





pFLD and pFLD α

Pichia pastoris expression vectors for inducible expression with methylamine and selection on Zeocin[™]

Catalog number V230-20

Revision date: 29 March 2012 Publication Part number 25-0547

MAN0000304





Contents

Kit Contents and Storage	V
Introduction	1
Product Overview	1
Methods	2
General Cloning Considerations	2
Cloning into pFLD	
Cloning into pFLDα	5
Transforming <i>E. coli</i>	7
Pichia Transformation	9
Expression in <i>Pichia</i>	
Purification	
Appendix	
Recipes	
Zeocin [™]	
Map of pFLD	
Map of pFLDα	
Vector Features	
Map of pFLD/CAT	
Lithium Chloride Transformation Method	
Accessory Products	
Technical Support	
Purchaser Notification	
References	

Kit Contents and Storage

Shipping/Storage	Vectors are shipped at room temperature. Upon receipt, store at -20°C.
Kit Contents	 Catalog number V230-20 includes the following vectors, supplied at a concentration of 0.5 μg/μL in 10 mM Tris-HCl, 1 mM EDTA, pH 8.0 in a total volume of 40 μL: pFLD pFLDα pFLD/CAT (control plasmid)
Reference Sources	The pFLD vectors may be used with the EasySelect [™] <i>Pichia</i> Expression Kit (Catalog no. K1740-01) or the Original <i>Pichia</i> Expression Kit (Catalog no. K1710- 01), available for purchase. Additional general information about recombinant protein expression in <i>Pichia pastoris</i> is provided in the manuals for the EasySelect [™] <i>Pichia</i> Expression Kit and the Original <i>Pichia</i> Expression Kit. The manuals can be downloaded from <u>www.lifetechnologies.com/support</u> or obtained from Technical Support (see page 31). For more information about the EasySelect [™] <i>Pichia</i> Expression Kit or the Original <i>Pichia</i> Expression Kit, refer to <u>www.lifetechnologies.com/support</u> or call Technical Support. More detailed information and protocols dealing with <i>Pichia pastoris</i> may also be found in the following general reference: Higgins, D. R., and Cregg, J. M. (1998) <i>Pichia</i> Protocols. In <i>Methods in Molecular</i> <i>Biology</i> , Vol. 103. (J. M. Walker, ed. Humana Press, Totowa, NJ)
Product Use	For research use only. Not intended for any human or animal therapeutic or diagnostic use.

Kit Contents and Storage, Continued

Recommended <i>Pichia</i> Host Strain	We recommend using the X-33 <i>Pichia</i> strain as the host for expression of recombinant proteins from pFLD. Other <i>Pichia</i> strains including GS115, KM71H, and SMD1168H are suitable. The X-33 <i>Pichia</i> strain and other strains are available for purchase (see page 29 for ordering information). The X-33 <i>Pichia</i> strain has the following genotype and phenotype: Genotype: Wild-type Phenotype: Mut ⁺
Materials Supplied by the User	 Fquipment Microbiological equipment Electroporation device and 0.2 cm cuvettes or reagents for transformation 16°C and 37°C water baths or temperature blocks 30°C and 37°C shaking and non-shaking incubators Hemacytometer Microtiter plates (optional) Reagents <i>Pichia</i> host strain (e.g., X-33, GS115, SMD1168H, KM71H) Electrocompetent or chemically competent <i>E. coli</i> (must be <i>rec</i>A, <i>end</i>A) for transformation Restriction enzymes and appropriate buffers Agarose and low-melt agarose PureLink[®] HQ Mini Plasmid Purification Kit (see page 29 for ordering) or glass milk Sterile water CIAP (calf intestinal alkaline phosphatase, 1 unit/ μL) 10X CIAP Buffer Phenol/chloroform 3 M sodium acetate 100% ethanol 80% ethanol T4 Ligase (2.5 units/ μL) 10X Ligation Buffer (with ATP) Zeocin[™] selection agent (see page 29 for ordering information) YPDS plates containing the appropriate concentration of Zeocin[™] (see page 18 for recipe) 50 mL conical centrifuge tubes 15 mL polypropylene tubes ProBond[™] Purification System (optional, see page 30 for ordering)

Introduction

Product Overview

Introduction

pFLD (4.4 kb) and pFLD α (4.7 kb) are vectors used to express recombinant proteins in *Pichia pastoris*. Recombinant proteins are expressed as fusions to a C-terminal peptide containing the V5 epitope and a polyhistidine (6xHis) tag. The vector allows high-level, methanol- and methylamine-inducible expression of the gene of interest in *Pichia*, and can be used in any *Pichia* strain, including X-33, GS115, SMD1168H, and KM71H. The pFLD vectors contain the following elements:

- 5' fragment containing the *FLD1* promoter for methanol- or methylamineinduced expression of the gene of interest (Shen *et al*, 1998)
- Zeocin[™] resistance gene for selection in both *E. coli* and *Pichia* (Baron *et al.*, 1992; Drocourt *et al.*, 1990)
- Ampicillin resistance gene for selection in E. coli
- C-terminal peptide containing the V5 epitope and a polyhistidine (6xHis) tag for detection and purification of a recombinant fusion protein (if desired)

Experimental Overview

The following table describes the basic steps needed to clone and express your gene of interest in pFLD.

Step	Action	Page
1	Propagate pFLD by transformation into a <i>rec</i> A, <i>end</i> A1 <i>E</i> . <i>coli</i> strain such as TOP10, DH5 α^{TM} , or JM109.	2
2	Develop a cloning strategy and ligate your gene into one of the pFLD vectors in frame with the C-terminal tag.	3–6
3	Transform into <i>E. coli</i> and select transformants on select on LB with $50-100 \ \mu g/mL$ ampicillin (LB-Amp).	7
4	Analyze 10–20 transformants by restriction mapping or sequencing to confirm in-frame fusion of your gene with the C-terminal tag.	8
5	Purify and linearize the recombinant plasmid for transformation into <i>Pichia pastoris</i> .	7–10
6	Transform your <i>Pichia</i> strain and plate onto YPDS plates containing the appropriate concentration of Zeocin [™] .	9–12
7	Select for Zeocin [™] -resistant transformants.	9
8	Optimize expression of your gene.	13–14
9	Purify your fusion protein on metal-chelating resin (<i>i.e.</i> ProBond [™]).	16–17

Methods

General Cloning Considerations

Introduction	The multiple cloning site for pFLD is shown on page 4 and the multiple cloning site for pFLD α is shown on page 6. Use these diagrams to design a strategy to clone your gene of interest in frame with the C-terminal peptide. General considerations for cloning and transformation are discussed in this section.						
General Molecular Biology Techniques	For assistance with <i>E. coli</i> transformations, restriction enzyme analysis, DNA biochemistry, and plasmid preparation, refer to <i>Molecular Cloning: A Laboratory Manual</i> (Sambrook <i>et al.</i> , 1989) or <i>Current Protocols in Molecular Biology</i> (Ausubel <i>et al.</i> , 1994).						
<i>E. coli</i> Strain	Many <i>E. coli</i> strains are suitable for the propagation of the pFLD vectors, including TOP10, JM109, and DH5 α^{TM} . We recommend that you propagate the pFLD vectors in <i>E. coli</i> strains that are recombination deficient (<i>recA</i>) and endonuclease A deficient (<i>endA</i>).						
	For your convenience, TOP10 <i>E. coli</i> are available as chemically competent or electrocompetent cells (see page 29 for ordering).						
Transformation Method	You may use any method of choice for transformation. Chemical transformation is the most convenient for many researchers. Electroporation is the most efficient and the method of choice for large plasmids.						
Maintaining Plasmids	The pFLD vectors contain the ampicillin resistance gene and the Zeocin ^{TT} resistance gene to allow selection of the plasmid in <i>E. coli</i> using either ampicillin or Zeocin ^{TT} . The following procedure uses ampicillin to propagate and maintain pFLD plasmids:						
	1. Use the supplied stock solution to transform a <i>recA</i> , <i>end</i> A <i>E</i> . <i>coli</i> strain like TOP10, DH5α [™] , JM 109, or equivalent.						
	2. Select transformants on LB agar with 50–100 μ g/mL ampicillin (LB-Amp).						
	3. Prepare a glycerol stock from each transformant containing plasmid for long-term storage (see page 7).						

