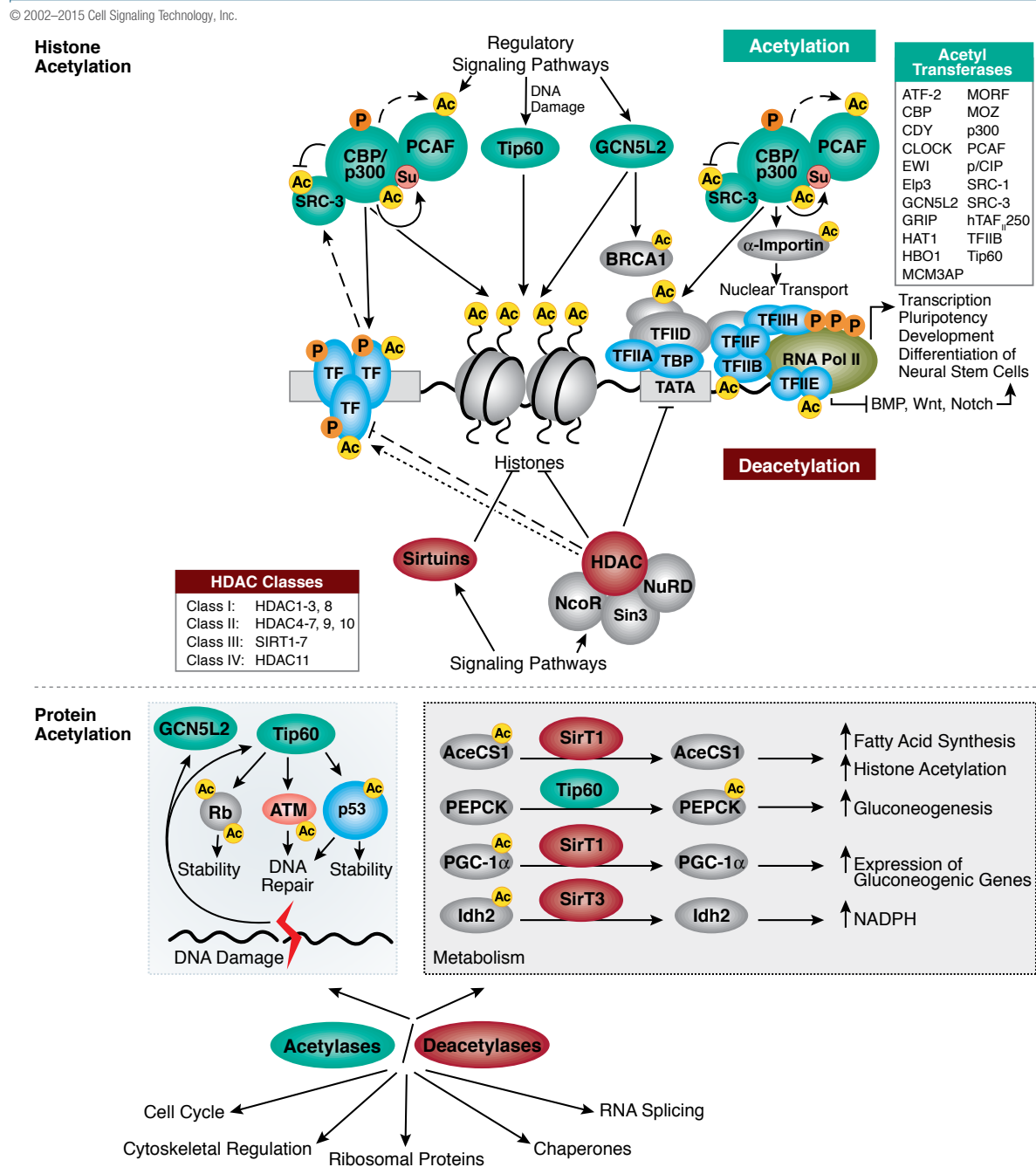
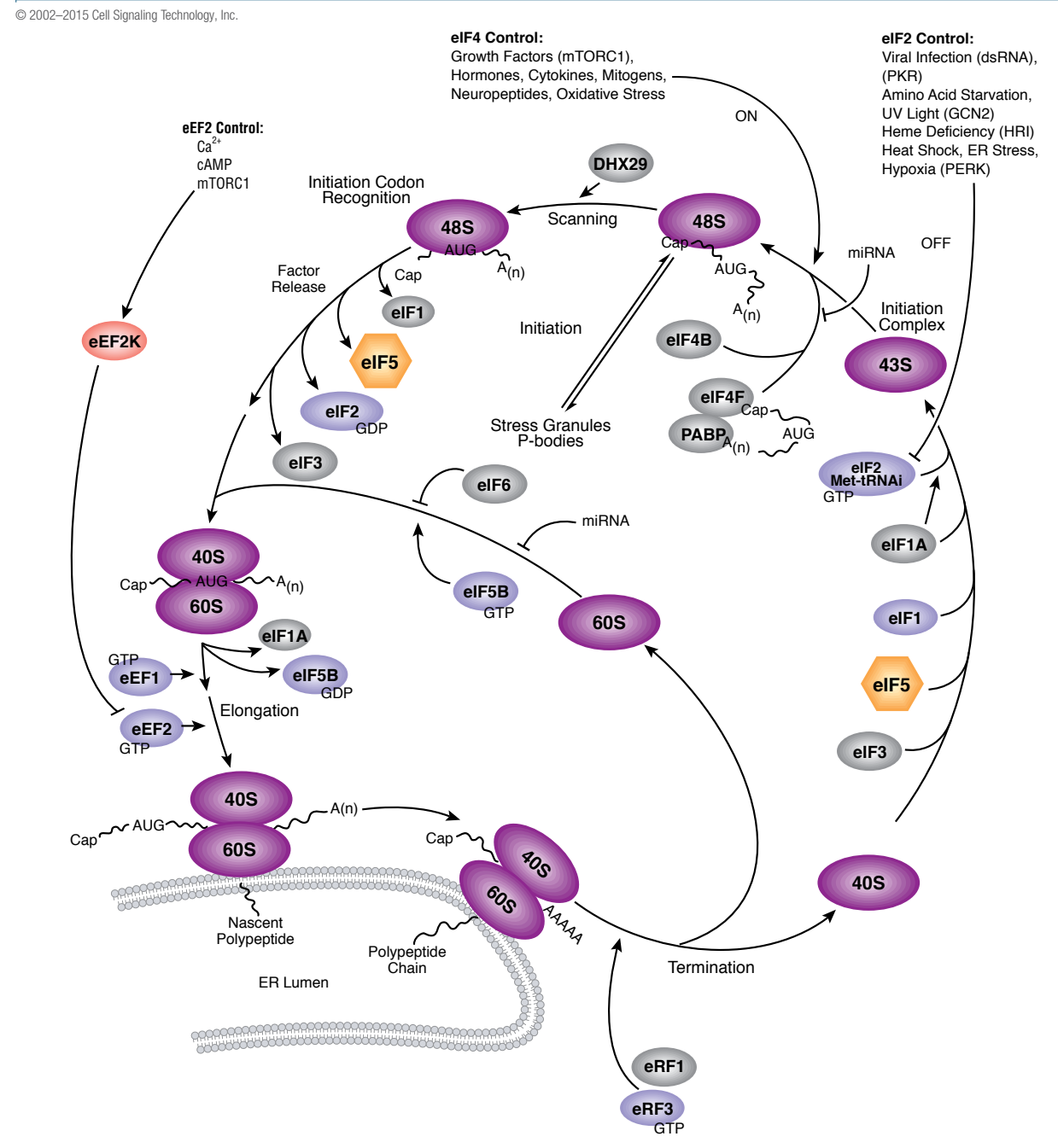


# Epigenetics & Translational Control

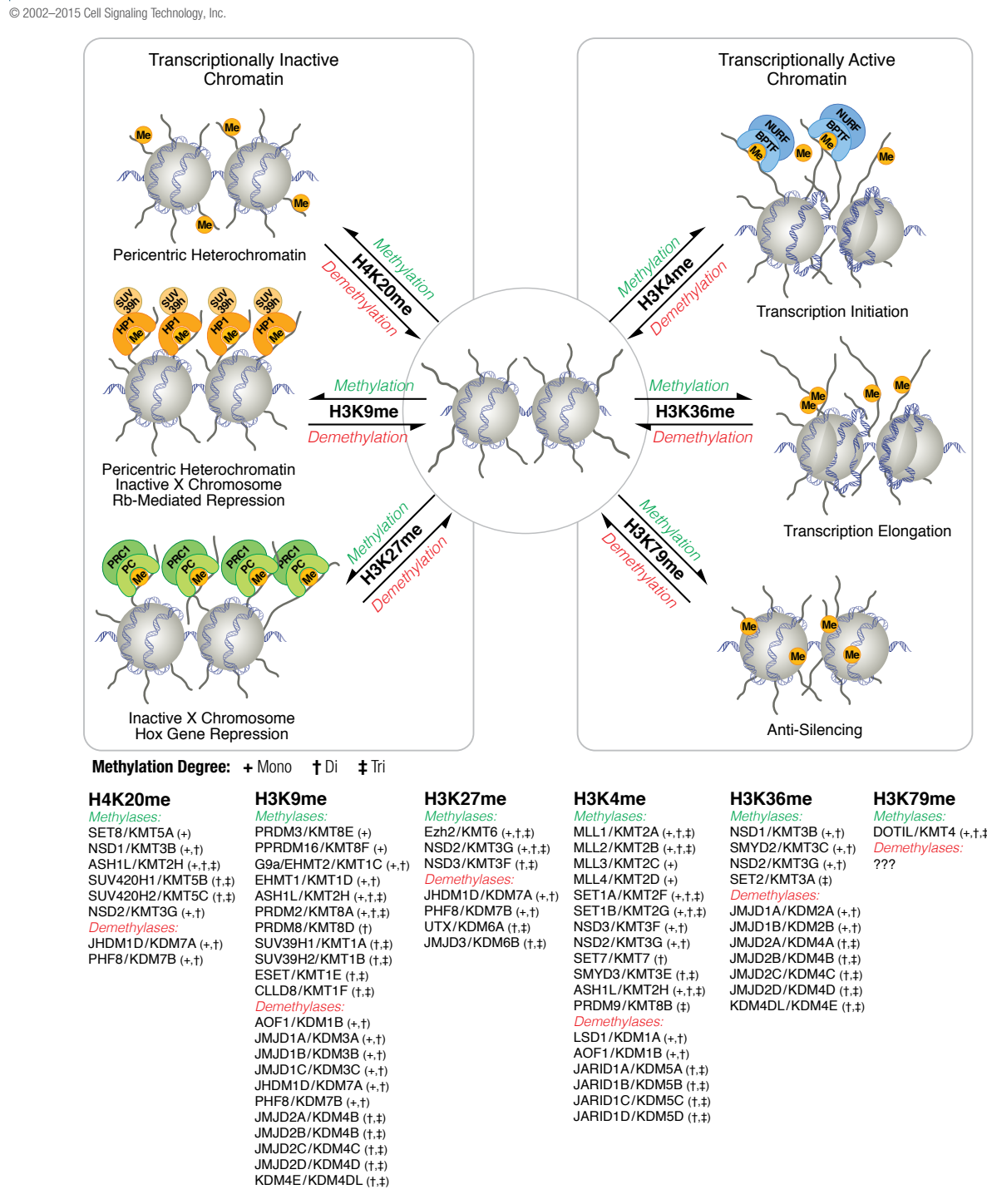
## Protein Acetylation



