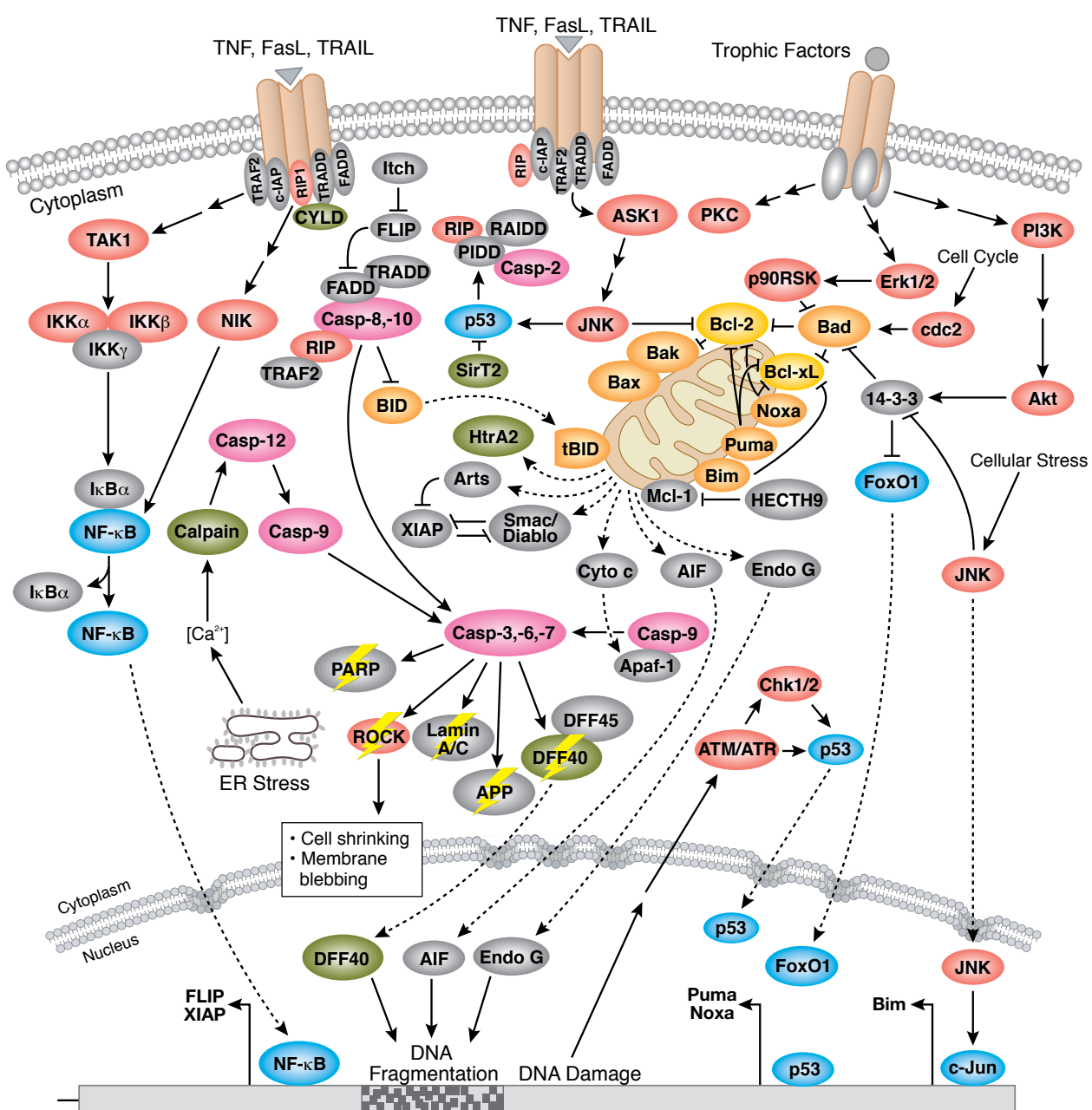


Cell Growth and Death

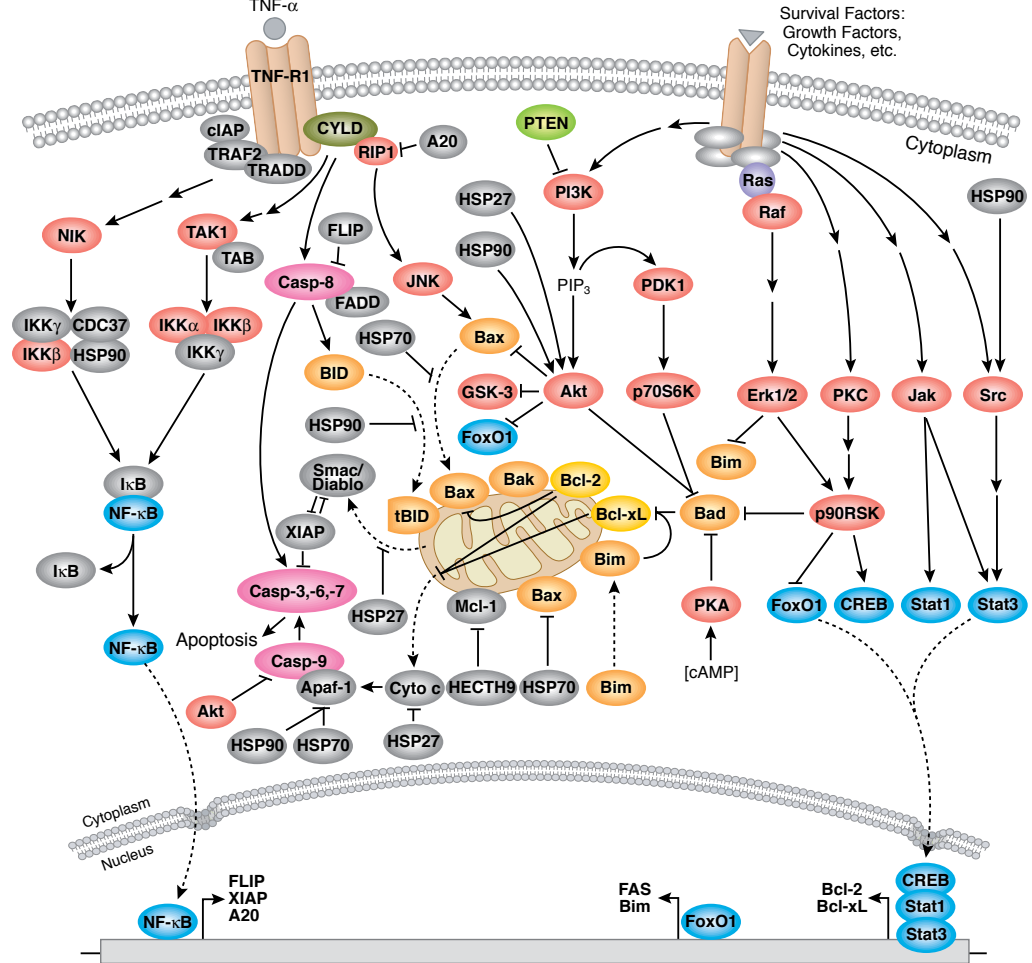
Regulation of Apoptosis Overview

© 2002-2015 Cell Signaling Technology, Inc.