General Cloning Considerations, Continued

General Considerations	The following are some general guidelines when using pFLD or pFLD α to express your gene of interest in <i>Pichia</i> :					
	• The codon usage in <i>Pichia</i> is believed to be similar to <i>Saccharomyces cerevisiae</i> .					
	• Many <i>Saccharomyces</i> genes have proven to be functional in <i>Pichia</i> .					
	• The premature termination of transcripts because of "AT rich regions" has been observed in <i>Pichia</i> and other eukaryotic systems (Henikoff and Cohen, 1984; Irniger <i>et al.</i> , 1991; Scorer <i>et al.</i> , 1993; Zaret and Sherman, 1984). If you have problems expressing your gene, check for premature termination by northern analysis and check your sequence for AT rich regions. It may be necessary to change the sequence in order to express your gene (Scorer <i>et al.</i> , 1993).					
Special	For pFLD only:					
Considerations for pFLD	• Your insert should contain an initiation ATG codon as part of a yeast consensus sequence (Romanos <i>et al.</i> , 1992). An example of a yeast consensus sequence is provided below. The ATG initiation codon is shown underlined.					
	(G/A)NN <u>ATG</u> G					
	• To express your gene as a recombinant fusion protein, you must clone your gene in frame with the C-terminal peptide containing the V5 epitope and the polyhistidine tag. Refer to the diagram on page 4 to develop a cloning strategy.					
	• To express your gene of interest without the C-terminal peptide, make sure that your gene contains a stop codon.					
Special	For pFLDa only:					
Considerations for $pFLD\alpha$	• pFLDα is an N-terminal fusion vector. To express your gene, you must clone your gene in frame with the N-terminal α-factor secretion signal. Refer to the diagram on page 6 to develop a cloning strategy.					
	• To express your gene of interest without the C-terminal peptide, make sure that your gene contains a stop codon.					
	• The predicted protease cleavage sites for the α-factor signal sequence are indicated in the figure on page 6.					

Cloning into pFLD

Multiple Site of p	•	indicate the cleavage sequencing and fur downloading from	n sites are labeled to s been confirmed by e of pFLD is available for Technical Support (see f pFLD, see page 24.		
		3´ end of FLD1 Mfe I		Sfil	Asp718 Kpn Xho
581	AATTCTTGA	AT ATTCACACAA TI	IGTTCACGT GGC	CCAGCCG GCCGTC	TCGG ATCGGTACCT
	Sac	• H	Apal Mfel		V5 epitope
641	CGAGCCGCG	GG CGGCCGCCAG CI	TTGGGCCCA ATT		CCT ATC CCT AAC Pro Ile Pro Asn
					Polyhistidine (6xHis) tag
697		CTC GGT CTC GAT Leu Gly Leu Asp		GTC GAC CAT CA Val Asp His Hi	T CAT CAT CAT CAT S His His His His
748	TGA GTTTO ***	GTAGCC TTAGACATO	GA CTGTTCCTCA	GTTCAAGTTG GG	CACTTACG AGAAGACCGG
			3´ <i>AOX1</i> pr	iming site	
811	TCTTGCTAG	ЗА ТТСТААТСАА GA	AGGATGTCA GAA	TGCCATT TGCCTG	AGAG ATGCAGGCTT

Cloning into $pFLD\alpha$

Introduction	pFLD α possesses the α -factor mating signal sequence for secretion of your protein. The information in this section is provided to assist you in designing a cloning strategy. Details of the multiple cloning site of the pFLD α can be found on page 6.				
Signal Sequence Processing	The processing of the α -factor mating signal sequence in pFLD α occurs in three steps:				
	1. Signal peptidase cleavage between Ala and Ala in the 19 and 20 positions.				
	2. Kex2 cleavage between Arg and Glu in the sequence Glu-Lys-Arg * Glu- Ala-Glu-Ala, where * is the site of cleavage.				
	3. The Glu-Ala repeats are further cleaved by the <i>STE13</i> gene product.				
Optimization of Signal Cleavage	In <i>Saccharomyces cerevisiae</i> , the Glu-Ala repeats are not necessary for cleavage by Kex2, but cleavage after Glu-Lys-Arg may be more efficient when followed by Glu-Ala repeats. A number of amino acids are tolerated at site X instead of Glu in the sequence Glu-Lys-Arg-X. These amino acids include the aromatic amino acids, small amino acids, and histidine. Proline, however, will inhibit Kex2 cleavage. For more information on Kex2 cleavage, see (Brake <i>et al.</i> , 1984). There are some cases where Ste13 cleavage of Glu-Ala repeats is not efficient, and Glu-Ala repeats are left on the N-terminus of the expressed protein of interest. This is generally dependent on the protein of interest.				
Expression of Recombinant Protein with Native N-terminus	To have your protein expressed with a native N-terminus, you can use the <i>Xho</i> I site at bp 865–870 to clone your gene flush with the Kex2 cleavage site. Use PCR to rebuild the sequence from the <i>Xho</i> I site to the arginine codon at nucleotides 744–746. Remember to include the first amino acid(s) of your protein, if necessary, for correct fusion to the Kex2 cleavage site.				

Cloning into pFLDα, Continued

Multiple Site of pl		-	ine see for		the c ing a nload	leava nd fui ling fo	ge site nctior or <u>ww</u>	e. The nal tes <u>vw.lif</u> e	e mult sting. etechi	iple c The v nolog	clonin rector <mark>ies.co</mark>	g site seque <u>m</u> or :	has b ence c from '	een co of pFL Techr	onfirr .Dα is nical S	ned b avail suppo	у
				3´	end of	FLD1											_
581	AAT'	ICTT	GAT 2	ATTC	ACAC	AA T'	TCAA	CAAC!	r at:	FTCG	AAAC		t Ar	g Ph	e Pr	o Se	r
													Signa	peptid	lase cle	eavage	
637				GCT Ala													
							α-fac	tor sign	al sequ	ence							
688	AAC Asn			ACA Thr													
739	GGT Gly			GAT Asp										Leu	Pro	Phe	Ser
														α-fa	ctor pri	ning sit	e
790				AAT Asn													
									Xho I		Kex	2 signa	al cleav	age			Mfe I
841				GAA Glu										Ala	Glu	Ala	
			Pml I		5	Sfi I			Asp	718 J	Kpn I		Sac I		13 sigr		vage
891	ATT		ACG :	rggco	CCAG	c'c Go	GCCG:	TCTCO	G GAT				scco	GCC (GCCAC	GCTTO	GG
951	Apa I GCC	1	rgg (GGT (_			-		AAC	V5 epi CCT (Pro I	CTC (_			
1003	ACG Thr	GGT Gly	GTC Val	GAC Asp	CAT His		•	ine (6x CAT His			TGA ***	GTT	IGTA (GCC 1	[TAG2	ACATO	SA
1056	CTG	TTCC:	ICA (GTTCI	AGT	rg go	GCAC!	TACO	G AGZ	AAGA	CCGG	TCT	IGCT	AGA 1	FTCT	ATC	AA
				<1 prim	-												
1116	GAG	GATG'	FCA (GAATO	GCCA?	CT TC	JCCT	GAGA	G ATC	GCAG	GCTT						

