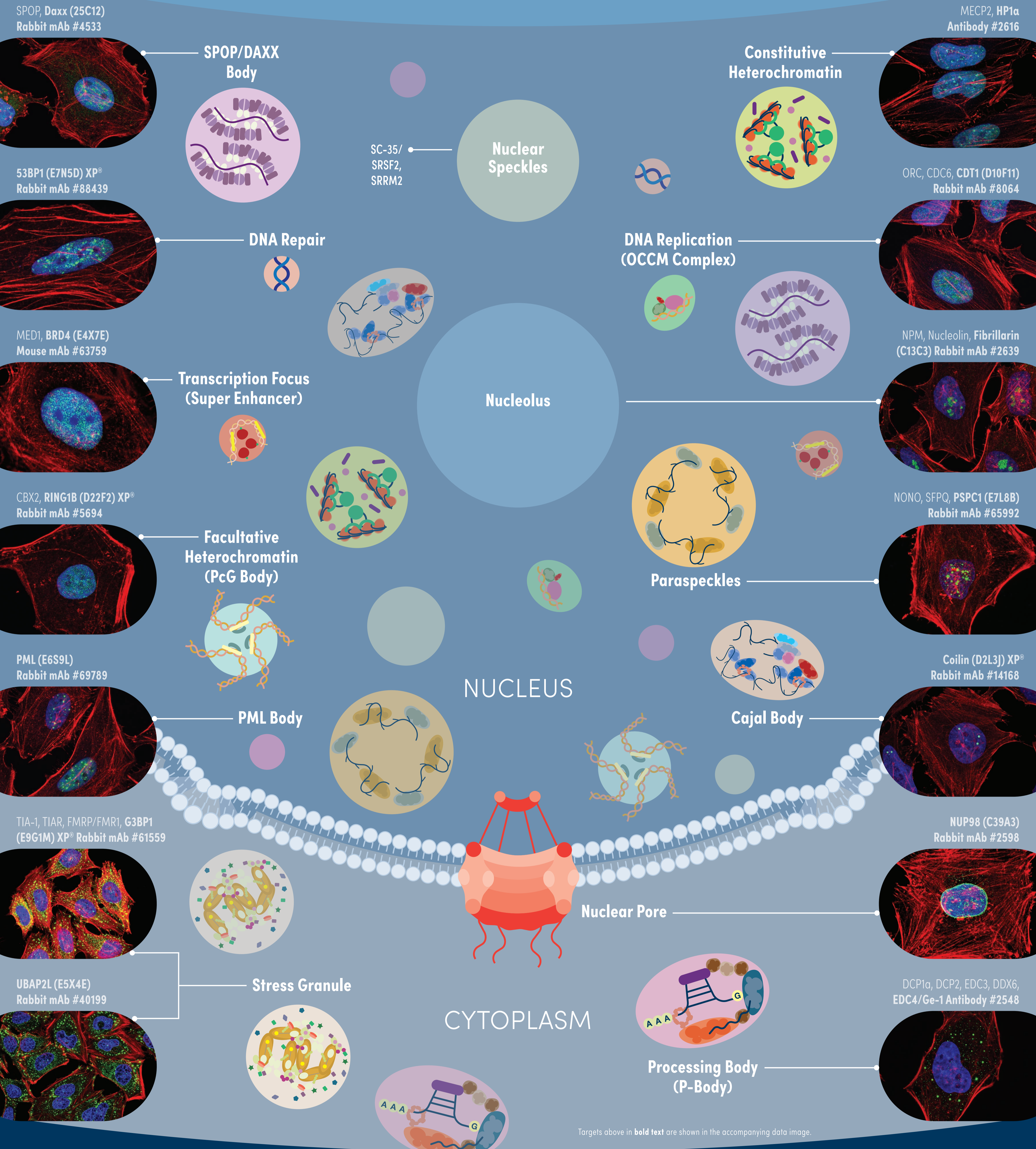


Biomolecular Condensate Markers

Biomolecular condensates, or membraneless organelles (MLOs), are involved in a myriad of cellular processes. MLOs are characterized by compartmentalization and concentration of proteins and nucleic acids, but without a lipid membrane. MLOs function across numerous cellular pathways, such as transcription and translation, DNA damage response, organelle biogenesis, and signal transduction; as such, they are highly varied in composition. This provides new opportunities

for deeper study; MLO proteins act as markers allowing researchers to study MLOs and their dynamics in more detail, while simultaneously enabling researchers to determine if other proteins of interest function in these MLOs. As these MLOs and their components are still being identified and characterized, presented here are routinely used markers for MLOs and antibodies available from Cell Signaling Technology (CST) for their study.



Targets above in **bold text** are shown in the accompanying data image.