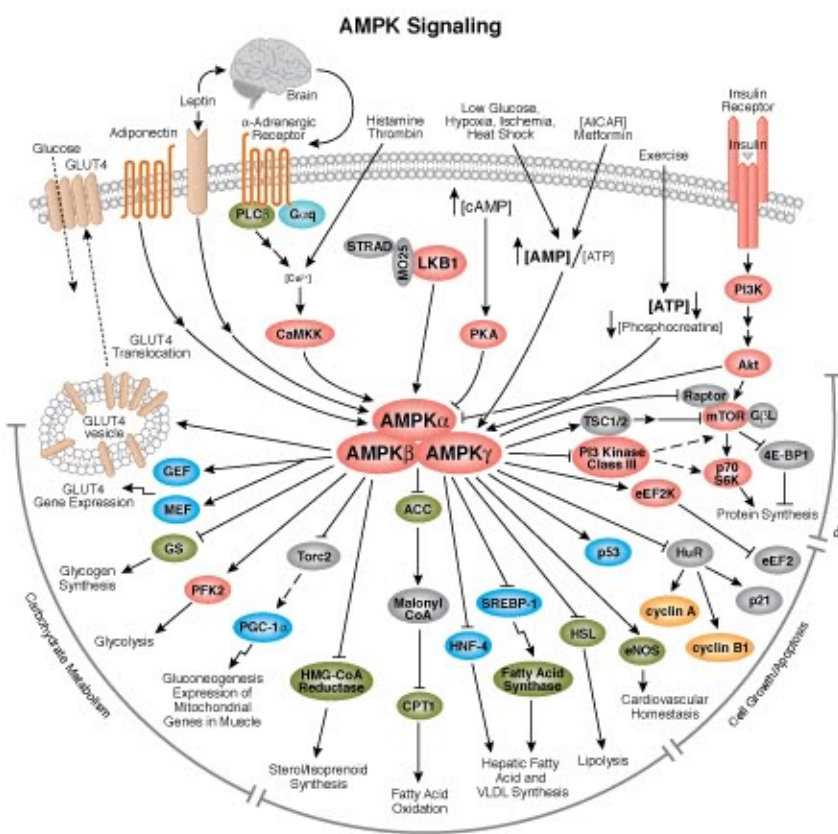
MITOGEN ACTIVATED PROTEIN KINASE



IMMUNOLOGY