Transforming E. coli

Transforming <i>E. coli</i>	Transform your ligation mixtures into a competent <i>recA</i> , <i>endA E</i> . <i>coli</i> strain (e.g. TOP10, DH5 α^{TM} , JM109) and select on LB agar plates containing 50-100 µg/mL ampicillin. Once you have obtained ampicillin-resistant colonies, pick 10 transformants and screen for the presence and orientation of your insert. Note that there is no blue/white screening for the presence of insert with the pFLD vectors.
	We recommend that you sequence your construct to confirm that your gene is in the correct orientation for expression and cloned in frame with the C-terminal peptide (if desired). Refer to the multiple cloning site diagrams on the previous pages for the sequences and location of the priming sites. To order custom-synthesized primers, visit <u>www.lifetechnologies.com</u> and select Custom Primers.
Preparing a Glycerol Stock	Once you have identified the correct clone, purify the colony and make a glycerol stock for long-term storage. We also recommend keeping a DNA stock of your plasmid at –20°C.
	 Streak the original colony out on an LB plate containing 50–100 μg/mL ampicillin. Incubate the plate at 37°C overnight.
	 Isolate a single colony and inoculate into 1–2 mL of LB containing 50-100 µg/mL ampicillin.
	3. Grow the culture to mid-log phase ($OD_{600} = 0.5-0.7$).
	4. Mix 0.85 mL of culture with 0.15 mL of sterile glycerol and transfer to a cryovial.
	5. Store at -80° C.
Plasmid Preparation	Once you have cloned and sequenced your insert, generate enough plasmid DNA to transform <i>Pichia</i> (5–10 µg of each plasmid per transformation). We recommend isolating plasmid DNA using the We recommend isolating plasmid DNA using the PureLink [®] HiPure Miniprep Kit (up to 30 µg) or the PureLink [®] HiPure Midiprep Kit (up to 150 µg), or CsCl gradient centrifugation. Once you have purified plasmid DNA, proceed to <i>Pichia</i> Transformation on page 9.

Transformation into E. coli, Continued

Sequencing Recombinant Clones

We recommend that you sequence your construct before transforming into *Pichia* to confirm that your gene is in frame with the α -factor secretion signal and/or the C-terminal tag. To sequence your construct in pFLD or pFLD α , we recommend using *FLD1* Forward and 3' *AOX1* primers. The 3' *AOX1* primer is available separately for purchase (see page 29 for ordering information). The *FLD1* Forward primer can be custom-ordered from us; visit www.lifetechnologies.com and select Custom Primers.

Sequencing Primer	Sequence
α-Factor	5'-TACTATTGCCAGCATTGCTGC-3'
3´ AOX1	5'-GCAAATGGCATTCTGACATCC-3'

For sequencing protocols, refer to Unit 7 in *Current Protocols in Molecular Biology* (Ausubel *et al.*, 1994) or Chapter 13 in *Molecular Cloning: A Laboratory Manual* (Sambrook *et al.*, 1989).

Pichia Transformation

Introduction	After your gene has been correctly cloned into one of the pFLD vectors, you are ready to transform the vector into your <i>Pichia</i> strain. This section provides general guidelines for preparing plasmid DNA, transformation, and selecting for Zeocin [™] -resistant clones.
Zeocin [™] Selection	We typically use 100 µg/mL Zeocin TM to select for transformants when using the X-33 <i>Pichia</i> strain. If you are transforming your pFLD or pFLD α construct into another <i>Pichia</i> strain, note that selection conditions may vary. We recommend performing a dose response curve to determine the appropriate concentration of Zeocin TM to use for selection of transformants in your strain.
Method of Transformation	We do not recommend spheroplasting for transformation of <i>Pichia</i> with plasmids containing the Zeocin [™] resistance marker. Spheroplasting involves removal of the cell wall to allow DNA to enter the cell. Cells must first regenerate the cell wall before they are able to express the Zeocin [™] resistance gene. For this reason, plating spheroplasts directly onto selective medium containing Zeocin [™] does not yield any transformants.
	We recommend electroporation for transformation of <i>Pichia</i> with pFLD. Electroporation yields 10 ³ to 10 ⁴ transformants per µg of linearized DNA and does not destroy the cell wall of <i>Pichia</i> . If you do not have access to an electroporation device, use the LiCl protocol on page 27 or the <i>Pichia</i> EasyComp [™] Transformation Kit available for purchase (see below).
<i>Pichia</i> EasyComp [™] Transformation Kit	For chemical transformation of <i>Pichia</i> strains with pFLD vectors, the <i>Pichia</i> EasyComp [™] Transformation Kit is available for purchase (see page 29 for ordering information). The <i>Pichia</i> EasyComp [™] Transformation Kit provides reagents to prepare six preparations of competent cells. Each preparation will yield enough competent cells for 20 transformations. Competent cells may be used immediately or frozen and stored for future use. For more information, refer to <u>www.lifetechnologies.com/support</u> or call Technical Support (see page 31).
Q Important	Since the pFLD vectors contain DNA sequences from both the <i>FLD1</i> and <i>AOX1</i> loci (promoter and polyA addition, respectively), homologous recombination in the host can occur at either of these locations. The restriction enzyme site used to linearize the vector will determine the predominant integration locus. Cutting the vector within the <i>FLD1</i> promoter sequence (see Linearizing Your pFLD Construct on page 10) will target the integration event to the <i>FLD1</i> locus. If the vector is completely linearized within the <i>FLD1</i> promoter—with no uncut, circular vector remaining—homologous recombination at the <i>AOX1</i> locus will not occur.

Pichia Transformation, Continued



pFLD vectors do not contain a yeast origin of replication. Transformants can only be isolated if recombination occurs between the plasmid and the *Pichia* genome.

transformants on Ze	eocin [™] . Note : Inclusion o	f sorbitol in YPD plates stabilizes
• 5–10 μg pure pF	FLD or pFLD α containing	g your insert
• YPD Medium		
• 50-mL conical polypropylene tubes		
• 1 liter cold (4°C)) sterile water (place on i	ce the day of the experiment)
• 25 mL cold (4°C) sterile 1 M sorbitol (pla	ice on ice the day of the experiment)
• 30°C incubator		
Electroporation	device and 0.2-cm cuvet	tes
• YPDS plates cor for recipe)	ntaining the appropriate	concentration of Zeocin [™] (see page 18
within the 5' <i>FLD1</i> r linearize pFLD prior sure that your insert	egion. The table below li to transformation. Othe does not contain the res	sts unique sites that may be used to er restriction sites are possible. Be
	 transformants on Ze electroporated cells, 5–10 µg pure pF YPD Medium 50-mL conical p 1 liter cold (4°C) 25 mL cold (4°C) 30°C incubator Electroporation YPDS plates corfor recipe) To promote integrate within the 5' <i>FLD1</i> r linearize pFLD prior sure that your inserted set of the	 50-mL conical polypropylene tubes 1 liter cold (4°C) sterile water (place on i 25 mL cold (4°C) sterile 1 M sorbitol (pla 30°C incubator Electroporation device and 0.2-cm cuvet YPDS plates containing the appropriate

Restriction Digest

Nde I

Cla I

1. Digest \sim 5–10 µg of plasmid DNA with one of the enzymes listed above.

2. Check a small aliquot of your digest by agarose gel electrophoresis for complete linearization.

406

493

- 3. If the vector is completely linearized, heat inactivate or add EDTA to stop the reaction, phenol/chloroform extract once, and ethanol precipitate using 1/10 volume 3 M sodium acetate and 2.5 volumes of 100% ethanol.
- 4. Centrifuge the solution to pellet the DNA, wash the pellet with 80% ethanol, air-dry, and resuspend in 10 μ L sterile, deionized water. Use immediately or store at -20° C.

