

pGEM[®]-T and pGEM[®]-T Easy Vector Systems

INSTRUCTIONS FOR USE OF PRODUCTS A1360, A1380, A3600 AND A3610.

Cloning PCR Products with pGEM[®]-T and pGEM[®]-T Easy Vectors

Ligation Using 2X Rapid Ligation Buffer

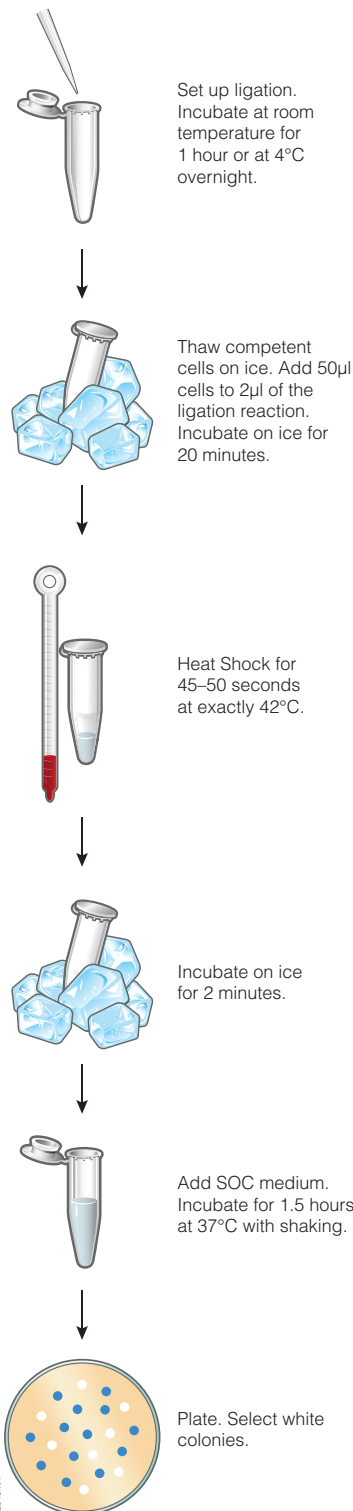
1. Briefly centrifuge the pGEM[®]-T or pGEM[®]-T Easy Vector and Control Insert DNA tubes to collect contents at the bottom of the tube.
2. Set up ligation reactions as described below. Vortex the 2X Rapid Ligation Buffer vigorously before each use. Use 0.5ml tubes known to have low DNA-binding capacity.

| Reagents | Standard Reaction | Positive Control | Background Control |
|---|-------------------|------------------|--------------------|
| 2X Rapid Ligation Buffer, T4 DNA Ligase | 5µl | 5µl | 5µl |
| pGEM [®] -T or pGEM [®] -T Easy Vector (50ng) | 1µl | 1µl | 1µl |
| PCR product | Xµl | – | – |
| Control Insert DNA | – | 2µl | – |
| T4 DNA Ligase (3 Weiss units/µl) | 1µl | 1µl | 1µl |
| Deionized water to a final volume of | 10µl | 10µl | 10µl |

3. Mix the reactions by pipetting. Incubate the reactions 1 hour at room temperature. Alternatively, incubate the reactions overnight at 4°C for the maximum number of transformants.

