# Pichia Spheroplast Kit

A Manual for the Preparation and Transformation of *Pichia pastoris* Spheroplasts

Catalog no. K1720-01

Version B

030502 25-0055



www.invitrogen.com tech\_service@invitrogen.com

### 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.

#### **Product User Registration Card**

Please complete and return the enclosed Product User Registration Card for each *Pichia* Expression Kit that you purchase. This will serve as a record of your purchase and registration and will allow Invitrogen to provide you with technical support and manual updates. It will also allow Invitrogen to update you on future developments of and improvements to the *Pichia* Expression Kit. The agreement outlined above becomes effective upon our receipt of your User Registration Card or 10 days following the sale of the *Pichia* Expression Kit to you. Use of the kit at any time results in immediate obligation to the terms and conditions stated in this Agreement.

#### **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.

United States Headquarters:	Japanese Headquarters:	European Headquarters:
Invitrogen Corporation	Invitrogen Japan K.K.	Invitrogen Ltd
1600 Faraday Avenue	Nihonbashi Hama-Cho Park Bldg. 4F	3 Fountain Drive
Carlsbad, CA 92008 USA	2-35-4, Hama-Cho, Nihonbashi	Inchinnan Business Park
Tel: 1 760 603 7200	Tel: 81 3 3663 7972	Paisley PA4 9RF, UK
Tel (Toll Free): 1 800 955 6288	Fax: 81 3 3663 8242	Tel (Free Phone Orders): 0800 269 210
Fax: 1 760 602 6500	E-mail: jpinfo@invitrogen.com	Tel (General Enquiries): 0800 5345 5345
E-mail: tech_service@invitrogen.com		Fax: +44 (0) 141 814 6287
		E-mail: eurotech@invitrogen.com

### **Table of Contents**

Table of Contents	v
Important Information	vi
Methods	1
Overview	1
Preparing the Transforming DNA and Pichia Cells	2
Preparing Spheroplasts	4
Transformation	6
Appendix	7
Recipes	7
References	10

### **Important Information**

Kit ContentsThe Pichia Sphreoplast Kit contains 2 boxes. Store each box as detailed below:Store Box 1 at +4°C.Store Box 2 at -20°C.

#### **Box 1 Reagents** Store Box 1 at $+4^{\circ}$ C.

Item	Components	Amount
SOS media	1 M Sorbitol, 0.3X YPD, 10 mM CaCl <sub>2</sub>	20 ml
Sterile Water		2 x 125 ml
SE	1 M Sorbitol; 25 mM EDTA, pH 8.0	2 x 125 ml
SCE	1 M Sorbitol; 1 mM EDTA; 10 mM sodium citrate buffer, pH 5.8	2 x 125 ml
1 M Sorbitol		2 x 125 ml
CaS	1 M Sorbitol; 10 mM Tris-HCl, pH 7.5; 10 mM CaCl <sub>2</sub>	2 x 60 ml
40% PEG	40% (w/v) PEG 3350 (Fisher Scientific grade) in water	25 ml
CaT	20 mM Tris-HCl, pH 7.5; 20 mM CaCl <sub>2</sub>	25 ml

#### **Box 2 Reagents** Store Box 2 at -20°C.

Item	Components	Amount
Zymolyase	3 mg/ml Zymolyase in water	10 x 20 μl
1 M DTT	1 M dithiothreitol in water	10 x 1 ml

#### Materials Supplied by User

**Materials Supplied** You will need the following reagents and equipment.

- 30°C rotary shaking incubator
  - Water baths capable of 37°C and 45°C
  - Centrifuge suitable for 50 ml conical tubes (floor or table-top)
  - Culture flask (500 ml)
  - Conical tubes, 50 ml, sterile
  - Snap-top tubes, 6 ml and 15 ml, sterile (Falcon 2059 or similar)
  - UV-VIS Spectrophotometer
  - Mini agarose gel apparatus
  - Media for transformation, growth, screening and expression
  - 5% SDS solution (10 ml required per transformation)

