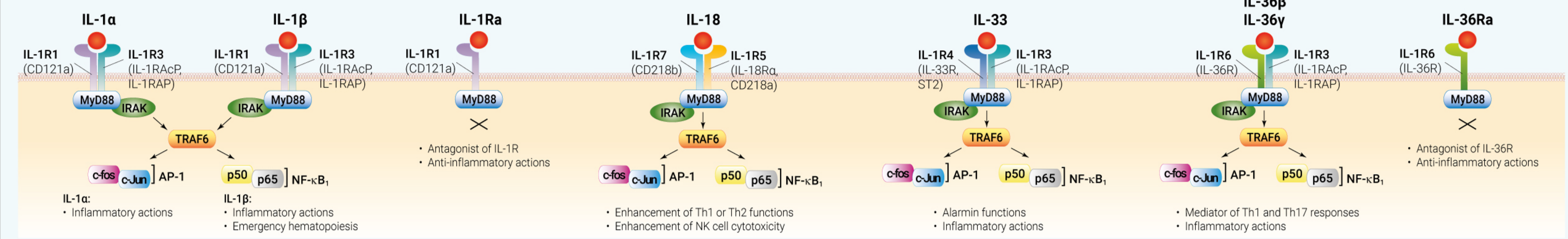
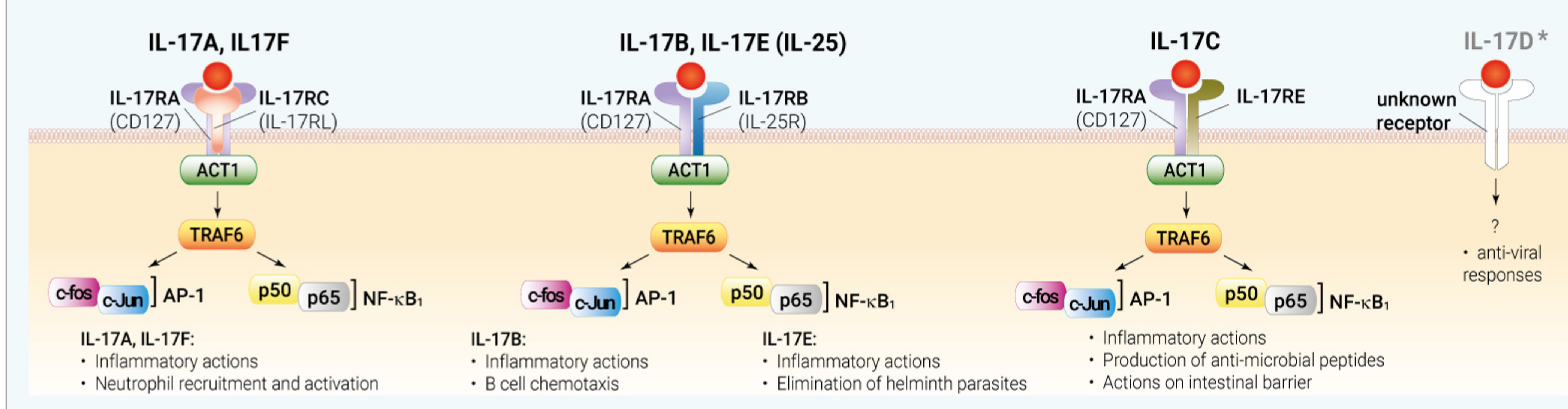