Pichia Transformation, Continued

Preparing Pichia		low the procedure below to prepare your Pichia pastoris strain for
for	elec	ctroporation.
Electroporation	1.	Grow 5 mL of your <i>Pichia pastoris</i> strain in YPD in a 50 mL conical tube at 30°C overnight.
	2.	Inoculate 500 mL of fresh medium in a 2 liter flask with 0.1–0.5 mL of the overnight culture. Grow overnight again to an $OD_{600} = 1.3-1.5$.
	3.	Centrifuge the cells at $1500 \times g$ for 5 minutes at 4°C. Resuspend the pellet with 500 mL of ice-cold (0°C), sterile water.
	4.	Centrifuge the cells as in Step 3, then resuspend the pellet with 250 mL of ice-cold (0°C), sterile water.
	5.	Centrifuge the cells as in Step 3, then resuspend the pellet in 20 mL of ice- cold (0°C) 1 M sorbitol.
	6.	Centrifuge the cells as in Step 3, then resuspend the pellet in 1 mL of ice- cold (0°C) 1 M sorbitol for a final volume of approximately 1.5 mL. Keep the cells on ice and use that day. Do not store cells.
Transformation by Electroporation	1.	Mix 80 μ L of the cells from Step 6 (above) with 5–10 μ g of linearized pFLD or pFLD α DNA (in 5–10 μ L sterile water) and transfer them to an ice-cold (0°C) 0.2-cm electroporation cuvette.
	2.	Incubate the cuvette with the cells on ice for 5 minutes.
	3.	Pulse the cells according to the parameters for yeast (<i>Saccharomyces cerevisiae</i>) as suggested by the manufacturer of the specific electroporation device being used.
	4.	Immediately add 1 mL of ice-cold 1 M sorbitol to the cuvette. Transfer the cuvette contents to a sterile 15 mL tube.
	5.	Let the tube incubate at 30°C without shaking for 1 to 2 hours.
	6.	Spread 50–200 μ L each on separate, labeled YPDS plates containing the appropriate concentration of Zeocin [™] .
	7.	Incubate plates for 2 to 3 days at 30°C until colonies form.
	8.	Pick 10–20 colonies and purify (streak for single colonies) on fresh YPD or YPDS plates containing the appropriate concentration of Zeocin [™] .

Pichia Transformation, Continued



Generally, several hundred Zeocin[™]-resistant colonies are generated using the protocol on page 11. If more colonies are needed, the protocol may be modified as described below. Note that you will need ~20 150-mm plates with YPDS agar containing the appropriate concentration of Zeocin[™].

- 1. Set up two transformations per construct and follow Steps 1 through 5 of the **Transformation by Electroporation** protocol, page 11.
- 2. After 1 hour in 1 M sorbitol at 30°C (Step 5, page 11), add 1 mL of YPD medium to each tube.
- 3. Shake (~200 rpm) the cultures at 30°C.
- After 1 hour, take one of the tubes and plate out all of the cells by spreading 200 μL on 150-mm plates containing the appropriate concentration of Zeocin[™].
- (Optional) Continue incubating the other culture for three more hours (for a total of four hours) and then plate out all of the cells by spreading 200 µL on 150-mm plates containing the appropriate concentration of Zeocin[™].
- 6. Incubate plates for 2 to 4 days at 30°C until colonies form.

Mut Phenotype If you used a *Pichia* strain containing a native *AOX1* gene (e.g., X-33, GS115, SMD1168H) as the host for your pFLD construct, your Zeocin[™]-resistant transformants will be Mut⁺. If you used a strain containing a deletion in the *AOX1* gene (e.g., KM71H), your transformants will be Mut^S.

To verify the Mut phenotype of your Zeocin[™]-resistant transformants, refer to the general guidelines provided in the EasySelect[™] *Pichia* Expression Kit manual or the Original *Pichia* Expression Kit manual or to published reference sources (Higgins and Cregg, 1998).

You are now ready to test your transformants for expression of your gene of interest. See **Expression in** *Pichia*, page 13.

Expression in Pichia

Introduction	The primary purpose of small-scale expression is to identify/confirm a recombinant <i>Pichia</i> clone that is expressing the correct protein. Small-scale expression conditions may not be optimal for your protein. For this reason, the method you choose for detection (e.g., SDS-PAGE, Western, or functional assay) may be an important factor in determining the success of expression. If your method of detection does not reveal any expression, you may want to consider using a more sensitive method.
	Once a positive clone has been identified, large-scale expression can be carried out in shake flask or fermentation, and expression conditions can be optimized.
Note	Note that once you have obtained Zeocin [™] -resistant transformants, it is not necessary to maintain your recombinant <i>Pichia</i> clone in medium containing Zeocin [™] for expression studies. Zeocin [™] is only required for initial screening and selection of recombinant clones.
Dotocting	We recommend that you use the following techniques to assay expression of

Detecting Recombinant Proteins in *Pichia*

We recommend that you use the following techniques to assay expression of your protein. Be sure to account for any additional amino acids that are in between the end of your protein and the C-terminal tag.

Technique	Method of Detection	Sensitivity
SDS-PAGE (Coomassie-stained)	Visualization by eye	Can detect as little as 100 ng in a single band
SDS-PAGE (Silver-stained)	Visualization by eye	Can detect as little as 2 ng in a single band
Western Analysis	Antibody to your particular protein Anti- <i>myc</i> antibodies (see the next page) Anti-His(C-term) antibodies (see the next page)	Can detect as little as 1-10 pg, depending on detection method (alkaline phosphatase, horseradish peroxidase, radiolabeled antibody)
Functional assay	Varies depending on assay.	Varies depending on assay Used to compare relative amounts of protein.



Reminder: Because the pFLD vector does not contain the *HIS4* gene, *his4 Pichia* strains containing the integrated plasmid must be grown in medium containing 0.004% histidine. If histidine is not present in the medium the cells will not grow. If you use X-33, SMD1168H, or KM71H as the host strain, supplementation of the medium with histidine is not required.

Expression in Pichia, Continued

Expression Guidelines	Information and guidelines for performing small-scale expression, optimizing expression, and performing scale-up of expression are provided in the EasySelect [™] <i>Pichia</i> Expression Kit manual and the Original <i>Pichia</i> Expression Kit manual. See below for information on inducing expression of the <i>FLD1</i> gene with methanol, methylamine, or methanol plus methylamine.
Inducing with Methanol, Methylamine, or Methanol Plus Methylamine	 Expression of the <i>FLD1</i> gene may be induced with methanol, methylamine, or methanol plus methylamine. The following guidelines for induction have been developed using the wild-type X-33 strain; other strains may be suitable. For media recipes, see pages 19–21. 1. For all types of induction, grow cultures in MGAs medium at 30°C in a shaking incubator (250–300 rpm) until the culture reaches an OD₆₀₀ of 4. 2. Harvest the cells by centrifuging at 1500–3000 × g for 5 minutes at room temperature. Then:
	 For induction with methanol, decant supernatant and resuspend cell pellet in 1/5 of the original culture volume of MMAs medium.
	 For induction with methylamine, decant supernatant and resuspend cell pellet in 1/5 of the original culture volume of MGMa medium.
	• For induction with methanol plus methylamine , decant supernatant and resuspend cell pellet in 1/5 of the original culture volume of MMMa medium.
	3. Incubate the resulting culture at 30°C for five days. For media containing methanol or methanol plus methylamine, add 100% methanol to a final concentration of 0.5% methanol every 24 hours to maintain induction.
	Time points: At each of the times indicated below, transfer 1 mL of the expression culture to a 1.5-mL microcentrifuge tube, and centrifuge at maximum speed in a tabletop microcentrifuge for 2–3 minutes at room temperature. These samples will be used to analyze expression levels and determine the optimal time post-induction to harvest. Time points (hours): 0, 24 (1 day), 48 (2 days), 72 (3 days), 96 (4 days), and 120 (5 days).
Polyacrylamide Gel Electrophoresis	To facilitate separation and visualization of your recombinant protein by polyacrylamide gel electrophoresis, a wide range of pre-cast NuPAGE [®] and Tris-Glycine polyacrylamide gels are available for purchase. The NuPAGE [®] Gel System avoids the protein modifications associated with Laemmli-type SDS-PAGE, ensuring optimal separation for protein analysis. In addition, we also carry a large selection of molecular weight protein standards and staining kits. For more information about the appropriate gels, standards, and stains to use to visualize your recombinant protein, refer to <u>www.lifetechnologies.com</u> or call Technical Support (see page 31).