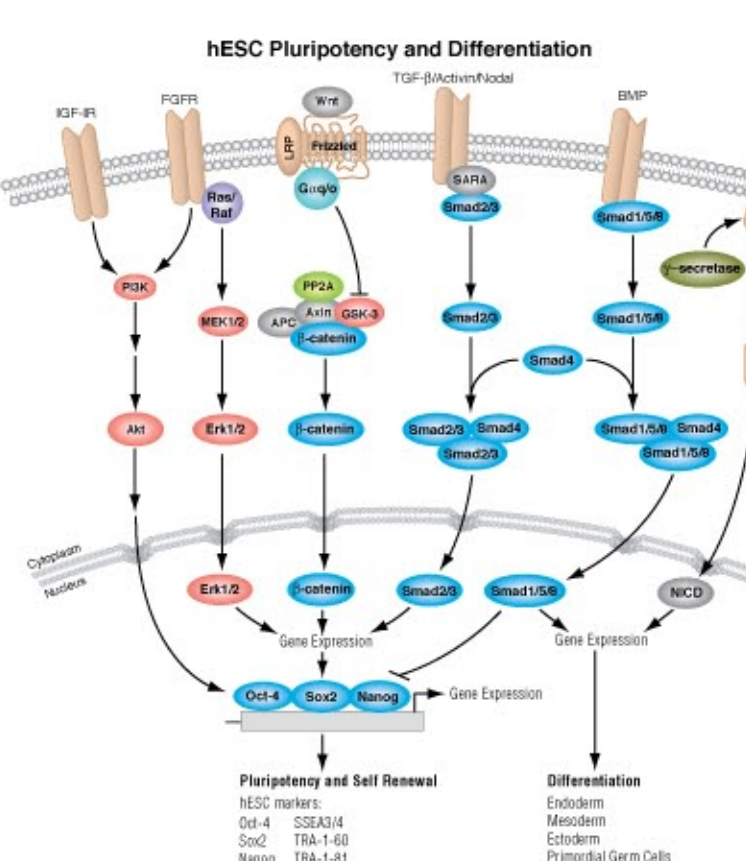
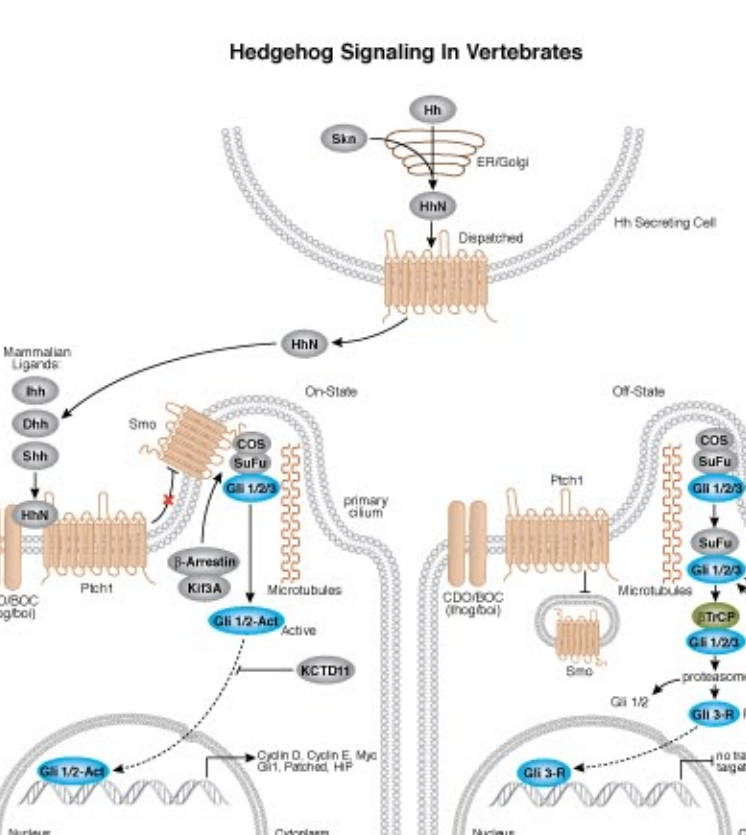
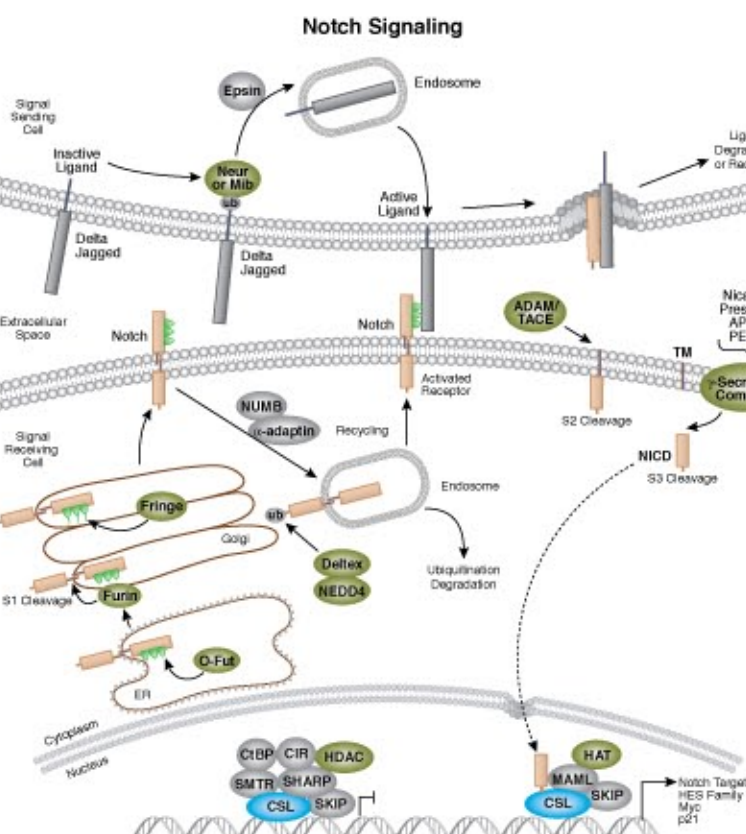
