



Instruction Manual

Gateway[®] pENTR[™] Vectors

**Catalog nos. 11813-011, 11816-014, 11817-012,
11818-010, 11819-018**

Version B
23 November 2004
25-0521

A Limited Label License covers this product (see Purchaser Notification). By use of this product, you accept the terms and conditions of the Limited Label License.

Table of Contents

Table of Contents.....	iii
Important Information	v
Accessory Products.....	vi
Methods.....	1
Overview	1
Using the pENTR™ Vectors	3
Cloning PCR Products.....	6
Multiple Cloning Site for pENTR™1A	7
Multiple Cloning Site for pENTR™2B	8
Multiple Cloning Site for pENTR™3C.....	9
Multiple Cloning Site for pENTR™4	10
Multiple Cloning Site for pENTR™11	11
Transforming and Analyzing Entry Clones	12
Transforming and Analyzing Entry Clones	12
Guidelines to Perform the LR Recombination Reaction.....	14
Appendix.....	16
Blunt Cloning of PCR Products	16
Map and Features of the pENTR™ Vectors	18
Technical Service	20
Purchaser Notification.....	22
Gateway® Clone Distribution Policy.....	24
References.....	25

Important Information

pENTR™ Vectors

This manual is supplied with the following products.

Product	Catalog no.
pENTR™1A Vector	11813-011
pENTR™2B Vector	11816-014
pENTR™3C Vector	11817-012
pENTR™4 Vector	11818-010
pENTR™11 Vector	11819-018

Shipping and Storage

pENTR™ vectors are shipped at room temperature. Upon receipt, store at -20°C. Products are guaranteed for six months from date of shipment when stored properly.

Contents

10 µg pENTR™ vector, lyophilized in TE, pH 8.0.

Quality Control

pENTR™ vectors are qualified by restriction enzyme digestion, and in a recombination assay using Gateway® LR Clonase™ II enzyme mix. The *ccdB* gene is assayed by transformation using an appropriate *E. coli* strain.

Accessory Products

Additional Products

Additional products that may be used with the pENTR™ vectors are available from Invitrogen. Ordering information is provided below.

Item	Quantity	Catalog no.
LR Clonase™ II Enzyme Mix	20 reactions	11791-020
	100 reactions	11791-100
Library Efficiency® DB3.1™ Competent Cells	5 x 0.2 ml	11782-018
One Shot® TOP10 Chemically Competent Cells	10 reactions	C4040-10
	20 reactions	C4040-03
One Shot® TOP10 Electrocompetent Cells	10 reactions	C4040-50
	20 reactions	C4040-52
One Shot® MAX Efficiency® DH5α™ -T1® Chemically Competent Cells	20 reactions	12297-016
Kanamycin Sulfate	1 g	11815-016

Gateway® Destination Vectors

A large selection of Gateway® destination vectors is available from Invitrogen to facilitate expression of your gene of interest in virtually any protein expression system. For more information about the vectors available, refer to the Gateway® Technology Central application portal on our Web site at www.invitrogen.com/gateway or contact Technical Service (see page 20).

Methods

Overview

Introduction

The pENTR™ vectors allow restriction cloning of a gene of interest into a vector for entry into the Gateway® System available from Invitrogen. A choice of pENTR™ vectors is available (see table below) for optimal expression of your gene of interest after recombination with the Gateway® destination vector of choice. For more information about the Gateway® Technology, see the next page.

Product	Benefit
pENTR™1A Vector	<ul style="list-style-type: none">• Three reading frames available• Kozak sequence for efficient initiation of translation in eukaryotic cells• <i>E. coli</i> ribosome binding site for efficient initiation of translation in prokaryotic cells (pENTR™1A and pENTR™3C only)
pENTR™2B Vector	
pENTR™3C Vector	
pENTR™4 Vector	<ul style="list-style-type: none">• Same multiple cloning site as pENTR™1A except that first restriction enzyme site is <i>Nco</i> I• Kozak sequence for efficient initiation of translation in eukaryotic cells
pENTR™11 Vector	<ul style="list-style-type: none">• Kozak sequence for efficient initiation of translation in eukaryotic cells• Two <i>E. coli</i> ribosome binding sites for efficient initiation of translation in prokaryotic cells

continued on next page

Overview, continued

Features of the pENTR™ Vectors

The pENTR™ vectors contain the following elements:

- *rrnB* transcription termination sequences to prevent basal expression of the PCR product of interest in *E. coli*
 - *attL1* and *attL2* sites for site-specific recombination of the entry clone with a Gateway® destination vector (for more information, refer to the Gateway® Technology with Clonase™ II manual or Landy, 1989)
 - Kozak consensus sequence for efficient translation initiation in eukaryotic systems
 - Ribosome binding site for efficient translation initiation in prokaryotic systems (**pENTR™1A, pENTR™3C, and pENTR™11 only**)
 - The *ccdB* gene located between the two *attL* sites for negative selection
 - Kanamycin resistance gene for selection in *E. coli*
 - pUC origin for high-copy replication and maintenance of the plasmid in *E. coli*
-

The Gateway® Technology

The Gateway® Technology is a universal cloning method that takes advantage of the site-specific recombination properties of bacteriophage lambda (Landy, 1989) to provide a rapid and highly efficient way to move your gene of interest into multiple vector systems. To express your gene of interest using the Gateway® Technology, simply:

1. Clone your gene of interest into one of the pENTR™ vectors to generate an entry clone.
2. Generate an expression clone by performing a recombination reaction between the entry clone and a Gateway® destination vector of choice.
3. Introduce your expression clone into the appropriate host (*e.g.* bacterial, mammalian, yeast, insect) and express your recombinant protein.

