

# EasySelect<sup>™</sup> Echo<sup>™</sup>-Adapted *Pichia* Expression Kit

For expression of the gene of interest in *Pichia pastoris* using pPICZ-E and pPICZα-E with the Echo<sup>™</sup> Cloning System

Catalog nos. ET230-xx, ET231-xx, ET232-xx

**Version E** 07 September 2010 *25-0388* 

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

**User Manual** 

### INDIVIDUAL PICHIA EXPRESSION KIT LICENSE AGREEMENT

The *Pichia* Expression Kit is based on the yeast *Pichia pastoris*. *Pichia pastoris* was developed into an expression system by scientists at Salk Institute Biotechnology/Industry Associates (SIBIA) for high-level expression of recombinant proteins. All patents for *Pichia pastoris* and licenses for its use as an expression system are owned by Research Corporation Technologies, Inc. Tucson, Arizona. Invitrogen has an exclusive license to sell the *Pichia* Expression Kit to scientists for research purposes only, under the terms described below. Use of *Pichia pastoris* by commercial corporations requires the user to obtain a commercial license as detailed below. Before using the *Pichia* Expression Kit, please read the following license a greement. If you do not agree to be bound by its terms, contact Invitrogen within 10 days for authorization to return the unused *Pichia* Expression Kit and to receive a full credit. If you do agree to the terms of this Agreement, please complete the User Registration Card and return it to Invitrogen before using the kit.

#### INDIVIDUAL PICHIA EXPRESSION KIT LICENSE AGREEMENT

Invitrogen Corporation (INVITROGEN) grants you a non-exclusive license to use the enclosed *Pichia* Expression Kit (EXPRESSION KIT) for academic research or for evaluation purposes only. The EXPRESSION KIT is being transferred to you in furtherance of, and reliance on, such license. You may not use the EXPRESSION KIT, or the materials contained therein, for any commercial purpose without a license for such purpose from RESEARCH CORPORATION TECHNOLOGIES, INC., Tucson, Arizona. Commercial purposes include the use in or sale of expressed proteins as a commercial product, or use to facilitate or advance research or development of a commercial product. Commercial entities may conduct their evaluation for one year at which time this license automatically terminates. Commercial entities will be contacted by Research Corporation Technologies during the evaluation period regarding the purchase of a commercial license.

Access to the EXPRESSION KIT must be limited solely to those officers, employees and students of your institution who need access thereto in order to perform the above-described research or evaluation. You must inform each of such officer, employee and student of the provisions of this Agreement and require them to agree, in writing, to be bound by the provisions of this Agreement. You may not distribute the EXPRESSION KIT to others, even those within your own institution. You may transfer modified, altered or original material from the EXPRESSION KIT to a third party following notification of INVITROGEN such that the recipient can be licensed. You may not assign, sub-license, rent lease or otherwise transfer this License or any of the rights or obligation hereunder, except as expressly permitted.

This License is effective until terminated. You may terminate it at any time by destroying all *Pichia* expression products in your control. It will also terminate automatically if you fail to comply with the terms and conditions of the Agreement. You shall, upon termination of the License, destroy all *Pichia* Expression Kits in your control, and so notify INVITROGEN in writing.

This License Shall be governed in its interpretation and enforcement by the laws of the State of California..

#### **Technical Services**

Invitrogen provides Technical Services to all of our registered *Pichia* Expression Kit users. Please contact us if you need assistance with the *Pichia* Expression Kit.

#### **Corporate Headquarters:** Invitrogen Corporation 1600 Faraday Avenue Carlsbad, CA 92008 USA Tel: 1 760 603 7200 Tel (Toll Free): 1 800 955 6288 Fax: 1 760 602 6500 E-mail: tech\_service@invitrogen.com

#### Japanese Headquarters

Invitrogen Japan LOOP-X Bldg. 6F 3-9-15, Kaigan Minato-ku, Tokyo 108-0022 Tel: 81 3 5730 6509 Fax: 81 3 5730 6519 E-mail: jpinfo@invitrogen.com

#### **European Headquarters:**

Invitrogen Ltd 3 Fountain Drive Inchinnan Business Park 3 Fountain Drive Paisley PA4 9RF, UK Tel: +44 (0) 141 814 6100 Fax: +44 (0) 141 814 6287 E-mail: eurotech@invitrogen.com

## **Table of Contents**

Table of Contents	V
Kit Contents and Storage	vi
Product Qualification	xi
Accessory Products	xiii
Purchaser Notification	xv
Using This Manual	xviii
Introduction	1
Overview of the Echo <sup>™</sup> Cloning System	1
Overview of Pichia pastoris Expression System	4
Methods	8
Recombining Your Gene into pPICZ-E or pPICZα-E	
Transforming the Recombination Reaction	9
Pichia Strains	
Pichia Transformation	
EasyComp <sup>™</sup> Transformation	
Determining the Mut Phenotype	
Expression in Pichia	
Analysis by SDS-Polyacrylamide Gel Electrophoresis	
Optimization of Pichia Protein Expression	
Scale-up of Expression	
Purification	
Pichia Media Recipes	
Appendix	45
Recipes	45
Maps of pPICZ-E and pPICZα-E	46
Features of pPICZ-E and pPICZ $\alpha$ -E	
Map of pPICZ-E/Uni-lacZ	
Map of pPICZα-E/Uni-HSA	
Recombination and Integration in Pichia	51
Zeocin <sup>™</sup>	53
Direct PCR Screening of Pichia Clones	
Total DNA Isolation from <i>Pichia</i>	
Determination of Copy Number of Multiple Integrants	
Procedure for Total RNA Isolation from Pichia	60
Technical Service	61
References	

## Kit Contents and Storage

### **Types of Kits**

Several EasySelect<sup>TM</sup> Echo<sup>TM</sup>-adapted *Pichia* Expression Kits, pPICZ-E and pPICZ $\alpha$ -E Echo<sup>TM</sup>-adapted Expression Vector Kits are available (see table below). This manual is included with all of these kits. Note that not all kits contain the reagents discussed in this manual (see page xviii for more details).

Kit	Reagents Supplied	Catalog nos.
EasySelect <sup>TM</sup> Echo <sup>TM</sup> -adapted <i>Pichia</i>	pPICZ-E Echo <sup>™</sup> -adapted Expression Vector Kit	ET230-02
Expression Kit	pPICZ $\alpha$ -E Echo <sup>TM</sup> -adapted Expression Vector Kit	
	EasySelect <sup>™</sup> <i>Pichia</i> Strains Kit	
	<i>Pichia</i> EasyComp <sup>™</sup> Kit	
	Pichia Media Kit	
EasySelect <sup>™</sup> Echo <sup>™</sup> -adapted <i>Pichia</i>	pUni/V5-His-TOPO® TA Cloning Kit	ET230-10C
Expression Kit with a choice of	pUniBlunt/V5-His-TOPO <sup>®</sup> Cloning Kit	ET230-20C
Donor Vector Kit and One Shot <sup>®</sup> TOP10 Chemically Competent	pUni/V5-His A, B and C	ET230-30C
<i>E. coli</i> (see page xiii for more information on donor vectors)	pUniD/V5-His-TOPO <sup>®</sup> Cloning Kit	ET230-40C
pPICZ-E Echo <sup>™</sup> -adapted Expression	pPICZ-E Vector	ET231-01
Vector Kit	Expression Control Vector	
	Cre Recombinase and 10X Buffer	
	5' AOX1 Sequencing Primer	
	Zeocin <sup>TM</sup>	
pPICZα-E Echo <sup>™</sup> -adapted	pPICZ $\alpha$ -E Vector	ET232-01
Expression Vector Kit	Expression Control Vector	
	Cre Recombinase and 10X Buffer	
	α-Factor Sequencing Primer	
	Zeocin <sup>™</sup>	
pPICZ-E Echo <sup>™</sup> -adapted Expression	pUni/V5-His-TOPO® TA Cloning Kit	ET231-10C
Vector Kit with a choice of Donor Vector Kit and One Shot <sup>®</sup> TOP10	pUniBlunt/V5-His-TOPO® Cloning Kit	ET231-20C
Chemically Competent E. coli (see	pUni/V5-His A, B and C	ET231-30C
page xiii for more information on donor vectors)	pUniD/V5-His-TOPO <sup>®</sup> Cloning Kit	ET231-40C
pPICZα-E Echo <sup>™</sup> -adapted	pUni/V5-His-TOPO® TA Cloning Kit	ET232-10C
Expression Vector Kit with a choice	pUniBlunt/V5-His-TOPO <sup>®</sup> Cloning Kit	ET232-20C
of Donor Vector Kit and One Shot <sup>®</sup> TOP10 Chemically Competent	pUni/V5-His A, B and C	ET232-30C
<i>E. coli</i> (see page xiii for more information on donor vectors)	pUniD/V5-His-TOPO <sup>®</sup> Cloning Kit	ET232-40C

#### The EasySelect<sup>™</sup> Echo<sup>™</sup>-adapted *Pichia* Expression Kit is shipped on dry ice except for Shipping and the EasySelect<sup>TM</sup>*Pichia* strains and the *Pichia* Media Kit, which are shipped at room Storage temperature. Each EasySelect<sup>TM</sup> Echo<sup>TM</sup>-adapted *Pichia* Expression Kit contains two vector kits (1 box for each vector), EasySelect<sup>TM</sup>Pichia Strains (Box 2), EasyComp<sup>TM</sup> Pichia Kit (Box 3), and Pichia Media Kit.

The pPICZ-E and pPICZ $\alpha$ -E Echo<sup>TM</sup>-adapted Expression Kits are shipped on dry ice. Each of pPICZ-E and pPICZ $\alpha$ -E Echo<sup>TM</sup>-Adapted Expression Kit contains the Echo<sup>TM</sup>adapted expression vector, an expression control vector, a sequencing primer and Zeocin<sup>™</sup>.

Reagents	Storage
pPICZ-E Echo <sup>™</sup> -adapted Expression Kit	-20°C
pPICZα-E Echo <sup>™</sup> -adapted Expression Kit	-20°C
EasySelect <sup>™</sup> <i>Pichia</i> Strains Kit	+4°C
Pichia EasyComp <sup>™</sup> Kit	+4°C
Pichia Media Kit	Room temperature
One Shot <sup>®</sup> TOP10 Chemically Competent <i>E. coli</i> (Optional)	-80°C

# pPICZ-E Echo<sup>™</sup>

The following items are supplied in the pPICZ-E Echo<sup>™</sup>-adapted Expression Kit.

# Reagents

Item	Composition	Amount
pPICZ-E	Supercoiled, lyophilized in TE, pH 8.0	20 µg
pPICZ-E/Uni- <i>lacZ</i> expression control	Supercoiled, lyophilized in TE, pH 8.0	20 µg
5' AOX1 sequencing primer	Lyophilized in water	2 µg
Cre Recombinase	Check the label on the tube for exact concentration of the enzyme.	15 µl
	Enzyme is supplied in:	
	50 mM Tris-HCl, pH 8.0	
	5 mM EDTA	
	1 mM EGTA	
	10 mM β-mercaptoethanol	
	20% Glycerol	
10X Recombinase Buffer	500 mM Tris-HCl, pH 7.5	25 μl
	100 mM MgCl <sub>2</sub>	
	300 mM NaCl	
	1.0 mg/ml BSA	
Zeocin <sup>™</sup>	100 mg/ml	2 x 1.25 ml

### pPICZα-E Echo<sup>™</sup> Reagents

The following items are supplied in the pPICZ $\alpha$ -E Echo<sup>TM</sup>-adapted Expression Kit

Item	Composition	Amount
pPICZa-E	Supercoiled, lyophilized in TE, pH 8.0	20 µg
pPICZα-E/Uni-HSA expression control	Supercoiled, lyophilized in TE, pH 8.0	20 µg
$\alpha$ -Factor sequencing primer	Lyophilized in water	2 µg
Cre Recombinase	Check the label on the tube for exact concentration of the enzyme.	15 μl
	Enzyme is supplied in:	
	50 mM Tris-HCl, pH 8.0	
	5 mM EDTA	
	1 mM EGTA	
	10 mM $\beta$ -mercaptoethanol	
	20% Glycerol	
10X Recombinase Buffer	500 mM Tris-HCl, pH 7.5	25 μl
	100 mM MgCl <sub>2</sub>	
	300 mM NaCl	
	1.0 mg/ml BSA	
Zeocin <sup>TM</sup>	100 mg/ml	2 x 1.25 ml

### The *Pichia* EasyComp<sup>™</sup> Kit

This kit contains sufficient reagents for 6 preparations of competent cells. Each competent cell preparation yields enough cells for 20 transformations.

Upon receipt, store the kit at +4°C. Solution II can also be stored at room temperature.

Component	Description	Quantity
Solution I	Sorbitol solution containing ethylene glycol and DMSO for the preparation of competent cells	75 ml
Solution II	PEG solution for the transformation of competent cells	150 ml (2 x 75 ml)
Solution III	Salt solution for washing and plating transformed cells	150 ml (2 x 75 ml)

# EasySelect<sup>™</sup>*Pichia* Strains Kit

The table below lists the genotype and phenotype of the different *Pichia* stabs. Store the stabs at  $+4^{\circ}$ C.

Strain	Genotype	Phenotype (Pichia only)
X-33	wild-type	Mut <sup>+</sup>
GS115	his4	His <sup>-</sup> , Mut <sup>+</sup>
KM71H	arg4 aox1::ARG4	Mut <sup>S</sup> , Arg <sup>+</sup>
GS115/pPICZ/lacZ	his4	His <sup>-</sup> , Mut <sup>+</sup>
GS115/Albumin	his4	His <sup>-</sup> , Mut <sup>s</sup>
TOP10F	F´ { <i>proAB</i> , <i>lacI</i> <sup>q</sup> , <i>lacZ</i> ΔM15, Tn10 (Tet <sup>R</sup> )} <i>mcrA</i> , Δ( <i>mrr-hsd</i> RMS- <i>mcrBC</i> ), $\phi$ 80 <i>lacZ</i> ΔM15, Δ <i>lacX</i> 74, , <i>recA</i> 1, λ- <i>ara</i> D139, Δ( <i>ara-leu</i> )7697, <i>gal</i> U, <i>gal</i> K, <i>rpsL</i> (Str <sup>R</sup> ), <i>end</i> A1, <i>nup</i> G	

#### Media

The following prepackaged media is included for your convenience. Instructions for use are provided on the package. Keep the media dry and store at room temperature.

Media	Amount	Yield
YP Base Medium	2 pouches	2 liters of YP medium
YP Base Agar Medium	2 pouches	2 liters of YP agar medium
Yeast Nitrogen Base	1 pouch	500 ml of 10X YNB

#### Sequence of Primers

The table below lists the sequence and pmoles of the primers included in this kit.

Primer	Sequence	pmoles Supplied
5' AOX1 sequencing primer	5'-GACTGGTTCCAATTGACAAGC-3'	312 pmoles
α-Factor sequencing primer	5'-TACTATTGCCAGCAATTGCTGC-3'	315 pmoles

#### One Shot<sup>®</sup> TOP10 Reagents (Optional)

The table below describes the items included in the One Shot<sup>®</sup> TOP10 Chemically Competent *E. coli* kit.

Store at -80°C.

Item	Concentration	Amount
SOC Medium	2% Tryptone	6 ml
(may be stored at room	0.5% Yeast Extract	
temperature or at $+4^{\circ}C$ )	10 mM NaCl	
+4 C)	2.5 mM KCl	
	10 mM MgCl <sub>2</sub>	
	10 mM MgSO <sub>4</sub>	
	20 mM glucose	
TOP10 E. coli		11 x 50 µl
pUC19 Control DNA	10 pg/µl in 5 mM Tris-HCl, 0.5 mM EDTA, pH 8	50 µl

## Genotype of TOP10

**TOP10:** Use this strain for general cloning of your gene of interest. Note that this strain cannot be used for transformation and growth of the donor vectors.

F- mcrA  $\Delta$ (mrr-hsdRMS-mcrBC)  $\Phi$ 80lacZ $\Delta$ M15  $\Delta$ lacX74 recA1 araD139  $\Delta$ (ara-leu)7697 galU galK rpsL (Str<sup>R</sup>) endA1 nupG

## **Product Qualification**

## **Vectors** pPICZ-E, $pPICZ\alpha-E$ and the control plasmids are qualified by restriction digest. The table below lists the restriction enzymes and the expected fragments.

Restriction pPICZ-E pPICZ-E/Uni-lacZ pPICZa-E/Uni-HSA pPICZa-E Enzyme 1403 bp, 2042 bp Bgl I 1777 bp, 1403 bp Not tested Not tested EcoR I (linearizes) 3180 bp Not tested 3445 bp Not tested Hind III Not tested 3404 bp, 5148 bp Not tested 3404 bp, 4108 bp Not I Not tested 2133 bp, 6419 bp Not tested 2133 bp, 5379 bp

Primers	The sequencing primers are lot-qualified by DNA sequencing experiments using the dideoxy chain termination technique.
Cre Recombinase	Purity: >95% homogeneity
	Endonuclease activity: Negative
	Exonuclease activity: Negative
	Functional Assay: Cre recombinase is qualified using the assay on page 8 of this manual. The donor vector used is pUni/ <i>lacZ</i> and the acceptor vector is pcDNA3.1-E. Five microliters of the recombination reaction is transformed into 50 $\mu$ l One Shot <sup>®</sup> TOP10 Chemically Competent <i>E. coli</i> using the protocol on page 9. Twenty-five $\mu$ l of the transformation reaction is plated on LB plates containing 50 $\mu$ g/ml kanamycin (performed in duplicate). One microliter of Cre recombinase should yield > 500 blue, kanamycin-resistant transformants.
<i>Pichia</i> EasyComp <sup>™</sup> Kit	The <i>Pichia</i> EasyComp <sup>TM</sup> Kit is qualified by preparation of competent GS115 <i>Pichia</i> cells. 50 $\mu$ l of competent GS115 cells are transformed with 3 $\mu$ g of linearized pPICZ $\alpha$ A DNA.
	Transformation efficiency must be greater than 50 cfu/ $\mu$ g DNA.
	All buffers and solutions are tested for sterility.
<i>Pichia</i> strains	Each <i>Pichia</i> strain is qualified by recovery on YPD medium.

## Product Qualification, Continued

Zeocin <sup>™</sup>	Zeocin <sup>TM</sup> is lot-qualified by demonstrating that Low Salt LB medium containing 25 $\mu$ g/ml Zeocin <sup>TM</sup> prevents growth of the <i>E. coli</i> strain, TOP10.
Growth and Expression Media	<i>Pichia</i> growth and expression media are qualified by the ability to support growth of the GS115 <i>Pichia</i> strain.
One Shot Competent <i>E. coli</i>	<ul> <li>All competent cells are qualified as follows:</li> <li>Cells are tested for transformation efficiency using the control plasmid included in the kit. Transformed cultures are plated on LB plates containing 100 µg/ml ampicillin and the transformation efficiency is calculated. Test transformations are performed in duplicate. Transformation efficiency should be ~1 x 10<sup>9</sup> cfu/µg DNA for chemically competent cells and &gt;1 x 10<sup>9</sup> for electrocompetent cells.</li> <li>To verify the absence of phage contamination, 0.5-1 ml of competent cells are added to LB top agar and poured onto LB plates. After overnight incubation, no</li> </ul>
	<ul> <li>plaques should be detected.</li> <li>Untransformed cells are plated on LB plates 100 μg/ml ampicillin, 25 μg/ml streptomycin, 50 μg/ml kanamycin, or 15 μg/ml chloramphenicol to verify the absence of antibiotic-resistant contamination.</li> </ul>

## **Accessory Products**

#### Additional Products

Many of the reagents in the EasySelect<sup>TM</sup> Echo<sup>TM</sup>-adapted *Pichia* Expression Kit and pPICZ-E or pPICZ $\alpha$ -E Echo<sup>TM</sup>-adapted Expression Vector Kits, as well as additional reagents that may be used with these kits, are available separately from Invitrogen. Ordering information is provided below.

Product	Amount	Catalog no.
One Shot <sup>®</sup> PIR1 Chemically Competent E.coli	11 x 50 μl	C1010-10
One Shot <sup>®</sup> PIR2 Chemically Competent <i>E</i> .coli	11 x 50 μl	C1111-10
One Shot <sup>®</sup> TOP10 Chemically Competent <i>E. coli</i>	11 x 50 μl	C4040-10
Cre Recombinase	10 reactions	R100-10
Kanamycin	5 g	11815-024
	25 g	11815-032
Zeocin <sup>™</sup>	1 g	R250-01
	5 g	R250-05
Pichia EasyComp <sup>™</sup> Transformation Kit	1 kit	K1730-01
Pichia Protocols	1 book	G100-01
X-33 Pichia strain	1 stab	C180-00
KM71H Pichia strain	1 stab	C182-00
SMD1168H Pichia strain	1 stab	C184-00
GS115 Pichia strain	1 stab	C181-00

### **Donor Vectors**

The table below lists a variety of donor vectors currently available from Invitrogen to facilitate cloning of your gene of interest for use with Echo<sup>™</sup> Cloning System.

Product	Application	Quantity	Catalog no.
pUniD/V5-His-TOPO <sup>®</sup> Cloning Kit	Directional cloning of blunt PCR products	10 reactions	ET004-10
pUni/V5-His-TOPO <sup>®</sup> TA Cloning Kit	Cloning A-tailed PCR products	10 reactions	ET001-10
pUniBlunt/V5-His-TOPO <sup>®</sup> Cloning Kit	Cloning blunt end products	10 reactions	ET002-10
pUni/V5-His A, B, and C	Cloning DNA fragments using restriction enzymes	10 reactions	ET003-10

## **Accessory Products, Continued**

#### Detection of Fusion Protein

A number of antibodies and immunodetection kits are available from Invitrogen to detect expression of your fusion protein from the pPICZ-E or pPICZ $\alpha$ -E vectors. Horseradish peroxidase (HRP)-or alkaline phosphatase (AP)-conjugated antibodies allow one-step detection in western blots using colorimetric or chemiluminescent detection methods. Sufficient antibody is provided for 25 westerns. The WesternBreeze<sup>TM</sup> kit contains enough reagents for 20 blots.

