

OmicLink™ Expression Clone (EF1α Promoter)

