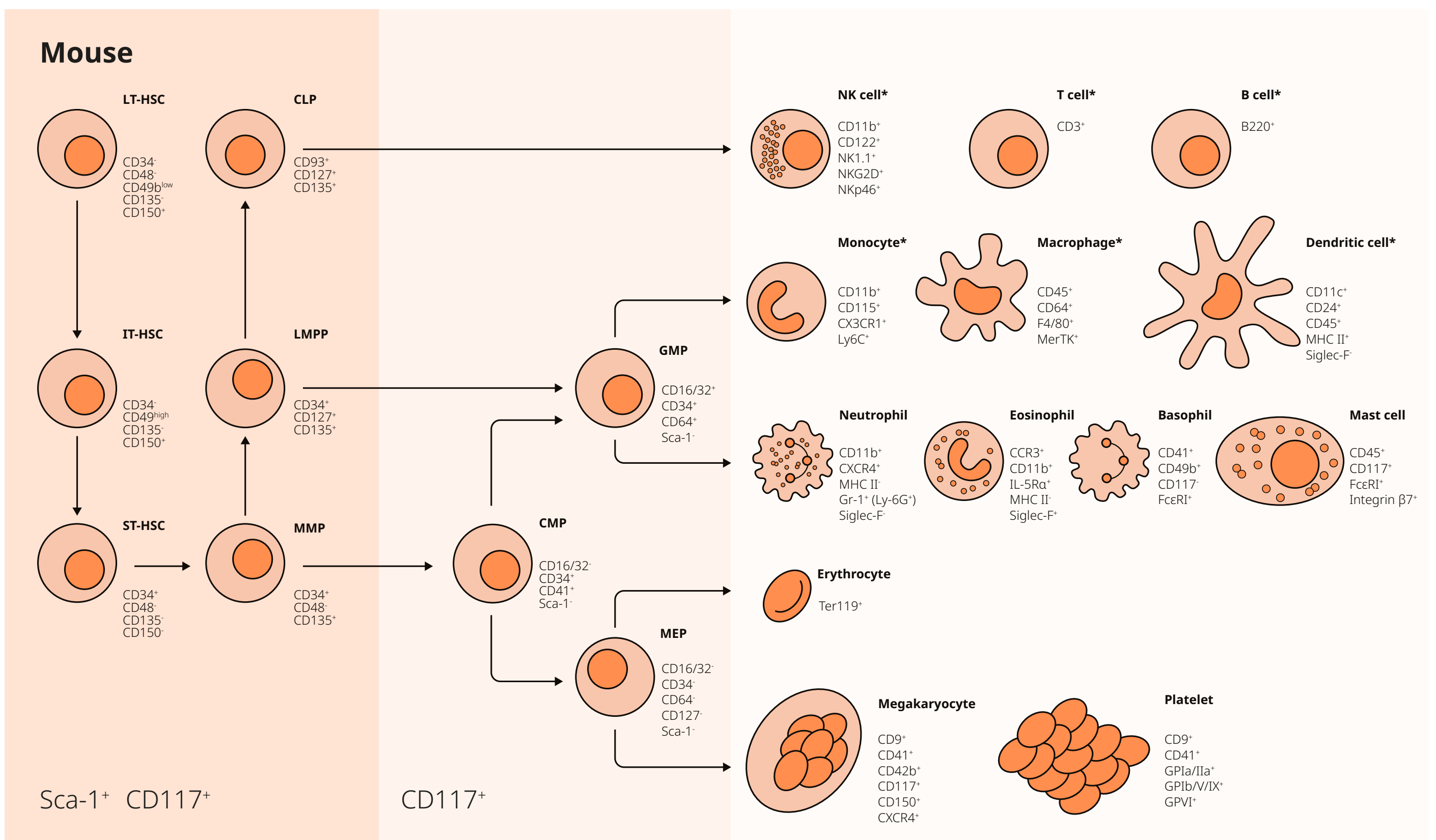
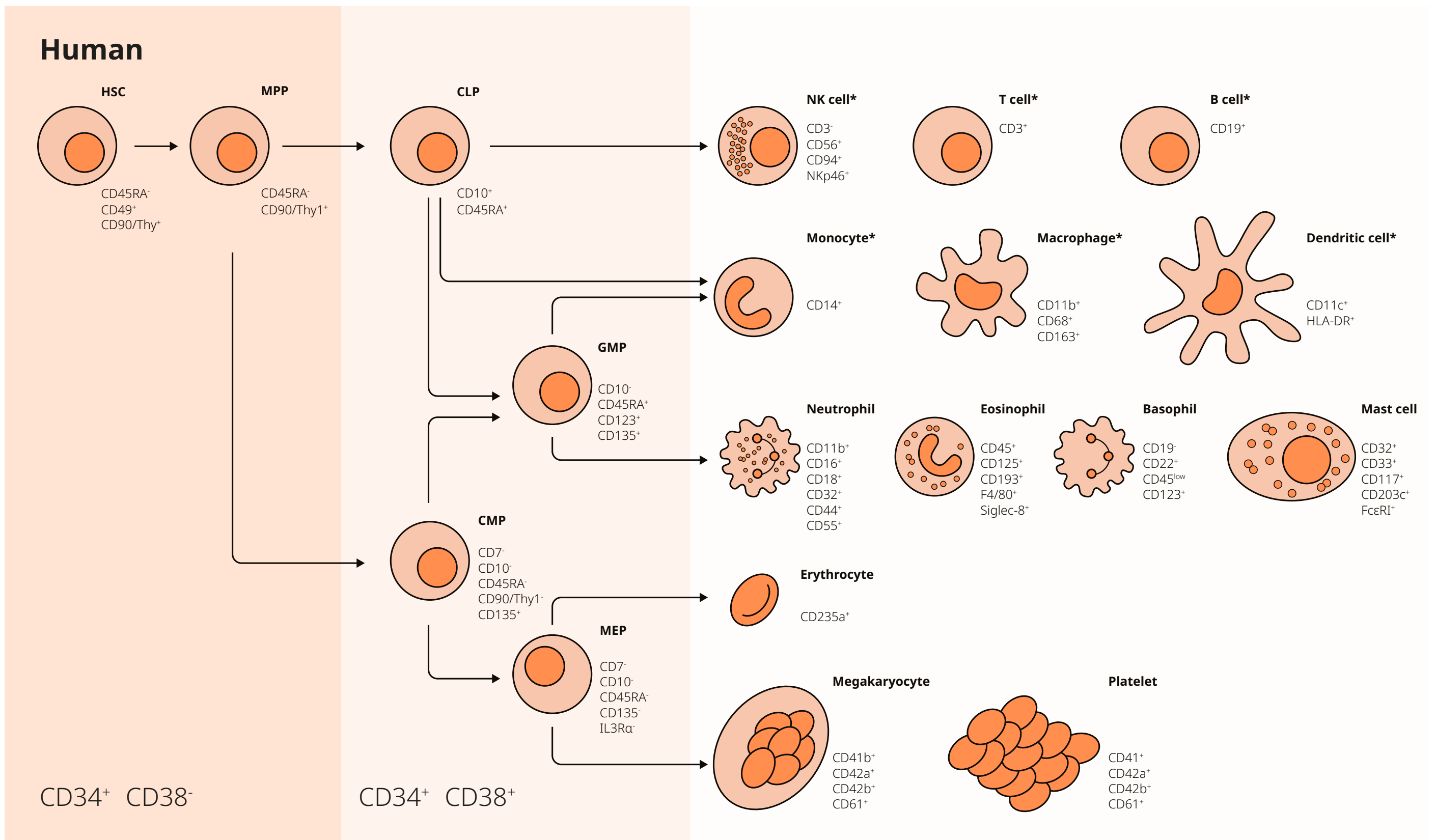


Immune cell markers



abcam.com/immune-cell-markers-poster

Markers are accompanied with a "+" or a "-" symbol to indicate whether the specific cell type expresses or lacks the antigen (eg CD10⁺ indicates a cell type expressing CD10). In some cases, markers are accompanied with "high" or "low", indicating different degrees of expression.
*The markers selected for these cell types are common to all subsets (eg T cell markers are common to killer, helper, and regulatory T cells).

progress happens together
abcam