

# OmicsLink™ Expression Clone (CMV Promoter)

