



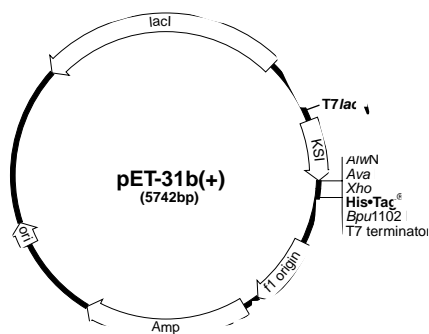
Table of Contents

I. Description	1
II. System Components	3
III. Construction of Target Peptide _n -His•Tag® Genes	3
IV. Expression of KSI-Target Peptide _n -His•Tag Fusion Protein	6
V. Purification by His•Bind® Chromatography and CNBr Cleavage	6
VI. Preparation of Peptide Conjugates and Affinity Supports	7
VII. References	8
Appendix A: Polyacrylamide Gel Purification of Oligonucleotides	9
Appendix B: pET-31b(+) Vector Maps and Sequence	10

Copyright © 1994 by Novagen, Inc. All rights reserved.

I. Description

pET Peptide Expression System 31	69953-3
pET Peptide Expression System 31 plus Competent Cells	69958-3
pET-31b(+) DNA, <i>Alw</i> N I digested, dephosphorylated	69954-3



The pET Peptide Expression System 31 is designed for the production of peptides in *E. coli*. Based on the pET-31b(+) vector, the system has several features that facilitate the expression and purification of target peptides (1). Peptide coding sequences are placed downstream of a 125aa ketosteroid isomerase (KSI) gene and upstream of a His•Tag sequence. Unique *Alw*N I and *Xho* I cloning sites are available. Use of the *Alw*N I site allows unidirectional insertion of coding sequences immediately adjacent to a methionine residue. For high-yield production, small (10-25aa) peptide coding sequences can be cloned as tandem repeats interspersed with single methionine residues. Long peptides (25-75aa) can be expressed in high yield from constructs containing single copies of coding sequences. The KSI-peptide-His•Tag fusion protein is expressed at high levels in the presence of a source of T7 RNA polymerase in the *recA*⁻, protease-deficient BLR(DE3)pLysS host. The fusion protein is purified in one step under denaturing conditions using Novagen's His•Bind metal chelation resin and then cleaved with CNBr to release monomer peptide units each terminating with a homoserine lactone, along with insoluble KSI and the His•Tag peptide. The HS-lactone can easily be converted to an amide or homoserine, or reacted with other compounds such as analogs of fluorescein and biotin to produce C-terminal derivatized peptides. The system has been used to make biologically active peptides of 14-67aa at yields of >50mg purified peptide per liter of culture (1).

The figure on p. 2 shows the cloning strategy for insertion of tandem repeats of peptide coding regions into the *Alw*N I site of pET-31b(+). Complementary oligonucleotides corresponding to the coding region are designed such that they lack internal methionine codons and produce the three-base 3' overhangs shown in the annealed molecule. The oligonucleotides are 5' phosphorylated to allow the formation of multimeric copies upon ligation. Products corresponding to the desired number of tandem copies (usually 3-6) are isolated by agarose gel electrophoresis, and then ligated into the *Alw*N I digested, dephosphorylated vector.

The resulting construct carries several tandem copies of the target peptide coding sequence interspersed with and flanked by single methionine codons. The initial vector + insert ligation is transformed into a non-expression host such as NovaBlue to allow the establishment of stable plasmid recombinants. The plasmid is then transferred into the BLR(DE3)pLysS expression host for target protein production upon the addition of IPTG to the culture.



pET Peptide Expression System 31

The KSI moiety of the fusion protein is extremely insoluble and causes the formation of large peptide-containing inclusion bodies in the host cell. The inclusion bodies are isolated by centrifugation, washed, and the fusion protein is purified by His•Bind[®] metal chelation chromatography under denaturing conditions. Elution is achieved by raising the imidazole concentration in the buffer. The subsequent steps vary somewhat depending on the characteristics of the target peptide. In general, the purified fusion protein is cleaved with CNBr under standard conditions, followed by differential precipitation of the extremely hydrophobic KSI protein to yield pure peptide. Two alternative protocols, suitable for purification of hydrophobic peptides and hydrophilic peptides, are provided in Section V.

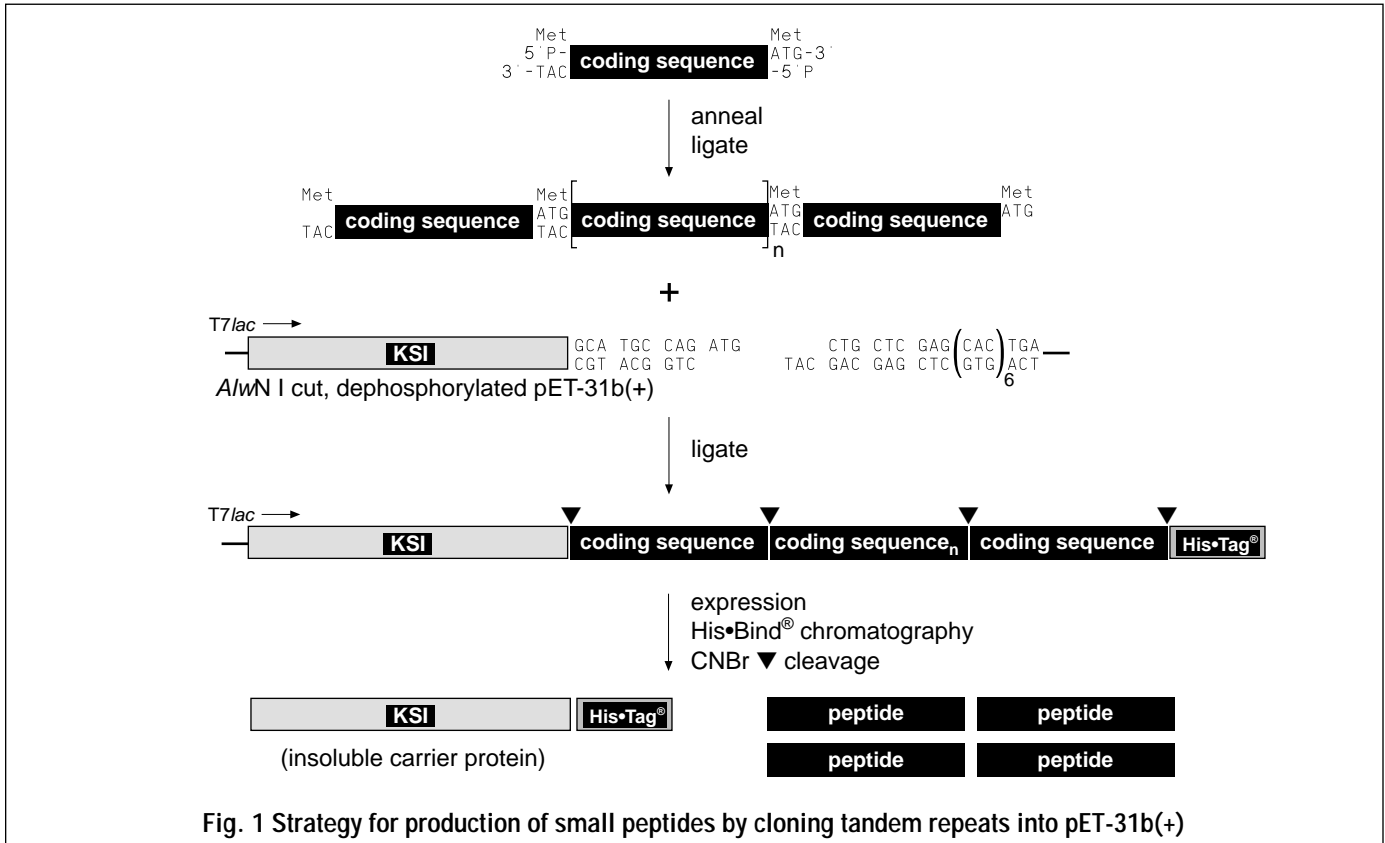


Fig. 1 Strategy for production of small peptides by cloning tandem repeats into pET-31b(+)



