

# GLuc-ON™ Promoter Reporter Vector

