

OmicsLink[™] Expression Clone (T7 Promoter)

