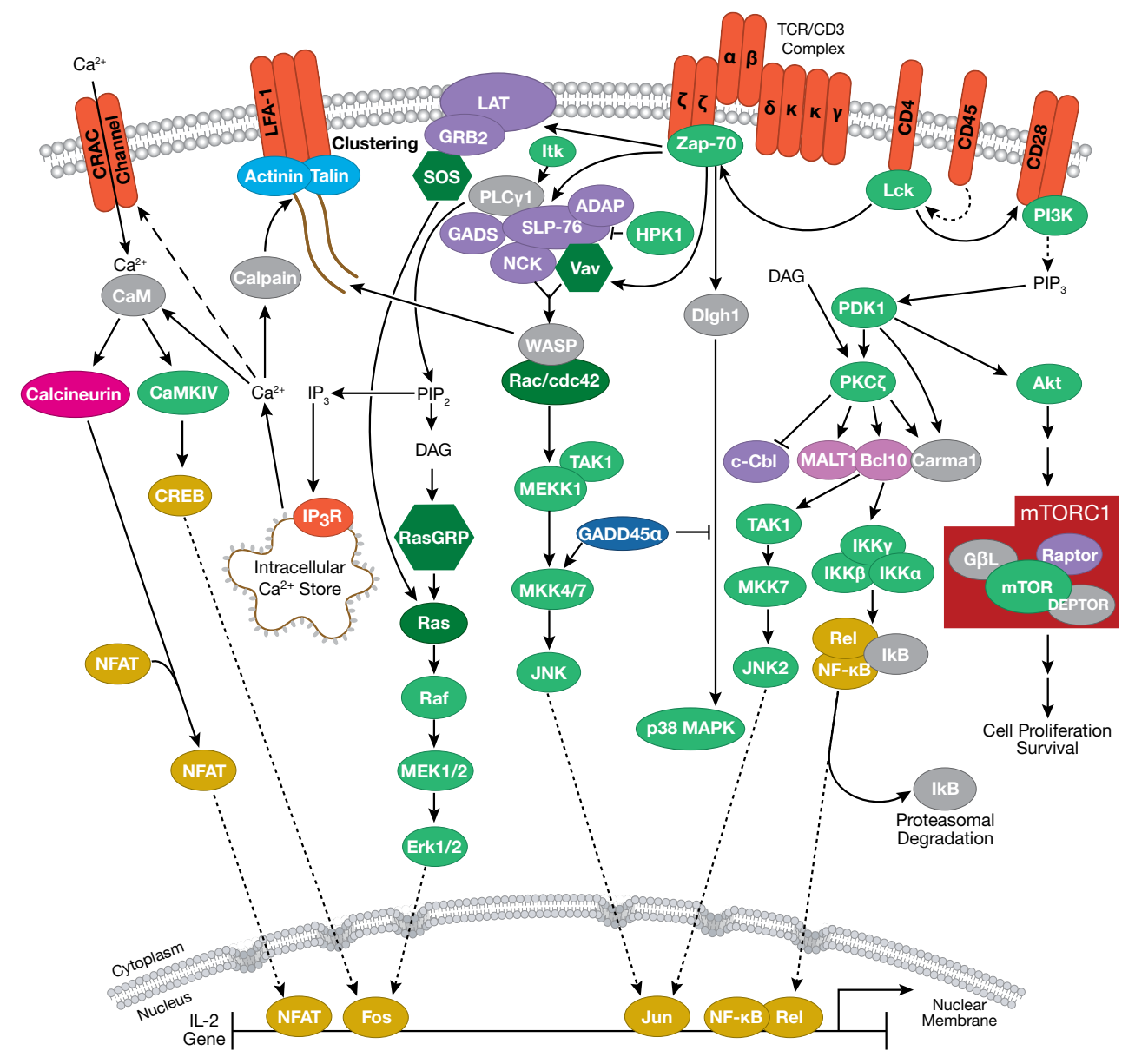


CST SIGNALING PATHWAY DIAGRAMS FOR Immunology