For more information about the Gateway® Technology, refer to the Gateway® Technology with Clonase™ II manual. You may download the manual from www.invitrogen.com or contact Technical Service (page 20).

Using the pENTR™ Vectors

Introduction

This section provides general guidelines for using the pENTR™ vectors. Diagrams are provided on pages 7-11 to help you ligate your gene of interest into the appropriate pENTR™ vector.

Propagating the pENTR™ Vectors

If you wish to propagate and maintain the pENTR™ vectors, we recommend using Library Efficiency® DB3.1™ Competent Cells (Catalog no. 11782-018) from Invitrogen for transformation. The DB3.1™ *E. coli* strain is resistant to CcdB effects and can support the propagation of plasmids containing the *ccdB* gene.

Note: Do not use general *E. coli* cloning strains including TOP10 or DH5α for propagation and maintenance as these strains are sensitive to CcdB effects.

Resuspension

Before using, resuspend your pENTR™ plasmid DNA in 100 µl of sterile water to a final concentration of 100 ng/µl.

General Molecular Biology Techniques

For help with DNA ligations, *E. coli* transformations, restriction enzyme analysis, DNA sequencing, and DNA biochemistry, refer to *Molecular Cloning: A Laboratory Manual* (Sambrook et al., 1989) or *Current Protocols in Molecular Biology* (Ausubel et al., 1994).



Important

Your gene of interest must replace the *ccdB* gene located between the two *attL* sites. Before cloning your gene of interest into a pENTR™ vector, we recommend that you:

- Digest the pENTR™ vector on each side of the *ccdB* gene
- Dephosphorylate and gel purify the pENTR™ vector

This will minimize the competition between the *ccdB* fragment and your gene of interest during the ligation process.

For more guidelines to help you develop your cloning strategy, see **Cloning Considerations** on page 5.

continued on next page

Using the pENTR™ Vectors, continued

Kozak Sequence for Mammalian Expression

If you will be recombining your entry clone with a destination vector for mammalian expression, your insert should contain a Kozak consensus sequence with an ATG initiation codon for proper initiation of translation (Kozak, 1987; Kozak, 1991; Kozak, 1990). An example of a Kozak consensus sequence is provided below. Other sequences are possible, but the G or A at position -3 and the G at position +4 are the most critical for function (shown in bold). The ATG initiation codon is shown underlined.

(G/A)NNATGG

Note: Cloning a blunt-ended fragment containing a 5' ATGG (where ATG is the initiation codon) into the *Xmn* I site of any of the pENTR™ vectors will constitute a Kozak consensus sequence (see diagrams on pages 7-11).

Ribosome Binding Site for Prokaryotic Expression

If you will be recombining your entry clone with a destination vector for prokaryotic expression, your insert should contain an *E. coli* ribosome binding site [AAGGA(A/G)] approximately 9-10 base pairs upstream of the ATG initiation codon (Gold, 1988; Miller, 1992). This will ensure the optimal spacing for proper translation.

Note: Ribosome binding sites are provided in pENTR™1A, pENTR™3C, and pENTR™11 (see diagrams on pages 7-11). If your insert will not be properly spaced from the vector-encoded ribosome binding site, you will need to include your own ribosome binding site for proper initiation of translation.

continued on next page

Using the pENTR™ Vectors, continued

Cloning

Considerations

Consider the following factors when cloning into the pENTR™ vectors.

If you wish to....	Then your insert...
express your native protein without an N-terminal or C-terminal tag	<ul style="list-style-type: none">• should contain a Kozak consensus sequence for mammalian expression or an <i>E. coli</i> ribosome binding site for prokaryotic expression (see previous page for more information)• should contain a stop codon if one is not provided in the destination vector
include an N-terminal tag (following recombination of the entry clone with a Gateway® destination vector)	<ul style="list-style-type: none">• does not need a Kozak consensus sequence, <i>E. coli</i> ribosome binding site, or an ATG initiation codon (these will be provided by the appropriate destination vector)• should be in frame with the tag after recombination (see diagrams on pages 7-11)• should contain a stop codon if one is not provided in the destination vector
include a C-terminal tag (following recombination of the entry clone with a Gateway® destination vector)	<ul style="list-style-type: none">• should contain a Kozak consensus sequence for mammalian expression or an <i>E. coli</i> ribosome binding site for prokaryotic expression (see previous page for more information)• should be in frame with the tag after recombination (see diagrams on pages 7-11)• should not contain a stop codon
include an N-terminal and C-terminal tag (following recombination of the entry clone with a Gateway® destination vector)	<ul style="list-style-type: none">• does not need a Kozak consensus sequence, <i>E. coli</i> ribosome binding site, or an ATG initiation codon (these will be provided by the appropriate destination vector)• should be in frame with both the N-terminal and C-terminal tags after recombination (see diagrams on pages 7-11)• should not contain a stop codon

Cloning PCR Products



Note

If you include an N-terminal tag following recombination with a destination vector, and your insert contains an ATG initiation codon, note that translation initiation may also occur at this site. This may result in a small amount of native, untagged protein being expressed along with your tagged fusion protein.



If you wish to clone a PCR product made using primers containing restriction enzyme sites, we recommend the following to ensure efficient cloning:

- Inactivate or remove the DNA polymerase (*Taq* DNA polymerase can fill in sticky ends and add bases to blunt ends of PCR products) using phenol extraction or the S.N.A.P.[™] MiniPrep Kit available from Invitrogen (Catalog no. K1900-01).
- Remove small DNA fragments such as primers, primer-dimers, and excess dNTP's. Refer to the Gateway[®] Technology with Clonase[™] II manual for a purification protocol using PEG/MgCl₂ precipitation.

