



www.kinexus.ca



Death receptor/Trail complex [pdb 1DOG]: Death receptor 5 extracellular domains (yellow, orange & ochre); & Trail/Apo2 ligands (ice blue, purple, mauve).

KINEXUS