IL-1, IL-17, TGF, and TNF-related cytokine receptor families

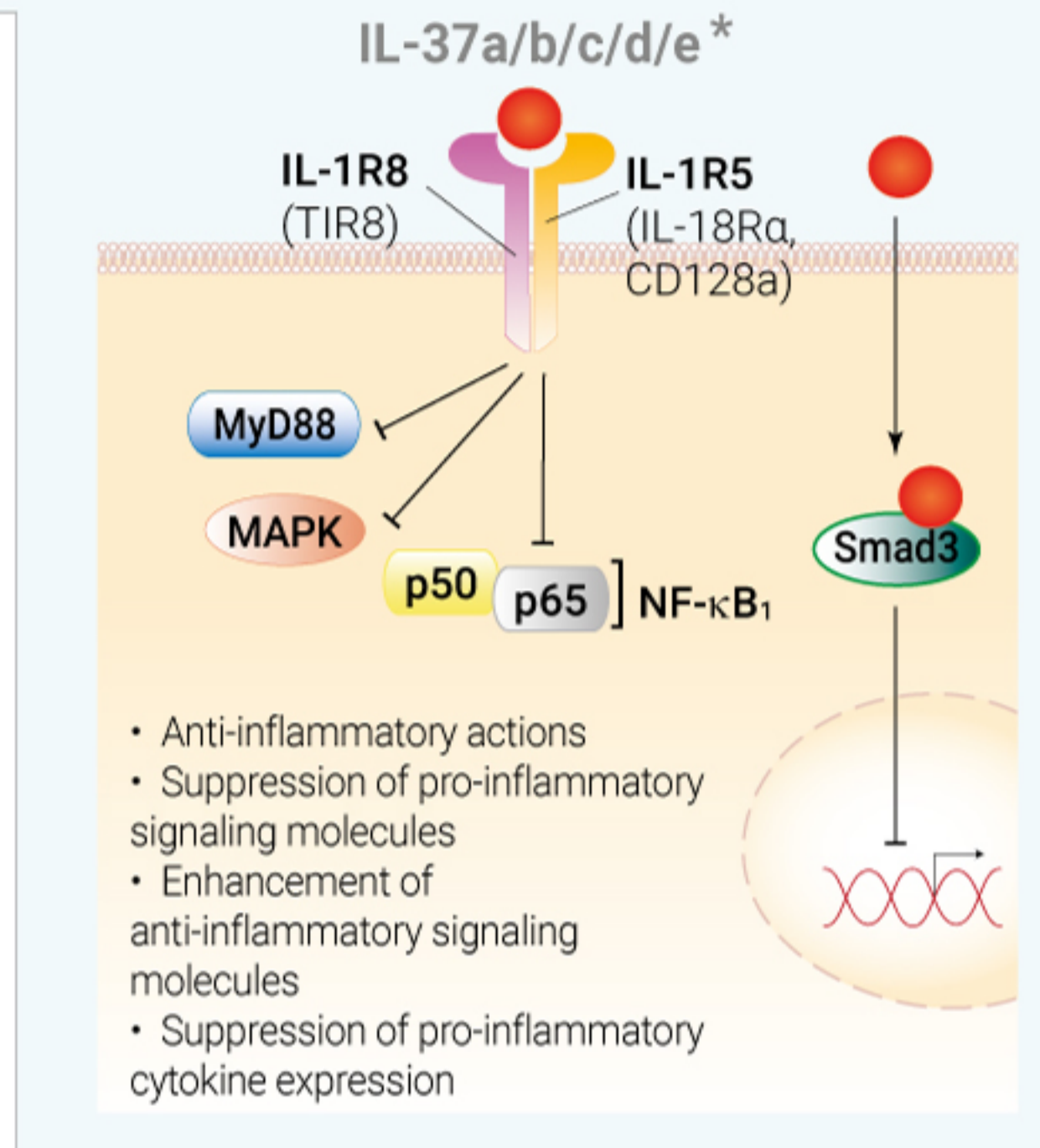
