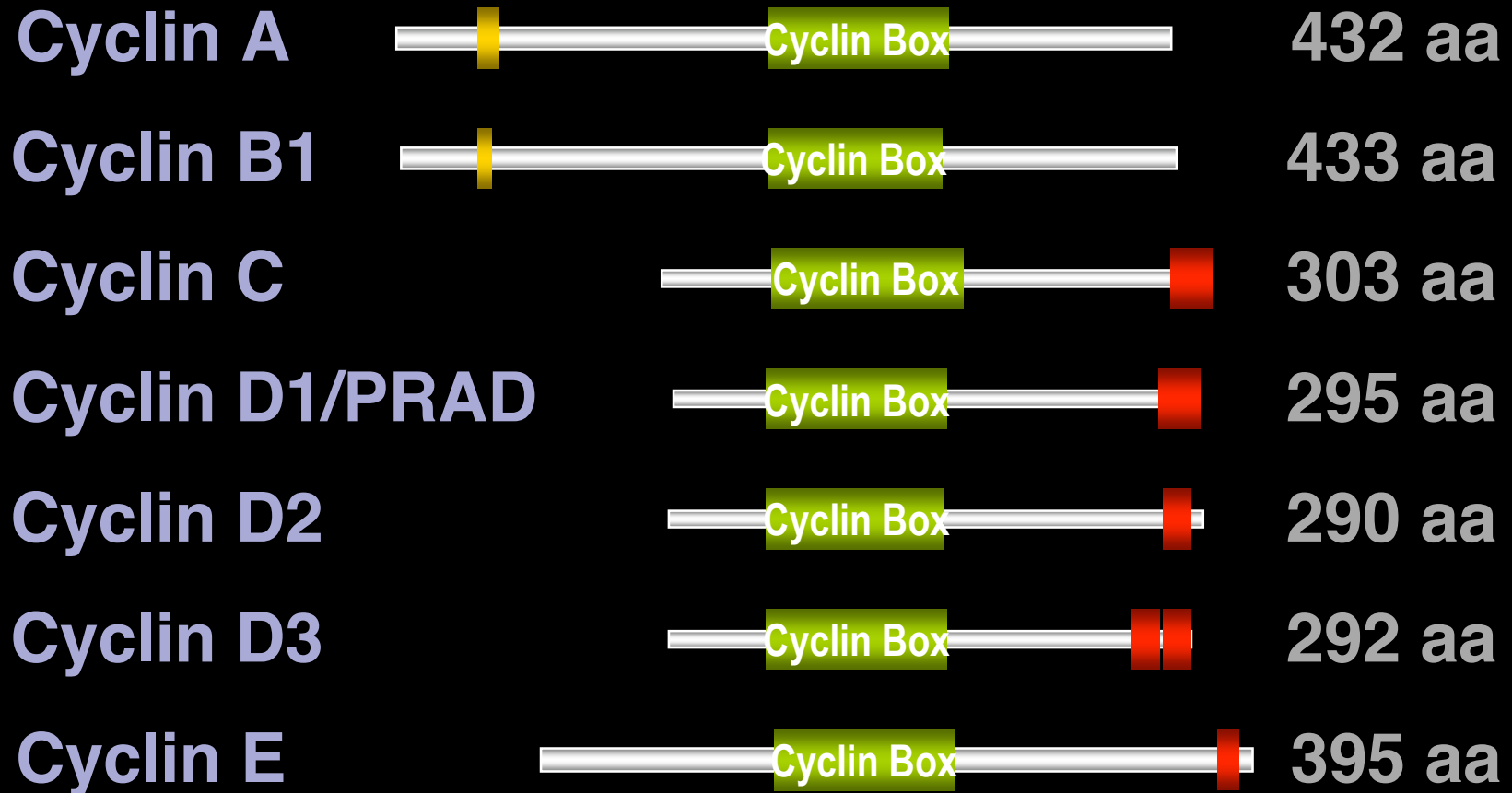


Domain Structures of Human Cyclins



PEST