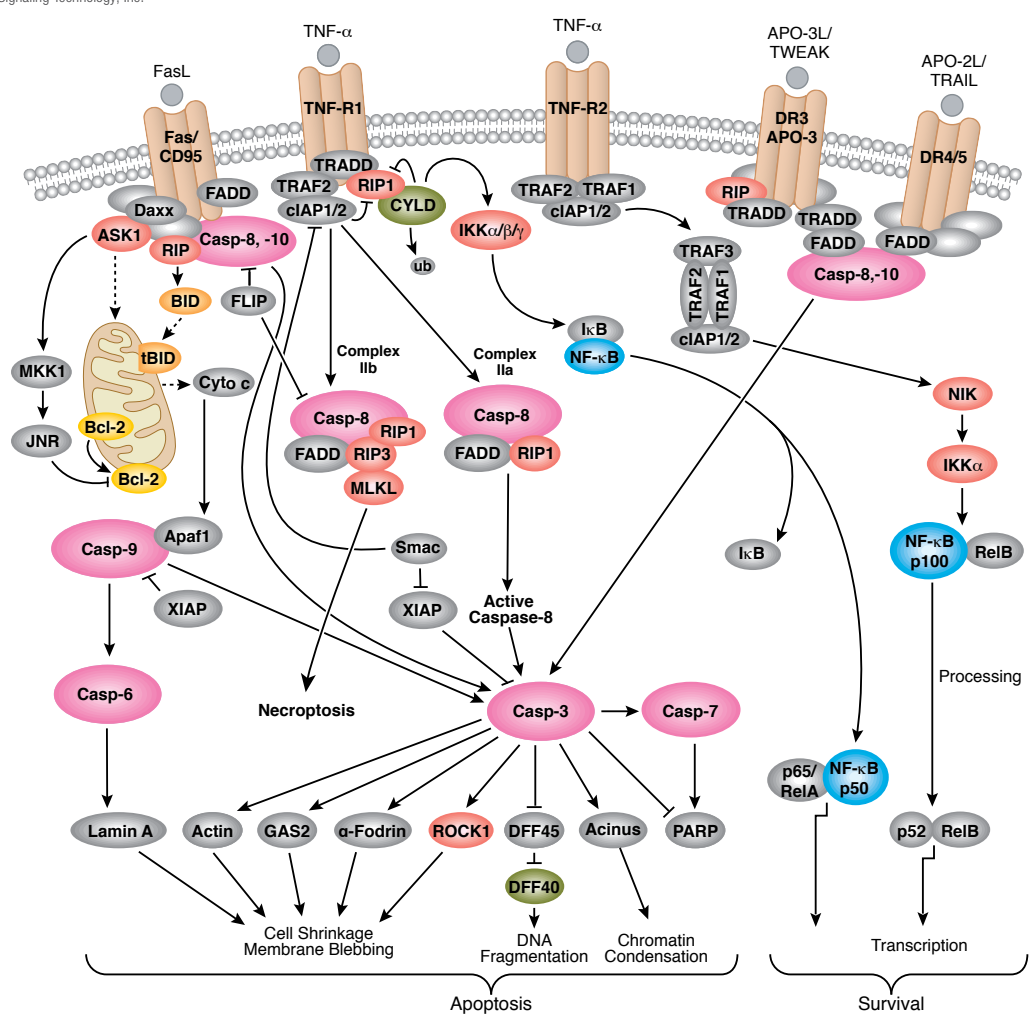
Inhibition of Apoptosis

© 2002-2015 Cell Signaling Technology, Inc.