Cloning Blunt PCR Products

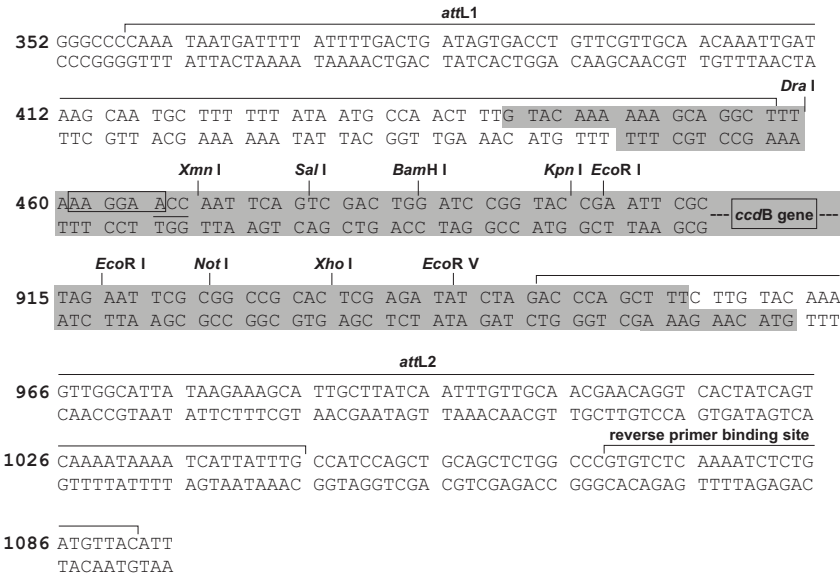
Because primers usually contain a 5' hydroxy group, PCR products generally do not have 5' phosphates and are not necessarily blunt. If you wish to clone a blunt PCR product into your entry vector, we recommend you perform the **Blunt Cloning of PCR Products** protocol provided in the **Appendix**, page 16.

Multiple Cloning Site for pENTR™ 1A

Multiple Cloning Site

Below is the multiple cloning site for pENTR™ 1A. Restriction sites are labeled to indicate the cleavage site. **Note:** Your gene of interest must replace the *ccdB* gene located between the two *attL* sites. Features are indicated as follows:

- The *attL* sites are properly spaced to indicate the correct reading frame for fusion of your gene of interest to an N-terminal or C-terminal tag following recombination with a destination vector.
- Shaded regions correspond to those DNA sequences transferred from the entry clone into the destination vector following recombination.
- Boxed and underlined sequences correspond to the *E. coli* ribosome binding site [AAGGA(A/G)] and the 5' end of the Kozak consensus sequence (ACC), respectively.



Multiple Cloning Site for pENTR™ 2B

Multiple Cloning Site

Below is the multiple cloning site for pENTR™2B. Restriction sites are labeled to indicate the cleavage site. **Note:** Your gene of interest must replace the *ccdB* gene located between the two *attL* sites. Features are indicated as follows:

- The *attL* sites are properly spaced to indicate the correct reading frame for fusion of your gene of interest to an N-terminal or C-terminal tag following recombination with a destination vector.
- Shaded regions correspond to those DNA sequences transferred from the entry clone into the destination vector following recombination.
- Underlined sequence corresponds to the 5' end of the Kozak consensus sequence (ACC).

attL1

352GGGCCCCAAA TAATGATTTT ATTTTGACTG ATAGTGACCT GTTCGTTGCA ACAAATTGAT
CCCGGGGTTT ATTACTAAAA TAAAACTGAC TATCACTGGA CAAGCAACGT TGTTTAACTA

412AAG CAA TGC TTT TTT ATA ATG CCA ACT TTG TAC AAA AAA GCA GGC TGG
TTC GTT ACG AAA AAA TAT TAC GGT TGA AAC ATG TTT TTT CGT CCG ACC

Ehe I Xmn I Sal I BamH I Kpn I EcoR I

460CGC CGG AAC CAA TTC AGT CGA CTG GAT CCG GTA CCG AAT TCG
GCG GCC TTG GTT AAG TCA GCT GAC CTA GGC CAT GGC TTA AGC

ccdB gene

EcoR I Not I Xho I EcoR V

916TAG AAT TCG CGG CCG CAC TCG AGA TAT CTA GAC CCA GCT TTC TTG TAC AAA
ATC TTA AGC GCC GGC GTG AGC TCT ATA GAT CTG GGT CGA AAG AAC ATG TTT

attL2

967GTTGGCATT TAAGAAAGCA TTGCTTATCA ATTTGTTGCA ACGAACAGGT CACTATCAGT
CAACCGTAAT ATTCTTTTCGT AACGAATAGT TAAACAACGT TGCTTGTTCCA GTGATAGTCA

reverse primer binding site

1027CAAAATAAAA TCATTATTTG CCATCCAGCT GCAGCTCTGG CCCGTGTCTC AAAATCTCTG
GTTTTATTTT AGTAATAAAC GGTAGGTCGA CGTCGAGACC GGGCACAGAG TTTTAGAGAC

1087ATGTTACATT
TACAATGTAA

Multiple Cloning Site for pENTR™ 3C

Multiple Cloning Site

Below is the multiple cloning site for pENTR™3C. Restriction sites are labeled to indicate the cleavage site. **Note:** Your gene of interest must replace the *ccdB* gene located between the two *attL* sites. Features are indicated as follows:

