

p65 post-translational modifications

p65 is modified by acetylation, phosphorylation and methylation. Find out where these modifications occur, and whether they enhance or repress p65 transcriptional activity.