### Methods

Overview	
Introduction	The yeast <i>Pichia pastoris</i> is a microbial eukaryote which has been developed into a recombinant protein expression system. To express recombinant proteins in <i>Pichia pastoris</i> , a <i>Pichia</i> expression vector containing the gene of interest must be introduced into the host. Following introduction, the gene of interest is integrated into the host genome by a method known as gene replacement. Spheroplast transformation is an efficient method for introducing DNA into <i>Pichia pastoris</i> cells. The <i>Pichia</i> Spheroplast Kit was designed for the production of <i>Pichia pastoris</i> spheroplasts for subsequent transformation with recombinant vector constructs.
Explanation of Spheroplasting	Yeast cells have a cell wall which limits the efficiency of transformation using chemical techniques common to bacterial cells. To enable the uptake of foreign DNA into the <i>Pichia pastoris</i> cells it is necessary to partially remove the cell wall. This is done by treating the cells with the enzyme zymolyase and the reducing agent DTT. Zymolyase is a $\beta$ -glucanase that hydrolyzes glucose polymers with $\beta$ -1,3 linkages. Enzyme treatment is followed by several isotonic washes to remove the enzyme. To transform the cells, DNA is incubated with the spheroplasts. To facilitate cell wall regeneration a solution containing polyethylene glycol (PEG) is added followed by a 10 minute incubation. Following removal of the PEG, the cells are resuspended in an isotonic regeneration by the spheroplast method may vary, though it is typically in the range of 10 <sup>3</sup> to 10 <sup>4</sup> colonies per microgram of DNA.
The <i>Pichia</i> Spheroplast Kit	The <i>Pichia</i> Spheroplast Kit contains buffers, media and reagents required for 10 spheroplast preparations. Each preparation of spheroplasts can be used for up to 5 transformations. The formation of spheroplasts is a carefully controlled procedure. All of the components in the kit have been tested to ensure the production of 70% spheroplasts in less than 30 minutes of zymolyase treatment. In addition, reagents, stock solutions, and final solutions are tested for sterility.
Guidelines for Growing <i>P. pastoris</i>	The growth temperature for <i>P. pastoris</i> is 30°C for liquid cultures, plates and slants. Log phase cultures of <i>Pichia pastoris</i> have a doubling time of approximately 2 hours in YPD media. When methanol utilization plus (Mut <sup>+</sup> ) recombinant <i>Pichia</i> strains are grown on minimal methanol (MM) plates or in liquid MM media, it is advisable to add methanol every two days to compensate for methanol lost from the medium due to evaporation. For plates, add 100 $\mu$ l of 100% methanol to the lid of the inverted plate. For liquid medium, add methanol directly to the liquid to a final concentration of 0.5%.
Guidelines for Storing <i>P. pastoris</i>	For medium-term storage of cells, use YPD medium or YPD agar slants. The cells can be stored in such media for several months at +4°C. For permanent storage, grow an overnight culture in YPD media. Add glycerol to the culture to a final concentration of 15%. Store the culture at -80°C.

### Preparing the Transforming DNA and Pichia Cells

Introduction	This section provides guidelines for preparing your transforming DNA and <i>Pichia</i> cells for the transformation procedure.
Preparing the Plasmid DNA	Your gene of interest should be cloned into a <i>Pichia</i> expression vector and the correct construction confirmed by sequencing. The DNA needed to generate His <sup>+</sup> <i>Pichia</i> transformants does not need to be ultra-pure, but must be pure enough to be successfully digested with the appropriate restriction enzyme prior to transformation. Cleaner DNA will result in better transformation efficiencies. Additional guidelines are provided below for preparing your transforming DNA.
Digesting the Transforming DNA	Digest <i>Pichia</i> constructions with the appropriate enzyme to yield His <sup>+</sup> Mut <sup>-</sup> (typically <i>Not</i> I or <i>Bgl</i> II) or His <sup>+</sup> Mut <sup>+</sup> (typically <i>Sal</i> I or <i>Stu</i> I) transformants.
	The digested plasmid DNA can be generated ahead of time and stored frozen. It is not necessary to purify the portion of the plasmid containing the <i>AOX1</i> sequences and the gene of interest away from the rest of the plasmid prior to transformation.
	We recommend that the success of the digestion of the plasmid be confirmed by agarose gel electrophoresis. Complete digestion is not critical, however, the number of His <sup>+</sup> transformants obtained will be reduced if the DNA is not completely digested.
	We also recommend that you include a plasmid-only (no insert) negative control in your transformation experiment to determine the protein background during the SDS-PAGE analysis.
Media	Prepare the following media several days in advance and store at +4°C (see <b>Recipes</b> , pages 7-9 for details).
	• YPD (Yeast extract Peptone Dextrose) medium, 1 liter
	<ul><li>YPD plates, 1 liter</li><li>RDB (Regeneration Dextrose Base) plates, 1 liter</li></ul>
	<ul> <li>RDHB (Regeneration Dextrose Histidine Base) plates, 1 liter</li> </ul>
	Prepare the following solution on the day of transformation and maintain at 45°C.
	• 5% SDS solution in water
	• RD (Regeneration Dextrose), molten agarose, 100 ml
	continued on next page