Product	Application	Catalog no.
Anti-V5 Antibody	Detects the 14 amino acid epitope derived	R960-25
Anti-V5-HRP Antibody	from the P and V proteins of the paramyxovirus, SV5 (Southern et al., 1991)	R961-25
Anti-V5-AP Antibody	GKPINPLLGLDST	R962-25
Anti-His(C-term) Antibody	Detects the C-terminal polyhistidine (6xHis)	R930-25
Anti-His(C-term)-HRP Antibody	detection) (Lindner et al., 1997):	R931-25
Anti-His(C-term)-AP Antibody		R932-25
WesternBreeze <sup>™</sup> Chromogenic kits-αMouse	Chromogenic detection of proteins following Western transfer with an alkaline phosphatase substrate	WB7103
WesternBreeze <sup>™</sup> Chemiluminescent kits- αMouse	Chemiluminescent detection of proteins following Western transfer with an alkaline phosphatase substrate	WB7104

### Purification of Fusion Protein

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

Product	Quantity	Catalog no.
ProBond <sup>™</sup> Purification System	6 purifications	K850-01
ProBond <sup>™</sup> Resin	50 ml	R801-01
	150 ml	R801-15
Purification Columns	50 polypropylene columns	R640-50

### **Purchaser Notification**

### Limited Use Label License No. 74: Pichia Pastoris

The Pichia Expression System is based on the yeast Pichia pastoris. Pichia pastoris was developed into an expression system by scientists at Salk Institute Biotechnology/ Industry Associates (SIBIA) and Phillips Petroleum for high-Expression System level expression of recombinant proteins. All patents for Pichia pastoris and licenses for its use as an expression system are owned by Research Corporation Technologies (RCT), Inc., Tucson, Arizona. Life Technologies has an exclusive license to sell Pichia expression kits and vectors to scientists for research purposes only, under the terms described below. Use of Pichia pastoris by commercial entities for any commercial purpose requires the user to obtain a commercial license as detailed below. Before using any Pichia expression product, please read the following license agreement. If you do not agree to be bound by its terms, contact Life Technologies within 10 days for authorization to return the unused Pichia expression products and to receive a full refund. If you do agree to the terms of this license agreement, please complete the User Registration Card and return it to Life Technologies before using the product. Life Technologies Corporation ("Life Technologies") grants you a nonexclusive license to use the enclosed Pichia expression vectors ("Expression Vector") for academic research or for evaluation purposes only. The Expression Vectors are being transferred to you in furtherance of, and reliance on, such license. You may not use the Expression Vectors for any commercial purpose without a license for such purpose from Research Corporation Technologies, Inc., Tucson, Arizona.

Commercial purposes include: any use of Expression Products or Expression Vectors in a Commercial Product; any use of Expression Products or Expression Vectors in the manufacture of a Commercial Product; any sale of Expression Products; any use of Expression Products or the Expression Kit to facilitate or advance research or development directed to a Commercial Product; and any use of Expression Products or the Expression Kit to facilitate or advance any research or development program the results of which will be directly applied to the development or manufacture of a Commercial Product. "Expression Products" means products expressed with the Expression Kit, or with the use of any Pichia expression vectors (including the Expression Vector) or host strains. "Commercial Product" means any product intended for sale or commercial use. Commercial entities may conduct their evaluation for one year at which time this license automatically terminates. Commercial entities will be contacted by Research Corporation Technologies during the evaluation period regarding their desire for a commercial license.

## **Purchaser Notification, Continued**

Limited Use Label License No. 74: Pichia Pastoris Expression System, continued	Commercial entities may conduct their evaluation for one year at which time this license automatically terminates. Commercial entities will be contacted by Research Corporation Technologies during the evaluation period regarding their desire for a commercial license. Access to the Expression Kit and Vector must be limited solely to those officers, employees and students of your institution who need access to per- form the above-described research or evaluation. You must inform each such officer, employee and student of the provisions of this license agreement and require them to agree, in writing, to be bound by the provisions of this license agreement. You may not distribute any Expression Vector or host strain contained herein or in the Expression Kit to others, even those within your own institution. You may only transfer modified, altered, or original material from the Expression Kit or Vector to a third party following written notification of, and written approval from, Life Technologies so that the recipient can be licensed. You may not assign, sub-license, rent, lease or otherwise transfer this license agreement or any of the rights or obligation there under, except as expressly permitted by Life Technologies and RCT. This license agreement is effective until terminated. You may terminate it at any time by destroying all Pichia Expression products in your control. It will also terminate automatically if you fail to comply with the terms and conditions of the license agreement. You shall, upon termination of the license agreement, destroy all Pichia Expression products in your control, and so notify Life Technologies in writing. You may contact Research Corporation Technologies, 101 North Wilmot Road, Suite 600, Tucson, Arizona 85711-3335. Tel: 520-748-4443, Fax: 520-748-0025.
Limited Use Label License No: 22 Vectors and Clones Encoding Histidine Hexamer	This product is licensed under U.S. Patent Nos. 5,284,933 and 5,310,663 and foreign equivalents from Hoffmann-LaRoche, Inc., Nutley, NJ and/or Hoffmann-LaRoche Ltd., Basel, Switzerland and is provided only for use in research. Information about licenses for commercial use is available from QIAGEN GmbH, Max-Volmer-Str. 4, D-40724 Hilden, Germany.

### **Purchaser Notification, Continued**

Limited Use Label License No: 119 Echo™ Cloning Products No license is conveyed to use this product with any recombination sites other than those purchased from Life Technologies Corporation or its authorized distributor. The buyer cannot modify the recombination sequence(s) contained in this product for any purpose.

#### Limited Use Label License No: 5 Invitrogen Technology

The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. The buyer may transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for any Commercial Purpose, and that such collaborator agrees in writing (a) not to transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes. Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. For products that are subject to multiple limited use label licenses, the terms of the most restrictive limited use label license shall control. Life Technologies Corporation will not assert a claim against the buyer of infringement of patents owned or controlled by Life Technologies Corporation which cover this product based upon the manufacture, use or sale of a therapeutic, clinical diagnostic, vaccine or prophylactic product developed in research by the buyer in which this product or its components was employed, provided that neither this product nor any of its components was used in the manufacture of such product. If the purchaser is not willing to accept the limitations of this limited use statement, Life Technologies is willing to accept return of the product with a full refund. For information about purchasing a license to use this product or the technology embedded in it for any use other than for research use please contact Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008; Phone (760) 603-7200 or e-mail: outlicensing@lifetech.com.

## **Using This Manual**



This manual is a comprehensive manual designed to support a variety of different *Pichia* kits. Some reagents discussed in this manual may not be included in the kit you purchased. In addition, you may have your own methods for transformation and expression in *Pichia*. Use the information from this manual according to your needs.

## Introduction

# Overview of the Echo<sup>™</sup> Cloning System

Introduction	The Echo <sup>TM</sup> Cloning System allows direct recombination of your gene of interest downstream of an appropriate promoter for expression in the host system of choice. pPICZ-E and pPICZ $\alpha$ E are members of the Echo <sup>TM</sup> Cloning System family of expression vectors and are specifically designed for expression in <i>Pichia pastoris</i> . The 5'AOX1 promoter controls high-level inducible recombinant protein expression in any <i>Pichia pastoris</i> strain of choice.
The Echo <sup>™</sup> Cloning System	The Echo <sup>TM</sup> Cloning System is based on the univector plasmid-fusion system (UPS) des- cribed by Elledge and coworkers to quickly and easily recombine a gene of interest into a series of recipient (acceptor) vectors (Liu et al., 1998; Liu et al., 1999). The system con- sists of the univector (donor) vector containing the gene of interest and recipient (acceptor) vector containing various regulatory sequences for expression in the host of choice. The Echo <sup>TM</sup> System utilizes the <i>cre-lox</i> site-specific recombination system of bacteriophage P1 (Abremski et al., 1983; Sternberg et al., 1981). The product of the <i>cre</i> gene is a site-specific recombinase that catalyzes conservative recombination between two 34 bp <i>loxP</i> sequences or, a <i>loxP</i> and a <i>loxH</i> sequence to resolve P1 dimers generated by replication of circular lysogens. It does not catalyze recombination between two <i>loxH</i> sequences.
Plasmid Fusion	The donor (pUni) vector and the acceptor vector (i.e. pPICZ-E or pPICZ $\alpha$ -E) each contain a single <i>lox</i> site. The donor vector and the acceptor vector each contain a <i>loxP</i> site. You may insert your gene of interest into the donor vector via the TOPO <sup>®</sup> Cloning method or traditional restriction enzyme-mediated cloning. pPICZ-E and pPICZ $\alpha$ -E contain the appropriate transcription regulatory sequences to control expression of your gene of interest in <i>Pichia</i> and a unique <i>loxP</i> site located downstream of these sequences. By mixing the donor vector containing the gene of interest with pPICZ-E or pPICZ $\alpha$ -E in the presence of Cre recombinase, a plasmid fusion is created that expresses the gene of interest in <i>Pichia pastoris</i> . A generic diagram is shown below.
	pUni (2.3 kb + gene) B <sup>fc</sup> on (xx <sup>2</sup> gene Recombinant

Cre

recombinase

pAcceptor (2.5 to 5.8 kb)

Amp<sup>ℝ</sup>

lox\* = loxP or loxH depending on acceptor vector

IOXP

Plasmid

(4.8 kb + gene to 8.1 kb + gene)

Amp<sup>R</sup>

# Overview of Echo<sup>™</sup> Cloning System, Continued

<i>loxP</i> or <i>loxH</i> Sites	The sequence of the <i>loxP</i> site is shown below. The <i>loxP</i> site consists of a 34 bp sequence containing two 13 bp inverted repeats (see underlined bases) separated by an 8 bp spacer (Hoess et al., 1982). The inverted repeats may form a stem and loop structure that may reduce expression of the gene of interest in some cases. A variation of the <i>loxP</i> site ( <i>loxH</i> , see below) was created to eliminate the formation of a stem and loop structure and improve expression. We have not observed any differences in expression levels in constructs containing a <i>loxP</i> or a <i>loxH</i> site. Mutated bases are shown in boldface. Cre-mediated recombination can still occur between a <i>loxP</i> and a <i>loxH</i> site although the efficiency may be slightly reduced.
	• <i>loxP</i> : <u>ATA ACT TCG TAT A</u> GC ATA CAT <u>TAT ACG AAG TTA T</u>
	• <i>loxH</i> : ATT ACC TCA TAT AGC ATA CAT TAT ACG AAG TTA T
Cre Recombinase	Cre recombinase (MW = 35 kDa) is a site-specific recombinase that binds to specific sequences ( <i>loxP</i> and <i>loxH</i> sites), brings together the target sites, cleaves them, and covalently attaches to the DNA. Recombination occurs following two pairs of strand exchanges and ligation of the DNAs in a novel (recombinant) form. A nucleophilic hydroxylated tyrosine initiates the DNA cleavage event by attack on a specific phosphodiester bond followed by the covalent attachment of the recombinase to the target sequence through a phosphoamino acid bond (Abremski & Hoess, 1992; Argos et al., 1986). The reaction does not require any host factors or ATP, but does require Mg <sup>2+</sup> or spermidine for activity (Abremski et al., 1983). Recombination between two supercoiled substrates, each containing a <i>loxP</i> or <i>loxH</i> site, results in a supercoiled dimer. The extent of the reaction is 10-20% and appears to be stoichiometric (Abremski & Hoess, 1984; Abremski et al., 1983).
Selection of Recombinants	By fusing the two plasmids, kanamycin resistance is now linked to the pUC origin of replication. The recombination reaction is transformed into TOP10 <i>E. coli</i> and recombinants are selected by plating the transformation reaction onto plates containing kanamycin. Because the donor plasmid carries the R6K $\gamma$ origin of replication, it will not propagate in <i>E. coli</i> strains such as TOP10 which do not carry the <i>pir</i> gene. In addition, the acceptor vector, which carries the ampicillin resistance gene will not be selected. Therefore every colony that is selected on kanamycin will represent a recombined fusion plasmid.

# Overview of Echo<sup>™</sup> Cloning System, Continued

pPICZ-E and pPICZα-E	pPICZ-E and pPICZ $\alpha$ -E (~ 3 kb) are derived from pPICZB and pPICZ $\alpha$ -B, respectively. They are designed for high-level recombinant protein expression in <i>Pichia pastoris</i> . The vectors contain the following elements:
	• 5' fragment containing the <i>AOX</i> 1 promoter for tightly regulated, methanol-induced expression of the gene of interest (Ellis <i>et al.</i> , 1985: Koutz <i>et al.</i> , 1989: Tschopp <i>et al.</i> , 1987a)
	• A <i>loxP</i> site for plasmid fusion
	• <i>Saccharomycese cerevisiae</i> α-factor secretion signal sequence (pPICZα-E) for secretion of recombinant proteins in the medium
	• Zeocin <sup>™</sup> resistance gene for selection in both <i>E. coli</i> and <i>Pichia</i> (Baron <i>et al.</i> , 1992: Drocourt <i>et al.</i> , 1990)
	• The pUC origin for high copy replication and maintenance of the plasmid in <i>E. coli</i>
	For a map and a description of the features of pPICZ-E or pPICZ $\alpha$ -E, refer to the <b>Appendix</b> , pages 46-48.
	Other Echo <sup>™</sup> -adapted acceptor vectors are available separately and are provided with their own manuals. For more information on other available acceptor vectors, visit our Web site (www.invitrogen.com) or call Technical Service (see page 61).
Selection of Vector	To utilize the strong, highly-inducible $P_{AOXI}$ promoter (see next page) for expression of your protein, there are two expression vectors included in this kit. One vector, pPICZ-E, is for intracellular expression while the other vector, pPICZ $\alpha$ -E, is for secreted expression. All vectors contain the Zeocin <sup>TM</sup> resistance gene for positive selection in <i>E. coli</i> and <i>Pichia</i> . See pages 46-48 for more information on these vectors.

## Overview of Pichia pastoris Expression System

Review Articles	The information presented here is designed to give you a concise overview of the <i>Pichia pastoris</i> expression system. It is by no means exhaustive. For further information, read the articles cited in the text along with recent review articles (Higgins & Cregg, 1998), (Buckholz & Gleeson, 1991; Cregg & Higgins, 1995; Cregg et al., 1993; Nico-Farber et al., 1995; Sreekrishna et al., 1988; Wegner, 1990). A general review of foreign gene expression in yeast is also available (Romanos et al., 1992)
General Characteristics of <i>Pichia pastoris</i>	As a eukaryote, <i>Pichia pastoris</i> has many of the advantages of higher eukaryotic expression systems such as protein processing, protein folding, and posttranslational modification, while being as easy to manipulate as <i>E. coli</i> or <i>Saccharomyces cerevisiae</i> . It is faster, easier, and less expensive to use than other eukaryotic expression systems such as baculovirus or mammalian tissue culture, and generally gives higher expression levels. As a yeast, it shares the advantages of molecular and genetic manipulations with <i>Saccharomyces</i> and has the added advantage of 10- to 100-fold higher heterologous protein expression levels. These features make <i>Pichia</i> very useful as a protein expression system.
Similarity to Saccharomyces	Many of the techniques developed for <i>Saccharomyces</i> may be applied to <i>Pichia</i> including transformation by complementation, gene disruption, and gene replacement. In addition, the genetic nomenclature used for <i>Saccharomyces</i> has been applied to <i>Pichia</i> . For example, histidinol dehydrogenase is encoded by the <i>HIS4</i> gene in both <i>Saccharomyces</i> and <i>Pichia</i> . There is also cross-complementation between gene products in both <i>Saccharomyces</i> and <i>Pichia</i> . Several wild-type genes from <i>Saccharomyces</i> complement comparable mutant genes in <i>Pichia</i> . Genes such as <i>HIS4</i> , <i>LEU2</i> , <i>ARG4</i> , <i>TRP1</i> , and <i>URA3</i> all complement their respective mutant genes in <i>Pichia</i> .
<i>Pichia pastoris</i> as a Methylotrophic Yeast	<i>Pichia pastoris</i> is a methylotrophic yeast, capable of utilizing methanol as its sole carbon source. Alcohol oxidase catalyses the oxidation of methanol to formaldehyde and hydrogen peroxide using molecular oxygen. To avoid hydrogen peroxide toxicity, methanol metabolism takes place within a specialized cell organelle, the peroxisome, which sequesters toxic by-products away from the rest of the cell. Alcohol oxidase has a poor affinity for $O_2$ , and <i>Pichia</i> compensates by generating large amounts of the enzyme. The promoter regulating the production of alcohol oxidase is used to drive recombinant protein expression in <i>Pichia</i> .
Two Alcohol Oxidase Proteins	Two genes in <i>Pichia pastoris</i> code for alcohol oxidase- <i>AOX1</i> and <i>AOX2</i> . The majority of alcohol oxidase activity in the cell is attributable to the product of the <i>AOX1</i> gene. Expression of the <i>AOX1</i> gene is tightly regulated and induced by methanol to very high levels, typically $\geq$ 30% of the total soluble protein in cells grown with methanol. The <i>AOX1</i> gene has been isolated and a plasmid-borne version of the <i>AOX1</i> promoter is used to drive expression of the gene of interest (Ellis et al., 1985; Koutz et al., 1989; Tschopp et al., 1987a). While <i>AOX2</i> is about 97% homologous to <i>AOX1</i> , growth on methanol is much slower than with <i>AOX1</i> . This slow growth on methanol allows isolation of Mut <sup>S</sup> strains (see below) ( <i>aox1</i> ) (Cregg <i>et al.</i> , 1989; Koutz <i>et al.</i> , 1989).

## Overview of Pichia pastoris Expression System, Continued

Expression of the <i>AOX1</i> gene is controlled at the level of transcription. In methanol- grown cells approximately 5% of the total polyA <sup>+</sup> RNA is from the <i>AOX1</i> gene. The regulation of the <i>AOX1</i> gene is a two step process of repression/derepression and induction mechanism. Briefly, growth on glucose represses transcription, even in the presence of the inducer methanol. For this reason, growth on glycerol is recommended for optimal induction with methanol. <b>Note:</b> Growth on glycerol alone (derepression) is not sufficient to generate even minute levels of expression from the <i>AOX1</i> gene. The inducer, methanol, is necessary for detectable levels of <i>AOX1</i> expression (Ellis <i>et al.</i> , 1985; Koutz <i>et al.</i> , 1989; Tschopp <i>et al.</i> , 1987a).
Loss of the <i>AOX1</i> gene results in a strain that is phenotypically Mut <sup>S</sup> (Methanol <u>utilization</u> slow, also referred to as Mut <sup>-</sup> in the past). The Mut <sup>S</sup> designation is chosen to accurately describe the phenotype of these mutants. This results in a reduction in the cells ability to metabolize methanol and they exhibit poor growth on methanol medium. Mut <sup>+</sup> (Methanol <u>utilization</u> plus) refers to the wild type ability of strains to metabolize methanol as the sole carbon source. These two phenotypes are used when evaluating <i>Pichia</i> transformants for integration of your gene ( <b>Experimental Outline</b> , page 7).
Heterologous expression in <i>Pichia pastoris</i> can be either intracellular or secreted. Secretion requires the presence of a signal sequence on the expressed protein to target it to the secretory pathway. The native secretion signal present on some heterologous proteins and several different secretion signal sequences have been used with varied success. <i>Saccharomyces cerevisiae</i> $\alpha$ factor secretion signal sequence has been used most successfully (Cregg et al., 1993; Scorer et al., 1993) The major advantages of expressing recombinant proteins as secreted proteins are: • <i>Pichia pastoris</i> secretes very low levels of native proteins
• Very low amount of protein is present in the minimal <i>Pichia</i> growth medium The secreted recombinant protein usually comprises the vast majority of the total protein in the medium and serves as the first step in purification of the protein (Barr et al., 1992). <b>Note</b> : If there are recognized glycosylation sites (Asn-X-Ser/Thr) in your protein's primary sequence, glycosylation may occur at these sites.
Unlike <i>Saccharomyces cerevisiae</i> , <i>Pichia</i> does not hyperglycosylate the secreted proteins. Both <i>Saccharomyces cerevisiae</i> and <i>Pichia pastoris</i> have a majority of N-linked glycosylation of the high-mannose type; however, the length of the oligosaccharide chains added posttranslationally to proteins in <i>Pichia</i> (average 8-14 mannose residues per side chain) is much shorter than those in <i>Saccharomyces cerevisiae</i> (50-150 mannose residues) (Grinna & Tschopp, 1989; Tschopp et al., 1987b) Very little O-linked glycosylation has been observed in <i>Pichia</i> . In addition, <i>Saccharomyces cerevisiae</i> core oligosaccharides have terminal $\alpha$ 1,3 glycan linkages whereas <i>Pichia pastoris</i> does not. It is believed that the $\alpha$ 1,3 glycan linkages in glycosylated proteins produced from <i>Saccharomyces cerevisiae</i> are primarily responsible for the hyper-antigenic nature of these proteins making them particularly unsuitable for therapeutic use. Although not yet proven, this is predicted to be less of a problem for glycoproteins generated in <i>Pichia pastoris</i> , because it may resemble the glycoprotein structure of higher eukaryotes (Cregg <i>et al.</i> , 1993).