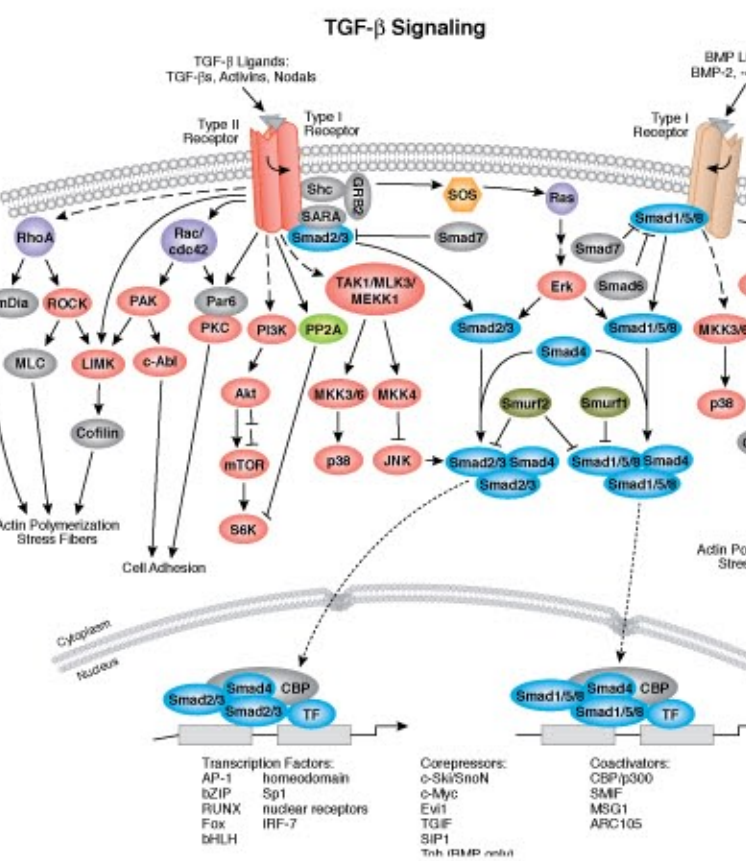
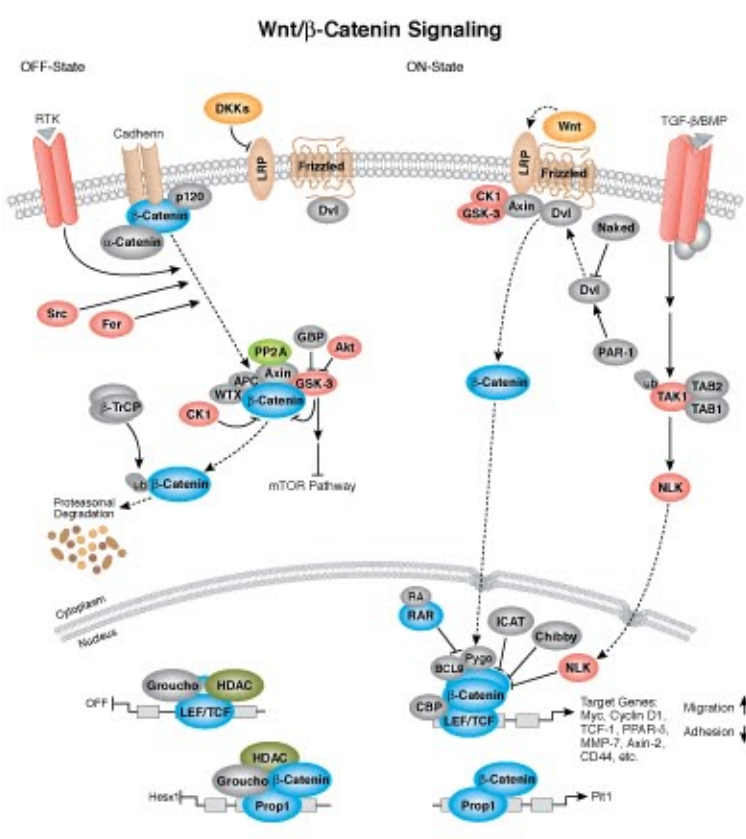
CELL CYCLE/DNA DAMAGE



