



T Cell Subsets

T helper 1 (Th1) Cells	Th2 Cells	Th9 Cells	Th17 Cells	Th22 Cells	Follicular Helper T (Tfh) Cells
Th1 cells promote cell-mediated immune responses and are required for host defense against microbial pathogens. Th1 cell cytokines promote macrophage activation, nitric oxide production, and cytotoxic T lymphocyte proliferation. When unregulated, Th1 responses may be involved in inflammatory disorders and autoimmunity.	Th2 cells are required for humoral immunity and play an important role in coordinating the immune response to large extracellular pathogens. Th2 cytokines also stimulate eosinophil activation and survival, mast cell activation, and have been implicated in the development of chronic allergic inflammation and asthma.	Th9 cells have a putative physiological role in host defense against parasitic helminth infections, but may also be involved in the development of chronic allergic inflammation, airway remodeling, and autoimmune diseases. They are closely related to Th2 cells, and there is considerable plasticity between the two cell types.	Th17 cells are involved in the immune response to specific fungi and extracellular bacteria. They stimulate chemokine secretion by resident cells, leading to the recruitment of neutrophils and macrophages to sites of inflammation. Recruited cells produce additional cytokines and proteases that can further promote the immune response.	Th22 cells express many of the same cell surface antigens as Th17 cells and primarily secrete IL-22. However, unlike Th17 cells, Th22 cells do not secrete IL-17, and they express the skin homing chemokine receptor CCR10. In the skin they may contribute to host defense against microbial pathogens and promote tissue repair or remodeling. Th22 cells may be involved in the pathogenesis of inflammatory skin disorders.	Tfh cells regulate the development of antigen-specific B cell immunity. Upon exposure to a foreign antigen, Tfh cells help centrocytes generate antibody-producing plasma cells and long-lived memory B cells. Tfh cells may be involved in the pathogenesis of immune cell-related cancers and autoimmunity.
Regulatory T (Treg) Cells	CD8 ⁺ Cytotoxic T Lymphocytes (CTL)	$\gamma\delta$ T Cells	Natural Killer T (NKT) Cells		
Treg cells play a crucial role in suppressing physiological immune responses and preventing sustained or exaggerated responses that can underlie inflammatory disorders and autoimmunity. They are best known for secreting cytokines that inhibit the activities of effector T cells and antigen-presenting cells.	CTLs are characterized by a TCR complex that includes an $\alpha\beta$ TCR and the co-receptor CD8 that recognizes peptide antigens presented by MHC I. CTLs are classically known for their direct interaction with cellular targets and the release of cytolytic granules containing a range of factors that induce apoptosis.	$\gamma\delta$ T cells are characterized by a TCR that consists of γ and δ subunits. Although traditional antigen presentation does not appear to be necessary for activation, phospho-antigens have been described as ligands for the $\gamma\delta$ TCR. Evidence suggests there are a wide range of subtypes that differ in their cytokine expression patterns and localization. Consequently, $\gamma\delta$ T cell activities are thought to be varied and context-dependent.	NKT cells express a semi-invariant $\alpha\beta$ TCR and surface antigens typically associated with natural killer cells. The TCR recognizes glycolipid antigens presented by the MHC I-like molecule CD1d. NKT cells can have divergent effects on immune function due to their ability to produce cytokines that promote either inflammation or tolerance. NKT cells may also be cytotoxic, but this is generally not thought to be their primary function.		

For more detailed information about the T cell subsets presented, please see our website at www.RnDSystems.com/TCellPoster

R&D
SYSTEMS®

For research use only. Not for use in diagnostic procedures. Note: The information in this poster should neither be considered comprehensive nor definitive. The particulars involved are understood to be subject to interpretation. © 2012 R&D Systems, Inc.