



Tumour suppressor protein p53 [pdb 1TUP] (yellow, lime, orange) complexed with DNA. DNA backbone (purple) and nucleotides (blue).