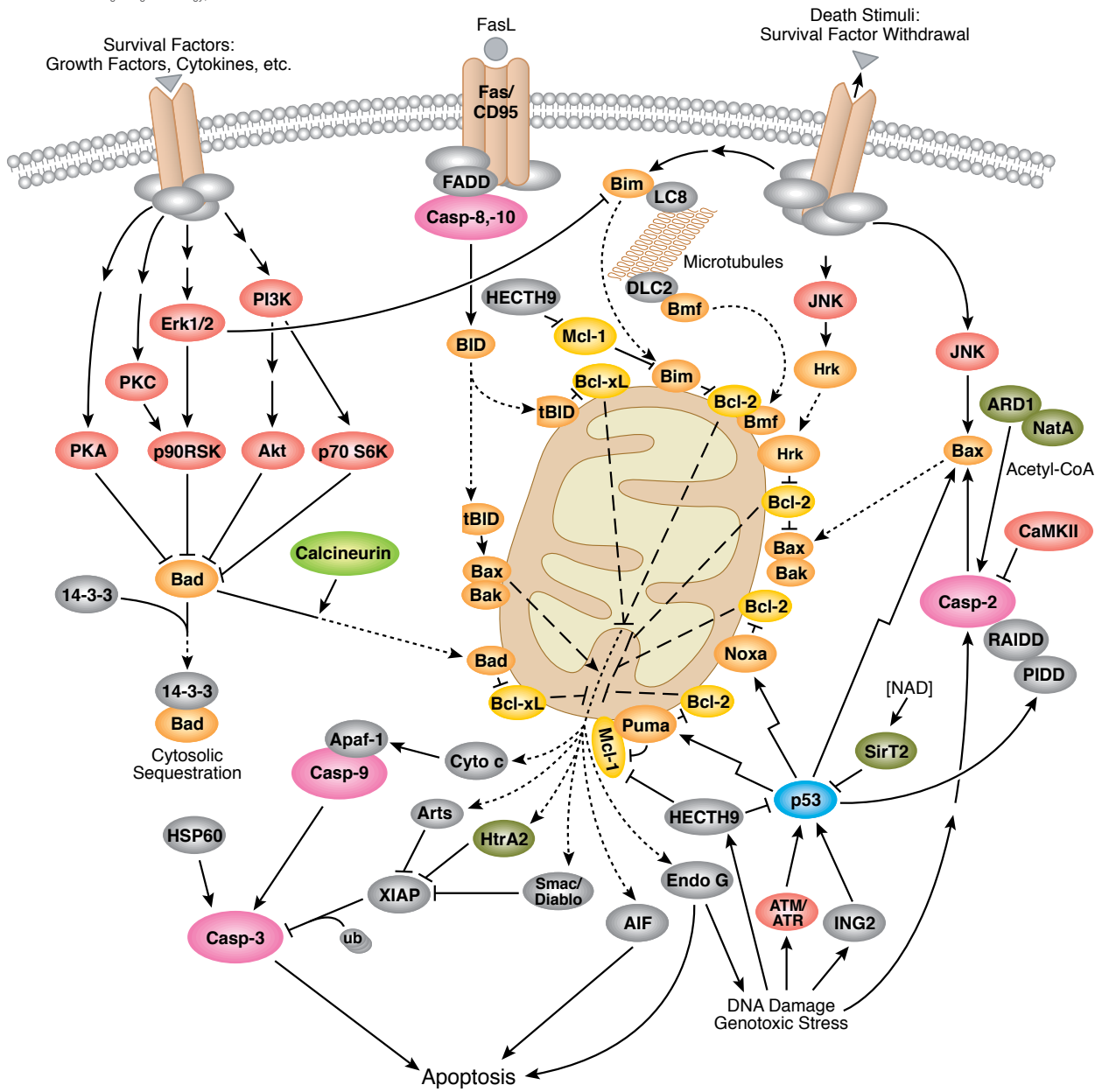
Death Receptor Signaling

© 2002-2015 Cell Signaling Technology, Inc.



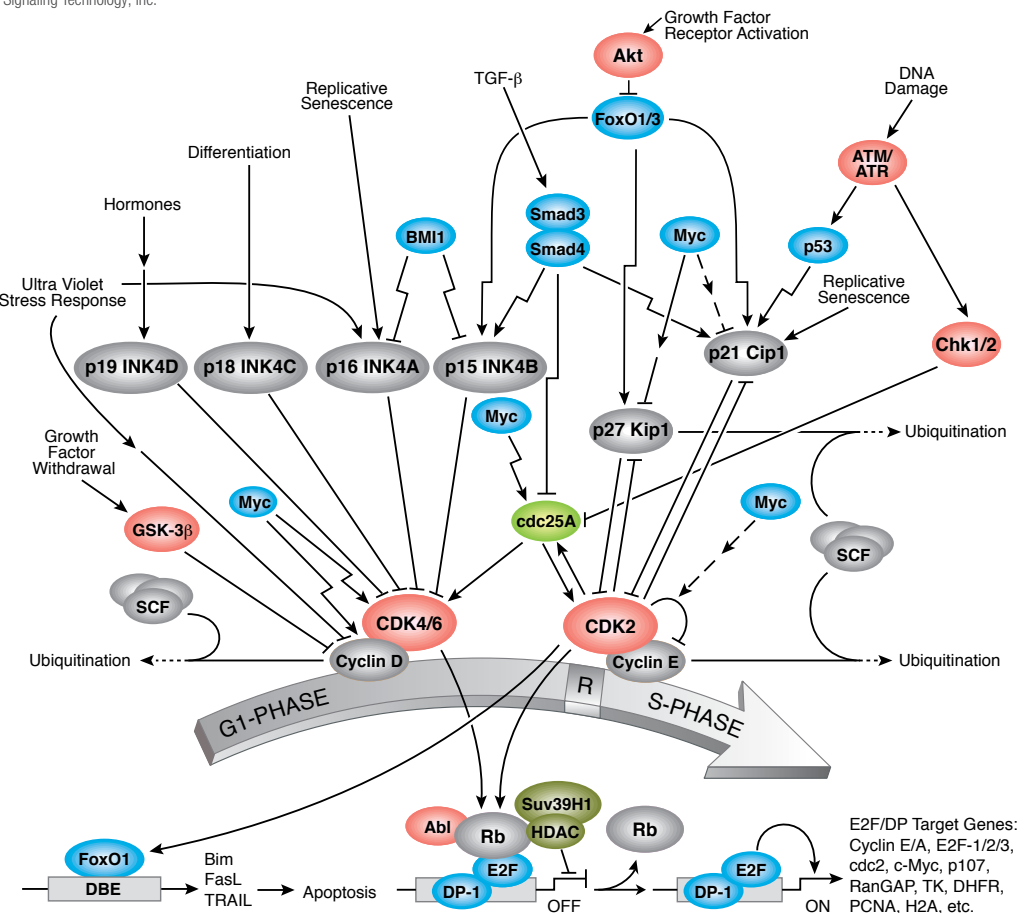
Mitochondrial Control of Apoptosis

© 2003-2015 Cell Signaling Technology, Inc.