II. System Components

The **pET Peptide Expression System 31** is available as a choice of kits with individual components available separately. The basic system contains:

- pET-31b(+) DNA, 10 μ g
- Host bacterial strains, BLR and BLR(DE3)pLysS, glycerol stocks
- Induction control clone H, glycerol stock
- His•Bind[®] Resin, 10ml
- 4 polypropylene chromatography columns
- Vector map, pET Manual and pET Peptide Expression System 31 protocols
- Novagen Vector Diskette containing all Novagen plasmid sequences (Macintosh or DOS format)

The **pET Peptide Expression System 31 plus Competent Cells** contains two host strains ready for high-efficiency transformation. One 0.2ml aliquot each of the initial cloning host NovaBlue and the expression host BLR(DE3)pLysS, plus SOC medium, are included.

A kit containing 2 μ g ***AIwN I* digested dephosphorylated pET-31b(+) DNA** and a Control Insert is also available (Cat. No. 69954-1).

Other Supplemental Products

- | | | |
|---------------------------------|------------|---------|
| • pET-31b(+) DNA (uncut) | 10 μ g | 69952-3 |
| • BLR(DE3)pLysS Competent Cells | 0.4ml | 69956-3 |
| | 1ml | 69956-4 |
| • His•Bind Resin | 10ml | 69670-3 |
| | 50ml | 69670-4 |
| • His•Bind Buffer Kit | | 69755-3 |

III. Construction of Target Peptide_n-His•Tag Genes

The procedure used to create the tandem repeat of the target peptide fused to the C-terminus of the 125 codons for the bacterial ketosteroid isomerase (KSI) gene is shown in Figure 1. The sense and anti-sense oligonucleotide strands encoding the target peptide should be designed using the preferred codon usage of *E. coli* (2). In general, the tandem repeat strategy is recommended for peptides up to 25 amino acids (75bp). Larger peptides may be produced in suitable yield from monomer constructs. Internal methionine residues must be mutated to other conserved residues in order to avoid cyanogen bromide cleavage at non-junctional methionines. In several cases, we have found that mutation of these internal methionines to isoleucine has no measurable effect on bioactivity (1). Tandem peptide construction requires that the 3' end of the coding strand include a 3-base ATG extension, and the 3' end of the non-coding strand include a 3-base TAC extension to create suitable overhangs in the duplex for unidirectional end-to-end self ligation.

The following protocols call for phosphorylation/ligation of the oligos on a relatively large scale. Sufficient quantities (25 μ g, or about 1 A₂₆₀ unit) of highly pure oligonucleotides help to ensure that the ligations proceed efficiently and that sufficient material will be recovered from gel fractionation. HPLC or gel purified oligonucleotides are suitable. A procedure for gel purification is provided in Appendix A. Note that it is also necessary to use 5' phosphorylated oligonucleotides to allow tandem ligation, and for ligation into the *AIwN I* digested dephosphorylated pET-31b(+) vector available from Novagen.

Phosphorylation

1. If the purified oligos have already been phosphorylated proceed to the annealing reaction below. If they are not phosphorylated assemble two reactions, each containing one oligo strand, as follows. We recommend Novagen's Polynucleotide Kinase (Cat. No. 69248-1, which includes 10X buffer and 10mM ATP).

25 μ g (~1500-2000pmol) purified oligo in TE buffer (10mM Tris-HCl pH 8.0, 1mM EDTA)



10 μ l 10X Kinase buffer
10 μ l 10mM ATP
125U T4 polynucleotide kinase
x μ l Nuclease-free water
100 μ l Total volume

2. Add the enzyme last, stir gently to mix and incubate at 37°C for 30 minutes.

Annealing

1. Combine the two phosphorylated oligos into one tube (total volume 200 μ l). Annealing can be performed in the phosphorylation reaction buffer. If pre-phosphorylated oligos are used, dissolve them in TE buffer or water at ~100pmol/ μ l and add equimolar amounts (1500-2000pmol) of each to one tube. Add 20 μ l of 10X annealing buffer (10X = 400mM Tris-HCl pH 8.0, 100mM MgCl₂, 500mM NaCl) and water to a total volume of 200 μ l.
2. Heat the combined oligos at 99°C for 10 minutes, then cool to 30°C over a 15 minute period. This procedure is conveniently performed in a thermal cycler.
3. Extract the reaction once with an equal volume of TE-buffered phenol:CIAA (1:1; CIAA is 24 parts chloroform and 1 part isoamyl alcohol). Extract the aqueous phase with 1 volume of CIAA. Transfer the final aqueous phase to a fresh tube. Add MgCl₂ to 10mM, 0.1 volume 3M Na acetate pH ~6 and 2 volumes of 100% ethanol. Vortex and place at -20°C for > 1 hour. Centrifuge at 12,000 \times g for 5 minutes, drain, rinse the pellet with 100% ethanol and air dry. Resuspend the annealed oligos in TE buffer at 100pmol/ μ l. Remember that 1pmol of each strand when combined equals 1pmol annealed oligo!

Ligation and purification of tandem repeats

1. Assemble the ligation reaction as follows (qualified ligation components are included in the DNA Ligation Kit, Cat. No. 69838-3):
2.5 μ l phosphorylated, annealed oligos (250pmol)
2.5 μ l 10X Ligase buffer (200mM Tris-HCl pH 7.6, 50mM MgCl₂)
2.5 μ l 100mM DTT
0.5 μ l 10mM ATP
0.5 μ l T4 DNA ligase (2 Weiss units)
16.5 μ l Nuclease-free water
25 μ l Total volume
2. Add the enzyme last and mix gently by stirring with the pipet tip. Incubate the reaction at 16°C for 2 hours. The extent of ligation can be checked by gel analysis of 2 μ l (~0.5 μ g) sample of the reaction.
3. Add concentrated gel loading buffer to 1X. Load the entire sample in a 1cm \times 1.5mm well of a 3% low melting point agarose gel containing 0.5 μ g/ml ethidium bromide. Run appropriate size markers (e.g. Novagen's PCR Markers, Cat. No. 69278-1) in an adjacent lane. Run the gel at 70V for 3 hr at 4°C (voltage and time given for standard 10cm long minigel apparatus).
4. Visualize the DNA bands under UV illumination (~300-360nm). Cut out bands containing the desired multimer species (several repeat sizes can be tested) and recover the DNA by standard melting/phenol extraction procedures (e.g. Sambrook et al. (1989) in "Molecular Cloning, A Laboratory Manual", 2nd edition, pp. 6.30-6.31).
5. Resuspend the final purified DNA in ~10 μ l TE buffer and determine its concentration spectrophotometrically or by gel analysis of a sample vs. known amounts of marker DNA. The DNA should now be ready for ligation with the pET-31b(+) vector. Calculate the molar concentration of the DNA as follows:
$$A_{260} \times \text{dilution factor} \times 50\mu\text{g/ml} = x \mu\text{g/ml}$$
$$x \mu\text{g/ml} \div (660\mu\text{g}/\mu\text{mol}/\text{bp} \times \#\text{bp}) = y \mu\text{mol/ml}$$
$$y \mu\text{mol/ml} \times 1/1000 = z \text{pmol}/\mu\text{l}$$

