

pBAD-DEST49 Gateway[®] Destination Vector

A destination vector for cloning and regulated expression of HP-thioredoxin N-terminal fusions in E. coli

Catalog no. 12283-016

Version F 29 December 2010 25-0427

User Manual

Table of Contents

Kit Contents and Storage	iv
Accessory Products	v
Methods	1
Overview	1
Using pBAD-DEST49	4
Expressing the Gene of Interest	7
Analyzing Samples	10
Appendix	13
Recipes	
Regulation by Arabinose	
Map of pBAD-DEST49	16
Features of pBAD-DEST49	17
Technical Support	
Purchaser Notification	20
References	22

Kit Contents and Storage

Shipping and	pBAD-DEST49 is shipped on wet ice. Upon receipt, store at –20°C.
Storage	pBAD-DEST49 is guaranteed for six months from the date of shipment when stored at –20°C.
Contents	6 μg of pBAD-DEST49 vector at 150 ng/μl in TE buffer, pH 8.0 (10 mM Tris-HCl, 1 mM EDTA, pH 8.0). Volume: 40 μl.

Accessory Products

Additional T Products r

The products below are also available from Invitrogen. For more details visit our Web site at <u>www.invitrogen.com</u>, or contact Technical Support (page 18).

Product	Quantity	Catalog no.
Gateway [®] LR Clonase [™] Enzyme Mix	20 reactions	11791-019
One Shot [®] TOP10 Chemically Competent E. Coli	10 reactions	C4040-10
Cells	20 reactions	C4040-03
One Shot [®] TOP10 Electrocompetent E. Coli Cells	10 reactions	C4040-50
	20 reactions	C4040-52
One Shot [®] ccdB Survival [™] 2 T1 Phage-Resistant	10 reactions	A10460
Cells		
pENTR Directional TOPO [®] Cloning Kit	20 reactions	K2400-20
PureLink™ HiPure Plasmid Midiprep Kit	25 preps	K2100-04
EKMax [™] Enterokinase	250 units	E180-01
	1000 units	E180-02
EK-Away [™] Resin	7.5 ml	R180-01

Detection of Recombinant Proteins

Expression of your recombinant fusion protein can be detected using an antibody to the appropriate epitope. Horseradish peroxidase (HRP) or alkaline phosphatase (AP)conjugated antibodies allow one-step detection using colorimetric or chemiluminescent detection methods. The antibody volume supplied is sufficient for 25 Western blots.

Product	Quantity	Catalog no.	Epitope
Anti-Thio™ Antibody	50 µl	R920-25	Detects His-Patch thioredoxin fusion proteins Note: The exact epitope detected by this antibody has not been mapped.
Anti-V5 Antibody Anti-V5-HRP Antibody	50 μl 50 μl	R960-25 R961-25	Detects 14 amino acid epitope derived from the P and V proteins of the paramyxovirus,
Anti-V5-AP Antibody	125 µl	R962-25	SV5 (Southern <i>et al.,</i> 1991) GKPIPNPLLGLDST
Anti-His (C-term) Antibody	50 µl	R930-25	Detects the C-terminal polyhistidine (6xHis) tag,
Anti-His (C-term)- HRP Antibody	50 µl	R931-25	requires the free carboxyl group for detection (Lindner <i>et</i>
Anti-His (C-term)- AP Antibody	125 µl	R932-25	al., 1997) НННННН-СООН

Accessory Products, Continued

Purification of Recombinant Fusion Protein	If your gene of interest is in frame with the C-terminal peptide containing the V5 epitope and the polyhistidine (6xHis) tag, you may use Immobilized Metal Affinity Chromatography (IMAC) to purify your recombinant fusion protein.
	The ProBond [™] Purification System or bulk ProBond [™] resin are available separately from Invitrogen. For more details

or contact Technical Support (page 18).

Product	Quantity	Catalog no.
ProBond [™] Nickel-chelating Resin	50 ml	R801-01
_	150 ml	R801-15
Purification Columns	50 each	R640-50
(10 ml polypropylene columns, empty)		
ProBond [™] Purification System	6 purifications	K850-01
Positope [™] Control Protein	5 µg	R900-50

GelA wide range of pre-cast NuPAGE® and Novex® Tris-ElectrophoresisGlycine polyacrylamide gels and electrophoresis
apparatuses are available to facilitate separation and
visualization of your recombinant fusion protein.

