



Instruction Manual

ProQuest™ Pre-made cDNA Libraries

For detecting protein-protein interactions

Version B
December 12, 2002
25-0617

Table of Contents

General Information	2
Overview	5
Using ProQuest™ Libraries.....	8
pPC86	13
pEXP-AD502.....	15
Recipe.....	17
Accessory Products	18
Purchaser Notification	19
Technical Service.....	22
References.....	24

General Information

Contents and Storage 2 x 0.5 ml aliquots
Each cDNA library is supplied in 80% SOB medium, 20% (v/v) glycerol.
Store the library at -80°C. The cDNA library is stable for six months when stored properly.

Titer of the Libraries Each library has greater than 5×10^9 cfu (colony forming units) derived from $> 10^7$ primary clones to ensure complete representation of rare sequences.

ProQuest™ Pre-made cDNA Libraries The following ProQuest™ Pre-made cDNA Libraries are available from Invitrogen. For more information on preparing the library, see page 7.

Product	Avg. Insert Size	Vector	Catalog no.
Human			
Brain (27 years, male)	1.9 kb	pEXP-AD502	11376-027
Brain (Fetal)	2.3 kb	pEXP-AD502	11386-026
Brain (Fetal)	2.1 kb	pPC86	11386-018
Heart (26 years, male)	1.4 kb	pPC86	11378-015
Heart (23 years, male)	1.7 kb	pEXP-AD502	11378-023
HeLa cells	1.5 kb	pPC86	11287-018
Liver (28 years, male)	2.0 kb	pPC86	11381-019
Skeletal muscle (24 years, male)	1.6 kb	pPC86	11367-018
Prostate (25 years, male)	1.3 kb	pPC86	11579-018
Prostate, poorly differentiated adenocarcinoma, (63 years, male)	1.3 kb	pPC86	11582-012
Prostate, moderately, differentiated leiomyosarcoma (29 years, male)	1.3 kb	pPC86	11583-010

Continued on next page

General Information, Continued

ProQuest™ Pre-made cDNA Libraries, continued

Product	Avg. Insert Size	Vector	Catalog no.
Mouse			
Brain	1.5 kb	pPC86	11298-015
Embryo (8.5 day)	1.2 kb	pPC86	11291-010
Embryo (10.5 day)	1.4 kb	pPC86	11292-018
Liver	1.5 kb	pPC86	11297-017
Lymph nodes (<i>in vivo</i> activated, see RNA Source , next page)	1.7 kb	pPC86	11289-014
Rat			
Brain	1.2 kb	pPC86	11290-012
Liver	1.3 kb	pPC86	11296-019
<i>Caenorhabditis elegans</i> (see RNA Source , next page)	1.4 kb	pPC86	11288-016

Continued on next page

General Information, Continued

RNA Source

Mouse Lymph Nodes

The RNA was obtained from mesenteric lymph node tissue at day 8 after oral inoculation of BALB/c mice with third stage larvae of the nematode, *Heligmosomoides polygyrus*. *H. polygyrus* triggers a strong Th2 mucosal response at day 8 after immunization (Greenwald *et al.*, 1997; Svetic *et al.*, 1993). The mesenteric lymph node at day 8 is composed of about 60% T cells, 30% B cells, 10% dendritic cells, macrophages, and other tissue-associated cells.

C. elegans

C. elegans (Bristol N2) RNA (supplied by Monique A Kreutzer and Sander van den Heuvel of the Massachusetts General Hospital Cancer Center, Charlestown, Massachusetts) isolated from populations of non-synchronously staged L1 to adult nematodes (including dauer larvae and embryos) was combined to generate approximately equal representation of RNA expressed at different developmental stages. The nematodes were mated to ensure representation of sex-specific RNAs.

Quality Control

The ProQuest™ Pre-made cDNA Libraries are qualified by PCR amplification of 23 randomly selected clones using primers flanking the multiple cloning site. Twenty out of twenty-three clones must contain inserts.

Overview

Introduction

ProQuest™ Pre-made cDNA Libraries are suitable for use with the ProQuest™ Two-Hybrid System (see page 11). Each library is constructed using the SuperScript™ II Plasmid System for cDNA Synthesis and Plasmid Cloning to generate full-length and high-yield cDNA.

