

GLuc-ON™ Promoter Reporter Clones