For more details on these products, visit our Web site at <u>www.invitrogen.com</u> or contact Technical Support (page 18).

on these products, visit our Web site at www.invitrogen.com

Introduction

Overview	
Description	pBAD-DEST49 is a 6.2 kb vector derived from pBAD/Thio and adapted for use with the Gateway [®] Technology. pBAD- DEST49 is designed for regulated expression of N-terminal HP-thioredoxin fusion proteins in <i>E. coli</i> .
	For a map of pBAD-DEST49, see page 16.
Features	pBAD-DEST49 contains the following features:
	 araBAD (pBAD) promoter: provides tight, dose- dependent regulation of heterologous gene expression
	• His-Patch Thioredoxin (HP-thioredoxin) fusion partner for efficient translation of the protein of interest and, in some cases, increased solubility
	• Two recombination sites, <i>att</i> R1 and <i>att</i> R2, downstream of the HP-thioredoxin gene for recombinational cloning of the gene of interest from an entry clone
	• Chloramphenicol resistance gene located between the two <i>att</i> R sites for counterselection
	• The <i>ccd</i> B gene located between the two <i>att</i> R sites for negative selection
	• The V5 epitope and 6xHis tag for detection and purification (optional)
	• <i>rrn</i> B transcription terminator for strong transcription termination
	• The pUC origin for replication and maintenance of the plasmid in <i>E. coli</i> .
	Note: Although pBAD-DEST49 contains a pUC origin, it acts as a low-copy number plasmid, resulting in lower yields of the vector.
	• The ampicillin (<i>bla</i>) resistance gene for selection in <i>E. coli</i>
	• The <i>ara</i> C gene for regulation of the <i>ara</i> BAD promoter (Lee, 1980; Schleif, 1992)

Overview, Continued

The Gateway [®] Technology	Gateway [®] is a universal cloning technology that takes advantage of the site-specific recombination properties of bacteriophage lambda (Landy, 1989) to provide a rapid and highly efficient way to move your gene of interest into multiple vector systems. To express your gene of interest using Gateway [®] Technology, simply:
	 Clone your gene of interest into a Gateway[®] entry vector to create an entry clone.
	2. Generate an expression clone by performing an LR recombination reaction between the entry clone and a Gateway [®] destination vector (<i>e.g.</i> pBAD-DEST49).
	3. Transform your expression clone into <i>E. coli</i> . Induce expression of your gene of interest using L-arabinose.
	For more information on the Gateway [®] Technology, refer to the Gateway [®] Technology Manual. This manual is available for downloading from our Web site (<u>www.invitrogen.com</u>) or by contacting Technical Support (page 18).
Regulation of Expression by L-Arabinose	Expression of pBAD is induced in the presence of L-arabinose. In the absence of L-arabinose, only very low levels of transcription are observed from pBAD (Lee, 1980; Lee <i>et al.</i> , 1987). Uninduced levels are further repressed by growth in the presence of glucose. Glucose reduces the levels of 3', 5'-cyclic AMP, lowering expression from the catabolite- repressed pBAD promoter (Miyada <i>et al.</i> , 1984). By varying the concentration of L-arabinose, protein expression levels can be optimized to ensure maximum protein expression. In addition, the tight regulation of pBAD by AraC is useful for expression of potentially toxic or essential genes (Carson <i>et al.</i> , 1991; Dalbey and Wickner, 1985; Guzman <i>et al.</i> , 1992; Kuhn and Wickner, 1985; Russell <i>et al.</i> , 1989; San Millan <i>et al.</i> , 1989).
	For additional information on the mechanism of expression and repression of the <i>ara</i> regulon, see page 15.
	Continued on next page

Overview, Continued

Thioredoxin	The 11.7 kDa thioredoxin protein is found in yeast, plants, and mammals, as well as in bacteria. It was originally isolated from <i>E. coli</i> as a hydrogen donor for ribonuclease reductase (see Holmgren, 1985 for a review). The gene has been completely sequenced (Wallace and Kushner, 1984), and the protein has been crystallized and its three-dimensional structure determined (Katti <i>et al.</i> , 1990). When overexpressed in <i>E. coli</i> , thioredoxin is able to accumulate to approximately 40% of the total cellular protein and still remain soluble. When used as a fusion partner, thioredoxin can increase translation efficiency and, in some cases, solubility of eukaryotic proteins expressed in <i>E. coli</i> .
	Examples of eukaryotic proteins that have been produced as soluble C-terminal fusions to the thioredoxin protein in <i>E. coli</i> (LaVallie <i>et al.,</i> 1993) include:
	Murine interleukin-2
	• Human interleukin-3
	Murine interleukin-4
	• Murine interleukin-5
	Human macrophage colony stimulating factor
	Murine steel factor
	Murine leukemia inhibitory factor
	• Human bone morphogenetic protein-2
His-Patch Thioredoxin	To create a metal binding domain in the thioredoxin protein, the glutamate residue at position 32 and the glutamine residue at position 64 were mutated to create histidine residues. When His-Patch thioredoxin folds, the histidines at positions 32 and 64 interact with a native histidine at position 8 to form a "patch". This histidine patch has high affinity for divalent cations (Lu <i>et al.</i> , 1996). His-Patch thioredoxin (HP-thioredoxin) proteins can therefore be purified on metal-chelating resins (<i>e.g.</i> ProBond TM , page vi).

