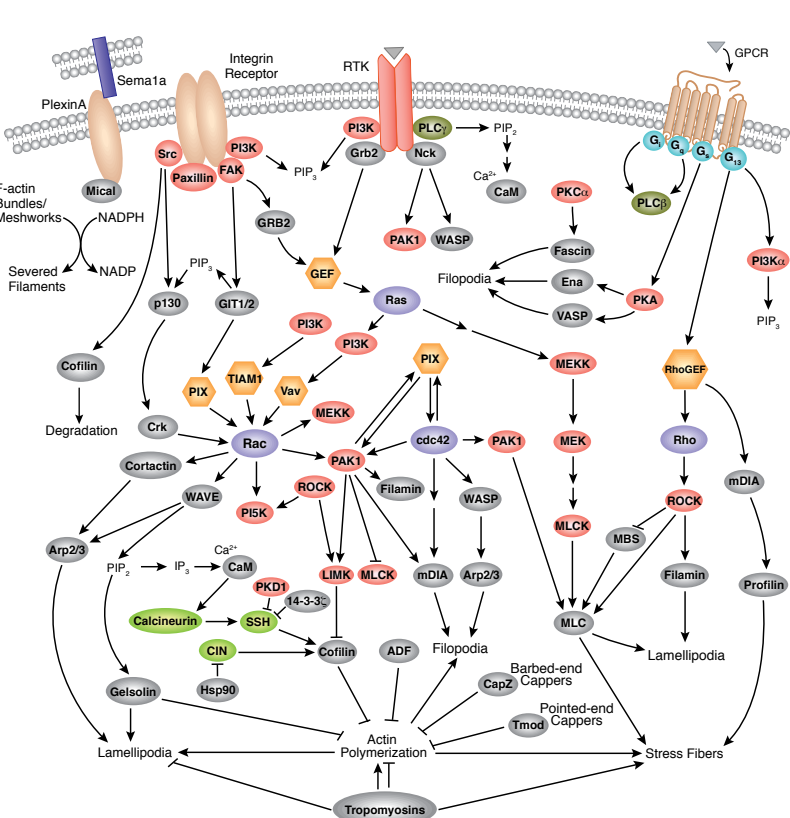
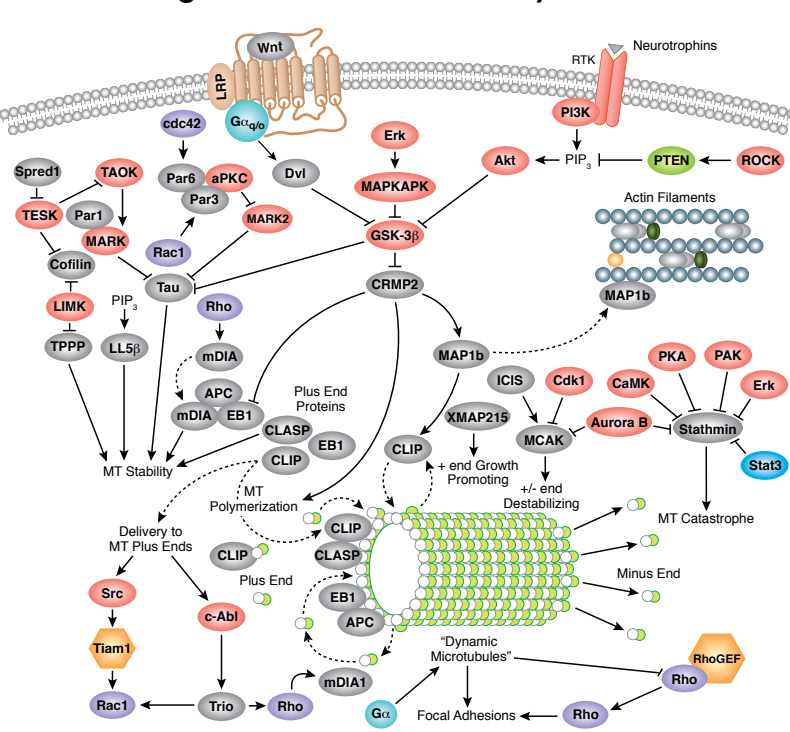