- The *attL* sites are properly spaced to indicate the correct reading frame for fusion of your gene of interest to an N-terminal or C-terminal tag following recombination with a destination vector.
- Shaded regions correspond to those DNA sequences transferred from the entry clone into the destination vector following recombination.
- Boxed and underlined sequences correspond to the *E. coli* ribosome binding site [AAGGA(A/G)] and the 5' end of the Kozak consensus sequence (ACC), respectively.

```

                                     attL1
352  GGGCCCCAAA TAATGATTTT ATTTTGACTG ATAGTGACCT GTTCGTTGCA ACAAATTGAT
    CCCGGGGTTT ATTACTAAAA TAAAACTGAC TATCACTGGA CAAGCAACGT TGTTTAACTA

412  AAG CAA TGC TTT TTT ATA ATG CCA ACT TTG TAC AAA AAA GCA GGC TCT
    TTC GTT ACG AAA AAA TAT TAC GGT TGA AAC ATG TTT TTT CGT CCG AGA

    Dra I      Xmn I      Sal I      BamH I      Kpn I EcoR I
460  TTA AAG GAA CCA ATT CAG TCG ACT GGA TCC GGT ACC GAA TTC --- ccdB gene ---
    AAT TTC CTT GGT TAA GTC AGC TGA CCT AGG CCA TGG CTT AAG

    EcoR I      Not I      Xho I      EcoR V
921  TAG AAT TCG CGG CCG CAC TCG AGA TAT CTA GAC CCA GCT TTC TTG TAC AAA
    ATC TTA AGC GCC GGC GTG AGC TCT ATA GAT CTG GGT CGA AAG AAC ATG TTT

                                     attL2
972  GTTGGCATT TAAGAAAGCA TTGCTTATCA ATTTGTTGCA ACGAACAGGT CACTATCAGT
    CAACCGTAAT ATTCTTTTCGT AACGAATAGT TAAACAACGT TGCTTGTTCCA GTGATAGTCA

                                     reverse primer binding site
1032 CAAAATAAAA TCATTATTTG CCATCCAGCT GCAGCTCTGG CCCGTGTCTC AAAATCTCTG
    GTTTTATTTT AGTAATAAAC GGTAGGTCGA CGTCGAGACC GGGCACAGAG TTTTAGAGAC

1092 ATGTTACATT
    TACAATGTAA
```

Multiple Cloning Site for pENTR™ 4

Multiple Cloning Site

Below is the multiple cloning site for pENTR™4. Restriction sites are labeled to indicate the cleavage site. **Note:** Your gene of interest must replace the *ccdB* gene located between the two *attL* sites. Features are indicated as follows:

- The *attL* sites are properly spaced to indicate the correct reading frame for fusion of your gene of interest to an N-terminal or C-terminal tag following recombination with a destination vector.
- Shaded regions correspond to those DNA sequences transferred from the entry clone into the destination vector following recombination.
- Underlined sequence corresponds to the 5' end of the Kozak consensus sequence (ACC).

attL1

```
352  GGGCCCCAAA TAATGATTTT ATTTTGACTG ATAGTGACCT GTTCGTTGCA ACAAATTGAT
    CCCGGGGTTC ATTACTAAAA TAAAACTGAC TATCACTGGA CAAGCAACGT TGTTTAACTA

412  AAG CAA TGC TTT TTT ATA ATG CCA ACT TTG TAC AAA AAA GCA GGC TCC
    TTC GTT ACG AAA AAA TAT TAC GGT TGA AAC ATG TTT TTT CGT CCG AGG

    Nco I      Xmn I      Sal I      BamH I      Kpn I      EcoR I
460  ACC ATG GGA ACC AAT TCA GTC GAC TGG ATC CGG TAC CGA ATT CGC
    TGG TAC CCT TGG TTA AGT CAG CTG ACC TAG GCC ATG GCT TAA GCG --- ccdB gene ---

    EcoR I      Not I      Xho I      EcoR V
918  TAG AAT TCG CGG CCG CAC TCG AGA TAT CTA GAC CCA GCT TTC TTG TAC AAA
    ATC TTA AGC GCC GGC GTG AGC TCT ATA GAT CTG GGT CGA AAG AAC ATG TTT

                                attL2

969  GTTGGCATT TAAGAAAGCA TTGCTTATCA ATTTGTTGCA ACGAACAGGT CACTATCAGT
    CAACCGTAAT ATTCTTTCGT AACGAATAGT TAAACAACGT TGCTTGTCCTA GTGATAGTCA

                                reverse primer binding site
1029  CAAAATAAAA TCATTATTTG CCATCCAGCT GCAGCTCTGG CCCGTGTCTC AAAATCTCTG
    GTTTTATTTT AGTAATAAAC GGTAGGTCGA CGTCGAGACC GGGCACAGAG TTTTAGAGAC

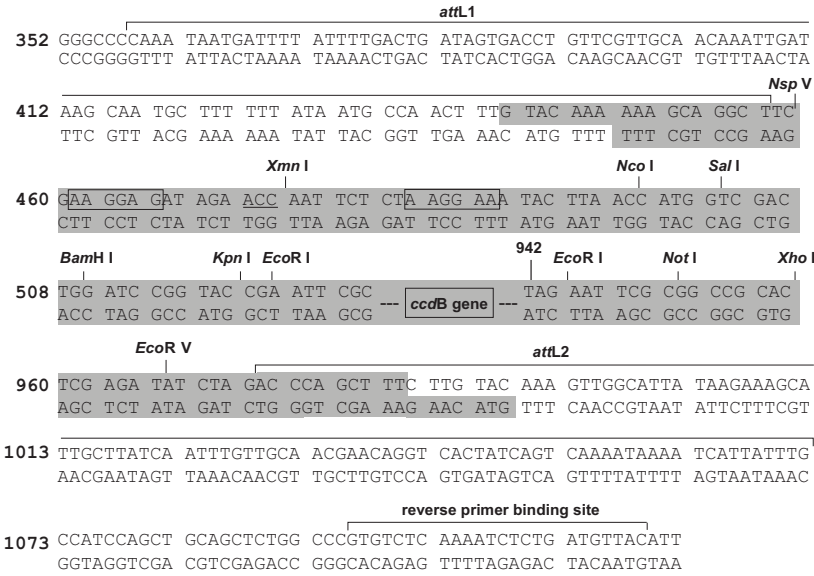
1089  ATGTTACATT
    TACAATGTAA
```