## Overview of Pichia pastoris Expression System, Continued

Transformation and Integration	Two different phenotypic classes of recombinant strains can be generated: Mut <sup>+</sup> and Mut <sup>S</sup> . Mut <sup>S</sup> refers to the " <u>M</u> ethanol <u>utilization slow</u> " phenotype caused by the loss of alcohol oxidase activity encoded by the <i>AOX1</i> gene. A strain with a Mut <sup>S</sup> phenotype has a mutant <i>aox1</i> locus, but is wild type for <i>AOX2</i> . This results in a slow growth phenotype on methanol medium. <i>Pichia</i> strains X-33 and GS115 are Mut <sup>+</sup> while KM71H is Mut <sup>S</sup> . Transformation of X-33 or GS115 with plasmid DNA linearized in the 5' <i>AOX1</i> region will yield Mut <sup>+</sup> transformants, while KM71H will yield only Mut <sup>S</sup> transformants. Both Mut <sup>+</sup> and Mut <sup>S</sup> recombinants are useful to have as one phenotype may favor better expression of your protein than the other. You should test between 6-10 recombinants per phenotype because the site of recombination may affect expression. There is no way to predict beforehand which construct or isolate will better express your protein. For more information on recombination in <i>Pichia</i> , please see page 51. Once you have successfully cloned your gene behind the <i>AOX1</i> promoter, you will then
	linearize your plasmid to permit recombination when the plasmid is transformed into <i>Pichia</i> .
Expression and Scale-up	After isolating your <i>Pichia</i> recombinants, you will then test expression of both Mut <sup>+</sup> and Mut <sup>S</sup> recombinants. This will involve growing a small culture of each recombinant, inducing with methanol, and taking time points. If looking for intracellular expression, analyze the cell pellet from each time point by SDS polyacrylamide gel electrophoresis (SDS-PAGE). If looking for secreted expression, analyze both the cell pellet and supernatant from each time point. We recommend that you analyze your SDS-PAGE gels by both Coomassie staining and western blot (for proteins expressed at low levels). We also suggest checking for protein by functional assay if one is available.
	Choose the <i>Pichia</i> recombinant strain that best expresses your protein and optimize induction based on the suggestions on pages 32. Once expression is optimized, scale-up your expression protocol to produce more protein for purification.
Purification	The donor vector contains a C-terminal tag consisting of a V5 epitope and a polyhistidine tag. If the donor vector is recombined with the acceptor vector (pPICZ-E or pPICZ $\alpha$ -E) such that the C-terminal tag is maintained, then both pPICZ-E or pPICZ $\alpha$ -E will contain the V5 epitope and the polyhistidine tag. The polyhistidine tag facilitates purification by binding to divalent cations like Ni <sup>2+</sup> usually present in metal-binding resins such as ProBond <sup>TM</sup> . We recommend that you use the ProBond <sup>TM</sup> Purification System (Catalog no. K850-01) to purify fusion proteins expressed using pPICZ-E or pPICZ $\alpha$ -E. Preliminary preparation steps are described on pages 36-37.
	If you are using a metal-chelating resin other than ProBond <sup>™</sup> , follow the manufacturer's recommendations for fusion proteins expressed in yeast.

## Overview of Pichia pastoris Expression System, Continued

### Experimental Outline

The table below describes the general steps needed to recombine, transform, and express your protein of interest.

Step	Action	Page
1	Perform the recombination reaction using your donor vector and pPICZ-E or pPICZ $\alpha$ -E.	8
2	Transform the recombination reaction into competent TOP10 E. coli.	9
3	Select transformants on LB plates containing 50 µg/ml kanamycin.	9
4	Analyze transformants by restriction digestion.	10
5	Select the correct clone and linearize the construct with appropriate restriction enzyme.	17
6	Transform your construct into appropriate competent <i>Pichia</i> host strain (X-33 or GS115 for Mut <sup>+</sup> , and KM71H for Mut <sup>s</sup> ) using your transformation method of choice. Select transformants on medium containing Zeocin <sup>TM</sup> .	18-21
7	Select 6-10 clones for small scale expression and analyze the expression of your recombinant protein by western blot analysis or functional assay.	25
8	Choose the highest expressers for large-scale expression in shake flask or fermenter.	34
9	Purify your protein using metal-chelating resin (Probond <sup><math>TM</math></sup> ) or any other method of choice.	36

## Methods

## Recombining Your Gene into pPICZ-E or pPICZ $\alpha$ -E

Introduction	You will need a plasmid preparation of your donor vector containing the gene of interest in addition to the pPICZ-E or pPICZ $\alpha$ -E vector. Review the information below and on th next page before performing the recombination reaction.					
Preparation and Maintenance of pPICZ-E or pPICZα-E	To prepare pPICZ-E or pPICZ $\alpha$ -E for use, add 20 µl sterile water to prepare a 1 µg/µl stock solution. You can further dilute a small aliquot of plasmid or use the stock solution as is. Store the stock solution at -20°C when you are finished.					
	If you wish to propagate the pPICZ-E or pPICZ $\alpha$ -E plasmid or prepare plasmid DNA, you may transform the plasmid into One Shot <sup>®</sup> TOP10 Chemically Competent <i>E. coli</i> as described on page 9 or TOP10F' provided in the kit. Use10-100 ng of plasmid DNA for transformation and select transformants on low salt LB plates containing 25 µg/ml Zeocin <sup>™</sup> . Be sure to prepare a glycerol stock of your plasmid-containing TOP10 strain for long-term storage (see page 12).					
Before Starting	You will need the following reagents and equipment.					
	٠	• 100 ng of your donor vector construct				
	٠	• 100 ng of pPICZ-E or pPICZ $\alpha$ -E (included in the kit)				
	•	Microcentrifuge tubes				
	•	• Heat blocks set at 37°C and 65°C				
	• Ice bucket with ice					
	•	• Cre recombinase (included in the kit)				
	•	10X Recombinase Buffer (				
Recombination	1.	Set up each 20 µl recombi	nation reaction on ice as follows	S:		
Reaction		Donor vector (100 ng)	x µl			
		pPICZ-E or pPICZ $\alpha$ -E (10	00 ng) y μl			
		10X Recombinase Buffer	2 µl			
		Deionized water	add to a total volume of 19 µl			
		Cre Recombinase	1 µl			
		Final Volume	20 µl			
	2.	Incubate at 37°C for 20 m				
	3.	Incubate at 65°C for 5 minutes to inactivate the recombinase.				
	4.		proceed to <b>Transformation</b> , nex combination reaction at +4°C or en tested.			

## **Transforming the Recombination Reaction**

Introduction	Once you have performed the recombination reaction, you are ready to transform your <i>E. coli</i> host. We recommend using TOP10 <i>E. coli</i> for transformation, but other strains are suitable. <i>E. coli</i> strains should be endonuclease A deficient ( <i>end</i> A) and recombination deficient ( <i>rec</i> A) to ensure quality plasmid preparations and reduce the chances of non-specific recombination, respectively.				
Materials Supplied by the User	In addition to general microbiological supplies (i.e. plates, spreaders), you will need the following reagents and equipment.				
	• 42°C water bath				
	• LB plates containing 50 µg/ml kanamycin (see <b>Important</b> , below)				
	• LB medium and SOB medium (see page 45 for recipe)				
	• 37°C shaking and non-shaking incubator				
<b>Q</b> Important	It is important to select the fusion plasmid using kanamycin. The donor vector contains the R6K $\gamma$ origin that can only be maintained in <i>E. coli</i> strains containing the <i>pir</i> gene. After recombination between the donor vector and acceptor vector, the kanamycin resistance gene (from the donor vector) in the fusion plasmid is linked to the pUC origin (from pPICZ-E or pPICZ $\alpha$ -E). The fusion plasmid can be maintained in <i>E. coli</i> strains that do not contain the <i>pir</i> gene (i.e. TOP10). By selecting for kanamycin resistance, you ensure that only colonies containing the fusion plasmid are selected.				
Preparing for Transformation	This transformation protocol is for use with the One Shot <sup>®</sup> TOP10 Chemically Competent <i>E. coli</i> available with the kit. Follow the manufacturer's protocol if you are using other competent cells.				
	For each transformation, you will need one vial of One Shot <sup>®</sup> TOP10 Chemically Competent <i>E. coli</i> and two selective plates. Perform the following steps before beginning.				
	1. Equilibrate a water bath to 42°C.				
	2. Thaw the vial of SOC medium from the kit and bring to room temperature.				
	3. Warm LB plates containing 50 µg/ml kanamycin at 37°C for 30 minutes.				
	4. Thaw <u>on ice</u> 1 vial of One Shot <sup>®</sup> cells for each transformation.				
One Shot <sup>®</sup> Transformation	<ol> <li>Add 5 μl of the recombination reaction (step 4, page 8) to a vial of One Shot<sup>®</sup> TOP10 E. coli and mix by stirring with a pipette tip. Do not mix by pipetting up and down.</li> </ol>				
Reaction	2. Heat-shock the cells for 30 seconds at 42°C without shaking.				
	3. Immediately transfer the tubes to ice.				
	4. Add 500 μl of room temperature SOC medium.				
	5. Cap the tube tightly and shake the tube horizontally at 37°C for 45 minutes.				
	<ol> <li>Spread 50 μl from each transformation onto a prewarmed LB plate containing 50 μg/ml kanamycin. Pellet the remaining cells, resuspend the cell pellet in 50 μl SOC and plate. Incubate plates overnight at 37°C.</li> </ol>				
	7. An efficient recombination reaction will produce hundreds of colonies. Pick five colonies for analysis.				

## Transforming the Recombination Reaction, Continued

Analyzing Pos Clones	itive		Culture the 5 colo ontaining 50 μg/i			vernight in 2	2-5 ml LB o	r SOB medi	um
		E K	<ol> <li>Isolate plasmid DNA using your method of choice. If you need ultra-pure plasmid DNA for automated or manual sequencing, we recommend the S.N.A.P.<sup>™</sup> MiniPrep Kit (10-15 µg DNA, Catalog no. K1900-01) or the S.N.A.P.<sup>™</sup> MidiPrep Kit (10- 200 µg DNA, Catalog no. K1910-01).</li> </ol>						
		0	nalyze the plasm nce in the donor re distinguishable	vector and one	ce in the acce	ptor vector	to yield two	o fragments	
		5' ne	AOX1 primer ar ext page for the s	onal) To sequence the fusion plasmid to confirm the fusion junctions, use the $DXI$ primer and the Uni1 Forward primer. Refer to the diagrams below and on the page for the sequence around the pPICZ-E or pPICZ $\alpha$ -E <i>loxP</i> site. Refer to the rector manual for the sequence around the donor vector <i>loxP</i> site.				on the	
		If you need help with setting up restriction enzyme digests or DNA sequencing, refer to general molecular biology texts (Ausubel et al., 1994; Sambrook et al., 1989)				to			
Sequencing Yo Construct in pPICZ-E		to indic	quence surroundicate the cleavage bading from our V	site. Note that	t the complet	e sequence	of pPICZ-E	E is available	for
			5′	end of AOX1 mF	NA		5	5' AOX1 priming	g site
801	CCCTG	ICTTA	AACCTTTTTT	TTTATCATC	Α ΤΤΑΤΤΑ(	GCTT ACT	ТТСАТАА	TTGCGACT	'GG
861	TTCCA	ATTGA	CAAGCTTTTG	ATTTTAACG	a ctttta	ACGA CAA	CTTGAGA	AGATCAAA	AA
			Sfu I	EcoR I			loxP site		
921	ACAAC	FAATT	ATTCGAAACG	AGGAATTC	ATA ACT	ICG TAT	AGC ATA	CAT TAT	ACG
						U	ni1 Forward I	priming site	
- 976	AAG TI	TA T	donor Gen vector inte		minal tag otional)	do	nor vector	→	
	loxP s	Age ite AC		AGATTC TAA	TCAAGAG (	GATGTCAG	AA TGCC <i>i</i>	ATTTGC	

### **Transforming the Recombination Reaction, Continued**

**Sequencing Your** The sequence surrounding your insert is shown below. Unique restriction sites are labeled **Construct** in to indicate the cleavage site. Note that the complete sequence of pPICZ $\alpha$ -E is available for downloading from our Web site: (www.invitrogen.com) or from Technical Service (see ρΡΙCΖα-Ε page 61). 5' end of AOX1 mRNA 5' AOX1 priming site 811 AACCTTTTTT TTTATCATCA TTATTAGCTT ACTTTCATAA TTGCGACTGG TTCCAATTGA 871 CAAGCTTTTG ATTTTAACGA CTTTTAACGA CAACTTGAGA AGATCAAAAA ACAACTAATT 931 ATTCGAAACG ATG AGA TTT CCT TCA ATT TTT ACT GCT GTT TTA TTC GCA GCA Met Arg Phe Pro Ser Ile Phe Thr Ala Val Leu Phe Ala Ala 983 TCC TCC GCA TTA GCT GCT CCA GTC AAC ACT ACA ACA GAA GAT GAA ACG Ser Ser Ala Leu Ala Ala Pro Val Asn Thr Thr Glu Asp Glu Thr 1031 GCA CAA ATT CCG GCT GAA GCT GTC ATC GGT TAC TCA GAT TTA GAA GGG Ala Gln Ile Pro Ala Glu Ala Val Ile Gly Tyr Ser Asp Leu Glu Gly  $\alpha$ -factor signal sequence 1079 GAT TTC GAT GTT GCT GTT TTG CCA TTT TCC AAC AGC ACA AAT AAC GGG Asp Phe Asp Val Ala Val Leu Pro Phe Ser Asn Ser Thr Asn Asn Gly α-factor priming site 1127 TTA TTG TTT ATA AAT ACT ACT ATT GCC AGC ATT GCT GCT AAA GAA GAA Leu Leu Phe Ile Asn Thr Thr Ile Ala Ser Ile Ala Ala Lys Glu Glu EcoR I loxP site Kex2 signal cleavage 1175 GGG GTA TCT CTC GAG AAA AGAVGAG GCT GAA GCT GAATTC ATA ACT TCG TAT Gly Val Ser Leu Glu Lys Arg Glu Ala Glu Ala Uni1 Forward priming site Ste13 signal cleavage donor C-terminal tag donor Gene of 1226 AGC ATA CAT TAT ACG AAG TTA T (optional) vector interest vector Age I ACCGGTCTTG CTAGATTCTA ATCAAGAGGA TGTCAGAATG CCATTTGCCT loxP site

## Transforming the Recombination Reaction, Continued

Fusion Plasmid Analysis	It should be clear from restriction analysis that you have a dimer plasmid consisting of the donor vector and pPICZ-E or pPICZ $\alpha$ -E. Occasionally, trimers will result. Trimers usually consist of two donor vector molecules and one acceptor molecule, but they usually express as well as the dimer product.				
	In theory, trimers may result from two sequential fusion events or a single fusion event between a pre-existing monomeric substrate and a dimeric substrate. The production of trimers can be eliminated if gel-purified monomeric supercoiled DNA is used in the recombination reaction.				
Preparing Plasmid DNA and Glycerol	Once you have identified the correct clone, isolate plasmid DNA using your method of choice for transformation of your construct into <i>Pichia</i> (see page 15-21).				
Stock for Long-	We recommend to make a glycerol stock for long term storage.				
Term Storage	1. Streak out the original colony on LB plates containing 50 $\mu$ g/ml kanamycin to isolate single colonies.				
	2. Select a single colony and inoculate into 1-2 ml of LB containing 50 $\mu$ g/ml kanamycin.				
	3. Grow overnight until culture is saturated.				
	4. Mix 0.85 ml of culture with 0.15 ml of sterile glycerol and transfer to a cryovial.				
	5. Store at -80°C. (You may also want to store a stock of plasmid DNA at -20°C.)				

Introduction	After isolating the plasmid DNA of your fusion construct, you are ready to transfom this construct into <i>Pichia</i> for expression of the recombinant protein. This section provides information on <i>Pichia</i> strains, initiation of <i>Pichia</i> culture, growth characteristics and storage of <i>Pichia pastoris</i> . You should be familiar with basic sterile microbiological, molecular biology and protein chemistry techniques before attempting to grow and manipulate <i>Pichia</i> . Some general references to consult are <i>Guide to Yeast Genetics and Molecular Biology</i> , (Guthrie & Fink, 1991), <i>Current Protocols in Molecular Biology</i> , (Ausubel <i>et al.</i> , 1994), <i>Molecular Cloning: A Laboratory Manual</i> , (Sambrook <i>et al.</i> , 1989), <i>Protein Methods</i> , (Bollag & Edelstein, 1991) and <i>Guide to Protein Purification</i> , (Deutscher, 1990).
Genotypes of <i>Pichia</i> Strains	X-33 is a wild-type <i>Pichia</i> strain that is useful for selection on Zeocin <sup>TM</sup> and large-scale growth. It will grow in YPD and in minimal media.
	The <i>Pichia</i> host strain GS115 has a mutation in the histidinol dehydrogenase gene ( <i>his4</i> ) that prevents it from synthesizing histidine. GS115 will grow on complex medium such as YPD (also known as YEPD) and on minimal media supplemented with histidine.
	The parent strain of KM71H has a mutation in the argininosuccinate lyase gene ( <i>arg4</i> ) that prevents the strain from growing in the absence of arginine. The wild-type <i>ARG4</i> gene was used to disrupt <i>AOX1</i> , creating KM71H, a Mut <sup>S</sup> , Arg <sup>+</sup> strain.
Construction of KM71H	The ARG4 gene (~2 kb) was inserted into the cloned, wild-type AOX1 gene between the BamH I site (codons 15/16 of AOX1) and the Sal I site (codons 227/228 of AOX1). ARG4 replaces codons 16 through 227 of AOX1. This construct was transformed into the parent strain of KM71 (arg4 his4) and Arg <sup>+</sup> transformants were isolated and analyzed for the Mut <sup>S</sup> phenotype. Genetic analysis of Arg <sup>+</sup> transformants showed that the wild-type AOX1 gene was replaced by the aox1::ARG4 construct. To create KM71H, KM71 was transformed with a gene fragment encoding the HIS4 gene and a His <sup>+</sup> convertant was isolated.
<b>O</b> Important	The advantage of using KM71H is that there is no need to screen for the Mut phenotype on methanol minimal medium. All transformants will be Mut <sup>S</sup> . Secondly, since the <i>AOX1</i> locus was not completely deleted, it is theoretically possible to replace $aox1::ARG4$ with your construct by gene replacement. The phenotype of this strain would be Mut <sup>S</sup> Arg <sup>-</sup> . This means the recombinant strain would require arginine in the medium to grow. Unfortunately, simple inclusion of arginine does not totally alleviate the effects of the <i>arg4</i> mutation, and <i>arg4</i> strains do not grow well on minimal medium supplemented with arginine. Therefore, we do not recommend that you generate transformants in KM71H by replacing the <i>aox1::ARG4</i> construct.

# Pichia Strains, Continued

Growth of <i>Pichia</i> Strains	The growth temperature of <i>Pichia pastoris</i> is 28-30°C for liquid cultures, plates, and stabs. Growth above 32°C during induction can be detrimental to protein expression and can even lead to cell death. Other important facts:				
	• Doubling time of log phase Mut <sup>+</sup> or Mut <sup>S</sup> <i>Pichia</i> in YPD is ~2 hours				
	• Mut <sup>+</sup> and Mut <sup>S</sup> strains do not differ in growth rates unless grown on methanol				
	• Doubling time of log phase Mut <sup>+</sup> <i>Pichia</i> in methanol medium (MM) is 4-6 hours				
	• Doubling time of log phase Mut <sup>S</sup> <i>Pichia</i> in MM is ~18 hours				
	• One $OD_{600} = ~5 \times 10^7 \text{ cells/ml}$				
	Note that growth characteristics may vary depending on the recombinant protein expressed				
Growth on Methanol	When plates or medium containing methanol are used as growth medium, it is advisable to add methanol every day to compensate for loss because of evaporation or consumption.				
	• For plates add 100 µl of 100% methanol to the lid of the inverted plate.				
	• For liquid medium add 100% methanol to a final concentration of 0.5%.				
	Some researchers have had success adding methanol to 1% every day for Mut <sup>S</sup> strains and up to 3% for Mut <sup>+</sup> without any negative effect to their liquid culture				
Initiating <i>Pichia</i> culture	To initiate a culture of <i>Pichia</i> from the stab provided with the kit, streak a small amount from the stab on YPD plate (see page 41) and incubate at 30°C. Once growth is established, you may pick a single colony and grow it in the appropriate medium for downstream applications. Remember to prepare frozen stocks for long-term storage of all <i>Pichia</i> strains included in this kit (see below).				
Storage of <i>Pichia</i> Strains	To store cells for weeks to months, use YPD medium and YPD agar slants (see page 41).				
	<ul> <li>Streak each strain for single colonies on YPD.</li> </ul>				
	• Transfer one colony to a YPD stab and grow for 2 days at 30°C.				
	• The cells can be stored on YPD for several weeks at $+4^{\circ}$ C.				
	<ul> <li>To store cells for months to years, store frozen at -80°C.</li> <li>Culture a single colony of each strain overnight in YPD.</li> </ul>				
	<ul> <li>Harvest the cells and suspend in YPD containing 15% glycerol at a final OD<sub>600</sub> of 50-100 (approximately 2.5-5.0 x 10<sup>9</sup> cells/ml).</li> </ul>				
	• Cells are frozen in liquid nitrogen or a dry ice/ethanol bath and then stored at -80°C.				
Note	After extended storage at $+4^{\circ}$ C or $-80^{\circ}$ C, it is recommended that Zeo <sup>R</sup> transformants be checked for correct phenotype and protein expression.				

