

# Matchmaker™ Mammalian Assay Kit 2 User Manual



Clontech

**United States/Canada**  
800.662.2566

**Asia Pacific**  
+1.650.919.7300

**Europe**  
+33.(0)1.3904.6880

**Japan**  
+81.(0)77543.6116

**Cat. No. 630305  
(122811)**

Clontech Laboratories, Inc.  
A Takara Bio Company  
1290 Terra Bella Ave.  
Mountain View, CA 94043  
Technical Support (US)  
E-mail: [tech@clontech.com](mailto:tech@clontech.com)  
[www.clontech.com](http://www.clontech.com)

---

## Table of Contents

---

<b>I. Introduction</b>	<b>3</b>
<b>II. List of Components</b>	<b>6</b>
<b>III. Additional Materials Required</b>	<b>7</b>
<b>IV. Matchmaker Mammalian Assay Kit 2 Protocols</b>	<b>8</b>
A. Transformation of Plasmids into <i>E. coli</i> and Plasmid Isolation	8
B. Construction of Plasmids Expressing Hybrid Proteins	8
C. Guidelines for Transfection of Mammalian Cells	9
D. Positive Control Experiment	11
E. Recommendations for Performing SEAP Assays	11
F. Further Experiments	12
<b>V. References</b>	<b>13</b>
<b>Appendix A: SEAP Protocol</b>	<b>14</b>
<b>Appendix B: Plasmid Maps &amp; Multiple Cloning Sites</b>	<b>16</b>

### List of Tables

<b>Table I.</b> Examples of Protein-Protein Interactions Characterized with Mammalian Two-Hybrid Assays	5
<b>Table II.</b> Recommended Set-up for Mammalian Two-Hybrid Assays	10
<b>Table III.</b> Positive Control Mammalian Two-Hybrid Experiment	11

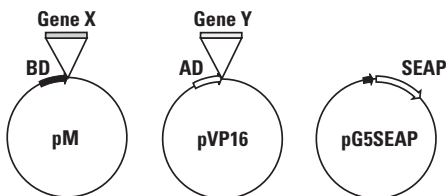
### List of Figures

<b>Figure 1.</b> Testing protein-protein interactions with the Matchmaker Mammalian Assay Kit 2	3
<b>Figure 2.</b> Map of pM GAL4 DNA-Cloning Vector and MCS	16
<b>Figure 3.</b> Map of pVP16 AD Cloning Vector and MCS	17
<b>Figure 4.</b> Map of pG5SEAP Reporter Vector	18

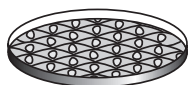
## I. Introduction

The Matchmaker Mammalian Assay Kit 2 provides a complete set of vectors for performing two-hybrid assays in mammalian cells. Two-hybrid assays are based on the fact that many eukaryotic transcriptional activators consist of two physically and functionally separable domains: a DNA-binding domain (DNA-BD) that specifically binds to a promoter or other *cis*-regulatory element, and an activation domain (AD) that directs RNA polymerase II to transcribe the gene downstream of the DNA-binding site. While these domains may be part of the same protein (as in the case of the native yeast GAL4 protein), they can also function as two separate proteins—as long as the AD is tethered to a DNA-BD bound to the promoter. In two-hybrid assays, that tether is the interaction between two additional proteins (X and Y) that are expressed as protein fusions to the AD and DNA-BD peptides, respectively.

- Insert genes for possible interacting proteins X & Y into pM & pVP16



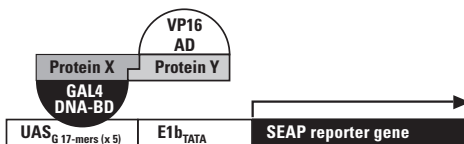
- Cotransfect mammalian host cells



- Incubate 48–72 hr
- Assay SEAP protein levels or activity

### Possible Results

- **Interaction between X & Y**— results in SEAP gene expression



- **No interaction between X & Y**— results in no significant SEAP gene expression

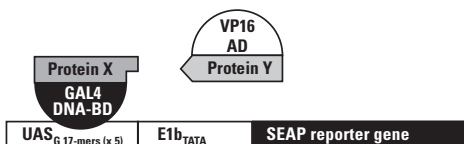


Figure 1. Testing protein-protein interactions with the Matchmaker Mammalian Assay Kit 2.

## I. Introduction *continued*

---

The Matchmaker Mammalian Assay Kit 2 is an improved two-hybrid assay using secreted alkaline phosphatase (SEAP) reporter, for testing protein-protein interactions in mammalian cells in addition to mapping interacting domains (*Clontechiques* 1996). The assay is an important follow-up to yeast screens because it tests interactions under conditions that allow for post-translational changes to hybrid proteins (i.e., phosphorylation, acetylation, proteolysis that cannot be replicated in yeast). Mammalian two-hybrid assays are often used to confirm protein-protein interactions identified through two-hybrid screening in yeast. SEAP, an enzyme that can be assayed simply by sampling the culture medium, avoids the need for cell lysis. If desired, sample collection can be automated using cultures grown in 96-well plates. The use of chemiluminescent substrates in this kit makes the assay extremely sensitive, with relatively few manipulations required, in order to obtain results.