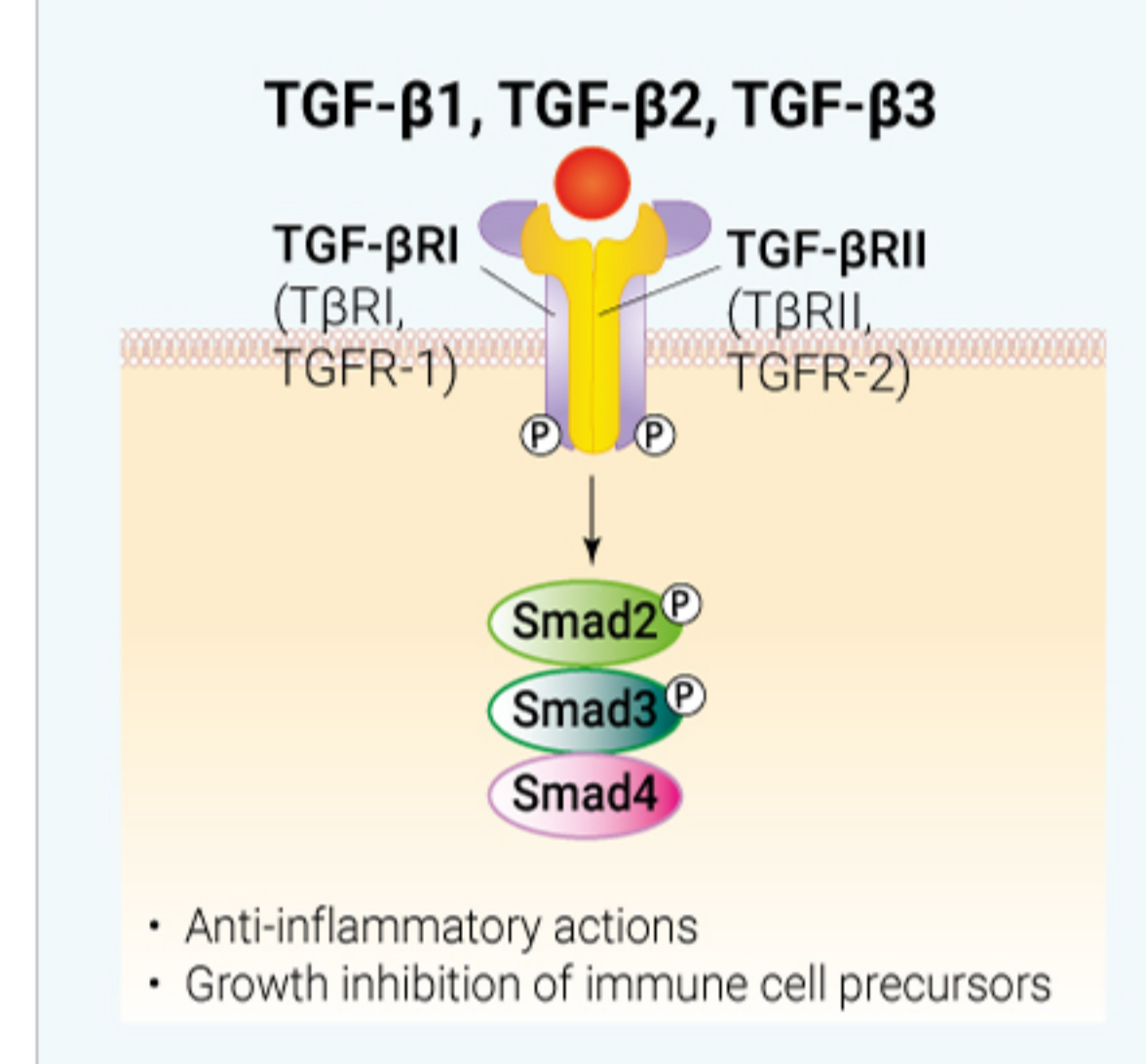
IL-1 / Toll-Like receptor cytokine family



