

pYESTrp2

**Expression Vector Designed for use with the
Hybrid Hunter™ Yeast Two-Hybrid Systems**

Catalog no. V615-20

Version C
December 9, 2002
25-0301



www.invitrogen.com
tech_service@invitrogen.com

Table of Contents

Table of Contents.....	iii
Important Information.....	iv
Methods	1
Overview.....	1
Accessory Products.....	3
Cloning into pYESTrp2	6
Expressing the Prey Protein	11
Appendix.....	13
Recipes.....	13
pYESTrp2 Vector	16
Small-Scale Yeast Transformation	18
Preparation of Denatured Salmon Sperm DNA	19
Technical Service.....	20
Purchaser Notification	22
References.....	24

Important Information

Kit Contents 20 µg pYESTrp2, lyophilized in TE, pH 8.0

Shipping/Storage Lyophilized plasmid is shipped at room temperature and should be stored at -20°C.

Product Qualification The pYESTrp2 vector is qualified by restriction enzyme digestion with *Hind* III, *Bam*H I, and *Hind* III/*Pvu* II. Restriction digests must demonstrate the correct banding pattern when electrophoresed on an agarose gel. The table below lists the restriction enzymes used to digest the vector and the expected fragments.

Vector	Restriction Enzyme	Expected Fragments (bp)
pYESTrp2	<i>Hind</i> III	5822
	<i>Bam</i> H I	5822
	<i>Hind</i> III/ <i>Pvu</i> II	372, 5450

Methods

Overview

Introduction

pYESTrp2 is a 5.8 kb expression vector designed for use with the following Hybrid Hunter™ Systems available from Invitrogen:

Kit	Catalog no.
Hybrid Hunter™ System	K5000-01
Dual Bait Hybrid Hunter™ Yeast Two-Hybrid System	K5200-01

The vector allows expression of a prey protein of interest in *Saccharomyces cerevisiae* as a fusion to the B42 acidic activation domain. The prey may be a known protein or a cDNA library. Transformation of a prey library (in pYESTrp2) into the appropriate *Saccharomyces cerevisiae* bait strain allows identification of novel interactors with a bait (or baits) of interest while transformation of a known prey into the bait strain allows verification of an interaction between the prey and a known bait (or baits). The vector contains the following elements:

- Yeast *GAL1* promoter for galactose-inducible expression of the prey fusion protein in *S. cerevisiae*
- V5 epitope for detection of purified prey fusion protein with Anti-V5 Antibodies
- Nuclear localization signal (NLS) for localization of prey fusion protein to the nucleus
- B42 acidic activation domain for reconstituted transcriptional activation of LexA or cI operator-based reporter genes when brought into proximity with the LexA or cI fusion bait proteins
- A multiple cloning site with 9 unique restriction sites for simplified cloning of the prey protein of interest
- *CYCI* transcriptional terminator for efficient termination of mRNA
- *TRP1* gene for selection of transformants in yeast host strains with a *trp1* genotype
- Ampicillin resistance gene for selection of transformants in *E. coli*

For more information about the various yeast two-hybrid systems available from Invitrogen, refer to the manual for each system, our Web site (www.invitrogen.com), or call Technical Service (see page 20). For more information about the general principles of interaction trap systems, refer to the system manuals, published reviews (Bartel *et al.*, 1993a; Fields and Sternglanz, 1994), or general reference sources (Golemis *et al.*, 1996; Golemis *et al.*, 1998). To order components of the various Hybrid Hunter™ Systems, see pages 3-5.

Compatibility

pYESTrp2 contains the *TRP1* gene for selection of transformants in yeast, therefore, the vector may be used with any yeast host strain with a *trp1* genotype (see page 3 for more information). In principle, pYESTrp2 may be used with any LexA, lambda cI, or GAL4-based bait plasmid in a two-hybrid screen. The prey protein or library expressed from pYESTrp2 will be fused to the B42 activation domain, and should activate transcription of specific reporter genes when an interaction with an appropriate fusion bait has occurred.

Continued on next page

Overview, continued

Experimental Outline

The table below outlines the major steps required to clone and express your prey protein of interest in pYESTrp2.

Step	Action
1	Consult the multiple cloning site diagram on page 8 to design a strategy to clone your prey gene in frame with sequence encoding the V5 epitope-NLS-B42 fusion in pYESTrp2.
2	Ligate your insert into pYESTrp2 and transform into <i>E. coli</i> . Select transformants on LB plates containing 50 µg/ml ampicillin.
3	Analyze your transformants for the presence of insert by restriction digestion.
4	Select a transformant with the correct restriction pattern and sequence to confirm that your gene is cloned in the proper orientation and in frame with the V5 epitope-NLS-B42 fusion protein.
5	Use a small-scale transformation protocol to transform your pYESTrp2 prey construct into the appropriate yeast host strain and select for tryptophan prototrophy.
6	Test for expression of your prey fusion protein by Western blot analysis, if desired.
7	Use a small-scale transformation protocol to transform your pYESTrp2 prey construct into the appropriate yeast bait strain and select for tryptophan prototrophy. Test for an interaction between your bait and prey by assaying for the appropriate reporter activity. If you are performing an interactor hunt, use a small-scale or large-scale library transformation protocol to transform your pYESTrp2 library into the appropriate yeast bait strain and select as above.

Accessory Products

Introduction

The products listed below are designed for use with the pYESTrp2 vector and the various Hybrid Hunter™ Systems available from Invitrogen. For more information about the components specific for each Hybrid Hunter™ system, refer to the manual for the particular system.

Yeast Strains

For your convenience, Invitrogen has available many yeast host strains that may be used with the pYESTrp2 vector in an interactor hunt. All of the strains carry the *trp1* genotype to allow selection of pYESTrp2 transformants. The strains have been specially designed for use with a specific Hybrid Hunter™ System. This information as well as the genotype of each strain is listed in the table below. In some cases, a particular yeast strain has been transformed with a LexA and/or cI-based reporter plasmid to facilitate detection of molecular interactions. For more information about a particular yeast strain, refer to the specific system manual or call Technical Service (see page 20).

Each of the yeast strains is supplied as a 20% glycerol stock in a 0.5 ml volume. The pSH18-34 and pLacGUS reporter plasmids may be ordered separately from Invitrogen (see page 5). Other reporter plasmids are also available (see page 5 for ordering information).