### Preparing the Transforming DNA and Pichia Cells, continued

Growing the <i>Pichia</i> Cells	1.	Streak <i>Pichia</i> strain onto a YPD plate such that isolated, single colonies will grow. Incubate the plate at 30°C for 2 days.
	2.	Inoculate 10 ml of YPD in a 50 ml conical tube or a 100 ml flask with a single colony of <i>Pichia</i> cells from the YPD plate and grow overnight at 30°C with vigorous shaking (> 200 rpm). This culture will be used as the inoculum in Step 3 and may be stored at $+4^{\circ}$ C for 3-4 days.
	3.	Place 200 ml of YPD in each of three 500 ml culture flasks. Inoculate the flasks with 5, 10, and 20 $\mu$ l of cells from Step 2 and incubate them overnight with vigorous shaking at 30°C.
	4.	The next morning, equilibrate the transformation solutions (SE, SCE, Sterile Water, SOS, PEG, CaS, CaT, 1 M sorbitol) provided in the kit, the RDB plates (for plating transformants), and the RDHB plates (for viability control) by bringing them to room temperature.
	5.	Check the $OD_{600}$ of each of the three culture flasks. Harvest the cells from the culture which has an $OD_{600}$ between 0.2 and 0.3 (the optical density of the cells is important for the successful formation of spheroplasts). The cells are harvested by centrifugation at room temperature for 5-10 minutes at 1500 x g. Use these cells for preparing the spheroplasts for transformation. Decant the supernatant and discard the other cultures.
		<b>Note</b> : If the cells have grown beyond an $OD_{600}$ of 0.3, dilute the culture with fresh media to an $OD_{600}$ less than 0.1 and allow the cells to grow back up to the correct density. During log phase, the doubling time of <i>Pichia</i> cells

is approximately 2 hours.

### **Preparing Spheroplasts**

Before Starting	<ul> <li>You should have a cell pellet from Step 5, previous page.</li> <li>Prepare 100 ml of molten RD agarose and keep at 45°C (see <b>Recipes</b>, page 9)</li> </ul>
	• Thaw one tube of 1 M DTT (provided)
	Prepare fresh SED for one batch of spheroplasts as follows:
	Using sterile technique, transfer 19 ml of SE (provided) to an appropriate sterile container ( <i>e.g.</i> 50 ml conical tube). Add 1 ml of 1 M DTT and mix well. For best results this solution of SED should be made and used immediately.
Important Important	The quality and freshness of DTT is critical for a successful spheroplast preparation. The 1 M DTT provided is analytical reagent grade and must be stored at -20°C.
Washing the Cells	<ol> <li>Wash the cells from Step 5, previous page by resuspending the pellet in 20 ml of sterile water (provided). Resuspend the pellet by swirling the tube. Transfer to a sterile, 50 ml conical tube.</li> </ol>
	<ol> <li>Pellet the cells by centrifugation at 1500 x g for 5 minutes at room temperature. Decant and discard the supernatant. Use the cell pellet to prepare spheroplasts.</li> </ol>
	3. Wash the cell pellet once by resuspending in 20 ml of fresh SED, prepared above and centrifuge at 1500 x g for 5 minutes at room temperature.
	4. Wash the cells once with 20 ml of 1 M sorbitol and centrifuge as described in Step 2.
	5. Resuspend the cells by swirling in 20 ml of SCE buffer and divide the suspension into two 50 ml conical tubes (~10 ml each).
	6. Remove one tube of Zymolyase from -20°C and place it on ice. Mix well by flicking the tube several times. Zymolyase is provided as a slurry and does not go into solution. It is important to mix the slurry thoroughly before each use to ensure addition of a consistent amount of Zymolyase.
Adding Zymolyase	You will use one tube of cells prepared above to determine the optimal time of digestion with Zymolyase to make spheroplasts. Once the optimal time is determined, use the other tube of cells to make spheroplasts.
	Zymolyase digests the cell wall and makes the cells extremely fragile. Handle the sample <b>gently</b> . The moment after addition of Zymolyase, digestion of the cell wall begins.
	• Prepare at least 20 ml of a 5% SDS solution (not provided) for.
	- Set your UV-Vis spectrophotometer to 800 nm and blank with 800 $\mu l$ 5% SDS and 200 $\mu l$ SCE.
	• Set up 17 sterile microcentrifuge tubes and label them 0, 2, 4, 5, 6, 7, 8, 9, 10, 15, 20, 25, 30, 35, 40, 45, and 50. Add 800 μl of 5% SDS to each tube.

