



Foreground [pbd IJSU]: Cyclin-dependent kinase (CDK2) (pink); Cyclin A (blue); & p27 KIP1 CDK inhibitor (yellow).
Background [pdb IFQ1: CDK phosphatase [KAP] (lime); & CDK2 (pink).