

OmicsLink™ Expression Clone (CMV Promoter)

