



Left [pdb 1PNY]: bacterial 30S ribosome with 19 protein subunits. Right [pdb 1PNX]: bacterial 50S ribosome with 31 protein subunits. RNA backbone (purple) and nucleotides (blue).