Note: If difficulties are encountered in producing the desired tandem species (as monitored in Step 2) or in recovering sufficient material for cloning, use more annealed oligos (up to 50 μ g, or 2000pmol) and more ligase (15U) in the 25 μ l



ligation reaction. In addition, increase the incubation time to 16 hours and reduce the temperature to 8°C. The extent of ligation may be further increased by adding more ligase (10U) and ATP (2µl 10mM) at this point and allowing the reaction to proceed for another 24 hr at 8°C.

Ligation with *A/wN* I digested dephosphorylated pET-31b(+)

The pET-31b(+) vector is prepared by *A/wN* I digestion and dephosphorylation. Qualified vector ready for ligation of tandem inserts and a positive control insert are available as a kit (Cat. No. 69954-1). Each different size tandem DNA multimer is ligated with the vector in separate reactions to create an array of pET-31(target peptide)_nHis•Tag[®] constructs.

1. Assemble the following ligation reaction:

2µl (0.026pmol) *A/wN* I digested, dephosphorylated pET-31b(+)
1µl 10X Ligase buffer (200mM Tris-HCl pH 7.6, 50mM MgCl₂)
1µl 100mM DTT
0.5µl 10mM ATP
xµl (0.13-0.26pmol) purified tandem DNA insert
0.5µl (2 Weiss units) T4 DNA ligase
yµl Nuclease-free water
10µl Total volume

In addition, set up a negative control reaction containing all of the same components except for the target insert (add an equivalent amount of TE buffer instead). This reaction will allow the establishment of vector background. A positive control reaction may also be set up, which should contain a known amount of target DNA that has been successfully used for cloning (see below for information on the Positive Control insert supplied with the *A/wN* I digested dephosphorylated vector).

2. Add the enzyme last and mix by stirring with the pipet tip. Incubate at 16°C overnight.
3. Transform competent NovaBlue cells and screen colonies for inserts.

Positive Control

A Positive Control insert is provided with the *A/wN* I digested, dephosphorylated vector to verify the cloning steps. The insert encodes a 15aa peptide corresponding to the S•Tag[™] peptide. It is provided as a 5' phosphorylated monomer with appropriate overhangs for *A/wN* I site ligation. Use 2µl (0.2pmol) of the insert for ligation. Note that this insert is provided for testing purposes only and may not be used for peptide production or distributed outside your laboratory.

Protocols for transformation and colony screening are found in the accompanying pET System Manual. Screen potential recombinants by colony PCR with the KSI primer (Cat. No. 69947-3) and T7 terminator primer (Cat. No. 69337-3) combination. The presence of inserts and number of target peptide encoding units can be determined by agarose gel electrophoresis of the PCR products. The vector without an insert produces a 158bp amplification product with the indicated primers. Candidate recombinants should also be sequenced to verify the construction. The appropriate constructs are then transformed into BLR(DE3)pLysS for protein overproduction.

The ligations containing the target oligo and Positive Control insert should produce many more colonies than the negative control ligation. If the target insert ligation produces nearly the same number of colonies as the negative control, cloning efficiency is low and screening will likely reveal a large percentage of clones containing only the religated vector. Instead of preparing more insert, it may be possible to obtain the desired clone by digesting the remainder of the ligation with *A/wN* I prior to transformation. This treatment will reduce the background of religated vector and should therefore increase the likelihood of finding the desired recombinant.



IV. Expression of KSI-Target Peptide_nHis•Tag[®] Fusion

A detailed discussion of induction strategies, culture maintenance and storage is found in the pET System Manual. The protease deficient strain BLR(DE3)pLysS should be used for protein production. The BLR strain is a *recA*⁻ derivative of BL21 and is potentially superior for stabilizing tandem repeats.

1. The day before induction, inoculate 25-50ml of LB media supplemented with 50µg/ml carbenicillin (Carb) or ampicillin and 34µg/ml chloramphenicol (Cam) in a 250ml flask with a single colony of the recombinant (grown on LB agar plus antibiotics). Shake at 250rpm at 37°C overnight.
2. The next morning, dilute the starter culture 40-fold into 500ml of fresh LB/Carb/Cam in a 2.8 liter Fernbach flask prewarmed to 37°C, using as many flasks as needed to suit the scale of the production.
3. Shake at 250rpm at 37°C until the OD₆₀₀ reaches 0.3-0.5. Induce the production of the target protein by adding 1mM IPTG. Harvest the cells (by centrifugation) at an OD ~ 2.0, which is generally reached 3-6 hr after induction. The drained cell pellet (~2.5g/l) can be stored at -20°C for at least one week without any loss in yield of fusion protein.
4. Resuspend the cell pellet in 0.1X the culture volume of 1X Binding Buffer (5mM imidazole, 40mM Tris-HCl pH 7.9, 500mM NaCl), sonicate on ice until no longer viscous, and centrifuge at 12,000 × g for 10 min at 4°C. If desired, remove a small sample prior to centrifugation for gel analysis of total cell protein.
5. Carefully remove the supernatant (soluble) fraction to a fresh tube. If desired, the pellet (insoluble) fraction can be washed by brief sonication in 1X Binding Buffer, recentrifuged as above, and drained. Resuspend the pellet in 0.1 culture volume of 1X Binding Buffer containing 6M guanidine-HCl. After the pellet has dissolved completely, centrifuge again at 12,000 × g for 10 min at 4°C to remove any residual particulate matter.
6. Since the KSI fusion partner is very insoluble, it is likely that the target fusion protein will be found in the insoluble fraction, which is the pellet from the centrifugation in Step 4. The supernatant, pellet and pre-centrifugation sample (if saved) can be quickly analyzed for the target protein by SDS-gel electrophoresis and staining with Coomassie blue. The following formula can be applied to determine an appropriate amount of sample to load on a standard mini-gel apparatus. Note that the presence of salt and guanidine may affect the mobility of the target protein; if this is crucial follow the protocol in the pET Manual for processing a portion of the culture in a low-salt buffer.
$$(30/\text{concentration factor}) \times (3/\text{OD}_{600} \text{ at harvest}) = x$$

Add xµl of sample to an equal volume of 2X SDS sample buffer.
Use the same amount of soluble, insoluble and total samples described above.

V. Purification by His•Bind[®] Chromatography and CNBr Cleavage

The following protocols are based on processing 200ml of *E. coli* culture. Scale all steps in proportion to the amount of culture used. All steps are performed at room temperature unless otherwise noted. **Use appropriate safety precautions when handling CNBr and other hazardous chemicals during this procedure.**

Protocol 1: Production and purification of hydrophobic peptides

The insoluble cellular protein pellet will generally contain the majority of the fusion protein as dense inclusion bodies. If gel analysis of the soluble and insoluble fractions shows that the target protein is in the redissolved inclusion bodies, His•Bind chromatography is carried out under denaturing conditions.

