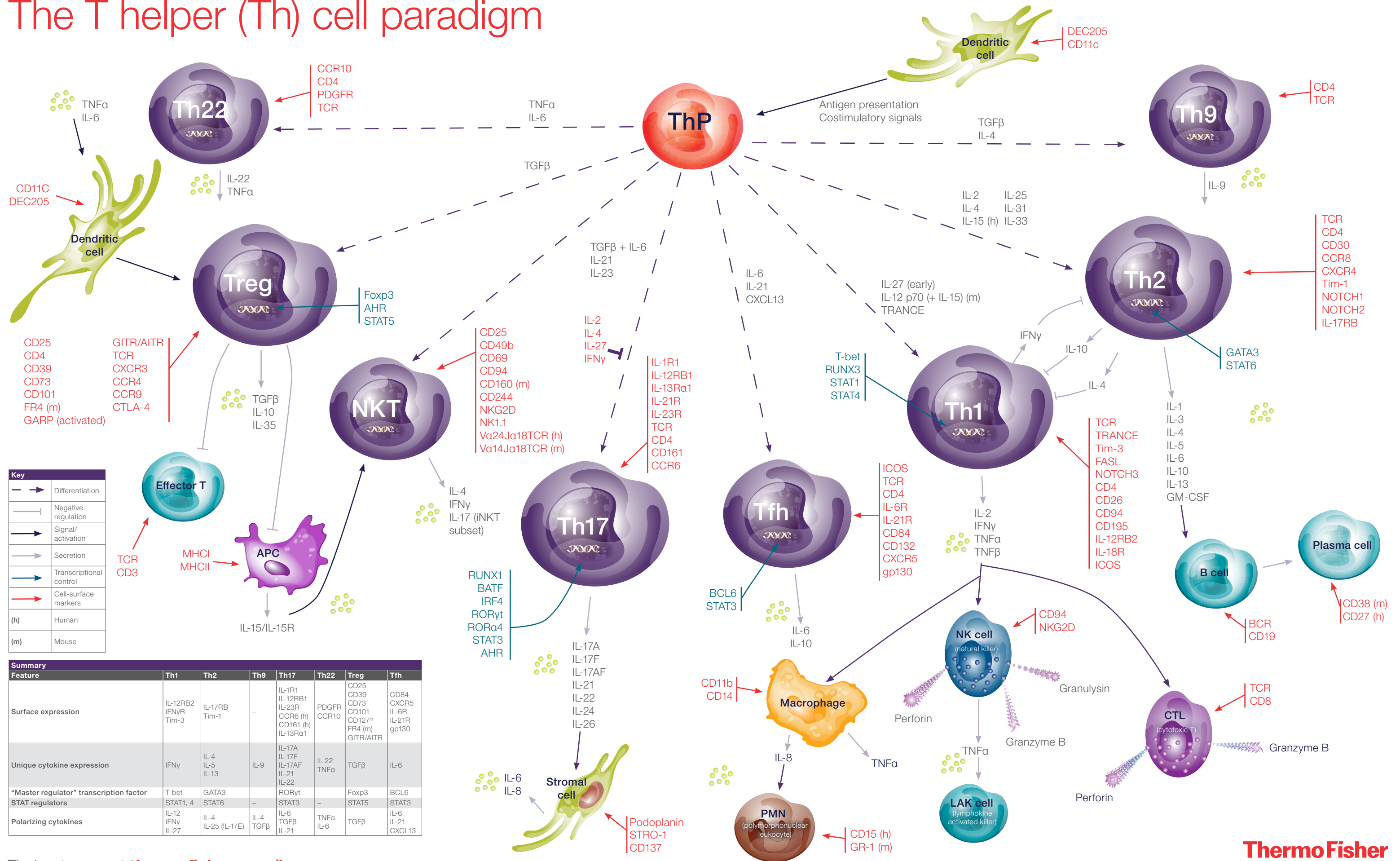


# The T helper (Th) cell paradigm

Naive Th cells play a central role in modulating the immune response. They are activated by recognition of a peptide antigen bound to the class II major histocompatibility complex (MHC) on antigen-presenting cells (APCs) through the interaction with the T cell receptor. After activation, Th precursor (ThP) cells begin to divide and give rise to effector cells. These effector Th cells are CD4<sup>+</sup> and can be divided into three main types with distinct cytokine secretions and functions: Th type 1 (Th1), Th type 2 (Th2), and most recently Th type 17 (Th17), expanding the group to what is now referred to as the Th1/Th2/Th17 paradigm.

# The T helper (Th) cell paradigm



Find out more at [thermofisher.com/immunoassay](https://thermofisher.com/immunoassay)

# invitrogen

Find out more at [thermofisher.com/immunoassay](https://www.thermofisher.com/immunoassay)

**For Research Use Only. Not for use in diagnostic procedures.** © 2020 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. **COL23856 0220**

**ThermoFisher**  
SCIENTIFIC