

OmicsLink[™] Expression Clone (CAG Promoter)