Libraries are constructed in a standard vector pPC86 or the Gateway® vector, pEXP-AD502 (see next page). For information on library preparation, see page 7.

ProQuest™ Libraries

ProQuest™ Pre-Made cDNA Libraries may be screened for interacting proteins using the ProQuest™ Two-Hybrid System (see page 11), PCR or plate screening procedures, further amplified (Sambrook *et al.*, 1989), or used for the production of single-stranded plasmid DNA *in vivo* (Li *et al.*, 1994). For more information on the ProQuest™ Two-Hybrid System, refer to the manual. This manual is available for downloading from our Web site at www.invitrogen.com or by contacting Technical Service (see page 22).

Libraries prepared in the Gateway® vector pEXP-AD502 allow you to shuttle DNA sequences of interest into a variety of Gateway®-compatible expression and analysis vectors using the Gateway® recombinational cloning Technology (see page 12).

Continued on next page

Overview, Continued

pPC86 and pEXP-AD502

The vectors pPC86 and pEXP-AD502 are activation domain expression vectors. pEXP-AD502 is similar to pPC86 (Chevray and Nathans, 1992) except pEXP-AD502 contains the Gateway[®] recombination sites *attB1* and *attB2*. The major features of the two vectors are listed below.

- The constitutive, moderate-strength yeast alcohol dehydrogenase (*ADH1*) promoter for expression of *GAL4* fusions
- The SV40 large T antigen nuclear localization sequence (NLS)
- The *GAL4* activation domain (AD) allowing expression of the reporter gene which is activated when brought into proximity with the DNA binding domain by interacting bait and prey proteins
- The recombination sites, *attB1* and *attB2* (only in pEXP-AD502) for transfer of cDNA into Gateway[®]-compatible vectors
- The *ADH1* transcription termination (TT) for efficient transcription termination and stabilization of the mRNA
- An *f1* origin for ss DNA production
- The *TRP1* promoter for expression of the *TRP1* gene
- The *TRP1* gene for auxotrophic selection of the plasmid in Trp⁻ yeast hosts
- An *ARSH4/CEN6* sequence for replication and low-copy number maintenance of plasmid in yeast
- Ampicillin resistance gene for selection of transformants in *E. coli*
- The pUC origin for high copy replication and maintenance of the plasmid in *E. coli*

For maps of vectors, see pages 13-15.

Continued on next page

Overview, Continued

Preparing ProQuest™ Libraries

ProQuest™ Pre-made cDNA Libraries are prepared as follows:

- mRNA is isolated using two steps. First, total RNA is isolated from tissues or cells using the TRIzol® Reagent. Second, mRNA is isolated from total RNA using oligo (dT) in a filter syringe.
- First-strand cDNA is synthesized using SuperScript™ Plasmid System for cDNA Synthesis and Plasmid Cloning with *Not* I primer-adaptor (D'Alessio, J. M., Gruber, C. E., Cain, C., and Noon, M. C. 1990) *Focus*® 12: 47)
- Second-strand cDNA is synthesized using *E. coli* RNase H, *E. coli* DNA polymerase I, and *E. coli* DNA ligase
- Blunt-end cDNA is created using T4 DNA polymerase
- cDNA is adapted with *Sal* I adaptor and digested with *Not* I
- cDNA is size-selected using column chromatography
- Size-selected cDNA is directionally cloned downstream of the *GAL4* activation domain into the *Sal* I-*Not* I region of the vector (pPC86 or pEXP-AD502) to generate *GAL4* AD-cDNA fusion proteins
- Ligation mixture is transformed into competent ElectroMAX™ DH10B™ *E. coli* and the number of primary recombinants is determined
- cDNA library is amplified once using a semi-solid procedure (Kriegler, 1990) to minimize representational biases.

Genotype of DH10B™

F' *mcrA* Δ(*mrr-hsdRMS-mcrBC*)φ80*lacZ*ΔM15 Δ*lacX74* *deoR recA1 endA1 araD139* Δ(*ara,leu*)7697 *galU galK rpsL nupG*

Using ProQuest™ Libraries

Introduction

ProQuest™ Pre-made cDNA Libraries may be screened for interacting proteins using the ProQuest™ Two-Hybrid System (see page 11), PCR or plate screening procedures.