Transformation of JM109 High Efficiency Competent Cells

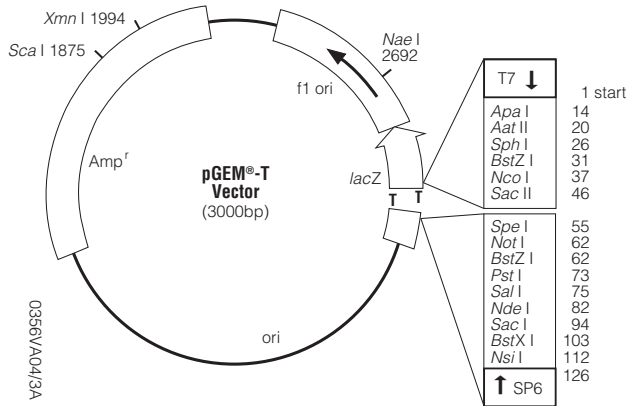
1. Prepare LB/ampicillin/IPTG/X-Gal plates.
2. Centrifuge the ligation reactions briefly. Add 2µl of each ligation reaction to a sterile 1.5ml tube on ice. Prepare a control tube with 0.1ng of uncut plasmid.
3. Place the JM109 High Efficiency Competent Cells in an ice bath until just thawed (5 minutes). Mix cells by gently flicking the tube.
4. Carefully transfer 50µl of cells to the ligation reaction tubes from Step 2. Use 100µl of cells for the uncut DNA control tube. Gently flick the tubes and incubate on ice for 20 minutes.
5. Heat-shock the cells for 45–50 seconds in a water bath at exactly 42°C. DO NOT SHAKE. Immediately return the tubes to ice for 2 minutes.
6. Add 950µl room temperature SOC medium to the ligation reaction transformations and 900µl to the uncut DNA control tube. Incubate for 1.5 hours at 37°C with shaking (~150rpm).
7. Plate 100µl of each transformation culture onto duplicate LB/ampicillin/IPTG/X-Gal plates. For the uncut DNA control, a 1:10 dilution with SOC is recommended.
8. Incubate plates overnight at 37°C. Select white colonies.



pGEM[®]-T and pGEM[®]-T Easy Vector Systems

INSTRUCTIONS FOR USE OF PRODUCTS A1360, A1380, A3600 AND A3610.

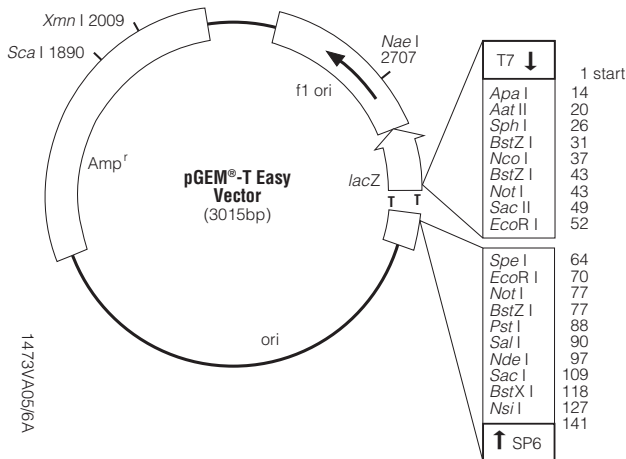
pGEM[®]-T Vector Circle Map and Sequence Reference Points



pGEM[®]-T Vector Sequence reference points:

| | |
|--|--------------------|
| T7 RNA Polymerase transcription initiation site | 1 |
| multiple cloning region | 10–113 |
| SP6 RNA Polymerase promoter (–17 to +3) | 124–143 |
| SP6 RNA Polymerase transcription initiation site | 126 |
| pUC/M13 Reverse Sequencing Primer binding site | 161–177 |
| <i>lacZ</i> start codon | 165 |
| <i>lac</i> operator | 185–201 |
| β-lactamase coding region | 1322–2182 |
| phage f1 region | 2365–2820 |
| <i>lac</i> operon sequences | 2821–2981, 151–380 |
| pUC/M13 Forward Sequencing Primer binding site | 2941–2957 |
| T7 RNA Polymerase promoter (–17 to +3) | 2984–3 |

pGEM[®]-T Easy Vector Circle Map and Sequence Reference Points



pGEM[®]-T Easy Vector Sequence reference points:

| | |
|--|--------------------|
| T7 RNA Polymerase transcription initiation site | 1 |
| multiple cloning region | 10–128 |
| SP6 RNA Polymerase promoter (–17 to +3) | 139–158 |
| SP6 RNA Polymerase transcription initiation site | 141 |
| pUC/M13 Reverse Sequencing Primer binding site | 176–197 |
| <i>lacZ</i> start codon | 180 |
| <i>lac</i> operator | 200–216 |
| β-lactamase coding region | 1337–2197 |
| phage f1 region | 2380–2835 |
| <i>lac</i> operon sequences | 2836–2996, 166–395 |
| pUC/M13 Forward Sequencing Primer binding site | 2949–2972 |
| T7 RNA Polymerase promoter (–17 to +3) | 2999-3 |

Ordering and Technical Information

www.promega.com • Phone 608-274-4330
Toll free in USA 800-356-9526 • Fax 608-277-2601