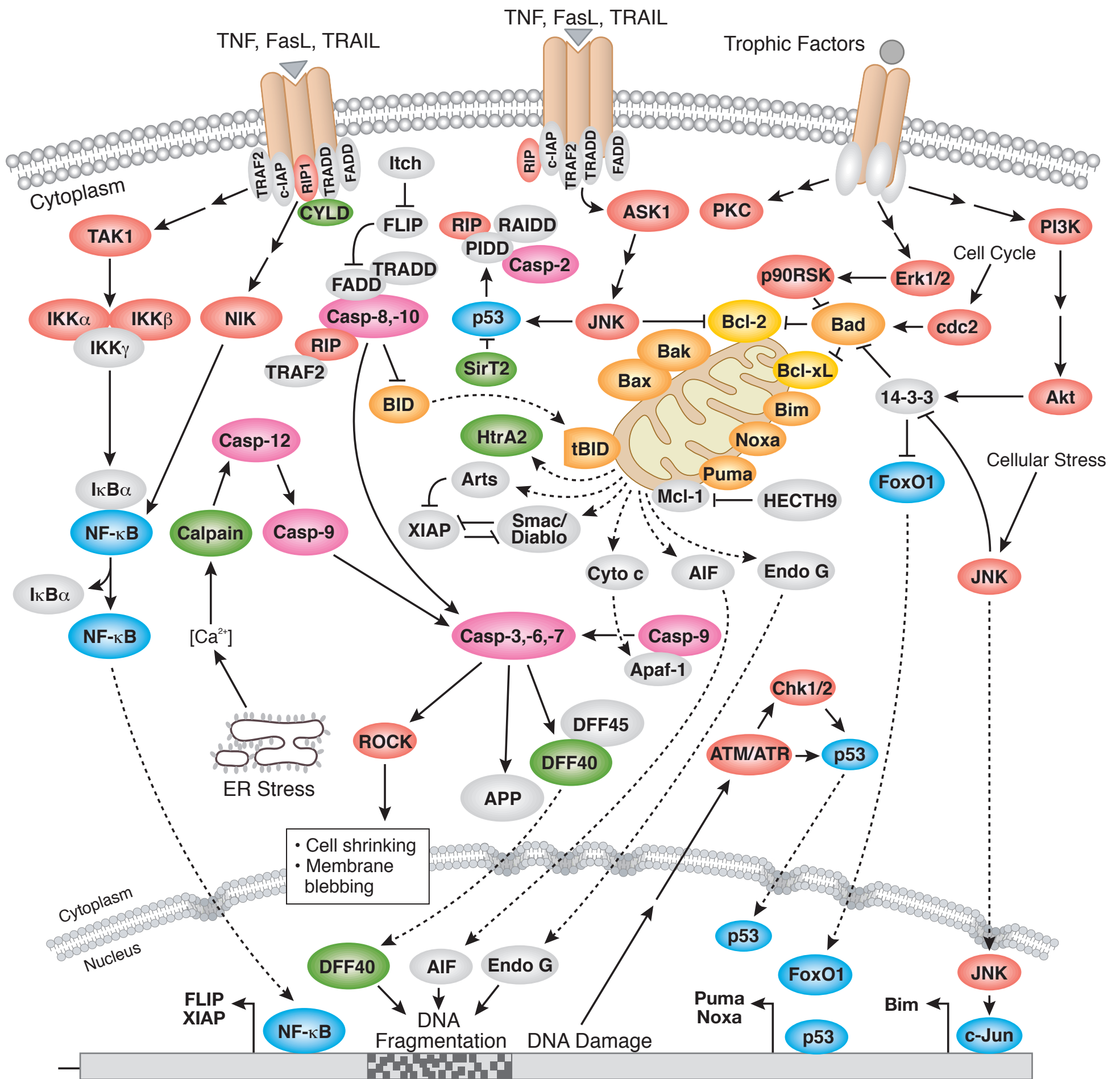
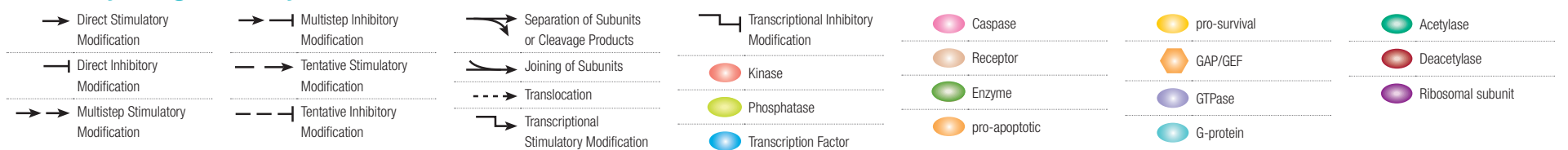


Regulation of Apoptosis



Pathway Diagram Key



Select Reviews

This pathway diagram has been assembled by Key Opinion Leaders in the field of apoptosis in collaboration with Cell Signaling Technology (CST) scientists. CST Pathways are reviewed and updated frequently to provide succinct and current overviews of selected signaling pathways. We would like to thank Prof. Junying Yuan, Harvard Medical School, Boston, MA, for reviewing this diagram.

Please visit www.cellsignal.com/pathways to find out more.

Degterev, A. and Yuan, J. (2008) *Nat. Rev. Mol. Cell Biol.* 9, 378–390.

Fuchs, Y. and Steller, H. (2011) *Cell* 147, 742–758.

Indran, I.R., Tufo, G., Pervaiz, S., and Brenner, C. (2011) *Biochim. Biophys. Acta.* 1807, 735–745.

Kaufmann, T., Strasser, A., and Jost, P.J. (2012) *Cell Death Differ.* 19, 42–50.

Kurokawa, M. and Kornbluth, S. (2009) *Cell* 138, 838–854.

Pradelli, L.A., Bénétteau, M., and Ricci, J.E. (2010) *Cell. Mol. Life Sci.* 67, 1589–1597.

Van Herreweghe, F., Festjens, N., Declercq, W., and Vandennebe, P. (2010) *Cell. Mol. Life Sci.* 67, 1567–1579.