Methods

Using pBAD-DEST49

Q Important	The pBAD-DEST49 vector is supplied as a supercoiled plasmid. Although Invitrogen has previously recommended using a linearized destination vector for more efficient recombination, further testing has found that linearization of this vector is not required to obtain optimal results for any downstream application.
Propagating pBAD-DEST49	If you wish to propagate and maintain pBAD-DEST49, we recommend using One Shot [®] <i>ccdB</i> Survival [™] 2 T1 Phage-Resistant Cells for transformation (page v). The <i>ccdB</i> Survival [™] 2 T1 Phage-Resistant <i>E. coli</i> strain is resistant to CcdB effects and can support the propagation of plasmids containing the <i>ccdB</i> gene.
	Note: Do not use general <i>E. coli</i> cloning strains including TOP10 or DH5 α for propagation and maintenance as these strains are sensitive to CcdB effects.
Entry Clone	To recombine your gene of interest into pBAD-DEST49, you should have an entry clone containing your gene of interest. For your convenience, Invitrogen offers the pENTR Directional TOPO [®] Cloning Kit (page v) for 5-minute cloning of your gene of interest into an entry vector.
	For detailed information on constructing an entry clone, refer to the specific entry vector manual. See the next page for factors to consider when designing your entry clone.
Specific Features	pBAD-DEST49 is designed with the following features to facilitate expression:
	• Initiation ATG is correctly spaced from the optimized ribosome binding site to ensure optimal translation of HP-thioredoxin fusions
	• HP-thioredoxin acts as a translation leader to facilitate high-level expression and in some cases, solubility
	• HP-thioredoxin can be removed after protein purification using enterokinase (<i>e.g.</i> EKMax [™] , page v).

Using pBAD-DEST49, Continued

Points to	It is very important to have a properly designed entry clone
Consider	before recombining with pBAD-DEST49. Refer to the table
Before	below and the Recombination Region on page 6.
Recombining	

If you wish to	Then your gene in the entry clone
clone in frame with HP-thioredoxin	must be in frame with the HP- thioredoxin gene after recombination.
include the V5 epitope and 6xHis tag	must not contain a stop codon and must be in frame with the C-terminal tag after recombination.
exclude the V5 epitope and 6xHis tag	must contain a stop codon.

Recombining Your Gene of Interest	Each entry clone contains <i>att</i> L sites flanking the gene of interest. Genes in an entry clone are transferred to the destination vector backbone by mixing the DNAs with the Gateway [®] LR Clonase [™] enzyme mix. The resulting recombination reaction is then transformed into <i>E. coli</i> and the expression clone selected. Recombination between the <i>att</i> R sites on the destination vector and the <i>att</i> L sites on the entry clone replaces the <i>ccd</i> B gene and the chloramphenico. (Cm ^R) gene with the gene of interest and results in the formation of <i>att</i> B sites in the expression clone.	
	Follow the instructions in the Gateway [®] Technology Manual to set up the LR Clonase TM reaction, transform a <i>recA endA E. coli</i> strain (<i>e.g.</i> TOP10 or DH5 α), and select for the expression clone.	
Resuspending pBAD-DEST49	Before you perform the LR Clonase [™] reaction, resuspend pBAD-DEST49 to 50–150 ng/µl in sterile water.	

Using pBAD-DEST49, Continued

Confirming the Expression Clone			ve wi Tr ge To LE ex	The <i>ccd</i> B gene mutates at a very low frequency, resulting in a very low number of false positives. True expression clones will be ampicillin-resistant and chloramphenicol-sensitive. Transformants containing a plasmid with a mutated <i>ccd</i> B gene will be both ampicillin- and chloramphenicol-resistant. To check your putative expression clone, test for growth on LB plates containing 30 μ g/ml chloramphenicol. A true expression clone will not grow in the presence of chloramphenicol.													
	omi jion	bina	tion	fro	om pl	BAD	-DES	ST49	×ent	ry cl	one i	s sho	own			ulting	5
				Fe •	Sha trar reco	ded sfer	regio red f natio	ons cons rom on. N	orres the e on-s	hade	d to t clone	hose e inte	DN. pBA	AD-E	DEST	ces 49 by rom 1	
				•	The cor	e und respo	lerlir ond t	ied n o bas	ucleo ses 72		d 24	08, re				egior the	l
				His	-Patch												
640				lys AAA	glu	phe	leu										
	CCA	GTC	AAC	TTT	CTC	AAG	GAG	CTG	CGA	TTG	GAC	CGG	CCG	AGA	CCT	AGG	CCA
	Ento	rokinas		ognitior	Sito	Entero	kinase	cleava	ge site			725					
				asp		1011	alv	ilo	ilo	thr	cor	1.1	tur	1.75	1.70	ala	alv
691	GAT	GAC	GAT	GAC CTG	AAG	CTG	GGA	ATT	ATC	ACA	AGT	T <u>T</u> G	TAC	AAA	AAA	GCA	GGC
										L				ttB1			_
						-	240				-	-		-		-	
742	 TNN	· · ·	· · · ·	NAC	pro	ala GCT	pne TTC	TTC	tyr Tac	lys aaa	Val GTG	Val GTG	11e ATC	TÀR VYC	Leu	GIU	
	ANN	_ <u>G</u> E	NE_	NAC NTG	GGT	CGA	AAG	AAC	ATG	TTT	CAC	CAC	TAG	TTC	GAA	CTT	
				L			att	B2									
							V5 epi										
	gly	lys	pro	ile	pro	asn	pro	leu	leu	gly	leu	asp	ser	thr	arg	thr	
2436				ATC TAG													
				6xHis	taq												
	gly	his	his	his		his	his	٦***									
2484	GGT	CAT	CAT	CAC	CAT	CAC	CAT	TGA									
	CCA	TTG	CCA	GTA	GTA	GTG	GTA	GTG	GTA	ACT	CAA	ATT	TGC	CAG	AGG	TCG	