CYTOSKELETAL REGULATION AND ADHESION

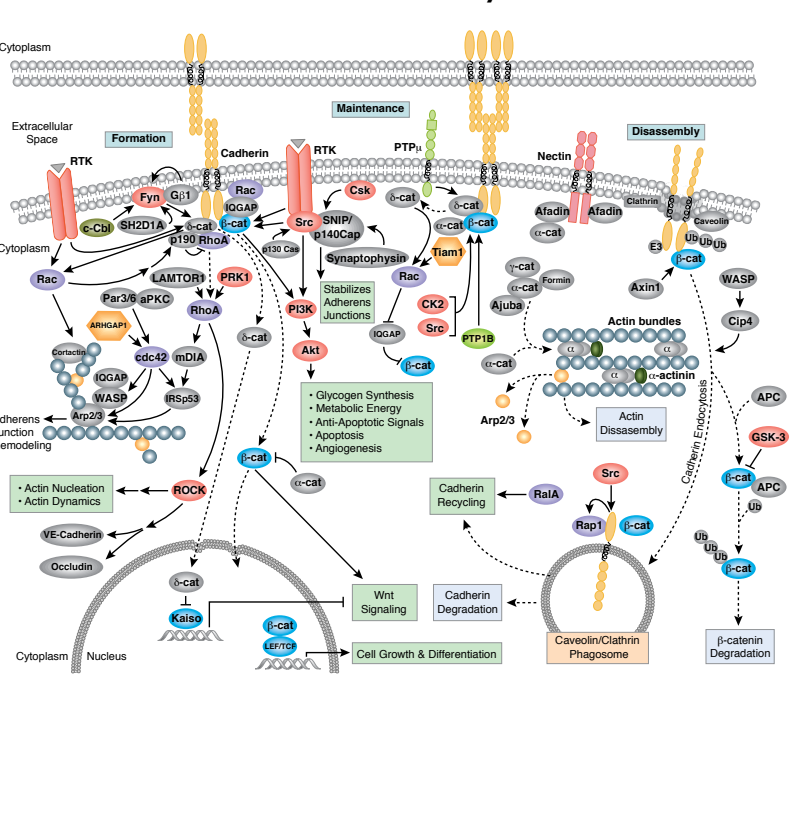
Regulation of Actin Dynamics



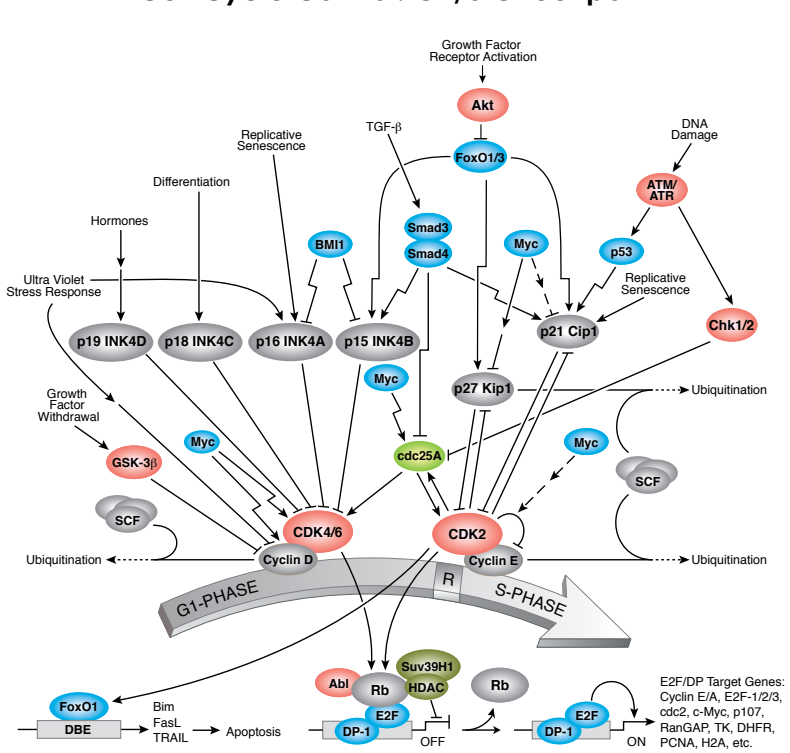
Regulation of Microtubule Dynamics



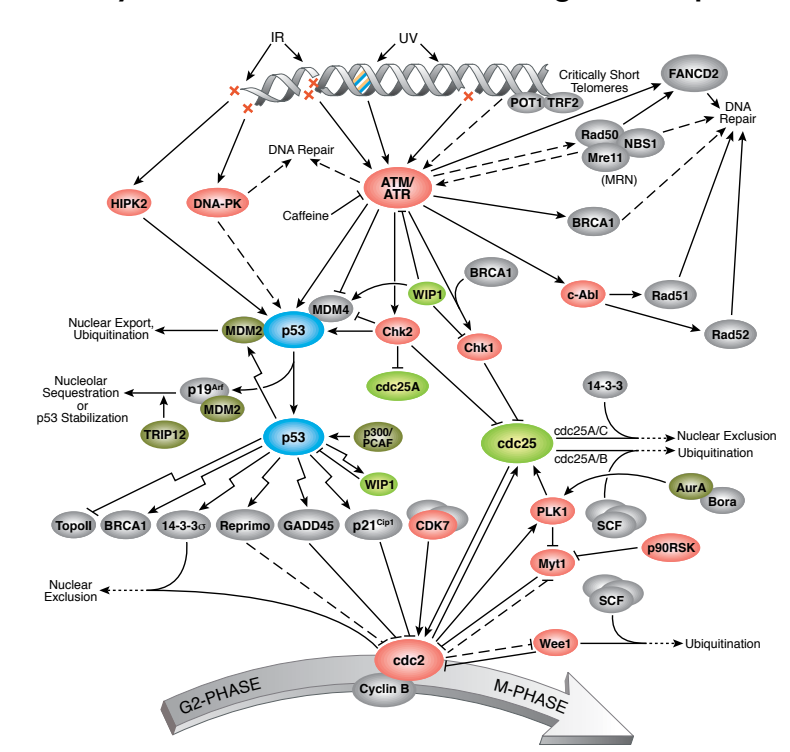