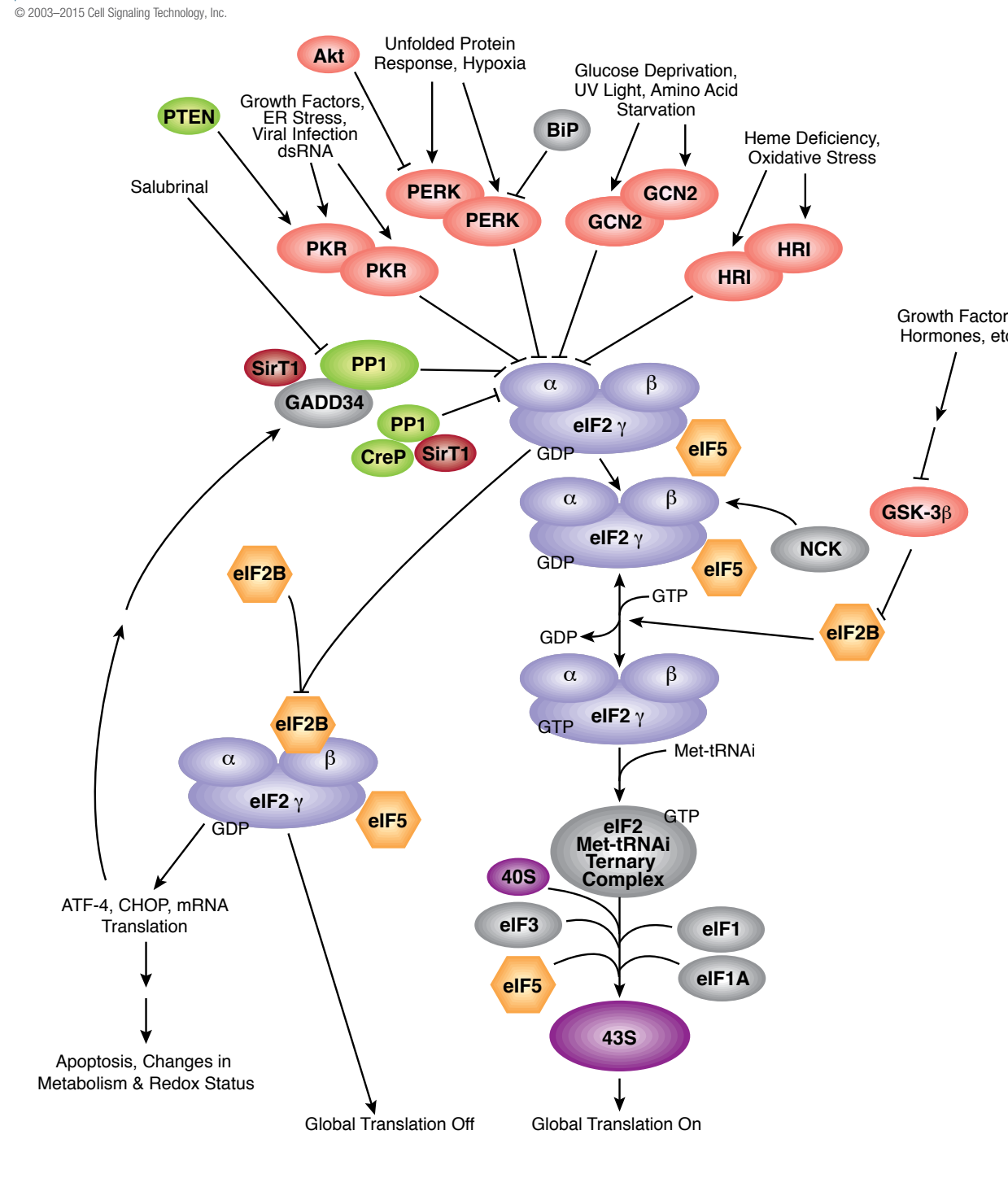
## Translational Control: Overview



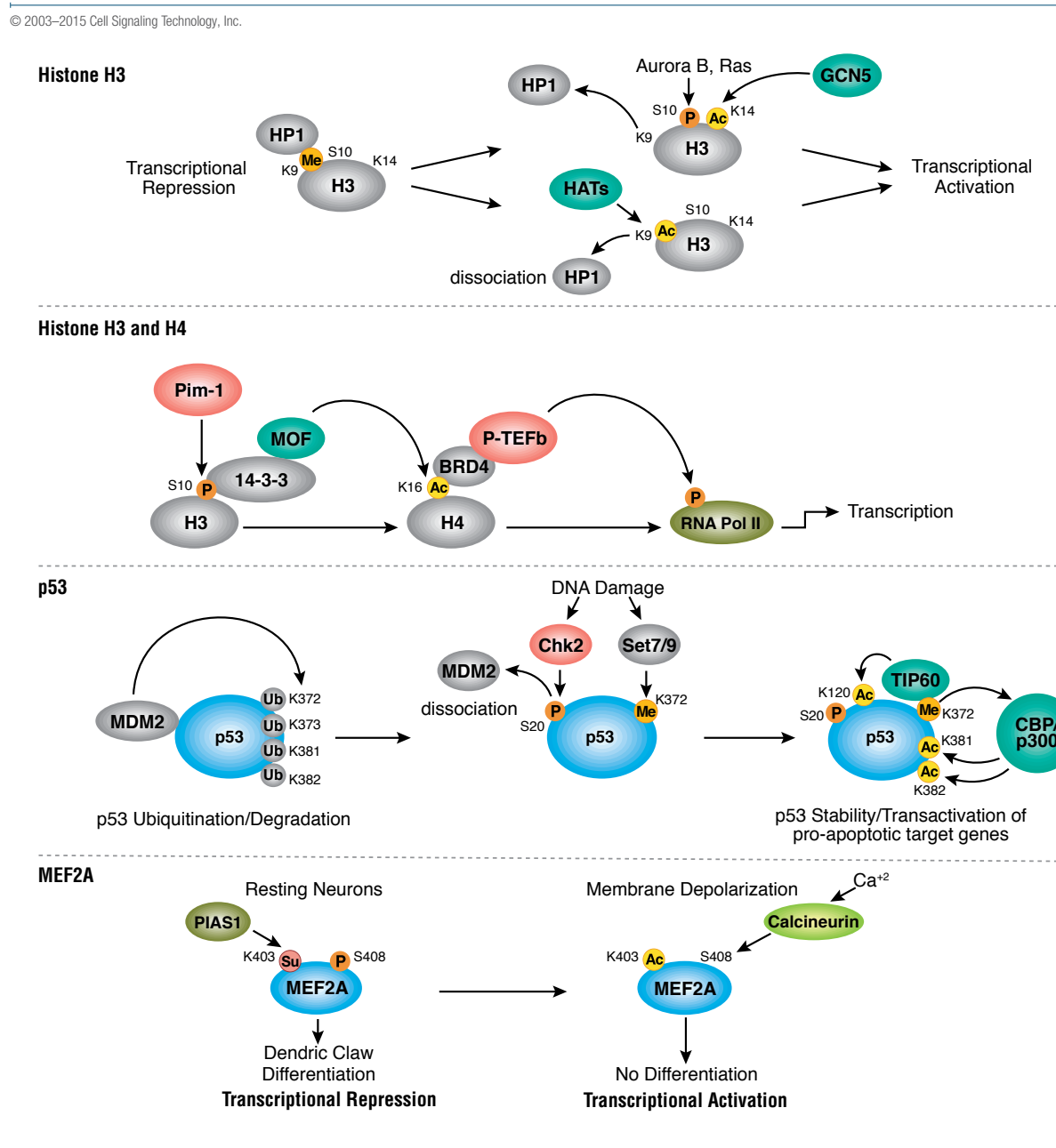
## Histone Lysine Methylation



## Translational Control: Regulation of eIF2



## Examples of Crosstalk Between Post-Translational Modifications



## Translational Control: Regulation of eIF4E and p70 S6K

