



RNA polymerase II complex [pdb 1NT9]: Rpb2 subunits 1-12.