Adherens Junction Dynamics



Cell Cycle Control: G1/S Checkpoint

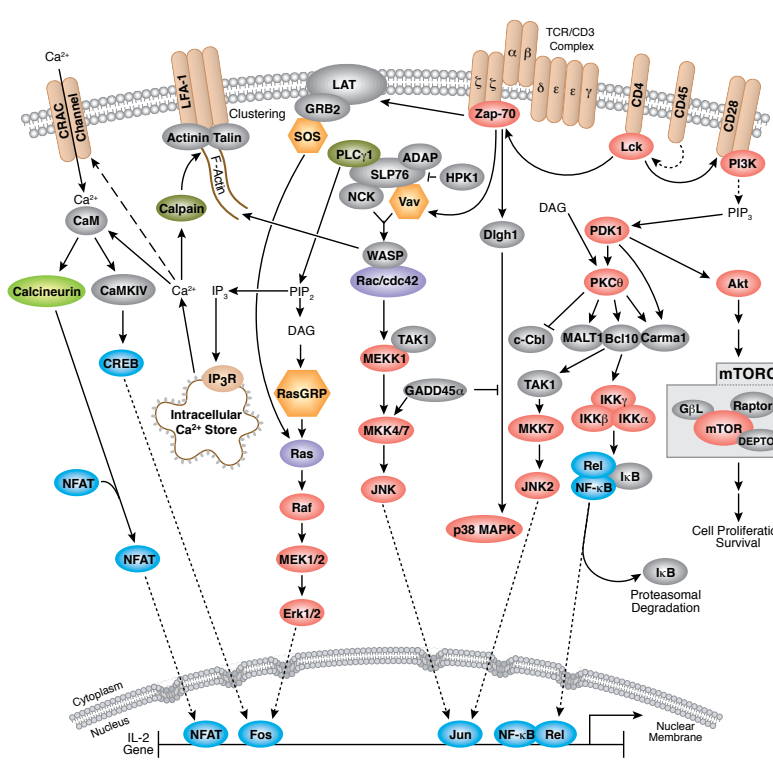


Cell Cycle Control: G2/M DNA Damage Checkpoint

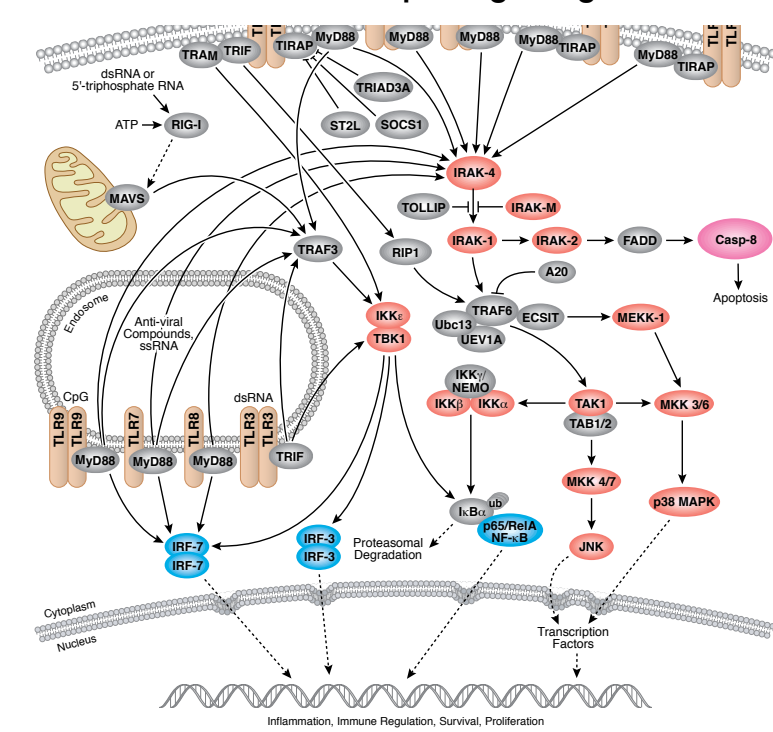


