

X-ChIP

Map proteins/histone modifications to genomic loci