Yeast Strain	Genotype	For Use With...	Catalog no.
L40	<i>MATa his3Δ200 trp1-901 leu2-3112 ade2 LYS2::(4lexAop-HIS3) URA3::(8lexAop-lacZ) GAL4</i>	Hybrid Hunter™	C830-00
EGY48	<i>MATα ura3 trp1 his3 6lexAop-LEU2</i>	Hybrid Hunter™	C835-00
EGY48/pSH18-34	<i>MATα ura3 trp1 his3 6lexAop-LEU2 pSH18-34(URA3)</i>	Hybrid Hunter™	C836-00
EGY191	<i>MATα ura3 trp1 his3 2lexAop-LEU2</i>	Hybrid Hunter™	C837-00
EGY191/pSH18-34	<i>MATα ura3 trp1 his3 2lexAop-LEU2 pSH19-34(URA3)</i>	Hybrid Hunter™	C838-00
SKY48	<i>MATα ura3 trp1 his3 6lexAop-LEU2 3cI-op-LYS2</i>	Dual Bait Hybrid Hunter™	C833-00
SKY48/pLacGUS	<i>MATα ura3 trp1 his3 6lexAop-LEU2 3cI-op-LYS2 pLacGUS (URA3)</i>	Dual Bait Hybrid Hunter™	C832-00
SKY191	<i>MATα ura3 trp1 his3 2lexAop-LEU2 3cI-op-LYS2</i>	Dual Bait Hybrid Hunter™	C834-00

Continued on next page

Accessory Products, continued

Hybrid Hunter™ Libraries

The following cDNA libraries in pYESTrp and pYESTrp2 are available from Invitrogen. All of the libraries are compatible for use with any of the Hybrid Hunter™ Systems. All libraries are amplified once on plates. For more information, see our Web site (www.invitrogen.com) or call Technical Service (see page 20).

Source	Plasmid	Primary Clones	Size Selection	Catalog no.
Cell Lines				
HeLa cells (Human cervical carcinoma)	pYESTrp	3.66 x 10 ⁶	0.3 to 1.2 kb	A201-01
BeWo cells (Human fetal placental choriocarcinoma)	pYESTrp	5.35 x 10 ⁶	0.3 to 0.8 kb	A208-01
BeWo cells (Human fetal placental choriocarcinoma)	pYESTrp	5.5 x 10 ⁶	0.3 to 1.2 kb	A208-02
Jurkat cells (Human T cell leukemia)	pYESTrp	3.2 x 10 ⁶	0.3 to 1.2 kb	A209-01
A20 cells (Mouse B cell lymphoma)	pYESTrp	3.11 x 10 ⁶	0.3 to 1.2 kb	A210-01
Human Adult Tissue				
Bladder	pYESTrp2	17.6 x 10 ⁶	0.4 to 1.2 kb	A225-01
Brain	pYESTrp2	10.8 x 10 ⁶	0.4 to 1.2 kb	A204-01
Breast	pYESTrp2	9.00 x 10 ⁶	0.4 to 1.2 kb	A217-01
Breast Tumor	pYESTrp2	8.84 x 10 ⁶	0.4 to 1.2 kb	A216-01
Colon Tumor	pYESTrp2	7.98 x 10 ⁶	0.4 to 1.2 kb	A222-01
Kidney	pYESTrp2	6.96 x 10 ⁶	0.4 to 1.2 kb	A223-01
Liver	pYESTrp	2.21 x 10 ⁶	0.3 to 1.2 kb	A203-01
Lung	pYESTrp2	5.95 x 10 ⁶	0.4 to 1.2 kb	A213-01
Lung Tumor	pYESTrp2	1.85 x 10 ⁶	0.4 to 1.2 kb	A215-01
Ovary	pYESTrp	4.54 x 10 ⁶	0.3 to 1.2 kb	A206-01
Placenta	pYESTrp	4.75 x 10 ⁶	0.3 to 1.2 kb	A207-01
Prostate	pYESTrp2	5.46 x 10 ⁶	0.4 to 1.2 kb	A218-01
Spleen	pYESTrp2	11.4 x 10 ⁶	0.4 to 1.2 kb	A214-01
Testes	pYESTrp	6.4 x 10 ⁶	0.3 to 1.2 kb	A205-01
Human Fetal Tissue				
Fetal Liver	pYESTrp	2.37 x 10 ⁶	0.3 to 1.2 kb	A202-01

Continued on next page

Accessory Products, continued

Additional Reagents

Additional reagents that may be used with pYESTrp2 and the various Hybrid Hunter™ Systems are available from Invitrogen. Ordering information is provided below. The quantity of antibody supplied is sufficient for 25 Westerns.

Item	Amount	Catalog no.
<i>S.c.</i> EasyComp™ Kit	20 reactions	K5050-01
pHybLex/Zeo (Bait Vector)	20 µg	V610-20
pHybcl/HK (Bait Vector)	20 µg	V614-20
pSH18-34 Reporter Plasmid	20 µg	V611-20
pRB1840 Reporter Plasmid	20 µg	V612-20
pJK103 Reporter Plasmid	20 µg	V613-20
pLacGUS Reporter Plasmid	20 µg	V616-20
Anti-V5 Antibody	50 µl	R960-25
Anti-V5-HRP Antibody	50 µl	R961-25

Cloning into pYESTrp2

Introduction

The pYESTrp2 vector (5822 bp) can be used to make two-hybrid cDNA libraries or to clone genes encoding known proteins. Use the diagram on page 8 to help you clone your gene of interest into pYESTrp2. General considerations for cloning and transformation are discussed below.

General Molecular Biology Techniques

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

E. coli Strain

Many *E. coli* strains are suitable for the propagation of the pYESTrp2 vector including TOP10, TOP10F', DH5 α or equivalent. We recommend that you propagate the vector in *E. coli* strains that are recombination deficient (*recA*) and endonuclease A deficient (*endA*). For your convenience, TOP10 *E. coli* are available as chemically competent or electrocompetent cells from Invitrogen.

Item	Quantity	Catalog no.
One Shot™ TOP10 (chemically competent cells)	21 x 50 μ l	C4040-03
Electrocomp™ TOP10 (electrocompetent cells)	5 x 80 μ l	C664-55

Transformation Method

You may use any method of your choice for transformation. Chemical transformation is the most convenient for most researchers. Electroporation is the most efficient and the method of choice for large plasmids.