Multiple Cloning Site for pENTR™ 11

Multiple Cloning Site

Below is the multiple cloning site for pENTR™11. Restriction sites are labeled to indicate the cleavage site. **Note:** Your gene of interest must replace the *ccdB* gene located between the two *attL* sites. Features are indicated as follows:

- The *attL* sites are properly spaced to indicate the correct reading frame for fusion of your gene of interest to an N-terminal or C-terminal tag following recombination with a destination vector.
- Shaded regions correspond to those DNA sequences transferred from the entry clone into the destination vector following recombination.
- Boxed and underlined sequences correspond to the two available *E. coli* ribosome binding site [AAGGA(A/G)] and the 5' end of the Kozak consensus sequence (ACC), respectively.



Transforming and Analyzing Entry Clones

Introduction

Once you have restriction cloned your gene of interest into your entry vector, you will transform the ligation reaction into competent *E. coli* and select for positive transformants. See below for general guidelines to transform and analyze your entry clones.

General Molecular Biology Techniques

For help with DNA ligations, *E. coli* transformations, restriction enzyme analysis, DNA sequencing, and DNA biochemistry, refer to *Molecular Cloning: A Laboratory Manual* (Sambrook et al., 1989) or *Current Protocols in Molecular Biology* (Ausubel et al., 1994).

E. coli Transformation

Transform your ligation mixture into a competent *recA*, *endA* *E. coli* strain (e.g. TOP10, DH5 α) and select on LB plates containing 50 μ g/ml kanamycin. For your convenience, competent TOP10 and DH5 α *E. coli* are available from Invitrogen in a One Shot[®] format (see page vi for ordering information).

Analyzing Positive Clones

1. Pick 5 colonies and culture them overnight in LB or SOB medium containing 50 μ g/ml kanamycin.
 2. Isolate plasmid DNA using your method of choice. We recommend using the S.N.A.P.[™] MidiPrep Kit (Catalog no. K1910-01) or the PureLink[™] HQ Mini Plasmid Purification Kit (Catalog no. K2100-01).
 3. Analyze the entry clones by restriction analysis to confirm the presence and correct orientation of the insert. Use a restriction enzyme or a combination of enzymes that cut once in the vector and once in the insert.
-

continued on next page

Transforming and Analyzing Entry Clones, continued

Analyzing Transformants by PCR

You may also analyze positive transformants using PCR. Use a primer that hybridizes within the pENTR™ vector and one that hybridizes within your insert. You will have to determine the amplification conditions. If you are using this technique for the first time, perform restriction analysis in parallel. Artifacts may be obtained because of mispriming or contaminating template. The protocol below is provided for your convenience. Other protocols are suitable.

Materials Needed:

PCR SuperMix High Fidelity (Catalog no. 10790-020)

Appropriate forward and reverse PCR primers, 20 μM each

Protocol:

1. For each sample, aliquot 48 μl of PCR SuperMix High Fidelity into a 0.5 ml microcentrifuge tube. Add 1 μl each of the forward and reverse PCR primer.
2. Pick 5 colonies and resuspend them individually in 50 μl of the PCR SuperMix containing primers (make a patch plate to preserve the colonies for further analysis).
3. Incubate reaction for 10 minutes at 94°C to lyse cells and inactivate nucleases.
4. Amplify for 20 to 30 cycles.
5. For the final extension, incubate at 72°C for 10 minutes. Store at +4°C.
6. Visualize by agarose gel electrophoresis.

Sequencing

You may sequence your entry clone using the recommended primer (see table below and the diagrams on pages 7-11) to confirm the presence and orientation of the insert. For your convenience, Invitrogen offers a custom primer synthesis service. See www.invitrogen.com or contact Technical Service (page 20) for more information.

Primer	Sequence
Reverse	5'-GTAACATCAGAGATTTTGAGACAC-3'

Guidelines to Perform the LR Recombination Reaction

Introduction

Once you have obtained an entry clone containing your gene of interest, you may perform an LR recombination reaction between the entry clone and a destination vector of choice. General guidelines are provided below.



Important

For most applications, we recommend performing the LR recombination reaction using a:

- Supercoiled *attL*-containing entry clone
- Supercoiled *attR*-containing destination vector

Note: If your destination vector or entry clone is large (>10 kb), you may linearize either vector to increase recombinational efficiency. You may also relax the destination vector using topoisomerase I to increase efficiency. For more details, refer to the Gateway® Technology with Clonase™ II manual.

Destination Vectors

A large selection of Gateway® destination vectors is available from Invitrogen to facilitate expression of your gene of interest in virtually any protein expression system. For more information about the vectors available, refer to our Web site (www.invitrogen.com) or contact Technical Service (page 20).

E. coli Host

Once you have performed the LR recombination reaction, you will transform the reaction mixture into competent *E. coli* and select for expression clones. You may use any *recA*, *endA* *E. coli* strain including TOP10, DH5α™, DH10B™ or equivalent for transformation. **DO NOT** transform the LR reaction mixture into *E. coli* strains that contain the F' episome (e.g. TOP10F'). These strains contain the *ccdA* gene and will prevent negative selection with the *ccdB* gene.

continued on next page

Guidelines to Perform the LR Recombination Reaction, continued

Performing the LR Reaction

To perform the Gateway® LR recombination reaction, you will need:

- Purified plasmid DNA of your entry clone
- A destination vector of choice
- LR Clonase™ II enzyme mix (Catalog no. 11791-020; see below)
- 2 µg/µl proteinase K solution (supplied with the LR Clonase™ II enzyme mix)
- TE Buffer, pH 8.0 (10 mM Tris-HCl, pH 8.0, 1 mM EDTA)
- Appropriate chemically competent *E. coli* host and growth media for expression
- Appropriate selective plates

For instructions to perform the LR recombination reaction, refer to the LR Clonase™ II enzyme mix manual or to the manual for the destination vector you are using.

