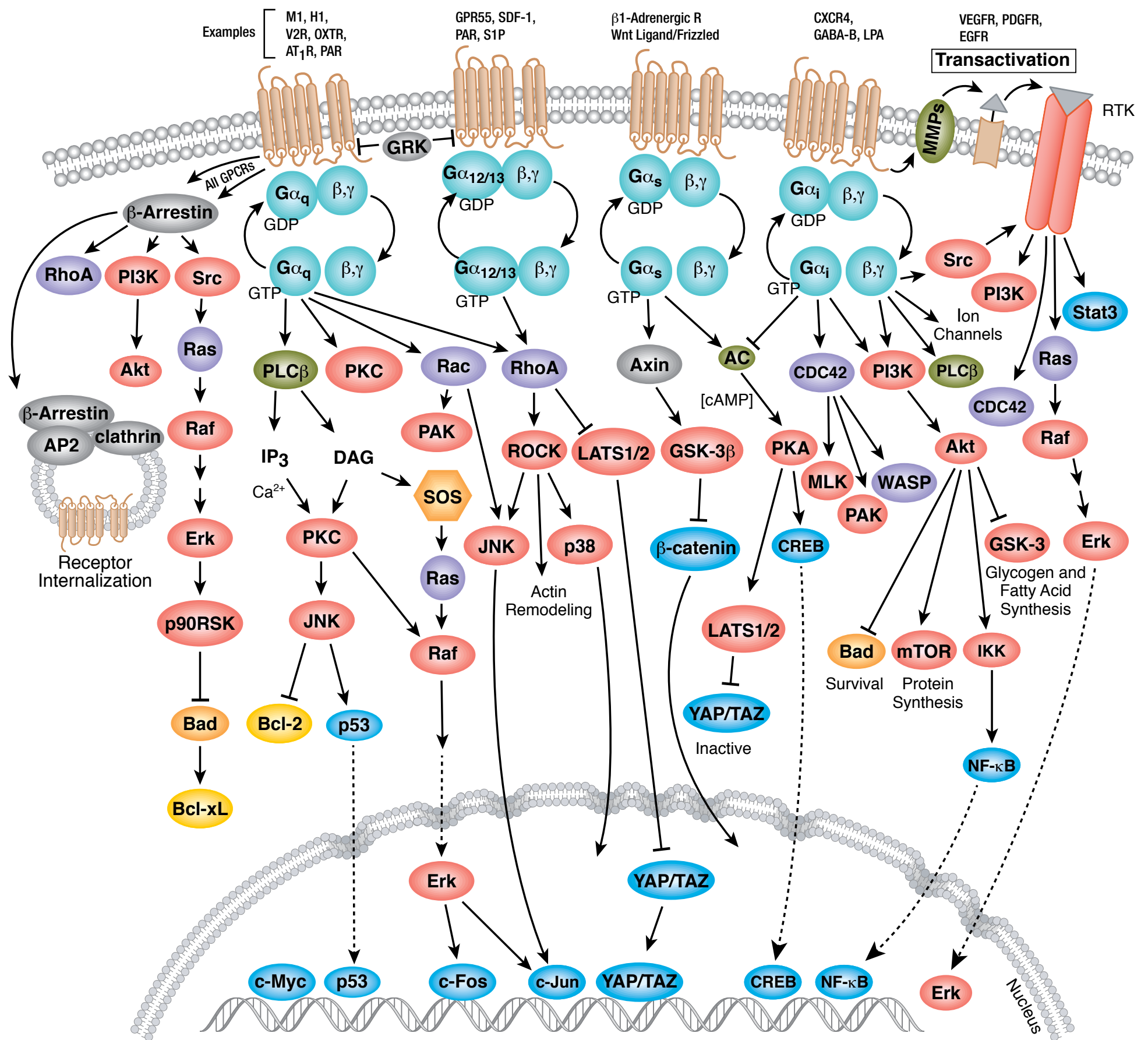
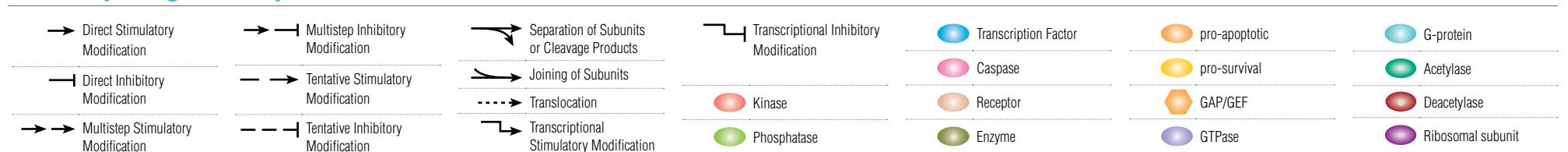


# G-Protein Coupled Receptor Signaling: Overview



## Pathway Diagram Key



## Select Reviews

This pathway diagram has been assembled by key opinion leaders in the field of G-Protein Coupled Receptor Signaling in collaboration with Cell Signaling Technology (CST) scientists. CST Pathways are reviewed and updated frequently to provide succinct and current overviews of the subject area. We would like to thank Dr. Jonathan Violin, Trevana Inc., King of Prussia, PA for reviewing this diagram.

Entschladen, F., Zänker, K.S., and Powe, D.G. (2011) *Cell Cycle* 10, 1086–1091. | Audet, M. and Bouvier, M. (2012) *Cell* 151, 14–23. | Lappano, R. and Maggiolini, M. (2012) *Acta. Pharmacol. Sin.* 33, 351–362. | O'Havre, M., Degese, M.S., and Gutkind, J.S. (2014) *Curr. Opin. Cell Biol.* 27C, 126–135. | Lappano, R. and Maggiolini, M. (2011) *Nat. Rev. Drug Discov.* 10, 47–60. | Yu, F.X. and Guan, K.L. (2013) *Genes Dev.* 27, 355–371. | Smrcka, A.V. (2008) *Cell. Mol. Life Sci.* 65, 2191–2214. | Kahn, S.M., Sleno, R., Gora, S., Zylbergold, P., Laverdure, J.P., Labbé, J.C., Miller, G.J., and Hébert, T.E. (2013) *Pharmacol. Rev.* 65, 545–577. | Shukla, A.K., Xiao, K., and Lefkowitz, R.J. (2011) *Trends Biochem. Sci.* 36, 457–469. | Raiagopal, S., Raiagopal, K., and Lefkowitz, R.J. (2010) *Nat. Rev. Drug Discov.* 9, 373–386.