



Bone morphogenetic protein (BMP) and its receptor [pbd 1ES7]: BMP-2 (lime, green); and BMP receptor 1A extracellular domains (ochre, tan).