## Pichia Transformation

Introduction	Once you have created the pPICZ-E or pPICZ $\alpha$ -E fusion plasmid and have verified its integrity, you are ready to transform the fusion plasmid into a <i>Pichia</i> host strain of your choice.
Zeocin <sup>™</sup> Selection	We generally use 100 $\mu$ g/ml Zeocin <sup>TM</sup> to select for transformants when using the X-33 <i>Pichia</i> strain. If you are transforming the fusion plasmid 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 <sup>TM</sup> 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 <sup>TM</sup> 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 <sup>TM</sup> resistance gene. For this reason, plating spheroplasts directly onto selective medium containing Zeocin <sup>TM</sup> does not yield any transformants.
	We recommend electroporation for transformation of <i>Pichia</i> with pPICZ-E or pPICZ $\alpha$ -E. Electroporation yields 10 <sup>3</sup> to 10 <sup>4</sup> 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 <i>Pichia</i> EasyComp <sup>TM</sup> Transformation Kit included in the kit (see page 19). The <i>Pichia</i> EasyComp <sup>TM</sup> Transformation Kit provides reagents to prepare 6 preparations of competent cells. Each preparation will yield enough competent cells for 20 transformations
<b>Q</b> Important	Since pPICZ-E or pPICZ $\alpha$ -E do not contain the <i>HIS4</i> gene, integration can only occur at the <i>AOX1</i> locus. Vector linearized within the 5' <i>AOX1</i> region will integrate by gene insertion into the host 5' <i>AOX1</i> region. Therefore, the <i>Pichia</i> host that you use will determine whether the recombinant strain will be Mut <sup>+</sup> or Mut <sup>S</sup> . Use a <i>Pichia</i> host that contains the native <i>AOX1</i> gene (e.g. X-33, GS115, SMD1168H) to generate a Mut <sup>+</sup> recombinant strain and a <i>Pichia</i> host that has a disrupted <i>AOX1</i> gene (i.e. KM71H) to generate a Mut <sup>S</sup> recombinant strain. Information on <i>Pichia</i> recombination is provided on page 51.
<i>His4</i> Host Strains	Host strains containing the <i>his4</i> allele (e.g. GS115) and transformed with the pPICZ-E or pPICZ $\alpha$ -E vectors require histidine when grown in minimal media. Add histidine to a final concentration of 0.004% to ensure growth of your transformants.

## Pichia Transformation, Continued



The pPICZ-E or pPICZ $\alpha$ -E vectors do not contain a yeast origin of replication. Transformants can only be isolated if recombination occurs between the plasmid and the *Pichia* genome.

**Pichia Controls** pPICZ-E/Uni-lacZ and pPICZ\alpha-E/Uni-HSA are control plasmids included in the kit to check for transformation and expression in Pichia pastoris. The lacZ gene was cloned into a donor vector, pUni/V5-His-Gene-TOPO®. The resulting vector was recombined with pPICZ-E to create pPICZ-E/Uni-lacZ, which can be used as an intracellular expression control after transformation into a *Pichia* host. The human serum albumin (HSA) gene was cloned into a donor vector, pUni/V5-His-Gene-TOPO<sup>®</sup>. The resulting vector was recombined with pPICZ $\alpha$ -E to create pPICZ $\alpha$ -E/Uni-HSA/V5-His, which can be used as a secreted expression control. For a map and more details on pPICZ-E/Uni-*lacZ* and pPICZ $\alpha$ -E/Uni-HSA controls, see pages 49-50. Use the control plasmids pPICZ-E/Uni-lacZ or pPICZ $\alpha$ -E/Uni-HSA to perform a control transformation using any method of choice in a *Pichia* host strain. We have used X-33 as the host strain. Linearize the plasmid using *Pme* I to promote integration (see next page). If you wish to propagate the supercoiled plasmids, you may transform the plasmid into TOP10 E. coli (see page 8). GS115/pPICZ/lacZ (Mut<sup>+</sup>) and GS115/Albumin (Mut<sup>s</sup>) are control strains included in the kit to check the Mut phenotype. Cloning the lacZ gene into the pPICZ vector and transforming the resulting construct into GS115 creating the GS115/pPICZ/lacZ (Mut<sup>+</sup>) strain. This strain expresses  $\beta$ -galactosidase and can also be used as a control for intracellular expression. GS115/Albumin is Mut<sup>s</sup> and is also a control for secreted expression. The gene for serum albumin was cloned with its native secretion signal and then transformed into GS115 to create GS115/Albumin. If you decide to use GS115/pPICZ/lacZ or GS115/Albumin as a control, then you need to initiate a culture from the stab (see page 14). We recommend you use pPICZ-E/Uni-*lacZ* and pPICZ $\alpha$ -E/Uni-HSA as transformation and expression controls, and GS115/pPICZ/lacZ and GS115/Albumin as controls for determining the Mut phenotype (see page 22) and confirming expression conditions. **Before Starting** The following reagents are needed for linearizing your construct, transforming *Pichia* and selecting transformants on  $\text{Zeocin}^{\text{TM}}$ . 10-50  $\mu$ g pPICZ-E or pPICZ $\alpha$ -E fusion plasmid and appropriate expression control plasmids Appropriate restriction enzymes and buffer Phenol/chloroform 3M sodium acetate Agarose gel apparatus and Electroporation device with 0.2 cm cuvettes 80% and 100% ethanol YPD Medium and YPDS plates containing the appropriate concentration of Zeocin<sup>™</sup> 50 ml conical polypropylene tubes 1 liter cold (+4°C) sterile water (place on ice the day of the experiment) 25 ml cold (+4°C) sterile 1 M sorbitol (place on ice the day of the experiment) Continued on next page

### Pichia Transformation, Continued

1.

#### Linearizing Your pPICZ-E or pPICZα-E Construct

To promote integration, we recommend that you linearize your pPICZ-E or pPICZ $\alpha$ -E fusion plasmid within the 5' *AOX1* region. The table below lists unique sites that may be used to linearize pPICZ-E or pPICZ $\alpha$ -E prior to transformation. **Other restriction sites are possible**. Be sure that your insert does not contain the restriction site you wish to use to linearize your vector.

Enzyme	<b>Restriction Site (bp)</b>	Supplier
Sac I	209	New England Biolabs
Pme I	414	New England Biolabs
BstX I	707	New England Biolabs

#### **Restriction Digest**

Digest ~10-50  $\mu$ g of plasmid DNA with one of the enzymes listed above. We recommend to use more DNA as you will lose some DNA during purification after the restriction digestion and will be helpful if you decide to do multiple transformations with varying concentrations of Zeocin<sup>TM</sup>.

- 2. Check a small aliquot of your digest by agarose gel electrophoresis for complete linearization.
- 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. You can also use any resin based purification methods to purify the linearized vector.
- 4. Centrifuge the solution to pellet the DNA, wash the pellet with 80% ethanol, air-dry, and resuspend in 10  $\mu$ l sterile, deionized water. Use immediately or store at -20°C.

Preparing *Pichia* for Electroporation Follow the procedure below to prepare your Pichia pastoris strain for electroporation.

- 1. Grow 5 ml of your *Pichia pastoris* 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 x g for 5 minutes at  $+4^{\circ}$ 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  $(0^{\circ}C)$ , 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.
- 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.

## Pichia Transformation, Continued

Transformation by	1. Mix 80 µl of the cells from Step 6 (above) with 5-10 µg of linearized DNA (in 5-		
Electroporation	$10 \mu l$ sterile water) and transfer them to an ice-cold 0.2 cm electroporation cuvette.		
	. 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.		
	<ol> <li>Immediately add 500 μl of ice-cold 1 M sorbitol to the cuvette. Transfer the cuvette contents to a sterile 15 ml tube. Rinse the cuvette with additional 500 μl of ice-cold 1 M sorbitol and transfer the contents to the 15 ml tube as above.</li> </ol>		
	5. Let the tube incubate at 30°C without shaking for 1 to 2 hours.		
	<ol> <li>Spread 50-200 µl each on separate, labeled YPDS plates containing the appropriate concentration of Zeocin<sup>™</sup>. Note: Inclusion of sorbitol in YPD plates stabilizes electroporated cells as they appear to be somewhat osmotically sensitive</li> </ol>		
	7. Incubate plates for 2 to 7 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 <sup>™</sup> .		
Isolation of Multi- copy Recombi-	A quick, direct way to select multi-copy recombinants is to plate the transformation mix on increasing concentrations of Zeocin <sup><math>TM</math></sup> .		
nants <i>in vivo</i>	1. Prepare YPDS plates containing 100, 500, 1000, and 2000 $\mu$ g/ml Zeocin <sup>TM</sup>		
	2. Plate 100 to 200 $\mu$ l of the transformation mix on each plate and incubate at 30°C for 2 to 7 days.		
	3. Test transformants for the Mut phenotype (page 22) and expression of your protein (page 25)		
Note	Generally, several hundred Zeocin <sup>TM</sup> -resistant colonies are generated using the protocol on the previous page. If more colonies are needed, the protocol may be modified as described below. Note that you will need twenty 150 mm plates with YPDS agar containing the appropriate concentration of Zeocin <sup>TM</sup> .		
	1. Set up two transformations per construct and follow Steps 1 through 5 of the <b>Transformation by Electroporation</b> protocol, see above.		
	2. After 1 hour in 1 M sorbitol at 30°C (Step 5, see above), add 1 ml YPD medium to each tube.		
	3. Shake (~200 rpm) the cultures at $30^{\circ}$ C.		
	<ol> <li>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<sup>™</sup>.</li> </ol>		
	<ol> <li>(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<sup>™</sup>.</li> </ol>		
	5. Incubate plates for 2 to 7 days at 30°C until colonies form.		
Analyzing <i>Pichia</i> Transformants	Select 6-10 of your Zeocin <sup>™</sup> -resistant <i>Pichia</i> transformants and confirm the Mut phenotype as described on page 22. You can also analyze for the presence of insert using PCR (page 55), or for copy number using Southern analysis (page 60).		

# EasyComp<sup>™</sup> Transformation

Introduction	The Pichia EasyComp <sup>™</sup> Kit produces chemically competent Pichia cells and is inclusion to provide a rapid and convenient alternative to electroporation. However, because low transformation efficiency (3 µg plasmid DNA yields about 50 colonies), it is very difficult to isolate multi-copy integrants. If multi-copy integrants are desired, use electroporation (page 16) for best results. Do not use cells prepared using the EasyComp <sup>™</sup> protocol for electroporation as cells are prepared differently for emethod.		
Required	• 30°C rotary shaking incubator		
Reagents and	• YPD (Yeast Extract Peptone Dextrose) medium (see <b>Recipes</b> , page 41)		
Equipment	• 50 ml, sterile conical tubes		
	• Centrifuge suitable for 50 ml conical tubes (floor or table-top)		
	• 1.5 ml sterile screw-cap microcentrifuge tubes		
	• Styrofoam box or paper towels		
Before Beginning	Streak a YPD plate with your <i>Pichia pastoris</i> strain such that isolated, single colonies will grow. Incubate the plate at 28-30°C for 2 days.		
	Equilibrate Solution I to room temperature.		
Preparing Competent Cells	<ol> <li>Inoculate 10 ml of YPD with a single colony of your <i>Pichia</i> strain. Grow overnight at 28-30°C in a shaking incubator (250-300 rpm).</li> </ol>		
	2. Dilute cells from the overnight culture to an $OD_{600}$ of 0.1-0.2 in 10 ml of YPD. Grow the cells at 28-30°C in a shaking incubator until the $OD_{600}$ reached 0.6-1.0. This will take approximately 4 to 6 hours.		
	<ol> <li>Pellet the cells by centrifugation at 500 x g for 5 minutes at room temperature. Discard the supernatant.</li> </ol>		
	4. Resuspend the cell pellet in 10 ml of Solution I. No incubation time is required.		
	5. Pellet the cells by centrifugation at 500 x g for 5 minutes at room temperature. Discard the supernatant.		
	6. Resuspend the cell pellet in 1 ml of Solution I. The cells are now competent.		
	<ol> <li>Aliquot 50 to 200 μl of competent cells into labeled 1.5 ml sterile screw-cap microcentrifuge tubes.</li> </ol>		
	<b>Note</b> : 50 $\mu$ l of cells are used for each transformation. Cells can be thawed and refrozen several times without significant loss in transformation efficiency.		
	8. At this point, the cells may be kept at room temperature, used directly for transformation or frozen for future use. To slowly freeze the cells, place tubes in a Styrofoam box or wrap in several layers of paper towels and place in a -80°C freezer. <b>Do not snap-freeze the cells in liquid nitrogen.</b>		
	9. Proceed to the transformation procedure.		
Note	Higher transformation efficiencies are obtained with frozen versus freshly prepared cells. You may choose to use some of the cells immediately following preparation and freeze the remaining cells in small aliquots.		

## EasyComp<sup>™</sup> Transformation, Continued

Transformation	The following protocol can be used to transform either freshly prepared or frozen compe- tent <i>Pichia</i> cells. Transformation efficiency may vary with each strain and vector used.	
Required Reagents and Equipment	•	30°C incubator Water baths or heat blocks at 30°C and 42°C Microcentrifuge at room temperature YPDS with 100 µg/ml Zeocin <sup>™</sup> plates (see <b>Recipes</b> , page 41)
Before Beginning	diss Equ Equ need You	bu see a precipitate in Solution II, warm the solution at 37°C until the precipitate olves. To prevent formation of a precipitate, store Solution II at room temperature. ilibrate Solution III to room temperature. ilibrate the appropriate number and type of plates to room temperature. You will d one plate for each transformation.
	cont	rol.
Transformation Protocol	1.	For each transformation, thaw one tube of competent cells at room temperature and aliquot 50 $\mu$ l into a sterile microcentrifuge tube. If transforming fresh cells, use 50 $\mu$ l of cells from <b>Preparation of Competent Cells</b> , Step 7, previous page.
	2.	Add 3 µg of linearized Pichia expression vector DNA to the competent cells.
		<b>Note</b> : Using greater than 3 $\mu$ g of DNA may increase transformation efficiencies in some cases. The volume of DNA should not exceed 5 $\mu$ l. Linearized DNA can be used directly from a restriction digest reaction without affecting transformation efficiency. Phenol chloroform extraction and ethanol precipitation are not necessary.
	3.	Add 1 ml of Solution II to DNA/cell mixture. Mix by vortexing or flicking the tube.
	4.	Incubate the transformation reactions for 1 hour at 30°C in a water bath or incubator. Mix the transformation reaction every 15 minutes by vortexing or flicking the tube. Failure to mix the transformation reaction every 15 minutes will result in decreased transformation efficiency.
	5.	Heat shock the cells in a 42°C heat block or water bath for 10 minutes.
	6.	Split the cells into 2 microcentrifuge tubes (approximately 525 $\mu$ l per tube) and add 1 ml of YPD medium to each tube.
	7.	Incubate the cells at 30°C for 1 hour to allow expression of Zeocin <sup>™</sup> resistance.
	8.	Pellet the cells by centrifugation at $3000 \text{ x}$ g for 5 minutes at room temperature. Discard the supernatant.
	9.	Resuspend each tube of cells in 500 $\mu l$ of Solution III and combine the cells into one tube.
	10.	Pellet the cells by centrifugation at $3000 \text{ x}$ g for 5 minutes at room temperature. Discard the supernatant.
	11.	Resuspend the cell pellet in 100 to 150 $\mu$ l of Solution III.
	12.	Plate the entire transformation on appropriate selection plates using a sterile spreader. Incubate the plates for 2 to 4 days at 30°C. Each transformation should yield approximately 50 colonies.

#### EasyComp<sup>™</sup> Transformation, Continued

#### Analyzing *Pichia* Transformants

Select 6-10 of your Zeocin<sup>™</sup>-resistant *Pichia* transformants and confirm the Mut phenotype as described on page 22. You may also wish to analyze for the presence of insert using PCR (page 55).

#### Troubleshooting

The table below provides solutions to possible problems you may encounter when preparing and transforming competent *Pichia pastoris* cells.

Problem	Probable Cause	Possible Solution
Low efficiency of transformation	The pH of Solution I or Solution III may have drifted. The pH of both solutions should be 8.0	Check the pH of Solutions I and III. If the pH is low, increase it by adding NaOH. If the pH is high, decrease it by adding HCl. Store solutions at +4°C in order to minimize drift in pH.
	Transformation reaction not mixed during incubation	Be sure to mix the transformation reaction every 15 minutes throughout the 1 hour incubation at 30°C. Vortexing works best.
	Incubation time is too short or temperature is too low	<i>Pichia pastoris</i> transformations may be incubated for longer periods of time (up to 3 hours) and at higher temperature (35-37°C). This may, in some instances, result in higher transformation efficiencies.
	Cell density is too low $(OD_{600} < 0.6)$	Resuspend cells from Preparation of Competent Cells, Step 6, page 19, in a smaller volume (i.e. 500 µl).

## **Determining the Mut Phenotype**

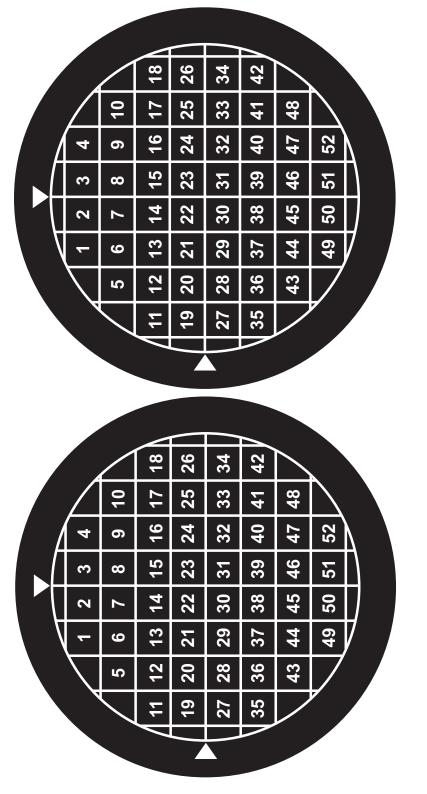
Introduction	To verify the Mut phenotype of your Zeocin <sup><math>TM</math></sup> -resistant transformants, refer to the guidelines below. <b>Note</b> : KM71H recombinants do not need to be screened for their Mut phenotype as they all will be Mut <sup>S</sup> .		
Screening for Mut <sup>+</sup> in X-33 and GS115	Transformation of X-33 or GS115 with linearized constructs favor single crossover recombination at the <i>AOX1</i> locus and result in Mut <sup>+</sup> transformants. However, with the presence of the <i>AOX1</i> sequences in the plasmid, there is a chance that recombination will also occur in the 3'AOX1 region, disrupting the wild-type $AOX1$ gene and creating Mut <sup>S</sup> transformants. Testing on MDH and MMH plates will allow you to confirm the Mut <sup>+</sup> phenotype (see below).		
Mut <sup>S</sup> in KM71H	All Zeo <sup>R</sup> transformants in KM71H will be Mut <sup>S</sup> because of the disruption of the $AOXI$ gene ( $aox1::ARG4$ ). There is no need to test recombinants for the Mut phenotype. Transformants need to be purified on minimal plates to ensure pure clonal isolates before either testing for expression (see page 28) or confirming integration by PCR (page 55).		
	Host strains containing the <i>his4</i> allele (e.g. GS115) and transformed with the pPICZ-E or pPICZ $\alpha$ -E vectors require histidine when grown in minimal media. Add histidine to a final concentration of 0.004% to ensure growth of your transformant.		
•	We recommend purifying your Zeo <sup>R</sup> transformants to ensure isolation of a pure clonal isolates. This is done by streaking for single colonies on YPD or minimal plates with histidine. You may do this before or after testing for the Mut phenotype.		
Preparation	The following media (see page 42-43) and materials can be prepared several days in advance and stored at $+4^{\circ}C$ :		
	<ul> <li>Minimal Dextrose with histidine (MDH) agar plates, 1 liter</li> <li>Minimal Methanol with histidine (MMH) agar plates, 1 liter</li> <li>Sterile toothpicks and Scoring Templates (see page 24).</li> </ul>		
Mut+ in GS115 or X-33	Use the plates containing the Zeo <sup>R</sup> transformants and confirm the Mut <sup>+</sup> phenotype as described below. <b>Note</b> : Instructions are for GS115 strains. These strains require histidine for growth. X-33 <b>does not</b> require histidine for growth, so you may leave it out of the medium.		
	1. Pick one colony using a sterile toothpick and streak/patch one Zeo <sup>R</sup> transformant in a regular pattern first on an MMH plate and then on an MDH plate.		
	3. Use a new toothpick for each transformant and continue until 10 transformants have been patched (1 plate).		
	4. To differentiate Mut <sup>+</sup> from Mut <sup>s</sup> , make one patch for each of the controls (GS115/ Albumin Mut <sup>s</sup> and GS115/pPICZ/ <i>lacZ</i> Mut <sup>+</sup> ) onto the MDH and MMH plates.		
	5. Incubate the plates at 30°C for 2 days.		
	6. After >2 days at 30°C, score the plates (see page 24). Mut <sup>+</sup> strains will grow normally on both plates, while Mut <sup>S</sup> strains will grow normally on the MDH plate but show little or no growth on the MMH plate.		

## Determining the Mut Phenotype, Continued

Replica-Plating Procedure	<ul> <li>This procedure gives a lower rate of misclassifications, but it increases the overall Mut<sup>+</sup>/Mut<sup>S</sup> screening procedure by 2 days. You will need equipment to replica-plate.</li> <li>1. Using sterile toothpicks, patch 10 Zeo<sup>R</sup> transformants on an MDH plate.</li> <li>2. Incubate the plate at 28-30°C for 2 days.</li> <li>3. After 2 days, replica-plate the patches from the MDH plate onto fresh MMH and MDH plates to screen for Mut<sup>S</sup> transformants.</li> <li>4. Incubate the replica plates at 28-30°C for 2 days.</li> <li>5. After 2 days at 28-30°C, score the replica plates. Look for patches that grow normally on the MDH replica plate but show little or no growth on the MMH replica plate.</li> </ul>	
Screening by Functional Assay	Some researchers have used a functional assay to directly screen for high expressing <i>Pichia</i> recombinant clones without first screening for Mut <sup>S</sup> or Mut <sup>+</sup> phenotypes. If you elect to screen directly for high-expressing recombinants, be sure to also check the Mut phenotype. This will help you optimize expression of your recombinant clone.	
Multiple Integration Events	<i>Pichia pastoris</i> is capable of integrating multiple copies of transforming DNA via recombination into the genome at sites of sequence homology (see page 52 for figure). Although the exact mechanism of multiple integration events is not fully understood, such events are reasonably common among selected transformants.	
	Successful expression of the gene of interest to useful levels may depend upon the generation of a recombinant strain that contains multiple copies integrated at the <i>AOX1</i> locus. In addition to simply screening expression levels among several Mut <sup>S</sup> or Mut <sup>+</sup> recombinants via SDS-PAGE analysis, it may be desirable to determine the existence of strains that have multiple integrants in the Mut <sup>S</sup> or Mut <sup>+</sup> recombinant strain.	
	See page 58 for methods to determine copy number.	
The Next Step	After confirming the Mut phenotype, you may proceed to small-scale expression (page 25) to test for expression of your gene.	