In the Matchmaker Mammalian Assay 2 (Figure 1), the pM GAL4-BD Cloning Vector is used to generate fusions of some protein X to the GAL4 DNA-BD. Similarly, pVP16 AD Cloning Vector is used to construct fusions of some protein Y to an AD derived from the VP16 protein of herpes simplex virus. pG5SEAP is a reporter vector which contains the secreted alkaline phosphatase (SEAP) reporter gene downstream and under the control of a GAL4-responsive element in the form of five consensus GAL4 binding sites and the minimal promoter of the adenovirus E1b gene (Figure 1). Once you have constructed your two hybrid plasmids, all three plasmids are cotransfected into a suitable mammalian host cell line using standard methods. 48–72 hr later, the interaction between proteins X and Y is assayed by measuring SEAP gene expression. In the absence of activation from the GAL4 sites (i.e., no protein-protein interaction), the minimal E1b promoter will not express significant levels of SEAP, resulting in low background. To measure SEAP activity we recommend using the Great EscAPe™ SEAP Chemiluminescence Detection Kit (Cat. No. 631701), a kit for the detection of secreted alkaline phosphatase (SEAP) using the chemiluminescent substrate CSPD.

Once an interaction between two proteins has been detected in mammalian cells, the Matchmaker Mammalian Assay Kit 2 is a powerful tool for functionally analyzing that interaction using deletional or site-directed mutagenesis (Table I). This assay is also useful for confirming the relevance of protein-protein interactions identified via yeast two-hybrid screens using yeast-based two-hybrid systems such as our Matchmaker Two-Hybrid Systems and cDNA and Genomic Libraries. Such confirmation eliminates the possibility of a false positive that is an artifact of working in yeast cells. For more information on yeast two-hybrid technology, see Fields & Sternglanz (1994), Bartel *et al.* (1993) or the Matchmaker Two-Hybrid System 3 User Manual (PT3247-1) available at [www.clontech.com/matchmaker](http://www.clontech.com/matchmaker).

## I. Introduction *continued*

**TABLE I. EXAMPLES OF PROTEIN-PROTEIN INTERACTIONS CHARACTERIZED WITH MAMMALIAN TWO-HYBRID ASSAYS\***

<b>GAL4 DNA-BD Hybrid</b>	<b>VP16 AD Hybrid</b>	<b>Cell Type</b>	<b>Reference</b>
E2F-1	Rb	Saos-2	Fagan, R., <i>et al.</i> (1994) <i>Cell</i> <b>78</b> :799–811
E2F-1	Adenovirus E4 protein	Saos-2	Fagan, R., <i>et al.</i> (1994) <i>Cell</i> <b>78</b> :799–811
FOS	JUN	CHO	Dang, C. V., <i>et al.</i> (1991) <i>Mol. Cell. Biol.</i> <b>11</b> :954–962
LCK	CD4	CHO	Fearon, E. R., <i>et al.</i> (1992) <i>Proc. Natl. Acad. Sci. USA</i> <b>89</b> :7958–7962
R6-K	R6-K	CV1	Vasavada, H. A., <i>et al.</i> (1991) <i>Proc. Natl. Acad. Sci. USA</i> <b>88</b> :10686–10690
TAL1	E47	Jurkat	Hsu, H., <i>et al.</i> (1994) <i>Proc. Natl. Acad. Sci. USA</i> <b>91</b> :3181–3185
VSV protein N	VSV protein PI	CHO	Takacs, A. M., <i>et al.</i> (1993) <i>Proc. Natl. Acad. Sci. USA</i> <b>90</b> :10375–10379

\* Detected with previous version of the Matchmaker Mammalian Two-Hybrid Assay Kit (Cat. No. 630301); CAT reporter used

## II. List of Components

---

**Store all components at –20°C.** Refer to Appendix B for maps and detailed descriptions of the cloning and reporter vectors.

- 20 µl **pM GAL4 DNA-BD Cloning Vector** (500 ng/µl)  
3.5 kb DNA-binding domain cloning vector used to express a fusion of a test protein with the GAL4 DNA-BD
- 20 µl **pVP16 AD Cloning Vector** (500 ng/µl)  
3.3 kb activation domain cloning vector used to express a fusion of a test protein with VP16 AD, a herpes virus protein that acts as a transcriptional activator in mammalian cells
- 20 µl **pG5SEAP Reporter Vector** (500 ng/µl)  
4.5 kb SEAP mammalian reporter plasmid for cotransfection into mammalian cells with recombinant plasmids derived from pM and pVP16
- 20 µl **pM3-VP16 Positive Control Vector** (500 ng/µl)  
4.4 kb positive control plasmid that expresses a fusion of the GAL4 DNA-BD to the VP16 AD
- 20 µl **pM-53 Positive Control Vector** (500 ng/µl)  
4.6 kb positive control plasmid that expresses a fusion of the GAL4 DNA-BD to the mouse p53 protein
- 20 µl **pVP16-T Positive Control Vector** (500 ng/µl)  
5.3 kb positive control plasmid that expresses a fusion of the VP16 AD to the SV40 large T-antigen, which is known to interact with p53
- 20 µl **pVP16-CP Negative Control Vector** (500 ng/µl)  
4.5 kb negative control plasmid that expresses a fusion of the VP16AD to a viral coat protein, which does not interact with p53

### III. Additional Materials Required

The following reagents are required but not supplied.

