

OmicLink™ Expression Clone (EF1a Promoter)