Maintenance of pYESTrp2

To propagate and maintain the pYESTrp2 vector, we recommend resuspending the vector in 20 μ l sterile water to prepare a 1 μ g/ μ l stock solution. Store the stock solution at -20°C. Use this stock solution to transform a *recA*, *endA* *E. coli* strain like TOP10, TOP10F' or equivalent. Select transformants on LB plates containing 50 μ g/ml ampicillin. Be sure to prepare a glycerol stock of the plasmid for long-term storage (see page 9 for a protocol).

B42 Activation Domain

Your prey protein of interest will be fused to the *E. coli* B42 acidic activation domain in pYESTrp2 (Ruden *et al.*, 1991). The B42 activation domain acts as a transcriptional activator when brought into proximity with a DNA binding protein such as LexA (Horii *et al.*, 1981; Markham *et al.*, 1981) or cI (Nilsson *et al.*, 1983). Transcriptional activation then occurs via binding of the LexA or cI protein to LexA or cI operators upstream of heterologous genes (i.e. reporter genes).

Continued on next page

Cloning into pYESTrp2, continued

Cloning Considerations

When designing your cloning strategy, remember that you must clone your gene in frame with the sequence encoding the V5 epitope-NLS-B42 fusion protein in order to create a prey fusion protein with a nuclear localization signal, activation domain, and an epitope for detection.



Note

The N-terminal peptide contains a V5 epitope to allow detection of your expressed prey fusion protein by immunoblot (Western analysis). Anti-V5 antibodies are available from Invitrogen to facilitate detection (see page 12 for more information).

Continued on next page

Cloning into pYESTrp2, continued

Multiple Cloning Site of pYESTrp2

Below is the multiple cloning site for pYESTrp2. Restriction sites are labeled to indicate the cleavage site. The multiple cloning site has been confirmed by sequencing and functional testing. **The complete sequence of pYESTrp2 is available for downloading from our Web site (www.invitrogen.com) or from Technical Service (see page 20).** For a map and a description of the features of pYESTrp2, see pages 16-17.

```

5' end of GALI promoter
1  CGCGCTTAAT GGGGCGCTAC AGGGCGCGTG GGGATGATCC ACTAGTACGG ATTAGAAGCC

61  GCCGAGCGGG TGACAGCCCT CCGAAGGAAG ACTCTCTCC GTGCGTCCTC GTCTTCACCG

121 GTCGCGTTCC TGAACGCAG ATGTGCCTCG CGCCGCACTG CTCCGAACAA TAAAGATTCT

181 ACAATACTAG CTTTTATGGT TATGAAGAGG AAAAATTGGC AGTAACCTGG CCCACAAAC

241 CTTCAAATGA ACGAATCAAA TTAACAACCA TAGGATGATA ATGCGATTAG TTTTTAGCC

301 TTATTTCTGG GGTAATTAAT CAGCGAAGCG ATGATTTTTG ATCTATTAAC AGATATATAA
                                     TATA Box

361 ATGCAAAAAC TGCATAACCA CTTTAACTAA TACTTTCAAC ATTTTCGGTT TGTATTACTT

421 CTTATTCAA  TGTAATAAAA GTATCAACAA AAAATTGTTA ATATACCTCT ATACTTTAAC
                                     transcriptional start →

481 GTCAAGGAGA AAAAACCCCG GATCGGACTA CTAGCAGCTG TAATACGACT CACTATAGGG
                                     T7 promoter/priming site

541 AATATTAAGC TCACC ATG GGT AAG CCT ATC CCT AAC CCT CTC CTC GGT CTC
    Met Gly Lys Pro Ile Pro Asn Pro Leu Leu Gly Leu
                                     V5 epitope

592 GAT TCT ACA CAA GCT ATG GGT GCT CCT CCA AAA AAG AAG AGA AAG GTA GCT
    Asp Ser Thr Gln Ala Met Gly Ala Pro Pro Lys Lys Lys Arg Lys Val Ala
                                     SV40 NLS

643 GGT ATC AAT AAA GAT ATC GAG GAG TGC AAT GCC ATC ATT GAG CAG TTT ATC
    Gly Ile Asn Lys Asp Ile Glu Glu Cys Asn Ala Ile Ile Glu Gln Phe Ile

694 GAC TAC CTG CGC ACC GGA CAG GAG ATG CCG ATG GAA ATG GCG GAT CAG GCG
    Asp Tyr Leu Arg Thr Gly Gln Glu Met Pro Met Glu Met Ala Asp Gln Ala
                                     B42 activation domain

745 ATT AAC GTG GTG CCG GGC ATG ACG CCG AAA ACC ATT CTT CAC GCC GGG CCG
    Ile Asn Val Val Pro Gly Met Thr Pro Lys Thr Ile Leu His Ala Gly Pro

796 CCG ATC CAG CCT GAC TGG CTG AAA TCG AAT GGT TTT CAT GAA ATT GAA GCG
    Pro Ile Gln Pro Asp Trp Leu Lys Ser Asn Gly Phe His Glu Ile Glu Ala

847 GAT GTT AAC GAT ACC AGC CTC TTG CTG AGT GGA GAT GCC TCC AAG CTT GGT
    Asp Val Asn Asp Thr Ser Leu Leu Leu Ser Gly Asp Ala Ser Lys Leu Gly
                                     Hind III

898 ACC GAG CTC GGA TCC ACT AGT AAC GGC CGC CAG TGT GCT GGA ATT CTG CAG
    Thr Glu Leu Gly Ser Thr Ser Asn Gly Arg Gln Cys Ala Gly Ile Leu Gln
    Kpn I Sac I BamH I BstX I* EcoR I

949 ATA TCC ATC ACA CTG GCG GCC GCT CGA GGC ATG CAT CTA GAG GGC CGC ATC
    Ile Ser Ile Thr Leu Ala Ala Ala Arg Gly Met His Leu Glu Gly Arg Ile
    BstX I* Not I Xho I Sph I

1000 ATG TAA TTAGTTA TGTACGCTT ACATTCACGC CCTCCCCCA
    Met ***
  
```

*Please note that there are two *BstX I* sites in the polylinker.