LR Clonase™ II Enzyme Mix

To catalyze the LR recombination reaction, we recommend using Gateway® LR Clonase™ II enzyme mix. The LR Clonase™ II enzyme mix combines the proprietary enzyme formulation and 5X LR Reaction Buffer previously supplied by Invitrogen as separate components in LR Clonase™ enzyme mix (Catalog no. 11791-019) into an optimized single-tube format for easier set-up of the LR recombination reaction.

Note: You may perform the LR recombination reaction using LR Clonase™ enzyme mix, if desired. To use LR Clonase™ enzyme mix, follow the instructions included with the product. **Do not** use the protocol for LR Clonase™ II enzyme mix as reaction conditions differ.

Appendix

Blunt Cloning of PCR Products

Introduction

Use this protocol to clone blunt-end PCR products into your pENTR™ vector.

Materials Needed

You should have the following materials on hand before beginning:

- PCR product (~40 ng as judged from an agarose gel)
 - 3 M sodium acetate
 - 100% ethanol
 - 10 mM ATP
 - 2 mM dNTP's
 - 5X T4 forward reaction buffer (350 mM Tris-HCl, pH 7.6; 50 mM MgCl₂; 500 mM KCl; 5 mM 2-mercaptoethanol)
 - T4 polynucleotide kinase and buffer (10 units/μl) (Catalog no. 18004-010)
 - T4 DNA polymerase (5 units/μl) (Catalog no. 18005-017)
 - 30% PEG 8000/30 mM MgCl₂
 - T4 DNA ligase and buffer (1 unit/μl) (Catalog no. 15224-017)
 - Entry vector (blunt, dephosphorylated, ~50 ng)
-

continued on next page

Blunt Cloning of PCR Products, continued

Protocol

1. In a 0.5 ml tube, precipitate approximately 40 ng of PCR product by adding 0.1 volume of 3 M sodium acetate followed by 2.5 volumes of 100% ethanol.
2. Pellet the DNA by centrifugation. Wash the pellet twice with 70% ethanol.
3. Add the following reagents to the DNA:

Distilled H ₂ O	4 µl
10 mM ATP	1 µl
2 mM of each dNTP (i.e. dATP, dCTP, dTTP, dGTP)	1 µl
5X T4 Forward Reaction Buffer	2 µl
T4 polynucleotide kinase	1 µl
<u>T4 DNA polymerase</u>	<u>1 µl</u>
Total Volume	10 µl
4. Incubate at 37°C for 10 minutes, then at 65°C for 15 minutes. Cool on ice for 5 minutes. Centrifuge briefly to bring any condensate to the bottom of the tube.
5. Add 5 µl of 30% PEG 8000/30 mM MgCl₂. Mix and centrifuge immediately at room temperature for 10 minutes.
6. Carefully remove and discard supernatant.
7. Dissolve the invisible pellet in a 10 µl cocktail containing:

2 µl of 5X T4 DNA ligase buffer
0.5 units T4 DNA ligase
~ 50 ng of blunt, dephosphorylated entry vector
Sterile water up to 10 µl
8. Incubate at 25°C for 1 hour, then at 65°C for 10 minutes. Add 40 µl TE.
9. Transform competent *E. coli* using your method of choice.

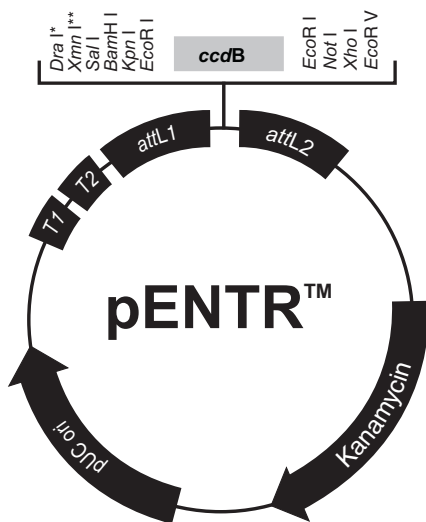
Transformation

Refer to the Gateway® Technology with Clonase™ II manual for instructions to transform the appropriate competent *E. coli* host. Make sure to digest isolated DNA from positive clones with the appropriate restriction enzymes to determine the orientation of the PCR fragment.

Map and Features of the pENTR™ Vectors

Map of the pENTR™ Vectors

The map below shows the features of the pENTR™ vectors. Maps and a complete sequence for each pENTR™ vector are available for downloading from our Web site (www.invitrogen.com) or by contacting Technical Service (page 20).



Comments for pENTR™1A 2717 nucleotides

rrnB T1 transcription termination sequence: bases 106-149

rrnB T2 transcription termination sequence: bases 281-308

attL1: bases 358-457 (complementary strand)

ccdB gene: bases 612-917

attL2: bases 946-1045

Kanamycin resistance gene: bases 1168-1977

pUC origin: bases 2041-2714

* There is a unique *Ehe* I site but no *Dra* I site in pENTR™2B.
There is a unique *Nco* I site but no *Dra* I site in pENTR™4.
There is a unique *Nsp* V site but no *Dra* I site in pENTR™11.