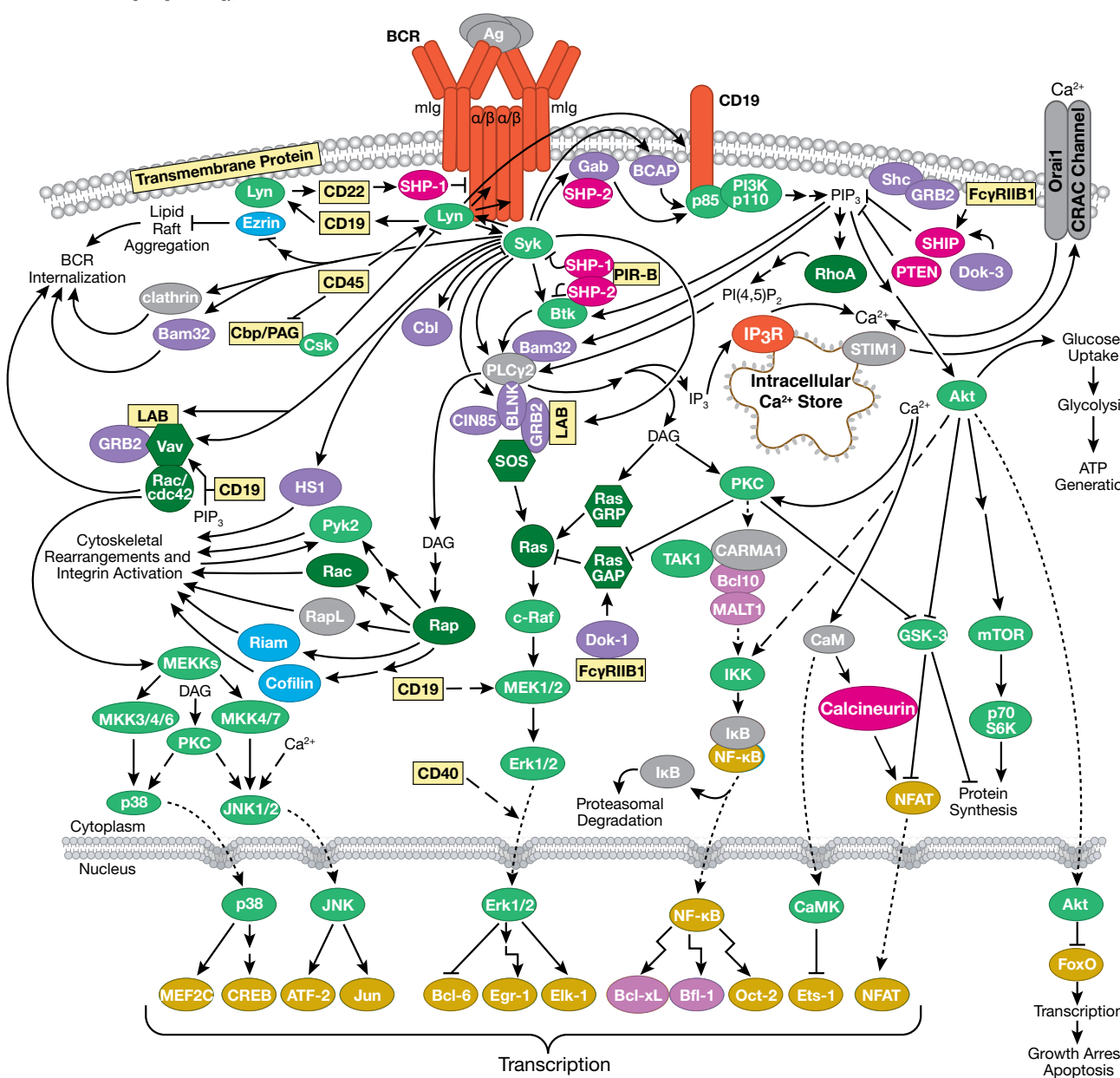
T Cell Receptor Signaling

© 2014-2018 Cell Signaling Technology, Inc.