Continued on next page

Cloning into pYESTrp2, continued

E. coli **Transformation**

Once you have completed your ligation reaction, transform your ligation mixture into a competent *recA*, *endA* *E. coli* strain (e.g. TOP10, TOP10F'). Select for transformants on LB plates containing 50 µg/ml ampicillin. Select 10-20 clones and analyze by restriction digest or sequencing for the presence and orientation of your insert.



We recommend that you sequence your construct to confirm that your gene is fused in frame with the V5 epitope, nuclear localization signal, and B42 activation domain. For your convenience, Invitrogen offers a Custom Primer service and can construct primers according to your specifications. For more information on this service, visit our Web site (www.invitrogen.com) or call Technical Service (see page 20).

Preparing a Glycerol Stock

Once you have identified the correct clone, be sure to purify the colony and make a glycerol stock for long-term storage. It is also a good idea to keep a DNA stock of your plasmid at -20°C in case you lose the glycerol stock.

1. Streak the original colony out on an LB plate containing 50 µg/ml ampicillin. Incubate the plate at 37°C overnight.
 2. Isolate a single colony and inoculate into 1-2 ml of LB containing 50 µg/ml ampicillin.
 3. Grow the culture to mid-log phase (OD₆₀₀ = 0.5-0.7).
 4. Mix 0.85 ml of culture with 0.15 ml of sterile glycerol and transfer to a cryovial.
 5. Store at -80°C.
-

Constructing a cDNA Library

You may use pYESTrp2 to construct your own two-hybrid cDNA library. General guidelines are provided below to generate a unidirectional cDNA library in pYESTrp2. Refer to *Current Protocols in Molecular Biology*, Unit 5 (Ausubel *et al.*, 1994) for details on cDNA library construction.

1. Isolate mRNA from the source of interest.
 2. Prepare first strand cDNA using random primers, Oligo dT(*Not* I) primer (Catalog no. N430-01), or an Oligo dT(*Xho* I) primer.
Alternatively, the Copy™ Kit (Catalog no. L1311-03) is available from Invitrogen for efficient production of double-stranded blunt-ended cDNA for either bidirectional or unidirectional cloning. Call Technical Service (see page 20 for more information).
 3. After second strand synthesis, be sure the ends are blunt prior to adding *Bst*X I/*Eco*R I adaptors. *Bst*X I/*Eco*R I adaptors (Catalog no. N418-18) are available from Invitrogen.
 4. Digest with *Not* I or *Xho* I and electrophorese on an agarose gel for size selection.
 5. Isolate cDNA for ligation into pYESTrp2.
 6. Digest pYESTrp2 with either *Bst*X I (or *Eco*R I) and *Not* I (or *Xho* I) to complement the ends on the cDNA.
 7. Ligate cDNA into digested vector and transform into *E. coli*.
 8. Determine the number of primary recombinants. You may wish to amplify the library prior to large-scale isolation of plasmid DNA for the library screen.
-

Continued on next page

Cloning into pYESTrp2, continued

Small-Scale Plasmid Preparation

You may use any method of your choice to prepare purified plasmid DNA for small-scale yeast transformation. Standard protocols can be found in *Current Protocols in Molecular Biology* (Ausubel *et al.*, 1994) or *Molecular Cloning: A Laboratory Manual* (Sambrook *et al.*, 1989).

If you need ultrapure DNA for sequencing, we recommend isolating plasmid DNA using the S.N.A.P.[™] MiniPrep Kit (10-15 µg DNA, Catalog no. K1900-01) or the S.N.A.P.[™] MidiPrep Kit (10-200 µg DNA, Catalog no. K1910-01).

Large-Scale Plasmid Preparation

If you have constructed a cDNA library in pYESTrp2, you will need to isolate purified plasmid DNA before you can proceed with library transformation. You will need 30 µg of library plasmid DNA to perform a small-scale library transformation and 500 µg of library plasmid DNA to perform a large-scale library transformation. To isolate plasmid DNA, follow the procedure below. Other methods are suitable. If you are using a Hybrid Hunter[™] Premade Library, follow the directions supplied with your library.

1. Inoculate 1-2 liters of LB medium containing 50 µg/ml ampicillin with sufficient bacterial library stock to insure 2-3 times the number of independent clones in the library.
 2. Incubate at 37°C overnight with shaking.
 3. After incubation, pellet the cells and proceed with large- or mega-scale isolation of plasmid DNA. Any standard method is suitable. You may have to adjust the plasmid preparation protocol to account for the density of the culture.
 4. Store the plasmid at -20°C until ready for use.
-

Expressing the Prey Protein

Introduction

If you wish to test for expression of your prey fusion protein before proceeding with an interactor hunt, you may use a small-scale transformation protocol to transform your pYESTrp2 prey construct into a suitable yeast host strain. Once you have selected transformants for tryptophan prototrophy, expression of the prey fusion protein may be induced by addition of galactose to the medium. A procedure to prepare lysates for immunoblot (Western) analysis is provided below.

Basic Yeast Molecular Biology

The user should be familiar with basic yeast molecular biology and microbiological techniques. Refer to *Current Protocols in Molecular Biology* (1996) *Saccharomyces cerevisiae*, pp. 13.01 to 13.2.12 for information on preparing yeast media and handling yeast.

Reagents for Yeast Transformation

The *S. c.* EasyComp™ Kit (Catalog no. K5050-01) provides a quick and easy method to prepare competent yeast cells that can be used immediately or stored frozen for future use. Transformation efficiency is guaranteed at $>10^3$ transformants per μg DNA.

For your convenience, a small-scale transformation protocol is included in the **Appendix**, page 18. Alternatively, there are published references for other small-scale transformation methods (Gietz *et al.*, 1992; Gietz and Schiestl, 1996; Hill *et al.*, 1991; Schiestl and Gietz, 1989).

Materials Required

To assay for expression of your prey fusion protein by immunoblot (Western blot) analysis, be sure to have the following reagents and equipment on hand before proceeding:

- 30°C incubator and shaking incubator
 - 60°C and 70°C water baths or temperature blocks and a boiling water bath
 - Clinical centrifuge and low-speed centrifuge
 - Selective medium and plates
 - Cracking buffer (see page 15 for a recipe), prewarmed to 60°C
 - Acid washed glass beads (Sigma-Aldrich G-8772, 425-600 microns)
 - Reagents for SDS-PAGE and immunoblotting
 - Antibody to your prey protein or Anti-V5 Antibody (see the next page)
-

Continued on next page

Expressing the Prey Protein, continued

Expression of the Prey Fusion Protein

Use the protocol below to prepare lysates from your Trp⁺ transformants and untransformed yeast cells for Western blot analysis. Test several transformants in case of heterogeneity in prey expression levels.