IMMUNOLOGY AND INFLAMMATION

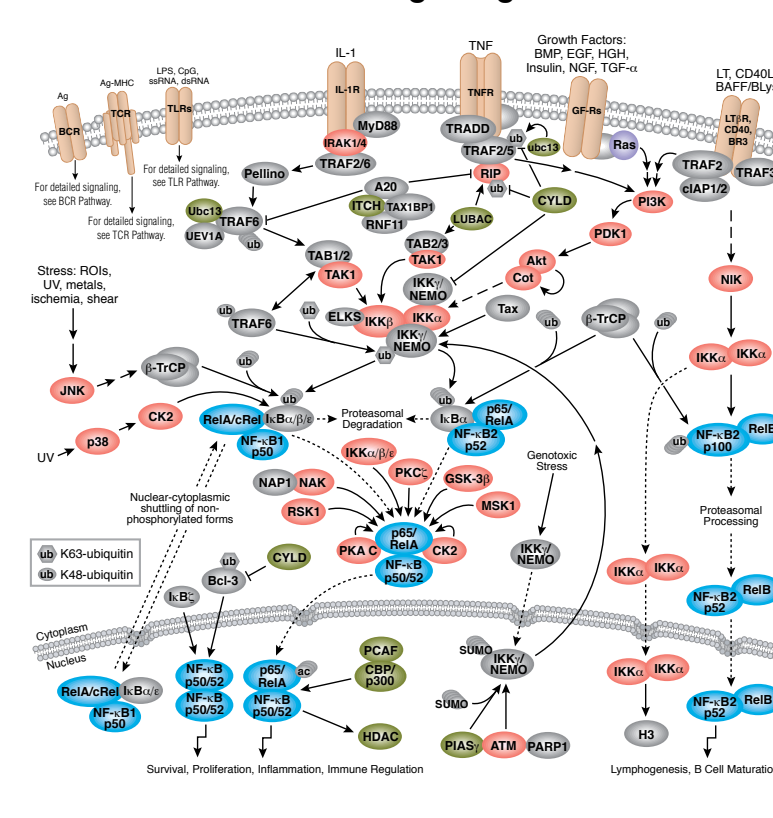
T Cell Receptor Signaling



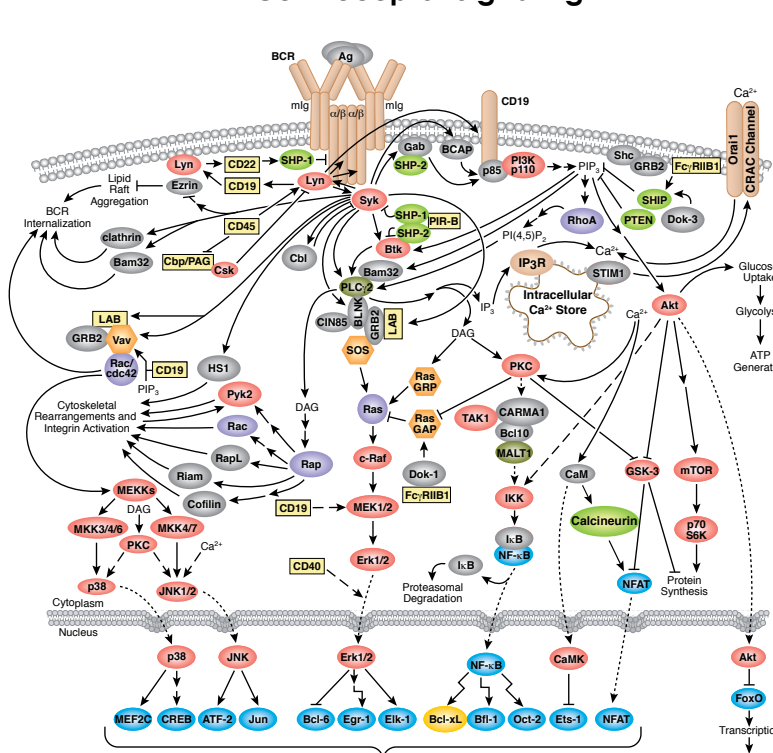
Toll-like Receptor Signaling



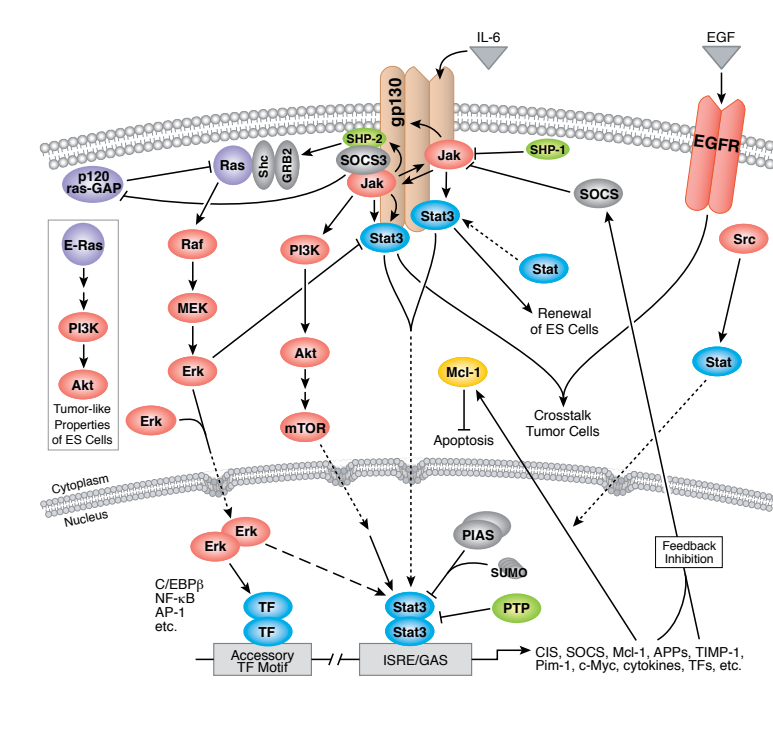