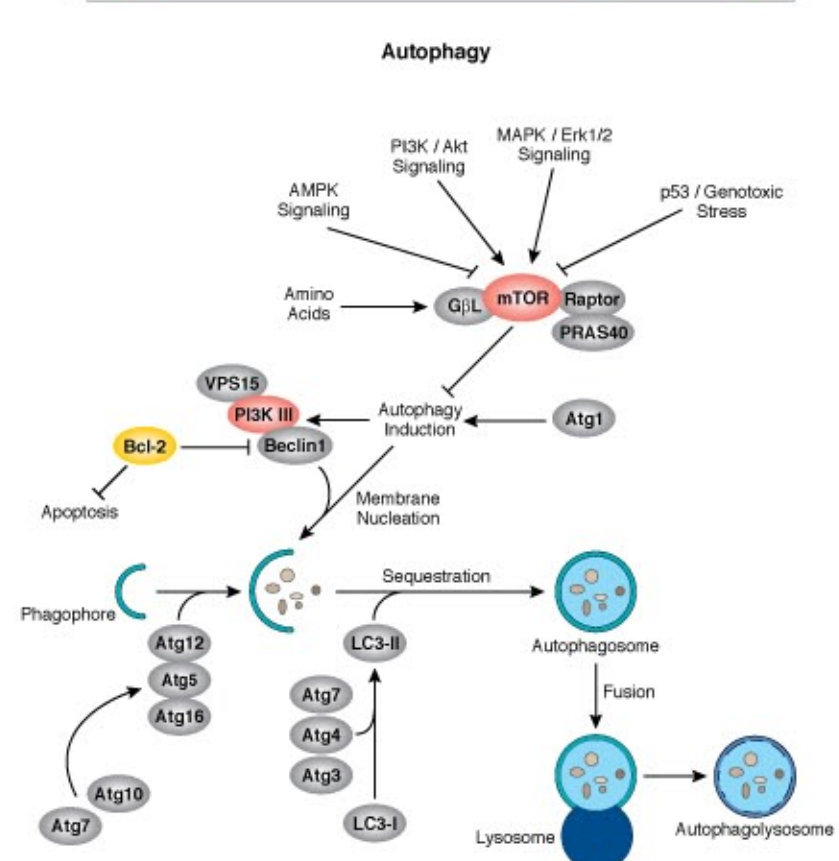
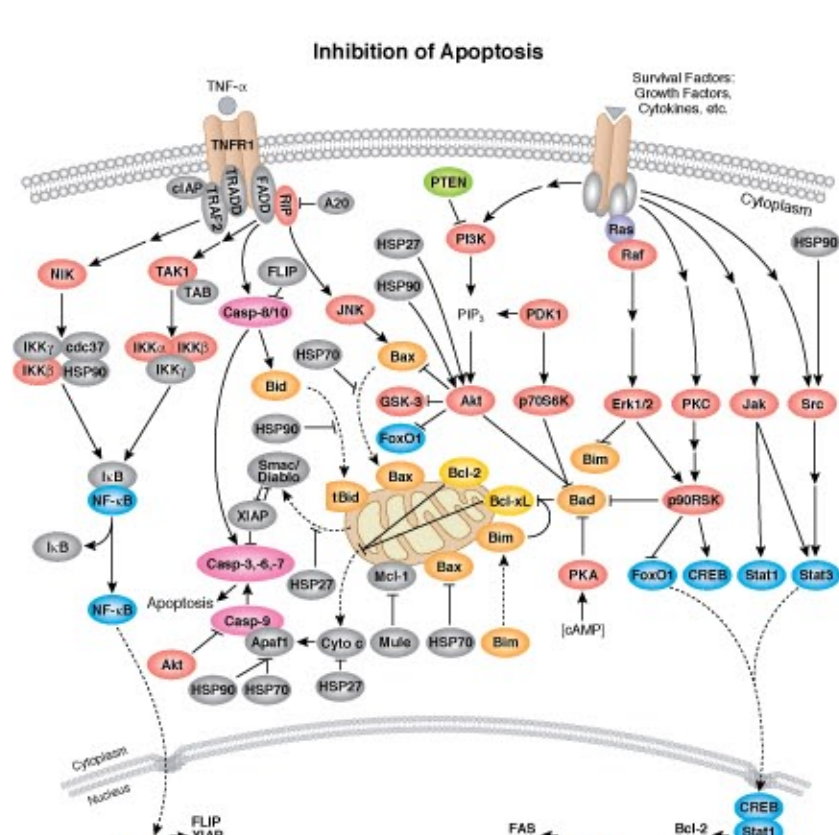
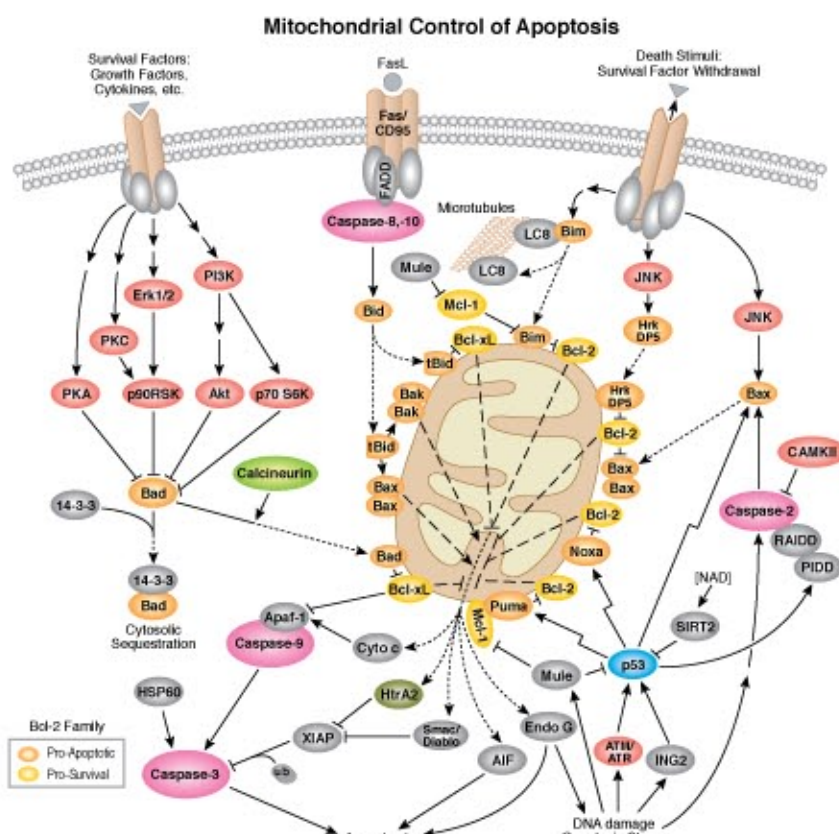