#### Mammalian Cell Culture and Harvesting

- **Cell culture medium** (e.g., Dulbecco's Modified Eagle Medium [DMEM, Sigma Cat. No. D5671] or another appropriate growth medium for mammalian cells in culture)
- **Fetal bovine serum, newborn calf serum, or equivalent** (to supplement growth medium; see Related Products, Sec. VI)
- **Phosphate buffered saline (PBS) (pH 7.4)**

<u>Component</u>	<u>Final Conc.</u>	<u>Amount for 2 L</u>
Na <sub>2</sub> HPO <sub>4</sub>	58 mM	16.5 g
NaH <sub>2</sub> PO <sub>4</sub>	17 mM	4.1 g
NaCl	68 mM	8.0 g

Dissolve the above components in 1.8 L of deionized H<sub>2</sub>O. Adjust to pH 7.4 with 0.1N NaOH. Add deionized H<sub>2</sub>O to final volume of 2 L. Store at room temperature.

- **1X Trypsin/EDTA**

#### Mammalian Cell Transfection

Reagents will depend on which transfection method you use. We generally use calcium-phosphate-mediated transfection for Matchmaker Mammalian Two-Hybrid experiments and recommend using the CalPhos Mammalian Transfection Kit (Cat. No. 631312). For liposome-mediated transfection, we recommend CLONfectin Transfection Reagent (Cat. No. 631301). See Appendix A for recipes and protocols. Protocols for other methods can be found in Freshney (1993) and Ausubel *et al.* (1994).

#### SEAP Assay

There are several nonradioactive alternatives that are easy to perform and provide quantitative data. At Clontech, we generally measure SEAP protein directly in Matchmaker Mammalian Two-Hybrid experiments using the Great EscAPe SEAP Chemiluminescence Detection Kit (Cat. Nos. 631701 & 631702), a simple, nonradioactive assay for the detection of secreted alkaline phosphatase (SEAP) using the chemiluminescent substrate CSPD.

## IV. Matchmaker Mammalian Assay Kit 2 Protocols

---

PLEASE READ THROUGH THE ENTIRE PROTOCOL BEFORE BEGINNING.

### A. Transformation of Plasmids into *E. coli* and Plasmid Isolation

1. Transform each of the plasmids provided in this kit into a suitable *E. coli* host strain such as Stellar™ Competent Cells (Cat. No. 636763) to ensure that you have a renewable source of DNA.
2. You will need to perform large-scale plasmid preparations of any plasmid that will be introduced into mammalian cells (e.g., your hybrid gene constructs derived from pM and pVP16, pG5SEAP, and the control plasmids). To ensure the purity of the DNA, isolate all plasmids for transfection by banding on CsCl gradient (Sambrook & Russell, 2001), or by an equivalent method. For optimal results, you may wish to purify twice on CsCl gradients. We recommend Nucleo-Bond Plasmid Maxi Kit (Cat. Nos. 635933 & 635934) for this purpose.

### B. Construction of Plasmids Expressing Hybrid Proteins

Construct fusion genes using standard molecular biology techniques. A brief outline of the protocol is given below (for more detailed information, see Sambrook & Russell, 2001). The gene of test protein X is fused to the GAL4 DNA-BD in the pM vector. The gene of test protein Y is fused to the VP16 AD in the pVP16 vector. The orientation and reading frame of both fusions must be correct for hybrid proteins to be expressed.

In some cases, the fusion gene can be generated using compatible restriction sites that are present in the test gene and the cloning vector. Many of the sites in pM are in the same reading frame as in the yeast two-hybrid vectors pGBKT7 (see map, Appendix B). If no such sites are present, the gene fragment can be generated by PCR with useful restriction sites incorporated into the primers (Scharf, 1990). Often a restriction site at the end of the gene of interest can be changed into a different site or put into a different reading frame by using a PCR primer which incorporates the desired restriction site at the desired place.

1. Purify the gene fragment, whether generated by PCR or cut out of a plasmid, by any standard method (Sambrook & Russell, 2001).
2. Digest pM (or pVP16) with the appropriate restriction enzyme(s), treat with phosphatase, and purify.
3. Ligate pM (or pVP16) and the insert encoding protein X (or protein Y).
4. Transform ligation mixtures into *E. coli* to maintain a working stock of the recombinant plasmid.
5. Identify insert-containing plasmid by restriction analysis.
6. Check orientation and reading frame by sequencing fusion junctions.