** There is a unique *Nco* I site between the *Xmn* I site and the *Sal* I site in pENTR™11 only.

continued on next page

Map and Features of the pENTR™ Vectors, continued

Features of the pENTR™ Vectors

pENTR™1A (2717 bp), pENTR™2B (2718 bp), pENTR™3C (2723 bp), pENTR™4 (2720 bp), and pENTR™11 (2744 bp) contain the following elements. All features have been functionally tested.

Feature	Benefit
<i>rrnB</i> T1 and T2 transcription termination sequences	Protects the cloned gene from expression by vector-encoded promoters, thereby reducing possible toxicity (Orosz <i>et al.</i> , 1991)
<i>attL</i> 1 and <i>attL</i> 2 sites	Allows site-specific recombination of the entry clone with a Gateway® destination vector (Landy, 1989)
<i>ccdB</i> gene	Allows negative selection of expression clones
Kanamycin resistance gene	Allows selection of the plasmid in <i>E. coli</i>
pUC origin	Allows high-copy number replication and growth in <i>E. coli</i>

Technical Service

World Wide Web



Visit the Invitrogen Web Resource using your World Wide Web browser. At the site, you can:

- Get the scoop on our hot new products and special product offers
- View and download vector maps and sequences
- Download manuals in Adobe® Acrobat® (PDF) format
- Explore our catalog with full color graphics
- Obtain citations for Invitrogen products
- Request catalog and product literature

Once connected to the Internet, launch your Web browser (Internet Explorer 5.0 or newer or Netscape 4.0 or newer), then enter the following location (or URL):

<http://www.invitrogen.com>

...and the program will connect directly. Click on underlined text or outlined graphics to explore. Don't forget to put a bookmark at our site for easy reference!

Contact Us

For more information or technical assistance, please call, write, fax, or email. Additional international offices are listed on our Web page (www.invitrogen.com).

Corporate Headquarters:

Invitrogen Corporation
1600 Faraday Avenue
Carlsbad, CA 92008 USA
Tel: 1 760 603 7200
Tel (Toll Free): 1 800 955 6288
Fax: 1 760 602 6500
E-mail:
tech_service@invitrogen.com

European Headquarters:

Invitrogen Ltd
Inchinnan Business Park
3 Fountain Drive
Paisley PA4 9RF, UK
Tel: +44 (0) 141 814 6100
Fax: +44 (0) 141 814 6117
E-mail:
eurotech@invitrogen.com

continued on next page

Technical Service, continued

MSDS Requests

To request an MSDS, visit our Web site (www.invitrogen.com). On the home page, go to 'Technical Resources', select 'MSDS', and follow instructions on the page.

Limited Warranty

Invitrogen is committed to providing our customers with high-quality goods and services. Our goal is to ensure that every customer is 100% satisfied with our products and our service. If you should have any questions or concerns about an Invitrogen product or service, please contact our Technical Service Representatives.

Invitrogen warrants that all of its products will perform according to the specifications stated on the certificate of analysis. The company will replace, free of charge, any product that does not meet those specifications. This warranty limits Invitrogen Corporation's liability only to the cost of the product. No warranty is granted for products beyond their listed expiration date. No warranty is applicable unless all product components are stored in accordance with instructions. Invitrogen reserves the right to select the method(s) used to analyze a product unless Invitrogen agrees to a specified method in writing prior to acceptance of the order.

Invitrogen makes every effort to ensure the accuracy of its publications, but realizes that the occasional typographical or other error is inevitable. Therefore Invitrogen makes no warranty of any kind regarding the contents of any publications or documentation. If you discover an error in any of our publications, please report it to our Technical Service Representatives.

Invitrogen assumes no responsibility or liability for any special, incidental, indirect or consequential loss or damage whatsoever. The above limited warranty is sole and exclusive. No other warranty is made, whether expressed or implied, including any warranty of merchantability or fitness for a particular purpose.

Purchaser Notification

**Limited Use
Label License
No. 19:
Gateway®
Cloning
Products**

This product and its use is the subject of one or more of U.S. Patent Nos. 5,888,732, 6,143,557, 6,171,861, 6,270,969, and 6,277,608 and/or other pending U.S. and foreign patent applications owned by Invitrogen Corporation. The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for profit entity). The purchase of this product does not convey a license under any method claims in the foregoing patents or patent applications, or to use this product with any recombination sites other than those purchased from Invitrogen Corporation or its authorized distributor. The right to use methods claimed in the foregoing patents or patent applications with this product for research purposes only can only be acquired by the use of Clonase™ purchased from Invitrogen Corporation or its authorized distributors. The buyer cannot modify the recombination sequence(s) contained in this product for any purpose. The buyer cannot sell or otherwise transfer (a) this product, (b) its components, or (c) materials made by the employment of this product or its components to a third party or otherwise use this product or its components or materials made by the employment of this product or its components for Commercial Purposes. The buyer may transfer information or materials made through the employment of this product to a scientific collaborator, provided that such transfer is not for any Commercial Purpose, and that such collaborator agrees in writing (a) not to transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes. Notwithstanding the preceding, any buyer who is employed in an academic or government institution may transfer materials made with this product to a third party who has a license from Invitrogen under the patents identified above to distribute such materials. Transfer of such materials and/or information to collaborators does not convey rights to practice any methods claimed in the foregoing patents or patent applications.

continued on next page

Purchaser Notification, continued

Limited Use Label License No. 19: Gateway® Cloning Products, continued

Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. Invitrogen Corporation will not assert a claim against the buyer of infringement of the above patents based upon the manufacture, use or sale of a therapeutic, clinical diagnostic, vaccine or prophylactic product developed in research by the buyer in which this product or its components was employed, provided that none of (i) this product, (ii) any of its components, or (iii) a method claim of the foregoing patents, was used in the manufacture of such product. Invitrogen Corporation will not assert a claim against the buyer of infringement of the above patents based upon the use of this product to manufacture a protein for sale, provided that no method claim in the above patents was used in the manufacture of such protein. If the purchaser is not willing to accept the limitations of this limited use statement, Invitrogen is willing to accept return of the product with a full refund. For information on purchasing a license to use this product for purposes other than those permitted above, contact Licensing Department, Invitrogen Corporation, 1600 Faraday Avenue, Carlsbad, California 92008. Phone (760) 603-7200.

