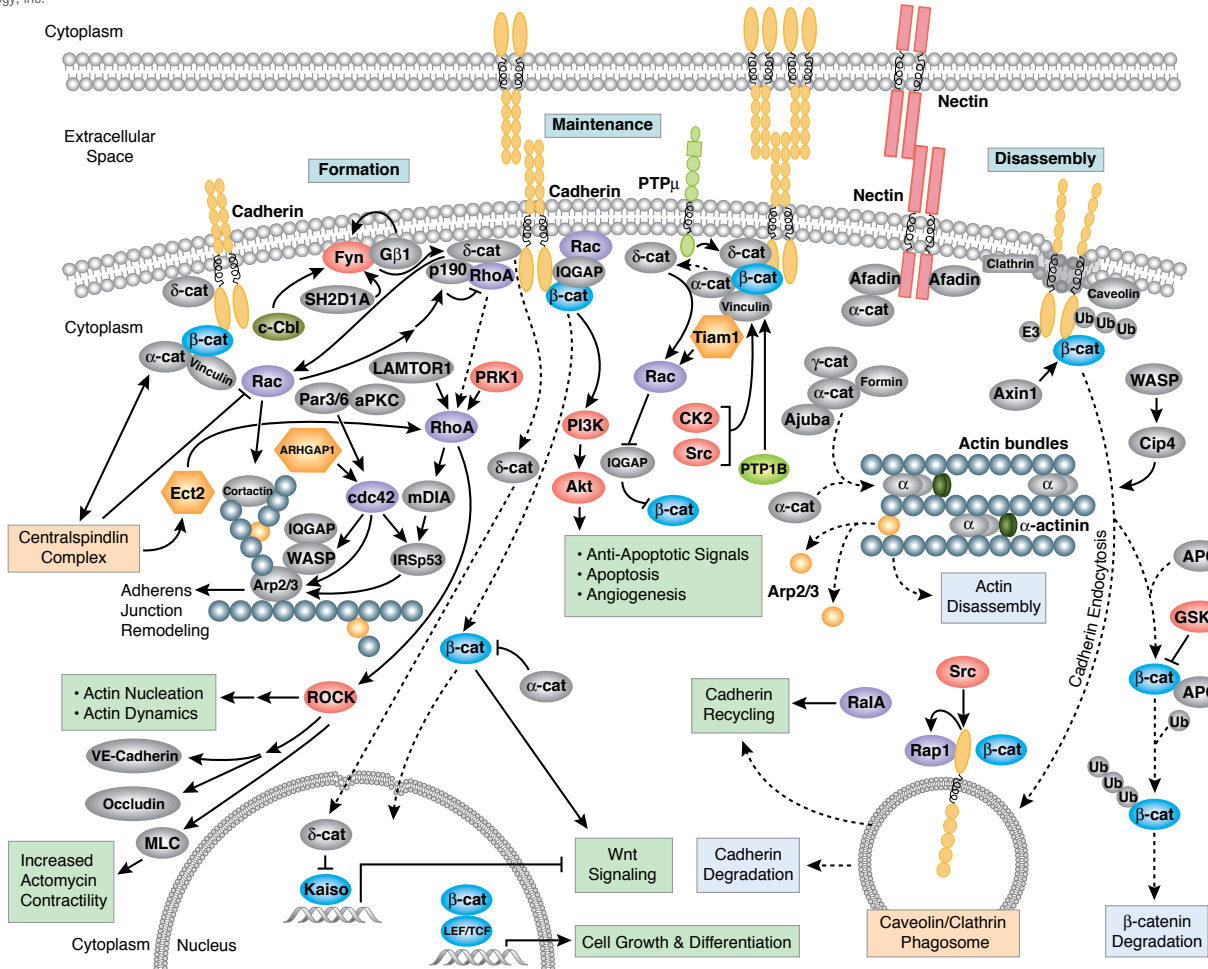


Cell Biology

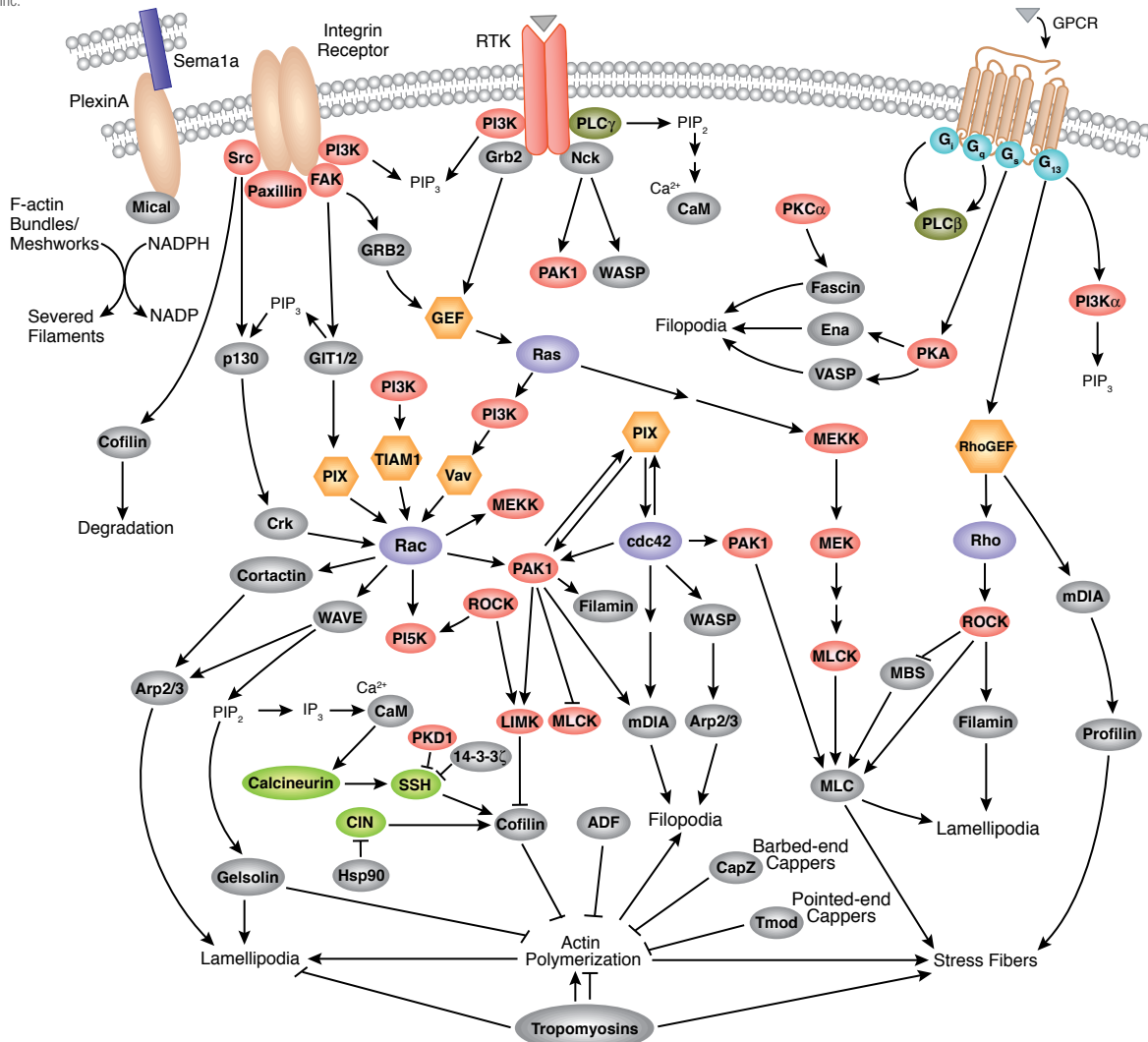
Adherens Junction Dynamics

© 2008–2015 Cell Signaling Technology, Inc.



