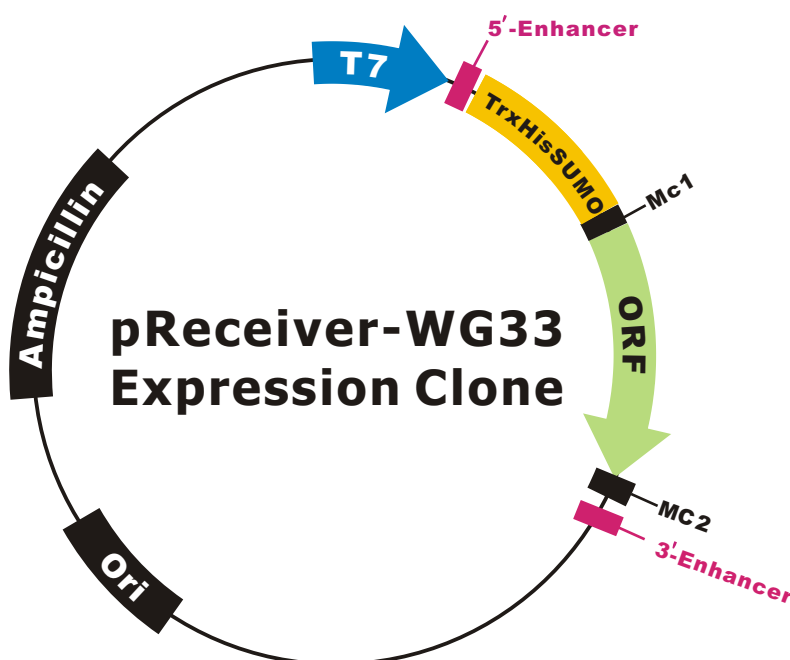


OmicsLink™ ORF Cell Free Expression Clone (T7 Promoter)



A detailed datasheet will be provided together with delivered OmicsLink™ ORF cell free expression clones, which contains restriction enzyme digestion information for the whole plasmid and linker sequence information at both ends of the ORF