General procedures for preparing DNA from the library and colony PCR are provided in this section. For detailed information on library screening refer to published references (Ausubel *et al.*, 1994; Sambrook *et al.*, 1989).

Preparing dsDNA from a Plasmid cDNA Library

You will need the following items.

- Terrific Broth (see page 17 for a recipe)
 - Buffer I with RNase (15 mM Tris-HCl, pH 8.0; 10 mM EDTA; 100 µg/ml RNase A; 1200 U/ml RNase T1)
 - Buffer II (0.2 M NaOH, 1% SDS)
1. Inoculate 100 ml Terrific Broth containing 100 µg/ml ampicillin with 2.5×10^9 cells from the library in a 500-ml flask.
 2. Incubate the culture for 16 hours at 30 °C with shaking at 275 rpm. **Clones are prone to deletions when the culture is grown at > 30°C.**
 3. Read the A_{590} of the culture. For accurate A_{590} determination, dilute the cells 1:10-1:20, so that the observed value is between 0.2 and 0.8. In two 50-ml centrifuge tubes, process ~ 500 OD_{590} units.
 4. Centrifuge the tubes at 4800 x g for 15 minutes at 4°C. Discard the supernatant.
 5. Resuspend the cell pellets in a total volume of 10 ml Buffer I with RNase. The cells must be < 50 OD/ml.
 6. Add 10 ml of Buffer II to the resuspended cells. Invert the tubes to mix the cells and incubate for 5 minutes at room temperature.
-

Continued on next page

Using ProQuest™ Libraries, Continued

Preparing dsDNA from a Plasmid cDNA Library, continued

7. Add 10 ml cold 7.5 M ammonium acetate to the cell mixture. Invert the tubes to mix the cells and incubate for 10 minutes on ice.
8. Centrifuge the tubes at 3,000 x g for 15 minutes at 4°C. Pour the supernatant through cheesecloth or a clean, DNase-free, porous filter into a fresh 50-ml centrifuge tube. Avoid the white flocculant material.
9. Add an equal volume of cold isopropanol to the tube, mix well, and centrifuge the tubes at 3,000 x g for 15 minutes at 4°C. Discard the supernatant.
10. Resuspend the cell pellet in 1 ml of Buffer I with RNase and transfer it to a microcentrifuge tube.
11. Centrifuge the tubes at 14,000 x g for 1 minute at 4°C. Transfer the supernatant to a fresh microcentrifuge tube. Incubate the tube at 37°C for 10 minutes.
12. Incubate the tubes at 65 °C for 5 minutes. Split each sample into two equal parts (~ 500 µl each) in 1.5-ml microcentrifuge tubes.
13. Add an equal volume of phenol:chloroform:isoamyl alcohol (25:24:1) to each sample and vortex the tubes. Centrifuge the tubes at 14,000 x g for 5 minutes at room temperature.
14. Avoiding the interface, transfer 450 µl of the upper (aqueous) phase to a fresh microcentrifuge tube.
15. Repeat the phenol : chloroform : isoamyl alcohol extraction at least twice. If an interface remains, repeat steps 13 and 14 until the supernatant is clear.
16. Add an equal volume of isopropanol (4 °C) to each tube. Centrifuge the tubes at 14,000 x g for 15 minutes at 4°C. Discard the supernatant.
17. Add 500 µl 70 % ethanol to each tube. Centrifuge the tubes at 14,000 x g for 5 minutes at 4°C. Discard the supernatant.
18. Dry the pellet for 10 minutes at room temperature.
19. Completely dissolve the two pellets in 200 µl TE buffer. The plasmid DNA library concentration must be approximately 1 µg/µl. Store the DNA at -20°C.

Continued on next page

Using ProQuest™ Libraries, Continued

Colony PCR Screening

A colony PCR procedure to screen for the presence of specific cDNA is described below. You may also use this method to identify desired cDNA clones.

The recommended primers for PCR and their sequences are shown on page 14 for pPC86 and on page 16 for pEXP-AD502.