---

## IV. Matchmaker Mammalian Assay Kit 2 *continued*

---

### C. Guidelines for Transfection of Mammalian Cells

Plasmids may be cotransfected into mammalian cells by any standard transfection method. At Clontech, we typically use calcium phosphate (Chen & Okayama, 1988) and recommend the CalPhos Mammalian Transfection Kit (Cat. No. 631312) for mammalian transfections. A protocol is provided in Appendix A. For liposome-mediated transfection, we recommend CLONfectin Transfection Reagent (Cat. No. 631301). DEAE-dextran (Rosenthal, 1987) also works well in our hands. Useful general references on cell culture techniques are:

- *Current Protocols in Molecular Biology*, ed. by F. M. Ausubel *et al.* (1994, Greene Publishing Associates and Wiley & Sons)

The efficiency of transfection for different cell lines may vary by several orders of magnitude. A method that works well for one host cell line may be inferior for another. Therefore, when working with a cell line for the first time, you may want to compare the efficiencies of several transfection protocols. This can be done by cotransfecting the pG5SEAP and pM3-VP16 plasmids and assaying for SEAP activity.

After a method of transfection is chosen, it may be necessary to optimize parameters such as cell density, the amount and purity of the DNA, media conditions, and transfection time. Once optimized, these parameters should be kept constant to obtain reproducible results. With each method, SEAP activity may be detected 48–72 hrs after transfection, depending on the host cell line used.

For rapid, qualitative results, we generally obtain reproducible results using single transfections. For quantitative data, we recommend that you perform duplicate or triplicate transfections and average the results. You may also wish to normalize for transfection efficiency by cotransfecting a constant amount of a second reporter under the control of a constitutive promoter. The values obtained in each sample for the primary reporter (i.e., SEAP) are then normalized to the values obtained for the second reporter in the same sample. Useful vectors for this purpose are Clontech pCMV $\beta$  (Cat. No. 631719), p $\beta$ gal-Control (Cat. No. 631709), or pSEAP2-Control (Cat. No. 631717).

## IV. Matchmaker Mammalian Assay Kit 2 *continued*

Table II describes the basic set-up for Matchmaker Mammalian Assay Kit 2 experiments, including positive and negative controls.

**TABLE II. RECOMMENDED SET-UP FOR MAMMALIAN TWO-HYBRID ASSAYS**

<b>Transfection</b>	<b>GAL4 DNA-BD Plasmid (10 µg<sup>a</sup>)</b>	<b>VP16 AD Plasmid (10 µg<sup>a</sup>)</b>	<b>Reporter Plasmid (2 µg<sup>a</sup>)</b>	<b>SEAP Protein or SEAP Activity</b>
1 (experiment)	pM-insert X	pVP16-insertY	pG5SEAP	to be determined
2 (untransfected control)	none	none	none	– <sup>b</sup>
3 (basal control)	pM	pVP16	pG5SEAP	– <sup>c</sup>
4 (X control)	pM-insert X	pVP16	pG5SEAP	to be determined <sup>d</sup>
5 (Y control)	pM	pVP16-insertY	pG5SEAP	to be determined <sup>e</sup>
6 (positive control)	pM3-VP16 <sup>f</sup>	(pM3-VP16)	pG5SEAP	+++

<sup>a</sup> These amounts of DNA are optimized for calcium-phosphate mediated transfection of HeLa cells. When using other transfection methods, adjust the total amount of DNA accordingly, but maintain the ratio of the different plasmids. The optimal amount of DNA may also vary with different cell types.

<sup>b</sup> This value is the background SEAP signal in your cells. For quantitative SEAP assays, subtract this value from all other experimental values.

<sup>c</sup> This control (or a control using just pM and pG5SEAP) provides the basal expression level of SEAP protein or SEAP activity in your experiments. To determine the fold-induction caused by an interaction between protein X and protein Y, divide your experimental SEAP values by this number.

<sup>d</sup> Critical control to determine whether or not your protein X functions autonomously as a transcriptional activator. The level of SEAP should be similar to the background value obtained in transfection 3.

<sup>e</sup> Critical control to determine whether or not your protein Y functions autonomously as a DNA-BD or binds directly to the DNA-BD encoded by pM. The level of SEAP should be similar to the background value obtained in transfection 3.

<sup>f</sup> Use 20 µg when transfecting *only* a DNA-BD or AD vector with pG5SEAP. The pM3-VP16 positive control plasmid encodes a fusion of the GAL4 DNA-BD and the VP16 AD and therefore gives very strong SEAP expression when cotransfected with pG5SEAP.