Gateway® Clone Distri- bution Policy

For additional information about Invitrogen's policy for the use and distribution of Gateway® clones, see the section entitled **Gateway® Clone Distribution Policy**, page 24.

Gateway[®] Clone Distribution Policy

Introduction

The information supplied in this section is intended to provide clarity concerning Invitrogen's policy for the use and distribution of cloned nucleic acid fragments, including open reading frames, created using Invitrogen's commercially available Gateway[®] Technology.

Gateway[®] Entry Clones

Invitrogen understands that Gateway[®] entry clones, containing *attL1* and *attL2* sites, may be generated by academic and government researchers for the purpose of scientific research. Invitrogen agrees that such clones may be distributed for scientific research by non-profit organizations and by for-profit organizations without royalty payment to Invitrogen.

Gateway[®] Expression Clones

Invitrogen also understands that Gateway[®] expression clones, containing *attB1* and *attB2* sites, may be generated by academic and government researchers for the purpose of scientific research. Invitrogen agrees that such clones may be distributed for scientific research by academic and government organizations without royalty payment to Invitrogen. Organizations other than academia and government may also distribute such Gateway[®] expression clones for a nominal fee (\$10 per clone) payable to Invitrogen.

Additional Terms and Conditions

We would ask that such distributors of Gateway[®] entry and expression clones indicate that such clones may be used only for research purposes, that such clones incorporate the Gateway[®] Technology, and that the purchase of Gateway[®] Clonase[™] from Invitrogen is required for carrying out the Gateway[®] recombinational cloning reaction. This should allow researchers to readily identify Gateway[®] containing clones and facilitate their use of this powerful technology in their research. Use of Invitrogen's Gateway[®] Technology, including Gateway[®] clones, for purposes other than scientific research may require a license and questions concerning such commercial use should be directed to Invitrogen's licensing department at 760-603-7200.

References

Ausubel, F. M., Brent, R., Kingston, R. E., Moore, D. D., Seidman, J. G., Smith, J. A., and Struhl, K. (1994). *Current Protocols in Molecular Biology* (New York: Greene Publishing Associates and Wiley-Interscience).

Gold, L. (1988). Posttranscriptional Regulatory Mechanisms in *Escherichia coli*. *Ann. Rev. Biochem.* 57, 199-233.

Kozak, M. (1987). An Analysis of 5'-Noncoding Sequences from 699 Vertebrate Messenger RNAs. *Nucleic Acids Res.* 15, 8125-8148.

Kozak, M. (1991). An Analysis of Vertebrate mRNA Sequences: Intimations of Translational Control. *J. Cell Biology* 115, 887-903.

Kozak, M. (1990). Downstream Secondary Structure Facilitates Recognition of Initiator Codons by Eukaryotic Ribosomes. *Proc. Natl. Acad. Sci. USA* 87, 8301-8305.

Landy, A. (1989). Dynamic, Structural, and Regulatory Aspects of Lambda Site-specific Recombination. *Ann. Rev. Biochem.* 58, 913-949.

Miller, J. H. (1992). *A Short Course in Bacterial Genetics: A Laboratory Manual and Handbook for Escherichia coli and Related Bacteria* (Plainview, New York: Cold Spring Harbor Laboratory Press).

Orosz, A., Boros, I., and Venetianer, P. (1991). Analysis of the Complex Transcription Termination Region of the *Escherichia coli* *rrnB* Gene. *Eur. J. Biochem.* 201, 653-659.

Sambrook, J., Fritsch, E. F., and Maniatis, T. (1989). *Molecular Cloning: A Laboratory Manual*, Second Edition (Plainview, New York: Cold Spring Harbor Laboratory Press).

©1999-2004 Invitrogen Corporation. All rights reserved.

For research use only. Not intended for any animal or human therapeutic or diagnostic use.

Notes

Notes

Notes



Corporate Headquarters:

*Invitrogen Corporation
1600 Faraday Avenue
Carlsbad, California 92008
Tel: 1 760 603 7200
Tel (Toll Free): 1 800 955 6288
Fax: 1 760 603 7229
Email: tech_service@invitrogen.com*

European Headquarters:

*Invitrogen Ltd
3 Fountain Drive
Inchinnan Business Park
Paisley PA4 9RF, UK
Tel (Free Phone Orders): 0800 269 210
Tel (General Enquiries): 0800 5345 5345
Fax: +44 (0) 141 814 6287
Email: eurotech@invitrogen.com*

International Offices:

*Argentina 5411 4556 0844
Australia 1 800 331 627
Austria 0800 20 1087
Belgium 0800 14894
Brazil 0800 11 0575
Canada 800 263 6236
China 10 6849 2578
Denmark 80 30 17 40*

*France 0800 23 20 79
Germany 0800 083 0902
Hong Kong 2407 8450
India 11 577 3282
Italy 02 98 22 201
Japan 03 3663 7974
The Netherlands 0800 099 3310
New Zealand 0800 600 200
Norway 00800 5456 5456*

*Spain & Portugal 900 181 461
Sweden 020 26 34 52
Switzerland 0800 848 800
Taiwan 2 2651 6156
UK 0800 838 380
For other countries see our Web site*

www.invitrogen.com

