



Instruction Manual

Expressway™ *In Vitro* Protein Synthesis System

For high-yield, *in vitro* protein synthesis from
plasmid or linear templates

Catalog nos. K9600-01, K9600-02, K9605-01, K9605-02

Version D

November 9, 2010

25-0455

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

Table of Contents

Table of Contents	iii
Procedure for Experienced Users	v
Kit Contents and Storage	vii
Accessory Products.....	viii
Introduction	1
Overview.....	1
Methods	2
Generating the DNA Template.....	2
Performing the Protein Synthesis Reaction.....	6
Analyzing Samples.....	9
Troubleshooting	13
Appendix	15
Technical Service.....	16
Purchaser Notification	18
References	20

Procedure for Experienced Users

Introduction

This quick reference sheet is provided for experienced users of the Expressway™ *In Vitro* Protein Synthesis System. If you are a first time user of the Expressway™ System, we recommend following the detailed protocols in the manual.

Step	Action								
Generate the DNA Template	For optimal expression in the Expressway™ System, clone your gene of interest into a pEXP-DEST or other T7 expression vector or use PCR to generate a DNA template containing the elements discussed on page 3.								
Purify the DNA Template	Use your method of choice or a commercially available kit to purify your plasmid or linear DNA template. Resuspend purified DNA in 1X TE Buffer or water such that the final concentration is greater than 500 ng/μl. Take care to avoid contamination of your DNA template with ethanol, salt, or RNases.								
Perform the Protein Synthesis Reaction	<ol style="list-style-type: none"> Prepare the following IVPS Reaction Cocktail on ice for each reaction to be performed. Remember to include a positive control (supplied with the kit) in your experiment. <table style="margin-left: 20px;"> <tr> <td>DNase/RNase-free water</td> <td style="text-align: right;">4 μl</td> </tr> <tr> <td>2.5X IVPS <i>E. coli</i> Reaction Buffer</td> <td style="text-align: right;">20 μl</td> </tr> <tr> <td><u>T7 RNA polymerase</u></td> <td style="text-align: right;"><u>1 μl</u></td> </tr> <tr> <td>Total Volume</td> <td style="text-align: right;">25 μl</td> </tr> </table> <p>Note: To generate radiolabeled proteins, add 2 μl ³⁵S Methionine to the IVPS Reaction Cocktail and decrease the volume of DNase/RNase-free water to 2 μl.</p> For each reaction, thaw one 20 μl tube of IVPS <i>E. coli</i> Extract on ice and add the 25 μl IVPS Reaction Cocktail from Step 1. Add DNA (1 μg of plasmid DNA or 2-3 μg of linear DNA) to the reaction tube. Bring the final reaction volume to 50 μl with DNase/RNase-free water. Vortex for 3 seconds to mix. Place tubes in a microcentrifuge and briefly centrifuge. Shake tubes at 1,400 rpm in a thermomixer at 37°C for 2 hours. Alternative: Transfer the 50 μl reaction to a 2 ml screw-cap tube (supplied with the kit) and incubate at 37°C for 2 hours in a standard shaking incubator (275-325 rpm) or in a water bath. Add 5 μl of RNase A to the 50 μl reaction and vortex briefly. Place tubes in a microcentrifuge and briefly centrifuge. Shake tubes at 1,400 rpm in a thermomixer at 37°C for 15 minutes. Alternative: Incubate tubes at 37°C for 15 minutes in a standard shaking incubator (275-325 rpm) or in a water bath. Place tubes in a microcentrifuge and briefly centrifuge. Place the reaction on ice and proceed to analyze samples. 	DNase/RNase-free water	4 μl	2.5X IVPS <i>E. coli</i> Reaction Buffer	20 μl	<u>T7 RNA polymerase</u>	<u>1 μl</u>	Total Volume	25 μl
DNase/RNase-free water	4 μl								
2.5X IVPS <i>E. coli</i> Reaction Buffer	20 μl								
<u>T7 RNA polymerase</u>	<u>1 μl</u>								
Total Volume	25 μl								
Analyze Sample	Use any method of choice including Coomassie® blue staining, Western blot, or activity assay to analyze your sample. If you plan to use Coomassie® blue staining or Western blot to analyze your sample, precipitate the proteins with acetone prior to performing polyacrylamide gel electrophoresis (see page 9).								

Kit Contents and Storage

Types of Kits

This manual is supplied with the following products.

Product	Amount	Catalog no.
Expressway™ <i>In Vitro</i> Protein Synthesis System <i>with pEXP1-DEST</i>	20 reactions	K9600-01
	8 reactions	K9605-01
<i>with pEXP2-DEST</i>	20 reactions	K9600-02
	8 reactions	K9605-02

Shipping/Storage

The Expressway™ *In Vitro* Protein Synthesis System is shipped on dry ice. Upon receipt, store the kit at -80°C.