#### **Determining the Mut Phenotype, Continued**

Scoring Templates



#### Expression in Pichia

Western Analysis

Functional assay

Pichia c not be o (e.g. SD determin expressi		Pichia cl not be op (e.g. SDS determin expressio	ary purpose of small-scale expression is t one that is expressing the correct protein. timal for your protein. For this reason, the S-PAGE, western, or functional assay) ma- ing the success of expression. If your met on, you may want to consider using a more ositive clone has been identified, large-sc	Small-scale expression conditions may e method you choose for detection y be an important factor in hod of detection does not reveal any e sensitive method.
			sk or fermentation, and expression condit	
Control	Strains	into a <i>Pia</i> expression maximum Blue stain Kit, Cata cannot be	positive control strains, pPICZ-E/Uni-lacZ chia host, which you made in the previous on. Expression in shake flasks is detectabl n at 96 hours (4 days). $\beta$ -galactosidase (11 ning of the SDS-PAGE gel or performing log no. K1455-01). <b>Note</b> : $\beta$ -galactosidase e detected using either anti-V5 or anti-His ow, then use anti- $\beta$ -galactosidase antibody	section (see page 16), to test for e after 48 hours for <i>lacZ</i> and reaches a l6 kDa) can be detected by Coomassie the $\beta$ -galactosidase assay ( $\beta$ -Gal Assay does not contain a C-terminal tag and (C-term) antibodies. If the expression
		maximur	kDa) expression starts as early as 24 hou n by 96 hours as determined by Coomassi ed to do a western analysis for detection,	e <sup>®</sup> Blue staining of the SDS/PAGE gel.
		for intrac β-galacto	also use the strains, GS115/pPICZ/lacZ of cellular and secreted expression, respective sidase (121 kDa) fused at the C-terminus 15/Albumin secretes albumin (67 kDa) in	ely. GS115/pPICZ/ <i>lacZ</i> expresses to the <i>myc</i> epitope and the polyhistidine
	Note	your reco	a have obtained Zeocin <sup>™</sup> -resistant transfor ombinant <i>Pichia</i> clone in medium contain is only required for initial screening and s	ng Zeocin <sup><math>^{TM}</math></sup> for expression studies.
Detectio	n	We recon protein.	nmend that you use the following techniq	ues to assay expression of your
	Techniq	lne	Method of Detection	Sensitivity
SDS-PAGE (Coomassie-stain		ained)	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

Antibody to your particular protein

Anti-V5 antibodies (see the next page)

Anti-His(C-term) antibodies (see the

Varies depending on assay

next page)

Can detect as little as 1-10 pg,

Varies depending on assay

depending on detection method

(alkaline phosphatase, horseradish

peroxidase, radiolabeled antibody)

Important	Host strains containing the <i>his4</i> allele (e.g. GS115) and transformed with the pPICZ-E or pPICZα-E vectors require histidine when grown in minimal media. Add histidine to a final concentration of 0.004% to ensure growth of your transformant. Complex media such as YPD and BMGY already contain histidine.
Media	You will need either BMGY/BMMY (buffered complex glycerol or methanol medium), BMGH/BMMH (buffered minimal glycerol or methanol medium containing histidine) or MGYH/MMH (minimal glycerol or minimal methanol medium containing histidine) for expression (see <b>Recipes</b> , pages 43-44). BMGH, BMMH, BMGY, and BMMY are usually used for the expression of secreted proteins, particularly if pH is important for the activity of your protein. Unlike MGYH and MMH, they are all buffered media. Because these media are buffered with phosphate buffer, a wide range of pH values may be used to optimize production of your protein. BMGY/BMMY contain yeast extract and peptone to stabilize secreted proteins and prevent or decrease proteolysis of secreted proteins. Inclusion of yeast extract and peptone allow better growth and biomass accumulation.
Proteases	There are some proteins specifically susceptible to proteases that have optimal activity at neutral pH. If this is the case, use MGYH or MMH media for expression. As <i>Pichia</i> expression progresses in an unbuffered medium such as MMH, the pH drops to 3 or below, inactivating many neutral pH proteases (Brierley et al., 1994). <i>Pichia</i> is resistant to low pH, so the low pH will not affect growth. In contrast, it has been reported that by including 1% Casamino acids (Difco) and buffering the medium at pH 6.0, extracellular proteases were inhibited, increasing the yield of mouse epidermal growth factor (Clare et al., 1991b).
	If your secreted protein of interest is susceptible to neutral pH proteases you may prefer to do your expressions in an unbuffered medium (MMH). If your secreted protein of interest is not susceptible to proteases at neutral pH, then you can use BMMY for initial expression. Try expression in an unbuffered medium if the expressed protein is degraded.
Aeration	The most important parameter for efficient expression in <i>Pichia</i> is adequate aeration during methanol induction. As a general rule when inducing expression, never allow cultures to be more than 10-30% of your total flask volume. It is strongly recommended that baffled flasks be used. See page 44 for suppliers of baffled flasks. Cover the flasks with cheesecloth (2-3 layers) or another loose fitting cover. Never use tight fitting covers. (Aeration is not as critical when generating biomass before induction.)
Kinetics of Growth	Note that while Mut <sup>+</sup> and Mut <sup>S</sup> strains will grow at essentially the same rate in YPD or glycerol media, Mut <sup>+</sup> will grow faster than Mut <sup>S</sup> when both are grown on methanol because of the presence of the <i>AOX1</i> gene product.
Temperature and Shaking	All expression is done at 30°C, in a shaking incubator. It is critical that the temperature does not exceed 30°C. If your incubator temperature fluctuates, set the temperature at 28°C. If using a floor shaking incubator, shake at 225-250 rpm. If using a table-top shaker that sits inside an incubator, shake at 250-300 rpm.

Signal Sequence	The processing of the $\alpha$ -factor signal sequence in pPICZ $\alpha$ -E	occurs in two steps:	
Processing	1. The preliminary cleavage of the signal sequence by the <i>K</i> final Kex2 cleavage occurring between arginine and glut. Lys-Arg * Glu-Ala-Glu-Ala, where * is the site of cleavage	amine in the sequence Glu-	
	2. The Glu-Ala repeats are further cleaved by the <i>STE13</i> ge	ne product.	
Optimizing Signal Cleavage	In <i>Saccharomyces cerevisiae</i> , it has been noted that 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 et al., 1984).		
	There are some cases where Ste13 cleavage of Glu-Ala repeat Ala repeats are left on the N-terminus of the expressed protein generally dependent on the protein of interest.		
Before Starting	When performing your expression, it is important to run the p will be able to interpret your expression results. The expression used are:		
	pPICZα-E/Uni-HSA in <i>Pichia</i> strain of your choice	Secretion control	
	pPICZ-E/Uni-lacZ in Pichia strain of your choice	Intracellular control	
	$GS115/pPICZ/lacZ Mut^+$	Mut <sup>+</sup> -Intracellular control	
	GS115/Albumin Mut <sup>s</sup>	Mut <sup>s</sup> -Secretion control	
	Appropriate Pichia strain /Vector (no insert)	Background control	
	Since recombination can occur in many different ways that ca variation), we recommend you screen 6-10 verified recombin levels. Start with fresh colonies as colony viability drops over the cultures with a small sample from a frozen glycerol stock single colony).	ant clones for expression r time. (You may also start	
Guidelines for Expression	The following steps should be viewed as guidelines and are p with expression. You may have to change the conditions to op particular protein. Use bottom or side baffled flasks whenever 2000 ml sizes). If you are analyzing a number of recombinant tubes. Be sure that the medium is well-aerated by increasing t placing the tubes at an angle in the shaker.	ptimize expression for your r possible (available in 50 - ts, you can try 50 ml conical	

Mut <sup>+</sup> Intracellular or Secreted	To evaluate expression, we recommend including the control strains discussed earlier (see page 16). Include X-33, GS115, or KM71H transformed with the parent vector as a control for background expression.
	1. Inoculate a single colony into 25 ml of MGYH, BMGH, or BMGY in a 250 ml baffled flask. Grow at 28-30°C in a shaking incubator (250-300 rpm) until culture reaches log phase ( $OD_{600} = 2$ -6) (approximately 16-18 hours).
	<ol> <li>Harvest the cells by centrifuging at 1500-3000 x g for 5 minutes at room temperature. Decant supernatant and resuspend cell pellet to an OD<sub>600</sub> of 1.0 in MMH, BMMH, or BMMY medium to induce expression (approximately 100- 200 ml).</li> </ol>
	3. Place culture in a 1 liter baffled flask. Cover the flask with 2 layers of sterile gauze or cheesecloth and return to incubator to continue growth.
	4. Add 100% methanol to a final concentration of 0.5% methanol every 24 hours to maintain induction.
	5. At each of the time points (below), transfer 1 ml of the expression culture to a 1.5 ml microcentrifuge tube. These samples will be used to analyze expression levels and determine the optimal time post-induction to harvest. Centrifuge at maximum speed in a tabletop microcentrifuge for 2-3 minutes at room temperature.
	Time points (hours): 0, 6, 12, 24 (1 day), 36, 48 (2 days), 60, 72 (3 days), 84, and 96 (4 days).
	6. For secreted expression, <b>transfer the supernatant to a separate tube.</b> Store the supernatant and the cell pellets at -80°C until ready to assay. Freeze quickly in liquid N <sub>2</sub> or a dry ice/alcohol bath.
	For intracellular expression, decant the supernatant and store just the cell pellets at $-80^{\circ}$ C until ready to assay. Freeze quickly in liquid N <sub>2</sub> or a dry ice/alcohol bath.
	7. Analyze the supernatants and cell pellets for protein expression by Coomassie- stained SDS-PAGE and western blot or functional assay (see page 30).
	Continued on next page

Mut <sup>s</sup> Intracellular or Secreted	To evaluate expression, we recommend including the control strains discussed earlier (see page 16). Include X-33, GS115 or KM71H transformed with the parent vector as a control for background expression.			
	1.	Using a single colony, inoculate 100 ml of MGYH, BMGH, or BMGY in a 1 liter baffled flask. Grow at 28-30°C in a shaking incubator (250-300 rpm) until the culture reaches an $OD_{600} = 2-6$ (approximately 16-18 hours.)		
	2.	Harvest the cells by centrifuging at 1500-3000 x g for 5 minutes at room temperature. To induce expression, decant the supernatant, and resuspend cell pellet in MMH, BMMH, or BMMY medium using 1/5 to 1/10 of the original culture volume (approximately 10-20 ml).		
	3.	Place in a 100 ml baffled flask. Cover the flask with 2 layers of sterile gauze or cheesecloth and return to incubator to continue to grow.		
	4.	Add 100% methanol to a final concentration of 0.5% every 24 hours to maintain induction.		
	5.	At each of the time points (below) transfer 1 ml of the expression culture to a 1.5 ml microcentrifuge tube. These samples will be used to analyze expression levels and determine the optimal time post-induction to harvest. Centrifuge at maximum speed in a tabletop microcentrifuge for 2-3 minutes at room temperature.		
		Time points (hours): 0, 24 (1 day), 48 (2 days), 72 (3 days), 96 (4 days), 120 (5 days), and 144 (6 days).		
	6.	For secreted expression, <b>transfer the supernatant to a separate tube.</b> Store the supernatant and the cell pellets at $-80^{\circ}$ C until ready to assay. Freeze quickly in liquid N <sub>2</sub> or a dry ice/alcohol bath.		
		For intracellular expression, decant the supernatant, and store just the cell pellets at -80°C until ready to assay. Freeze quickly in liquid N <sub>2</sub> or a dry ice/alcohol bath.		
	7.	Analyze the cell pellets for protein expression by Coomassie-stained SDS-PAGE and western blot or functional assay (see <b>Analysis by SDS-Polyacrylamide Gel</b> <b>Electrophoresis</b> , next page).		
	The	e C-terminal tag will add 2.5 kDa to the size of your protein and the $\alpha$ -factor signal		

Note

The C-terminal tag will add 2.5 kDa to the size of your protein and the  $\alpha$ -factor signal sequence will add 9.3 kDa to the size of the protein, if it is not processed. Be sure to account for any additional amino acids that are in between the signal sequence site and the N-terminus of your protein, and also the end of your protein and the C-terminal tag.

## Analysis by SDS-Polyacrylamide Gel Electrophoresis

Introduction	polyacrylamide gel electrophoresis.         To facilitate separation and visualization of your recombinant protein by polyacrylamide gel electrophoresis, a wide range of pre-cast NuPAGE <sup>®</sup> and Tris-Glycine polyacrylamide		
Polyacrylamide Gel Electrophoresis			
Preparing Samples	You will need to prepare Breaking Buffer (see page 44) and have acid-washed 0.5 mm glass beads on hand.		
	Preparation of cell pellets (Intracellular and Secreted Expression):		
	1. Thaw cell pellets quickly and place on ice.		
	2. For each 1 ml sample, add 100 µl Breaking Buffer to the cell pellet and resuspend.		
	3. Add an equal volume of acid-washed glass beads (size 0.5 mm). Estimate equal volume by displacement.		
	4. Vortex 30 seconds, then incubate on ice for 30 seconds. Repeat for a total of 8 cycles.		
	5. Centrifuge at maximum speed for 10 minutes at +4°C. Transfer the clear supernatant to a fresh microcentrifuge tube.		
	<ol> <li>Take 50 μl of supernatant (cell lysate) and mix with 50 μl 2X SDS-PAGE Gel Loading buffer (Sample Buffer).</li> </ol>		
	<ol> <li>Boil for 10 minutes and load 10-20 μl per well. Thickness of the gel and number of wells will determine volume loaded. Remaining sample may be stored at -20°C for western blots, if necessary. Cell lysates may be stored at -80°C for further analysis.</li> </ol>		
	Preparation of supernatant (Secreted Expression only):		
	1. Thaw supernatants and place on ice.		
	2. Mix 50 µl of the supernatant with 50 µl of SDS-PAGE Gel Loading buffer.		
	3. Boil 10 minutes, then load 10-30 μl onto the gel. Remaining sample may be stored at -20°C for western blots, if necessary. Supernatants may be stored at -80°C for further analysis.		
	4. If no protein is seen by Coomassie or by western blot, then concentrate the supernatant 5-10 fold and analyze samples again by western blot. Centricon and Centriprep filters (Amicon) are very useful for this purpose.		

#### Analysis by SDS-Polyacrylamide Gel Electrophoresis, Continued

Protein Concentration	Lowry, BCA (Pierce) or Bradford protein determinations can be performed to quantify the amounts of protein in the cell lysates and medium supernatants. In general, <i>Pichia</i> cell lysates contain 5-10 $\mu$ g/ $\mu$ l protein. <i>Pichia</i> medium supernatants will vary in protein concentration primarily due to the amount of your secreted protein. <i>Pichia</i> secretes very few native proteins. If the protein concentration of the medium is > 50 $\mu$ g/ml, 10 $\mu$ l of medium will give a faint band on a Coomassie-stained SDS-PAGE gel.
Controls	Include the following samples as controls on your SDS-PAGE:
	<ul> <li>Molecular weight standards appropriate for your desired protein</li> <li>A sample of your protein as a standard (if available)</li> <li>A sample of X-33, GS115, or KM71H with the parent plasmid transformed into it. Inclusion of this sample shows the background of native <i>Pichia</i> proteins that are present intracellularly and will help you differentiate your protein from background</li> <li>Analyze the <i>lac</i>Z and HSA controls also as they should indicate any problems with the media or expression conditions</li> </ul>
	In addition to Coomassie-stained SDS-PAGE, we strongly recommend performing a western blot or another more sensitive assay to detect your protein. Visualization of the expressed protein will depend on several factors, including its expression level, its solubility, its molecular weight, and whether it will be masked by an abundant cellular protein of the same size.
Western Blot 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 available from Invitrogen (see page xiii for ordering information) or an antibody to your protein of interest. The ready-to-use WesternBreeze <sup>™</sup> Chromogenic Kits and WesternBreeze <sup>™</sup> Chemiluminescent Kits are available from Invitrogen to facilitate detection of antibodies by colorimetric or chemiluminescent methods. The Positope <sup>™</sup> Control Protein from Invitrogen can be used a s a positive control for detection of fusion proteins containing a V5 epitope and a polyhistidine tag. For more information, refer to our Web site (www.invitrogen.com) or call Technical Service (see page 61).
Analysis of Protein Expression	The Coomassie-stained SDS-PAGE gel should reveal the induction over time of your recombinant protein co-migrating with your standard. If you are satisfied with the level of expression, try a test purification (page 36) or scale-up expression (page 34).
•	If there is no recombinant protein visible, then perform either a western blot or a functional assay if available.
	If you detect low expression of your recombinant protein, see <b>Optimization of</b> <i>Pichia</i> <b>Protein Expression</b> , page 31, for guidelines to optimize expression.
	Test your expression conditions with the control strains included in the kit (see page 25).
	If there is no indication of expression at all, use PCR to analyze your recombinants for the correctly sized PCR product (page 55). If you find that you have recombinants, perform a northern analysis to see if and how much full-length mRNA is induced (see page 60).

## Optimization of *Pichia* Protein Expression

Introduction	Based on available data, there is approximately a 50 to 75% chance of expressing your protein of interest in <i>Pichia pastoris</i> at reasonable levels. The biggest hurdle seems to be generating initial success-i.e. expression of your protein at <u>any</u> level. While there are relatively few examples of expression of >10 grams/liter, there are many examples of expression in the >1 gram/liter range, making the <i>Pichia pastoris</i> expression system one of the most productive eukaryotic expression systems available. Likewise, there are several examples of proteins that have been successfully expressed in <i>Pichia pastoris</i> that were completely unsuccessful in baculovirus or <i>Saccharomyces cerevisiae</i> , suggesting that the <i>Pichia pastoris</i> system is an important alternative to have available. If you obtain no or low protein expression in your initial expression experiment, use the following guidelines to optimize expression.
Proteolysis or Degradation	• Do a time course study of expression. Check to see if there is a time point that yields a larger percentage of full-length protein.
	• If secreting your protein, check to see if your protein is susceptible to neutral pH proteases by expressing in unbuffered medium (MMH). In addition, try 1% Casamino acids with buffered medium to inhibit extracellular proteases.
Low Secreted Expression Levels	• Check cell pellet to see if overall expression is low or if the protein did not secrete. If it did not secrete, try a different signal sequence (e. g. a native or α-factor signal sequence).
	• Concentrate your supernatant by ammonium sulfate precipitation or ultrafiltration (see page 35).
	• For Mut <sup>+</sup> , induce expression with a higher density culture.
Low Expression Levels	• Look for multi-copy recombinants (i.e. jackpot clones) by dot blot (see page 57). There are quite a few examples of increasing the expression levels of a particular protein by increasing the gene dosage. See Clare, 1991a; Clare <i>et al.</i> , 1991b; Romanos <i>et al.</i> , 1991).
	• Check both Mut <sup>+</sup> and Mut <sup>S</sup> recombinants for increased expression. Some proteins express better in one type of genetic background than another.
	• If secreting your protein, try intracellular expression. The protein may not be processed correctly and fail to secrete. Be sure you check your cell pellets for evidence of expression. If you are having problems with intracellular expression, try secreting your protein. It probably will glycosylate which may be desirable or not. If glycosylation is undesirable, oligosaccharides can be removed with Peptide:N-Glycosidase F (New England BioLabs).
	• Scale up to fermentation (page 35). <i>Pichia</i> is a yeast and is particularly well suited to fermentation.

## Optimization of *Pichia* Protein Expression, Continued

No Expression	Be sure to try some of the things listed above as no expression can be the same thing as very low expression. If none of these things improve protein expression, use PCR to check for insertion of your gene into the Pichia genome (page 55). If your gene is present, perform a northern blot analysis to check for transcription of your gene. There is a protocol in the <b>Appendix</b> for RNA isolation from <i>Pichia</i> (see page 60).
	If you see premature transcriptional termination, check the AT content of your gene. In <i>Saccharomyces</i> , there are a few consensus sequences which promote premature termination. One of these, TTTTTATA, resembles a sequence in HIV-1 gp120, ATTATTTTAT AAA, which when expressed in <i>Pichia</i> gave premature termination of the mRNA. When this sequence was changed, longer transcripts were found (Scorer <i>et al.</i> , 1993).
Hyperglycosy-	If your protein is hyperglycosylated:
lation	• Try intracellular expression as your protein will not go through the secretion pathway and therefore, not be modified.
	• Try deglycosylating the protein with Peptide:N-Glycosidase F or other enzymes.

