



Erythropoietin (Epo) and its receptor [pdb 1EBA]: Epo (yellow, lime) and Epo receptor extracellular domains (blue, ice blue).