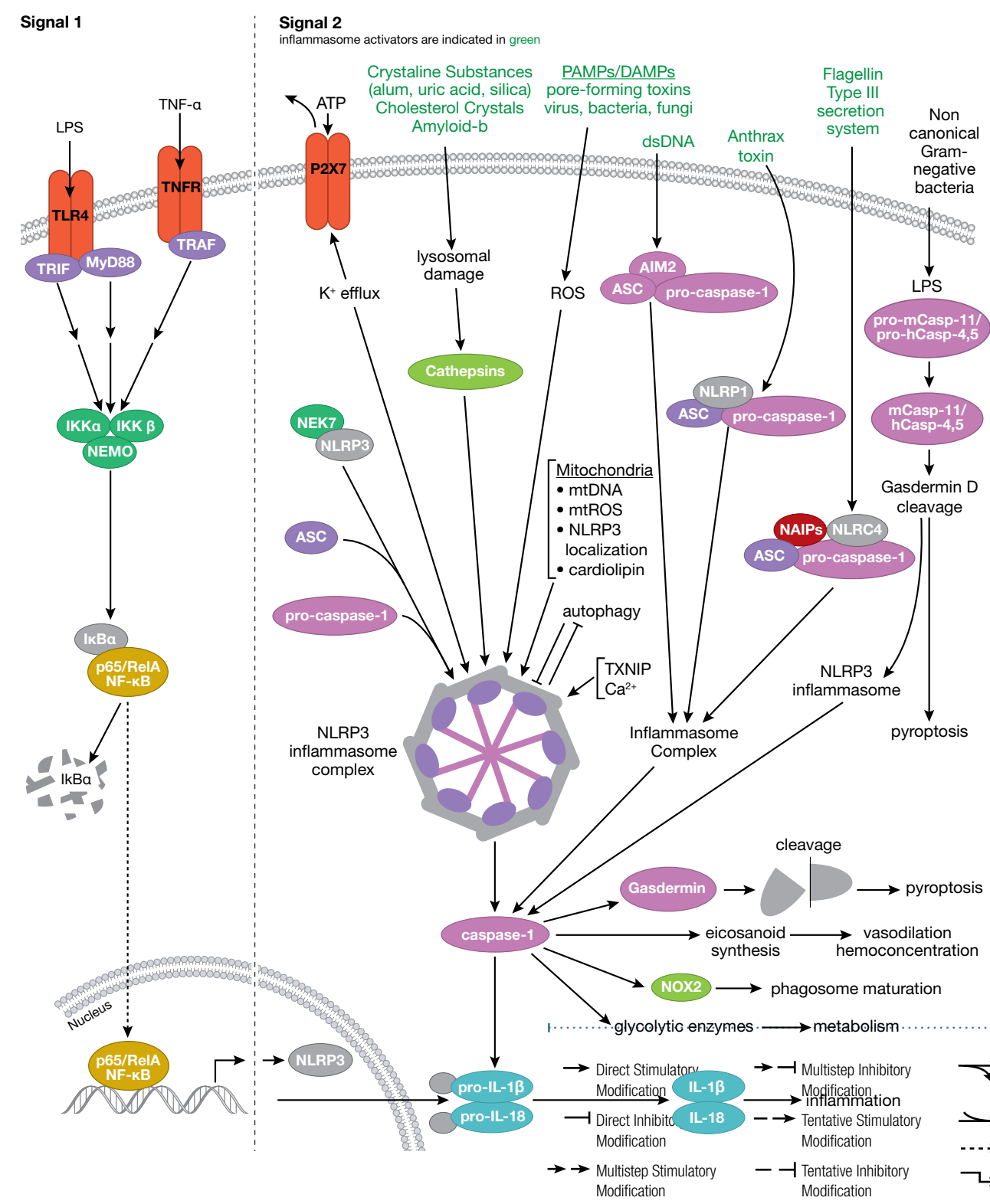
B Cell Receptor Signaling

© 2002-2018 Cell Signaling Technology, Inc.



