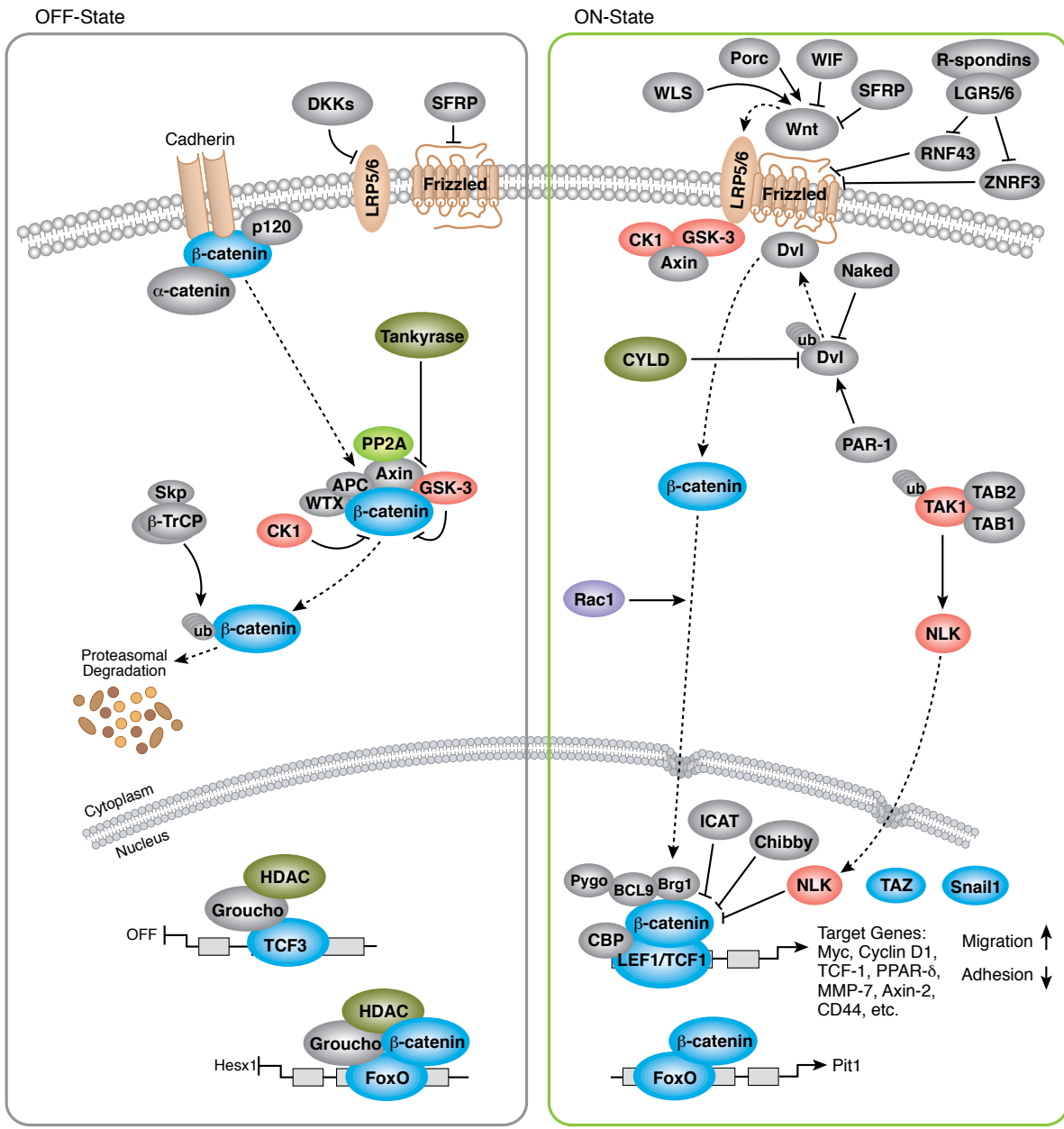


Development and Differentiation