1. Add 10 μ l TE to each labeled, 0.5 ml microcentrifuge tube.
2. Pick individual colonies using a pipette tip and place the colonies directly into separate tubes containing TE. Pipette up and down to mix.
3. Incubate the tubes in a pre-warmed thermal cycler at 99°C for 5 minutes.
4. Incubate the tubes on ice for 2 minutes.
5. Centrifuge briefly to collect the sample at the bottom of the tube. Replace the tubes on ice.
6. Prepare the appropriate amount of the following reaction mix and add 40 μ l of reaction mix to each tube.
 - 1X PCR Buffer (contains no $MgCl_2$)
 - 0.2 mM dNTP mix (see page 18)
 - 0.5 μ M primers
 - 2.4 mM $MgCl_2$
 - 2.5 units Platinum® *Taq* DNA polymerase (see page 18)
7. Bring the volume to 50 μ l with sterile water.

Continued on next page

Using ProQuest™ Libraries, Continued

Colony PCR Screening, continued

8. Perform PCR using the following cycling parameters:

Temperature	Time	Cycles
95°C	2 minutes	1
94°C	1 minute	40
55°C	1 minute	
72°C	1 minute	
72°C	5 minutes	1

9. Transfer 10 µl of each reaction to a new tube containing 2 µl 10X gel loading buffer.
10. Electrophorese the samples on a 1.5% agarose gel and analyze your results.
-

ProQuest™ Two-Hybrid System

The ProQuest™ Two-Hybrid System is an *in vivo* yeast-based system for identifying protein-protein interactions (Chevray and Nathans, 1992). The major features of the system are:

- Low copy (ARS/CEN) vectors for reduced toxicity
- Three reporter genes with independent promoters to reduce false positives due to non-specific interactions
- Large panel of control yeast strains
- Gateway®-compatible vectors for transferring your DNA sequences of interest into a variety of expression and analysis vectors using recombinational cloning

For more details on screening ProQuest™ Pre-made cDNA Libraries using the ProQuest™ Two-Hybrid System, refer to the ProQuest™ Two-Hybrid System manual. This manual is available for downloading from our Web site at www.invitrogen.com or by contacting Technical Service (see page 22).

Continued on next page

Using ProQuest™ Libraries, Continued

Gateway® Cloning

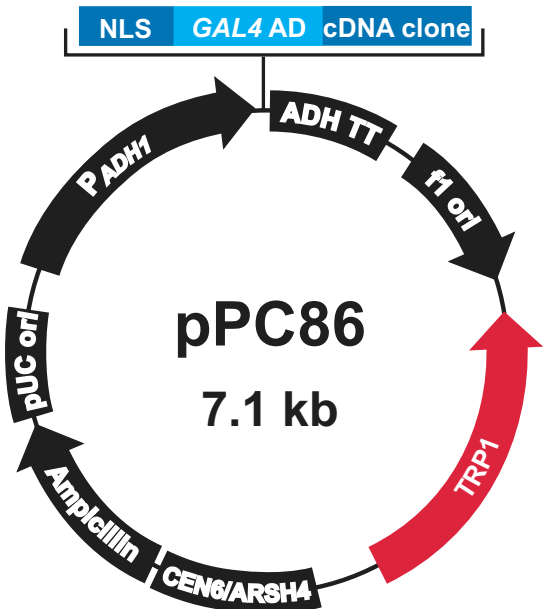
The vector, pEXP-AD502 contains *attB1* and *attB2* recombination sites flanking the cDNA cloning site. The cDNA insert can be transferred into other Gateway®-compatible vectors for expression by performing a BP recombination reaction with a pDONR™ vector (see page 18).

For details on Gateway® technology, refer to the Gateway® Technology Manual on our Web site at www.invitrogen.com or contact Technical Service (see page 22).

pPC86

Map of pPC86

The figure below shows the features of pPC86 vector. The complete sequence of pPC86 is available for downloading from our Web site (www.invitrogen.com) or by contacting Technical Service (see page 22).