1. Inoculate 10 ml of the appropriate selective medium with a single colony of your prey strain (above) and inoculate 10 ml of the appropriately-supplemented YC medium with your untransformed yeast strain as a negative control. Remember that the medium needs to contain galactose to induce expression of the prey protein. Grow overnight with shaking at 30°C.
 2. Pellet the cells in Step 1 by centrifuging at 2500 rpm for 5 minutes at room temperature. Decant the medium.
 3. Transfer the cell pellets to a -80°C freezer for 10 minutes.
 4. Thaw cell pellet in 100 µl of prewarmed (60°C) cracking buffer and resuspend by pipetting the cell pellet in the buffer.
 5. Transfer cell suspension to a 1.5 ml microcentrifuge tube containing 100 µl of glass beads.
 6. Incubate the solution at 70°C for 10 minutes.
 7. Vortex solution for 1 minute.
 8. Centrifuge in a microcentrifuge at 14,000 rpm for 5 minutes at room temperature and transfer supernatant to a new tube.
 9. Add SDS-PAGE sample buffer and boil sample for 5 minutes. Use 30 to 50 µl for immunoblot analysis. Detect the prey fusion protein using an antibody to your prey of interest or Anti-V5 antibodies available from Invitrogen (see below).
-

Antibodies for Detection

The prey fusion protein contains the N-terminal V5 epitope to allow detection of the expressed prey protein by Western blot analysis. The Anti-V5 Antibody (Catalog no. R960-25) and the Anti-V5-HRP Antibody (Catalog no. R961-25) are available from Invitrogen for detection of the expressed prey protein (see page 5 for ordering information). For the sequence of the V5 epitope, refer to the diagram on page 8.



Note

Note that the N-terminal peptide containing the V5 epitope, nuclear localization signal, and B42 activation domain will add approximately 12 kDa to the size of your prey protein.

Performing a Two-Hybrid Screen

Once you have verified expression of your prey fusion protein in yeast, you may proceed to assay for an interaction between your prey protein and the bait (or baits) of interest or to identify novel interactors for your bait of interest. For details, refer to the manual for the appropriate Hybrid Hunter™ System.

Retrieving Putative Interactors

There are several methods available to retrieve the prey plasmid or the gene encoding a putative interactor. We generally use PCR to amplify the prey gene of interest and clone it into a PCR cloning vector. Other methods allow retrieval of the original prey plasmid by isolation of yeast plasmid DNA or by plasmid segregation. Further details and protocols for each of these methods are provided in each Hybrid Hunter™ System manual.

Appendix

Recipes

YC Medium and Plates

YC is minimal defined medium for yeast.

0.12% yeast nitrogen base (**without either** amino acids or ammonium sulfate)

0.5% ammonium sulfate

1% succinic acid

0.6% NaOH

2% glucose

0.01% (adenine, arginine, cysteine, leucine, lysine, threonine, tryptophan, uracil)

0.005% (aspartic acid, histidine, isoleucine, methionine, phenylalanine, proline, serine, tyrosine, valine)

2% agar (for plates)

1. Dissolve the following reagents in 900 ml deionized water. **Note:** We make medium and plates as we need them and weigh out each amino acid. Many researchers prepare 100X solutions of each amino acid mix that they need.

1.2 g Yeast Nitrogen Base

0.1 g each

0.05 g each

5 g Ammonium sulfate

adenine

aspartic acid

10 g Succinic acid

arginine

histidine

6 g NaOH

cysteine

isoleucine

leucine

methionine

lysine

phenylalanine

threonine

proline

tryptophan (W)

serine

uracil

tyrosine

valine

Note: The amino acids with the one letter code are those you need to omit to make selective plates, depending on the genotype of the host, plasmid markers, and reporters.

2. If you are making plates, add the agar after dissolving the reagents above.
3. Autoclave at 15 psi, 121°C for 20 minutes.
4. Cool to 50°C and add 100 ml of filter-sterilized 20% glucose. **Note:** You may add the sugar before autoclaving; however, the medium will be darker in color because of heating the glucose.
For plates that contain galactose and raffinose, add 100 ml 20% galactose and 50 ml 20% raffinose instead of glucose.
5. Pour plates and allow to harden. Invert the plates and store at +4°C. Plates are stable for up to 6 months.



Note

The recipe for YC medium has been optimized for use with the Hybrid Hunter™ Systems and its components. Other recipes may be suitable, but should be tested with the host strain, plasmid markers, and reporters.

Continued on next page

Recipes, continued

YPD

Yeast Extract Peptone Dextrose Medium (1 liter)

1% yeast extract
2% peptone
2% dextrose (D-glucose)

1. Dissolve the following in 1000 ml of water:
 - 10 g yeast extract
 - 20 g peptone
 - 20 g dextrose (see note below if making plates)
2. Optional: Add 20 g agar, if making plates.
3. Autoclave for 20 minutes on liquid cycle.
4. Store medium at room temperature or cool the medium and pour plates. The shelf life is approximately one to two months.

Note: If making plates, omit dextrose from Step 1. Autoclaving agar and dextrose together will cause the dextrose to caramelize. Prepare a separate stock solution of 20% dextrose and autoclave or filter-sterilize. After the YPD broth has been autoclaved, add 100 ml of 20% dextrose to the medium.

10X TE

100 mM Tris, pH 7.5
10 mM EDTA

1. For 100 ml, dissolve 1.21 g of Tris base and 0.37 g of EDTA in 90 ml of deionized water.
2. Adjust the pH to 7.5 with concentrated HCl and bring the volume up to 100 ml.
3. Filter sterilize and store at room temperature.

Alternatively, you can make the solution using 1 M Tris-HCl, pH 7.5 and 0.5 M EDTA, pH 8.0.

1X TE

10 mM Tris, pH 7.5
1 mM EDTA

Dilute 10X TE 10-fold with sterile water.

