

OmicsLink[™] Expression Clone (CMV Promoter)