Comments for pPC86 (no insert) 7093 nucleotides

ADH1 promoter: bases 114-1567

Initiation ATG: bases 1573-1575

SV40 nuclear localization signal (NLS): bases 1576-1596

GAL4 activation domain: bases 1606-1953

ADH1 transcription termination region: bases 2042-2508

f1 origin: bases 2700-3155

TRP1 gene (c): bases 3252-3926

TRP1 promoter (c): bases 3947-4029

CEN6/ARSH4 (c): bases 4458-4976

bla promoter: bases 5009-5113

Ampicillin (*bla*) resistance gene: bases 5108-5968

pUC origin: bases 6113-6786

(c) = complementary strand

Continued on next page

pPC86, Continued

Cloning Site of pPC86

The cloning site for pPC86 is shown below. Restriction sites are labeled to indicate the cleavage site.

1481 TTCTCTGTC A TTGTTCTCGT TCCCTTTCTT CCTTGTTTCT TTTTCTGCAC AATATTTCAA GCTATACCAA

1551 GCATACAATC AACTCCAAGC TT ATG CCC AAG AAG AAG CGG AAG GTC TCG AGC GGC GCC AAT
Met Pro Lys Lys Lys Arg Lys Val Ser Ser Gly Ala Asn
SV40 nuclear localization signal (NLS)

1612 TTT AAT CAA AGT GGG AAT ATT GCT GAT AGC TCA TTG TCC TTC ACT TTC ACT AAC AGT
Phe Asn Gln Ser Gly Asn Ile Ala Asp Ser Ser Leu Ser Phe Thr Phe Thr Asn Ser
GAL4 DNA activation domain

1669 AGC AAC GGT CCG AAC CTC ATA ACA ACT CAA ACA AAT TCT CAA GCG CTT TCA CAA CCA
Ser Asn Gly Pro Asn Leu Ile Thr Thr Gln Thr Asn Ser Gln Ala Leu Ser Gln Pro

1726 ATT GCC TCC TCT AAC GTT CAT GAT AAC TTC ATG AAT AAT GAA ATC ACG GCT AGT AAA
Ile Ala Ser Ser Asn Val His Asp Asn Phe Met Asn Asn Glu Ile Thr Ala Ser Lys

1783 ATT GAT GAT GGT AAT AAT TCA AAA CCA CTG TCA CCT GGT TGG ACG GAC CAA ACT GCG
Ile Asp Asp Gly Asn Asn Ser Lys Pro Leu Ser Pro Gly Trp Thr Asp Gln Thr Ala

Suggested forward sequencing/PCR primer

1840 TAT AAC GCG TTT GGA ATC ACT ACA GGG ATG TTT AAT ACC ACT ACA ATG GAT GAT GTA
Tyr Asn Ala Phe Gly Ile Thr Thr Gly Met Phe Asn Thr Thr Thr Met Asp Asp Val

1897 TAT AAC TAT CTA TTC GAT GAT GAA GAT ACC CCA CCA AAC CCA AAA AAA GAG GGT GGG
Tyr Asn Tyr Leu Phe Asp Asp Glu Asp Thr Pro Pro Asn Pro Lys Lys Glu Gly Gly
Sal I

1954 TCG ACC CAC GCG TCC G cDNA insert GCGGCCGCTA AGTAAGTAAG ACGTCGAGCT CTAAGTAAGT
Ser Thr His Ala Ser Not I Sst I

2022 AACGGCCGCC ACCGCGGTGG AGCTTTGGAC TTCTTCGCCA GAGGTTTGGT CAAGICTCCA ATCAAGGTTG
Sst II

Suggested reverse sequencing/PCR primer

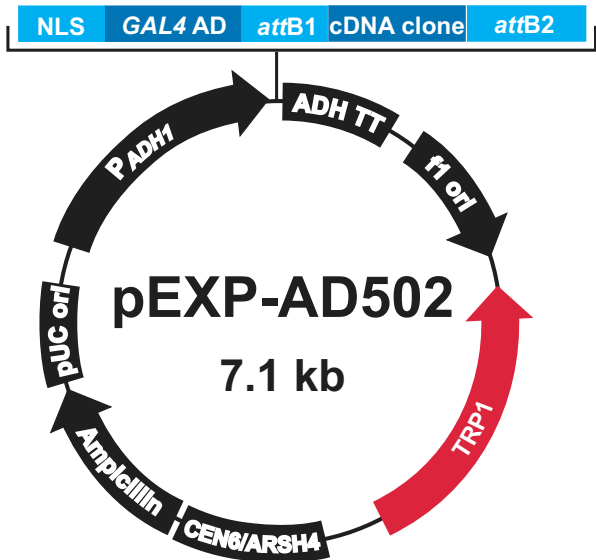
2092 TCGGCTTGTC TACCTTGCCA GAAATTTACG AAAAGAT

