

OmicLink™ Expression Clone (CAG Promoter)