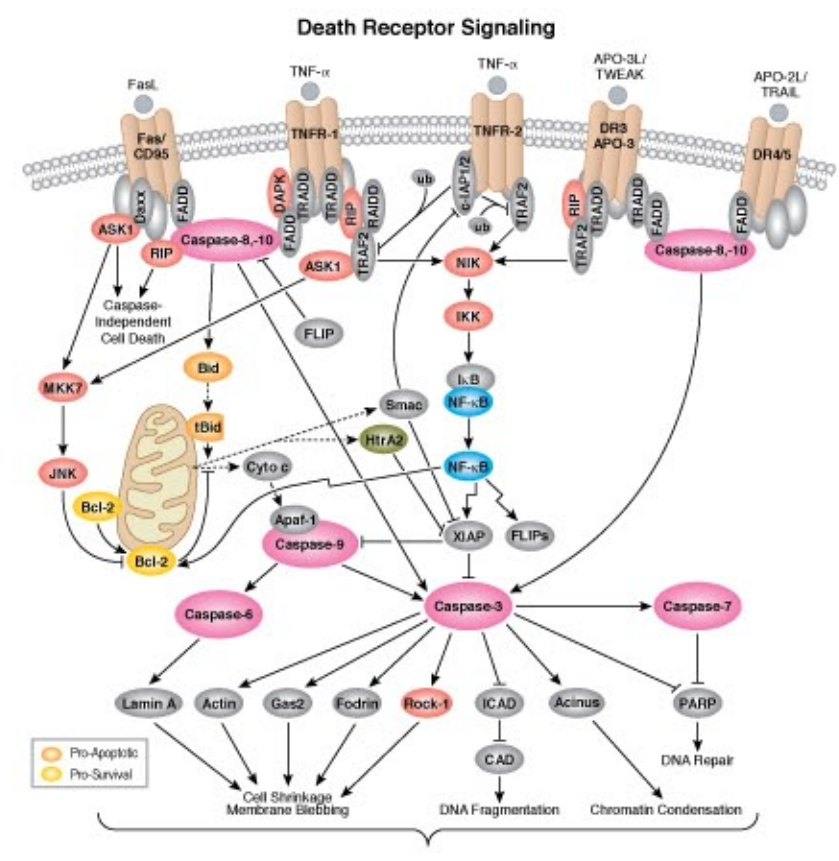
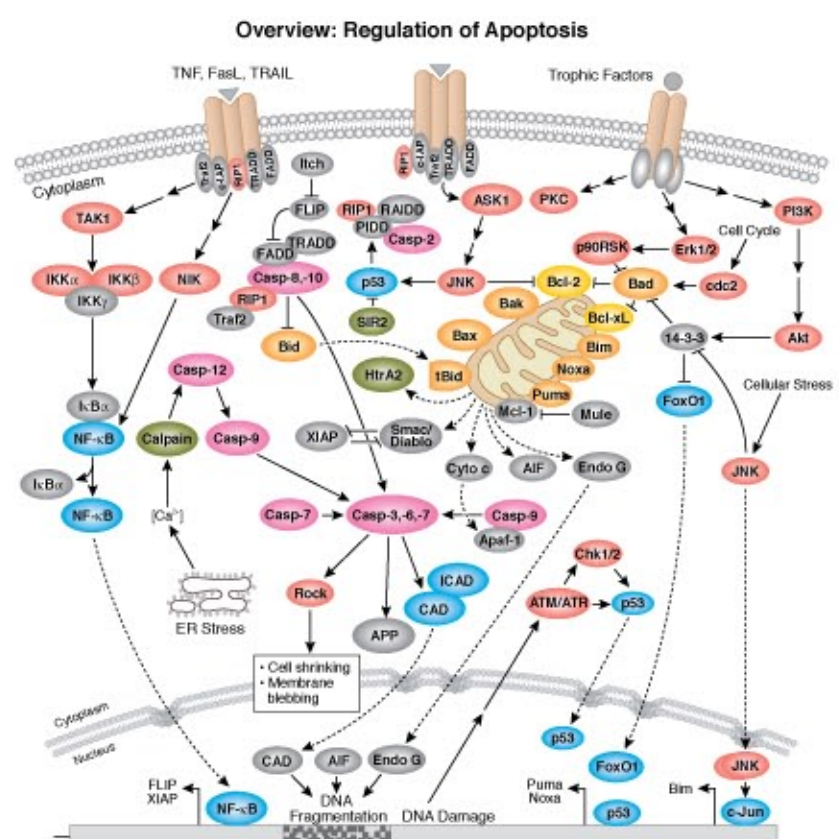
Signal Transduction Pathways