Expressing the Gene of Interest

Introduction	Once you have selected your expression clone, you are ready to transform it into <i>E. coli</i> and test for expression of your gene. Since each recombinant protein has different characteristics that may affect optimal expression, it is helpful to vary the L-arabinose concentration and/or run a time course of expression to determine the best conditions for optimal expression of your particular protein.
Plasmid Preparation	You may prepare plasmid DNA using your method of choice. We recommend using the PureLink [™] HiPure Plasmid Midiprep Kit (page v) for isolation of pure plasmid DNA. Note that since you are purifying a vector that acts as a low- copy number plasmid, you may need to increase the amount of bacterial culture that you use to prepare your plasmid construct.
<i>E. coli</i> Strain	We recommend that you transform your construct into an <i>E. coli</i> strain that is deficient for the <i>ara</i> operon. TOP10 competent cells are deleted for <i>ara</i> BADC and are available from Invitrogen as chemically competent or electrocompetent cells (page v).
L-Arabinose	Prepare a 20% solution of L-arabinose in deionized water and filter sterilize. Store at room temperature. To determine the optimal concentration for expression of
	your particular protein, see the next page. D-arabinose will not induce expression and should not be
	used.
	Note: L-arabinose can be ordered from Sigma (Catalog no. A3256).
	Continued on worth nace

Expressing the Gene of Interest, Continued

Basic Strategy	Once you have clones that you wish to characterize, we recommend the following strategy to determine the optimal expression level.
	1. Pilot Expression: In this expression experiment you will vary the amount of arabinose over a 10,000-fold range (0.00002% to 0.2%) to determine the approximate amount of arabinose needed for maximum expression of your protein. (See next page for protocol).
	2. To optimize expression of your protein, you may wish to try arabinose concentrations spanning the amount determined in Step 1, or you may perform a time course.
	Note : If your protein is insoluble, remember to analyze the supernatant and the pellet of lysed cells for expression of soluble protein (see page 10).
Materials Needed	Be sure to have the following solutions and equipment on hand before starting the experiment:
	 SOB or LB medium containing 50–100 µg/ml ampicillin (see pages 13-14 for recipes)
	• 37°C shaking incubator
	• 20% L-arabinose
	• 37°C heat block or water bath
	• 42°C water bath
	Liquid nitrogen
	• 1X and 2X SDS-PAGE sample buffer
	Reagents and apparatus for SDS-PAGE gel
	Boiling water bath
	• Lysis Buffer (see page 14 for recipe)
	Sterile water
	Continued on next page

Expressing the Gene of Interest, Continued

Pilot Expression	 In addition to testing your transformants, we recommend that you include cells without a vector as a negative control. 1. For each transformant or control, inoculate 2 ml of SOB or LB medium containing 50–100 µg/ml ampicillin with a single recombinant <i>E. coli</i> colony. 2. Grow cells overnight on a 37°C shaking incubator at 225–250 rpm. The cells should have an optical density at 600 (OD₆₀₀) of 1–2 3. Label five tubes 1 through 5 and add 10 ml of SOB or LB medium containing 50–100 µg/ml ampicillin. 4. Inoculate each tube with 0.1 ml of the overnight culture. 5. Grow the cultures in the tubes on a 37°C shaking incubator (use vigorous shaking) Grow cells to an OD₆₀₀ of ~0.5 (the cells should be in mid-log phase). 6. While the cells are growing, prepare four 10-fold serial dilutions of 20% L-arabinose with sterile water using aseptic technique (<i>e.g.</i> 2%, 0.2%, and 0.002%). 7. Take a 1 ml aliquot of grown cells from each tubes. Centrifuge at maximum speed in a microcentrifuge for 30 seconds, and aspirate the supernatant. 8. Freeze the cell pellet at -20°C. This is the zero time point-sample. 9. Using the solutions prepared in Step 6, add the following volumes of L-arabinose to the five 9 ml cultures. Note: It is only necessary to test the highest concentration of L-arabinose. 			
	L-aral	pinose.	-	
	Tube	Stock Solution		Final Concentration
	1	0.002%	0.09	0.00002%
	2	0.02%	0.09	0.0002%
	3	0.2%	0.09	0.002%
	4	2%	0.09	0.02%
	5	20%	0.09	0.2%