1. Load the resolubilized inclusion bodies from Step 5 above onto a 2.5ml His•Bind column which has been charged with NiSO₄ and equilibrated with 1X Binding Buffer (1X = 20mM Tris-HCl pH 7.9, 0.5M NaCl, 5mM imidazole) plus 6M guanidine-HCl, at a flow rate of 0.25ml/min.



2. Wash the column with 15ml of 1X Binding Buffer plus 6M guanidine-HCl and then with 20ml of the same buffer containing 16mM imidazole instead of 5mM. An 8X stock of this solution (without guanidine-HCl) can be prepared by adding 4 parts 8X Binding Buffer plus 1 part 8X Wash Buffer supplied in the His•Bind Buffer Kit. Collect and save the flow-through and wash fractions.
3. Elute the fusion protein with 15ml of 300mM imidazole in the same buffer. (A 4X stock of this elution buffer without guanidine-HCl can be prepared from His•Bind Buffer kit components as follows: first dilute the kit's 8X Binding Buffer with an equal volume of water to bring it to 4X, then add 2.35 parts of 4X Binding Buffer to 1 part of the kit's 4X Elute Buffer.) Collect 2ml fractions. Analyze samples of each fraction by SDS-polyacrylamide gel electrophoresis. Combine the peak fractions (ca. 10ml), dialyze overnight at 4°C against 2 × 1 liter of H₂O in 12-14 kDa cut-off dialysis bags. The majority of the protein should form a white precipitate which can be pelleted by centrifugation at 2000 × g for 10 min at 4°C.
4. Working in a fume hood, dissolve the dense white pellet in 6ml of 80% formic acid. Transfer to a 50ml round-bottomed flask and add 0.2g CNBr. Bubble in nitrogen, wrap the flask in aluminum foil and stir for 18-22 hr. Attach to a rotary evaporator and evaporate to dryness at 28°C. Resuspend the resultant transparent proteinaceous gel in a minimal volume of 40% CH₃CN/60% H₂O/0.1 % TFA (ca. 2-4ml) and stir for 1 hr. Centrifuge this suspension at 12,000 × g for 10 min at 4°C. The supernatant can be further clarified by passage through a 0.22µm filter and the purity of the recombinant target peptide analyzed by reverse-phase HPLC and UV spectroscopy.

Protocol 2: Production and purification of hydrophilic peptides

The purification and CNBr cleavage of the target peptide from the carrier fusion protein is the same as in Protocol 1 until the CNBr cleavage step. Here the two protocols diverge due to the differing solubility between the hydrophilic and the hydrophobic target peptides. After the CNBr cleaved peptide mixture is dried by rotary evaporation, the gelatinous material is resuspended in 20mM KH₂PO₄, 100mM NaCl, and the pH adjusted to 7.4 with 1M NaHCO₃. This mixture is stirred overnight under N₂ and wrapped in aluminum foil. The suspension is centrifuged at 5000 × g for 15 min at 4°C and the peptide in the supernatant analyzed by UV spectroscopy. The hydrophilic peptide in the supernatant can be either concentrated by rotary evaporation or dried by lyophilization.

VI. Preparation of Peptide Conjugates and Affinity Supports

CNBr cleavage of the tandem peptides at the junctional methionines will generate a C-terminal homoserine lactone. In most cases, the bioactivity should be unaffected by the presence of this extra amino acid. One may wish to take advantage of this uniquely reactive residue and couple a wide variety of primary amine-containing compounds such as biotin, fluorescein, and other fluorescent reagents to the C-terminus of the target peptide, or prepare C-terminally coupled peptide affinity supports.

Preparation of peptide-homoserine conjugates

In the simplest case, the lactone ring can be completely and immediately converted to the homoserine (HS) free acid form by dropwise addition of 0.2N NaOH, until a final pH of 12.5–13 is reached and then neutralized with phosphate-buffered saline, pH 7.4. For each of the following reactions, 0.3-0.8mg of peptide-lactone are placed in 1.7ml polypropylene tubes and completely lactonized by addition of 20µl of 100% trifluoroacetic acid (TFA) and immediately dried in a Speed Vac (Savant). The dried pellets are dissolved in 50µl anhydrous DMF delivered by a gas-tight syringe and 8µl of Et₃N are added. All reaction tubes are equipped with a micro-stir bar.

Synthesis of peptide-HS-amide: 50µl of anhydrous DMF saturated with NH₃ gas are added directly to the dried lactonized pellet prior to the addition of Et₃N. This reaction is stirred at 44°C and amidation is complete in 30 min.



Synthesis of peptide-HS-biotin: 100 equivalents of biotin ethylenediamine (Molecular Probes) are added to the lactonized peptide solution and the reaction allowed to proceed for 5 hr at 48°C.

Synthesis of peptide-HS-fluorescein: 28 equivalents of fluoresceinyl glycine amide (Molecular Probes) are dissolved in an additional 100µl DMF and added to the lactonized peptide solution and the reaction is allowed to proceed for 8 hr at 48°C.

Synthesis of peptide-HS-dansyl: 37 equivalents of dansyl ethylenediamine (Molecular Probes) are added to the lactonized peptide solution and the reaction is allowed to proceed for 7 hr at 48°C.

Synthesis of peptide-HS-ANS: 100 equivalents of 5((2-aminoethyl)amino)naphthalene-1-sulfonic acid (Molecular Probes) are dissolved in an additional 200µl DMF plus an additional 32µl Et₃N and added to the lactonized peptide solution and the reaction is allowed to proceed for 6 hr at 45°C.

Preparation of C-terminally coupled peptide-affinity supports

Four mg of lyophilized peptide-HS-lactone are placed into a 1.7ml polypropylene tube and solubilized with 765µl 0.88 M NaOAc, and the pH is adjusted to 6 with NaOH. 0.5ml of aminoalkyl Sepharose (Affigel 102, Bio Rad) is added and the mixture is gently rocked for 170 hr at 37°C. The extent of the coupling is monitored by removing small aliquots at various time points and determining the amount of peptide remaining in the supernatant by the Bradford Method.

VII. References

1. Kuliopulos, A. and Walsh, C.T. (1994) *J. Am. Chem. Soc.* **116**, 4599-4607.
2. Boer, H. A. d., and Kastelein, R. A. (W. Reznikoff and L. Gold, Ed.) in "Maximizing Gene Expression", Butterworths: Stoneham, MA, 1986; pp 225-285.