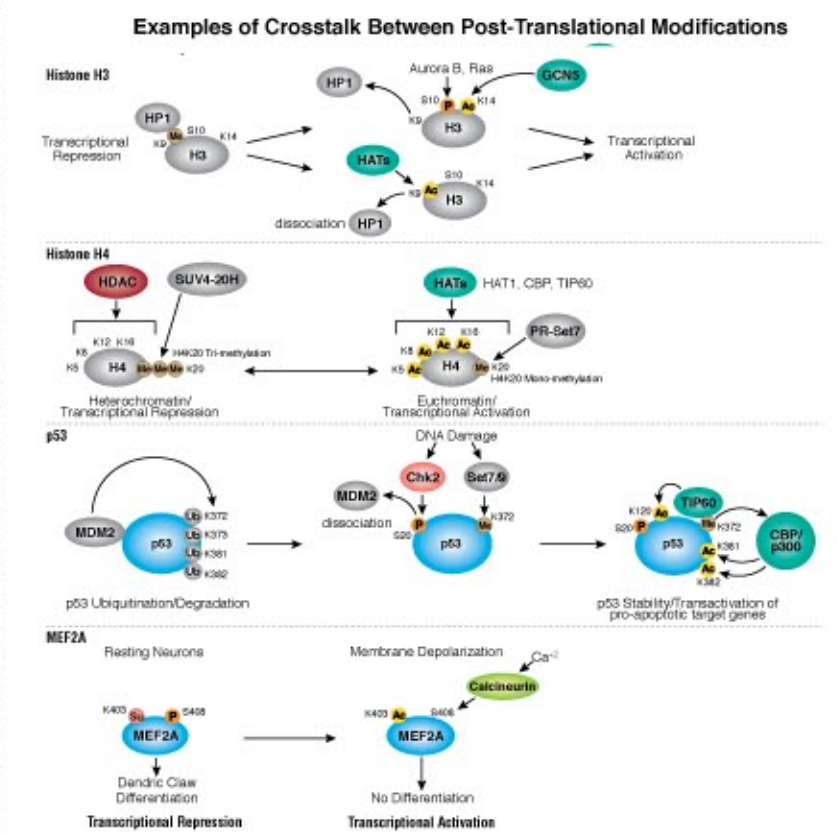
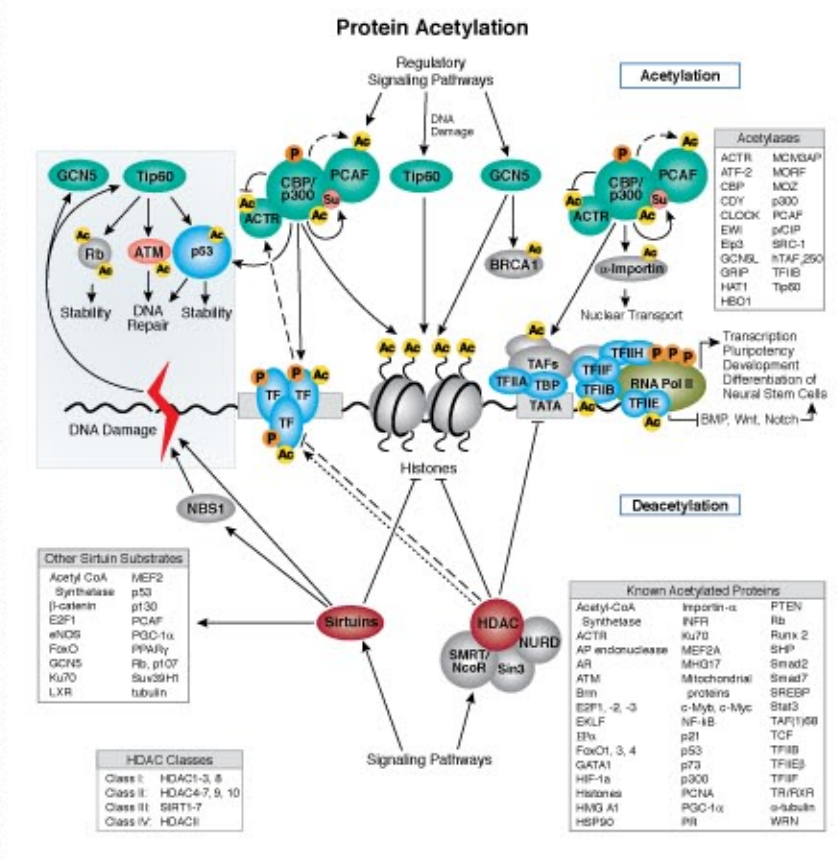
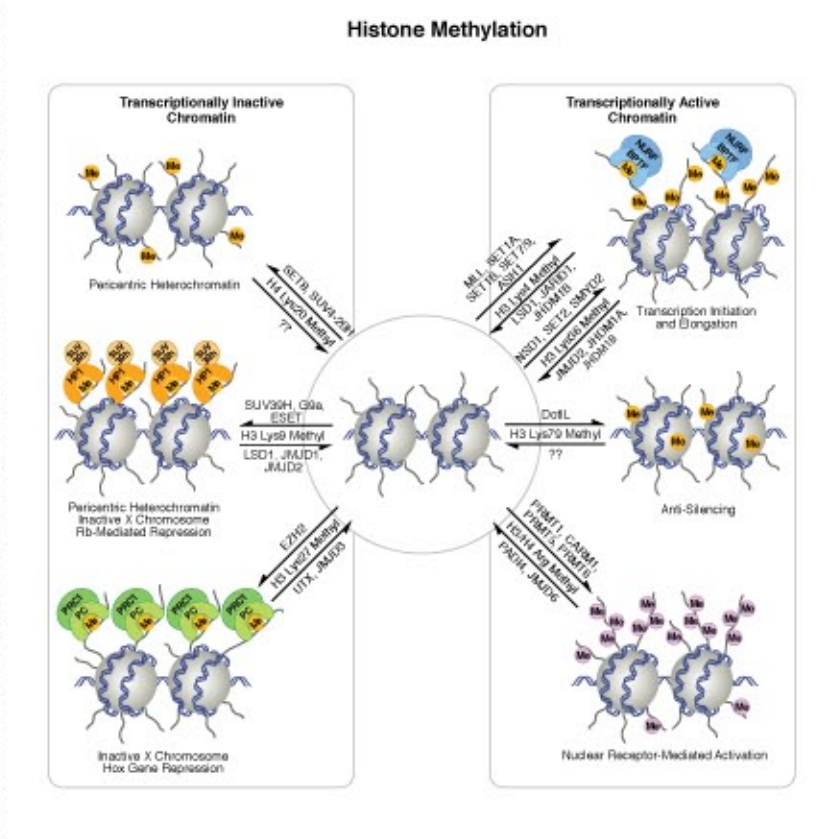