10. Grow at 37°C with shaking for 4 hours.

- 11. Take 1 ml samples at 4 hours and treat samples as in Steps 7 and 8. You will have a total of 10 samples for each transformant and two samples for the negative control.
- 12. Proceed to Analyzing Samples, next page.

Analyzing Samples

Preparation of Samples	Before starting, prepare SDS-PAGE gels or use one of the pre-cast polyacrylamide gels available from Invitrogen (page vi) to analyze all the samples you collected. Note : If you wish to analyze your samples for soluble protein, see the section below.
	 When all the samples have been collected from Steps 8 and 11 in the Pilot Expression, resuspend each pellet in 100 μl of 1X SDS-PAGE sample buffer.
	2. Heat for 5 minutes at 70°C and centrifuge briefly.
	3. Load 5–10 μl of each sample on an SDS-PAGE gel and electrophorese. Save your samples by storing at –20°C.
Preparing Samples for	 Thaw and resuspend each pellet in 500 µl of Lysis Buffer (see page 14 for recipe).
Soluble/ Insoluble Protein	2. Freeze sample in dry ice or liquid nitrogen and then thaw at 42°C. Repeat 2 to 3 times.
FIOLEIN	Note: To facilitate lysis, you may need to add lysozyme or sonicate the cells.
	3. Centrifuge samples at maximum speed in a microcentrifuge for 1 minute at +4°C to pellet insoluble proteins. Transfer supernatant to a fresh tube and store on ice.
	4. Mix together equivalent amounts of supernatant and 2X SDS-PAGE sample buffer and boil for 5 minutes.
	5. Add 500 μl of 1X SDS-PAGE sample buffer to the pellets from Step 3 and boil 5 minutes.
	6. Load 10 μl of the supernatant sample and 5μl of the pellet sample onto an SDS-PAGE gel and electrophorese.
	Continued on next page

Analyzing Samples, Continued

Analyzing Samples	To determine the success of your expression experiment, you may want to perform the following types of analyses:			
	1. Stain the polyacrylamide gel with Coomassie blue and look for a band of increasing intensity in the expected size range for the recombinant protein. Use the uninduced culture as a negative control.			
	2. Perform a Western blot to confirm that the overexpressed band is your desired protein (see below).			
	3. Determine the approximate arabinose concentration for maximum expression.			
Detection of Recombinant Fusion Proteins	To detect expression of your recombinant fusion protein by Western blot analysis, you may use antibodies against the appropriate epitope (page v) or an antibody to your protein of interest. In addition, the Positope [™] Control Protein (page vi) is available from Invitrogen for use as a positive control for detection of fusion proteins containing a thioredoxin, V5, or C-terminal 6xHis epitope.			
Note	Expression of your protein with the N-terminal HP-thioredoxin peptide and the C-terminal tag will increase the size of your protein by approximately 14 kDa and 4 kDa, respectively.			

Analyzing Samples, Continued

Optimization of Expression	Once you have detected expression of your protein of interest, you may wish to perform experiments to further optimize expression. Use the Pilot Expression protocol (page 9), but vary the L-arabinose concentration over a smaller range. For example, if you obtained the best expression at 0.002%, try 0.0004%, 0.0008%, 0.001%, 0.004%, and 0.008%.
	You may also perform a time course of induction to determine if varying the time increases expression. Take time points every hour, over a 5 to 6 hour period.
	If your protein is insoluble, you may wish to analyze the supernatant and pellet of lysed cells when you vary the L-arabinose concentration (see Preparing Samples for Soluble/Insoluble Protein , page 10).
	Remember to store your cell lysates at –20°C.
Expression of Toxic Proteins	To ensure low levels of expression, you may want to utilize glucose to further repress the <i>ara</i> BAD promoter. Follow the Pilot Expression protocol (page 9) using SOB or LB containing 50–100 μ g/ml ampicillin plus glucose at all steps (see pages 13-14 for recipes).
Purification of Recombinant Fusion Proteins	The presence of the C-terminal polyhistidine (6xHis) tag in your recombinant fusion protein allows use of a metal- chelating resin such as ProBond [™] (page vi) to purify your fusion protein. Note: Other metal-chelating resins and purification methods are suitable.
Removal of the N-terminal Leader by Enterokinase	The enterokinase (EK) recognition site can be used to remove the N-terminal leader from your recombinant fusion protein after purification. Note that after digestion with enterokinase, there will be twelve vector-encoded amino acids remaining at the N-terminus of the protein (see diagram on page 6). A recombinant preparation of the catalytic subunit of bovine enterokinase (EKMax [™]) is available from Invitrogen (see page v).