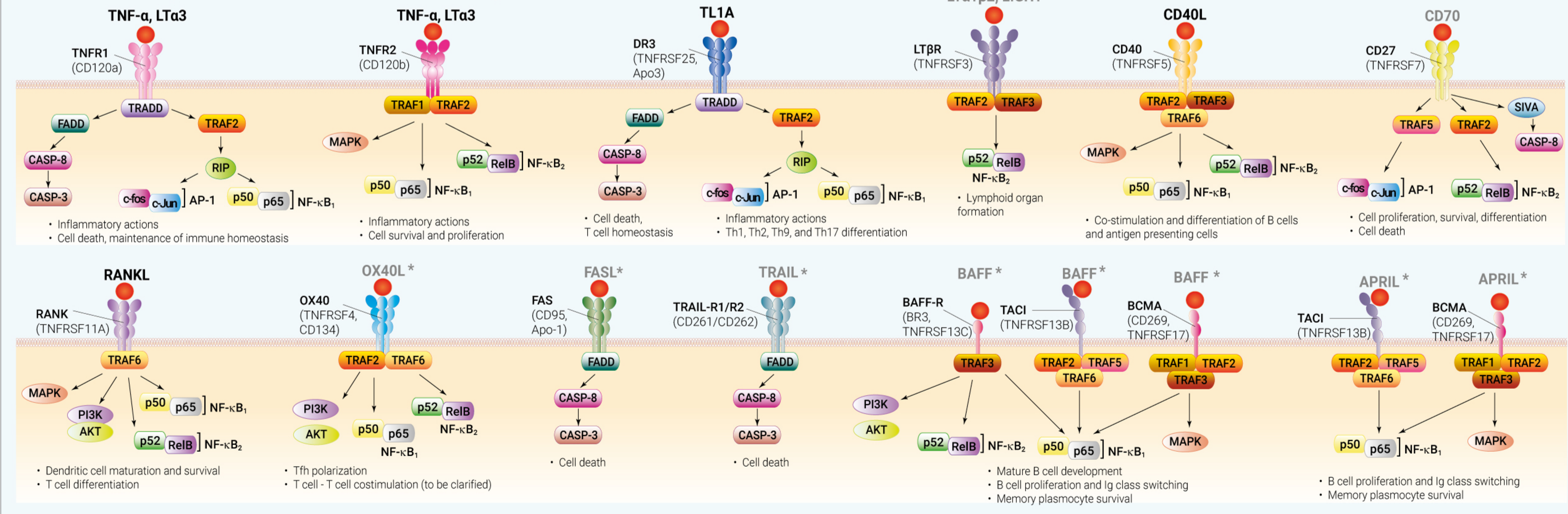
IL-17 cytokine receptor family



TGF cytokine family

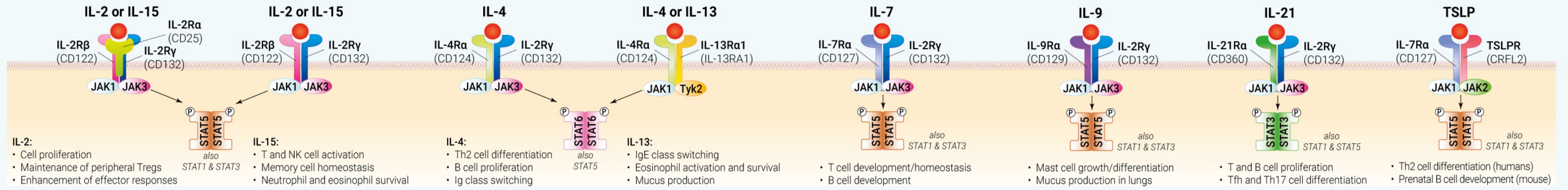


TNF-related cytokine receptor superfamily (Note: for clarity purposes, the transmembrane form of ligands is not represented)