## IV. Matchmaker Mammalian Assay Kit 2 *continued*

### D. Controls

Perform the control experiments described in Table III to confirm that the mammalian two-hybrid assay works in your hands and in your cells, and to optimize your transfection and SEAP assay protocols. These experiments compare the interaction between the p53 protein and two other proteins: the SV40 large T-antigen, which is known to interact with p53; and a polyoma virus coat protein (CP), which does not interact with p53. All the plasmids in Table III are included in this kit.

TABLE III. CONTROLS FOR MAMMALIAN TWO-HYBRID EXPERIMENT

	Transfection Control	GAL4 DNA-BD Vector (10 µg <sup>a</sup> )	VP16 AD Vector (10 µg <sup>a</sup> )	Reporter Vector (2 µg <sup>a</sup> )	SEAP Protein or SEAP Activity
1	Positive	pM-53	pVP16-T	pG5SEAP	+
2	Negative	pM-53	pVP16-CP	pG5SEAP	–
3	Untransfected	none	none	none	– <sup>b</sup>
4	Basal	pM	pVP16	pG5SEAP	– <sup>c</sup>
5	Negative	pM-53	pVP16	pG5SEAP	– <sup>d</sup>
6	Negative	pM	pVP16-T	pG5SEAP	– <sup>e</sup>
7	Positive	pM3-VP16 <sup>f</sup>	(pM3-VP16)	pG5SEAP	++

<sup>a-f</sup> See Table II (previous page).

### E. Recommendations for Performing SEAP Assays

At Clontech, we generally measure SEAP protein directly in Matchmaker Mammalian Assay Kit 2 experiments using the Great EscAPe SEAP Chemiluminescence Detection Kit (Cat. Nos. 631701 & 631702), a simple, nonradioactive assay for the detection of secreted alkaline phosphatase (SEAP) using the chemiluminescent substrate CSPD.

## IV. Matchmaker Mammalian Assay Kit 2 *continued*

---

### F. Further Experiments

#### 1. Verification of Hybrid Protein Expression [Optional]

Expression of the GAL4 DNA-BD/Protein X fusion in mammalian cells can be verified by preparing Western blots using soluble protein extracts from the cells. Probe the blots with the GAL4 DNA-BD Monoclonal Antibody (Cat. No. 630403) using standard Western blotting procedures (Harlow & Lane, 1988; Sambrook & Russell, 2001).

#### 2. Mapping Protein Structure

Once an interaction between two proteins has been detected in mammalian cells, the Matchmaker Mammalian Assay Kit 2 provides a powerful tool for functionally dissecting that interaction using deletional or site-directed mutagenesis. Transformer Site-Directed Mutagenesis Kit (Cat. No. 630702) is useful for generating mutant proteins which can then be tested in the Matchmaker Mammalian Assay Kit 2.

#### 3. Confirmation in Yeast Matchmaker Two-Hybrid System

Interactions that are detected in the Matchmaker Mammalian Two-Hybrid Assay may also be detectable in a yeast-based two-hybrid assay such as Matchmaker Two-Hybrid System 3 (Cat. No. 630303). Much less DNA is required for each two-hybrid experiment in yeast. Furthermore, yeast two-hybrid assays are much less expensive, since they do not require tissue-culture facilities and supplies for the growth and transfection of mammalian cells. For more information on yeast two-hybrid technology, see the references by Fields & Sternglanz (1994) and Bartel *et al.* (1993) or the references in the Matchmaker Two-Hybrid System 3 User Manual (PT3247-1) available at [www.clontech.com/matchmaker](http://www.clontech.com/matchmaker).