Appendix

Recipes

LB (Luria-Bertani) Medium and Plates

1.0% Tryptone 0.5% Yeast Extract 1.0% NaCl 0.1% glucose (optional) pH 7.0

- 1. Dissolve 10 g tryptone, 5 g yeast extract, and 10 g NaCl in 950 ml deionized water.
- 2. Adjust solution to pH 7.0 with NaOH and add deionized water to 1 liter.
- 3. Autoclave on liquid cycle for 20 minutes at 15 psi. Allow solution to cool to 55°C. Add antibiotic if needed.
- 4. Add 5 ml of a 20% sterile glucose solution (optional).
- 5. Store at room temperature or at $+4^{\circ}$ C.

LB agar plates

- 1. Prepare LB medium as above, but add 15 g/L agar before autoclaving.
- 2. Autoclave on liquid cycle for 20 minutes at 15 psi.
- 3. After autoclaving, cool to ~55°C. Add antibiotic if needed.
- 4. Add 5 ml of a 20% sterile glucose solution (optional).
- 5. Pour into 10 cm plates and let harden. Invert plates and store at +4°C in the dark.

Recipes, Continued

SOB Medium	2% Tryptone					
	0.5% Yeast Extract					
	0.05% NaCl					
	2.5 mM KCl					
	10 mM MgCl ₂					
	0.1% glucose (optional)					
	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.					
	 Autoclave this solution and cool to ~55°C. Add 10 ml of sterile 1 M MgCl₂. Add antibiotic if needed. 					
	5. Add 5 ml of a 20% sterile glucose solution (optional).					
	6. Store at +4°C. Medium is stable for only 1–2 weeks.					
Lysis Buffer	50 mM potassium phosphate, pH 7.8 400 mM NaCl 100 mM KCl 10% glycerol 0.5% Triton X-100 10 mM imidazole					
	 Prepare 1 M stock solutions of KH₂PO₄ and K₂HPO₄. For 100 ml, dissolve the following reagents in 90 ml of deionized water: anl KH₂PO₄ ml K₂HPO₄ g NaCl r5 g KCl ml glycerol ml Triton X-100 mg imidazole 					
	3. Mix thoroughly and adjust pH to 7.8 with HCl. Bring the volume to 100 ml.					

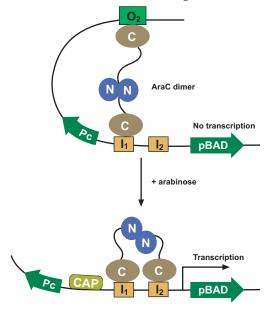
4. Store at +4°C.

Regulation by L-Arabinose

Regulation of the P_{BAD} Promoter

The *ara*BAD promoter used in pBAD-DEST49 is both positively and negatively regulated by the product of the *ara*C gene (Ogden *et al.*, 1980; Schleif, 1992). The AraC protein is a transcriptional regulator that forms a complex with L-arabinose. In the absence of L-arabinose the AraC dimer contacts the O_2 and I_1 half sites of the *ara*BAD operon, forming a 210 bp DNA loop (see figure below). For maximum transcriptional activation two events are required.

- L-arabinose binds to AraC and causes the protein to release the O₂ site and bind the I₂ site, which is adjacent to the I₁ site. This releases the DNA loop and allows transcription to begin.
- The cAMP activator protein (CAP)-cAMP complex binds to the DNA and stimulates binding of AraC to I₁ and I₂.

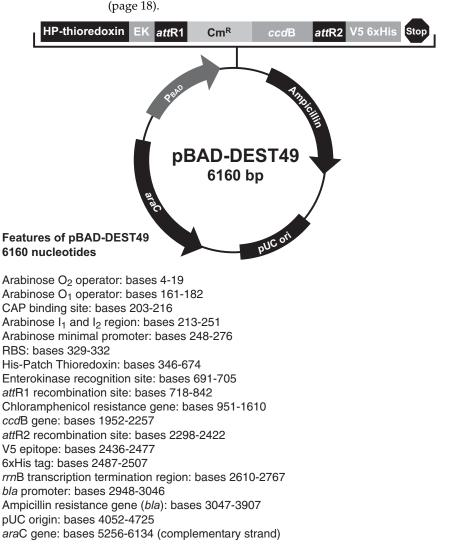


Glucose Repression

Glucose represses basal expression levels by lowering cAMP levels which in turn decrease the binding of CAP. As cAMP levels are lowered, transcriptional activation is decreased. Add glucose to the culture medium to repress basal expression levels.