## Preparing Spheroplasts, continued

a one tube of cells (Step 5, previous spage), withdraw 200 µl cells and to the tube marked "0". This is your zero time point. Set the tube aside e. 7.5 µl of Zymolyase to the same tube of cells, mix gently by inversion, ncubate the cells at 30°C. <b>Do not shake the sample.</b> This sample will be to establish the incubation time for optimal spheroplasting. Keep econd tube of cells at room temperature for use in Step 6. Keep the ining Zymolyase on ice. itor the formation of the spheroplasts as follows: At time 2 minutes, draw 200 µl of cells (from the suspension in Step 2) and add to the tube ted "2". Repeat at time t= 4, 5, 6, 7, 8, 9, 10, 15, 20, 25, 30, 35, 40, 45, and inutes after adding Zymolyase. Read the OD800 for all samples. rmine the percent of spheroplasting for each time point using the tion: heroplasting = 100 - [(OD800 at time t/OD800 at time 0) x 100] <b>xample:</b> At time t = 0, the OD800 = 0.256 At time t = 15, the OD800 = 0.032 allation: % spheroplasting = 100-[(0.032/0.256) x 100] = 100 - [(0.125)x100] = 100 - [(0.125)x100]
ncubate the cells at 30°C. <b>Do not shake the sample.</b> This sample will be to establish the incubation time for optimal spheroplasting. Keep econd tube of cells at room temperature for use in Step 6. Keep the ining Zymolyase on ice. itor the formation of the spheroplasts as follows: At time 2 minutes, draw 200 µl of cells (from the suspension in Step 2) and add to the tube eted "2". Repeat at time t= 4, 5, 6, 7, 8, 9, 10, 15, 20, 25, 30, 35, 40, 45, and inutes after adding Zymolyase. Read the OD800 for all samples. rmine the percent of spheroplasting for each time point using the tion: heroplasting = 100 - [(OD800 at time t/OD800 at time 0) x 100] <b>xample</b> : At time t = 0, the OD800 = 0.256 At time t = 15, the OD800 = 0.032 alation: % spheroplasting = 100-[(0.032/0.256) x 100] = 100 - [(0.125)x100]
draw 200 µl of cells (from the suspension in Step 2) and add to the tube ted "2". Repeat at time t= 4, 5, 6, 7, 8, 9, 10, 15, 20, 25, 30, 35, 40, 45, and inutes after adding Zymolyase. Read the OD800 for all samples. rmine the percent of spheroplasting for each time point using the tion: heroplasting = 100 - [(OD800 at time t/OD800 at time 0) x 100] <b>xample</b> : At time t = 0, the OD800 = 0.256 At time t = 15, the OD800 = 0.032 allation: % spheroplasting = 100-[(0.032/0.256) x 100] = 100 - [(0.125)x100]
tion: heroplasting = 100 - [(OD800 at time t/OD800 at time 0) x 100] xample: At time t = 0, the OD800 = 0.256 At time t = 15, the OD800 = 0.032 alation: % spheroplasting = 100-[(0.032/0.256) x 100] = 100 - [(0.125)x100]
xample: At time t = 0, the OD800 = 0.256 At time t = 15, the OD800 = 0.032 alation: % spheroplasting = $100 - [(0.032/0.256) \times 100]$ = $100 - [(0.125) \times 100]$
At time t = 15, the OD800 = $0.032$ alation: % spheroplasting = $100 - [(0.032/0.256) \times 100]$ = $100 - [(0.125) \times 100]$
$= 100 - [(0.125) \times 100]$
= 100 - 12.5 = 87.5%
rmine the time of incubation that results in approximately 70% roplasting. This time of incubation is variable due to differences in of Zymolyase. In Invitrogen labs, it takes approximately 15-40 minutes rmolyase treatment to achieve optimal spheroplasting.
: It is important to establish the minimum time required for the ed amount of spheroplasting. Prolonged incubation with Zymolyase is erious to spheroplasts and will result in lower transformation ency.
7.5 μl Zymolyase to the remaining tube of cells as described in Step 1. pate the tube at 30°C for the time that was established in Step 5 to n the optimal level (70%) of spheroplasting.
est the spheroplasts by centrifugation at 750 x g for 10 minutes at temperature. Decant and discard the supernatant.
Vash the spheroplasts once with 10 ml of 1 M sorbitol ( <b>gently</b> disperse ellet by tapping the tube, <b>do not vortex</b> ). Collect the spheroplasts by ifugation at 750 x g for 10 minutes at room temperature.
Vash the spheroplasts once with 10 ml of CaS and centrifuge as in Step

