



Miltenyi Biotec

Get reliable results in your T cell differentiation culture

Polarizing T_H cells from naive CD4⁺ T cells

T helper cells – important subsets regulating immune responses

CD4⁺ T helper (T_H) cells play a central role in the adaptive immune system by controlling a variety of cellular responses, defending against intra-vesicular agents (T_H1) and external pathogens (T_H2). Their cytokine secretion suppresses or

stimulates immune responses and leads to antibody production of B cells, immunoglobulin class switch, and macrophage activation. Their crucial impact on immune responses and distinct role in disease, makes T_H cells a focus of many researchers studying immune regulation. The various T_H cell subsets can be differentiated from naive CD4⁺ T cells *in vitro* using specific combinations of cytokines (fig. 1).

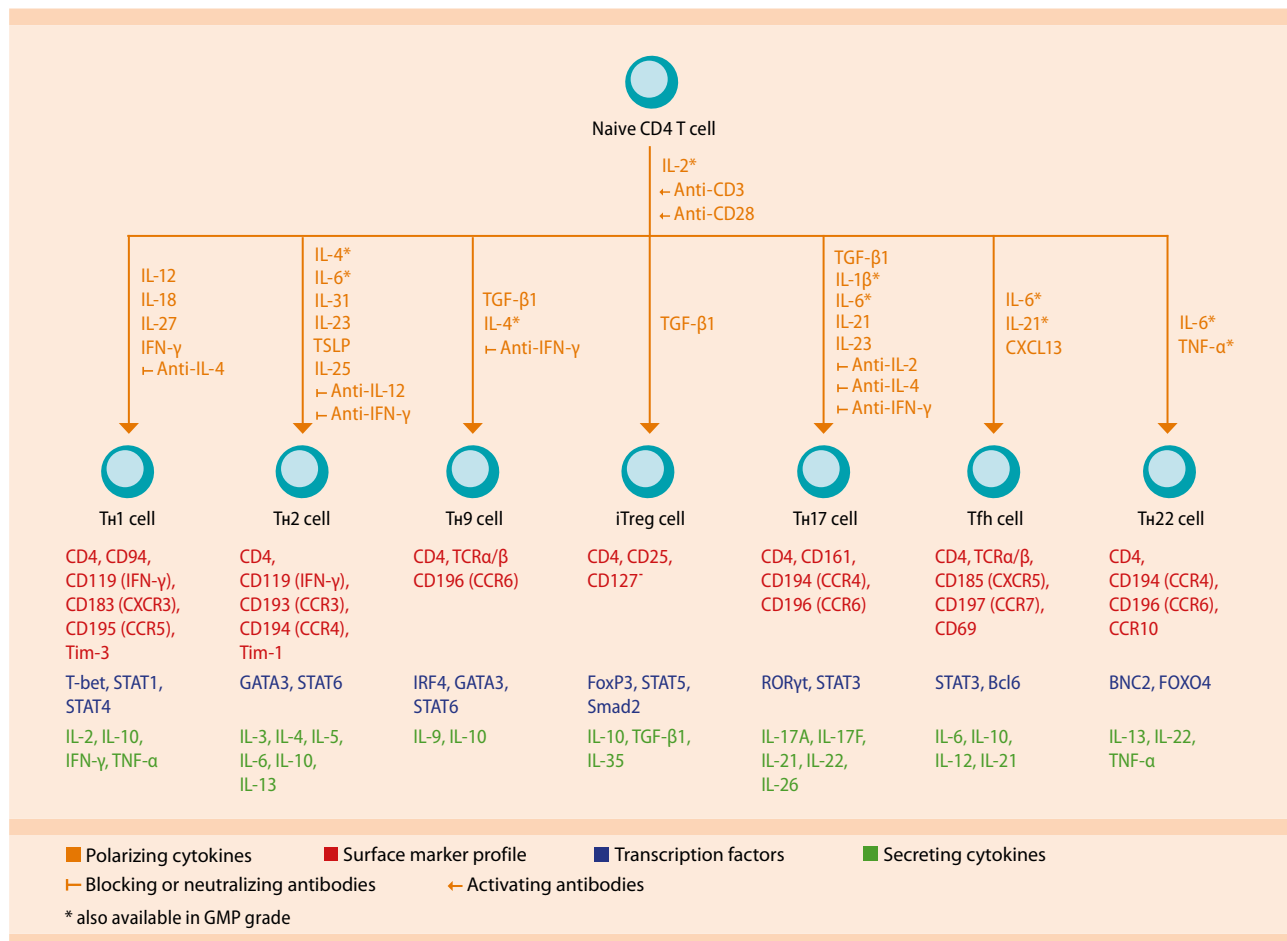


Figure 1: Polarization of naive CD4⁺ T cells into T_H cell subsets.