Kit Contents

The following reagents are supplied in the Expressway™ *In Vitro* Protein Synthesis System. **Store at -80°C.**

Item	Composition	Amount/ 8 Rxn Kit	Amount/ 20 Rxn Kit
IVPS <i>E. coli</i> Extract	Proprietary	8 x 20 µl	20 x 20 µl
2.5X IVPS <i>E. coli</i> Reaction Buffer	Proprietary	400 µl	
DNase/RNase-Free Distilled Water	---	400 µl	
RNase A	1 mg/ml in DNase/RNase-free water	100 µl	
T7 RNA Polymerase	Proprietary	20 µl	
2 ml screw-cap tubes	--	8	20

Product Qualification

Each lot of the Expressway™ *In Vitro* Protein Synthesis System is functionally tested for protein generation by incorporation of ³⁵S-methionine. Using a control plasmid in a 50 µl Expressway™ protein synthesis reaction, greater than 18 µg recombinant protein must be obtained.

Accessory Products

Accessory Products

DNase/RNase-free water and other reagents suitable for use with the Expressway™ *In Vitro* Protein Synthesis System are available separately from Invitrogen. Ordering information is provided below.

Item	Quantity	Catalog no.
DNase/RNase-Free Distilled Water	500 ml	10977-015
S.N.A.P.™ MiniPrep Kit	25 reactions	K1900-01
S.N.A.P.™ MidiPrep Kit	20 reactions	K1910-01
Coomassie Brilliant Blue R®-250 Protein Stain	10 g	15528-011

Coomassie Brilliant Blue R® is a registered trademark of Imperial Chemical Industries PLC

Introduction

Overview

Introduction

The Expressway™ *In Vitro* Protein Synthesis System (“Expressway™ System”) is designed for T7-based, *in vitro* transcription and translation of target DNA to protein in a single tube. The Expressway™ System uses extract from an *E. coli* strain to drive the reaction, and a transcription/translation mix that allows for optimal protein production (up to 50 µg protein per 50 µl reaction, depending on the protein of interest) in about 2 hours. For higher yields, the reaction can be extended up to 4 hours.

Components of the System

The major components of the Expressway™ System include:

- Single-use tubes of S30 *E. coli* extract (Zubay, 1973)
 - T7 RNA polymerase (Studier *et al.*, 1990)
 - A reaction buffer composed of all the required amino acids and an ATP regenerating system (Kim *et al.*, 1996; Lesley *et al.*, 1991; Pratt, 1984)
-

Applications

The Expressway™ System is suitable for use in the following applications:

- Characterizing proteins
 - Analyzing mutants
 - Verifying cloned gene products
 - Producing analytical quantities of protein or radiolabeled protein
 - Producing proteins which are toxic to cells
-

Experimental Outline

The table below describes the major steps necessary to synthesize your recombinant protein of interest using the Expressway™ *In Vitro* Protein Synthesis System. Refer to the specified pages for details to perform each step.

Step	Action	Pages
1	Generate the DNA template.	2-4
2	Purify your DNA template.	5
3	Perform the protein synthesis reaction.	6-8
4	Analyze your sample using polyacrylamide gel electrophoresis, Western blot analysis, or activity assay.	9-12

Methods

Generating the DNA Template

Introduction

Successful use of the Expressway™ *In Vitro* Protein Synthesis System requires only the addition of a DNA template containing the gene of interest placed within the proper context of transcription/translation regulatory elements including a bacteriophage T7 RNA polymerase promoter (“T7 promoter”), prokaryotic Shine-Dalgarno ribosome binding site (RBS), ATG initiation codon, stop codon, and T7 terminator. However, protein yield can be significantly enhanced if the DNA template is optimally configured. The Expressway™ System includes an optimally configured pEXP-DEST vector from which to express your gene of interest. If you wish to use a different T7 expression vector, guidelines are provided to produce your DNA template in an optimal configuration for protein expression in the Expressway™ System.

Factors Affecting Protein Yield

The yield of protein produced in *in vitro* transcription and translation systems is generally dependent on many factors, including:

- The size of the protein
- The sequence of the gene of interest
- The spacing of the T7 promoter and the gene of interest in the DNA template
- The quality of the DNA template
- Stability of mRNA

Recommendations and guidelines to generate a DNA template with the optimal configuration and to purify the DNA template are provided in this section. The size of the protein and its sequence will vary depending on your gene of interest. Any variability in protein yield due to these two factors will require empiric experimentation to optimize expression conditions.

DNA Templates

The following DNA templates may be used in the Expressway™ System.

- Supercoiled plasmid DNA (recommended for most applications for highest protein yields)
- Linear DNA
- PCR product

For proper expression in the Expressway™ System, all templates **must** contain the T7 promoter, an initiation codon, and a prokaryotic Shine-Dalgarno ribosome binding site (RBS) upstream of the gene of interest. See the next page for a discussion of template optimization.