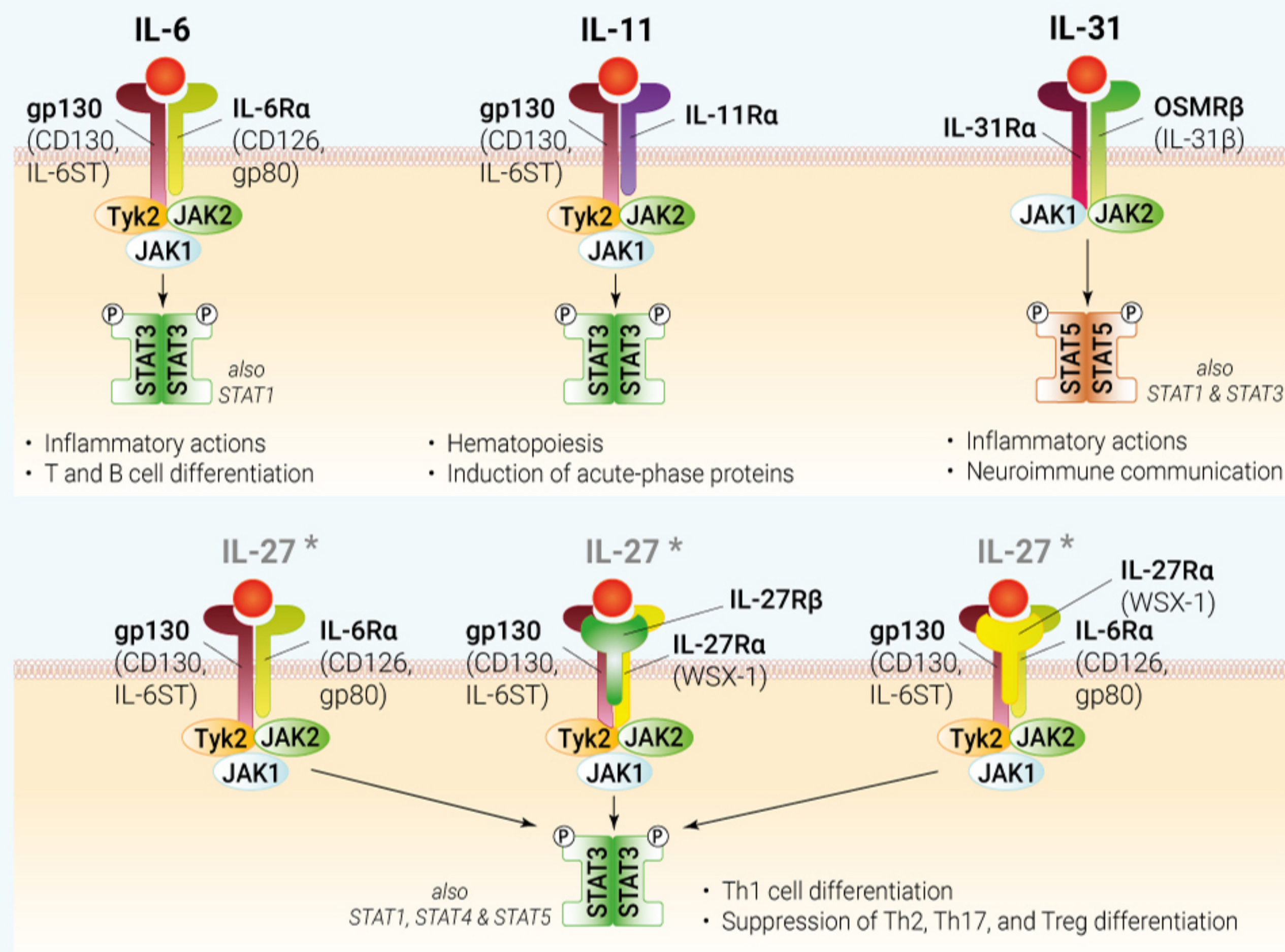


InvivoGen's Cytokine Reporter Cell Lines

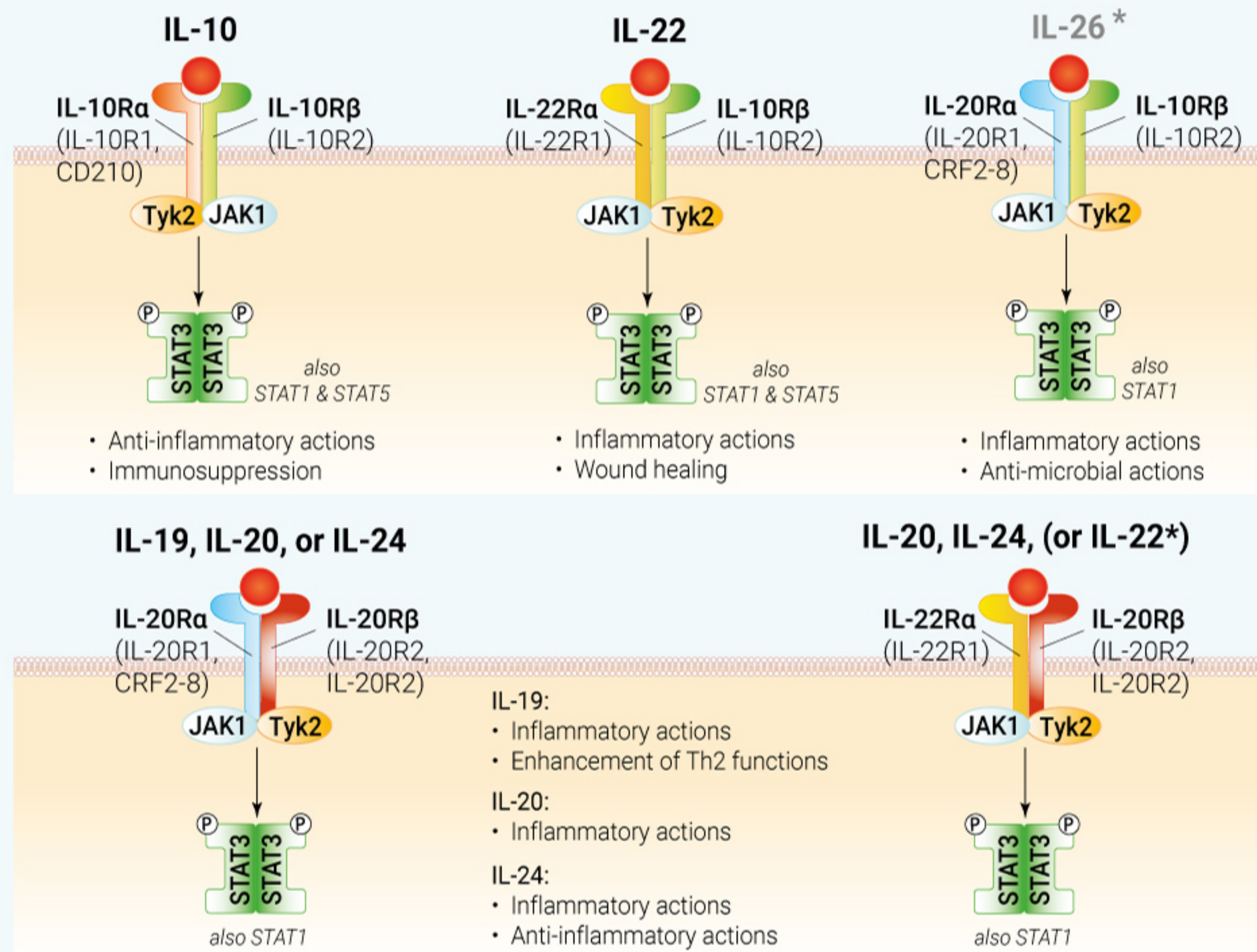