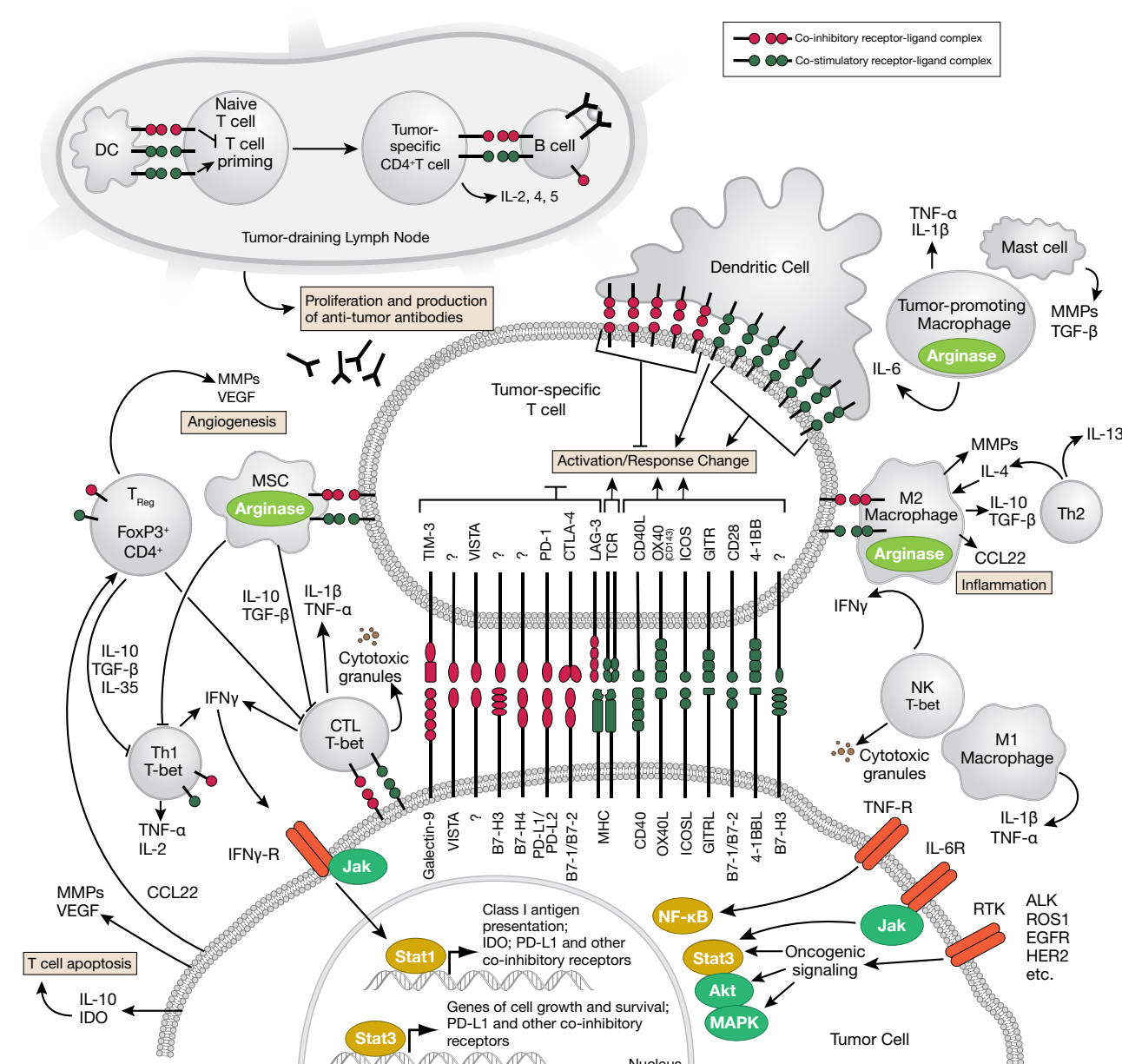
Inflammasome Signaling

© 2014-2018 Cell Signaling Technology, Inc.