## Scale-up of Expression

Introduction	prot	e expression is optimized, the expression protocol can be scaled up to produce more ein by increasing the culture volume using larger baffled flasks or fermentation. Use guidelines below to scale-up your expression protocol.
Mut <sup>+</sup> Intracellular or Secreted	1.	Using a single colony, inoculate 25 ml of MGYH, BMGH, or BMGY in a 250 ml baffled flask. Grow at 28-30°C in a shaking incubator (250-300 rpm) until culture reaches an $OD_{600} = 2-6$ (approximately 16-18 hours).
	2.	Use this 25 ml culture to inoculate 1 liter of MGYH, BMGH, or BMGY in a 3 or 4 liter baffled flask and grow at 28-30°C with vigorous shaking (250-300 rpm) until the culture reaches log phase growth (OD <sub>600</sub> = 2-6).
	3.	Harvest the cells using sterile centrifuge bottles by centrifuging at 1500-3000 x g for 5 minutes. To induce expression, decant the supernatant and resuspend cell pellet to an $OD_{600} = 1.0$ (2-6 liters) in MMH, BMMH, or BMMY medium to start induction.
	4.	Aliquot the culture between several 3 or 4 liter baffled flask. Cover the flasks with 2 layers of sterile gauze or cheesecloth and return to incubator. Continue to grow at 28-30°C with shaking.
	5.	Add 100% methanol to 0.5% every 24 hours until the optimal time of induction is reached as determined from the time course study.
	6.	Harvest cells by centrifuging at 1500-3000 x g for 5 minutes at room temperature.
		For intracellular expression, decant the supernatant. Cells can be processed immediately or stored at -80°C until ready for use.
		For secreted expression, <b>save the supernatant, chill to</b> +4°C, <b>and concentrate it if desired</b> (see next page). Proceed directly to purification (page 36) or store the supernatant at -80°C until ready to process further.
Mut <sup>s</sup> Intracellular or Secreted	1.	Using a single colony, inoculate 10 ml of MGYH, BMGH, or BMGY in a 100 ml baffled flask. Grow at 28-30°C in a shaking incubator (250-300 rpm) to an $OD_{600} = 2-6$ (approximately 16-18 hours).
	2.	Use this 10 ml culture to inoculate 1 liter of MGYH, BMGH, or BMGY in a 3 or 4 liter baffled flask and grow at 28-30°C with vigorous shaking (250-300 rpm) until the culture reaches log phase growth (OD <sub>600</sub> = 2-6).
	3.	Harvest the cells by centrifuging at 1500-3000 x g for 5 minutes. To induce expression, decant the supernatant and resuspend cell pellet in 1/5 to 1/10 of the original culture volume of MMH, BMMH, or BMMY medium (~ 100-200 ml).
	4.	Place the culture in a 1 liter baffled flask. Cover the flask with 2 layers of sterile gauze or cheesecloth and return to incubator. Continue to grow at 28-30°C with shaking.
	5.	Add 100% methanol to 0.5% every 24 hours until the optimal induction time is reached.
	6.	Harvest cells by centrifuging at 1500-3000 x g for 5 minutes at room temperature.
		For intracellular expression, decant the supernatant and store the cell pellets at -80°C until ready to process.
		For secreted expression, save the supernatant, chill to $+4^{\circ}C$ , and concentrate it if desired (see next page). Proceed directly to purification (page 36) or store the supernatant at -80°C until ready to process further.

#### Scale-up of Expression, Continued

To increase the amount of cells for Mut<sup>S</sup> recombinants, increase the number of flasks, put 200-300 ml in a 3 liter flask, or try fermentation.

Concentration of Proteins	Proteins secreted into the media are usually > 50% homogeneous and will require some additional purification (see page 35). It is optimal to concentrate the protein if the expression level is not particularly high. There are several general methods to concentrate proteins secreted from <i>Pichia</i> . A general guide to protein techniques is <i>Protein Methods</i> (Bollag & Edelstein, 1991). Some general methods include:
	Ammonium sulfate precipitation
	• Centrifuge concentrator for small volumes (e.g. Centricon or Centriprep devices available from Amicon)
	• Pressurized cell concentrators for large volumes (Amicon ultrafiltration devices)
	Lyophilization
Cell Lysis	A general procedure for cell lysis using glass beads is provided on the next page. There is also a cell lysis protocol in <i>Current Protocols in Molecular Biology</i> , page 13.13.4. (Ausubel <i>et al.</i> , 1994) and in <i>Guide to Protein Purification</i> (Deutscher, 1990). We also recommend lysis by French Press (follow the manufacturer's suggestions for yeast).
Fermentation	Basic guidelines are available for fermentation of <i>Pichia</i> from Invitrogen. We recommend that only those with fermentation experience or those who have access to people with experience attempt fermentation. Call Technical Service (see page 61) for more information or you may download the manual from our Web site at (www.invitrogen.com).

#### 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 <sup>TM</sup> . You may harvest the cells and store them at -80°C until you are ready to purify your fusion protein, or you may proceed directly with protein purification. <b>This section only describes the preparation of cell lysates and sample application onto ProBond<sup>TM</sup></b> . For instructions on how to prepare and use ProBond <sup>TM</sup> resin, refer to the ProBond <sup>TM</sup> Purification manual.
ProBond <sup>™</sup> Resin	We recommend that you use the ProBond <sup>TM</sup> Purification System (Catalog no. K850-01) to purify fusion proteins expressed from pPICZ-E or pPICZ $\alpha$ -E. Note that instructions for equilibration of and chromatography on ProBond <sup>TM</sup> resin are contained in the ProBond <sup>TM</sup> Purification Kit. You may download the manual from our Web site (www.invitrogen.com) or call Technical Service (see page 61).
	If you are using a metal-chelating resin other than $ProBond^{TM}$ , follow the manufacturer's recommendations to purify fusion proteins expressed in bacteria or yeast.
Binding Capacity of ProBond <sup>™</sup>	One milliliter of ProBond <sup>™</sup> resin binds at least 1 mg of recombinant protein. This amount can vary depending on the protein.
Important	Throughout the following protocol, be sure to keep the cell lysate and fractions on ice. Small-scale purifications using the 2 ml ProBond <sup>TM</sup> columns and buffers can be performed at room temperature on the bench top. For large scale purifications, all reagents must be kept at $+4^{\circ}$ C.
Ordering Information	ProBond <sup>™</sup> resin is available precharged and prepacked in 2 ml columns or in bulk quantities. For ordering information, refer to page xiv.

## **Purification**, Continued

Preparation of Cell Lysates	Express your protein using a small-scale culture (10-20 ml for Mut <sup>S</sup> strains; 100-200 ml for Mut <sup>+</sup> ) and the optimal conditions for expression (if determined). Once your protein is expressed, follow the protocol below to prepare a cell lysate for chromatography on ProBond <sup>TM</sup> .		
	Prepare Breaking Buffer (BB) as described in the <b>Recipes</b> , page 44. Do not add 1 mM EDTA to the breaking buffer as it will interfere with binding to Probond <sup><math>\mathbb{M}</math></sup> .		
	1. Wash cells once in BB by resuspending them and centrifuging 5-10 minutes at 3000 x g at +4°C.		
	2. Resuspend the cells to an $OD_{600}$ of 50-100 in BB.		
	3. Add an equal volume of acid-washed glass beads (0.5 mm). Estimate volume by displacement.		
	4. 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 $+4^{\circ}$ C for 5-10 minutes at 12,000 x g.		
	6. 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.		
Sample Application	For sample application onto ProBond <sup>TM</sup> , you will need Native Binding Buffer, pH 7.8 and a 2 ml ProBond <sup>TM</sup> column, pre-equilibrated using native conditions.		
(Native Conditions)	1. Combine 1 ml (2-3 mg/ml total protein) of <i>Pichia</i> lysate with 7 ml Native Binding Buffer.		
	2. Take a pre-equilibrated ProBond <sup>™</sup> column and resuspend the resin 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 x g) and carefully remove the supernatant. Save the supernatant to check for unbound protein.		
	<ol> <li>Repeat Steps 2 through 4 with the remaining 4 ml of diluted lysate. Proceed to Column Washing and Elution Under Native Conditions in the ProBond<sup>™</sup> Purification manual. Use the recommendations noted for bacterial cell lysates.</li> </ol>		
Sample Applica- tion (Denaturing Conditions)	Use the protocol above except pre-equilibrate the ProBond <sup><math>TM</math></sup> column using Denaturing Binding Buffer and combine 1 ml of the <i>Pichia</i> cell lysate with 7 ml of the Denaturing Binding Buffer.		

## **Purification**, Continued

Note	We have observed that some <i>Pichia</i> proteins may be retained on the ProBond <sup>TM</sup> column using native purification conditions. Optimization of the purification (see ProBond <sup>TM</sup> Purification manual) or using denaturing purification may remove these non-specific <i>Pichia</i> proteins.
Analysis of Purification	Be sure to 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 <sup>™</sup> Purification 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 with 1/10 volume of 10X Stock Solution B (ProBond <sup>™</sup> Purification Kit) 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. Using 10X Stock Solution B to adjust the pH and the ionic strength keeps the total volume small for sample application.

#### Pichia Media Recipes

#### Using *Pichia* Media

The table below is designed to help you decide which *Pichia* medium to use for a particular application.

Medium	Description	Application
YPD or YEPD	Rich, complex broth	General growth and storage
$YPDS + Zeocin^{TM}$	YPD with sorbitol and Zeocin <sup>™</sup>	Selection of Pichia Zeo <sup>R</sup> transformants
MGYH or MGY	Minimal medium containing glycerol and/or histidine	Intracellular Expression: Generation of biomass prior to methanol induction
MDH or MD	Minimal medium containing glucose and/or histidine	Determination of Mut phenotype
MMH or MM	Minimal medium containing methanol and/or histidine	Determination of Mut phenotype
		Intracellular expression of desired protein
BMGH or BMG	Buffered minimal medium containing glycerol and/or histidine	Secreted Expression: Used to control the pH of the medium and generate biomass
BMMH or BMM	Buffered minimal medium containing methanol and/or histidine	Secreted Expression: Used to control the pH of the medium and induce expression of the desired protein
BMGY	Buffered complex medium containing glycerol	Secreted Expression: Used to control the pH of the medium, decrease protease activity, and generate biomass.
BMMY	Buffered complex medium containing methanol	Secreted Expression: Used to control the pH of the medium, decrease protease activity, and induce expression.

#### **Stock Solutions**

## **10X YNB (13.4% Yeast Nitrogen Base with Ammonium Sulfate without amino acids)**

Dissolve 134 g of yeast nitrogen base (YNB) with 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. Alternatively, use 34 g of YNB without ammonium sulfate and amino acids and 100 g of ammonium sulfate. The shelf life of this solution is approximately one year. If you are using the YNB pouch included in the kit, follow the directions on the pouch. **Note:** *Pichia* cells exhibit optimal growth with higher YNB concentrations, therefore, the amount of YNB used in this kit is twice as concentrated as YNB formulations for *Saccharomyces*.

#### 500X B (0.02% Biotin)

Dissolve 20 mg biotin in 100 ml of water and filter sterilize. Store at  $+4^{\circ}$ C. The shelf life of this solution is approximately one year.

#### 100X H (0.4% Histidine)

Dissolve 400 mg of L-histidine in 100 ml of water. Heat the solution, if necessary, to no greater than  $50^{\circ}$ C in order to dissolve. Filter sterilize and store at +4°C. The shelf life of this solution is approximately one year.

#### 10X D (20% Dextrose)

Dissolve 200 g of D-glucose in 1000 ml of water. Autoclave for 15 minutes or filter sterilize. 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 GY (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.

#### 1 M potassium phosphate buffer, pH 6.0:

Combine 132 ml of 1 M K<sub>2</sub>HPO<sub>4</sub>, 868 ml of 1 M KH<sub>2</sub>PO<sub>4</sub> and confirm that the pH =  $6.0 \pm 0.1$  (if the pH needs to be adjusted, use phosphoric acid or KOH). Sterilize by autoclaving and store at room temperature. The shelf life of this solution is greater than one year.

YPD or YEPD	<ul> <li>Yeast Extract Peptone Dextrose Medium (1 liter)</li> <li>1% yeast extract</li> <li>2% peptone</li> <li>2% dextrose (glucose)</li> <li>Note: If you are using the YP Base Medium or the Y included with the Original Pichia Expression K</li> <li>Dissolve 10 g yeast extract and 20 g of peptone of agar if making YPD slants or plates.</li> <li>Autoclave for 20 minutes on liquid cycle.</li> <li>Add 100 ml of 10X D.</li> </ul>	it, follow the directions on the pouch.	
	The liquid medium is stored at room temperature. YPD slants or plates are stored at $+4^{\circ}$ C. The shelf life is several months.		
YPD (+ Zeocin™)	<ul> <li>Yeast Extract Peptone Dextrose Medium (1 liter)</li> <li>1% yeast extract</li> <li>2% peptone</li> <li>2% dextrose (glucose)</li> <li>± 2% agar</li> <li>± 100 µg/ml Zeocin<sup>™</sup></li> <li>1. Dissolve the following in 900 ml of water:</li> <li>2. Include 20 g of agar if making YPD slants or pl</li> <li>3. Autoclave for 20 minutes on liquid cycle.</li> <li>4. Cool solution to ~60°C and add 100 ml of 10X Zeocin<sup>™</sup>, if desired.</li> <li>Liquid medium without Zeocin<sup>™</sup> can be stored at root Zeocin<sup>™</sup> should be stored at +4°C in the dark. YPD store at +4°C in the dark. YPD store at the shelf life of medium is several months. Medium of one to two weeks.</li> </ul>	D. Add 1.0 ml of 100 mg/ml om temperature. Medium containing slants or plates are stored at $+4^{\circ}$ C.	

YPDS + Zeocin <sup>™</sup>	Yeast Extract Peptone Dextrose Medium (1 liter)			
Agar	<ul> <li>1% yeast extract</li> <li>2% peptone</li> <li>2% dextrose (glucose)</li> <li>1 M sorbitol</li> <li>2% agar</li> <li>100 µg/ml Zeocin<sup>™</sup></li> </ul>			
	1.	Dissolve the following in 900 ml of water:	10 g yeast extract 182.2 g sorbitol 20 g of peptone	
	2.	Add 20 g of agar.		
	3.	Autoclave for 20 minutes on liquid cycle.		
	4.	Add 100 ml of 10X D.		
	<ul> <li>5. Cool solution to ~60°C and add 1.0 ml of 100 mg/ml Zeocin<sup>™</sup>.</li> <li>Store YPDS slants or plates containing Zeocin<sup>™</sup> at +4°C in the dark. The shelf life is one to two weeks.</li> </ul>			
MGY and MGYH	<u>Minimal Glycerol Medium + Histidine</u> (1 liter) 1.34% YNB 1% glycerol 4 x 10 <sup>-5</sup> % biotin ± 0.004% histidine			
	<ol> <li>Combine aseptically 800 ml autoclaved water with 100 ml of 10X YNB, 2 ml of 500X B, and 100 ml of 10X GY.</li> </ol>			
	2. For growth of <i>his4</i> strains in this medium, a version can be made that contains histidine (called MGYH) by adding 10 ml of 100X H stock solution.			
	Store at +4°C. The shelf life of this solution is approximately two months.			
MD and MDH	<u>Minimal Dextrose Medium + H</u> istidine (1 liter) 1.34% YNB 4 x 10 <sup>-5</sup> % biotin 2% dextrose			
	1.	For medium, autoclave 800 ml of water for 2	0 minutes on liquid cycle.	
	2.	Cool to about 60°C and then add:		
		100 ml of 10X YNB 2 ml of 500X B 100 ml of 10X D		
	3.	To make MDH, add 10 ml of 100X H stock s	olution. Mix and store at +4°C.	
	4.	For plates, add 15 g agar to the water in Step	1 and proceed.	
	5.	If preparing plates, pour the plates immediate at $+4^{\circ}C$ .	ly. MD stores well for several months	

MM and MMH	<u>Minimal Methanol + H</u> istidine (1 liter) 1.34% YNB 4 x 10 <sup>-5</sup> % biotin 0.5% methanol			
	1. For medium, autoclave 800 ml of water for 20 minutes on liquid cycle.			
	2. Cool autoclaved water to 60°C and add:			
	100 ml of 10X YNB 2 ml of 500X B 100 ml of 10X M			
	3. To make MMH, add 10 ml of 100X H stock solution. Mix and store at $+4^{\circ}$ C.			
	4. For plates, add 15 g agar to the water in Step 1 and proceed.			
	5. After mixing, pour the plates immediately. MM and MMH stores well for several months at +4°C.			
BMGH and BMMH	Buffered Minimal Glycerol Buffered Minimal Methanol (1 liter)			
	100 mM potassium phosphate, pH 6.0			
	1.34% YNB 4 x 10 <sup>-5</sup> % biotin			
	1% glycerol or 0.5% methanol			
	1. Autoclave 690 ml water for 20 minutes on liquid cycle.			
	2. Cool to room temperature, then add the following and mix well:			
	100 ml 1 M potassium phosphate buffer, pH 6.0 100 ml 10X YNB 2 ml 500X B 100 ml 10X GY			
	3. For BMMH, add 100 ml 10X M instead of glycerol.			
	4. To add histidine, add 10 ml of 100X H stock solution. Mix and store at +4°C.			
	5. Store media at $+4^{\circ}$ C. The shelf life of this solution is approximately two months.			

BMGY and BMMY	<u>B</u> uffered <u>G</u> lycerol-complex Medium <u>B</u> uffered <u>M</u> ethanol-complex Medium (1 liter)				
	<ul> <li>1% yeast extract</li> <li>2% peptone</li> <li>100 mM potassium phosphate, pH 6.0</li> <li>1.34% YNB</li> <li>4 x 10<sup>-5</sup>% biotin</li> <li>1% glycerol or 0.5% methanol</li> </ul>				
	<ol> <li>Dissolve 10 g of yeast extract, 20 g peptone in 700 ml water.</li> </ol>				
	<ol> <li>Autoclave 20 minutes on liquid cycle.</li> </ol>				
	3. Cool to room temperature, then add the following and mix well:				
	100 ml 1 M potassium phosphate buffer, pH 6.0 100 ml 10X YNB 2 ml 500X B 100 ml 10X GY				
	4. For BMMY, add 100 ml 10X M instead of glycerol.				
	5. Store media at $+4^{\circ}$ C. The shelf life of this solution is approximately two months.				
Breaking Buffer	50 mM sodium phosphate, pH 7.4 1 mM PMSF (phenylmethylsulfonyl fluoride or other protease inhibitors) 1 mM EDTA (Omit EDTA if using this buffer for purification on metal chelating resins 5% glycerol				
	1. Prepare a stock solution of your desired protease inhibitors and store appropriately. Follow manufacturer's recommendations.				
	2. For 1 liter, dissolve the following in 900 ml deionized water:				
	6 g sodium phosphate (monobasic) 372 mg EDTA 50 ml glycerol				
	3. Use NaOH to adjust pH and bring up the volume to 1 liter. Store at $+4^{\circ}$ C.				
	4. Right before use, add the protease inhibitors.				
Vendors for Baffled Flasks	Bellco (1-800-257-7043) has a wide variety of baffled flasks from 50 to 2000 ml. Wheaton (1-609-825-1100) only sells side baffle flasks.				

## Appendix

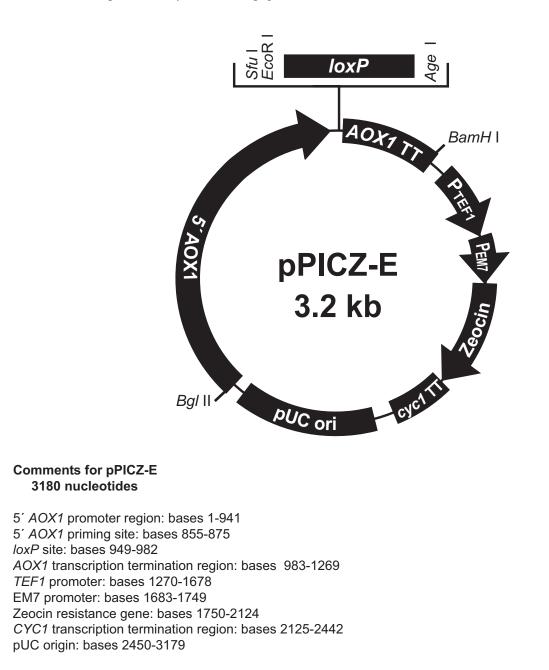
#### Recipes

Low Salt LB Medium with Zeocin <sup>™</sup>	10 g Tryptone <b>5 g NaCl</b> 5 g Yeast Extract
	<ol> <li>Combine the dry reagents above and add deionized, distilled water to 950 ml. Adjust pH to 7.5 with 1N NaOH. Bring the volume up to 1 liter. For plates, add 15 g/L agar before autoclaving.</li> </ol>
	2. Autoclave on liquid cycle at 15 psi and 121°C for 20 minutes.
	3. Allow the medium to cool to at least 55°C before adding the Zeocin <sup>™</sup> to 25 μg/ml final concentration.
	<ol> <li>Store plates at +4°C in the dark. Plates containing Zeocin<sup>™</sup> are stable for up to 2 weeks.</li> </ol>
LB (Luria-Bertani)	Composition:
Medium	1.0% Tryptone 0.5% Yeast Extract 1.0% NaCl pH 7.0
	1. For 1 liter, dissolve 10 g tryptone, 5 g yeast extract, and 10 g NaCl in 950 ml deionized water.
	2. Adjust the pH of the solution to 7.0 with NaOH and bring the volume up to 1 liter.
	3. Autoclave on liquid cycle for 20 minutes at 15 psi. Allow solution to cool to 55°C and add antibiotic if needed.
	4. Store at room temperature or at $+4^{\circ}$ C.
SOB Medium (with Kanamycin)	SOB (per liter) 2% Tryptone 0.5% Yeast Extract 0.05% NaCl 2.5 mM KCl 10 mM MgCl <sub>2</sub>
	1. Dissolve 20 g tryptone, 5 g yeast extract, and 0.5 g NaCl in 950 ml deionized water.
	2. Make a 250 mM KCl solution by dissolving 1.86 g of KCl in 100 ml of deionized water. Add 10 ml of this stock KCl solution to the solution in Step 1.
	3. Adjust pH to 7.5 with 5 M NaOH and add deionized water to 1 liter.
	<ol> <li>Autoclave this solution, cool to ~55°C, and add 10 ml of sterile 1 M MgCl<sub>2</sub>. You may also add kanamycin to 50 μg/ml.</li> </ol>
	5. Store at $+4^{\circ}$ C. Medium is stable for only $\sim 1$ month.
	·

#### Maps of pPICZ-E and pPICZ $\alpha$ -E

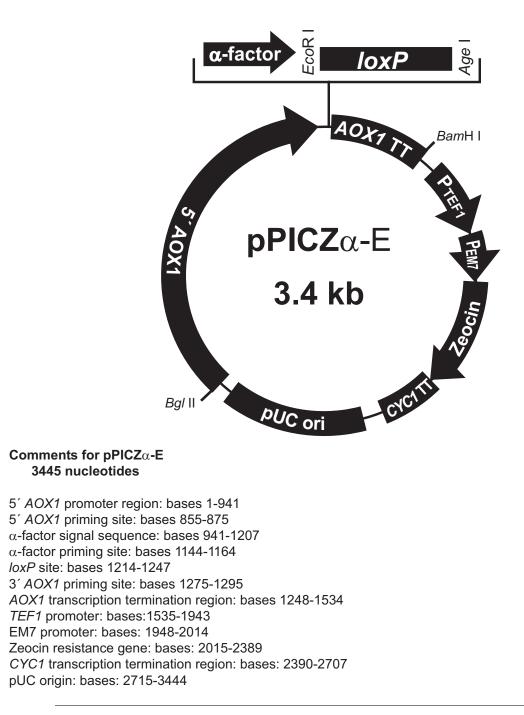
Map of pPICZ-E

The figure below summarizes the features of the pPICZ-E vector. **The complete sequence for pPICZ-E is available for downloading from our Web site (www.invitrogen.com) or from Technical Service (see page 61)**. Details of the sequences surrounding the *loxP* site in pPICZ-E may be found on page 10.