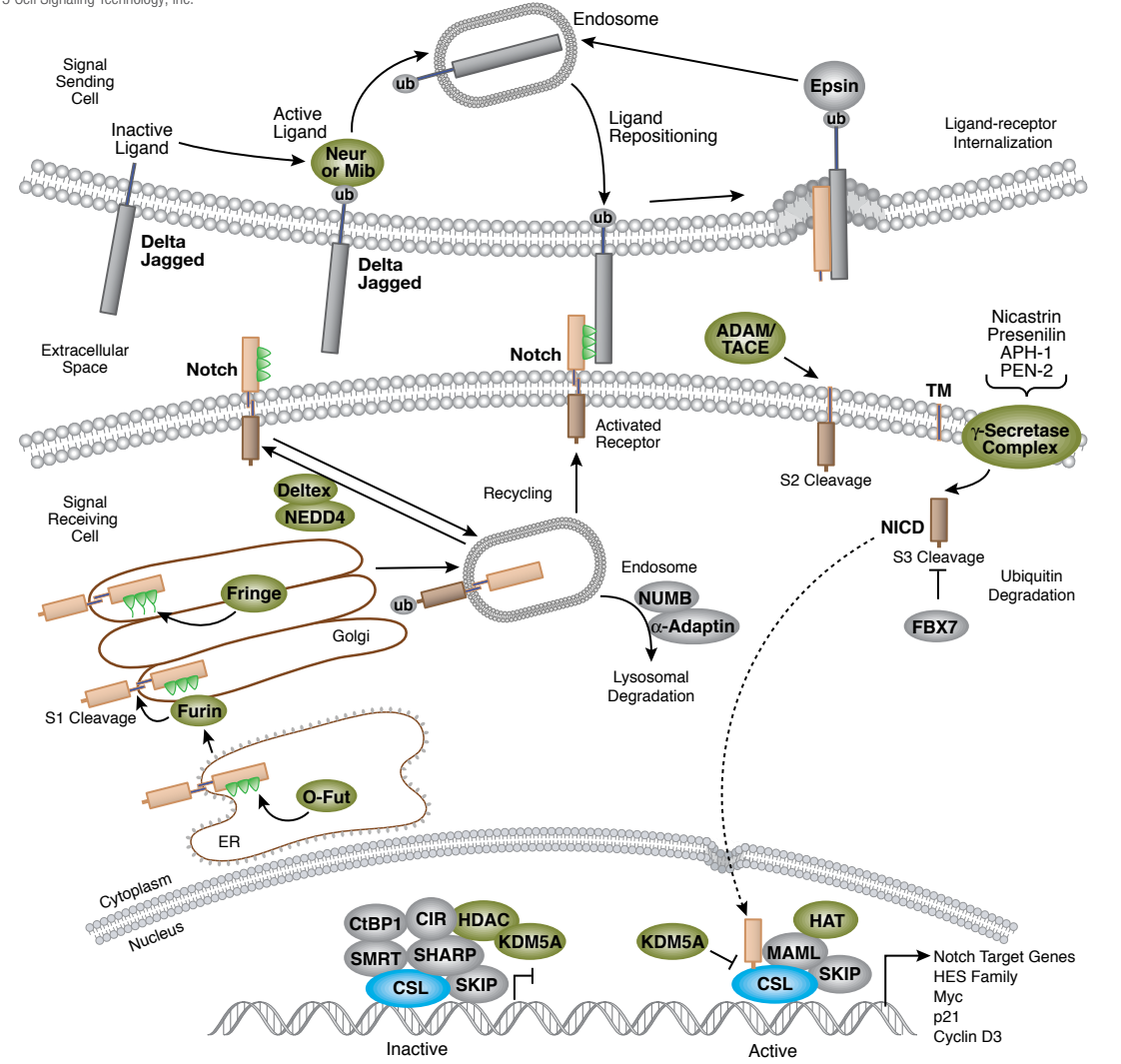
Wnt/ β -Catenin Signaling

© 2003–2015 Cell Signaling Technology, Inc.