## V. References

---

- Ausubel, F. M., Brent, R., Kingston, R. E., Moore, D. M., Seidman, J. G., Smith, J. A., & Struhl, K. (1994) *Current Protocols in Molecular Biology*. (Greene Publishing Associates, Inc. & John Wiley & Sons, Inc.).
- Bartel, P. L., Chien, C.-T., Sternglanz, R. & Fields, S. (1993) Elimination of false positives that arise in using the two-hybrid system. *BioTechniques* **14**:920–924.
- Chen, C. & Okayama, H. (1988) Calcium phosphate-mediated gene transfer: A highly efficient transfection system for stably transforming cells with plasmid DNA. *BioTechniques* **6**:632.
- Dang, C. V., Barrett, J., Villa-Garcia, M., Resar, L. M. S., Kato, G. J. & Fearon, E. R. (1991) Intracellular leucine zipper interactions suggest c-myc hetero-oligomerization. *Mol. Cell. Biol.* **11**:954–962.
- Fagan, R., Flint, K. J. & Jones, N (1994) Phosphorylation of E2F-1 modulates its interaction with the retinoblastoma gene product and the adenoviral E4 19-kDa protein. *Cell* **78**:799–811.
- Fearon, E. R., Finkel, T., Gillison, M. L., Kennedy, S. P., Casella, J. F., Tomaselli, G. F., Morrow, J. S. & Ding, C. V. (1992) Karyoplasmic interaction selection strategy: a general strategy to detect protein-protein interaction in mammalian cells. *Proc. Natl. Acad. Sci. USA* **89**:7958–7962.
- Fields, S. (1993) The two-hybrid system to detect protein-protein interactions. *Methods: A Companion to Meth. Enzymol.* **5**:116–124.
- Fields, S. & Sternglanz, R. (1994) The two-hybrid system: an assay for protein-protein interactions. *Trends Genet.* **10**: 286–292.
- Freshney, R. I. (1993) *Culture of Animal Cells*, Third Edition (Wiley-Liss, NY).
- Harlow, E. & Lane, E. (1988) *Antibodies: A Laboratory Manual* (Cold Spring Harbor Laboratory, Cold Spring Harbor, NY).
- Hsu, H., Wadman, I. & Baer, R. (1994) Formation of *in vivo* complexes between the TAL1 and E2A polypeptides of leukemic T-cells. *Proc. Natl. Acad. Sci. USA* **91**:3181–3185.
- Sadowski, I., Bell, B., Broad, P. & Hollis, M. (1992) GAL4 fusion vectors for expression in yeast or mammalian cells. *Gene* **118**:137–141.
- Sambrook, J. & Russell, D. W. (2001) *Molecular Cloning: A Laboratory Manual, Third Edition* (Cold Spring Harbor Laboratory Press, Cold Spring Harbor, NY).
- Scharf, S. J. (1990) Cloning with PCR. In *PCR Protocols: A Guide to Methods and Applications*, eds. Innis, M. A., Gelfand, D. H., Sninsky, J. J. & White, T. J. (Academic Press, Inc., San Diego), pp. 84–91.
- Silver, P. A., Keegan, L. P. & Ptashne, M. (1984) Amino terminus of the yeast GAL4 gene product is sufficient for nuclear localization. *Proc. Natl. Acad. Sci. USA* **81**:5951–5955.
- Takacs, A. M., Das, T. & Banerjee, A. K. (1993) Mapping of interacting domain between the nucleocapsid protein and the phosphoprotein of vesicular stomatitis virus by using a two-hybrid system. *Proc. Natl. Acad. Sci. USA* **90**:10375–10379.
- Vasavada, H. A., Ganguly, S., Germino, F. J., Wang, Z. X. & Weissman, S. M. (1991) A contingent replication assay for the detection of protein-protein interactions in animal cells. *Proc. Natl. Acad. Sci. USA* **88**:10686–10690.
- Matchmaker Mammalian Two-Hybrid Assay Kit (1996) *Clontechiques* **XI**(1):10–12.

## Appendix A: SEAP Protocol

### A. Materials Required

We recommend using the CalPhos Mammalian Transfection Kit (Cat. No. 631312) for mammalian transfection. Follow the protocol in the User Manual provided.

The recipes and calcium phosphate transfection protocol below are adapted from *Current Protocols in Molecular Biology*, Supplement 14, Section 9.1.3.

- **2X BES-Buffered Solution (BBS, pH 6.95)**

<u>Component</u>	<u>Final Conc.</u>
<i>N, N</i> -bis (2-hydroxyethyl)-2-aminoethane-sulfonic acid (BES; CALBIOCHEM Cat. No. 391334)	50 mM
NaCl	280 mM
Na <sub>2</sub> HPO <sub>4</sub>	1.5 mM

*It is critical that the pH of this solution be between pH 6.95 and 6.98.*

We recommend that you check each new batch of 2X BES buffer against a reference stock prepared (and tested) earlier.

Filter sterilize through a 0.45 µm nitrocellulose filter (Nalgene).

Store in aliquots at -20°C (can be frozen and thawed repeatedly).

- **2.5 mM CaCl<sub>2</sub>**

Add 183.7 g of CaCl<sub>2</sub> dihydrate (Sigma; tissue culture grade) to 500 ml of H<sub>2</sub>O.

Filter sterilize through a 0.45 µm nitrocellulose filter (Nalgene).

Store at -20°C in 10 ml aliquots (can be frozen and thawed repeatedly).

### B. Protocol for Calcium Phosphate Transfection

All plasmids should be CsCl-banded and diluted to a concentration of 1.0 µg/µl. Store the DNA solution at 4°C. For initial experiments, we recommend that each transfection be performed with a total of 22 µg of DNA (as shown in Tables II and III). However, the optimal concentration of DNA may vary with different cell types.

#### Day 1

1. For each transfection, seed 5 x 10<sup>5</sup> exponentially growing cells in a 10 cm tissue culture plate in 10 ml of complete medium (with serum). There should be < 10<sup>6</sup> cells/plate (~30% confluency) just prior to transfection. This provides enough surface area on the plate for at least two more doublings.

