

# GLuc-ON™ Promoter Reporter Clones