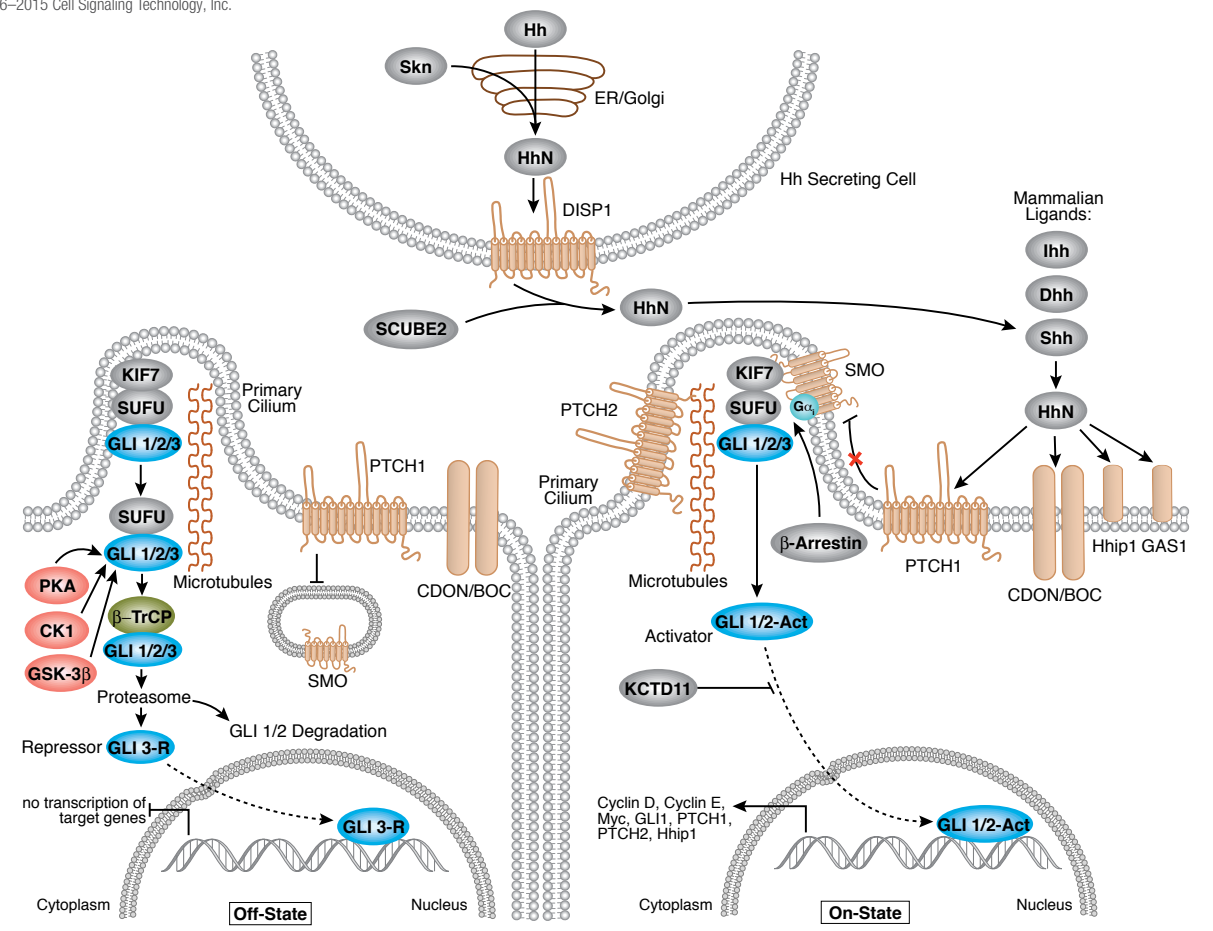
Notch Signaling

© 2006–2015 Cell Signaling Technology, Inc.