Polarizing T helper cells from naive CD4⁺ T Cells

Miltenyi Biotec provides the tools and support you need to differentiate T helper cells – from sample preparation, cell enrichment and cell culture, to analysis.

- Eliminate variability in cell culture with premium MACS® Cytokines
- Serum-free T cell culture medium for standardized cell culture conditions
- Comprehensive technical support from the experts in T cells

Start with gentle sample preparation

Get high yields of viable cells with preserved epitopes for your downstream applications. The gentleMACS™ Dissociators with MACS Tissue Dissociation Kits enable gentle and effective dissociation of many tissue types, including tumor and skin.



Figure 2: The gentleMACS Family for standardized tissue dissociation or homogenization.

Isolate highly pure naive CD4⁺ T cells

In about 18 minutes, you can have pure, viable naive CD4⁺ T cells for high-quality results. Automate the process with the autoMACS® Pro Separator for greater convenience and reproducibility. Need to process multiple samples in parallel? Choose the MultiMACS™ Cell24 Separator Plus.



Figure 3: The MultiMACS Cell24 Separator Plus and autoMACS Pro Separator enable automated multisample cell separation for reproducible results.

Standardize your cell culture

Increase reproducibility in your polarization experiments with TexMACS™ Medium, an optimized serum-free T cell culture medium available in research and GMP grade. Get consistent cell growth and eliminate lot-to-lot variability with premium MACS Cytokines to ensure identical unit dosage.



Figure 4: TexMACS Medium and MACS Cytokines enable consistent polarization of naive CD4⁺ T cells into Th subsets.

Sensitive flow cytometry

Detect and enrich viable cytokine-secreting cells without fixation of dead cells by using MACS Cytokine Secretion Assays. For fast cytokine detection, check out the Rapid Cytokine Inspector Kits. Also, detect multiple cytokines in a single sample with MACSplex Cytokine Kits and the MACSQuant® Analyzer 10. Our flow cytometer's pre-installed Express Mode enables full automation of acquisition, gating strategy, and analysis.

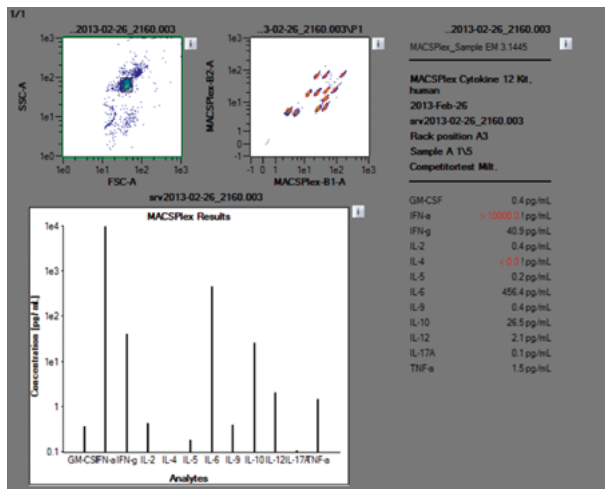


Figure 5: MACSplex Cytokine Kit analysis template using the MACSQuant Analyzer's Express Mode configuration.

See the whole solution for polarizing T helper cells from naive CD4⁺ T cells at

► miltenyibiotec.com/polarizing



Miltenyi Biotec Inc. | Phone 800 FOR MACS | Phone +1 530 888 8871 | Fax +1 877 591 1060 | macs@miltenyibiotec.com | www.miltenyibiotec.com
Miltenyi Biotec provides products and services worldwide. Visit www.miltenyibiotec.com/local to find your nearest Miltenyi Biotec contact.

Unless otherwise specifically indicated, Miltenyi Biotec products and services are for research use only and not for therapeutic or diagnostic use. autoMACS, gentleMACS, MACS, MACSQuant, MultiMACS and TexMACS are either registered trademarks or trademarks of Miltenyi Biotec GmbH. Copyright © 2015 Miltenyi Biotec GmbH. All rights reserved.