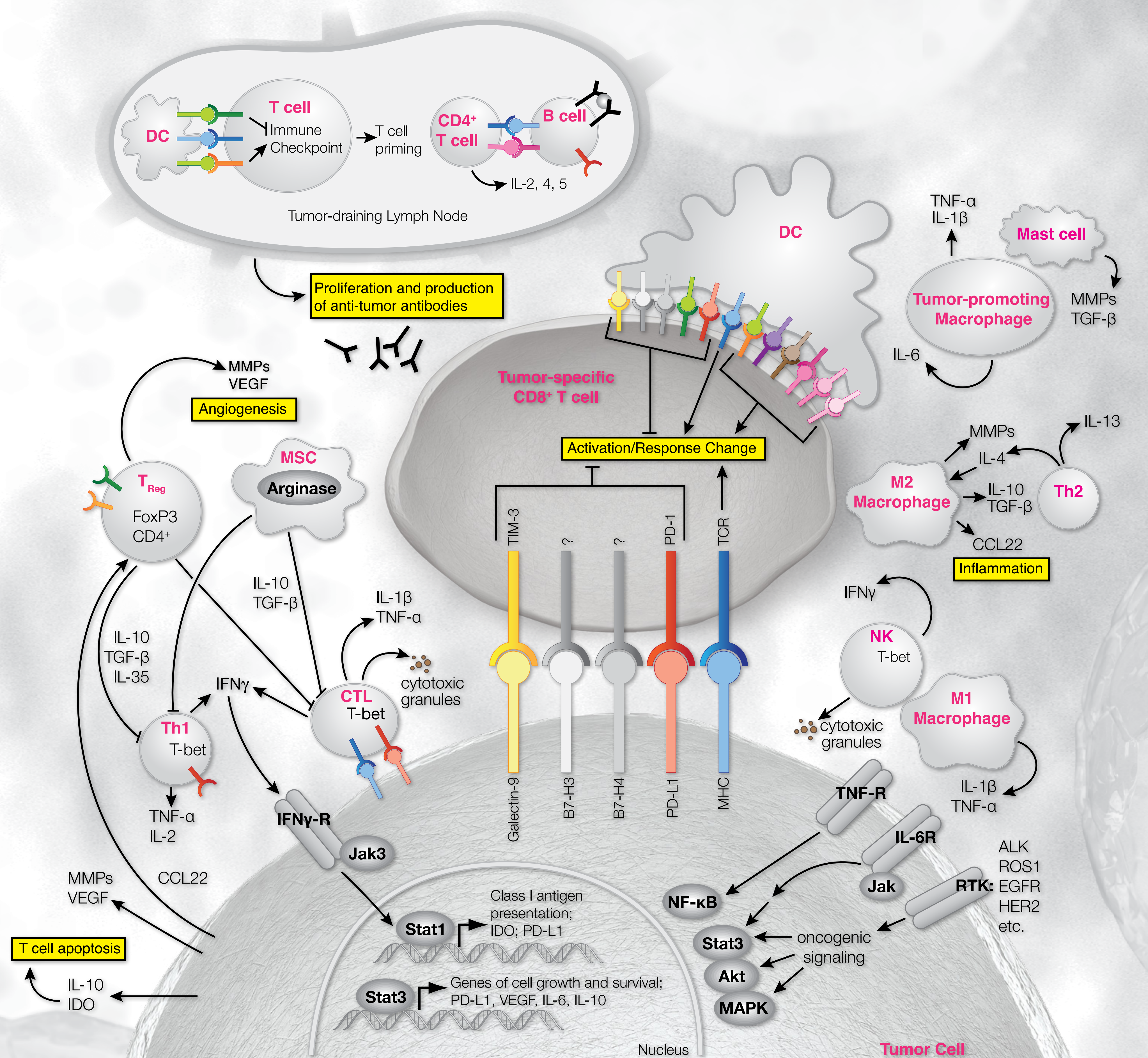
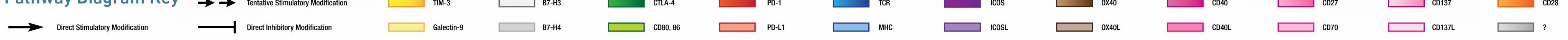


# Tumor Immunology



© 2014 Cell Signaling Technology, Inc.

## Pathway Diagram Key



Tumor cells employ multiple defense strategies to evade detection by the immune system. One common strategy, upregulation of immune checkpoint proteins and ligands, takes advantage of a natural immune mechanism for self-tolerance and prevention of collateral tissue damage (1,2). Immune checkpoint proteins, such as PD-1, CTLA-4, and many others, are located on T cells and engage with their corresponding ligand on tumor cells or dendritic cells, sending inhibitory signals that repress T cell activation or response (1,2). One of the first discovered checkpoint proteins, CTLA-4, plays a role at the stage of T cell priming by binding to the CD28 ligands CD80 or CD86 to prevent co-stimulatory signals necessary for T cell activation (1). In contrast, the PD-1/PD-L1 checkpoint acts later in the process, inhibiting anti-tumor immune responses by effector T cells such as CD4<sup>+</sup> T helper 1 (Th1) cells and CD8<sup>+</sup> cytotoxic T lympho-

cytes (CTLs), leading to decreases in IFN $\gamma$  production and cytolytic activity (1). Upregulation of PD-L1 expression on the tumor cell surface is mediated by IFN $\gamma$  signaling to Stat1, as well as oncogenic signaling through several receptor tyrosine kinases (EGFR, ALK, ROS, HER2, and others) to activate the MAPK, Akt, and Stat3 pathways (1-3).

Cells in the tumor microenvironment can also influence tumor progression. FoxP3<sup>+</sup>/CD4<sup>+</sup> T regulatory cells (T<sub>Reg</sub>) and myeloid-derived suppressor cells (MSCs) secrete immunosuppressive cytokines IL-10 and TGF- $\beta$  to inhibit the activity of Th1 cells and CTLs (1,2,4). Natural killer (NK) cells release cytotoxic granules against the tumor cell and secrete IFN $\gamma$ , which stimulates surrounding pro-inflammatory M1 macrophages (5). Pro-tumorigenic M2 macrophages sup-

press anti-tumor immune responses via production of IL-10 and TGF- $\beta$  and promote metastasis through release of MMPs (4,6). MMPs and TGF- $\beta$  are also released by surrounding mast cells (7).

### Select Reviews:

1. Pardoll, D.M. (2012) *Nat. Rev. Cancer* 12, 252-264.
2. Vanneman, M. and Dranoff, G. (2012) *Nat. Rev. Cancer* 12, 237-251.
3. Kawakami, Y., Yaguchi, T., and Park, J.H., et al. (2013) *Front. Oncol.* 3, 136.
4. Elinav, E., Nowarski, R., and Thaiss, C.A., et al. (2013) *Nat. Rev. Cancer* 13, 759-771.

5. Mentlik, J.A., Cohen, A.D., and Campbell, K.S. (2013) *Front. Immunol.* 4, 481.
6. Gajewski, T.F., Schreiber, H., and Fu, Y.X. (2013) *Nat. Immunol.* 14, 1014-1022.
7. Krstic, J. and Santibanez, J.F. (2014) *Scientific World Journal* 521754.

We would like to thank Girija Goyal, Susanne Baumeister, M.D., Karrie Wong, Ph.D., and Glenn Dranoff, M.D. (Dana-Farber Cancer Institute and Harvard Medical School) for reviewing this poster.