

OmicLink™ Expression Clone (T7 Promoter)