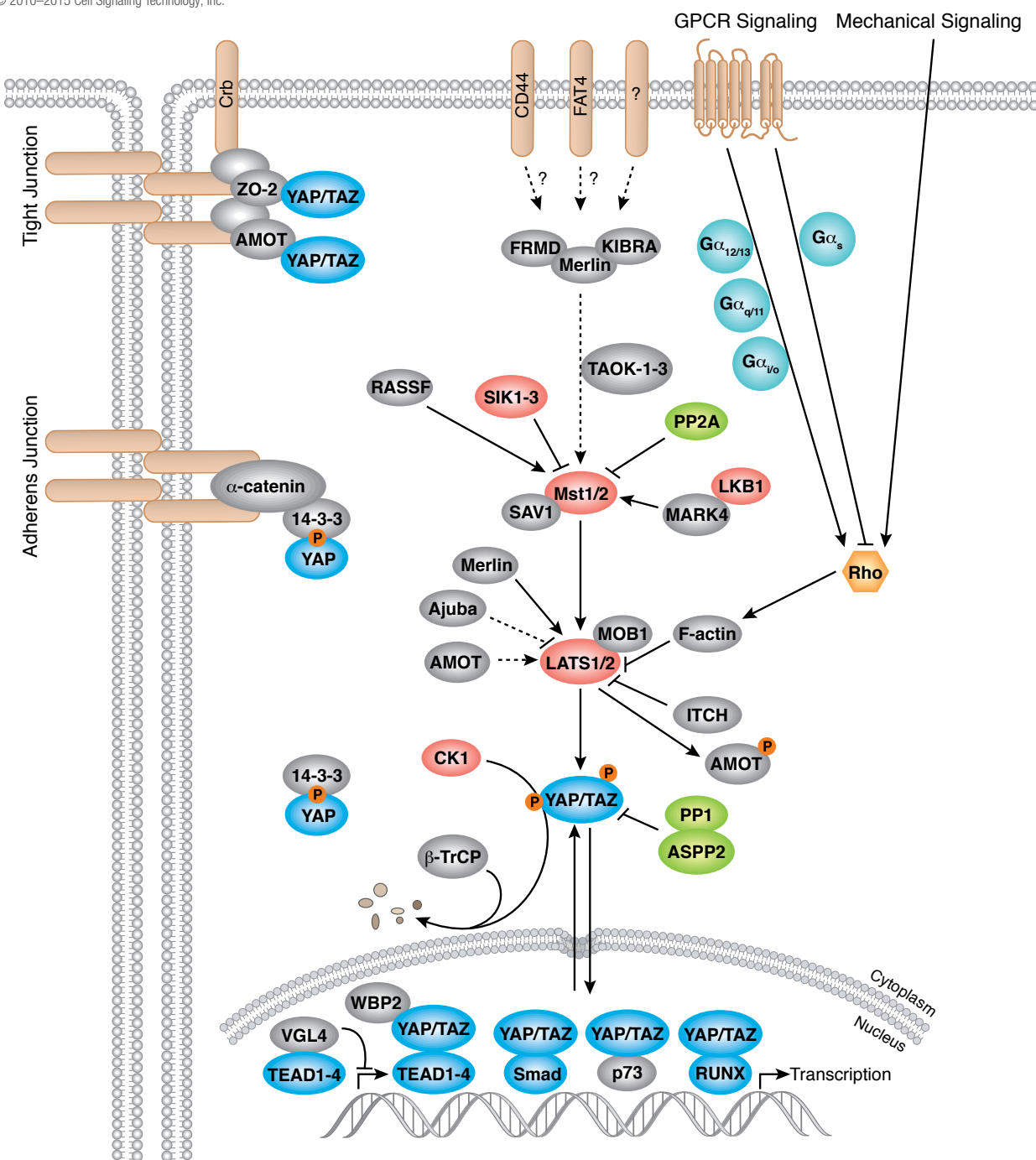
Hedgehog Signaling

© 2006–2015 Cell Signaling Technology, Inc.



