



Epidermal growth factor (EGF) and its receptor [pdb 1M14 and 1IVO]: EGF (orange); and EGF receptor intracellular and extracellular domains (pink).