

OmicLink™ Expression Clone (PGK Promoter)