*This *Mlu* I site is contained within the *Sal* I adapter introduced into the vector upon ligation of the cDNA insert.

pEXP-AD502

Map of pEXP-AD502

The figure below shows the features of pEXP-AD502 vector. The complete sequence of pEXP-AD502 is available for downloading from our Web site (www.invitrogen.com) or by contacting Technical Service (see page 22).



Comments for pEXP-AD502 (no insert) 7146 nucleotides

ADH1 promoter: bases 64-1517

Initiation ATG: bases 1523-1525

SV40 nuclear localization signal (NLS): bases 1526-1548

GAL4 activation domain: bases 1556-1903

attB1: bases 1910-1934

attB2: bases 1979-2003

ADH1 transcription termination region: bases 2045-2511

f1 origin: bases 2703-3158

TRP1 gene (c): bases 3255-3929

TRP1 promoter (c): bases 3930-4031

CEN6/ARSH4 (c): bases 4461-4979

bla promoter: bases 5012-5116

Ampicillin (*bla*) resistance gene: bases 5111-5971

pUC origin: bases 6116-6789

(c) = complementary strand

Continued on next page

pEXP-AD502, Continued

Cloning Site of pEXP- AD502

The cloning site for pEXP-AD502 is shown below. Restriction sites are labeled to indicate the cleavage site.

```
1451 TCCCTTCTCT CCTTGTTTCT TTTTCTGCAC AATATTTCAA GCTATACCAA GCATACAATC AACTCCAAGC

1521 TT ATG CCC AAG AAG AAG CGG AAG GTC TCG AGC GGC GCC AAT TTT AAT CAA AGT GGG AAT
    Met Pro Lys Lys Lys Arg Lys Val Ser Ser Gly Ala Asn Phe Asn Gln Ser Gly Asn
    SV40 nuclear localization signal (NLS)

1580 ATT GCT GAT AGC TCA TTG TCC TTC ACT TTC ACT AAC AGT AGC AAC GGT CCG AAC CTC
    Ile Ala Asp Ser Ser Leu Ser Phe Thr Phe Thr Asn Ser Ser Asn Gly Pro Asn Leu
    GAL4 DNA activation domain

1637 ATA ACA ACT CAA ACA AAT TCT CAA GCG CTT TCA CAA CCA ATT GCC TCC TCT AAC GTT
    Ile Thr Thr Gln Thr Asn Ser Gln Ala Leu Ser Gln Pro Ile Ala Ser Ser Asn Val

1694 CAT GAT AAC TTC ATG AAT AAT GAA ATC ACG GCT AGT AAA ATT GAT GAT GGT AAT AAT
    His Asp Asn Phe Met Asn Asn Glu Ile Thr Ala Ser Lys Ile Asp Asp Gly Asn Asn
    Suggested forward sequencing/PCR primer

1751 TCA AAA CCA CTG TCA CCT GGT TGG ACG GAC CAA ACT GCG TAT AAC GCG TTT GGA ATC
    Ser Lys Pro Leu Ser Pro Gly Trp Thr Asp Gln Thr Ala Tyr Asn Ala Phe Gly Ile

1808 ACT ACA GGG ATG TTT AAT ACC ACT ACA ATG GAT GAT GTA TAT AAC TAT CTA TTC GAT
    Thr Thr Gly Met Phe Asn Thr Thr Thr Met Asp Asp Val Tyr Asn Tyr Leu Phe Asp

1865 GAT GAA GAT ACC CCA CCA AAC CCA AAA AAA GAG GGT GGG TCG ATC ACA AGT TTG TAC
    Asp Glu Asp Thr Pro Pro Asn Pro Lys Lys Glu Gly Gly Ser Ile Thr Ser Leu Tyr
    attB1

1922 AAA AAA GCA GGC TTG TCG ACC CAC GCG TCC G GGGCGGGCGC ACGCGTACCC
    Lys Lys Ala Gly Leu Ser Thr His Ala Ser
    Sal I Mu I Not I

1983 AGCTTCTTGG TACAAAGTGG TGACGTGCGAG CTCTAAGTAA GTAACGGCCG CCACCGCGGT GGAGCTTTGG
    Sst I Sst II
    attB2

2053 ACTTCTTCGC CAGAGGTTTG GTCAAGTCTC CAATCAAGGT TGTCGGCTTG TCTACCTTGC CAGAAATTTA
    Suggested reverse sequencing/PCR primer

2123 CGAAAAGATG GAAAAGGG
```