## Appendix A: SEAP Protocol *continued*

---

### Day 2

2. Two hrs before transfection, replace the medium with fresh medium (with serum) to stimulate cell growth.
3. For each transfection, combine the DNAs in a sterile 1.5 ml microcentrifuge tube.
4. Bring the total volume to 450  $\mu$ l by adding TE or deionized H<sub>2</sub>O.
5. Add 50  $\mu$ l of 2.5 M CaCl<sub>2</sub> and vortex thoroughly. (The final concentration of CaCl<sub>2</sub> is 0.25 M.)
6. Add 500  $\mu$ l of 2X BBS, mix well, and incubate at room temperature for 10–20 min to allow a precipitate to form.
7. Add the calcium phosphate-DNA solution dropwise to the medium while swirling the plate.
8. Incubate overnight in a 37°C, 5% CO<sub>2</sub> incubator.

**Note:** The level of carbon dioxide is critical. We recommend using a Fyrite gas analyzer to measure percent CO<sub>2</sub> prior to incubation.

### Day 3

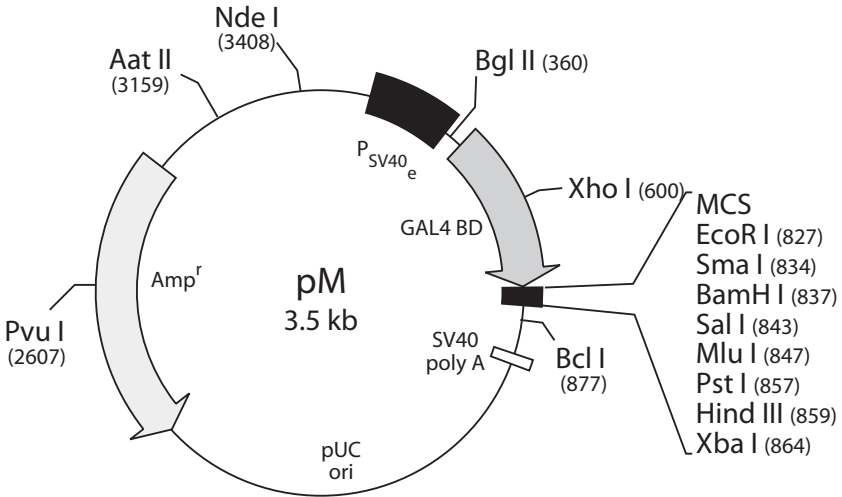
9. Carefully wash the cells twice with 5 ml of PBS, then add 10 ml of complete medium (with serum).
10. Incubate the cells for an additional 24–48 hrs.

**Troubleshooting note:** The amounts of DNA in Tables II and III have been optimized for calcium phosphate transfection of HeLa cells; however, the optimal amount of total DNA may vary with different cell types. (The ratio of input DNAs should stay the same.) To determine the optimum amount of plasmid, transfect three plates of cells with 10, 20, and 30  $\mu$ g of plasmid DNA and incubate overnight. The next day examine the plates with a microscope at 100X. A coarse, clumpy precipitate will form at DNA concentrations that are too low, a fine (almost invisible) precipitate will form at concentrations that are higher than optimal, and an even, granular precipitate will form with optimal DNA concentrations.

### C. SEAP Assay

There are several nonradioactive alternatives that are easy to perform and provide quantitative data. At Clontech, we generally measure SEAP protein directly in Matchmaker Mammalian Assay Kit 2 experiments using the Great EscAPe SEAP Chemiluminescence Detection Kit (Cat. Nos. 631701 & 631702).

## Appendix B: Plasmid Maps & Multiple Cloning Sites



775

•  
 TCA TCG GAA GAG AGT AGT AAC AAA GGT CAA AGA CAG TTG ACT GTA TCG CCG  
 a.a. 147

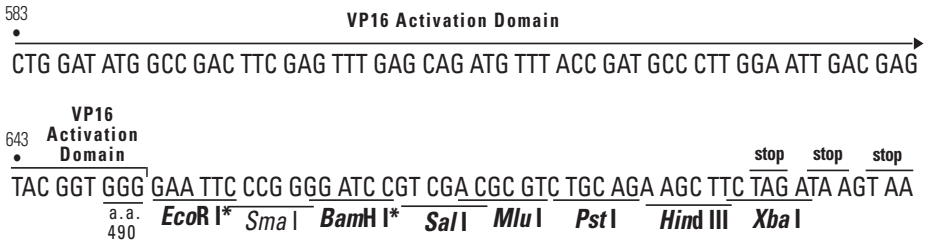
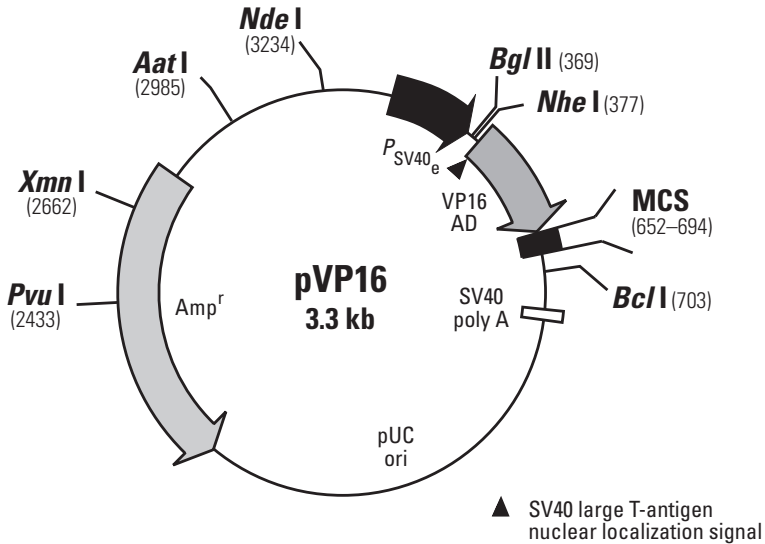
826