NF-κB Signaling



B Cell Receptor Signaling

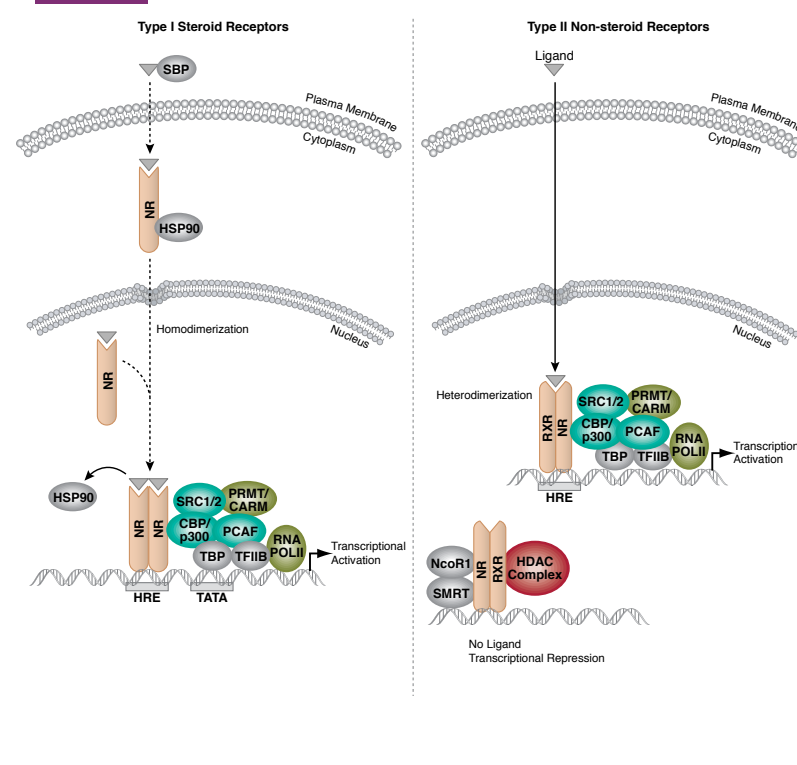


Jak/Stat Signaling: IL-6 Receptor Family

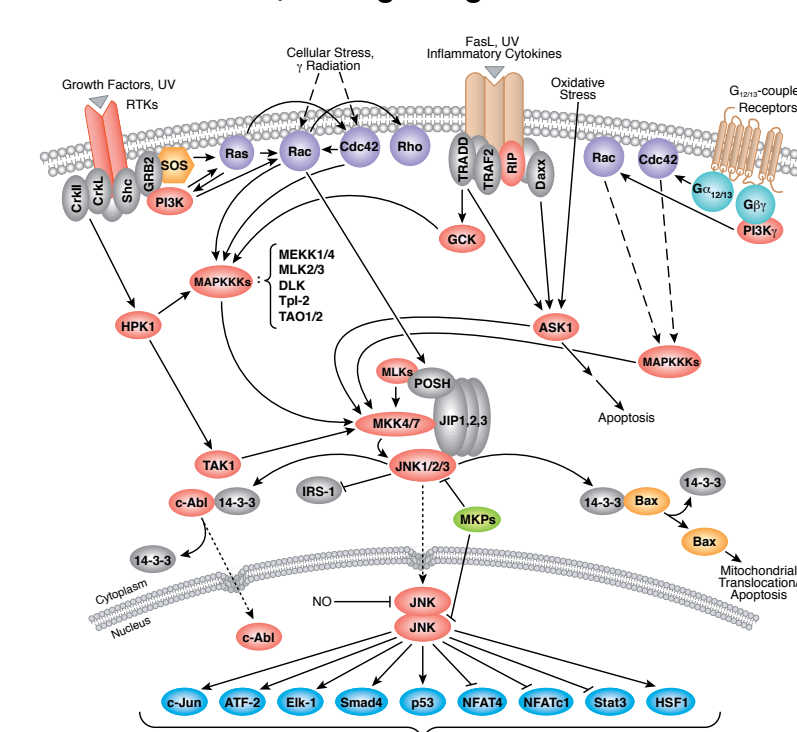


MAP KINASE SIGNALING

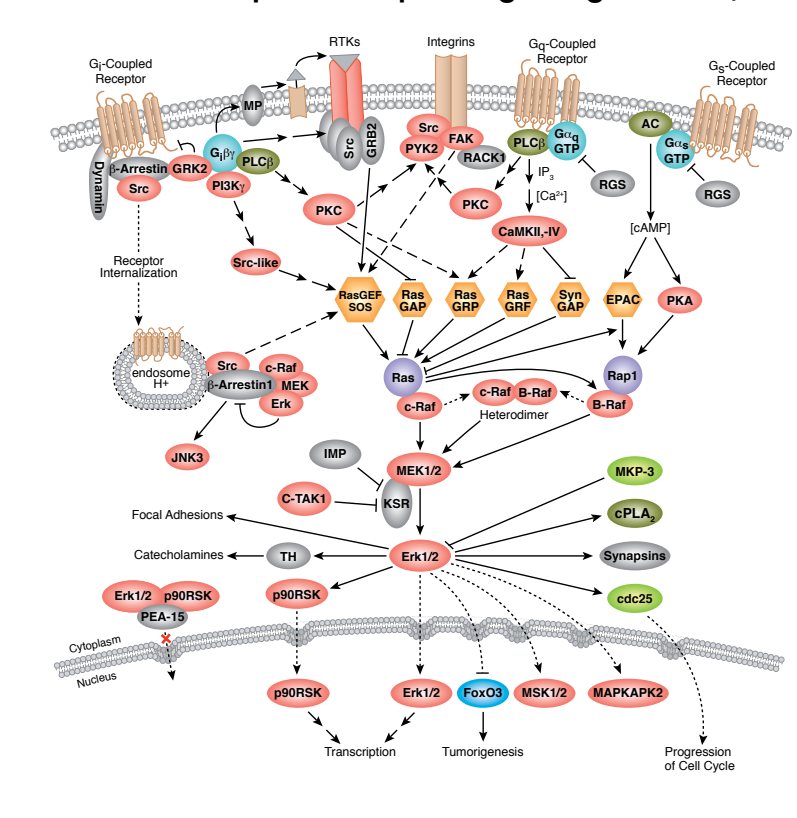