Western Analysis	To detect expression of your recombinant fusion protein by Western blot
	analysis, you may use the Anti-V5 antibodies or the Anti-His(C-term) antibodies
	(see page 30 for ordering information) or an antibody to your protein of interest.
	In addition, the Positope [™] Control Protein available for use as a positive control
	for detection of fusion proteins containing a V5 epitope or a polyhistidine (6xHis)
	tag (see page Error! Bookmark not defined. for ordering). WesternBreeze®
	Chromogenic Kits and WesternBreeze® Chemiluminescent Kits are available for
	detection of antibodies by colorimetric or chemiluminescent methods. For more
	information, refer to <u>www.lifetechnologies.com</u> or call Technical Support (see
	page 31).

Purification

Introduction	In this section, you will grow and induce a 10–200 mL culture of your <i>Pichia</i> transformant for trial purification on a metal-chelating resin such as ProBond [™] . You can harvest the cells and store them at -80°C until you are ready to purify your fusion protein, or you can proceed directly with protein purification. Note that this section only describes preparation of cell lysates and sample application onto ProBond[™] . For instructions on how to prepare and use ProBond [™] resin, refer to the ProBond [™] Purification System manual.
ProBond [™] Resin	We recommend the ProBond [™] Purification System to purify fusion proteins expressed from pFLD or pFLDα (see page Error! Bookmark not defined.). Note that instructions for equilibration of and chromatography on ProBond [™] resin are contained in the ProBond [™] Purification Kit.
	If you are using a metal-chelating resin other than ProBond [™] , follow the manufacturer's recommendations to purify fusion proteins expressed in bacteria or yeast.
Binding Capacity of ProBond [™]	One milliliter of ProBond [™] resin binds at least 1 mg of recombinant protein. This amount can vary depending on the protein.
O Important	Throughout the following protocol, be sure to keep the cell lysate and fractions on ice. Small-scale purifications using the 2-mL ProBond [™] columns and buffers can be performed at room temperature on the bench top. For large scale purifications, all reagents must be kept at 4°C.
Preparing Cell Lysates	Express your protein using a small-scale culture (10–20 mL for Mut ^S strains; 100-200 mL for Mut ⁺) and the optimal conditions for expression (if determined). Refer to the <i>Pichia</i> Expression Kit manual for details. Once your protein is expressed, follow the protocol below to prepare a cell lysate for chromatography on ProBond [™] .
	Prepare Breaking Buffer (BB) as described in the Recipes , page 21.
	1. Wash cells once in BB by resuspending them and centrifuging 5–10 minutes at $3000 \times g$ at 4°C.
	2. Resuspend the cells to an OD_{600} of 50–100 in BB.
	 Add an equal volume of acid-washed glass beads (0.5 mm). Estimate volume by displacement.
	 Vortex the mixture for 30 seconds, then incubate on ice for 30 seconds. Repeat 7 more times. Alternating vortexing with cooling keeps the cell extracts cold and reduces denaturation of your protein.
	5. Centrifuge the sample at 4°C for 5–10 minutes at 12,000 × g .
	 Transfer the clear supernatant to a fresh container and analyze for your protein. The total protein concentration should be around 2–3 mg/mL.
	7. Save the pellet and extract with 6 M urea or 1% Triton X-100 to check for insoluble protein.

Purification, Continued

Sample Application	For sample application onto ProBond [™] , you will need Native Binding Buffer, pH 7.8 and a 2-mL ProBond [™] column, pre-equilibrated using native conditions.
(Native Conditions)	 Combine 1 mL (2–3 mg/mL total protein) of <i>Pichia</i> lysate with 7 mL Native Binding Buffer.
	 Prepare a ProBond[™] column with resin as described in the ProBond[™] manual, and resuspend the resin in the column in 4 mL of the diluted lysate from Step 1.
	3. Seal the column and batch-bind by rocking gently at room temperature for 10 minutes.
	4. Let the resin settle by gravity or low speed centrifugation $(800 \times g)$ and carefully remove the supernatant. Save the supernatant to check for unbound protein.
	5. Repeat Steps 2 through 4 with the remaining 4 mL of diluted lysate. Proceed with column washing and elution under native conditions as described in the ProBond [™] manual. Use the recommendations noted for bacterial cell lysates.
Sample Applica- tion (Denaturing Conditions)	Use the protocol above except pre-equilibrate the ProBond [™] column using Denaturing Binding Buffer and combine 1 mL of the <i>Pichia</i> cell lysate with 7 mL of the Denaturing Binding Buffer.
Note	We have observed that some <i>Pichia</i> proteins may be retained on the ProBond [™] column using native purification conditions. Optimization of the purification (see the ProBond [™] manual) or using denaturing purification may remove these non-specific <i>Pichia</i> proteins.
Analysis of Purification	Save all fractions, washes, and flow-through for analysis by SDS-PAGE. You may need to use western blot analysis to detect your protein if expression is low or not enough protein was loaded onto the column. Refer to the ProBond [™] System manual for a guide to troubleshoot chromatography.
Scale-up	You may find it necessary to scale-up your purification to obtain sufficient amounts of purified protein. Adjust the pH and NaCl concentration of your lysate as indicated in the ProBond [™] manual before adding it to the column. The pH should be greater than or equal to 7.5 and the NaCl concentration should be ~500 mM.

Appendix

Recipes	
YPD (+ Zeocin [™])	Yeast Extract Peptone Dextrose Medium (1 liter)
	1% yeast extract
	2% peptone
	2% dextrose (glucose)
	<u>+</u> 2% agar
	<u>+</u> appropriate concentration of Zeocin [™]
	1. Dissolve 10 g yeast extract and 20 g of peptone in 900 mL of water.
	2. Include 20 g of agar if making YPD slants or plates.
	3. Autoclave for 20 minutes on liquid cycle.
	4. Add 100 mL of 20% dextrose (filter-sterilize dextrose before use).
	5. Cool solution to ~60°C and add the appropriate amount of Zeocin [™] from a 100 mg/mL stock solution. Note: It is necessary to include Zeocin [™] in the medium for selection of <i>Pichia</i> transformants only. Zeocin [™] may be omitted from the medium when performing expression studies.
	Store YPD slants or plates containing Zeocin [™] at 4°C. The shelf life is 1–2 weeks.
YPDS (+ Zeocin [™])	<u>Y</u>east Extract <u>Peptone</u> <u>Dextrose</u> Medium with <u>S</u> orbitol (1 liter)
	1% yeast extract
	2% peptone
	2% dextrose (glucose)
	1 M sorbitol
	<u>+</u> 2% agar
	<u>+</u> appropriate concentration of Zeocin [™]
	1. Dissolve: 10 g yeast extract
	182.2 g sorbitol
	20 g of peptone
	in 900 mL of water.
	2. Add 20 g of agar.
	3. Autoclave for 20 minutes on liquid cycle.
	4. Add 100 mL of 20% dextrose (filter-sterilize dextrose before use).
	5. Cool solution to ~60°C and add the appropriate amount of Zeocin [™] from a 100 mg/mL stock solution. Note: It is necessary to include Zeocin [™] in the medium for selection of <i>Pichia</i> transformants only. Zeocin [™] may be omitted from the medium when performing expression studies.
	Store YPDS slants or plates containing Zeocin [™] at 4°C. The shelf life is 1–2 weeks.

