

OmicLink™ Expression Clone (UBC Promoter)