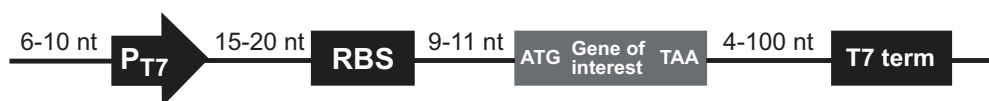
continued on next page

Generating the DNA Template, continued

Optimal Configuration of DNA Template

Optimized expression constructs are provided with the Expressway™ System, however, other expression constructs may be used. If you are designing your own expression construct, we recommend generating a DNA template that contains the following elements (see the figure below for reference). See the next page for an example.

- Gene of interest placed downstream of a T7 promoter and a ribosome binding site (RBS). The gene of interest **must** contain an ATG initiation codon and a stop codon.
- Sequence upstream of the T7 promoter containing a minimum of 6-10 nucleotides (nt) for efficient promoter binding (**required** for linear PCR products). This sequence need not be specific.
- Sequence following the T7 promoter containing a **minimum** of 15-20 nt which forms a potential stem-and-loop structure as described by Studier *et al.*, 1990 (see **T7 Expression Vectors**, below for more information).
- Sequence of 9-11 nt between the RBS and the ATG initiation codon for optimal translation efficiency of the protein of interest. This sequence need not be specific.
- A T7 terminator located 4-100 nt downstream of the gene of interest for efficient transcription termination and message stability. For the sequence of the T7 terminator, see page 4.



T7 Expression Vectors

Many T7-based expression vectors contain a T7 promoter, RBS, and T7 terminator with the suitable spacing and sequence configuration for optimal expression of protein in the Expressway™ System. In general, T7-based expression vectors that contain the bacteriophage ϕ 10-s10 segment (*i.e.* ϕ 10 promoter and the translation initiation region for the gene 10 protein) are recommended for use (see Studier *et al.*, 1990 for a list). The ϕ 10-s10 segment contains a region that forms a hypothetical stem-and-loop structure as described by Studier *et al.*, 1990. Examples of T7-based expression vectors containing this configuration include:

- pEXP1-DEST or pEXP2-DEST (supplied with the kit, as specified)
- pCR®T7/CT-TOPO® (Invitrogen, Catalog no. K4210-01)
- pCR®T7/NT-TOPO® (Invitrogen, Catalog no. K4200-01)
- pRSET (Invitrogen, Catalog no. V351-20)

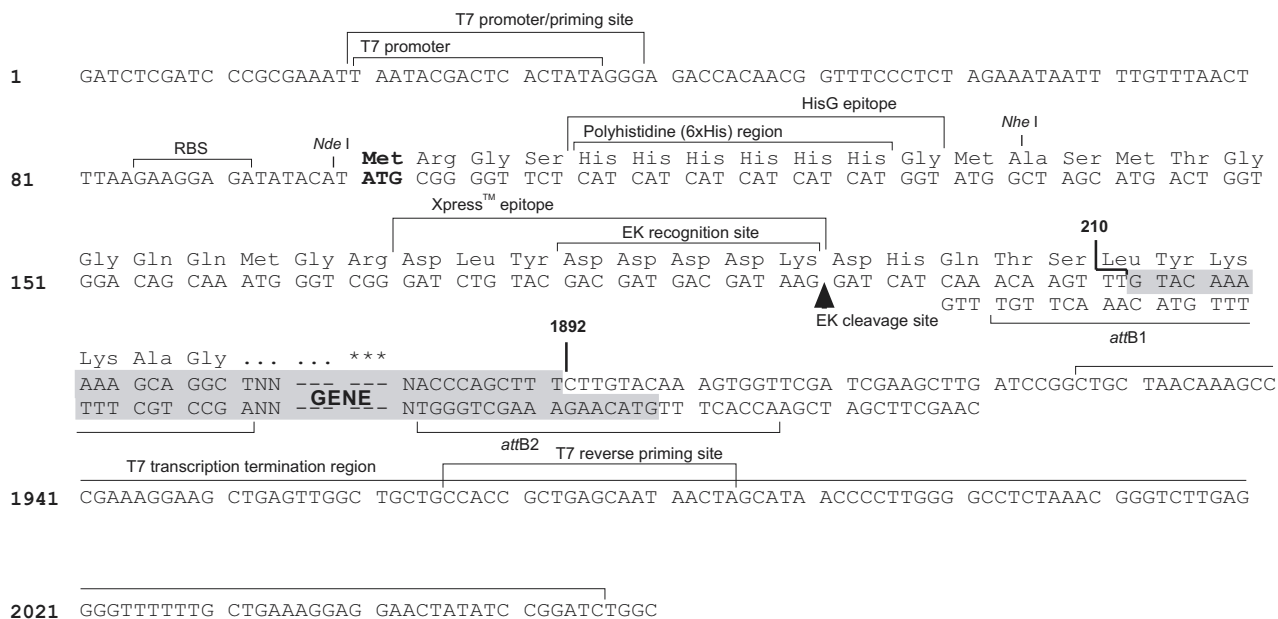
Note: T7 expression vectors that contain the T7_{lac} promoter (*e.g.* pET101/D-TOPO® available from Invitrogen™) are suitable for use in the Expressway™ System although protein yield may be reduced. In these vectors, the *lac* operator replaces the region containing the hypothetical stem-and-loop structure.

continued on next page

Generating the DNA Template, continued

Example

The T7-based pEXP1-DEST or pEXP2-DEST vector supplied with the kit allows you to generate an expression clone containing your gene of interest in an optimal configuration for expression in the Expressway™ System. To illustrate the points discussed on the previous page, the region surrounding the gene of interest after generating an expression clone in pEXP1-DEST is diagrammed below.



