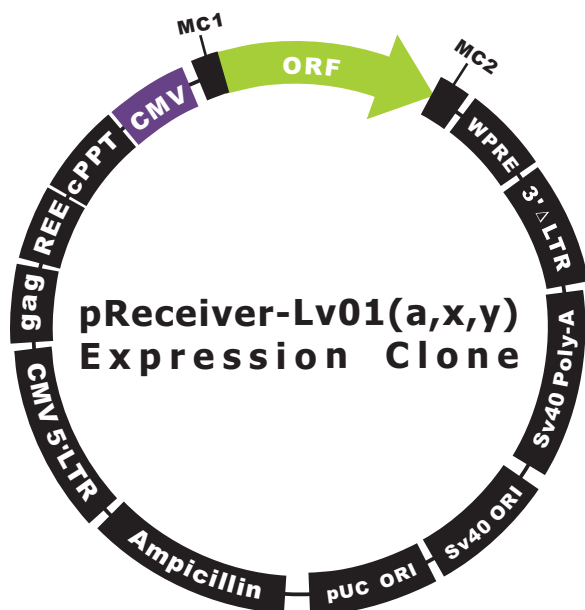


OmicLink™ Expression Clone (CMV Promoter)



Vector Features

| | |
|--------------------------------|------------|
| Promoter | CMV |
| Host Cell | Mammalian |
| Bacterial selection antibiotic | Ampicillin |
| Tag | n/a |