#### Maps of pPICZ-E and pPICZα-E, Continued

Map of pPICZα-E The figure below summarizes the features of the pPICZ $\alpha$ -E vector. The complete sequence for pPICZ $\alpha$ -E is available for downloading from our Web site (www.invitrogen.com) or from Technical Service (see page 61). Details of the sequences surrounding the *loxP* site in pPICZ $\alpha$ -E may be found on page 11.



#### Features of pPICZ-E and pPICZ $\alpha$ -E

#### Features

The important elements of pPICZ-E (3.2 kb) and pPICZ $\alpha$ -E (3.4 kb) are described in the table below. All features have been functionally tested.

Feature	Benefit
5' AOX1	A 942 bp fragment containing the <i>AOX1</i> promoter that allows methanol-inducible, high-level expression in <i>Pichia</i> .
	Targets plasmid integration to the AOX1 locus.
Native Saccharomyces cerevisiae $\alpha$ -factor secretion signal in pPICZ $\alpha$ -E	Allows for efficient secretion of most proteins from <i>Pichia</i> .
<i>loxP</i> site	Allows recombination between the donor vector and pPICZ-E or pPICZ $\alpha$ -E acceptor vector ((Hoess et al., 1982).
<i>AOX1</i> Transcription Termination (TT)	Native transcription termination and polyadenylation signal from AOX1 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 <i>Sh ble</i> gene in <i>Pichia</i> , conferring Zeocin <sup>TM</sup> resistance.
EM7 (synthetic prokaryotic promoter)	Constitutive promoter that drives expression of the <i>Sh ble</i> gene in <i>E. coli</i> , conferring Zeocin <sup>TM</sup> resistance.
Sh ble gene (Streptoalloteichus hindustanus ble gene)	Zeocin <sup>™</sup> resistance gene.
<i>CYC1</i> transcription termination region	3' end of the <i>Saccharomyces cerevisiae CYC1</i> gene that allows efficient 3' mRNA processing of the
(GenBank accession number M34014)	Sh ble gene for increased stability.
pUC origin	Allows high copy replication and maintenance of the plasmid in <i>E. coli</i> .
Sac I	Unique restriction sites that permit linearization of
Pme I	the vectors at the AOX1 locus for efficient
BstX I	integration into the <i>Pichia</i> genome.

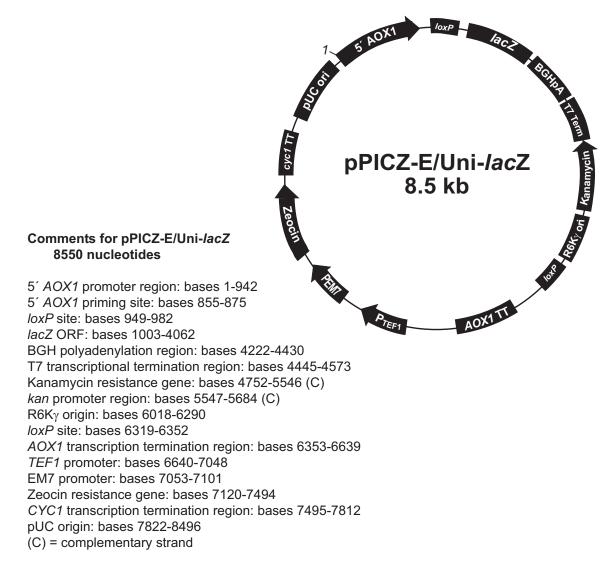
#### Map of pPICZ-E/Uni-IacZ

pPICZ-E/Uni-*lacZ* is a 8.5 kb control vector expressing a β-galactosidase protein (116 kDa). The *lacZ* gene was amplified and cloned into a donor vector, pUni/V5-His-Gene-TOPO<sup>®</sup>. The resulting vector was recombined with pPICZ-E using Cre recombinase to create pPICZ-E/Uni-*lacZ*. **Note**: pUniV5/Gene-TOPO<sup>®</sup> is similar to pUni/V5-His-TOPO<sup>®</sup> TA except that it contains additional DNA between the TOPO<sup>®</sup> Cloning site and the V5 epitope.

#### Map of pPICZ-E/Uni-*lacZ*

Description

The figure below summarizes the features of pPICZ-E/Uni-*lacZ* vector. The complete nucleotide sequence for pPICZ-E/Uni-*lacZ* is available for downloading from our Web site (www.invitrogen.com) or by contacting Technical Service (see page 61).

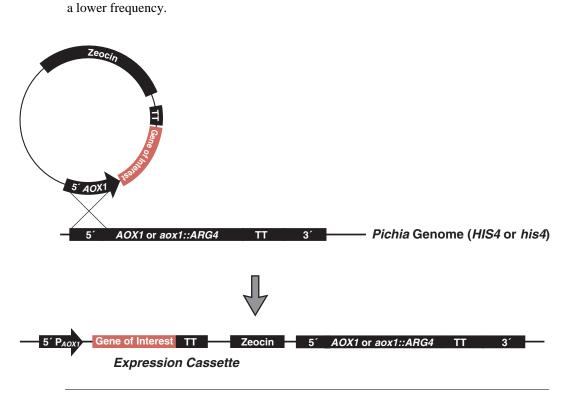


#### Map of pPICZ $\alpha$ -E/Uni-HSA

#### Description pPICZα-E/Uni-HSA is a 7.5 kb control vector expressing a human serum albumin (HSA) protein (69 kDa). The HSA gene was amplified and cloned into a donor vector, pUni/V5-His-Gene-TOPO<sup>®</sup>. The resulting vector was recombined with pPICZα-E using Cre recombinase to create pPICZα-E/Uni-HSA Note: pUniV5/Gene-TOPO<sup>®</sup> is similar to pUni/V5-His-TOPO® TA except that it contains additional DNA between the TOPO<sup>®</sup> Cloning site and the V5 epitope. Map of pPICZ $\alpha$ -The figure below summarizes the features of pPICZα-E/Uni-HSA vector. The E/Uni-HSA complete nucleotide sequence for pPICZα-E/Uni-HSA is available for downloading from our Web Site (www.invitrogen.com) or by contacting Technical Service (see page 61). VSHIST BCIDA pUC ori ni - Kanamycin pPICZα-E/Uni-HSA Comments for pPICZ<sub>α</sub>-E/Uni-HSA 7512 nucleotides 7.5 kb 5' AOX1 promoter region: bases 1-941 5' AOX1 priming site: bases 855-875 $\alpha$ -factor signal sequence: bases 942-1207 $\alpha$ -factor priming site: bases 1144-1164 *loxP* site: bases 1214-1247 HSA gene: bases 1268-3024 AOX1TI V5 epitope: bases 3035-3076 6xHis tag: bases 3086-3103 BGH polyadenylation region: bases 3184-3392 P<sub>TEF</sub>1 T7 transcriptional termination region: bases 3407-3535 Kanamycin resistance gene: bases 3714-4508 (C) kan promoter region: bases 4509-4646 (C) R6Ky origin: bases 4827-5266 *loxP* site: bases 5281-5314 AOX1 transcription termination region: bases 5315-5601 TEF1 promoter: bases 5602-6010 EM7 promoter: bases 6015-6081 Zeocin resistance gene: bases 6082-6456 CYC1 transcription termination region: bases 6457-6774 pUC origin: bases 6785-7511 (C) =complementary strand

#### Recombination and Integration in Pichia

Introduction	Like <i>Saccharomyces cerevisiae</i> , linear DNA can generate stable transformants of <i>Pichia pastoris</i> via homologous recombination between the transforming DNA and regions of homology within the genome (Cregg et al., 1985; Cregg et al., 1989). Such integrants show extreme stability in the absence of selective pressure even when present as multiple copies. Note that single crossover events (insertions) are much more likely to happen than double crossover events (replacements). Multiple insertion events occur spontaneously at about 1-10% of the single insertion events.
Gene Insertion at AOX1 or aox1::ARG4	Gene insertion events at the <i>AOX1</i> (X-33 or GS115) or <i>aox1::ARG4</i> (KM71H) loci arise from a single crossover event between the loci and either of the two <i>AOX1</i> regions on the pPICZ-E or pPICZ $\alpha$ -E vectors: the <i>AOX1</i> promoter or the <i>AOX1</i> transcription termination region (TT). This results in the insertion of one or more copies of the vector upstream or downstream of the <i>AOX1</i> or the <i>aox1::ARG4</i> genes. The phenotype of such a transformant is Mut <sup>+</sup> (X-33 or GS115) or Mut <sup>S</sup> (KM71H). By linearizing the recombinant vector at a restriction enzyme site located in the 5' <i>AOX1</i> regions, Mut <sup>+</sup> or Mut <sup>S</sup> recombinants can be conveniently generated depending on the host strain used.
	The figure below shows the result of an insertion of the plasmid 5' to the intact <i>AOX1</i> locus (Mut <sup>+</sup> ) and the gain of $P_{AOX1}$ , your gene of interest, and the Zeocin <sup>TM</sup> resistance gene. This also occurs with non-linearized plasmid and plasmid that religates, although at

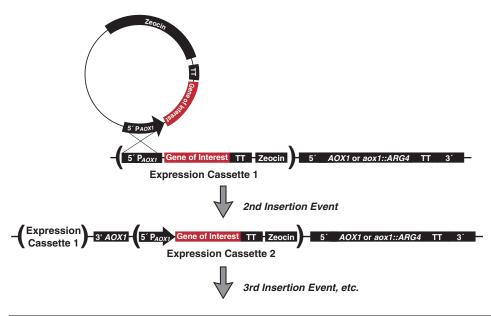


#### Recombination and Integration in Pichia, Continued

#### Multiple Gene Insertion Events

Multiple gene insertion events at a single locus in a cell do occur spontaneously with a low, but detectable frequency-between 1 and 10% of all selected Zeo<sup>R</sup> transformants. Because of the low frequency of multiple gene insertion events, you will need to screen hundreds to thousands of Zeocin<sup>™</sup>-resistant transformants to locate these "jack-pot" clones. We recommend that you use electroporation to generate Zeo<sup>R</sup> transformants for screening.

Multi-copy events can occur as gene insertions either at the *AOX1* or the *aox1::ARG4* loci. This results in a Mut<sup>+</sup> phenotype in X-33 or GS115 and a Mut<sup>S</sup> phenotype in KM71H. Multiple gene insertion events can be detected by quantitative dot blot analysis, Southern blot analysis, and differential hybridization. See page 18 for a protocol to screen for multiple inserts.



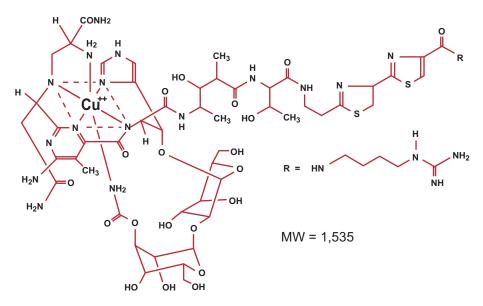
Zeocin<sup>™</sup>

Zeocin<sup>™</sup> is a member of the bleomycin/phleomycin family of antibiotics isolated from *Streptomyces*. Antibiotics in this family are broad spectrum antibiotics that act as strong anti-bacterial and anti-tumor drugs. They show strong toxicity against bacteria, fungi (including yeast), plants, and mammalian cells (Baron et al., 1992; Drocourt et al., 1990; Mulsant et al., 1988; Perez et al., 1989)

The Zeocin<sup>TM</sup> resistance protein has been isolated and characterized (Calmels et al., 1991; Drocourt et al., 1990). This protein, the product of the *Sh ble* gene (*Streptoalloteichus hindustanus* bleomycin gene), is a 13.7 kDa protein that binds Zeocin<sup>TM</sup> and inhibits its DNA strand cleavage activity. Expression of this protein in eukaryotic and prokaryotic hosts confers resistance to Zeocin<sup>TM</sup>.

#### Molecular Weight, Formula, and Structure

The formula for Zeocin<sup>TM</sup> is  $C_{60}H_{89}N_{21}O_{21}S_3$  and the molecular weight is 1,535. The diagram below shows the structure of Zeocin<sup>TM</sup>.



#### Applications of Zeocin<sup>™</sup>

Zeocin<sup>TM</sup> 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<sup>TM</sup> for selection in *Pichia* and *E. coli* are listed below:

Organism	Zeocin <sup>™</sup> Concentration and Selective Medium
E. coli	25-50 μg/ml in <b>Low Salt LB</b> medium <sup>*</sup> (see page 45 for a recipe)
Pichia	100-1000 $\mu$ 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<sup>™</sup>, Continued

# Handling Zeocin<sup>™</sup> High salt and acidity or basicity inactivate Zeocin<sup>™</sup>; therefore, we recommend that you reduce the salt in bacterial medium and adjust the pH to 7.5 to keep the drug active (see Low Salt LB Medium, page 45). Note that the salt concentration and pH do not need to be adjusted when preparing tissue culture medium containing Zeocin<sup>™</sup>. Store Zeocin<sup>™</sup> at -20°C and thaw on ice before use. Zeocin<sup>™</sup> is light sensitive. Store drug, plates, and medium containing drug in the dark.

- Wear gloves, a laboratory coat, and safety glasses or goggles when handling solutions containing Zeocin<sup>™</sup>.
- $\operatorname{Zeocin}^{\mathbb{M}}$  is toxic. Do not ingest or inhale solutions containing the drug.
- Store tissue culture medium containing Zeocin<sup>™</sup> at +4°C in the dark. Medium containing Zeocin<sup>™</sup> is stable for 1-2 months.

## Direct PCR Screening of Pichia Clones

Introduction	inse con	imple protocol has been ertion of your gene by I abination of enzyme, fr ectly as a PCR template	PCR (Linder et al., reezing, and heatin	1996) Briefly, the	
Before Starting	You	ı will need the followir	ng reagents and eq	uipment on hand:	
J	• • • •	A culture or single co 1.5 ml microcentrifug 5 U/µl solution of Lyt 30°C water bath or he Liquid nitrogen Reagents for PCR	e tube icase (Sigma)	ansformant	
Procedure	1.	Place 10 $\mu$ l of a <i>Pichia pastoris</i> culture into a 1.5 ml microcentrifuge tube. For relatively dense cultures, dilute 1 $\mu$ l of the culture into 9 $\mu$ l water. Alternatively, pick a single colony and resuspend in 10 $\mu$ l of water.			
	2.	Add 5 $\mu$ l of a 5 U/ $\mu$ l	solution of lyticas	e and incubate at 3	30°C for 10 minutes.
	3.	Freeze the sample at 1 minute.	-80°C for 10 minu	ites or immerse in	liquid nitrogen for
	4.	Set up a 50 µl PCR f	or a hot start:		
		10X Reaction Buffer		5 µl	l
		25 mM MgCl <sub>2</sub>		5 µl	l
		25 mM dNTPs		1 μl	l
		5' AOX1 primer (10)	pmol/µl)	1 μl	l
		3' AOX1 primer (10)	pmol/µl)	1 µl	l
		Sterile water		27 μl	l
		Cell lysate		5 µl	<u>l</u>
		Total Volume		45 μl	l
	5.	Mix the solution and	overlay with 20 $\mu$	l of mineral oil.	
	6.	Place the solution in	the thermocyler ar	nd incubate at 95°	C for 5 minutes.
	7.	Add 5 µl of a 0.16 U	/µl solution of <i>Taq</i>	polymerase (0.8	units).
	8.	Cycle 30 times using the following parameters:			
		Step	Temperature	Time	]
		Denaturation	95°C	1 minute	
		Annealing	54°C	1 minute	
		Extension	72°C	1 minute	
		Include a final extens	sion of 7 minutes	1	-

9. Analyze a 10 µl aliquot by agarose gel electrophoresis.

#### Total DNA Isolation from Pichia

<ul> <li>Solutions</li> <li>You will need to make the following solutions. There is not enough of some of the reagents in the kit to perform this experiment</li> <li>Minimal Medium (MD, MGYH)</li> <li>Sterile water</li> <li>SCED (1 M sorbitol, 10 mM sodium citrate, pH 7.5, 10 mM EDTA, 10 mM D</li> <li>Zymolyase, 3 mg/ml stock solution in water (Seikagaku America, Inc., 1-800-4512)</li> <li>1% SDS in water</li> <li>5 M potassium acetate, pH 8.9</li> <li>TE buffer, pH 7.4 (10 mM Tris-HCl, pH 7.4, 1 mM EDTA, pH 8.0)</li> <li>7.5 M ammonium acetate, pH 7.5</li> <li>Phenol:chloroform (1:1 v/v)</li> </ul>	e			
<ul> <li>Sterile water</li> <li>SCED (1 M sorbitol, 10 mM sodium citrate, pH 7.5, 10 mM EDTA, 10 mM D</li> <li>Zymolyase, 3 mg/ml stock solution in water (Seikagaku America, Inc., 1-800-4512)</li> <li>1% SDS in water</li> <li>5 M potassium acetate, pH 8.9</li> <li>TE buffer, pH 7.4 (10 mM Tris-HCl, pH 7.4, 1 mM EDTA, pH 8.0)</li> <li>7.5 M ammonium acetate, pH 7.5</li> </ul>	You will need to make the following solutions. There is not enough of some of these reagents in the kit to perform this experiment			
<ul> <li>SCED (1 M sorbitol, 10 mM sodium citrate, pH 7.5, 10 mM EDTA, 10 mM D</li> <li>Zymolyase, 3 mg/ml stock solution in water (Seikagaku America, Inc., 1-800-4512)</li> <li>1% SDS in water</li> <li>5 M potassium acetate, pH 8.9</li> <li>TE buffer, pH 7.4 (10 mM Tris-HCl, pH 7.4, 1 mM EDTA, pH 8.0)</li> <li>7.5 M ammonium acetate, pH 7.5</li> </ul>				
<ul> <li>Zymolyase, 3 mg/ml stock solution in water (Seikagaku America, Inc., 1-800-4512)</li> <li>1% SDS in water</li> <li>5 M potassium acetate, pH 8.9</li> <li>TE buffer, pH 7.4 (10 mM Tris-HCl, pH 7.4, 1 mM EDTA, pH 8.0)</li> <li>7.5 M ammonium acetate, pH 7.5</li> </ul>				
<ul> <li>4512)</li> <li>1% SDS in water</li> <li>5 M potassium acetate, pH 8.9</li> <li>TE buffer, pH 7.4 (10 mM Tris-HCl, pH 7.4, 1 mM EDTA, pH 8.0)</li> <li>7.5 M ammonium acetate, pH 7.5</li> </ul>	ΓT)			
<ul> <li>5 M potassium acetate, pH 8.9</li> <li>TE buffer, pH 7.4 (10 mM Tris-HCl, pH 7.4, 1 mM EDTA, pH 8.0)</li> <li>7.5 M ammonium acetate, pH 7.5</li> </ul>	237-			
<ul> <li>TE buffer, pH 7.4 (10 mM Tris-HCl, pH 7.4, 1 mM EDTA, pH 8.0)</li> <li>7.5 M ammonium acetate, pH 7.5</li> </ul>				
• 7.5 M ammonium acetate, pH 7.5				
-				
• Phenol:chloroform (1:1 v/v)				
<ul> <li>Preparation</li> <li>1. Grow at 30°C the recombinant strain and the parent strain to an OD<sub>600</sub> of 5-1 ml of minimal media such as MD or MDH. Note: <i>his4</i> strains require histidin growth.</li> </ul>				
2. Collect the cells by centrifugation at 1500 x g for 5-10 minutes at room temperature.				
3. Wash the cells with 10 ml sterile water by centrifugation as in Step 2.				
<b>Spheroplasting</b> 1. Resuspend the cells in 2 ml of SCED buffer, pH 7.5. Make this solution fresh				
and Lysis 2. Add 0.1-0.3 mg of Zymolyase (mix well before adding to the cells). Incubate 37°C for 50 minutes to achieve < 80% spheroplasting.	at			
3. Add 2 ml of 1% SDS, mix <b>gently</b> and set on ice (0 to $+4^{\circ}$ C) for 5 minutes.				
4. Add 1.5 ml of 5 M potassium acetate, pH 8.9, and mix gently.				
5. Centrifuge at 10,000 x g for 5-10 minutes at $+4^{\circ}$ C and save the supernatant.				

## Total DNA Isolation from Pichia, Continued

DNA Precipitation	1.	Transfer the supernatant from Step 5 above and add 2 volumes of ethanol to the supernatant. Incubate at room temperature for 15 minutes.
	2.	Centrifuge at 10,000 x g for 20 minutes at +4°C.
	3.	Resuspend the pellet <b>gently</b> in 0.7 ml of TE buffer, pH 7.4 and transfer to a microcentrifuge tube.
	4.	<b>Gently</b> extract with an equal volume of phenol:chloroform $(1:1 v/v)$ followed by an equal volume of chloroform:isoamyl alcohol (24:1). Split the aqueous layer into two microcentrifuge tubes.
	5.	Add 1/2 volume of 7.5 M ammonium acetate, pH 7.5, and 2 volumes of ethanol to each tube. Place on dry ice for 10 minutes or at -20°C for 60 minutes.
	6.	Centrifuge at 10,000 x g for 20 minutes at $+4^{\circ}$ C and wash the pellets once with 1 ml of 70% ethanol. Air dry the pellets and resuspend each in 50 µl of TE buffer, pH 7.5. Determine the concentration of the DNA sample. The two samples can be stored separately or combined and stored at $-20^{\circ}$ C until ready for use.