Generating an Expression Clone

You may generate an expression clone by recombination of an entry clone containing your gene of interest with the pEXP1-DEST or pEXP2-DEST vector using the Gateway® Technology. For more information about the pEXP1-DEST and pEXP2-DEST vectors, refer to the pEXP1-DEST and pEXP2-DEST Gateway® Vector Kits manual supplied with the vectors. For more information about the Gateway® Technology, creating an entry clone, and performing the LR recombination reaction, refer to the Gateway® Technology manual, which is available for downloading from our Web site (www.invitrogen.com) or by contacting Technical Service (see page 16).

continued on next page

Generating the DNA Template, continued

Purifying the DNA Template

Once you have generated the DNA template, you must purify the DNA before proceeding to the protein synthesis reaction. Depending on the type of DNA template (*i.e.* supercoiled plasmid, linear DNA, or PCR product), you may use any method of choice to purify your DNA template including commercial DNA purification kits (see below) or CsCl gradient centrifugation. For protocols to purify DNA, refer to published reference sources (Ausubel *et al.*, 1994; Sambrook *et al.*, 1989). When purifying your DNA template, keep the following in mind:

- **Do not** gel-purify your DNA template. Purified DNA solution obtained from agarose gels significantly inhibits the protein synthesis reaction. For PCR products or linear DNA, we recommend using a suitable purification kit.
- Ammonium acetate is not recommended for use in DNA precipitation as any residual contamination may inhibit translation. Use sodium acetate.
- Purified DNA **must** be free of RNases (wear gloves and use RNase-free reagents when preparing DNA).
- Purified DNA should be free of excess ethanol or salt as both can inhibit translation.

Note: Ethanol precipitated DNA should be carefully washed with 70% ethanol to remove excess salt and dried.

- Purified DNA should be resuspended in 1X TE Buffer or water such that the final concentration is at a minimum of 500 ng/ μ l.



Reminder: Do not gel-purify your DNA template. Purified DNA solution obtained from agarose gels significantly inhibits the protein synthesis reaction. For PCR products or linear DNA, we recommend using a suitable purification kit.



For rapid, resin-based isolation of purified plasmid DNA, we recommend using the S.N.A.P.[™] MiniPrep Kit (Catalog no. K1900-01) available from Invitrogen. Other commercial DNA purification kits are suitable.

Performing the Protein Synthesis Reaction

Introduction

Once you have obtained purified template DNA, you are ready to synthesize your recombinant protein using the Expressway™ System. General guidelines and instructions to produce your recombinant protein are provided in this section.



Important

RNase contamination may affect protein yield. To reduce the chances of RNase contamination, wear gloves and use RNase-free reagents (*i.e.* microcentrifuge tubes and pipette tips) when performing the protein synthesis reaction.

Incubation Conditions

To obtain the optimal yield of protein, we recommend using an Eppendorf Thermomixer (Fisher, Catalog no. 05-400-200) to shake your sample(s) at 37°C during the protein synthesis reaction (see Steps 8-11 in the protocol on page 8).

If a thermomixer is unavailable, you may use one of the following:

- Standard shaking incubator
- Standard shaking water bath
- Non-shaking water bath

We do not recommend using a non-shaking incubator because it produces a less stable and less consistent temperature environment.

2 ml Tubes

If you will not using a thermomixer, we recommend that you transfer the reaction mixture to a 2 ml screw-cap tube (supplied with the kit) before incubating your samples. Using one of the thermomixer alternatives listed above and incubating your samples in 2 ml tubes will result in protein yields that are comparable to that obtained using a thermomixer.

Positive Control

The pEXP1-GW/*lacZ* or pEXP2-GW/*lacZ* vector is provided in the kit as appropriate for use as a positive control for protein expression. pEXP1-GW/*lacZ* and pEXP2-GW/*lacZ* were constructed using the Gateway® LR recombination reaction between an entry clone containing the *lacZ* gene and pEXP1-DEST or pEXP2-DEST, respectively. For a map and a description of the features of each vector, see the pEXP1-DEST and pEXP2-DEST Gateway® Vector Kits manual.

continued on next page

Performing the Protein Synthesis Reaction, continued

Materials to Have on Hand

You should have the following materials on hand before beginning:

- DNA Template (purified; resuspended in TE or water at a concentration greater than 500 ng/ μ l)
 - IVPS *E. coli* Extract (supplied with the kit; keep on ice)
 - 2.5X IVPS *E. coli* Reaction Buffer (supplied with the kit; thaw on ice)
 - T7 RNA polymerase (supplied with the kit; keep on ice)
 - RNase A (supplied with the kit; do not thaw until needed)
 - DNase/RNase-free water (supplied with the kit)
 - ³⁵S Methionine (3,000 Ci/mmol; 15 μ Ci/ μ l; if needed)
 - Microcentrifuge tubes and microcentrifuge
 - Thermomixer (recommended), standard shaking incubator (set to 37°C), or water bath (set to 37°C)
 - 2 ml screw-cap tubes, optional (supplied with the kit)
 - Vortex mixer
-



Note

Upon thawing the 2.5X IVPS *E. coli* Reaction Buffer, you may notice some precipitate in the bottom of the tube. Gently flick the tube several times with your finger to mix and allow the precipitate to go back into solution. **Do not** pipette the buffer up and down or place the tube in a warm (greater than 37°C) water bath as this may result in loss of activity of the Reaction Buffer. Note that the solution may remain cloudy, but is suitable for use.



Important

The 2.5X IVPS *E. coli* Reaction Buffer is supplied in a single tube. Depending on the number of protein synthesis reactions performed, you may not use all of the 2.5X IVPS *E. coli* Reaction Buffer in a single experiment. In this case, make sure that you thaw the 2.5X IVPS *E. coli* Reaction Buffer on ice, mix (see **Note** above), remove the amount needed, and return the reaction buffer to a -80°C freezer. **Do not** store the 2.5X IVPS *E. coli* Reaction Buffer at -20°C or room temperature as this may result in loss of activity. To prevent contamination, use RNase-free, sterile pipette tips and wear gloves when removing the 2.5X IVPS *E. coli* Reaction Buffer from the tube.

continued on next page

Performing the Protein Synthesis Reaction, continued

Performing the Protein Synthesis Reaction

Use the protocol below to synthesize your protein from the DNA template.

1. For each sample, prepare the following IVPS Reaction Cocktail on ice. If you are performing multiple reactions, scale up the volume of each reagent accordingly.

DNase/RNase-free water 4 μ l

2.5X IVPS *E. coli* Reaction Buffer 20 μ l

T7 RNA polymerase 1 μ l

Total Volume 25 μ l

Note: If you wish to produce radiolabeled proteins, add 2 μ l 35 S Methionine to the IVPS Reaction Cocktail and decrease the volume of DNase/RNase-free water to 2 μ l. Include a negative control (no DNA) to determine background.

2. For each reaction to be performed, thaw one tube of IVPS *E. coli* Extract on ice.
3. To each tube of extract, add the 25 μ l IVPS Reaction Cocktail from Step 1.
4. Add the following amount of DNA to the reaction tube.
 - For plasmid DNA templates, add 1 μ g
 - For linear DNA templates or PCR products, add 2-3 μ g
 - For control DNA, add 1 μ g.
5. Bring the final reaction volume to 50 μ l with DNase/RNase-free water. For no-DNA controls, add 5 μ l of DNase/RNase-free water only.
6. Vortex for 3 seconds to mix.
7. Place tubes in a microcentrifuge and briefly centrifuge.
8. Shake tubes at 1,400 rpm in a thermomixer at 37°C for 2 hours.

Alternative: Transfer the 50 μ l reaction to a 2 ml screw-cap tube (supplied with the kit) and incubate at 37°C for 2 hours in a standard shaking incubator (275-325 rpm) or in a water bath.

Note: It is possible to shake tubes for up to 4 hours to obtain as much as 50% greater protein yield.

9. Add 5 μ l of RNase A to the 50 μ l reaction and vortex briefly.
 10. Place tubes in a microcentrifuge and briefly centrifuge.
 11. Shake tubes at 1,400 rpm in a thermomixer at 37°C for 15 minutes.

Alternative: Incubate tubes at 37°C for 15 minutes in a standard shaking incubator (275-325 rpm) or in a water bath.
 12. Place tubes in a microcentrifuge and briefly centrifuge. Place the reaction on ice and proceed to **Analyzing Samples**, next page.
-

Analyzing Samples

Introduction

Once you have performed the protein synthesis reaction, you may use any method of choice to analyze your sample. Generally, sufficient protein is produced for analysis on a Coomassie[®]-stained protein gel, by Western blot analysis, by enzymatic activity, or by affinity purification (if affinity tag is present), however, expression levels may vary depending on the nature of your protein and the configuration of the DNA template (see page 3 for more details). If you plan to analyze your sample using polyacrylamide gel electrophoresis, note that you should first precipitate the proteins with acetone to remove background smearing. A protocol for acetone precipitation and other general guidelines for gel electrophoresis are provided in this section.

If you have included radiolabeled methionine in the protein synthesis reaction, you may use TCA precipitation to determine the amount of radiolabeled methionine incorporated and to calculate the yield of protein (see pages 11-12).