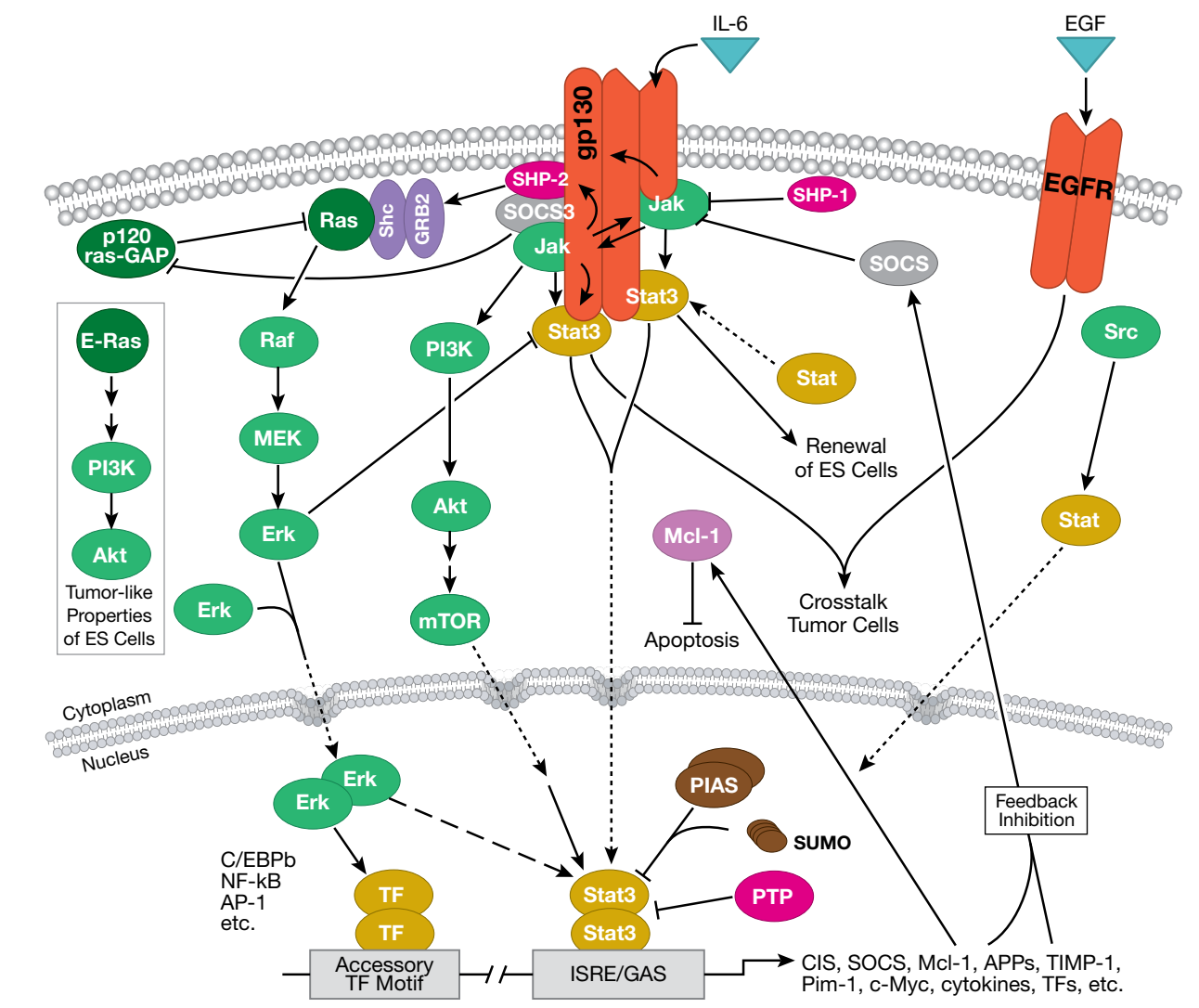
Immune Checkpoint Signaling in the Tumor Microenvironment

© 2014-2018 Cell Signaling Technology, Inc.



