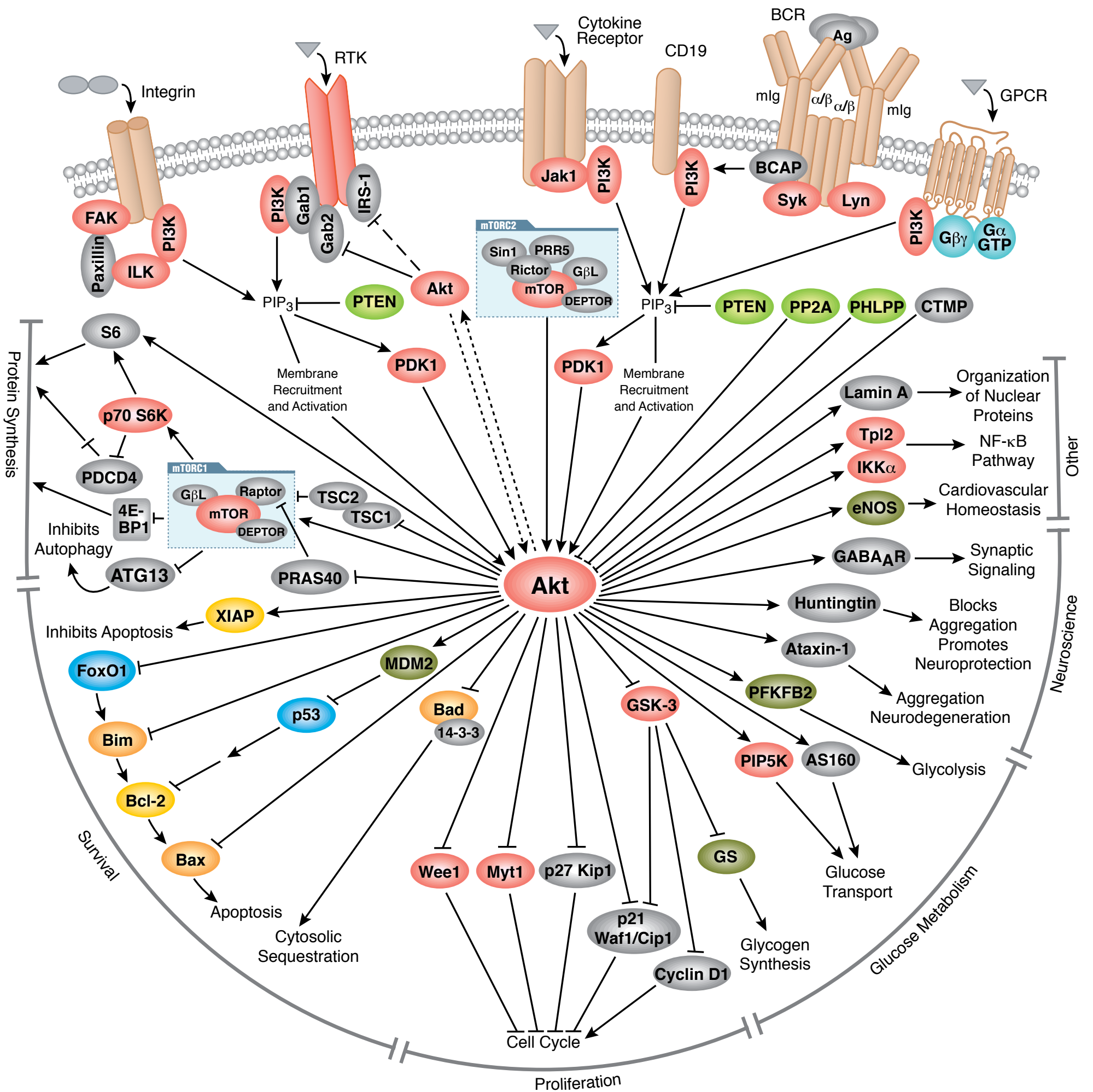
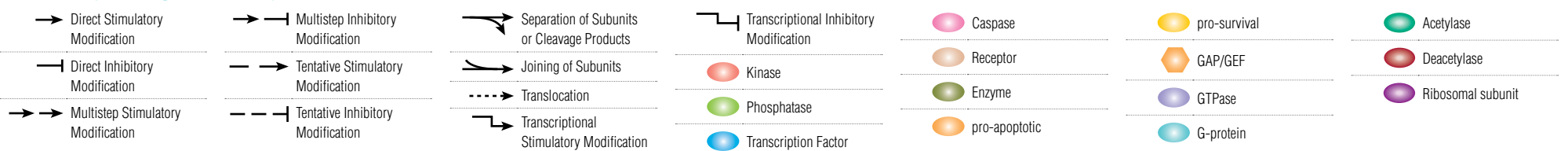


# PI3 Kinase/Akt Signaling



## Pathway Diagram Key



## Select Reviews

This pathway diagram and the Akt Substrates Guide have been assembled by Key Opinion Leaders in the field of Akt Signaling 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. Michael Scheid, York University of Toronto, Ontario, for reviewing this diagram.

Please visit [www.cellsignal.com/pathways](http://www.cellsignal.com/pathways) to find out more.

Carnero, A., Blanco-Aparicio, C., Renner, O., Link, W., and Leal, J.F. (2008) *Curr. Cancer Drug Targets* 8, 187–198.

Duronio, V. (2008) *Biochem. J.* 415, 333–344.

Hers, I., Vincent, E.E., and Tavaré, J.M. (2011) *Cell Signal.* 10, 1515–1527.

Hussain, K., Challis, B., Rocha, N., Payne, F., Minic, M., Thompson, A., Daly, A., Scott, C., Harris, J., Smillie, B.J., Savage, D.B., Ramaswami, U., De Lonlay, P., O'Rahilly, S., Barroso, I., and Semple, R.K. (2011) *Science* 334, 474.

Liu, P., Cheng, H., Roberts, T.M., and Zhao, J.J. (2009) *Nat. Rev. Drug Discov.* 8, 627–644.

Manning, B.D. and Cantley, L.C. (2007) *Cell* 129, 1261–1274.

Sheppard, K., Kinross, K.M., Solomon, B., Pearson, R.B., and Phillips, W.A. (2012) *Crit. Rev. Oncol.* 17, 69–95.

Salmena, L., Carracedo, A., and Pandolfi, P.P. (2008) *Cell* 133, 403–414.

Song, M.S., Salmena, L., and Pandolfi, P.P. (2012) *Nat. Rev. Mol. Cell Biol.* 13, 283–296.