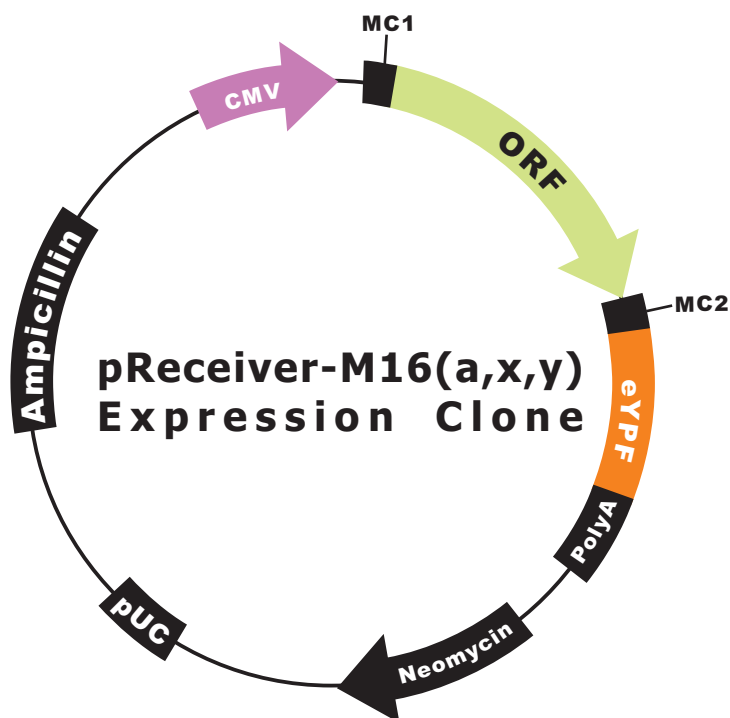


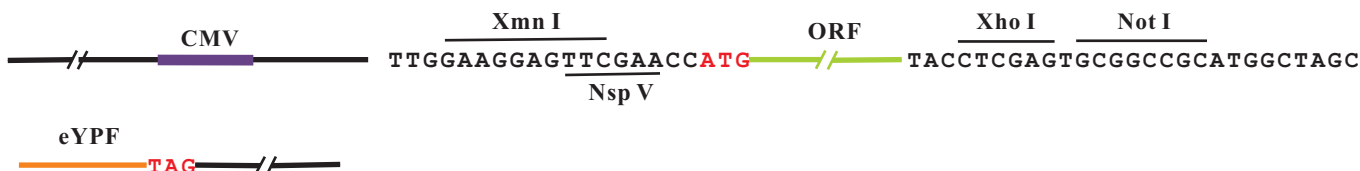
# OmicLink™ Expression Clone (CMV Promoter)



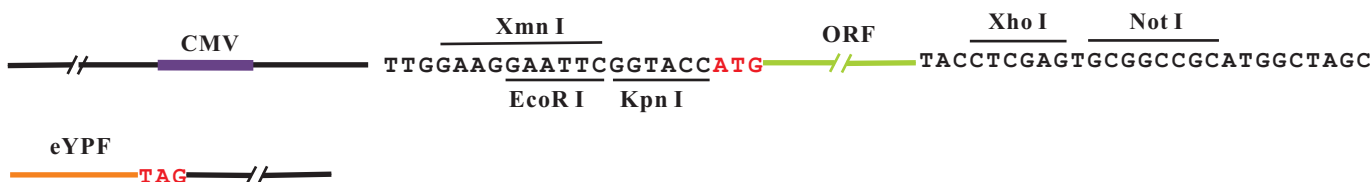
## Vector Features

|                                |            |
|--------------------------------|------------|
| Promoter                       | CMV        |
| Host Cell                      | Mammalian  |
| Bacterial selection antibiotic | Ampicillin |
| Mammalian selection marker     | Yes        |
| Tag                            | eYFP       |

## pReceiver-M16a



## pReceiver-M16x



## pReceiver-M16y