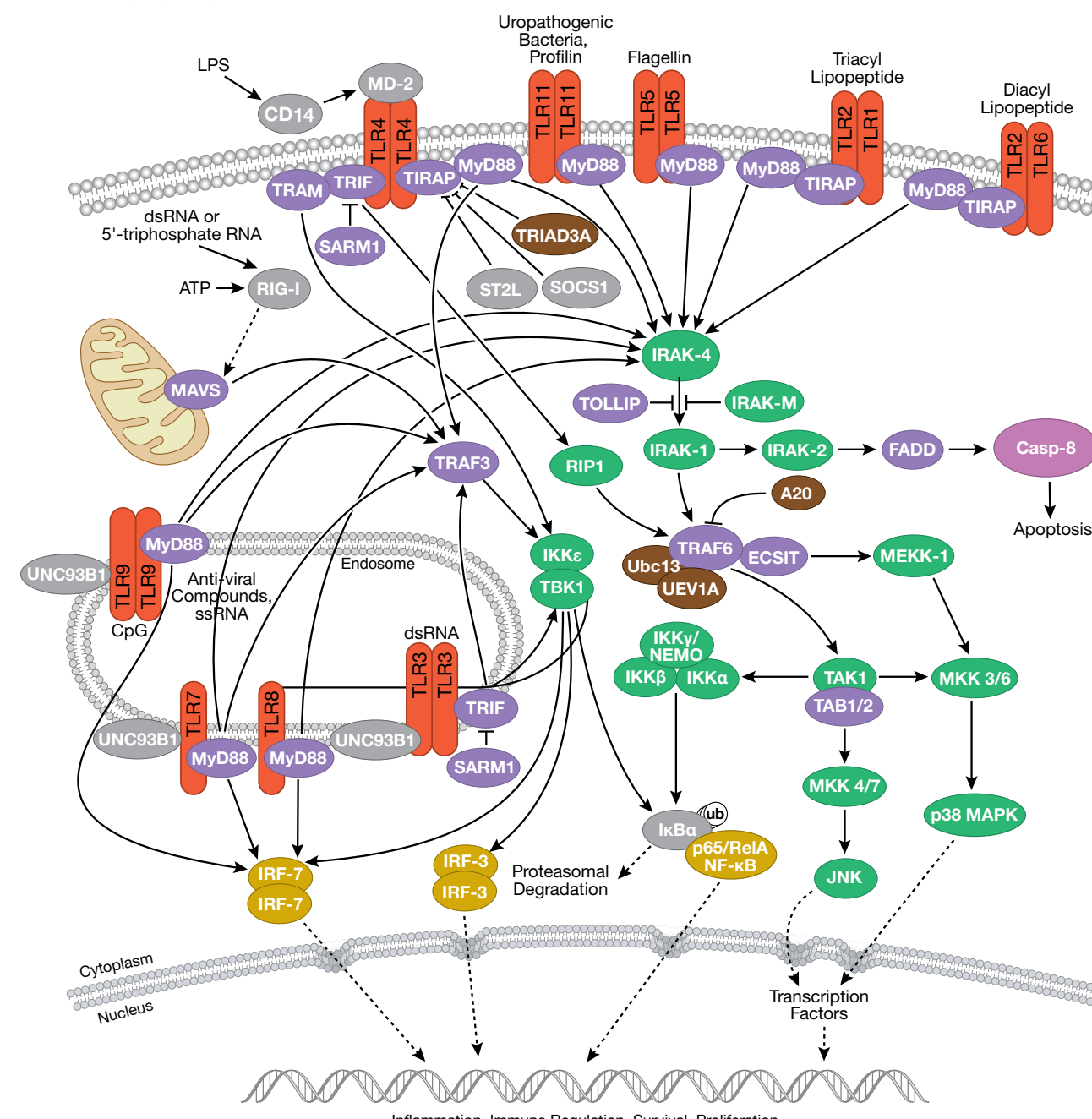
Jak/Stat Signaling: IL-6 Receptor Family

© 2002-2018 Cell Signaling Technology, Inc.