DEVELOPMENT



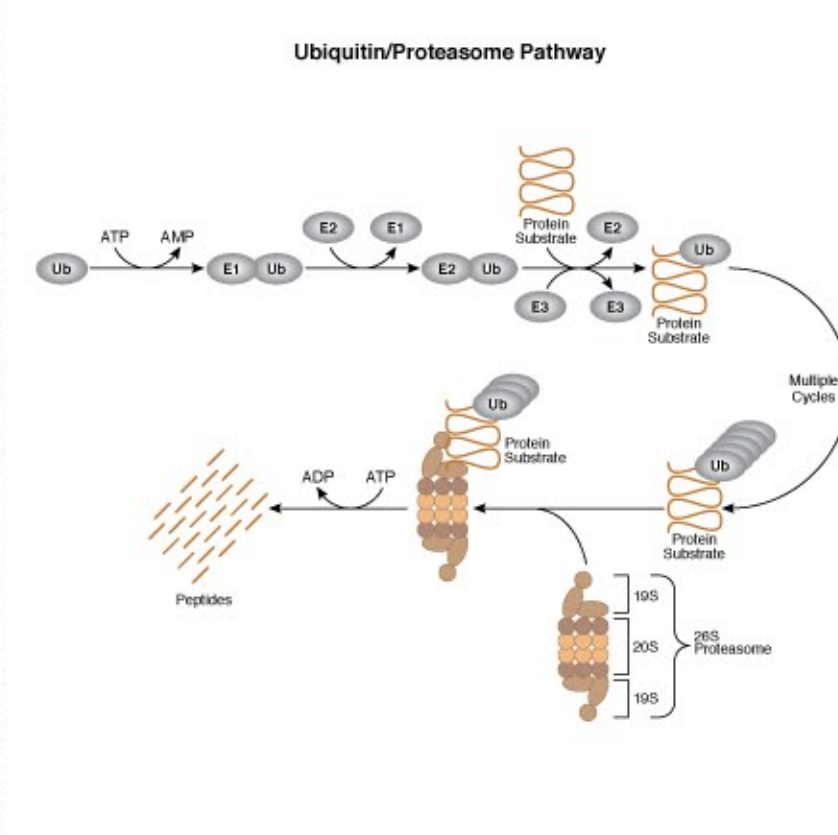
APOPTOSIS AND AUTOPHAGY