Nuclear Receptor Signaling



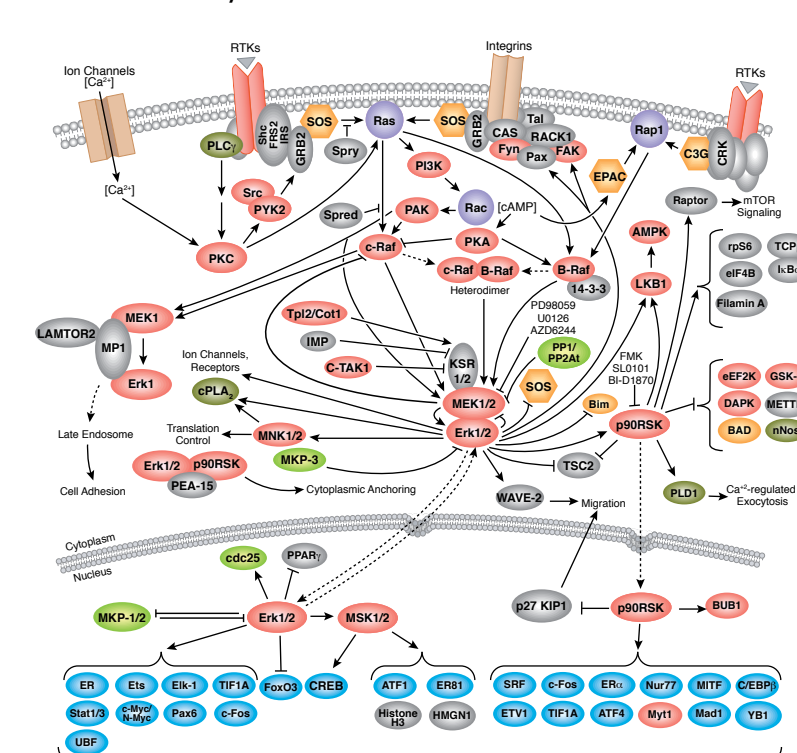
SAPK/JNK Signaling Cascades



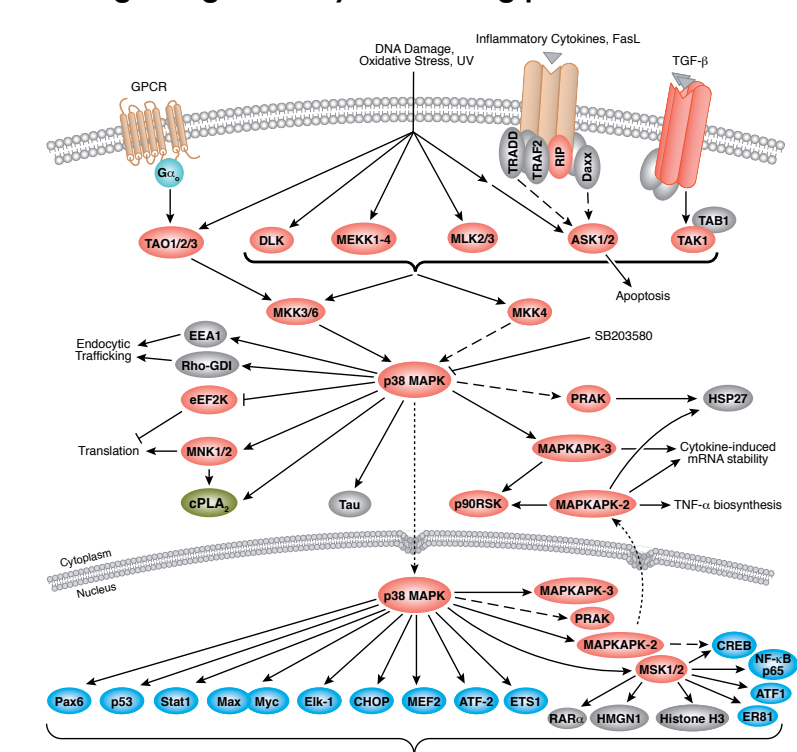
G Protein-coupled Receptors Signaling to MAPK/Erk