### Transformation

	ke sure your RDB plates are at room temperature and that you have molten top agarose available. Thaw your linearized DNA and keep on ice.
1.	For each transformation, dispense $100 \ \mu$ l of the spheroplast preparation from Step 9 (previous page) into a sterile 15 ml snap-top Falcon 2059 tube (or equivalent).
2.	Add 10 $\mu$ g of DNA and incubate the tube at room temperature for 10 minutes.
3.	During the 10 minute incubation, make a fresh PEG/CaT solution. Since each transformation requires 1.0 ml of the PEG/CaT solution, calculate the amount you need and prepare this volume by adding together equal volumes of 40% PEG and CaT (a 1:1 solution).
4.	Add 1.0 ml of fresh PEG/CaT solution to the cells and DNA, mix gently, and incubate at room temperature for 10 minutes.
5.	Centrifuge the tube at 750 x g for 10 minutes at room temperature and carefully aspirate the PEG/CaT solution. Invert the tube and tap it gently to drain the excess PEG/CaT solution.
6.	Resuspend the pellet of transformed cells in 150 $\mu$ l of SOS medium and incubate it at room temperature for 20 minutes.
7.	Add 850 µl of 1 M sorbitol. Proceed to <b>Plating</b> , below.
	<i>hia</i> spheroplasts need to be plated in top agarose or agar to protect them from is prior to selection.
1.	Mix together 100-300 $\mu$ l of each spheroplast-DNA solution from Step 7, above with 10 ml of molten RD agarose and pour on RDB plates. Allow the top agarose to harden. Note there is enough of the spheroplast-DNA solution to plate duplicate and triplicate plates.
2.	Invert plates and incubate at 28-30°C. Transformants should appear in 4-6 days.
3.	For cell viability: Mix 100 µl of spheroplasts with 900 µl of 1 M sorbitol.
4.	Mix 100 $\mu$ l of this diluted sample with 10 ml of molten RDH and pour on a RDHB plate. Allow top agarose to harden.
5.	Invert plates and incubate at 28-30°C. Appearance of colonies after 4-6 days demonstrates that the spheroplasts can regenerate into dividing cells.
	RD 1. 2. 3. 4. 5. 6. 7. <i>Pick</i> lysi 1. 2. 3. 4.