Map of pBAD-DEST49

Map of pBAD-
DEST49The map below shows the elements of pBAD-DEST49. DNA
from the entry clone replaces the region between bases 725
and 2408. The complete sequence of pBAD-DEST49 is
available for downloading from our Web site
(www.invitrogen.com) or by contacting Technical Support



Features of pBAD-DEST49

Features of
pBAD-DEST49pBAD-DEST49 (6160 bp) contains the following elements.Features have been functionally tested.

Feature	Benefit
araBAD promoter (PBAD)	Provides tight, dose-dependent regulation of heterologous gene expression (Guzman <i>et al.</i> , 1995)
O ₂ region	Binding site of AraC that represses transcription from PBAD
O1 region	Binding site of AraC that represses transcription of the <i>ara</i> C promoter (PC) (transcribed on the opposite strand)
CAP binding site	Site where CAP (cAMP binding protein) binds to activate transcription from $PBAD$ and PC
I_2 and I_1 regions	Binding sites of AraC that activate transcription from PBAD
Ribosome binding site	Increases efficiency of recombinant fusion protein expression
HP-thioredoxin	Provides a highly efficient fusion partner for translation of the fusion protein
Enterokinase recognition site	Allows removal of the N-terminal tag from the recombinant fusion protein
attR1 and attR2 sites	Allows recombinational cloning of the gene of interest from an entry clone
Chloramphenicol resistance gene	Allows counterselection of expression clones
ccdB gene	Allows negative selection of expression clones
V5 epitope	Allows detection of the recombinant protein by the Anti-V5 antibodies (Southern <i>et al.</i> , 1991)
C-terminal polyhistidine tag	Allows purification of the recombinant protein on metal-chelating resin such as ProBond [™]
	Allows detection of the recombinant protein by the Anti-His (C-term) antibodies (Lindner <i>et al.</i> , 1997)
<i>rrn</i> B transcription termination region	Strong transcription termination region
Ampicillin resistance gene	Allows selection of the plasmid in E. coli
pUC origin (see note on page 1)	Allows low-copy replication and growth in <i>E. coli</i>
araC gene	Encodes the regulatory protein for tight regulation of the PBAD promoter (Lee, 1980; Schleif, 1992)

Technical Support

Web Resources			the Invitrogen website at <u>www</u> Fechnical resources including					
		 Technical resources including manuals, vector maps and sequences, application notes, MSDSs, FAQs, formulations, citations, handbooks, etc. 						
		Complete technical support contact information.						
		• A	Access to the Invitrogen Onlin	e Catalog.				
		• A	Additional product informatio	on and special offers.				
	Contact Us		nore information or technical a ail. Additional international c					
				sinces are instea on our				
			ite (<u>www.invitrogen.com</u>).					
	Corporate Headquarter		ite (<u>www.invitrogen.com</u>). Japanese Headquarters:	European Headquarters:				
	Corporate Headquarter Invitrogen Corporation 5791 Van Allen Way Carlsbad, CA 92008 USA	S	Japanese Headquarters: Invitrogen Japan LOOP-X Bldg. 6F	Invitrogen Ltd Inchinnan Business Park 3 Fountain Drive				
	Invitrogen Corporation 5791 Van Allen Way	S	Japanese Headquarters: Invitrogen Japan LOOP-X Bldg. 6F 3-9-15, Kaigan	Invitrogen Ltd Inchinnan Business Park 3 Fountain Drive Paisley PA4 9RF, UK				
	Invitrogen Corporation 5791 Van Allen Way Carlsbad, CA 92008 USA	s	Japanese Headquarters: Invitrogen Japan LOOP-X Bldg. 6F	Invitrogen Ltd Inchinnan Business Park 3 Fountain Drive Paisley PA4 9RF, UK Tel: +44 (0) 141 814 6100				
	Invitrogen Corporation 5791 Van Allen Way Carlsbad, CA 92008 USA Tel: 1 760 603 7200	s	Japanese Headquarters: Invitrogen Japan LOOP-X Bldg. 6F 3-9-15, Kaigan Minato-ku, Tokyo 108-0022	Invitrogen Ltd Inchinnan Business Park 3 Fountain Drive Paisley PA4 9RF, UK				

tech_support@invitrogen.com

MSDS	MSDSs (Material Safety Data Sheets) are available on our website at <u>www.invitrogen.com/msds.</u>
Certificate of Analysis	The Certificate of Analysis (CofA) provides detailed quality control information for each product. CofAs are available on our website at <u>www.invitrogen.com/support</u> , and are searchable by product lot number, which is printed on each box.