Recipes, Continued

Stock Solutions	10X YNB (1.7% Yeast Nitrogen Base without Ammonium Sulfate)
	Dissolve 17 g of yeast nitrogen base (YNB) without ammonium sulfate and without amino acids in 1000 mL of water and filter sterilize. Heat the solution to dissolve YNB completely in water. Store at 4°C. The shelf life of this solution is approximately one year.
	10X As (5% Ammonium Sulfate)
	Dissolve 50 g of ammonium sulfate in 1000 mL of water and filter sterilize.
	500X B (0.02% Biotin)
	Dissolve 20 mg biotin in 100 mL of water and filter sterilize. Store at 4°C. The shelf life of this solution is approximately one year.
	10X M (5% Methanol)
	Mix 5 mL of methanol with 95 mL of water. Filter sterilize and store at 4°C. The shelf life of this solution is approximately two months.
	10X G (10% Glycerol)
	Mix 100 mL of glycerol with 900 mL of water. Sterilize either by filtering or autoclaving. Store at room temperature. The shelf life of this solution is greater than one year.
	10X Ma (2.5% Methylamine)
	Mix 2.5 mL of methylamine HCl with 97.5 mL of water. Filter sterilize and store at 4°C. The shelf life of this solution is approximately two months.
MGAs	Minimal Glycerol with Ammonium sulfate (1 liter)
	0.17% YNB
	0.5% ammonium sulfate
	1% glycerol $4 \times 10^{-5\%}$ biotin
	Combine aseptically 700 mL autoclaved water with 100 mL of 10X YNB, 100 mL of As, 2 mL of 500X B, and 100 mL of 10X G.
	Store at 4°C. The shelf life of this solution is approximately two months.

Recipes, Continued

MMAs	Minimal Methanol with Ammonium sulfate Medium (1 liter)
	 0.17% YNB 0.5% ammonium sulfate 4 × 10⁻⁵% biotin 0.5% methanol 1. For medium, autoclave 700 mL of water for 20 minutes on liquid cycle 2. Cool autoclaved water to 60°C and add: 100 mL of 10X YNB 100 mL of As 2 mL of 500X B 100 mL of 10X M
	 Mix and store at 4°C. For plates, add 20 g agar to the water in Step 1 and proceed. After mixing, pour the plates immediately. MMAs stores for several months at 4°C.
MGMa	 <u>Minimal Glycerol and Methylamine Medium</u> (1 liter) 0.17% YNB 4 × 10⁻⁵% biotin 1% glycerol 0.25% methylamine 1. For medium, autoclave 700 mL of water for 20 minutes on liquid cycle 2. Cool autoclaved water to room temperature and add: 100 mL of 10X YNB 2 mL of 500X B 100 mL 10X G 100 mL of 10X Ma
	 Mix and store at 4°C. For plates, add 20 g agar to the water in Step 1 and proceed. After mixing, pour the plates immediately. MGMa stores for several months

at 4°C.

Recipes, Continued

МММа	0.1 4 × 0.5 0.2 1.	inimal <u>M</u> ethanol and <u>M</u> ethy 7% YNB 10 ⁻⁵ % biotin % methanol 5% methylamine For medium, autoclave 700 Cool autoclaved water to 6 100 mL of 10X YNB 2 mL of 500X B 100 mL of 10X M 100 mL of 10X Ma	mL of water for 20 minutes on liquid cycle
	3. 4. 5.		the water in Step 1 and proceed. es immediately. MMMa stores for several months
Breaking Buffer	1 n inł 1 n	mM sodium phosphate, pH nM PMSF (phenylmethylsul: nibitors) nM EDTA 9 glycerol	7.4 fonyl fluoride. You may use other protease
	1.		your desired protease inhibitors and store ufacturer's recommendations.
	2.	For 1 liter, dissolve:	6 g sodium phosphate (monobasic) 372 mg EDTA 50 mL glycerol
		in 900 mL deionized water	
	3.	Use NaOH to adjust pH an	d bring up the volume to 1 liter. Store at 4°C.
	4.	Add protease inhibitors im	mediately before use.

Zeocin[™]

Zeocin[™] is a member of the bleomycin/phleomycin family of antibiotics isolated from *Streptomyces*. It shows strong toxicity against bacteria, fungi, plants and mammalian cell lines (Calmels *et al.*, 1991; Drocourt *et al.*, 1990; Gatignol *et al.*, 1987; Mulsant *et al.*, 1988; Perez *et al.*, 1989).

A Zeocin[™] resistance protein has been isolated and characterized (Calmels *et al.*, 1991; Drocourt *et al.*, 1990). This 13,665 Da protein, the product of the *Sh ble* gene (*Streptoalloteichus hindustanus* bleomycin gene), binds stoichiometrically to Zeocin[™] and inhibits its DNA strand cleavage activity. Expression of this protein in eukaryotic and prokaryotic hosts confers resistance to Zeocin[™].

Molecular Weight, Formula, and Structure

The formula for ZeocinTM is $C_{55}H_{86}O_{21}N_{20}S_2Cu$ -HCl and the molecular weight is 1527.5. The structure of ZeocinTM is shown below.



Applications of Zeocin[™]

Zeocin^M is used for selection in mammalian cells (Mulsant *et al.*, 1988); plants (Perez *et al.*, 1989); yeast (Baron *et al.*, 1992); and prokaryotes (Drocourt *et al.*, 1990). Suggested concentrations of Zeocin^M for selection in *Pichia* and *E. coli* are listed below:

Organism	Zeocin [™] Concentration and Selective Medium
E. coli	25–50 μg/mL in Low Salt LB medium [*]
Pichia	100–1000 μ g/mL (varies with strain and medium)
*Efficient selection requires that the concentration of NaCl be no more than 5 g/L (< 90 mM)	

Zeocin[™], Continued

Handling Zeocin [™]	•	High ionic strength and acidity or basicity inhibits the activity of Zeocin [™] . Therefore, we recommend that you reduce the salt in bacterial medium and adjust the pH to 7.5 to keep the drug active (see page 18 for a recipe).
	•	Store Zeocin ^{TM} at –20°C and thaw on ice before use.

- Zeocin[™] is light sensitive. Store the drug and plates or medium containing the drug in the dark.
- Wear gloves, a laboratory coat, and safety glasses when handling Zeocin[™]- containing solutions.
- Do not ingest or inhale solutions containing the drug.

Map of pFLDThe figure below summarizes the features of the pFLD vector. The vector
sequence for pFLD is available for downloading from www.lifetechnologies.com
or from Technical Support (see page 31). See page 26 for a description of the
features of the vector.



Map of pFLD α

Map of pFLDαThe figure below summarizes the features of the pFLDα vector. The complete
sequence for pFLDα is available for downloading from
www.lifetechnologies.com or from Technical Support (see page 31). See page 26
for a description of the features of the vector.



Zeocin resistance gene: bases 3983-4357

CYC1 transcription termination region: bases 4358-4675

Vector Features

Features

The vectors pFLD and pFLD α contain the following elements. All features have been functionally tested.

Feature	Benefit
FLD1 promoter	A 597-bp fragment containing the <i>FLD1</i> promoter that allows methanol- or methylamine-inducible, high-level expression of the gene of interest in <i>Pichia</i>
	Targets plasmid integration to the FLD1 locus.
α-Factor Secretion Signal (pFLDα only)	Encodes the native <i>Saccharomyces cerevisiae</i> α-factor secretion signal that allows for efficient secretion of most proteins from <i>Pichia</i> (Cregg <i>et al.</i> , 1993)
Multiple cloning site	Allows insertion of your gene into the expression vector
V5 epitope (Gly-Lys-Pro-Ile-Pro-Asn- Pro-Leu-Leu-Gly-Leu-Asp- Ser-Thr)	Permits detection of your recombinant fusion protein with the Anti-V5 Antibody or Anti-V5-HRP Antibody (Southern <i>et al.</i> , 1991) (see page 30 for ordering)
C-terminal polyhistidine (6xHis) tag	Permits purification of your recombinant fusion protein on metal-chelating resin such as $ProBond^{TM}$ In addition, the C-terminal polyhistidine tag is the epitope for the Anti-His(C-term) Antibody (Lindner <i>et al.</i> , 1997) and the Anti-His(C-term)-HRP Antibody (see page Error! Bookmark not defined. for ordering)
<i>AOX1</i> transcription termination (TT) region	Native transcription termination and polyadenylation signal from <i>AOX1</i> gene (~260 bp) that permits efficient 3´ mRNA processing, including polyadenylation, for increased mRNA stability
<i>TEF1</i> promoter (GenBank accession numbers D12478, D01130)	Transcription elongation factor 1 gene promoter from <i>Saccharomyces cerevisiae</i> that drives expression of the Zeocin [™] resistance gene in <i>Pichia</i>
T7 promoter	Synthetic prokaryotic promoter that drives constitutive expression of the Zeocin [™] resistance gene in <i>E. coli</i>
Zeocin TM resistance gene (<i>Sh ble</i>)	Allows selection of transformants in <i>E. coli</i> and <i>Pichia</i>
CYC1 transcription termination region (GenBank accession number M34014)	3´ end of the <i>Saccharomyces cerevisiae</i> CYC1 gene that allows efficient 3´ mRNA processing of the Zeocin [™] resistance gene for increased stability
pUC origin	Allows replication and maintenance of the plasmid in <i>E. coli</i>

Map of pFLD/CAT (5014 bp) is a control vector expressing chloramphenicol acetyltransferase (CAT). It was constructed by cloning the CAT gene into the pFLD vector. The figure below summarizes the features of the pFLD/CAT vector. The vector sequence for pFLD/CAT is available for downloading from www.lifetechnologies.com or from Technical Support (see page 31).