## **Determination of Copy Number of Multiple Integrants**

Introduction	You may wish to determine the actual number of gene copies in your <i>Pichia</i> recombinant using either quantitative dot blots or Southern hybridization. ((Brierley et al., 1994; Clare et al., 1991a; Romanos et al., 1991; Scorer et al., 1993; Scorer et al., 1994)). This requires isolation of genomic DNA from <i>Pichia</i> recombinants transformed with the parent vector (0 copies of your gene), pPICZ-E or pPICZ $\alpha$ -E containing 1 copy of your gene (single copy control), and the <i>Pichia</i> recombinants containing multiple copies of your gene. Use the protocol detailed on the previous page to isolate genomic DNA or you may use any method of your choice.
Quantitative Dot Blot Solutions	<ul> <li>You will need the following solutions, 10-15 ml of each for each dot blot.</li> <li>50 mM EDTA, 2.5% β-mercaptoethanol pH 9</li> <li>1 mg/ml Zymolyase 100T in water (Seikagaku America, Inc., 1-800-237-4512)</li> <li>0.1N NaOH, 1.5 M NaCl</li> <li>2 x SSC</li> <li>3MM paper</li> </ul>
Quantitative Dot Blot Procedure	<ul> <li>The following protocol is a summary of a rapid DNA dot blot technique to detect multiple integrants (Romanos, <i>et al.</i>, 1991). It is very important to spot equivalent numbers of cells onto filters in order to quantify copy number. Genomic DNA may also be isolated and spotted directly onto nitrocellulose or nylon, fixed, and analyzed.</li> <li>Grow Mut<sup>+</sup> or Mut<sup>§</sup> transformants in individual wells of a 96-well microtiter plate in 200 µl of YPD broth at 30°C until all wells have approximately the same density. Alternatively, individual transformants may be grown in culture tubes and the absorbance at 600 nm normalized with the addition of medium.</li> <li>Filter 50 µl of each sample onto a nitrocellulose or nylon filter placed into a dot (slot) blot apparatus using multi-channel pipettor. Air dry filters.</li> <li>Lyse the cells on the filter with four solutions as follows: soak two sheets of 3 MM paper in a tray with 10-15 ml of 50 mM EDTA, 2.5% β-mercaptoethanol pH 9. After the paper is uniformly soaked, place the nitrocellulose filter face down on the treated 3MM paper. Incubate for 15 minutes at room temperature.</li> <li>Remove the nitrocellulose filter from the 3MM paper and replace the 3MM paper with two new sheets. Soak with 10-15 ml of 1 mg/ml Zymolyase 100T as described in Step 3. Place the nitrocellulose filter face down on the 3MM paper and incubate for 4 hours at 37°C.</li> <li>Remove the nitrocellulose filter from the paper and replace the paper with two new sheets. Soak with 10-15 ml of 0.1 N NaOH, 1.5 M NaCl. Place the nitrocellulose filter face down on the 3MM paper and incubate for 5 minutes at room temperature.</li> <li>Remove the nitrocellulose filter and replace with two new 3MM sheets. Soak with 10-15 ml of 2 x SSC. Place the nitrocellulose filter face down on the 3MM paper and incubate for 5 minutes at room temperature. Repeat.</li> <li>Bake the nitrocellulose filters at 80°C or UV-crosslink DNA to nylon. The filters may be probed with a norradioactive-labeled or random-primed, <sup></sup></li></ul>

#### Determination of Copy Number of Multiple Integrants, Continued

Southern Blot Analysis	For a detailed description of this technique as applied to <i>Pichia pastoris</i> , see (Clare, <i>et al.</i> , 1991a). It is very important to digest your DNA with the right restriction enzyme(s) to generate a blot of digested and gel-separated genomic DNA. We recommend that you use a restriction enzyme that cuts <b>outside</b> of the expression cassette. This will generate a restriction fragment whose size reflects the number of multimers. For example, if a 1.2 kb fragment is cloned into pichia vector and transformed into <i>Pichia</i> , DNA from this recombinant can be digested with <i>Nhe</i> I (which does not cut within the vector or the insert). The size of the fragment will increase in additions of 4.7 kb over the size of the fragment containing 1 copy of the gene.
	Alternatively, digestion of DNA from <i>Pichia</i> recombinants containing multiple copies will produce a band that will vary in intensity depending on the number of copies of your gene. It is very important to include a control to show the intensity of a single copy gene. The band intensities can be relatively quantified using densitometry to estimate gene dosage.
Controls	It is very important to include DNA from the host strain alone (X-33, GS115 or KM71H), the host strain transformed with the parent vector, and the host strain transformed with a vector containing one copy of your gene.
General Guidelines	• Use standard procedures and solutions for Southern blotting as outlined in <i>Molecular Cloning: A Laboratory Manual</i> (Sambrook, <i>et al.</i> , 1989), pages 9.31-9.58.
	• Isolate genomic DNA and quantify using fluorometry. Be sure to eliminate RNA. It is very important to load the same amount of DNA into each lane in order to accurately determine copy number if you are using relative intensities.
	Probe your Southern blot with a fragment complementary to your gene

#### Procedure for Total RNA Isolation from Pichia

Introduction	This protocol is designed to isolate 60-300 $\mu$ g total RNA (Schmitt et al., 1990) from <i>Pichia</i> which is suitable for mRNA isolation using Invitrogen's FastTrack <sup>®</sup> 2.0 or Micro FastTrack <sup>M</sup> 2.0 mRNA Isolation Kit. If you wish to use another protocol, you should scale-up the reaction to yield about 2 mg of total RNA per time point. The mRNA is for northern blot analysis of <i>Pichia</i> recombinants to determine if the gene of interest is being induced and transcribed. RNA isolation should be done from induced cultures using an uninduced culture as a negative control.			
Solutions	You will need the following solutions. Remember to use DEPC-treated water and to us equipment free of RNase.	e		
	MGYH or BMGY mediumDEPC-treated waterAE buffer (50 mM sodium acetate, pH 5.3, 1 mM EDTA)Buffered phenol10% SDS in DEPC treated waterPhenol:chloroform (1:1)Chloroform:isoamyl alcohol (24:1)65°C water bath3 M sodium acetate, pH 5.3Solution acetate, pH 5.3	)		
Growth of Cells	<ol> <li>Grow up two cultures (100-200 ml in MGYH or BMGY), but induce only one of them. Use the same protocol for induction that you used in the Expression section.</li> </ol>			
	2. Take 10 ml time points at 1, 2, 3, 4, and 6 days.			
	3. Harvest the cells from each time point by centrifugation at 1500 x g for 10 minute at room temperature.	es		
	4. Resuspend cell pellet in 400 $\mu$ l AE buffer and transfer to a microcentrifuge tube.			
Lysis of Cells	1. Add 40 $\mu$ l 10% SDS and vortex for ~ 20 seconds.			
	<ol> <li>Add an equal volume (450-500 μl) of buffer saturated phenol and vortex for ~20 seconds.</li> </ol>			
	3. Incubate at 65°C for 4 minutes.			
	4. Incubate in a dry ice/ethanol bath until crystals show (~1 minute). Centrifuge at maximum speed for 2 minutes at +4°C.			
	<ol> <li>Transfer aqueous phase to new centrifuge tube and add an equal volume of phenol/chloroform and vortex for ~20 seconds. Centrifuge at maximum speed for minutes at +4°C.</li> </ol>	2		
	5. Remove upper phase to a new tube and add 40 $\mu$ l of 3 M sodium acetate, pH 5.3 and 2.5 volumes of 100% ethanol (-20°C). Centrifuge at maximum speed for 15 minutes at +4°C. Remove ethanol.			
	<ol> <li>Wash pellet with 80% ethanol and air dry briefly. Resuspend total RNA in 20 μl DEPC-treated water and store at -80°C. Yield is 60-300 μg total RNA.</li> </ol>			
mRNA Isolation and Northern Analysis	See (Ausubel <i>et al.</i> , 1994) for a protocol for mRNA isolation and northern analysis. The FastTrack <sup>®</sup> 2.0 mRNA Kit (Catalog no K1593-02; 6 reactions) is designed to isolate mRNA from 0.2 to 1 mg total RNA. The Micro-FastTrack <sup>™</sup> 2.0 Kit (Catalog no. K152) (2; 20 reactions) is designed to isolate mRNA from ~100 µg total RNA. You will need approximately 1-5 µg mRNA per time point.	0-		

#### **Technical Service**

Visit the Invitrogen Web site at <u>www.invitrogen.com</u> for:

## Web Resources

- Technical resources, including manuals, vector maps and sequences, application notes, MSDSs, FAQs, formulations, citations, handbooks, etc.
- Complete technical service contact information
- Access to the Invitrogen Online Catalog
- Additional product information and special offers

#### Contact Us

#### Corporate Headquarters: Invitrogen Corporation 1600 Faraday Avenue Carlsbad, CA 92008 USA Tel: 1 760 603 7200 Tel (Toll Free): 1 800 955 6288 Fax: 1 760 602 6500 E-mail: tech service@invitrogen.com

#### Japanese Headquarters: Invitrogen Japan LOOP-X Bldg. 6F 3-9-15, Kaigan Minato-ku, Tokyo 108-0022 Tel: 81 3 5730 6509 Fax: 81 3 5730 6519 E-mail: jpinfo@invitrogen.com

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

#### Material Data Safety Sheets (MSDSs)

Limited

Warranty

MSDSs are available on our Web site at <u>www.invitrogen.com</u>. On the home page, click on **Technical Resources** and follow instructions on the page to download the MSDS for your product.

For more information or technical assistance, call, write, fax, or email. Additional

international offices are listed on our Web page (www.invitrogen.com).

Invitrogen is committed to providing our customers with high-quality goods and services. Our goal is to ensure that every customer is 100% satisfied with our products and our service. If you should have any questions or concerns about an Invitrogen product or service, contact our Technical Service Representatives.

Invitrogen warrants that all of its products will 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 limits Invitrogen Corporation's liability</u> <u>only to the cost 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. Invitrogen reserves the right to select the method(s) used to analyze a product unless Invitrogen agrees to a specified method in writing prior to acceptance of the order.

Invitrogen makes every effort to ensure the accuracy of its publications, but realizes that the occasional typographical or other error is inevitable. Therefore Invitrogen makes no warranty of any kind regarding the contents of any publications or documentation. If you discover an error in any of our publications, please report it to our Technical Service Representatives.

Invitrogen assumes no responsibility or liability for any special, incidental, indirect or consequential loss or damage whatsoever. The above limited warranty is sole and exclusive. No other warranty is made, whether expressed or implied, including any warranty of merchantability or fitness for a particular purpose.

#### References

- Abremski, K. & Hoess, R. (1984). Bacteriophage P1 Site-Specific Recombination. Purification and Properties of the Cre Recombinase Protein. J. Biol. Chem. 259, 1509-1514.
- Abremski, K., Hoess, R. & Sternberg, N. (1983). Studies on the Properties of P1 Site-Specific Recombination: Evidence for Topologically Unlinked Products Following Recombination. *Cell* **32**, 1301-1311.
- Abremski, K. E. & Hoess, R. H. (1992). Evidence for a Second Conserved Arginine Residue in the Integrase Family of Recombination Proteins. *Protein Eng.* **5**, 87-91.
- Argos, P., Landy, A., Abremski, K., Egan, J. B., Haggard-Ljungquist, E., Hoess, R. H., Kahn, M. L., Kalionis, B., Narayana, S. V. L., Pierson III, L. S., Sternberg, N. & Leong, J. M. (1986). The Integrase Family of Site-Specific Recombinases: Regional Similarities and Global Diversity. *EMBO J.* 5, 433-440.
- Ausubel, F. M., Brent, R., Kingston, R. E., Moore, D. D., Seidman, J. G., Smith, J. A. & Struhl, K. (1994). *Current Protocols in Molecular Biology*, Greene Publishing Associates and Wiley-Interscience, New York.
- Baron, M., Reynes, J. P., Stassi, D. & Tiraby, G. (1992). A Selectable Bifunctional β-Galactosidase::Phleomycinresistance Fusion Protein as a Potential Marker for Eukaryotic Cells. *Gene* **114**, 239-243.
- Barr, K. A., Hopkins, S. A. & Sreekrishna, K. (1992). Protocol for Efficient Secretion of HSA Developed from *Pichia pastoris. Pharm. Eng.* 12(2), 48-51.
- Bollag, D. M. & Edelstein, S. J. (1991). Protein Methods, Wiley-Liss, New York.
- Brierley, R. A., Davis, G. R. & Holtz, G. C. (1994). Production of Insulin-Like Growth Factor-1 in Methylotrophic Yeast Cells. *United States Patent* **5**,324,639.
- Buckholz, R. G. & Gleeson, M. A. G. (1991). Yeast Systems for the Commercial Production of Heterologous Protein. *Bio/Technology* **9**, 1067-1072.
- Calmels, T., Parriche, M., Burand, H. & Tiraby, G. (1991). High Efficiency Transformation of *Tolypocladium* geodes Conidiospores to Phleomycin Resistance. *Curr. Genet.* **20**, 309-314.
- Clare, J. J., Rayment, F. B., Ballantine, S. P., Sreekrishna, K. & Romanos, M. A. (1991a). High-level Expression of Tetanus Toxin Fragment c in *Pichia pastoris* Strains Containing Multiple Tandem Integrations of the Gene. *Bio/Technology* 9, 455-460.
- Clare, J. J., Romanos, M. A., Rayment, F. B., Rowedder, J. E., Smith, M. A., Payne, M. M., Sreekrishna, K. & Henwood, C. A. (1991b). Production of Epidermal Growth Factor in Yeast: High-Level Secretion Using *Pichia pastoris* Strains Containing Multiple Gene Copies. *Gene* **105**, 205-212.
- Cregg, J. M., Barringer, K. J. & Hessler, A. Y. (1985). Pichia pastoris as a Host System for Transformations. Mol. Cell. Biol. 5, 3376-3385.
- Cregg, J. M. & Higgins, D. R. (1995). Production of Foreign Proteins in the Yeast *Pichia pastoris*. *Canadian J. Botany Supp.* **73**, 5981-5987.
- Cregg, J. M., Madden, K. R., Barringer, K. J., Thill, G. & Stillman, C. A. (1989). Functional Characterization of the Two Alcohol Oxidase Genes from the Yeast, *Pichia pastoris. Mol. Cell. Biol.* **9**, 1316-1323.
- Cregg, J. M., Vedvick, T. S. & Raschke, W. C. (1993). Recent Advances in the Expression of Foreign Genes in *Pichia pastoris. Bio/Technology* **11**, 905-910.
- Deutscher, M. P., Ed. (1990). Guide to Protein Purification. Vol. 182. Methods in Enzymology. Edited by Abelson, J. N. & Simon, M. I. San Diego, CA.: Academic Press.
- Drocourt, D., Calmels, T. P. G., Reynes, J. P., Baron, M. & 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. & Gingeras, T. R. (1985). Isolation of Alcohol Oxidase and Two other Methanol Regulatable Genes from the Yeast, *Pichia pastoris. Mol. Cell. Biol.* 5(5), 1111-1121.
- Grinna, L. S. & Tschopp, J. F. (1989). Size Distribution and General Structural Features of N-Linked Oligosaccharides from the Methylotrophic Yeast, *Pichia pastoris*. *Yeast* **5**, 107-115.
- Guthrie, C. & Fink, G. R., Eds. (1991). Guide to Yeast Genetics and Molecular Biology. Vol. 194. Methods in Enzymology. Edited by Abelson, J. N. & Simon, M. I. San Diego, CA: Academic Press.
- Higgins, D. R. & Cregg, J. M., Eds. (1998). *Pichia* Protocols. Vol. 103. Methods in Molecular Biology. Edited by Walker, J. M. Totowa, NJ: Humana Press.
- Hoess, R. H., Ziese, M. & Sternberg, N. (1982). P1 Site-Specific Recombination: Nucleotide Sequence of the Recombining Sites. *Proc. Natl. Acad. Sci USA* 79, 3398-3402.
- Holm, C., Meeks-Wagner, D. W., Fangman, W. L. & Botstein, D. (1986). A Rapid, Efficient Method for Isolating DNA from Yeast. *Gene* 42, 169-173.
- Koutz, P. J., Davis, G. R., Stillman, C., Barringer, K., Cregg, J. M. & Thill, G. (1989). Structural Comparison of the *Pichia pastoris* Alcohol Oxidase Genes. *Yeast* 5, 167-177.
- Linder, S., Schliwa, M. & Kube-Granderath, E. (1996). Direct PCR Screening of *Pichia pastoris* Clones. *BioTechniques* **20**, 980-982.
- Lindner, P., Bauer, K., Krebber, A., Nieba, L., Kremmer, E., Krebber, C., Honegger, A., Klinger, B., Mocikat, R. & Pluckthun, A. (1997). Specific Detection of His-tagged Proteins With Recombinant Anti-His Tag scFv-Phosphatase or scFv-Phage Fusions. *BioTechniques* 22, 140-149.
- Liu, Q., Li, M. Z., Leibham, D., Cortez, D. & Elledge, S. (1998). The Univector Plasmid-Fusion System, a Method for Rapid construction of Recombinant DNA Without Restriction Enzymes. *Current Biology* 8, 1300-1309.
- Liu, Q., Li, M. Z., Liu, D. & Elledge, S. J. (1999). Rapid Construction of Recombinant DNA by the Univector Plasmid-Fusion System. *Methods in Enzymology, in press.*
- Mulsant, P., Tiraby, G., Kallerhoff, J. & Perret, J. (1988). Phleomycin Resistance as a Dominant Selectable Marker in CHO Cells. *Somat. Cell Mol. Genet.* **14**, 243-252.
- Nico-Farber, K., Harder, W., Ab, G. & Veenhuis, M. (1995). Review: Methylotrophic Yeasts as Factories for the Production of Foreign Proteins. *Yeast* **11**, 1331-1344.
- Perez, P., Tiraby, G., Kallerhoff, J. & Perret, J. (1989). Phleomycin Resistance as a Dominant Selectable Marker for Plant Cell Transformation. *Plant Mol. Biol.* **13**, 365-373.
- Romanos, M. A., Clare, J. J., Beesley, K. M., Rayment, F. B., Ballantine, S. P., Makoff, A. J., Dougan, G., Fairweather, N. F. & Charles, I. G. (1991). Recombinant *Bordetella pertussis* Pertactin p69 from the Yeast *Pichia pastoris* High Level Production and Immunological Properties. *Vaccine* 9, 901-906.
- Romanos, M. A., Scorer, C. A. & Clare, J. J. (1992). Foreign Gene Expression in Yeast: A Review. Yeast 8, 423-488.
- Sambrook, J., Fritsch, E. F. & Maniatis, T. (1989). Molecular Cloning: A Laboratory Manual. Second edit, Cold Spring Harbor Laboratory Press, Plainview, New York.
- Schmitt, M. E., Brown, T. A. & Trumpower, B. L. (1990). A Rapid and Simple Method for Preparation of RNA from Saccharomyces cerevisiae. Nucleic Acids Res. 18(10), 3091.
- Scorer, C. A., Buckholz, R. G., Clare, J. J. & Romanos, M. A. (1993). The Intracellular Production and Secretion of HIV-1 Envelope Protein in the Methylotrophic Yeast *Pichia pastoris*. *Gene* 136, 111-119.

#### **References**, Continued

- Scorer, C. A., Clare, J. J., McCombie, W. R., Romanos, M. A. & Sreekrishna, K. (1994). Rapid Selection Using G418 of High Copy Number Transformants of *Pichia pastoris* for High-level Foreign Gene Expression. *Bio/Technology* 12, 181-184.
- Southern, J. A., Young, D. F., Heaney, F., Baumgartner, W. & 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.
- Sreekrishna, K., Potenz, R. H. B., Cruze, J. A., McCombie, W. R., Parker, K. A., Nelles, L., Mazzaferro, P. K., Holden, K. A., Harrison, R. G., Wood, P. J., Phelps, D. A., Hubbard, C. E. & Fuke, M. (1988). High Level Expression of Heterologous Proteins in Methylotrophic Yeast *Pichia pastoris*. J. Basic Microbiol. 28(4), 265-278.
- Sternberg, N., Hamilton, D., Austin, S., Yarmolinsky, M. & Hoess, R. (1981). Site-Specific Recombination and its Role in the Life Cycle of P1. CSH Symp. Quant. Biol. 45, 297-309.
- Strathern, J. N. & Higgins, D. R. (1991). Recovery of Plasmids from Yeast into *Escherichia coli*: Shuttle Vectors. In *Guide to Yeast Genetics and Molecular Biology* (Guthrie, C. & Fink, G. R., eds.), Vol. 194. Academic Press, San Diego, CA.
- Tschopp, J. F., Brust, P. F., Cregg, J. M., Stillman, C. & Gingeras, T. R. (1987a). Expression of the *lacZ* Gene from Two Methanol Regulated Promoters in *Pichia pastoris*. *Nucleic Acids Res.* **15**(9), 3859-3876.
- Tschopp, J. F., Sverlow, G., Kosson, R., Craig, W. & Grinna, L. (1987b). High Level Secretion of Glycosylated Invertase in the Methylotrophic Yeast *Pichia pastoris*. *Bio/Technology* **5**, 1305-1308.
- Wegner, G. H. (1990). Emerging Applications of the Methylotrophic Yeasts. *FEMS Microbiology Reviews* 87, 279-284.

©2000-2006 Invitrogen Corporation. All rights reserved.

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

## invitrogen®

Corporate Headquarters Invitrogen Corporation 1600 Faraday Avenue Carlsbad, CA 92008 T: 1 760 603 7200 F: 1 760 602 6500 E: tech.service@invitrogen.com

For country-specific contact information visit our web site at **www.invitrogen.com**