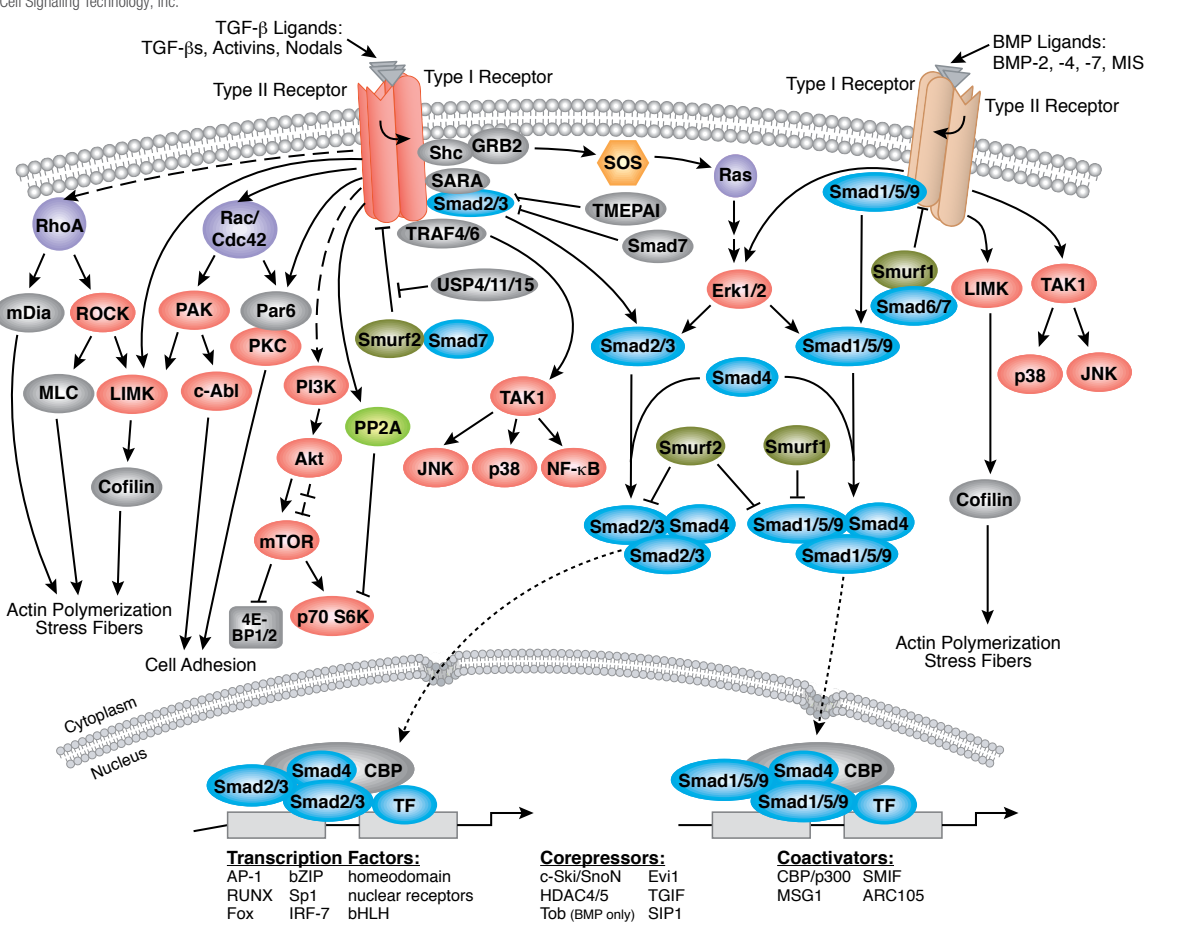
Hippo Signaling

© 2010–2015 Cell Signaling Technology, Inc.