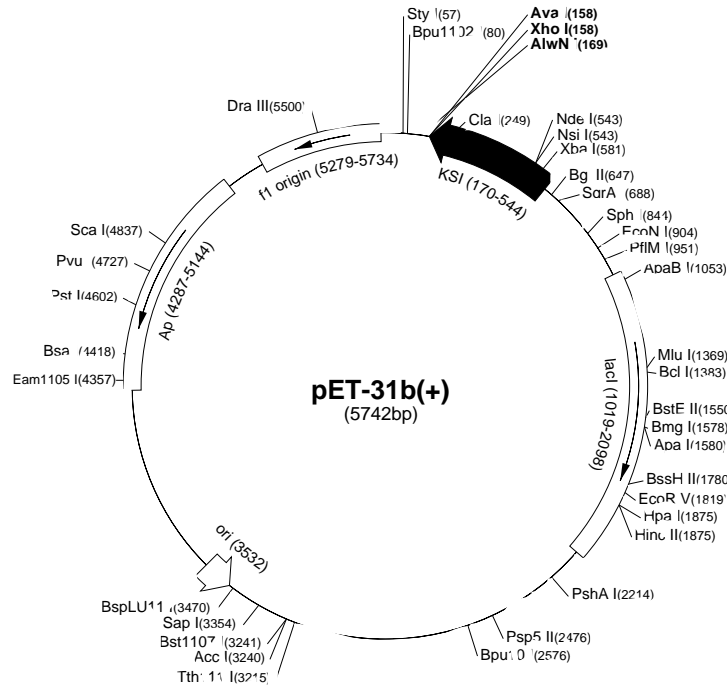
Appendix A: Polyacrylamide Gel Purification of Oligonucleotides

1. Synthesize oligonucleotide, determine concentration by reading the absorbance at 260nm. Use $20\mu\text{g/ml} = 1 A_{260}$
2. Pour 12% polyacrylamide minigel:
 - 4.2g urea
 - 1.0ml 10X TBE (Tris-Borate-EDTA)
 - 1.8ml water
 - 4.0ml 30% acrylamide/0.8% bisacrylamide Protogel (National Diagnostics) solution
 - 100 μl 10% ammonium persulfate
 - 5 μl TEMED
 - 10ml total volumeDegas gel, pour 0.75 mm gel, overlay with tert-amyl alcohol, let gel polymerize.
3. Rinse away urea prior to loading oligo. Load 200 μg of oligo in 1X DNA loading buffer with bromophenol blue dye, run at 200 volts \times ca. 30 min in minigel apparatus. Take apart gel, place on Saran wrap.
4. Using hand-held short-wave UV lamp with fluorescent-indicator TLC plate, excise dark band of interest (UV absorbing) with a clean razor blade. Using forceps, place gel slice into 15ml Falcon tube and mash-up gel slice with a glass rod. Add 10ml 100mM NH_4Ac , 1 mM EDTA, pH 6-6.5 and rock overnight at room temperature.
5. Pellet gel at $3000 \times g$ 1 min. Transfer supernatant to fresh Falcon tube.
6. Pre-equilibrate Waters C-18 Sep-Pak with 10ml MeOH, then 20ml H_2O . Slowly load 10ml of oligo-containing supernatant, avoid pushing bubbles through Sep-Pak at all times. Wash with 10ml H_2O . Elute very slowly with $2 \times 1\text{ml}$ 50% MeOH/50% H_2O into two 1.7ml polypropylene tubes labeled E1 and E2.
7. Take absorbance readings at 260nm of 50% MeOH/50% H_2O eluates E1 and E2, blanked against same solvent. Use $30\mu\text{g/ml} = 1 A_{260}$
8. Dry in Speed Vac.



pET Peptide Expression System 31

Appendix B: pET-31b(+) Maps and Sequence



<u>Bgl II</u>	<u>T7 promoter</u>	<u>lac operator</u>	<u>Xba I</u>	<u>rbs</u>
AGATCTCGATCCCGCAAATTAATACGACTC	ATAGGGGAATTGTGAGCGGATAACCAAT	TCCCCTCTAGAAATAATTTGTTAACTTTA	AAGAGGAGA	
<u>Nsi I</u>	<u>KSI primer #69947-3</u>	<u>KSI</u>	<u>AlwNI</u>	<u>Ava I</u>
TATACATATG...	105aa...GGCAAGGTGGTGAGCATCCGCGCCTT	GGCGAGAAGAATATTCACGCATGCCAGATGCTGCTCGAGCACCACCACCACCACTGA	Met...315bp...GlyLysValValSerIleArgAlaLeuPheGlyIuLysAsnIleHisAlaCysGlnMetLeuLeuGluHisHisHisHisHisEnd	
<u>Nde I</u>	<u>Bpu1102 I</u>	<u>T7 terminator</u>	<u>Xho I</u>	<u>His*Tag®</u>
GATCCGGCTGCTAACAAAGCCGAAAGGAAGCTGAGTTGGCTGCTGCCACCGCTGAGCAATAACTAGCATAACCCCTTGGGGCCTCTAAACGGGCTTTGAGGGGTTTTTTTG				
	<u>T7 terminator primer #69337-3</u>			

pET-31b(+) cloning/expression region

Unique sites are shown. Note that the sequence is numbered by the pBR322 convention, so the T7 expression region is reversed (*i.e.* the sequence below shows the reverse complement of the expressed strand; the above diagram shows the expressed strand). The f1 origin in pET-31b(+) is oriented such that the single stranded DNA produced upon helper phage infection will anneal with the T7 terminator primer.

pET-31b(+) Sequence

```

1  ATCCGGATAT AGTTCCTCCT TTCAGCAAAA AACCCCTCAA GACCCGTTTA
51  GAGGCCCAA GGGTTATGC TAGTTATTGC TCAGCGGTGG CAGCAGCCAA
101 CTCAGCTTCC TTTCCGGGCTT TGTTAGCAGC CGGATCTCAG TGGTGGTGGT
151 GGTGGTGCTC GAGCAGCATC TGGCATGCGT GAATATTCTT CTCGCCAAAC
201 AAGGCGCGGA TGCTCACCAC CTTGCCGGCG CCATTGAAGC GAAAGTGATC
251 GATGGGCGCA ACTACGGTCT TGCGGCCCTG AACTCTGAAG CTGACGGTGA
301 AAGCGAAGGC CGCTTCGTTG GCGACCGCGC GTACCTCCTG CGTCAGCTCC
351 ACCGCCAAAG GCAGTTTGAG CGAGTTGGCG TAAACTCAC GAATCGCAGC
401 CGTACCGGAC CTGGGCTCGG AACCCACGGG GTCTTCCACC GTGGCGTCAT
451 CGGCAAACAG CGCGACGATG CCGTCCAGAT CGCCGGCATT GAGCGCAGCC
501 ACAAAGCGCT GTACCACGGC GGTGATGTGT TCTGGGGTAT GCATATGTAT
551 ATCTCCTTCT TAAAGTTAAA CAAAATTATT TCTAGAGGGG AATTGTTATC
601 CGCTCACAAT TCCCTATAG TGAGTCGTAT TAATTCGCGG GGATCGAGAT
651 CTCGATCTC TACGCCGGAC GCATCGTGGC CGGCATCACC GGCGCCACAG
701 GTGCGGTTGC TGGCGCTAT ATCGCCGACA TCACCGATGG GGAAGATCGG
751 GCTCGCCACT TCGGGCTCAT GAGCGCTTGT TTCGGCGTGG GTATGGTGGC