*This *Mlu* I site is contained within the *Sal* I adapter introduced into the vector upon ligation of the cDNA insert.

Recipe

Terrific Broth Terrific Broth is available from Invitrogen (see page 18 for ordering information).

1. Dissolve the following reagents in 800 ml of distilled water:

Tryptone	12 g
Yeast Extract	24 g
Glycerol	4 ml
 2. Mix well and adjust the volume to 900 ml with distilled water.
 3. Autoclave on liquid cycle for 20 minutes. Allow solution to cool to ~55°C
 4. Dissolve the following reagents in 80 ml of distilled water:

KH_2PO_4 (monobasic)	2.3 g
K_2HPO_4 (dibasic)	12.5 g
 5. Mix well and adjust the volume to 100 ml with distilled water.
 6. Autoclave on liquid cycle for 20 minutes. Allow solution to cool to ~55°C. Mix this solution with the solution prepared in Step 3.
 7. After the media is cooled, add antibiotic to the desired concentration.
 8. Store at +4°C.
-

Accessory Products

Additional Products

The table below lists additional products that may be used with the ProQuest™ Pre-made cDNA Libraries.

Product	Quantity	Catalog no.
ProQuest™ Two-Hybrid System with Gateway™ Technology	5 reactions	10835-031
MaV203 Competent Yeast Cells; Library Scale	2 x 0.55 ml	11281-011
Terrific Broth	500 g	22711-022
2.5 mM dNTP Mix	1 ml	R725-01
Platinum® <i>Taq</i> DNA Polymerase	100 reactions	10966-018
pDONR™ 221	6 µg	12536-017
pDONR™ 207	6 µg	12213-013
pDONR™ 221	6 µg	12536-017
Phenol:Chloroform:Isoamyl Alcohol, (25:24:1, v/v/v)	100 ml	15593-031

Purchaser Notification

**Limited Use
Label License
No. 11:
ProQuest™
Reverse Two-
Hybrid System**

The ProQuest™ Reverse Two-Hybrid System is the subject of one or more patents licensed to Invitrogen Corporation by Massachusetts General Hospital and Johns Hopkins University (including U.S. Patent Nos. 5,955,280 and/or 5,965,368 and foreign equivalents) ("Reverse Two-Hybrid"). 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. 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 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, Invitrogen is willing to accept return of the product with a full refund. For information on purchasing a license to this product for purposes other than research, contact Licensing Department, Invitrogen Corporation, 1600 Faraday Avenue, Carlsbad, California 92008. Phone (760) 603-7200. Fax (760) 602-6500.

Continued on next page

Purchaser Notification, Continued

Limited Use Label License No. 67: Two Hybrid Technology

Practice of the two-hybrid system is covered by U.S. Patents Nos. 5,283,173, 5,468,614 and 5,667,973 assigned to The Research Foundation of the State University of New York (SUNY). Purchase of any Invitrogen two-hybrid reagents does not imply or convey a license to practice the two-hybrid system covered by these patents, beyond use of the enclosed kit for non-commercial research. Commercial entities purchasing these reagents must obtain a license from The Research Foundation of the State University of New York before using them. For license information, please contact: Barbara A. Sawitsky, The Research Foundation of SUNY at Stony Brook, Office of Technology Licensing, W5530 Melville Memorial Library, Stony Brook, NY 11794-3368, Phone: 516-632-9077; Fax: 516-632-9839; E-mail: bsawitsky@notes.cc.sunysb.edu. You may not distribute the System or the vectors or host strains contained in it to others. You may not transfer modified, altered, or original material from the System to a third party without written notification to, and written approval from Invitrogen. You may not assign, sub-license, rent, lease or otherwise transfer any of the rights or obligations set forth herein, except as expressly permitted by Invitrogen.