### Appendix

Recipes	
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. Store at $+4^{\circ}$ C. The shelf life of this solution is one year.
	500X B (0.02% Biotin)
	Dissolve 20 mg biotin in 100 ml of water and filter sterilize. Store at +4°C. The shelf life of this solution is approximately one year.
	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°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 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.
	100X AA (0.5% of each Amino Acid)
	Dissolve 500 mg each of L-glutamic acid, L-methionine, L-lysine, L-leucine, and L-isoleucine in 100 ml of water. Filter sterilize and store at +4°C. The shelf life of this solution is approximately one year.
YPD or YEPD	Yeast Extract Peptone Dextrose Medium (1 liter)
	1% yeast extract 2% peptone 2% dextrose (glucose)
	1. Dissolve 10 g yeast extract and 20 g of peptone in 900 ml of water. <b>Note</b> : Add 20 g of agar if making YPD slants or plates.
	2. Autoclave for 20 minutes on liquid cycle.
	3. Add 100 ml of 10X D.
	Store the liquid medium at room temperature. Store YPD slants or plates at +4°C. The shelf life is several months.
	continued on next page

### Recipes, continued

MGY and MGYH	<ul> <li>Minimal Glycerol Medium ± Histidine (1 liter)</li> <li>1.34% YNB</li> <li>1% glycerol</li> <li>4 x 10<sup>-5</sup>% biotin</li> <li>± 0.004% histidine</li> <li>1. Combine aseptically 800 ml autoclaved water with 100 ml of 10X YNB, 2 ml of 500X B, and 100 ml of 10X GY.</li> <li>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.</li> <li>Store at +4°C. The shelf life of this solution is approximately two months.</li> </ul>
RD and RDH Liquid Media	Regeneration Dextrose Medium $\pm$ Histidine (1 liter)1 M sorbitol2% dextrose1.34% YNB4 x 10 <sup>-5</sup> % biotin0.005% amino acids $\pm$ 0.004% histidine
	<ol> <li>Dissolve 186 g of sorbitol in 700 ml of water and proceed to Step 2.</li> <li>Autoclave 20 minutes on liquid cycle.</li> </ol>
	<ol> <li>Cool and maintain the liquid medium in a 45°C water bath.</li> </ol>
	<ol> <li>Prepare a prewarmed (45°C) mixture of the following stock solutions:</li> </ol>
	<ul> <li>100 ml of 10X D</li> <li>100 ml of 10X YNB</li> <li>2 ml of 500X B</li> <li>10 ml of 100X AA</li> <li>88 ml of sterile water</li> <li>Add to sorbitol solution.</li> </ul> 5. For growth of <i>his4</i> strains you must add histidine to the media. Add 10 ml of 100X H (histidine) to the prewarmed mixture in Step 4. Store liquid medium at +4°C.

continued on next page

### **Recipes**, continued

RDB and RDHB Agar Plates	<ol> <li>Dissolve 186 g of sorbitol in 700 ml of water and add 20 g of agar.</li> <li>Autoclave 20 minutes on liquid cycle.</li> <li>Place the autoclaved solution in a 60°C water bath prior to addition of prewarmed mixture of stock solutions. This will keep the medium from becoming too thick to mix reagents.</li> <li>Prepare the prewarmed (45°C) mixture from RD and RDH Liquid Media, Step 4, previous page. Add to sorbitol/agar solution. If you are selecting for His<sup>+</sup> transformants, do not add histidine.</li> </ol>
	<ol> <li>Pour the plates immediately after mixing the solutions in Step 4. The plates should be stored at +4°C and should last for several months.</li> </ol>
RD and RDH Top Agar	<ol> <li>Dissolve 186 g of sorbitol in 700 ml of water and add 10 g of agar or agarose.</li> <li>Autoclave 20 minutes on liquid cycle.</li> <li>Place the autoclaved solution in a 60°C water bath prior to addition of prewarmed mixture of stock solutions. This will keep the medium from becoming too thick to mix reagents.</li> <li>Prepare the prewarmed (45°C) mixture from RD and RDH Liquid Media, Step 4, previous page. Add to sorbitol/agar solution. If you are selecting for His<sup>+</sup> transformants, do not add histidine.</li> <li>Place the solution to 45°C after adding the solutions in Step 4. During transformation, use as a molten solution at 45°C.</li> </ol>

