

OmicsLink[™] Expression Clone (EF1a Promoter)