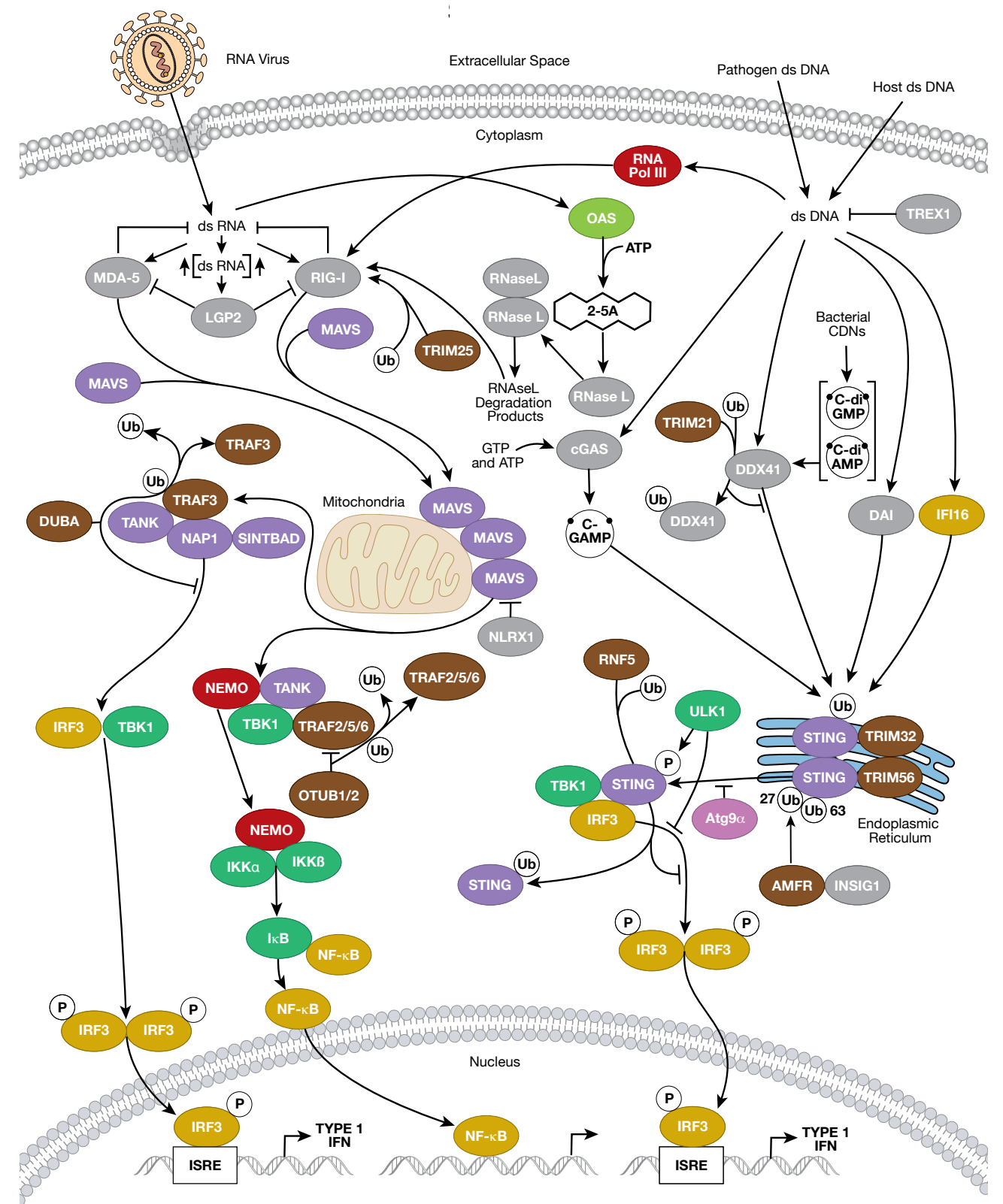
Toll-like Receptor Signaling

© 2009-2018 Cell Signaling Technology, Inc.



Cell Intrinsic Innate Immunity (STING) Signaling

© 2014-2018 Cell Signaling Technology, Inc.



NF-κB Signaling

© 2009-2018 Cell Signaling Technology, Inc.

