



Left [pdb 1I51]: cleaved Caspase-7 p20 (ice blue) and p10 (orange) chains interacting with the Xenopus inhibitor of apoptosis protein Xiap-Bir2 (lime). Right [pdb 1FIJ]: cleaved Caspase-7 p20 (ice blue) and p10 (orange) chains with the substrate peptide acetyl-Asp-Glu-Val-Asp-Cho.