IL-2 / Common γ c cytokine family



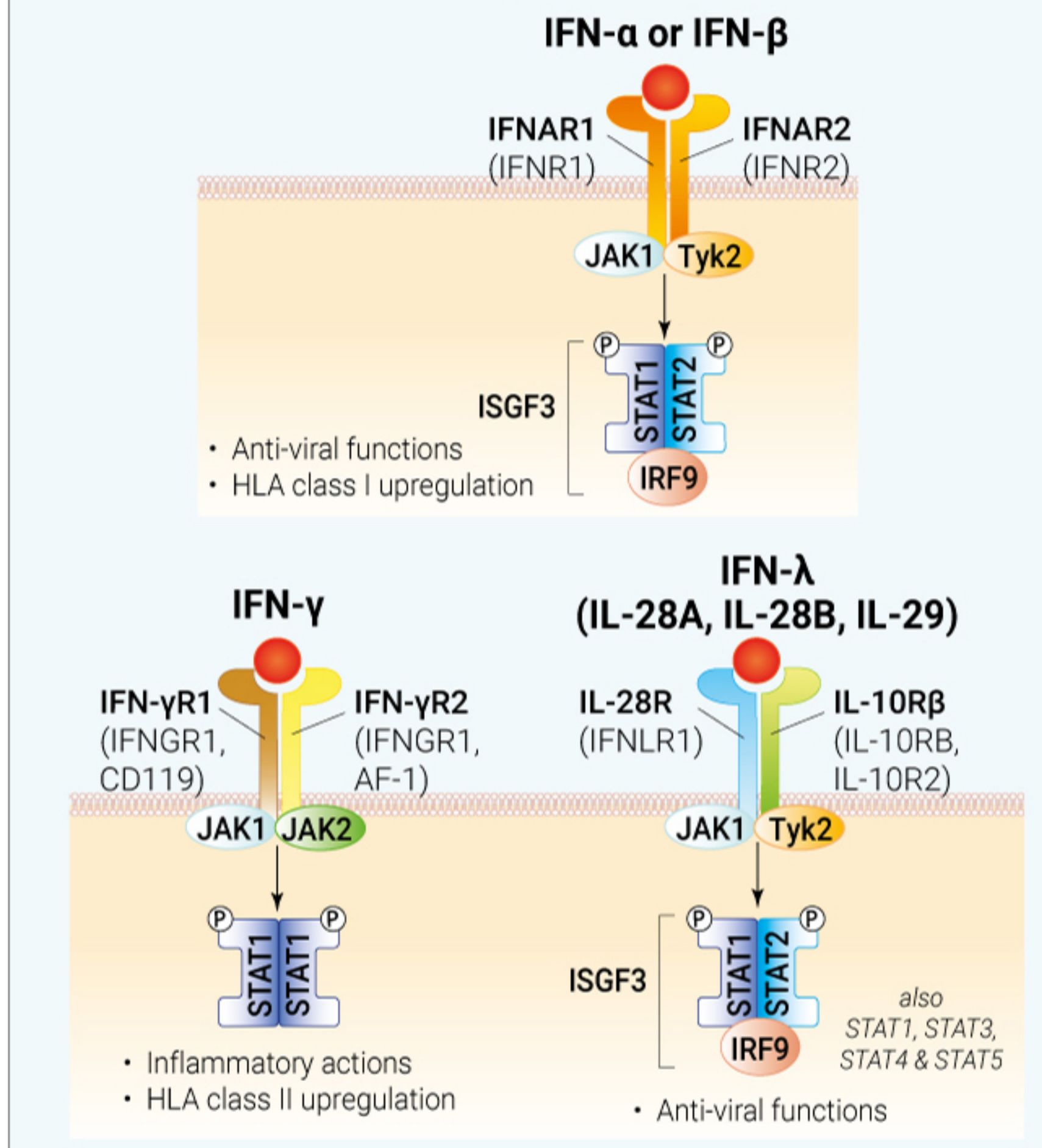
gp130 cytokine family



