

OmicsLink™ Expression Clone (CAG Promoter)

