

OmicsLinkTM Expression Clone (EF1 α Promoter)

