

OmicLink™ Expression Clone (CMV Promoter)