Comments for pFLD/CAT: 5014 nucleotides

FLD1 promoter region: bases 1-597 CATORF: bases 612-1268 V5 epitope: bases 1284-1325 Polyhistidine (6xHis) tag: bases 1335-1352 *AOX1* transcription termination region: bases 1363-1691 3' *AOX1* priming site: bases 1438-1458 *bla* promoter: bases 2109-2115 Ampicillin resistance gene: bases 2150-3010 pUC origin: bases 3155-3827 *TEF1* promoter: bases 3828-4230 T7 promoter: bases 4236-4302 Zeocin resistance gene: bases 4303-4677 *CYC1* transcription termination region: bases 4678-4995

Lithium Chloride Transformation Method

Introduction	This modified version of the procedure described for <i>S. cerevisiae</i> (Gietz and Schiestl, 1996) is provided as an alternative to transformation by electroporation. Transformation efficiency is between 10^2 and 10^3 cfu/µg linearized DNA.
Preparing Solutions	Lithium acetate does not work with <i>Pichia pastoris.</i> Use only lithium chloride. 1 M LiCl in distilled, deionized water. Filter-sterilize. Dilute as needed with sterile water.
	50% polyethylene glycol (PEG-3350) in distilled, deionized water. Filter-sterilize. Store in a tightly capped bottle.
	2 mg/mL denatured, sheared salmon sperm DNA in TE (10 mM Tris-HCl, pH 8.0, 1.0 mM EDTA). Store at -20° C.
Preparing Cells	1. Grow a 50 mL culture of <i>Pichia pastoris</i> in YPD at 30°C with shaking to an OD_{600} of 0.8 to 1.0 (approximately 10 ⁸ cells/mL).
	2. Harvest the cells, wash with 25 mL of sterile water, and centrifuge at $1500 \times g$ for 10 minutes at room temperature.
	3. Resuspend the cell pellet in 1 mL of 100 mM LiCl and transfer the suspension to a 1.5 mL microcentrifuge tube.
	 Pellet cells at maximum speed for 15 seconds and remove LiCl with a pipette.
	5. Resuspend the cells in 400 μ L of 100 mM LiCl.
	6. Dispense 50 μL of cell suspension into a 1.5-mL microcentrifuge tube for each transformation and use immediately. Do not store on ice or freeze at -20°C.
Transformation	 Boil a 1 mL sample of single-stranded DNA for 5 minutes, then quickly chill on ice. Keep on ice. Note: It is neither necessary nor desirable to boil the carrier DNA prior to each use. Store a small aliquot at -20°C and boil every 3-4 times the DNA is thawed.
	2. Centrifuge the cells from Step 6, above, and remove the LiCl with a pipette.
	3. For each transformation, add the following reagents IN THE ORDER GIVEN to the cells. PEG shields the cells from the detrimental effects of the high LiCl concentration.
	240 μL 50% PEG 36 μL 1 M LiCl
	$25 \ \mu\text{L} 2 \ \text{mg/mL}$ single-stranded DNA Plasmid DNA (5–10 \mug) in 50 \mu L sterile water
	 Vortex each tube vigorously until the cell pellet is completely mixed (~1 minute).
	5. Incubate the tube at 30°C for 30 minutes without shaking.
	6. Heat shock in a water bath at 42° C for 20–25 minutes.
	7. Centrifuge the cells at 6000 to 8000 rpm to pellet.
	 Resuspend the pellet in 1 mL of YPD and incubate at 30°C with shaking. After 1 hour and 4 hours, plate 25 to 100 µL on YPD plates containing the appropriate concentration of Zeocin[™]. Incubate the plates for 2–3 days at 30°C.

Accessory Products

Introduction	For more in		to <u>www.lifetechno</u>	l for use with the pFLD vectors. plogies.com/support or call	
Zeocin [™]	prepared i	Zeocin [™] may be available for purchase. For your convenience, the drug is prepared in autoclaved, deionized water and available in 1.25 mL aliquots at a concentration of 100 mg/mL.			
		Amount	Catalog no.		
		1 g	R250-01	_	
		5 g	R250-05		

Additional Products Many reagents that may be used with the pFLD vectors and for *Pichia* expression are available for purchase. Ordering information is provided below.

Item	Amount	Catalog no.
Ampicillin, Sodium Salt, lyophilized	20 mL	11593-019
X-33 Pichia strain	1 stab	C180-00
GS115 Pichia strain	1 stab	C181-00
KM71H Pichia strain	1 stab	C182-00
SMD1168H Pichia strain	1 stab	C184-00
3' AOX1 Pichia Primer	2 µg	N720-02
Original Pichia Expression Kit	1 kit	K1710-01
EasySelect [™] <i>Pichia</i> Expression Kit	1 kit	K1740-01
<i>Pichia</i> EasyComp [™] Transformation Kit	1 kit	K1730-01
Pichia Protocols	1 book	G100-01
CAT Antiserum (supplied amount is sufficient to perform 25 Western blots using 10 mL of working solution per reaction)	50 µL	R902-25
One Shot [®] TOP10 Chemically Competent <i>E. coli</i>	$20 \times 50 \ \mu L$	C4040-03
One Shot [®] TOP10 Electrocomp [™] (electrocompetent cells)	21 × 50 μL	C4040-52
Electrocomp [™] TOP10 (electrocompetent cells)	$5 \times 80 \ \mu L$	C664-55
PureLink [®] HQ Mini Plasmid Purification Kit	100 reactions	K2100-01
PureLink [®] HiPure Miniprep Kit	100 preps	K2100-03
PureLink [®] HiPure Midiprep Kit	25 preps	K2100-04

Accessory Products, Continued

Detecting Fusion Protein A number of antibodies are available for purchase to detect expression of your fusion protein from the pFLD vector. Horseradish peroxidase (HRP) or alkaline phosphatase (AP)-conjugated antibodies allow one-step detection using colorimetric or chemiluminescent detection methods. The fluorescein isothiocyanate (FITC)-conjugated antibody allows one-step detection in immunofluorescence experiments.

The amount of antibody supplied is sufficient for 25 Western blots or 25 immunostaining reactions (FITC-conjugated antibody only).

Antibody	Epitope	Catalog no.
Anti-V5	Detects 14 amino acid epitope derived	R960-25
Anti-V5-HRP	from the P and V proteins of the paramyxovirus, SV5 (Southern et al.,	R961-25
Anti-V5-AP Antibody	1991).	R962-25
Anti-V5-FITC Antibody	GKPIPNPLLGLDST	R963-25
Anti-His(C-term)	Detects the C-terminal polyhistidine	R930-25
Anti-His(C-term)-HRP	(6xHis) tag (requires the free carboxyl group for detection) (Lindner <i>et al.,</i> 1997):	R931-25
	ННННН-СООН	
Positope [™] Control Protein		R900-50

Purifying Fusion Protein

The polyhistidine (6xHis) tag allows purification of the recombinant fusion protein using metal-chelating resins such as ProBond[™]. Ordering information for ProBond[™] resin is provided below.