MAPK/Erk in Growth and Differentiation

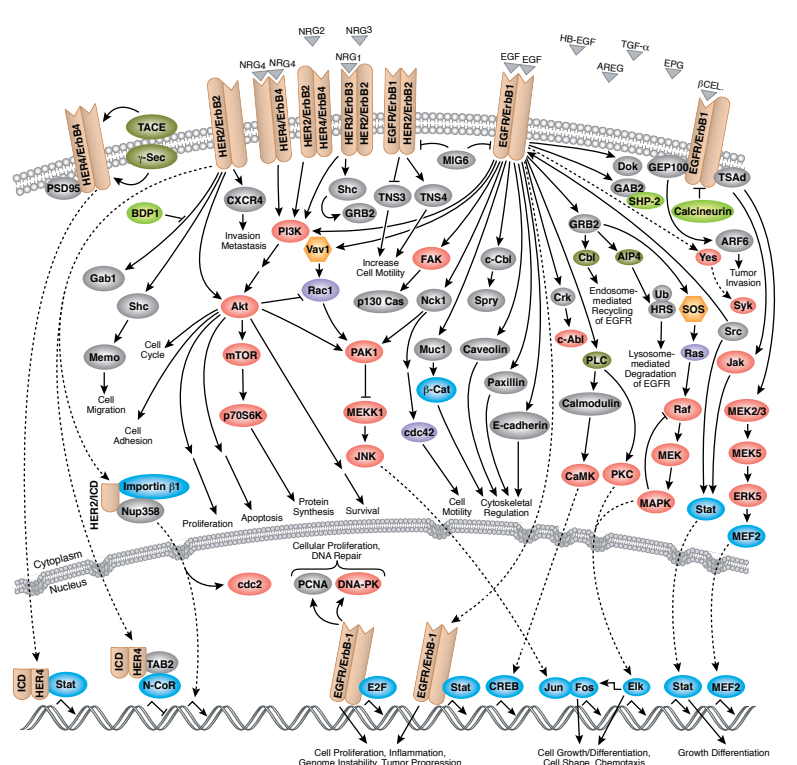


Signaling Pathways Activating p38 MAP Kinase



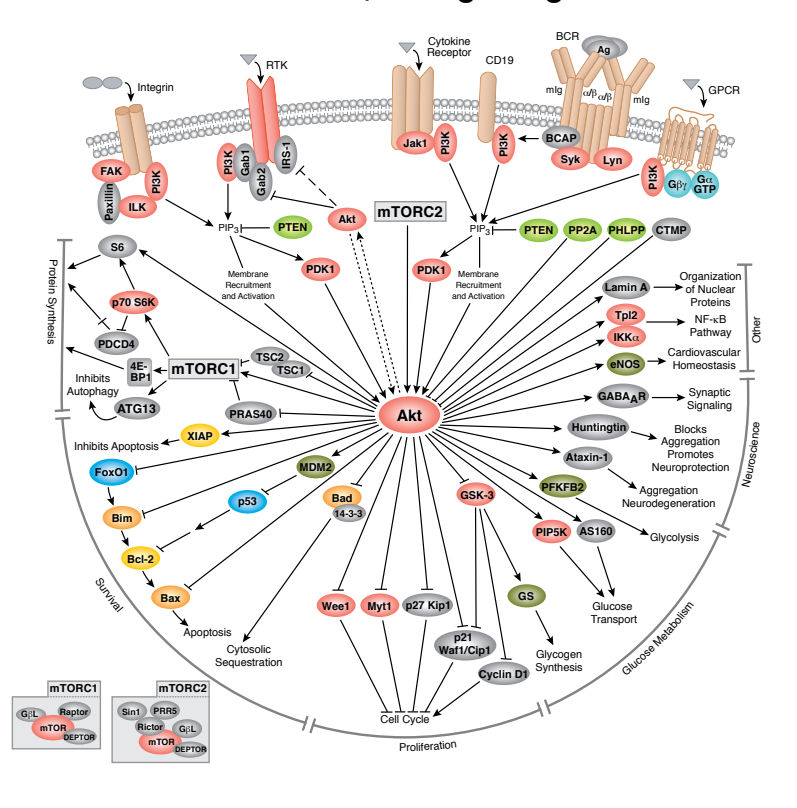
EGF RECEPTOR SIGNALING

Erb/HER Signaling



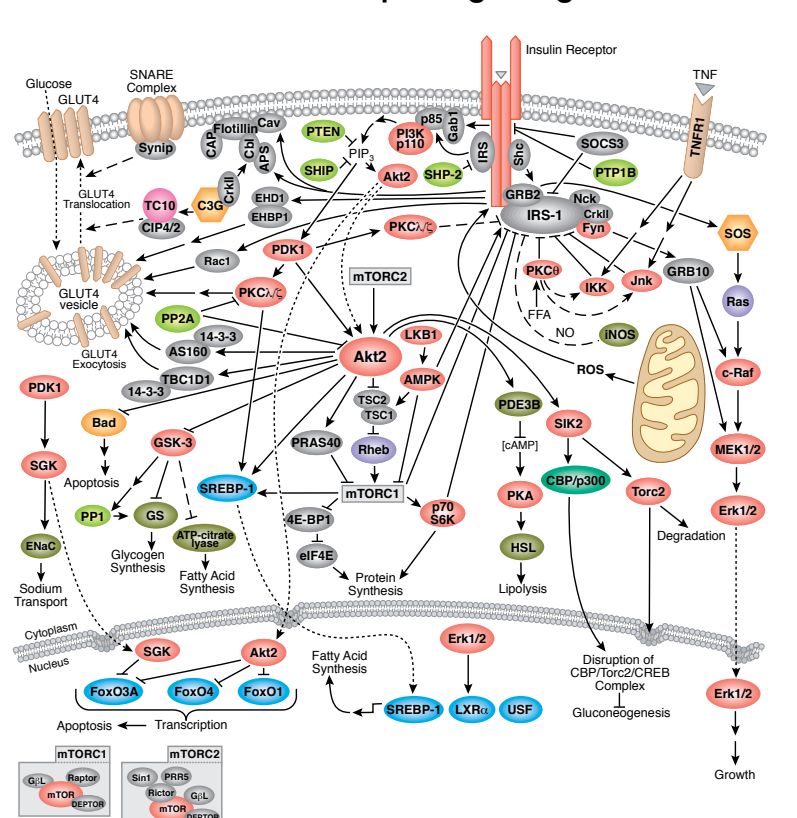
PI3 KINASE/AKT SIGNALING

PI3 Kinase/Akt Signaling

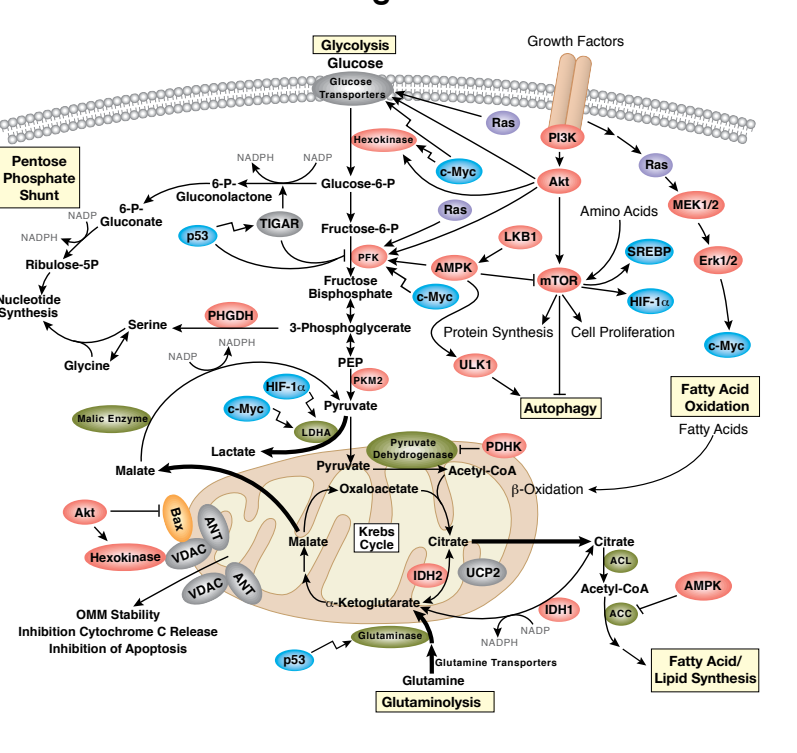


CELLULAR METABOLISM

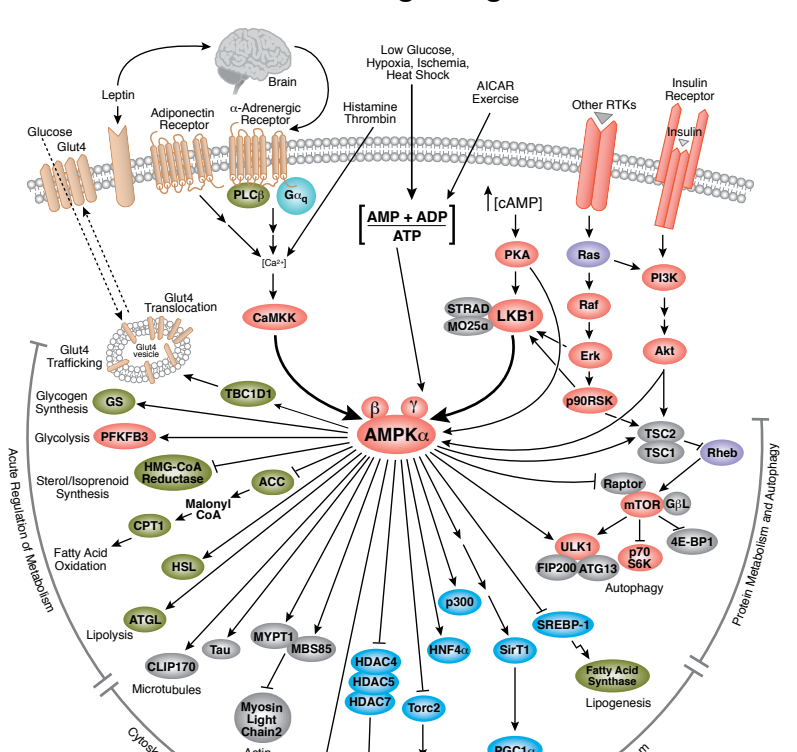
Insulin Receptor Signaling



Warburg Effect



AMPK Signaling



CELL CYCLE CONTROL/DNA DAMAGE

The above pathways were revised and updated by Cell Signaling Technology scientists and external opinion leaders in June 2012. For complete pathway descriptions and selected reviews please visit www.cellsignal.com/reference. © 09/2012 Printed in the USA on 55% recycled paper (35% post-consumer waste fiber) using soy inks and processed chlorine free.