Materials Needed

You should have the following materials on hand before proceeding:

- Acetone (room temperature)
 - Microcentrifuge
 - SpeedVac[®] concentrator (Thermo Savant)
 - 1X SDS-PAGE sample buffer (see page 15 for a recipe)
 - Appropriate polyacrylamide gel to resolve your protein of interest (see the next page)
 - Coomassie[®] blue stain
-

Acetone Precipitation

Before starting, prepare an SDS-PAGE gel or use one of the pre-cast polyacrylamide gels available from Invitrogen[™] (see the next page) to analyze your samples. Use the following protocol to precipitate your proteins prior to loading on the polyacrylamide gel.

1. Add 5 μ l of the protein reaction product from Step 12, page 8 to 20 μ l of acetone. Mix well.
 2. Centrifuge for 5 minutes at room temperature in a microcentrifuge at 12,000 rpm.
 3. Carefully remove the supernatant, taking care not to disturb the protein pellet.
 4. Dry the sample under vacuum for 15 minutes. **Note:** Alternatively, you may air dry the sample for 1 hour at room temperature.
 5. Resuspend pellet in 20 μ l of 1X SDS-PAGE sample buffer.
 6. Heat at 70-80°C for 10-15 minutes and centrifuge briefly. Proceed to **Polyacrylamide Gel Electrophoresis**, next page. **Note:** Alternatively, samples may be stored at -20°C until needed.
-

[®]SpeedVac is a registered trademark of Thermo Savant

continued on next page

Analyzing Samples, continued

Polyacrylamide Gels Available from Invitrogen™

To facilitate separation and visualization of your recombinant protein by polyacrylamide gel electrophoresis, a wide range of pre-cast NuPAGE® and Novex® Tris-Glycine polyacrylamide gels and electrophoresis apparatus are available from Invitrogen. In addition, Invitrogen carries a large selection of molecular weight protein standards and staining kits. For more information about the appropriate gels, standards, and stains to use to visualize your recombinant protein, refer to our Web site (www.invitrogen.com) or call Technical Service (see page 16).

Polyacrylamide Gel Electrophoresis

1. Load 5-10 µl of the sample from Step 6, previous page on an SDS-PAGE gel and electrophorese at 120V. You may save your sample by storing at -20°C, if desired.
2. Depending on your assay of choice, perform the following:

If you are...	Then...
Visualizing your protein using Coomassie® blue staining	Stain gel with Coomassie® blue stain. Note: For radiolabeled proteins, the signal may be enhanced by placing the gel in a commercially available reagent that enhances the signal. Dry the gel and expose to x-ray film for 1-4 hours.
Analyzing your protein by Western blot	Transfer proteins electrophoretically to a suitable membrane and use an appropriate antibody to detect the protein of choice.

continued on next page

Analyzing Samples, continued

Determining Total Counts Incorporated

For radiolabeled reactions, we recommend using TCA precipitation to determine the amount of ^{35}S methionine incorporated. A protocol is provided below.

Determining Total Counts

1. Mix and spot 5 μl of each radiolabeled reaction from Step 12, page 8 on an individual glass microfiber filter (Type GF/C; Whatman, Catalog no. 1822-021).
2. Set aside and let dry. **Do not** wash or TCA precipitate these filters.

Performing TCA Precipitation

1. Mix and spot 5 μl of each radiolabeled reaction from Step 12, page 8 on a separate set of individual glass fiber (GF/C) filters and allow to air dry for approximately 5-10 seconds.
 2. Place filter in a beaker and wash once with cold 10% TCA/1% sodium pyrophosphate for 10 minutes at room temperature while shaking gently (use approximately 10-20 ml per filter).
 3. Wash with 5% TCA for 5 minutes at room temperature while shaking gently. Repeat wash.
 4. Rinse filters with methanol to facilitate drying.
 5. Allow filters to dry, place in scintillation vials, and add scintillation fluid.
 6. Count samples in a scintillation counter.
 7. Proceed to **Calculating the Yield of Protein**, below to determine your protein yield.
-

Calculating the Yield of Protein

If you generate radiolabeled proteins using 2 μl of ^{35}S methionine (3,000 Ci/mmol, 15 $\mu\text{Ci}/\mu\text{l}$) in the protein synthesis reaction (see page 8), you may calculate the yield of protein as shown below. For a sample calculation, see the next page.

Before beginning, you will need to determine the total counts incorporated using TCA precipitation (see above).

Total pmoles of methionine (includes radiolabeled and non-radiolabeled methionine) = 1,240 pmoles/ μl (or 62,000 total pmoles for a 50 μl reaction)	
Total counts:	Total cpm per 5 μl spotted x total reaction volume/5
Specific activity:	$\frac{\text{Total counts}}{\text{pmoles of methionine}}$
pmoles methionine incorporated:	$\frac{(\text{TCA precipitable counts} - \text{background}) \times 50/5}{\text{specific activity}}$
pmoles of protein:	$\frac{\text{pmoles of methionine incorporated for protein}}{\text{number of methionines in protein}}$
Yield of protein:	pmoles of protein x molecular weight of protein

continued on next page

Analyzing Samples, continued

Sample Calculation for Protein Yield

The sample calculation below illustrates how to use the data obtained from the TCA precipitation procedure, page 11 to calculate the yield of protein.