10X LiAc

1 M Lithium Acetate, pH 7.5

1. For 100 ml, dissolve 10.2 g of lithium acetate in 90 ml of deionized water.
 2. Adjust pH to 7.5 with dilute glacial acetic acid and bring up the volume to 100 ml.
 3. Filter sterilize and store at room temperature.
-

1X LiAc

100 mM Lithium Acetate, pH 7.5

Dilute 10X LiAc solution 10-fold with sterile, deionized water.

Continued on next page

Recipes, continued

1X LiAc/0.5X TE

100 mM Lithium Acetate, pH 7.5
5 mM Tris-HCl, pH 7.5
0.5 mM EDTA

1. For 100 ml, mix together 10 ml of 10X LiAc and 5 ml of 10X TE.
 2. Add deionized water to 100 ml.
 3. Filter-sterilize and store at room temperature.
-

1X LiAc/40% PEG-3350/1X TE

100 mM Lithium Acetate, pH 7.5
40% PEG-3350
10 mM Tris-HCl, pH 7.5
1 mM EDTA

1. Prepare solution immediately prior to use. For 100 ml, mix together 10 ml of 10X LiAc, 10 ml of 10X TE, and 40 g of PEG-3350.
 2. Add deionized water to 100 ml and dissolve the PEG. You may have to heat the solution to fully dissolve the PEG.
 3. Autoclave at 121°C, 15 psi for 20 minutes. Store at room temperature.
-

Cracking Buffer

8 M urea
5% SDS
40 mM Tris-HCl pH 6.8
0.1 mM EDTA
1% β -mercaptoethanol
0.4 mg/ml bromophenol blue

1. Prepare a 1 M Tris-HCl, pH 6.8 stock. (12.11 g in 90 ml deionized water and adjust pH to 6.8. Bring the volume to 100 ml).

2. Mix together the following reagents:

Urea	48.0 g
SDS	5 g
1 M Tris-HCl, pH 6.8	4 ml
EDTA	3.72 mg (or 20 μ l of a 0.5 M stock)
β -mercaptoethanol	1 ml
Bromophenol blue	40 mg

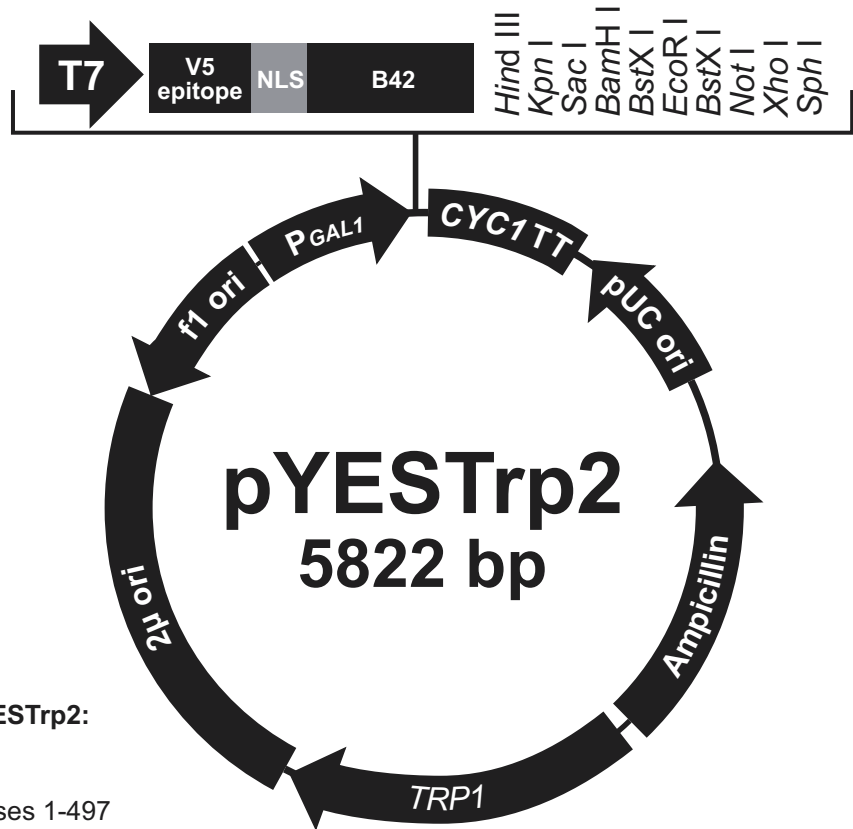
Bring up in 100 ml deionized water and dissolve reagents.

3. Store at +4°C or -20°C.
-

pYESTrp2 Vector

Map of pYESTrp2

pYESTrp2 is a 5822 bp prey vector that can be used to make two-hybrid cDNA libraries or to clone genes encoding known proteins. The figure below summarizes the features of the pYESTrp2 vector. **The complete nucleotide sequence for pYESTrp2 is available for downloading from our Web site (www.invitrogen.com) or by contacting Technical Service (see page 20).**



Comments for pYESTrp2: 5822 nucleotides

GAL1 promoter: bases 1-497

T7 promoter/priming site: bases 521-540

Initiation ATG: bases 556-558

V5 epitope: bases 559-600

Nuclear localization signal (NLS): bases 616-642

B42 activation domain: bases 646-883

Multiple cloning site: bases 889-982

CYC1 transcription termination region: bases 997-1245

pUC origin: bases 1427-2100 (complementary strand)

Ampicillin (*bla*) resistance gene: bases 2219-3105 (complementary strand)

bla promoter: bases 3106-3204 (complementary strand)

TRP1 promoter: bases 3313-3414

TRP1 gene: bases 3415-4089

2μ origin: bases 4493-5327

f1 origin: bases 5396-5768 (complementary strand)

Continued on next page

pYESTrp2 Vector, continued

Features of pYESTrp2

The table below describes the features of pYESTrp2. All elements have been functionally tested.