Item	Quantity	Catalog no.
ProBond [™] Purification System	Precharged ProBond [™] resin and buffers plus 6 × 2 mL columns for native and denaturing purification	K850-01
ProBond [™] Purification System with Anti-V5-HRP Antibody	1 Kit (same as above plus 50 μL of antibody) The amount of antibody supplied is sufficient for 25 Westerns	K854-01
ProBond [™] Purification System with Anti-His(C- term)-HRP Antibody	1 Kit (same as above plus 50 μL of antibody) The amount of antibody supplied is sufficient for 25 Westerns	K853-01
ProBond [™] Resin	50 mL	R801-01
	150 mL	R801-15
Purification Columns	50 polypropylene columns	R640-50

Technical Support

Obtaining support	 For the latest services and support information for all locations, go to <u>www.lifetechnologies.com/support</u>. At the website, you can: Access worldwide telephone and fax numbers to contact Technical Support and Sales facilities Search through frequently asked questions (FAQs) Submit a question directly to Technical Support (<u>techsupport@lifetech.com</u>) Search for user documents, SDSs, vector maps and sequences, application notes, formulations, handbooks, certificates of analysis, citations, and other product support documents Obtain information about customer training Download software updates and patches
Safety Data Sheets (SDS)	Safety Data Sheets (SDSs) are available at <u>www.lifetechnologies.com/support</u> .
Certificate of Analysis	The Certificate of Analysis provides detailed quality control and product qualification information for each product. Certificates of Analysis are available on our website. Go to <u>www.lifetechnologies.com/support</u> and search for the Certificate of Analysis by product lot number, which is printed on the box.
Limited warranty	Life Technologies Corporation 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 a Life Technologies product or service, contact our Technical Support Representatives. All Life Technologies products are warranted to perform according to specifications stated on the certificate of analysis. The Company will replace, free of charge, any product that does not meet those specifications. <u>This warranty</u> <u>limits the Company's liability to only the price of the product</u> . 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. The Company reserves the right to select the method(s) used to analyze a product unless the Company agrees to a specified method in writing prior to acceptance of the order. Life Technologies makes every effort to ensure the accuracy of its publications, but realizes that the occasional typographical or other error is inevitable. Therefore the Company makes no warranty of any kind regarding the contents of any publications or documentation. If you discover an error in any of our publications, report it to our Technical Support Representatives. Life Technologies Corporation shall have 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

License No: to Research Use cc Only pt w of ob	he purchase of this product conveys to the purchaser the limited, non-transferable right o use the purchased amount of the product only to perform internal research for the sole enefit of the purchaser. No right to resell this product or any of its components is ponveyed expressly, by implication, or by estoppel. This product is for internal research urposes only and is not for use in commercial applications of any kind, including, vithout limitation, quality control and commercial services such as reporting the results f purchaser's activities for a fee or other form of consideration. For information on btaining additional rights, please contact outlicensing@lifetech.com or Out Licensing, ife Technologies, 5791 Van Allen Way, Carlsbad, California 92008.
--	--

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).
- Baron, M., Reynes, J. P., Stassi, D., and Tiraby, G. (1992). A Selectable Bifunctional β-Galactosidase: Phleomycin-resistance Fusion Protein as a Potential Marker for Eukaryotic Cells. Gene 114, 239-243.
- Calmels, T., Parriche, M., Burand, H., and Tiraby, G. (1991). High Efficiency Transformation of *Tolypocladium geodes* Conidiospores to Phleomycin Resistance. Curr. Genet. 20, 309-314.
- Drocourt, D., Calmels, T. P. G., Reynes, J. P., Baron, M., and Tiraby, G. (1990). Cassettes of the *Streptoalloteichus hindustanus ble* Gene for Transformation of Lower and Higher Eukaryotes to Phleomycin Resistance. Nucleic Acids Res. *18*, 4009.
- Ellis, S. B., Brust, P. F., Koutz, P. J., Waters, A. F., Harpold, M. M., and Gingeras, T. R. (1985). Isolation of Alcohol Oxidase and Two other Methanol Regulatable Genes from the Yeast, *Pichia pastoris*. Mol. Cell. Biol. *5*, 1111-1121.
- Evans, G. I., Lewis, G. K., Ramsay, G., and Bishop, V. M. (1985). Isolation of Monoclonal Antibodies Specific for *c-myc* Proto-oncogene Product. Mol. Cell. Biol. 5, 3610-3616.
- Gietz, R. D., and Schiestl, R. H. (1996) Transforming Yeast with DNA. In *Methods in Molecular and Cellular Biology*: in press).
- Henikoff, S., and Cohen, E. H. (1984). Sequences Responsible for Transcription Termination on a Gene Segment in *Saccharomyces cerevisiae*. Mol. Cell. Biol. *4*, 1515-1520.
- Higgins, D. R., and Cregg, J. M. (1998) *Pichia* Protocols. In *Methods in Molecular Biology*, Vol. 103. (J. M. Walker, ed. Humana Press, Totowa, NJ.
- Irniger, S., Egli, C. M., and Braus, G. H. (1991). Different Classes of Polyadenylation Sites in the Yeast *Saccharomyces cerevisiae*. Mol. Cell. Bio. *11*, 3060-3069.
- Koutz, P. J., Davis, G. R., Stillman, C., Barringer, K., Cregg, J. M., and Thill, G. (1989). Structural Comparison of the *Pichia pastoris* Alcohol Oxidase Genes. Yeast 5, 167-177.
- Lindner, P., Bauer, K., Krebber, A., Nieba, L., Kremmer, E., Krebber, C., Honegger, A., Klinger, B., Mocikat, R., and Pluckthun, A. (1997). Specific Detection of His-tagged Proteins With Recombinant Anti-His Tag scFv-Phosphatase or scFv-Phage Fusions. BioTechniques 22, 140-149.
- Mulsant, P., Tiraby, G., Kallerhoff, J., and Perret, J. (1988). Phleomycin Resistance as a Dominant Selectable Marker in CHO Cells. Somat. Cell Mol. Genet. *14*, 243-252.
- Perez, P., Tiraby, G., Kallerhoff, J., and Perret, J. (1989). Phleomycin Resistance as a Dominant Selectable Marker for Plant Cell Transformation. Plant Mol. Biol. *13*, 365-373.
- Romanos, M. A., Scorer, C. A., and Clare, J. J. (1992). Foreign Gene Expression in Yeast: A Review. Yeast 8, 423-488.
- Sambrook, J., Fritsch, E. F., and Maniatis, T. (1989). Molecular Cloning: A Laboratory Manual, Second Edition (Plainview, New York: Cold Spring Harbor Laboratory Press).
- Scorer, C. A., Buckholz, R. G., Clare, J. J., and Romanos, M. A. (1993). The Intracellular Production and Secretion of HIV-1 Envelope Protein in the Methylotrophic Yeast *Pichia pastoris*. Gene *136*, 111-119.
- Shen, S., Sulter, G., Jeffries, T.W., and Cregg, J.M. (1998) A Strong Nitrogen Source-Regulated Promoter for Controlled Expression of Foreign Genes in the Yeast *Pichia pastoris*. Gene 216, 93-102
- 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.

References, Continued

Tschopp, J. F., Brust, P. F., Cregg, J. M., Stillman, C., and Gingeras, T. R. (1987a). Expression of the *lacZ* Gene from Two Methanol Regulated Promoters in *Pichia pastoris*. Nucleic Acids Res. 15, 3859-3876.

Zaret, K. S., and Sherman, F. (1984). Mutationally Altered 3' Ends of Yeast CYC1 mRNA Affect Transcript Stability and Translational Efficiency. J. Mol. Biol. 177, 107-136.

©2012 Life Technologies Corporation. All rights reserved.

The trademarks mentioned herein are the properties of Life Technologies Corporation or their respective owners.

Zeocin is a trademark of CAYLA, SA.

Notes

Notes





www.lifetechnologies.com