#### Pichia pastoris References

- Barr, K.A., Hopkins, S.A. and Sreekrishna, K. (1992). Protocol for Efficient Secretion of HSA Developed from *Pichia pastoris*. *Pharm. Eng.* **12(2)**: 48-51.
- Buckholz, R.G. and Gleeson, M.A.G. (1991). Yeast Systems for the Commercial Production of Heterologous Proteins. *Bio/Technology* **9**: 1067-1072.
- Clare, J.J., Rayment, F.B., Ballantine, S.P., Sreekrishna, K. and 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., and Henwood, C.A. (1991b). Production of Mouse 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. and Madden, K.R. (1985). *Pichia pastoris* as a Host System for Transformations. *Mol. Cell. Biol.* **5(12)**: 3376-3385.
- Cregg, J.M. (1987) Genetics of Methylotrophic Yeasts In: Proceedings of the Fifth International Symposium on Microbial Growth on C1 Compounds (Van Verseveld, H.W. and Duine, J.A., eds.), Nijhoff, Dordrecht.
- Cregg, J.M., Tschopp, J.F., Stillman, C., Siegel, R., Akong, M., Craig, W.S., Buckholz, R.G., Madden, K.R., Kellaris, P.A., Davis, G.R., Smiley, B.L., Cruze, J., Torregrossa, R., Velicelebi, G. and Thill, G.P. (1987). High-Level Expression and Efficient Assembly of Hepatitis B Surface Antigen in the Methylotrophic Yeast, *Pichia pastoris*. *Bio/Technology* 5: 479-485.
- Cregg, J.M. and Madden, K.R. (1989). Use of Site-specific Recombination to Regenerate Selectable Markers. *Mol. Gen. Genet.* **219**: 320-323.
- Cregg, J.M., Vedvick, T.S. and Raschke, W.C. (1993). Recent Advances in the Expression of Foreign Genes in *Pichia pastoris*. *Bio/Technology* **11**: 905-910.
- Digan, M.E., Lair, S.V., Brierley, R.A., Siegel, R.S., Williams, M.E., Ellis, S.B., Kellaris, P.A., Provow, S.A., Craig, W.S., Velicelebi, G., Harpold, M.M. and Thill, G.P. (1989). Continuous Production of a Novel Lysozyme via Secretion from the Yeast, *Pichia pastoris*. *Bio/Technology* **7**: 160-164.
- Dohmen, R.J., Strasser, A.W.M., Honer, C.B. and Hollenberg, C.P. (1991). An Efficient Transformation Procedure Enabling Long Term Storage of Competent Cells of Various Yeast Genera. *Yeast* 7: 691-692.
- Ellis, S.B., Brust, P.F., Koutz, P.J., Waters, A.F., Harpold, M.M. and Gingeras, T.R. (1985). Isolation of Alcohol Oxidase and Two Other Methanol Regulatable Genes from the Yeast *Pichia pastoris*. *Mol. Cell. Biol.* **5(5)**: 1111-1121.
- Gould, S.J., McCollum, D., Spong, A.P., Heyman, J.A. and Subramani, S. (1992). Development of the Yeast *Pichia pastoris* as a Model Organism for a Genetic and Molecular Analysis of Peroxisome Assembly. *Yeast* **8**: 613-628.
- Grinna, L.S. and Tschopp, J.F. (1989). Size Distribution and General Structural Features of N-linked Oligosaccharides from the Methylotrophic Yeast *Pichia pastoris*. *Yeast* **5**: 107-115.
- Hagenson, M.J., Holden, K.A., Parker, K.A., Wood, P.J., Cruze, J.A., Fuke, M., Hopkins, T.R. and Stroman, D.W. (1989). Expression of Streptokinase in *Pichia pastoris* Yeast. *Enzyme Microbiol. Technol.* **11**: 650-656.
- Ito, H., Fukuda, Y., Murata, K., and Kimura, A. (1983). Transformation of Intact Yeast Cells Treated with Alkali Cations. *J. Bacteriol.* **153**: 163-168.
- Koutz, P.J., Davis, G.R., Stillman, C., Barringer, K., Cregg, J.M. and Thill, G. (1989). Structural Comparison of the *Pichia pastoris* Alcohol Oxidase Genes. *Yeast* **5**: 167-177.