Feature	Benefit
<i>GAL1</i> promoter	Allows galactose-inducible expression of prey genes cloned into pYESTrp2
T7 promoter/priming site	Permits sequencing of insert or <i>in vitro</i> transcription of sense strand
V5 epitope	Allows detection of fusion protein(s) using the Anti-V5 Antibody (Catalog no. R960-25) or Anti-V5-HRP Antibody (Catalog no. R961-25) (Southern <i>et al.</i> , 1991)
SV40 large T antigen nuclear localization sequence (NLS)	Localizes prey fusions to the nucleus for potential interaction with bait fusions
B42 activation domain (AD) ORF	Transcriptional activation domain that allows expression of reporter genes when brought into proximity with an appropriate DNA binding domain (DBD) (i.e. LexA or lambda cI) by two interacting proteins (Ma and Ptashne, 1987)
Multiple cloning site with 8 unique sites, plus two <i>BstX</i> I sites.	Allows in-frame cloning of a cDNA library or a single gene with the B42 activation domain
<i>CYC1</i> transcription termination signal	Permits efficient termination and stabilization of mRNA
pUC origin	Maintenance and high-copy replication in <i>E. coli</i>
<i>bla</i> promoter	Allows expression of the <i>bla</i> resistance gene
Ampicillin (<i>bla</i>) resistance gene	Allows selection of transformants in <i>E. coli</i>
<i>TRP1</i> promoter	Allows expression of the <i>TRP1</i> gene
<i>TRP1</i> gene	Permits auxotrophic selection of the plasmid in Trp ⁻ yeast hosts (Tschumper and Carbon, 1980)
2 μ origin	Permits maintenance and high-copy replication in yeast
f1 origin	Allows rescue of single-stranded DNA

Small-Scale Yeast Transformation

Introduction

A small-scale yeast transformation protocol for routine transformations is provided below. Other protocols are suitable.

Materials Needed

Be sure to have the following reagents on hand before starting.

- YPD liquid medium (see recipe, page 14)
 - 1X TE (see recipe, page 14)
 - 1X LiAc/0.5X TE (see recipe, page 15)
 - Denatured salmon sperm DNA (see recipe, next page)
 - Plasmid DNA to be transformed
 - 1X LiAc/40% PEG-3350/1X TE (see recipe, page 15)
 - DMSO
 - Selective plates
-

Protocol

1. Inoculate 10 ml of YPD medium with a colony of the appropriate yeast strain and shake overnight at 30°C.
 2. Determine the OD₆₀₀ of your overnight culture. Dilute culture to an OD₆₀₀ of 0.4 in 50 ml of YPD medium and grow an additional 2-4 hours.
 3. Pellet the cells at 2500 rpm and resuspend the pellet in 40 ml 1X TE.
 4. Pellet the cells at 2500 rpm and resuspend the pellet in 2 ml of 1X LiAc/0.5X TE.
 5. Incubate the cells at room temperature for 10 minutes with shaking.
 6. For each transformation, mix together 1 µg plasmid DNA and 100 µg denatured sheared salmon sperm DNA with 100 µl of the yeast suspension from Step 5.
 7. Add 700 µl of 1X LiAc/40% PEG-3350/1X TE and mix well.
 8. Incubate solution at 30°C for 30 minutes with shaking.
 9. Add 88 µl DMSO, mix well, and heat shock at 42°C for 7 minutes.
 10. Centrifuge in a microcentrifuge for 10 seconds and remove supernatant.
 11. Resuspend the cell pellet in 1 ml 1X TE and re-pellet.
 12. Resuspend the cell pellet in 50-100 µl 1X TE and plate on a selective plate.
-



Note

To calculate the number of yeast cells, assume that 1 OD₆₀₀ unit = ~2.0 x 10⁷ yeast cells.

Preparation of Denatured Salmon Sperm DNA

Introduction

A convenient protocol to make denatured salmon sperm DNA (Schiestl and Gietz, 1989) is provided for your convenience. You may also purchase denatured salmon sperm DNA from Sigma-Aldrich (Catalog no. D9156). Alternatively, some researchers have found that using yeast transfer RNA (Sigma-Aldrich, Catalog no. R9001) as a carrier results in a cleaner transformation although there are fewer total colonies.

Materials Needed

Prepare or have the following reagents on hand before starting.

- Salmon Sperm DNA (Sigma-Aldrich, Catalog no. D1626)
 - 1X TE
 - Sonicator
 - 50 ml conical centrifuge tubes
 - TE-saturated phenol
 - TE-saturated phenol:chloroform:isoamyl alcohol (25:24:1)
 - Chloroform
 - Low-speed centrifuge
 - 3 M sodium acetate, pH 6.0
 - 95% ethanol (-20°C)
 - 250 ml centrifuge bottle
 - Boiling water bath
-

Protocol

1. In a 250 ml flask, dissolve 1 g salmon sperm DNA into 100 ml of TE (10 mg/ml). Pipet up and down with a 10 ml pipet to dissolve completely.
 2. Incubate overnight at +4°C on a rotating wheel.
 3. Using a sonicator with a large probe, sonicate the DNA twice for 30 seconds at 3/4 power. The resulting DNA will have an average size of 7 kb. You may verify the size of the DNA on a gel.
 4. Aliquot the sonicated DNA into four 50 ml conical centrifuge tubes (25 ml per tube).
 5. Extract with 25 ml of TE-saturated phenol. Centrifuge at 10,000 x g for 5 minutes at +4°C. Transfer the DNA (upper layer) to a fresh 50 ml conical centrifuge tube.
 6. Extract with 25 ml of TE-saturated pheno:chloroform:isoamyl alcohol (25:24:1). Centrifuge at 10,000 x g for 5 minutes at +4°C. Transfer the DNA (upper layer) to a fresh 50 ml conical centrifuge tube.
 7. Extract with 25 ml of chloroform. Centrifuge at 10,000 x g for 5 minutes at +4°C. Transfer the DNA (upper layer) to a 250 ml centrifuge bottle.
 8. Add 5 ml of 3 M sodium acetate, pH 6.0 (1/10 volume) and 125 ml ice-cold (-20°C) 95% ethanol (2.5 volume) to precipitate DNA.
 9. Pellet the DNA at 12,000 x g for 15 minutes at +4°C.
 10. Wash the DNA once with 200 ml 70% ethanol and centrifuge as described in step 9.
 11. Partially dry DNA by air or in a Speed-Vac (cover tubes with parafilm and poke holes in top) for 20 minutes.
 12. Transfer DNA to a 250 ml sterile flask and dissolve DNA in 100 ml sterile TE (10 mg/ml).
 13. Boil for 20 minutes to denature DNA. Immediately place on ice, aliquot in 1 ml samples, and freeze at -20°C.
-

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):

www.invitrogen.com

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

Contact Us

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

Corporate Headquarters:

Invitrogen Corporation
1600 Faraday Avenue
Carlsbad, CA 92008
USA

Tel: 1 760 603 7200

Tel (Toll Free): 1 800 955 6288

Fax: 1 760 602 6500

E-mail: tech_service@invitrogen.com

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, please visit our Web site (www.invitrogen.com) and follow the instructions below.