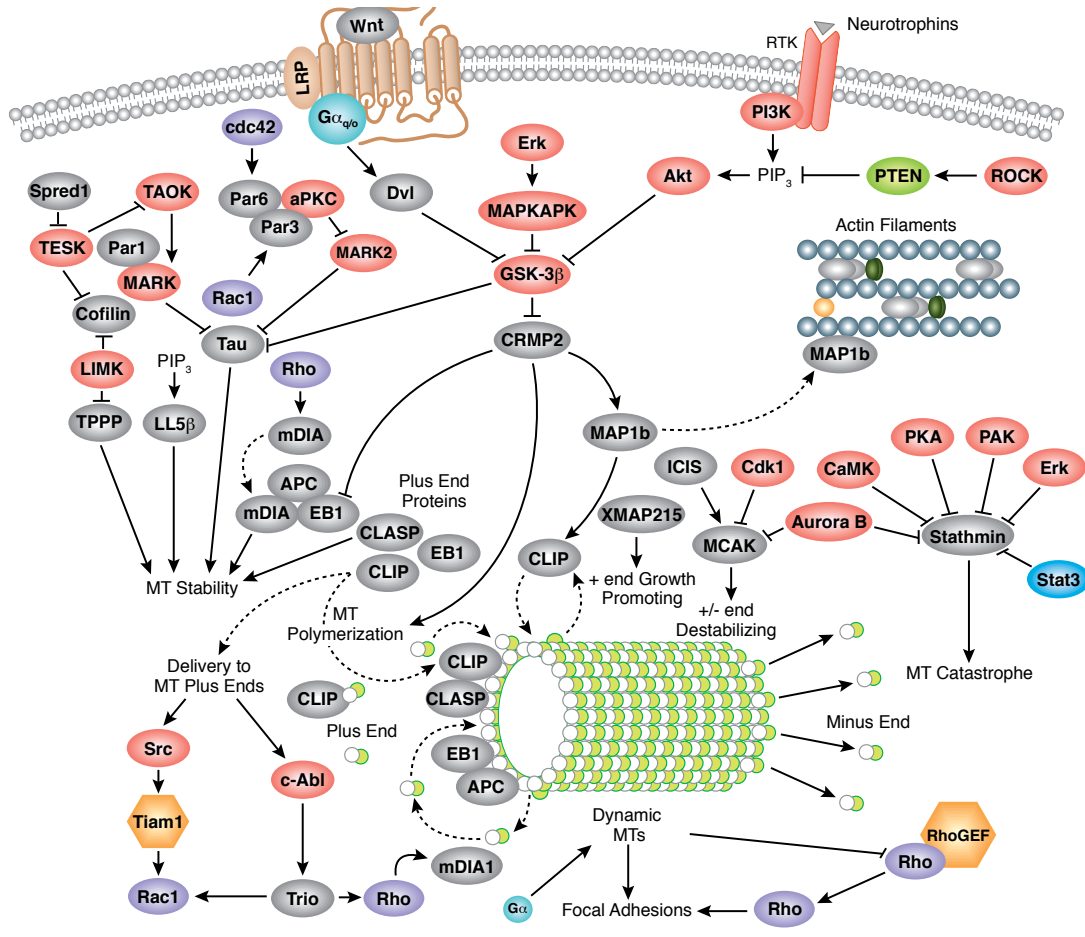
Regulation of Actin Dynamics

© 2008–2015 Cell Signaling Technology, Inc.



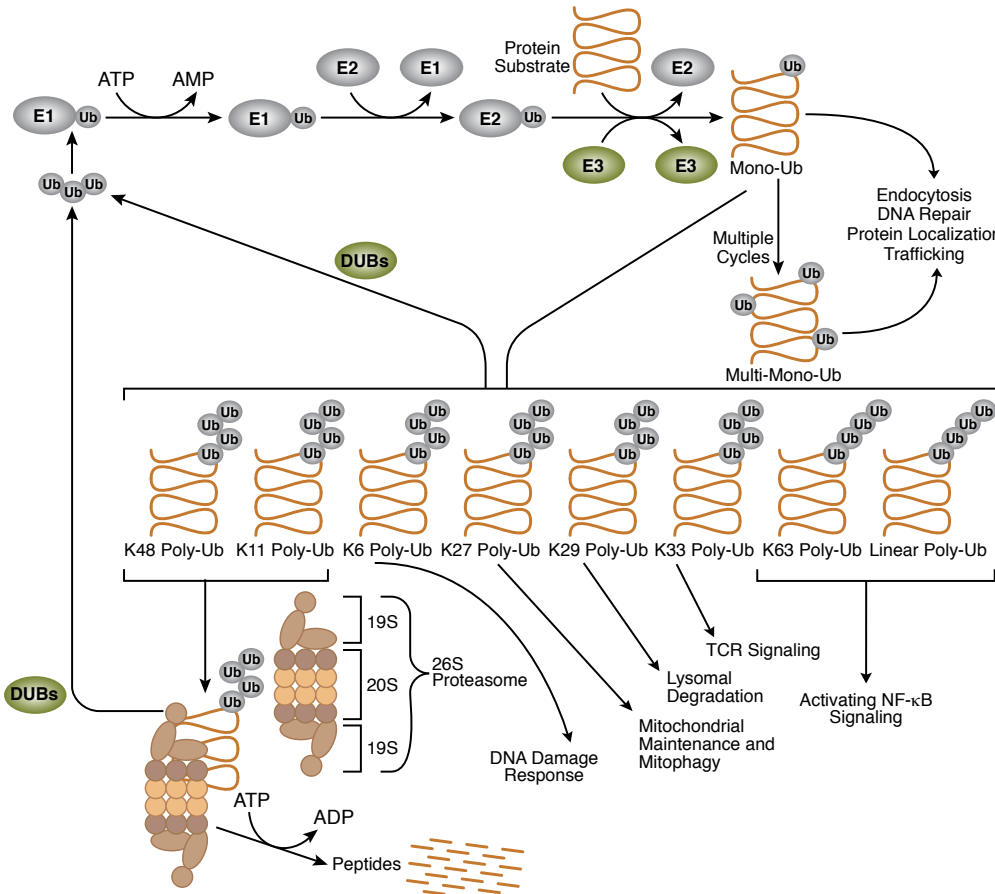
Regulation of Microtubule Dynamics

© 2008–2015 Cell Signaling Technology, Inc.



Ubiquitin/Proteasome

© 2004–2015 Cell Signaling Technology, Inc.



- | | | | | | |
|--------------------------------------|--------------------------------------|---|---|-----------------|---------------------|
| → Direct Stimulatory Modification | → Multistep Inhibitory Modification | → Separation of Subunits or Cleavage Products | → Transcriptional Inhibitory Modification | ● Receptor | ● GTPase |
| → Direct Inhibitory Modification | → Tentative Stimulatory Modification | → Joining of Subunits | ● Kinase | ● Enzyme | ● G-protein |
| → Multistep Stimulatory Modification | → Tentative Inhibitory Modification | → Translocation | ● Phosphatase | ● pro-apoptotic | ● Acetylase |
| | | → Transcriptional Stimulatory Modification | ● Transcription Factor | ● pro-survival | ● Deacetylase |
| | | | ● Caspase | ● GAP/GEF | ● Ribosomal subunit |