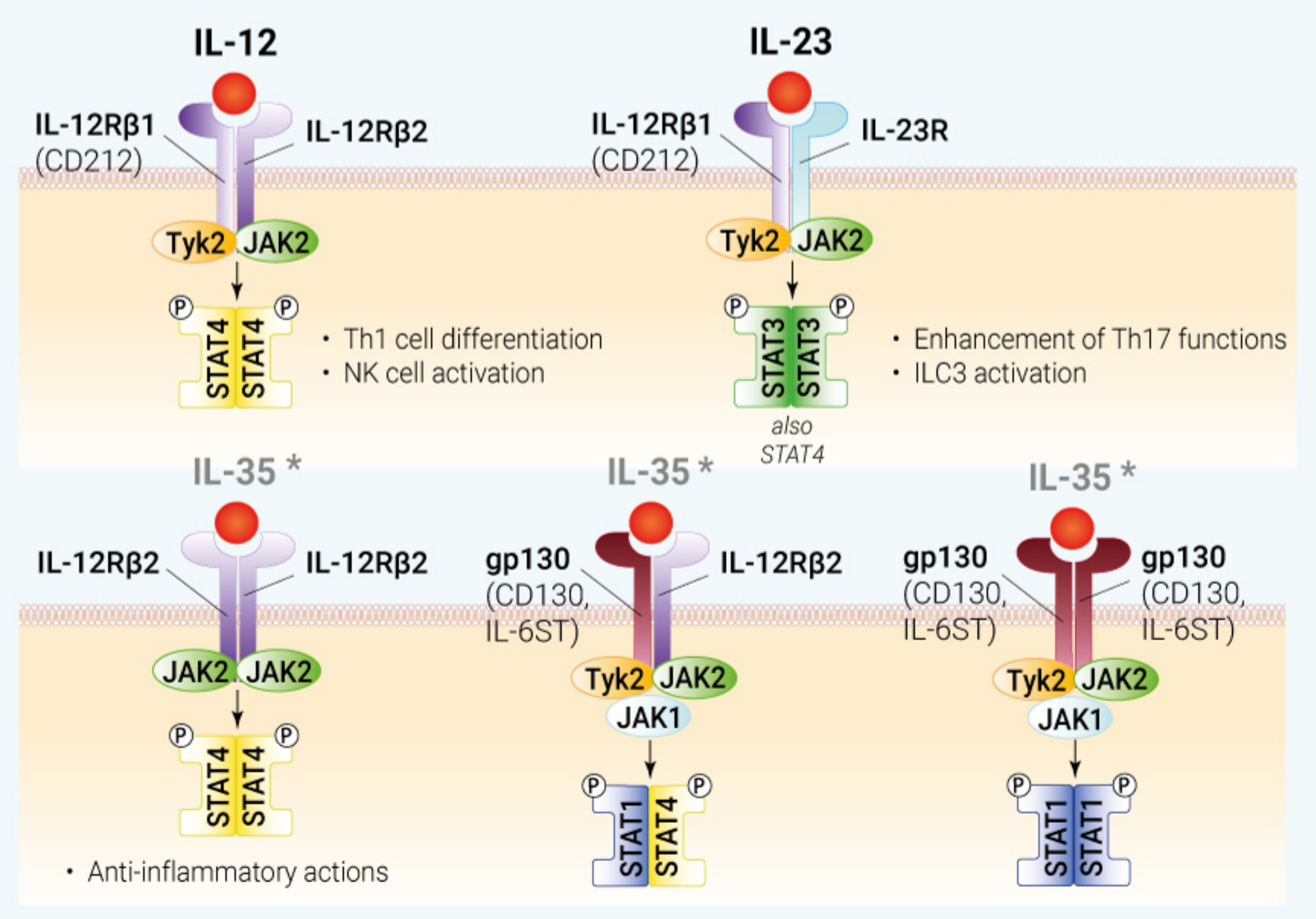
IL-10 cytokine family



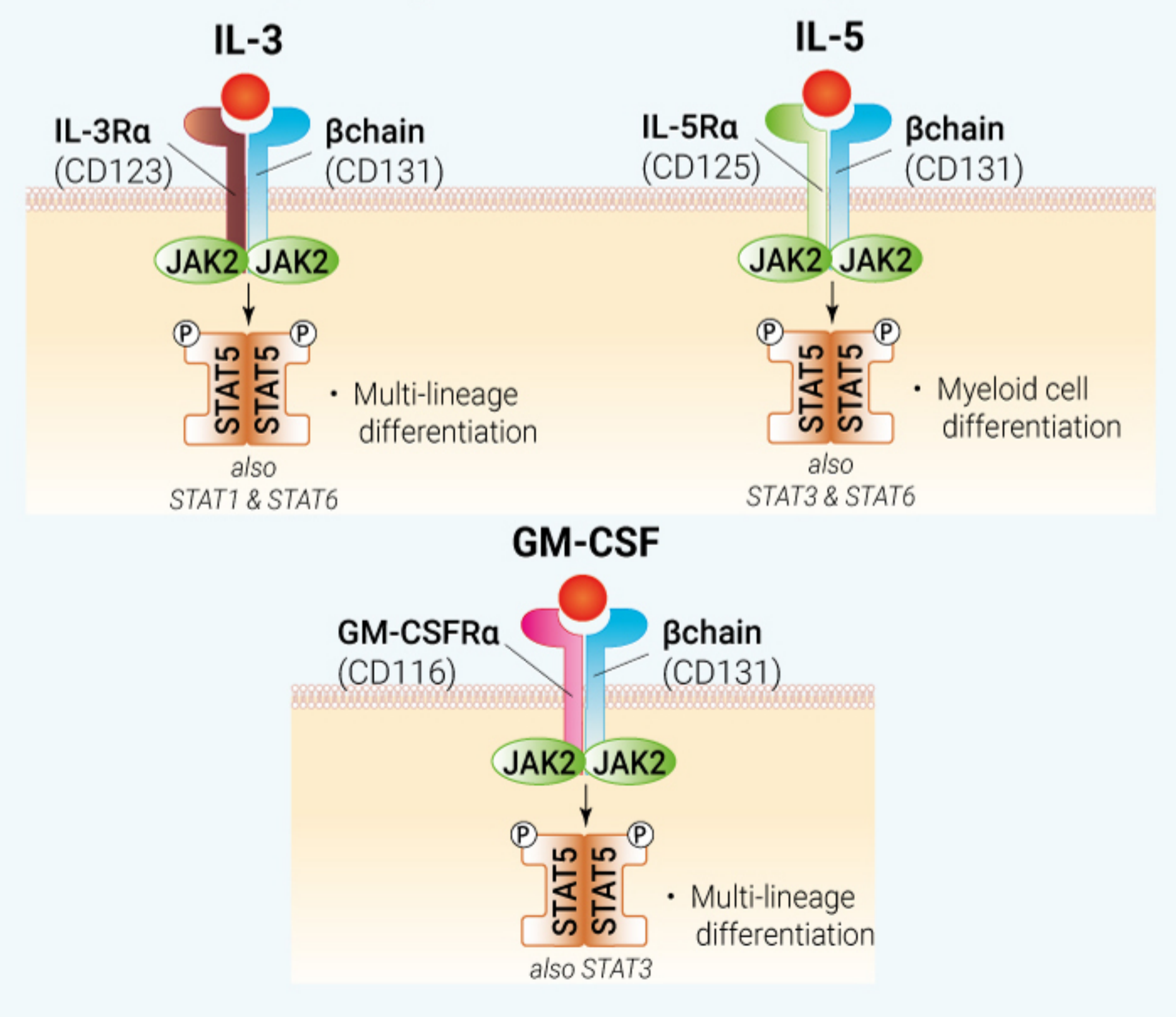
Interferon family



IL-12 cytokine family



Common β c cytokine family



Hormone-like cytokine family

