

OmicsLinkTM Expression Clone (CMV Promoter)

