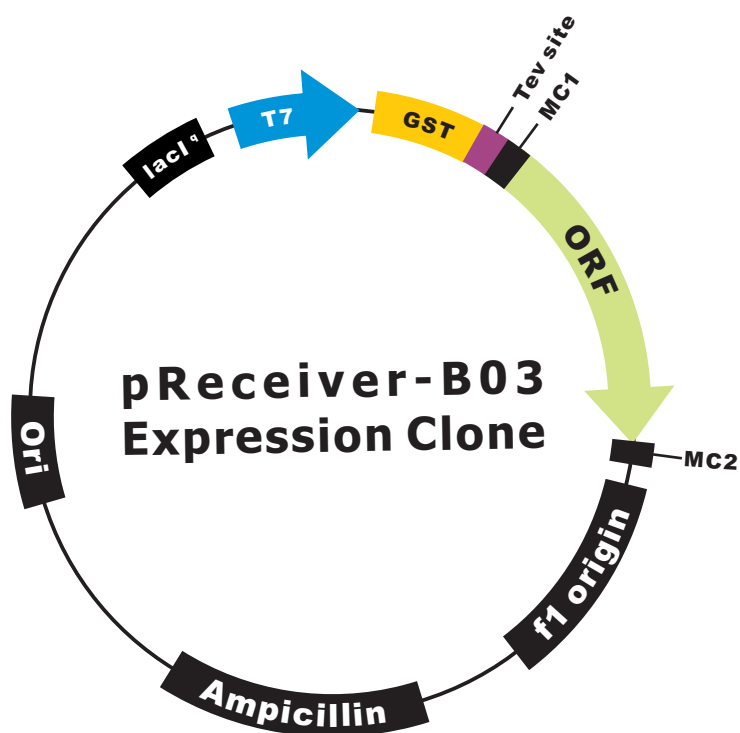


OmicLink™ Expression Clone (T7 Promoter)



Vector Features

| | |
|--------------------------------|----------------|
| Promoter | T7 |
| Host Cell | <i>E. Coli</i> |
| Bacterial selection antibiotic | Ampicillin |
| Tag | GST |