Example: We wish to calculate the yield of chloramphenicol acetyltransferase (CAT) protein obtained from a typical protein synthesis reaction using a T7-based plasmid containing the CAT gene.

1. We know the following:
 - The CAT protein has a molecular weight of 24 kDa and contains 9 methionines (0.024 $\mu\text{g}/\text{pmole}$)
 - There are 62,000 total pmoles of methionine in a 50 μl reaction
2. Using the TCA precipitation protocol on page 11, we obtain the following values after scintillation counting:

Sample	Radioactive counts (cpm)
Total counts (5 μl spotted):	4×10^6 cpm
Background TCA precipitable counts (No DNA; 5 μl spotted):	1.3×10^4 cpm
CAT TCA precipitable counts (5 μl spotted):	6.2×10^5 cpm

3. Determine total counts in the reaction:

$$4 \times 10^6 \text{ cpm in } 5 \mu\text{l} \times \frac{50}{5} = 4 \times 10^7 \text{ cpm in } 50 \mu\text{l reaction}$$

4. Calculate the specific activity:

$$\frac{4 \times 10^7 \text{ cpm}}{62000} = 661 \text{ cpm/pmoles methionine}$$

5. Calculate the pmoles methionine incorporated:

$$\frac{\left[(6.2 \times 10^5 \text{ cpm} - 1.3 \times 10^4 \text{ cpm}) \times \frac{50}{5} \right]}{661} = 9183 \text{ pmoles methionine}$$

6. Calculate the pmoles of protein obtained:

$$\frac{9183 \text{ pmoles}}{9} = 1020 \text{ pmoles protein}$$

7. Calculate the yield of protein obtained:

$$1020 \text{ pmoles} \times 0.024 \mu\text{g}/\text{pmole} = 24 \mu\text{g protein}$$

Troubleshooting

Introduction

The table below and on the next page lists some potential problems and possible solutions that you may use to help you troubleshoot your *in vitro* transcription and translation experiment.

Problem	Reason	Solution
Low or no yield of target protein (but control reaction produces protein)	DNA template not optimally configured	<ul style="list-style-type: none"> Use one of the pEXP-DEST vectors or follow the guidelines on pages 2-3 to clone your gene of interest into a T7 expression vector with the optimal configuration. Make sure that the ATG initiation codon is in the proper context for expression (<i>i.e.</i> check spacing and placement after the RBS). For linear DNA templates, make sure that at least 6-10 additional nucleotides are present upstream of the T7 promoter.
	DNA template not pure <ul style="list-style-type: none"> Contaminated with ethanol, sodium salt, or ammonium acetate Contaminated with RNases 	<ul style="list-style-type: none"> Prepare new DNA template taking care to remove excess ethanol and/or salt after precipitation. Do not use ammonium acetate to precipitate DNA. Use sodium acetate. Wear gloves and use RNase-free reagents when preparing DNA.
	DNA template purified from agarose gel	We recommend using a commercial DNA purification kit to purify your DNA template. See page 5 for additional guidelines.
	DNA concentration not optimal (generally more of a problem for linear DNA or PCR products)	<ul style="list-style-type: none"> Determine the concentration of your DNA template and adjust the concentration to at least 500 ng/μl. Titrate the amount of DNA template used in the protein synthesis reaction to determine the optimal yield.
	Sample incubated in a non-shaking incubator during protein synthesis reaction	<ul style="list-style-type: none"> Use a thermomixer and shake sample at 1,400 rpm (see protocol on page 8). Incubate samples in a standard shaking incubator (275-325 rpm) or in a water bath (see recommend alternatives on page 6).
	Expression time too short	Extend expression time up to 4 hours.