Limited Use Label License No. 141: Expression of Poly-peptides in Yeast

This product is the subject of one or more of U.S. Patent Nos. 5,618,676, 5,854,018, 5,856,013, 5,919,651 and foreign equivalents. Rights to use this product are limited to academic research use only. For-profit entities are required to obtain a separate license from Washington Research Foundation to utilize this product for any use. Washington Research Foundation, 2815 Eastlake Avenue East, Suite 300, Seattle, Washington 98102. Tel: 206-336-5600. Fax: 206-336-5615.

Product Use by European Customers

These cells are genetically modified and contain the pUC-derived plasmid pPC86 or pEXP-AD502. As a condition of sale, this product must be used only according to applicable local legislation and guidelines, including EC Directive 90/219/EEC on the contained use of genetically modified organisms.

Continued on next page

Purchaser Notification, Continued

**Limited Use
Label License
No. 19:
Gateway®
Cloning
Products
(for catalog
nos.
11376-027,
11378-023,
11386-026
only)**

The Gateway® Cloning Technology products and their use are 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). No license is conveyed under the foregoing patents to use this product with any recombination sites other than those purchased from Invitrogen Corporation or its authorized distributor. 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 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. 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 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, Invitrogen is willing to accept return of the product with a full refund. For information on purchasing a license to this product for purposes other than research, contact Licensing Department, Invitrogen Corporation, 1600 Faraday Avenue, Carlsbad, California 92008. Phone (760) 603-7200. Fax (760) 602-6500.

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 at www.invitrogen.com.

United States 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
Tech 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).

1. On the home page, go to the left-hand column under 'Technical Resources' and select 'MSDS Requests'.
 2. Follow instructions on the page and fill out all the required fields.
 3. To request additional MSDSs, click the 'Add Another' button.
 4. All requests will be faxed unless another method is selected.
 5. When you are finished entering information, click the 'Submit' button. The MSDS will be sent within 24 hours.
-

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.

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).
- Chevray, P. M., and Nathans, D. (1992). Protein Interaction Cloning in Yeast: Identification of Mammalian Proteins that React with the Leucine Zipper of Jun. *Proc. Nat. Acad. Sci.* *89*, 5789-5793.
- Greenwald, R. J., Lu, P., Halvorson, M. J., Zhou, X., Chen, S., Madden, K. B., Perrin, P. J., Morris, S. C., Finkelman, F. D., Peach, R., Linsley, P. S., Urban, J. F. J., and Gause, W. C. (1997). Effects of Blocking B7-1 and B7-2 Interactions During a Type 2 In Vivo Immune Response. *J. Immunol.* *158*, 4088-4096.
- Kriegler, M. (1990). *Gene Transfer and Expression: A Laboratory Manual* (New York: Stockton Press).
- Li, W. B., Gruber, C. E., Lin, J. J., Lim, R., D'Alessio, J. M., and Jessee, J. A. (1994). The Isolation of Differentially Expressed Genes in Fibroblast Growth Factor Stimulated BC3H1 Cells by Subtractive Hybridization. *Biotechniques* *16*, 722-729.
- Sambrook, J., Fritsch, E. F., and Maniatis, T. (1989). *Molecular Cloning: A Laboratory Manual, Second Edition* (Plainview, New York: Cold Spring Harbor Laboratory Press).
- Svetic, A., Madden, K. B., Zhou, X. D., Lu, P., Katona, I. M., Finkelman, F. D., Urban, J. F. J., and Gause, W. C. (1993). A Primary Intestinal Helminthic Infection Rapidly Induces a Gut-Associated Elevation of Th2-Associated Cytokines and IL-3. *J. Immunol.* *150*, 3434-3441.

©2001-2002 Invitrogen Corporation. All rights reserved.

TRIzol[®] is a registered trademark of Molecular Research Center, Inc.



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