- Kumagai, M.H., Sverlow, G.G., della-Cioppa, G. and Grill, L.K. (1993). Conversion of Starch to Ethanol in a Recombinant *Saccharomyces cerevisiae* Strain Expressing Rice Amylase from a Novel *Pichia pastoris* Alcohol Oxidase Promoter. *Bio/Technology* **11**: 606-610.
- Ratner, M. (1989). Protein Expression in Yeast. Bio/Technology 7: 1129-1133.
- Romanos, M.A., Clare, J.J., Beesley, K.M., Rayment, F.B., Ballantine, S.P., Makoff, A.J., Dougan, G., Fairweather, N.F. and Charles, I.G. (1991). Recombinant *Bordetella pertussis* Pertactin (P69) from the Yeast *Pichia pastoris*: High Level Production and Immunological Properties. *Vaccine* **9**: 901-906.
- Schiestl, R.H., and Gietz, R.D. (1989). High Efficiency Transformation of Intact Yeast Cells Using Singlestranded Nucleic Acids as a Carrier. *Curr. Genet.* **16**: 339-346.
- 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. and Fuke, M. (1988). High Level Expression of Heterologous Proteins in Methylotrophic Yeast *Pichia pastoris*. *J. Basic Microbiol*. 28(4): 265-278.
- Sreekrishna, K., Nelles, L., Potenz, R., Cruze, J., Mazzaferro, P., Fish, W., Fuke, M., Holden, K., Phelps, D., Wood, P. and Parker, K. (1989). High Level Expression, Purification, and Characterization of Recombinant Human Tumor Necrosis Factor Synthesized in the Methylotrophic Yeast *Pichia pastoris*. *Biochem*. **28(9)**: 4117-4125.
- Thill, G., Davis, G., Stillman, C., Tschopp, J.F., Graig, W.S., Velicelebi, G., Greff, J., Akong, M., Stroman, D., Torregrossa, R. and Siegel, R.S. (1987) The Methylotrophic Yeast *Pichia pastoris* as a Host for Heterologous Protein Production. In: Proceedings of the Fifth International Symposium on Microbial Growth on C1 Compounds (Van Verseveld, H.W. and Duine, J.A., eds.), Nijhoff, Dordrecht.
- Tschopp, J.F., Brust, P.F., Cregg, J.M., Stillman, C.A. and Gingeras, T.R. (1987). Expression of the *lacZ* Gene from Two Methanol-Regulated Promoters in *Pichia pastoris*. *Nuc. Acids Res.* **15(9)**: 3859-3876.
- Tschopp, J.F., Sverlow, G., Kosson, R., Craig, W. and Grinna, L. (1987). High Level Secretion of Glycosylated Invertase in the Methylotrophic Yeast, *Pichia pastoris*. *Bio/Technology* **5**: 1305-1308.
- Vedvick, T., Buckholz, R.G., Engel, M., Urcan, M., Kinney, J., Provow, S., Siegel, R.S., and Thill, G.P. (1991). High-Level Secretion of Biologically Active Aprotonin from the Yeast *Pichia pastoris*. *J. Ind. Microbiol.* **7**: 197-202.
- Wagner, S.L., Siegel, R.S., Vedvick, T.S., Raschke, W.C., and Van Nostrand, W.E. (1992). High-Level Expression, Purification, and Characterization of the Kunitz-type Protease Inhibitor Domain of Protease Nexin-2/amyloid -Protein Precursor. *Biochem. Biophys. Res. Commun.* 186(2): 1138-1145.
- Wegner, G.H. and Harder, W. (1987). Methylotrophic Yeasts -- 1986. In: Proceedings of the Fifth International Symposium on Microbial Growth on C1 Compounds (Van Verseveld, H.W. and Duine, J.A., eds.), Nijhoff, Dordrecht.
- Wegner, G.H. (1990). Emerging Applications of the Methylotrophic Yeasts. *FEMS Microbiol. Rev.* 87: 279-284.

#### General References

#### Yeast Molecular Biology

Guthrie, C. and Fink, G.R., eds. <u>Guide to Yeast Genetics and Molecular Biology</u>. (1991). Methods in Enzymology, vol 194, Academic Press, San Diego.

#### **Molecular Biology Techniques**

Sambrook, J., Fritsch, E.F., and Maniatis, T. <u>Molecular Cloning, A Laboratory Manual</u>, second edition. (1989). Cold Spring Harbor Laboratory Press, Cold Spring Harbor.