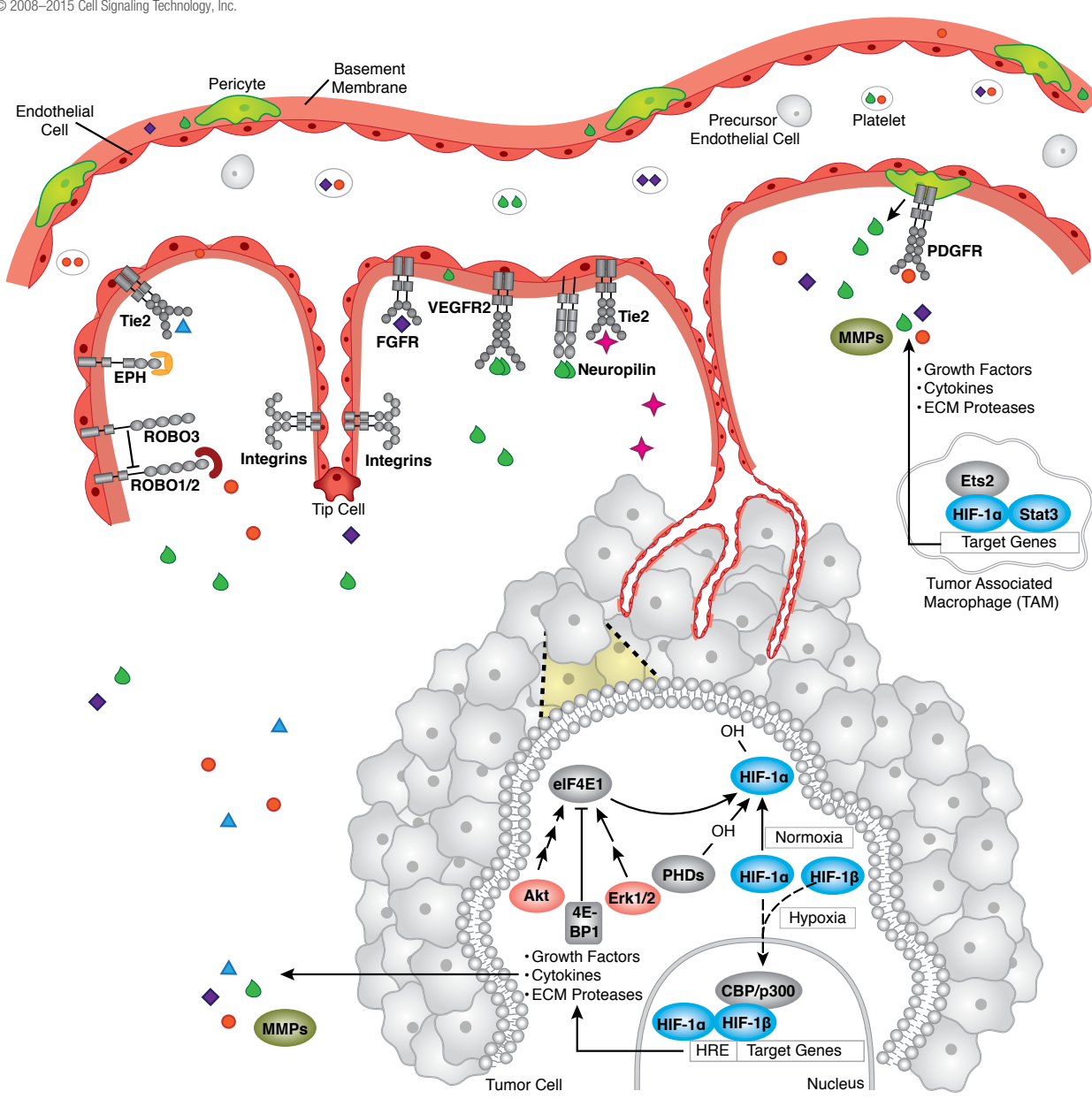
TGF- β Signaling

© 2003–2015 Cell Signaling Technology, Inc.



Angiogenesis Signaling in Tumor Neovascularization

© 2006–2015 Cell Signaling Technology, Inc.



ESC Pluripotency Differentiation

© 2009–2015 Cell Signaling Technology, Inc.