•  
 GAA TTC CCG GGG ATC CGT CGA CGC GTC TGC AGA AGC TTC TAG ATA AGT AA  
 EcoR I\* Sma I\* BamH I\* Sal I\* Mlu I Pst I Hind III Xba I

**Figure 2. Map of pM GAL4 DNA-BD Cloning Vector and MCS.** pM is used to generate a fusion of the GAL4 DNA-BD (amino acids 1–147) and a protein of interest. The hybrid protein is targeted to the cell’s nucleus by the GAL4 nuclear localization sequence (Silver *et al.*, 1984). Genes encoding test proteins should be cloned, *in the correct orientation and reading frame*, into one of the unique restriction sites in the MCS region at the 3’ end of the GAL4 DNA-BD. Restriction sites marked with an asterisk (\*) are in the same reading frame as pGBKT7. Transcription is initiated from the constitutive SV40 early promoter ( $P_{SV40e}$ ); transcription is terminated at the SV40 poly A transcription termination signal.

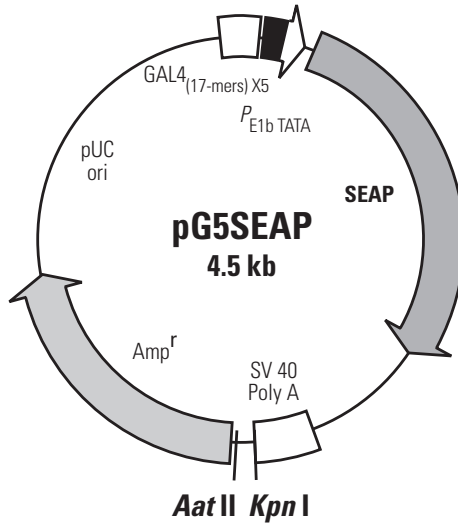


## Appendix B: Plasmid Maps & MCS *continued*



**Figure 3. Map of pVP16 AD Cloning Vector and MCS.** Unique sites are in bold. pVP16 is used to generate a fusion of the VP16 AD (amino acids 446–490) and a protein of interest. The hybrid protein is targeted to the cell's nucleus by the SV40 nuclear localization sequence. Genes encoding test proteins should be cloned, *in the correct orientation and reading frame*, into one of the unique restriction sites in the MCS region at the 3' end of the VP16 AD. Restriction sites marked with an asterisk (\*) are in the same reading frame as pGADT7. Transcription is initiated from the constitutive SV40 early promoter (*P<sub>SV40e</sub>*); transcription is terminated at the SV40 poly A transcription termination signal.

## Appendix B: Plasmid Maps & MCS *continued*



**Figure 4. Map of pG5SEAP Reporter Vector.** pG5SEAP contains five consensus GAL4 binding sites (UAS<sub>G 17-mer (x 5)</sub>) and an adenovirus E1b minimal promoter upstream of the secreted alkaline phosphatase (SEAP) gene. As seen in Figure 1 in the Introduction, the DNA-BD portion of a hybrid protein expressed from a pM-derived plasmid localizes to the GAL4 binding sites in pG5SEAP. If the hybrid test protein X interacts with test protein Y (expressed as a hybrid protein from a pVP16-derived plasmid), the SEAP gene will be transcribed.

## Notes

---

### Notice to Purchaser

Clontech products are to be used for research purposes only. They may not be used for any other purpose, including, but not limited to, use in drugs, in vitro diagnostic purposes, therapeutics, or in humans. Clontech products may not be transferred to third parties, resold, modified for resale, or used to manufacture commercial products or to provide a service to third parties without prior written approval of Clontech Laboratories, Inc.

Your use of this product is also subject to compliance with any applicable licensing requirements described on the product's web page at <http://www.clontech.com>. It is your responsibility to review, understand and adhere to any restrictions imposed by such statements.

Clontech, the Clontech logo, Great EscAPe, Matchmaker, and Stellar are trademarks of Clontech Laboratories, Inc. All other marks are the property of their respective owners. Certain trademarks may not be registered in all jurisdictions. Clontech is a Takara Bio Company. ©2011 Clontech Laboratories, Inc.

This document has been reviewed and approved by the Clontech Quality Assurance Department.