continued on next page

Troubleshooting, continued

Problem	Reason	Solution
	Size of protein	Protein yield may decrease as the size of the protein increases; optimize expression conditions.
	Sample not mixed before spotting on filter for TCA precipitation (radiolabeled samples only)	Mix sample before spotting on filter for TCA precipitation.
Control reaction produces no protein	Reagents have lost activity	<ul style="list-style-type: none"> • Store reagents at -80°C. • Use care when freeze thawing the 2.5X IVPS <i>E. coli</i> Reaction Buffer.
	2.5X IVPS <i>E. coli</i> Reaction Buffer contaminated	Wear gloves and use RNase-free reagents when working with 2.5X IVPS <i>E. coli</i> Reaction Buffer to prevent contamination.
Multiple bands on the polyacrylamide gel	Proteins denatured for too long	Add 1X SDS-PAGE sample buffer to sample and heat at 70-80°C for 10-15 minutes before loading on gel.
	Old ³⁵ S methionine	Use fresh ³⁵ S methionine.
	Not enough SDS in the 1X SDS-PAGE sample buffer	Prepare new 1X SDS-PAGE sample buffer according to standard instructions.
	Internal ATG codons in the context of RBS-like sequences	<ul style="list-style-type: none"> • Check the sequence of your gene and search for potential RBSs with the proper spacing from internal methionines. • Replace the methionine or change RBS sequence(s) using point mutation(s).
Smearing on the gel	Samples not precipitated with acetone	Precipitate the proteins with acetone to remove background smearing. Follow the protocol provided on page 9.
	Too much protein loaded	Reduce the amount of protein loaded on the gel.
	Gel not clean	<ul style="list-style-type: none"> • Rinse the gel briefly before exposing to film. • If you have stained the gel with Coomassie® blue, destain the gel in water or 50% methanol, 7.5% glacial acetic acid for 15-30 minutes before drying. If you have already destained the gel, repeat destaining procedure.
	Ethanol present in the protein synthesis reaction	Make sure that any residual ethanol is removed during DNA purification.
	Old pre-cast gels	Do not use pre-cast gels after the expiration date.

Appendix

Recipes

1X SDS-PAGE Sample Buffer

1. Combine the following reagents:

0.5 M Tris-HCl, pH 6.8	2.5 ml
Glycerol (100%)	2 ml
β -mercaptoethanol	0.4 ml
Bromophenol Blue	0.02 g
SDS	0.4 g
 2. Bring the volume to 20 ml with sterile water.
 3. Aliquot and freeze at -20°C until needed.
-

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, 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

Japanese Headquarters:

Invitrogen Japan K.K.
Nihonbashi Hama-Cho Park Bldg. 4F
2-35-4, Hama-Cho, Nihonbashi
Tel: 81 3 3663 7972
Fax: 81 3 3663 8242
E-mail: jpinfo@invitrogen.com

European Headquarters:

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

MSDS Requests

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

continued on next page

Technical Service, continued

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

Introduction

Use of the Expressway™ *In Vitro* Protein Synthesis System is covered under the licenses detailed below.

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

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 Life Technologies 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 Life Technologies 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 Life Technologies 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. 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. Life Technologies 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. Life Technologies 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, Life Technologies 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, Life Technologies Corporation, 5791 Van Allen Way, Carlsbad, California 92008. Phone (760) 603-7200.

continued on next page

Purchaser Notification, continued

**Limited Use Label
License No.5:
Invitrogen Techno-
logy**

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 buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. The buyer may transfer information or materials made through the use 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. 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. For products that are subject to multiple limited use label licenses, the terms of the most restrictive limited use label license shall control. Life Technologies Corporation will not assert a claim against the buyer of infringement of patents owned or controlled by Life Technologies Corporation which cover this product 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 neither this product nor any of its components was used in the manufacture of such product. If the purchaser is not willing to accept the limitations of this limited use statement, Life Technologies is willing to accept return of the product with a full refund. For information about purchasing a license to use this product or the technology embedded in it for any use other than for research use please contact Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008; Phone (760) 603-7200 or e-mail: outlicensing@lifetech.com.

**Limited Use Label
License No. 22:
Vectors & Clones
Encoding
Histidine Hexamer**

This product is licensed under U.S. Patent Nos. 5,284,933 and 5,310,663 and foreign equivalents from Hoffmann-LaRoche, Inc., Nutley, NJ and/or Hoffmann-LaRoche Ltd., Basel, Switzerland and is provided only for use in research. Information about licenses for commercial use is available from QIAGEN GmbH, Max-Volmer-Str. 4, D-40724 Hilden, Germany.

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*.

Kim, D. M., Kigawa, T., Choi, C. Y., and Yokoyama, S. (1996). A Highly Efficient Cell-free Protein Synthesis System from *E. coli*. *Eur. J. Biochem.* 239, 881-886.

Lesley, S. A., Brow, M. A., and Burgess, R. R. (1991). Use of *in vitro* Protein Synthesis from Polymerase Chain Reaction-generated Templates to Study Interaction of *Escherichia coli* Transcription Factors with Core RNA Polymerase and for Epitope Mapping of Monoclonal Antibodies. *J. Biol. Chem.* 266, 2632-2638.

Pratt, J. M. (1984). *Transcription and Translation* (Oxford: S.J. IRL Press).

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

Studier, F. W., Rosenberg, A. H., Dunn, J. J., and Dubendorff, J. W. (1990). Use of T7 RNA Polymerase to Direct Expression of Cloned Genes. *Meth. Enzymol.* 185, 60-89.

Zubay, G. (1973). *In vitro* Synthesis of Protein in Microbial Systems. *Annu. Rev. Genet.* 7, 267-287.

©2001-2003, 2010 Invitrogen Corporation. All rights reserved.

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

Notes

Notes



United States 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 website

www.invitrogen.com