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. Your MSDS will be sent within 24 hours.
-

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

**Limited Use Label
License No. 57:
Dual-Baid Hybrid
Hunter™
Technology**

Practice of “Improvements to the Yeast Interaction Trap” is covered by U.S. Patent No. 6,326,150 assigned to the Fox Chase Cancer Center (FCCC). Purchase of this product does not imply or convey a license to practice the “Improvements” covered by this patent application, beyond use of the enclosed kit for non-commercial research. If you are a commercial entity, your right from FCCC to use the System expires after one year. Any commercial entity that wishes to use the product beyond this one-year period must obtain a commercial license from FCCC. Commercial entities will be contacted during this one-year period by FCCC regarding their desire to obtain a commercial license. For license information, please contact: Fox Chase Cancer Center, Office of Business Development, 7701 Burholme Avenue, Philadelphia, PA 19111, Phone: 215-728-1113; Fax:215-728-2594.

**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, sublicense, 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. 76:
Prokaryotic
Control Elements
in Eukaryotes**

This product is licensed under U.S. Patent No. 4,833,080. It may not be used for gene expression in plants nor certain molecular modifications and screening. Inquiries about licensing for commercial manufacture or use outside the research kits and reagents market should be directed to: Office of Technology and Trademark Licensing, Harvard University, University Place, Suite 410 South, 124 Mt. Auburn Street, Cambridge, MA 02138, Phone: 617-495-3067, Fax: 617-495-9568.

Inquiries regarding commercial use in the research market should be directed to the Licensing Department, Invitrogen Corporation, 1600 Faraday Avenue, Carlsbad, CA 92008.

Continued on next page

Purchaser Notification, continued

**Limited Use Label
License No. 122:
Interaction Trap
System**

This product is subject to U.S. Patent No. 5,580,736 owned by Massachusetts General Hospital and licensed to Invitrogen for research and may not be administered to humans or animals or used for any therapeutic or diagnostic purposes.

**Limited Use Label
License No. 141:
Expression of
Polypeptides 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.

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).
- Bartel, P. L., Chien, C.-T., Sternglanz, R., and Fields, S. (1993a) Using the Two-Hybrid System to Detect Protein-Protein Interactions. In *Cellular Interactions in Development: A Practical Approach*, D. A. Hartley, ed. (Oxford: Oxford University Press), pp. 153-179.
- Fields, S., and Sternglanz, R. (1994). The Two-Hybrid System: An Assay For Protein-Protein Interactions. *Trends Genet.* *10*, 286-292.
- Gietz, D., Jean, A. S., Woods, R. A., and Schiestl, R. H. (1992). Improved Method for High-Efficiency Transformation of Intact Yeast Cells. *Nuc. Acids Res.* *20*, 1425.
- Gietz, R. D., and Schiestl, R. H. (1996) Transformation of Lithium-Treated Yeast Cells and the Selection of Auxotrophic and Dominant Markers. In *Methods in Molecular Biology*, I. H. Evans, ed. (Totowa, NJ: Humana Press).
- Golemis, E. A., Gyuris, J., and Brent, R. (1996) Interaction Trap/Two-Hybrid Systems to Identify Interacting Proteins. In *Current Protocols in Molecular Biology*, F. M. Ausubel, R. Brent, R. E. Kingston, D. D. Moore, J. G. Seidman, J. A. Smith and K. Struhl, eds. (New York: Greene Publishing Associates and Wiley-Interscience), pp. 20.1.1.-20.1.28.
- Golemis, E. A., Serebriiskii, I., R.L. Finley, J., Kolonin, M. G., Gyuris, J., and Brent, R. (1998) Interaction Trap/Two-Hybrid System to Identify Interacting Proteins. In *Current Protocols in Protein Science*, J. E. Coligan, B. M. Dunn, H. L. Ploegh, D. W. Speicher and P. T. Wingfield, eds. (New York: John Wiley), pp. 19.2.1-19.2.40.
- Hill, J., Donald, K. A., and Griffiths, D. E. (1991). DMSO-Enhanced Whole Cell Yeast Transformation. *Nuc. Acids Res.* *19*, 5791.
- Horii, T., Ogawa, T., and Ogawa, H. (1981). Nucleotide Sequence of the *lexA* Gene of *E. coli*. *Cell* *23*, 689-697.
- Ma, J., and Ptashne, M. (1987). A New Class of Yeast Transcriptional Activators. *Cell* *51*, 113-119.
- Markham, B. E., Little, J. W., and Mount, D. W. (1981). Nucleotide Sequence of the *lexA* Gene of *Escherichia coli* K-12. *Nuc. Acids Res.* *9*, 4149-4161.
- Nilsson, B., Uhlen, M., Josephson, S., Gatenbeck, S., and Philipson, L. (1983). An Improved Positive Selection Plasmid Vector Constructed by Oligonucleotide Mediated Mutagenesis. *Nuc. Acids Res.* *11*, 8019-8030.
- Ruden, D. M., Ma, J., Li, Y., Wood, K., and Ptashne, M. (1991). Generating Yeast Transcriptional Activators Containing No Yeast Protein Sequences. *Nature* *350*, 250-252.
- Sambrook, J., Fritsch, E. F., and Maniatis, T. (1989). *Molecular Cloning: A Laboratory Manual*, Second Edition (Plainview, New York: Cold Spring Harbor Laboratory Press).
- Schiestl, R. H., and Gietz, R. D. (1989). High Efficiency Transformation of Intact Cells Using Single Stranded Nucleic Acids as a Carrier. *Curr. Genet.* *16*, 339-346.
- Southern, J. A., Young, D. F., Heaney, F., Baumgartner, W., and Randall, R. E. (1991). Identification of an Epitope on the P and V Proteins of Simian Virus 5 That Distinguishes Between Two Isolates with Different Biological Characteristics. *J. Gen. Virol.* *72*, 1551-1557.
- Tschumper, G., and Carbon, J. (1980). Sequence of a Yeast DNA Fragment Containing a Chromosomal Replicator and the *TRP1* Gene. *Gene* *10*, 157-166.