

IL3 Receptor Alpha and Common Beta Subunits