```



```
801 AGGCCCCGTG GCCGGGGGAC TGTGGGGCGC CATCTCCTTG CATGCACCAT
851 TCCTTGCGGC GCGGGTGCTC AACGGCCTCA ACCTACTACT GGGCTGCTTC
901 CTAATGCAGG AGTCGCATAA GGGAGAGCGT CGAGATCCCG GACACCATCG
951 AATGGCGCAA AACCTTTCGC GGTATGGCAT GATAGCGCCC GGAAGAGAGT
1001 CAATTCAGGG TGGTGAATGT GAAACCAGTA ACGTTATACG ATGTGCGAGA
1051 GTATGCCGGT GTCTCTTATC AGACCGTTTC CCGCGTGGTG AACCAAGGCCA
1101 GCCACGTTTC TGCGAAAACG CGGGAAAAAG TGGAAAGCGC GATGGCGGAG
1151 CTGAATTACA TTCCCAACCG CGTGGCACAA CAACTGGCGG GCAAAACAGTC
1201 GTTGCTGATT GGCCTTGCCA CCTCCAGTCT GGCCCTGCAC GCGCCGTCGC
1251 AAATTGTCGC GCGGATTAAT TCTCGCGCCG ATCAACTGGG TGCCAGCGTG
1301 GTGGTGTGCGA TGGTAGAACG AAGCGGCGTC GAAGCCTGTA AAGCGGCGGT
1351 GCACAATCTT CTCGCGCAAC GCGTCAGTGG GCTGATCATT AACTATCCCG
1401 TGGATGACCA GGATGCCATT GCTGTGGAAG CTGCCTGCAC TAATGTTCCG
1451 GCGTTATTTT TTGATGTCTC TGACCAGACA CCCATCAACA GTATTATTTT
1501 CTCCCATGAA GACGGTACGC GACTGGGCGT GGAGCATCTG GTCGCATTGG
1551 GTCACCAGCA AATCGCGCTG TTAGCGGGCC CATTAAAGTT TGTCTCGGCG
1601 CGTCTGCGTC TGGCTGGCTG GCATAAATAT CTCACTCGCA ATCAAATTCA
1651 GCCGATAGCG GAACGGGAAG GCGACTGGAG TGCCATGTCC GGTTTTCAAC
1701 AAACCATGCA AATGCTGAAT GAGGGCATCG TTCCCACTGC GATGCTGGTT
1751 GCCAACGATC AGATGGCGCT GGGCGCAATG CGCGCCATTA CCGAGTCCGG
1801 GCTGCGCGTT GGTGCGGATA TCTCGGTAGT GGGATACGAC GATACCGAAG
1851 ACAGCTCATG TTATATCCCG CCGTTAACCA CCATCAAACA GGATTTTCGC
1901 CTGCTGGGGC AAACCAGCGT GGACCGCTTG CTGCAACTCT CTCAGGGCCA
1951 GGCGGTGAAG GGCAATCAGC TGTTCGCCGT CTCACTGGTG AAAAGAAAAA
2001 CCACCCTGGC GCCCAATACG CAAACCGCCT CTCCCCGCGC GTTGGCCGAT
2051 TCATTAATGC AGCTGGCACG ACAGGTTTCC GACTGGAAA GCGGGCAGTG
2101 AGCGCAACGC AATTAATGTA AGTTAGCTCA CTCATTAGGC ACCGGGATCT
2151 CGACCGATGC CTTGAGAGC CTTCAACCCA GTCAGCTCCT TCCGGTGGGC
2201 GCGGGGCATG ACTATCGTCG CCGCACTTAT GACTGTCTTC TTTATCATGC
2251 AACTCGTAGG ACAGGTGCCG GCAGCGCTCT GGGTCATTTT CCGCGAGGAC
2301 CGTTTTCGCT GGAGCGCGAC GATGATCGGC CTGTGCTTGG CGGTATTCGG
2351 AATCTTGCA CCCCCTGCTC AAGCCTTCGT CACTGGTCCC GCCACCAAAC
2401 GTTTCGGCGA GAAGCAGGCC ATTATCGCCG GCATGGCGGC CCCACGGGTG
2451 CGCATGATCG TGCTCCTGTC GTTGAGGACC CGGCTAGGCT GCGGGGGTTG
2501 CCTTACTGGT TAGCAGAATG AATCACCGAT ACGCGAGCGA ACGTGAAGCG
2551 ACTGCTGCTG CAAAACGTCT GCGACCTGAG CAACAACATG AATGGTCTTC
2601 GGTTTCCGTG TTTTCGTAAG TCTGGAACG CCGAAGTCAG CGCCCTGCAC
2651 CATTATGTTT CGGATCTGCA TCGCAGGATG CTGCTGGCTA CCCTGTGGAA
2701 CACCTACATC TGTATTAACG AAGCGCTGGC ATTGACCCTG AGTGATTTTT
2751 CTCTGGTCCC GCCGCATCCA TACCGCCAGT TGTTTACCCT CACAACGTTT
2801 CAGTAACCGG GCATGTTTCA CATCAGTAAC CCGTATCGTG AGCATCCTCT
2851 CTCGTTTCAT CGGTATCATT ACCCCATGTA ACAGAAATCC CCCTTACACG
2901 GAGGCATCAG TGACCAAACA GGA AAAAACC GCCCTTAACA TGGCCCGCTT
2951 TATCAGAAGC CAGACATTA CCGTTCTGGA GAAACTCAAC GAGCTGGACG
3001 CGGATGAACA GGCAGACATC TGTGAATCGC TTCACGACCA CGCTGATGAG
3051 CTTTACCGCA GCTGCCTCGC GCGTTTCGGT GATGACGGTG AAAACCTCTG
3101 ACACATGCAG CTCCCGGAGA CGGTCACAGC TTGTCTGTAA GCGGATGCCG
3151 GGAGCAGACA AGCCCGTCAG GCGCGCTCAG CGGGTGTTGG CGGGTGTCGG
3201 GGCGCAGCCA TGACCCAGTC ACGTAGCGAT AGCGGAGTGT ATACTGGCTT
```