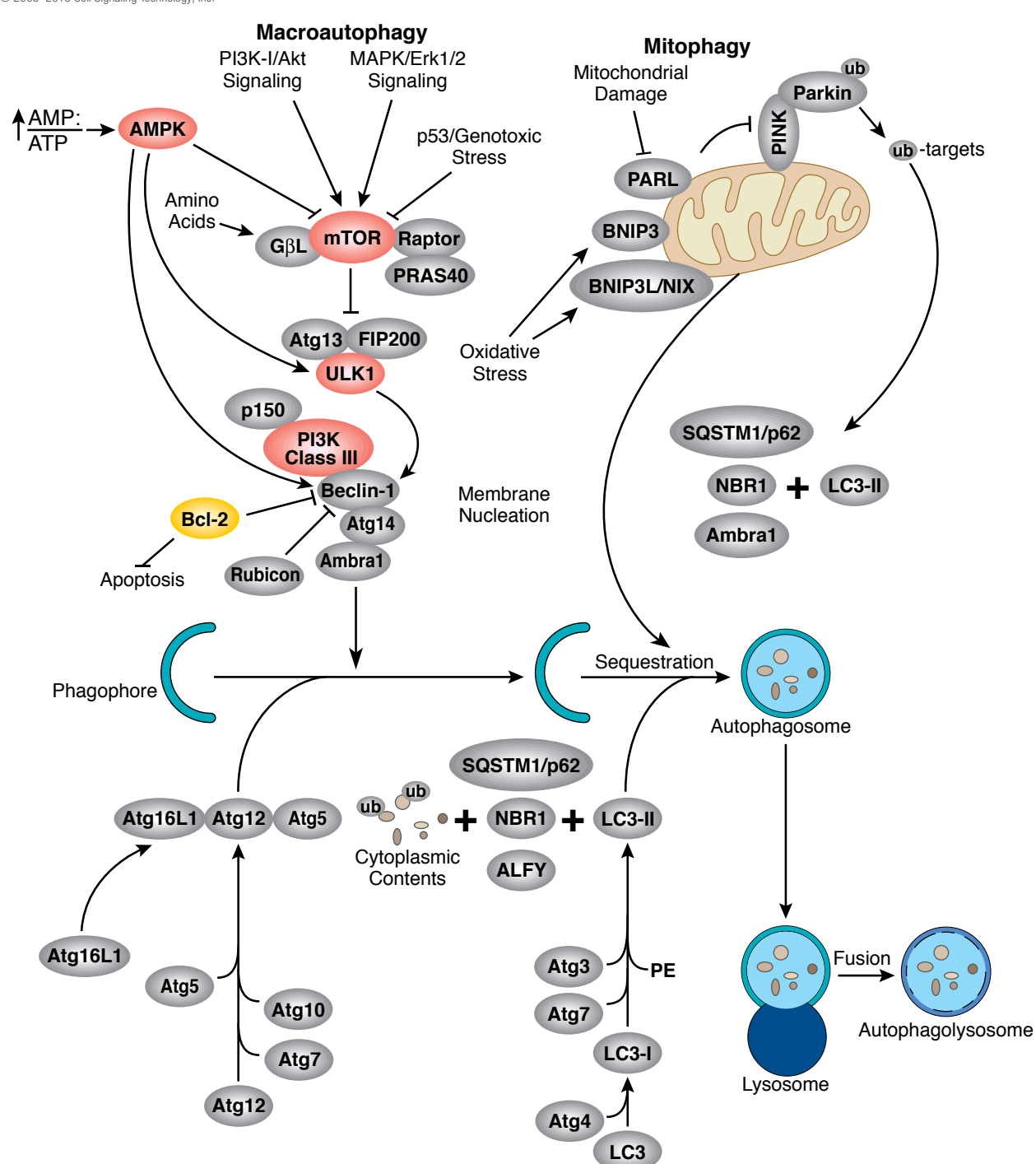
Cell Cycle Control: G1/S Checkpoint

© 2002-2015 Cell Signaling Technology, Inc.



Autophagy Signaling

© 2003-2015 Cell Signaling Technology, Inc.



Cell Cycle Control: G2/M DNA Damage Checkpoint

© 2002-2015 Cell Signaling Technology, Inc.