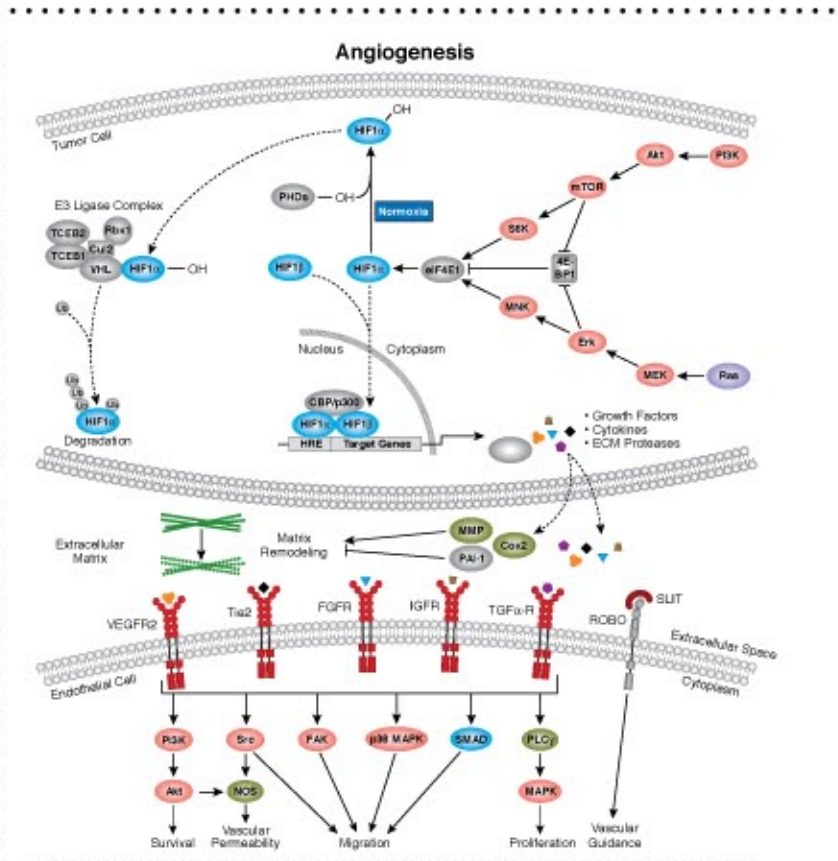
CHROMATIN MODIFICATION



UBIQUITIN/PROTEASOME



ANGIOGENESIS



TRANSLATIONAL