pET Peptide Expression System 31

```
3251 AACTATGCGG CATCAGAGCA GATTGTA CTG AGAGTGCACC ATATATGCGG
3301 TGTGAAATAC CGCACAGATG CGTAAGGAGA AAATACCGCA TCAGGCGCTC
3351 TTCCGCTTCC TCGCTCACTG ACTCGCTGCG CTCGGTCGTT CGGCTGCGGC
3401 GAGCGGTATC AGCTCACTCA AAGGCGGTAA TACGGTTATC CACAGAATCA
3451 GGGGATAACG CAGGAAAGAA CATGTGAGCA AAAGGCCAGC AAAAGGCCAG
3501 GAACCGTAAA AAGGCCGCGT TGCTGGCGTT TTTCCATAGG CTCCGCCCCC
3551 CTGACGAGCA TCACAAAAAT CGACGCTCAA GTCAGAGGTG GCGAAACCCG
3601 ACAGGACTAT AAAGATACCA GGC GTTTCC CCGTGAAGCT CCCTCGTGCG
3651 CTCTCCTGTT CCGACCTGCG CGCTTACCGG ATACCTGTCC GCCTTTCTCC
3701 CTTCCGGGAA CGTGGCGCTT TCTCATAGCT CACGCTGTAG GTATCTCAGT
3751 TCGGTGTAGG TCGTTCGCTC CAAGCTGGGC TGTGTGCACG AACCCCCCGT
3801 TCAGCCCGAC CGCTGCGCCT TATCCGGTAA CTATCGTCTT GAGTCCAACC
3851 CGGTAAGACA CGACTTATCG CCACTGGCAG CTGGTAACAG GATTAGCAGA
3901 GCGAGGTATG TAGGCGGTGC TACAGAGTTC TTGAAGTGGT GGCCTAACTA
3951 CGGCTACACT AGAAGGACAG TATTTGGTAT CTGCGCTCTG CTGAAGCCAG
4001 TTACCTTCGG AAAAAGAGTT GGTAGCTCTT GATCCGGCAA ACAAAACCACC
4051 GCTGGTAGCG GTGGTTTTTT TGTGTTGCAAG CAGCAGATTA CGCGCAGAAA
4101 AAAAGGATCT CAAGAAGATC CTTTGATCTT TTCTACGGGG TCTGACGCTC
4151 AGTGGAACGA AAACCTACGT TAAGGGATTT TGGTCATGAG ATTATCAAAA
4201 AGGATCTTCA CCTAGATCCT TTTAAATTAA AAATGAAGTT TTAATCAAT
4251 CTAAAGTATA TATGAGTAAA CTTGGTCTGA CAGTTACCAA TGCTTAATCA
4301 GTGAGGCACC TATCTCAGCG ATCTGTCTAT TTCGTTTCATC CATAGTTGCC
4351 TGA CTCCCG TCGTGTAGAT AACTACGATA CGGGAGGGCT TACCATCTGG
4401 CCCCAGTGCT GCAATGATAC CGCGAGACCC ACGCTCACCG GCTCCAGATT
4451 TATCAGCAAT AAACCAGCCA GCCGGAAGGG CCGAGCGCAG AAGTGGTCTT
4501 GCAACTTTAT CCGCTCCAT CCAGTCTATT AATTGTTGCC GGAAGCTAG
4551 AGTAAGTAGT TCGCCAGTTA ATAGTTTGGC CAACGTTGTT GCCATTGCTG
4601 CAGGCATCGT GGTGTCACGC TCGTCGTTTG GTATGGCTTC ATTCAGCTCC
4651 GGTTCCCAAC GATCAAGGCG AGTTACATGA TCCCCATGT TGTGCAAAAA
4701 AGCGGTTAGC TCCTTCGGTC CTCCGATCGT TGTCAGAAGT AAGTTGGCCG
4751 CAGTGTATC ACTCATGGTT ATGGCAGCAC TGCATAATC TCTTACTGTC
4801 ATGCCATCCG TAAGATGCTT TTCTGTGACT GGTGAGTACT CAACCAAGTC
4851 ATTCTGAGAA TAGTGTATGC GCGCAGCGAG TTGCTCTTGC CCGGCGTCAA
4901 TACGGGATAA TACCGCGCCA CATAGCAGAA CTTTAAAAGT GCTCATCATT
4951 GGAAAACGTT CTTCCGGGCG AAAACTCTCA AGGATCTTAC CGCTGTTGAG
5001 ATCCAGTTCG ATGTAACCCA CTCGTGCACC CAACTGATCT TCAGCATCTT
5051 TTA CTTT CAC CAGCGTTTCT GGGTGAGCAA AAACAGGAAG GCAAAATGCC
5101 GCAAAAAAGG GAATAAGGGC GACACGGAAG TGTTGAATAC TCATACTCTT
5151 CCTTTTTCAA TATTATTGAA GCATTTATCA GGGTTATTGT CTCATGAGCG
5201 GATACATATT TGAATGTATT TAGAAAAATA AACAAATAGG GGTCCGCGC
5251 ACATTTCCCG GAAAAGTGCC ACCTGAAATT GTAAACGTTA ATATTTTGTG
5301 AAAATTGCGG TTAATTTTTT GTTAAATCAG CTCATTTTTT AACCAATAGG
5351 CCGAAATCGG CAAAATCCCT TATAAATCAA AAGAATAGAC CGAGATAGGG
5401 TTGAGTGTG TTCCAGTTTG GAACAAGAGT CCACTATTAA AGAACGTGGA
5451 CTCCAACGTC AAAGGGCGAA AAACCGTCTA TCAGGGCGAT GGCCCACTAC
5501 GTGAACCATC ACCCTAATCA AGTTTTTTGG GGTGAGGTG CCGTAAAGCA
5551 CTAAATCGGA ACCCTAAAGG GAGCCCCGA TTTAGAGCTT GACGGGGAAA
5601 GCCGGCGAAC GTGGCGAGAA AGGAAGGGAA GAAAGCGAAA GGAGCGGGCG
```

pET Peptide Expression System 31



5651 CTAGGGCGCT GGCAAGTGTG GCGGTCACGC TCGCGTAAC CACCACACCC
 5701 GCCGCGCTTA ATGCGCCGCT ACAGGGCGCG TCCCATTCGC CA

pET-31b(+) sequence landmarks

T7 promoter*	616-632
T7 transcription start*	615
<i>lacI</i> coding sequence	1019-2098
KSI coding sequence*	170-544
<i>AlwNI</i> I site	169
His•Tag® coding seq.*	140-157
T7 terminator*	26-72
pBR322 origin	3532
<i>bla</i> coding sequence*	4287-5144
f1 origin*	5279-5734

* indicates sequence shown is reverse complement

pET-31b(+) Restriction Map

Enzyme	Sequence	# Sites	Locations
AccI	GT mk_AC	1	3240
AccII	C CGC	84	
AflIII	A CryG_T	2	1369 3470
AluI	AG CT	26	
AlwI	GGATCnnnn n_	14	
Alw2II	G_wGCw C	8	159 869 1353 2464 3288 3788 4943 5028
AlwNI	CAG_nnnn CTG	1	169
ApaI	G_GGCC C	1	1580
ApaBI	GCA_nnnnn TGC	1	1053
ApaLI	G TGCA_C	4	1349 3284 3784 5024
ApoI	r AATT_y	3	1644 5302 5313
AvaI	C yCGr_G	1	158
AvaII	G GwC_C	8	407 1921 2297 2385 2476 2755 4495 4717
BanI	G GyrC_C	10	
BanII	G_rGCy C	5	417 753 767 1580 5575
BbsI	GAAGACnn nnnn_	5	424 1515 1854 2228 2588
BbvI	GCAGC(n)8 (n)4n_	28	
BccI	CCATC	14	
Bce83I	CTTGAG(n)14_nn	7	21 2183 2353 3561 3859 4094 4962
Bcefl	ACGGC(n)12 n_	8	385 456 532 888 1229 1856 3966 5526
BcgI	CGA(n)6TGC(n)10_nn	8	1661 1695 2195 2229 3047 3081 4862 4896
BclI	T GATC_A	1	1383
BfaI	C TA_G	7	70 582 2484 3959 4212 4547 5651
BglI	GCCn_nnn nGGC	2	2433 4477
BglII	A GATC_T	1	647
BmgI	GkGCC	1	1578
BpmI	CTGGAGn(14)_nn	5	1207 1696 2330 2997 4427
Bpu10I	CC TnA_GC	1	2576
Bpu1102I	GC TnA_GC	1	80
BsaI	GGTCTCn nnnn_	1	4418
BsaAI	yAC GTr	2	3222 5500
BsaBI	GATnn nnATC	3	646 652 2667
BsaHI	Gr CG_yC	8	228 444 692 713 827 1326 2009 4894
BsaJI	C CnnG_G	10	
BsaWI	w CCGG_w	8	2 404 1688 2191 2659 3676 3823 4648
BsaXI	ACnnnnnCTCC	2	2028 5448
BsbI	CAACAC	2	3186 5407
BscGI	CCCGT	14	
BsiI	C TCGT_G	2	3643 5021
BsiEI	CG_ry CG	6	325 2154 3386 3810 4727 4876
BsII	CCnn_nnn nnGG	22	
BsmAI	GTCTCn nnnn_	7	1066 1471 1597 1984 3111 4418 5194
BsmBI	CGTCTCn nnnn_	2	1984 3111
BsmFI	GGGACn(10) nnnn_	4	830 2371 2741 5715
BsoFI	GC n_GC	48	
Bsp24I	GACn(6)TGGn(10)_nn	10	
Bsp1286I	G_dGCh C	13	
BspEI	T CCGG_A	2	2 2659
BspGI	CTGGAC	2	475 2996
BspLU11I	A CATG_T	1	3470
BsrI	ACTG_Gn	24	
BsrBI	GAG CGG	4	602 3403 5198 5644
BsrDI	GCAATG_nn	4	1416 1782 4418 4592
BsrFI	r CCGG_y	9	224 482 679 688 1055 2267 2427 4437 5601
BssHII	G CGCG_C	1	1780
Bst1107I	GTA TAC	1	3241
BstEII	G GTnAC_C	1	1550
BstXI	CCAn_nnnn nTGG	3	1171 1300 1423
BstYI	r GATC_y	11	
Cac8I	GCn nGC	40	



