



www.kinexus.ca



Nucleosome core particle [pdb 1KX3]: DNA backbone (purple) and nucleotides (blue); Histone H2A.1 (green, lime); Histone H2B.2 (yellow, tan); Histone H3 (pink, red); & Histone H4 (orange, ochre).