pET Peptide Expression System 31

Enzyme	Sequence	# Sites	Locations																				
CjeI	ACn(6)TGGn(9) n(6)_	28																					
CjePI	CCAn(7)TCn(7) n(6)_	20																					
Clai	AT CG_AT	1	249																				
CviJI	rG Cy	89																					
CviRI	TG CA	25																					
DdeI	C TnA_G	11																					
DpnI	GA TC	28																					
DraI	TTT AAA	3	4223	4242	4934																		
DraIII	CAC_ennn GTG	1	5500																				
DrdI	GACnn_nn nnGTC	4	470	3163	3578	5455																	
DrdII	GAACCA	2	1092	5505																			
DsaI	C CryG_G	5	424	439	514	806	2442																
EaeI	y GGCC_r	4	677	809	2043	4745																	
Eam1105I	GACnn_n nnGTC	1	4357																				
EorI	CTCTTCn nnn_	3	987	3354	5152																		
EciI	TCCGCC	4	1146	3544	3690	4512																	
Eco4711I	AGC GCT	4	507	774	2275	2724																	
Eco57I	CTGAAGn(14)_nn	2	4012	5024																			
EcoNI	CCTnn n_nnAGG	1	904																				
EcoO109I	rG GnC_Cy	3	53	802	2476																		
EcoRII	=CCwGG_	8	409	1092	1407	1947	2004	3496	3617	3620													
EcoRV	GAT ATC	1	1819																				
FauI	CCCGCnnnn nn_	17																					
FokI	GGATGnnnnnnnnn nnnn_11	1																					
FspI	TGC GCA	2	2451	4579																			
GdiII	y GGCCG	4	677	809	2043	4745																	
HaeI	wGG CCw	5	1097	2418	3485	3496	3942																
HaeII	r_GCCG y	16																					
HaeIII	GG CC	25																					
HgaI	GACGCnnnnn nnnn_	14																					
HgiEII	ACCnnnnnGGT	2	967	4050																			
HhaI	G_CG C	52																					
Hin4I	GAbnnnnvTC	4	412	1268	4356	4430																	
HincII	GTy rAC	1	1875																				
HinfI	G AnT_C	15																					
HpaI	GTT AAC	1	1875																				
HphI	GGTGAnnnnnn_n	19																					
MaeII	A CG_T	15																					
MaeIII	=GTnAC_	18																					
MboII	GAAGAnnnnnn_n	16																					
MluI	A CGCG_T	1	1369																				
MmeI	TCCrACn(18)_nn	3	3685	3869	5477																		
MnI	CCTCnnnnnn_n	27																					
MseI	T TA_A	28																					
MsII	CAynnnnrTG	10																					
MspI	C CG_G	34																					
MspAII	CmG CkG	10																					
MwoI	GCnn_nnn nnGC	46																					
NarI	GG CG_CC	5	228	692	713	827	2009																
NciI	CC s_GG	12																					
NdeI	CA TA_TG	1	543																				
NgoAIV	G CCGG_C	6	224	482	679	2267	2427	5601															
NlaIII	_CATG	25																					
NlaIV	GGn nCC	26																					
NsiI	A_TGCA T	1	543																				
NspI	r_CATG y	5	177	844	2815	3107	3474																
Pfi1108I	TCGTAG	1	951																				
PfIMI	CCAn_nnn nTGG	2	395	569																			
PleI	GAGTCnnnn n_	9	630	918	1005	1801	3364	3849	4346	5435	5443												
PshAI	GACnn nnGTC	1	2214																				
Psp5II	rG GwC_Cy	1	2476																				
Psp1406I	AA CG_TT	6	1031	2399	2795	4583	4956	5285															
PstI	C_TGCA G	1	4602																				
PvuI	CG_AT CG	1	4727																				
PvuII	CAG CTG	4	1969	2062	3061	3880																	
RcaI	T CATG_A	3	767	4184	5192																		
RsaI	GT AC	6	332	403	512	1516	3276	4837															
SapI	GCTCTTCn nnn_	1	3354																				
Sau96I	G GnC_C	20																					
Sau3AI	=GATC_	28																					
ScaI	AGT ACT	1	4837																				
ScrFI	CC n_GG	20																					
SfaNI	GCATCnnnnn nnnn_	23																					
SfcI	C TryA_G	5	615	3735	3920	4598	5719																
SgrAI	Cr CCGG_yG	1	688																				
SphI	G_CATG C	2	177	844																			
SspI	AAT ATT	3	184	5161	5292																		
StyI	C CwG_G	1	57																				
TaqI	T CG_A	13																					
TaqII	GACCGAnnnnnnnnn_n	9	1277	1495	2168	3372	4705	4890	5043	5060	5404												
TfiI	G AwT_C	6	391	2048	2350	2520	3024	3445															
ThoI	CG CG	38																					
TseI	CG CG	28																					
Tsp45I	=GTsAC_	8	1550	2378	2909	3122	3217	4613	4824	5673													
Tsp509I	=AATT_	15																					
Tth1111I	GACn n_GTC	1	3215																				
Tth11111I	CAArCAnnnnnnnnn_n	8	212	470	1208	1901	2931	4054	4061	4093													
UbaJI	GCsGC	20																					
VspI	AT TA_AT	4	630	2054	2113	4529																	
XbaI	T CTAG_A	1	58																				
XcmI	CCAnnnn_n nnnnTGG	3	1225	1741	1759																		
XhoI	C TCGA_G	1	158																				
XmnI	GAAnn nnTTC	2	3028	4956																			



Enzymes that do not cut pET-31b(+):

AatII	AflIII	AgeI	AscI	AvrII	BaeI	BamHI	BseRI
BsmI	BspMI	BsrGI	Bsu36I	EagI	EcoRI	FseI	HindIII
KpnI	MunI	NcoI	NheI	NotI	NruI	NspV	PacI
PmeI	PmlI	RleAI	RsrII	SacI	SacII	SalI	SexAI
SfiI	SgfI	SmaI	SnaBI	SpeI	SrfI	Sse8387I	StuI
SunI	SwaI						