Technical Support, Continued

Invitrogen is committed to providing our customers with Limited Warranty 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 Support Representatives. Invitrogen warrants that all of its products will perform according to the specifications stated on the certificate of analysis. The company will replace, free of charge, any product that does not meet those specifications. This warranty limits Invitrogen Corporation's liability only to the cost of the product. No warranty is granted for products beyond their listed expiration date. No warranty is applicable unless all product components are stored in accordance with instructions. Invitrogen reserves the right to select the method(s) used to analyze a product unless Invitrogen agrees to a specified method in writing prior to acceptance of the order. Invitrogen makes every effort to ensure the accuracy of its publications, but realizes that the occasional typographical or other error is inevitable. Therefore Invitrogen makes no warranty of any kind regarding the contents of any publications or documentation. If you discover an error in any of our publications, report it to our Technical Support Representatives. Invitrogen assumes no responsibility or liability for any

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

Purchaser Notification

Limited Use Label License No. 19: Gateway[®] Cloning Products

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 purchase of this product does not convey a license under any method claims in the foregoing patents or patent applications, or to use this product with any recombination sites other than those purchased from Invitrogen Corporation or its authorized distributor. The right to use methods claimed in the foregoing patents or patent applications with this product for research purposes only can only be acquired by the use of Clonase[™] purchased from Invitrogen Corporation or its authorized distributors. The buyer cannot modify the recombination sequence(s) contained in this product for any purpose. The buyer cannot sell or otherwise transfer (a) this product, (b) its components, or (c) materials made by the employment of this product or its components to a third party or otherwise use this product or its components or materials made by the employment of this product or its components for Commercial Purposes. The buyer may transfer information or materials made through the employment 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. Notwithstanding the preceding, any buyer who is employed in an academic or government institution may transfer materials made with this product to a third party who has a license from Invitrogen under the patents identified above to distribute such materials. Transfer of such materials and/or information to collaborators does not convey rights to practice any methods claimed in the foregoing patents or patent applications. 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. Invitrogen Corporation will not assert a claim against the buyer of infringement of the above patents 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 none of (i) this product, (ii) any of its components, or (iii) a method claim of the foregoing patents, was used in the manufacture of such product. Invitrogen Corporation will not assert a claim against the buyer of infringement of the above patents based upon the use of this product to manufacture a protein for sale, provided that no method claim in the above patents was used in the manufacture of such protein. If the purchaser is not willing to accept the limitations of this limited use statement, Invitrogen is willing to accept return of the product with a full refund. For information on purchasing a license to use this product for purposes other than those permitted above, contact Licensing Department, Invitrogen Corporation, 5791 Van Allen Way, Carlsbad, California 92008. Phone (760) 603-7200.

Purchaser Notification, Continued

Limited Use Label License No. 22: Vectors and Clones Encoding Histidine Hexamer	This product is licensed under U.S. and foreign patents 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.
Limited Use	The ThioFusion TM Expression System is licensed under U.S.

Label License No. 29: ThioFusion[™] Expression System The ThioFusionTM Expression System is licensed under U.S. patents from Genetics Institute, Inc. for research use only. Licenses for commercial manufacture or use may be obtained directly from Genetics Institute, Inc., 87 Cambridgepark Drive, Cambridge, MA 02140.

References

Carson, M. J., Barondess, J. J., and Beckwith, J. (1991). The FtsQ Protein of *Escherichia coli*: Membrane Topology, Abundance, and Cell Division Phenotypes Due to Overproduction and Insertion Mutations. J. Bacteriol. *173*, 2187-2195.

Dalbey, R. E., and Wickner, W. (1985). Leader Peptidase Catalyzes the Release of Exported Proteins from the Outer Surface of the *Escherichia coli* Plasma Membrane. J. Biol. Chem. *260*, 15925-15931.

Guzman, L.-M., Barondess, J. J., and Beckwith, J. (1992). FtsL, an Essential Cytoplasmic Membrane Protein Involved in Cell Division in *Escherichia coli*. J. Bacteriol. *174*, 7716-7728.

Holmgren, A. (1985). Thioredoxin. Ann. Rev. Biochem. 54, 237-271.

Katti, S. K., LeMaster, D. M., and Eklund, H. (1990). Crystal Structure of Thioredoxin from *E. coli* at 1.68 Angstroms Resolution. J. Mol. Biol. 212, 167-184.

Kuhn, A., and Wickner, W. (1985). Isolation of Mutants in M13 Coat Protein That Affect its Synthesis, Processing and Assembly into Phage. J. Biol. Chem. 260, 15907-15913.

Landy, A. (1989). Dynamic, Structural, and Regulatory Aspects of Lambda Sitespecific Recombination. Annu. Rev. Biochem. 58, 913-949.

LaVallie, E. R., DiBlasio, E. A., Kovacic, S., Grant, K. L., Schendel, P. F., and McCoy, J. M. (1993). A Thioredoxin Gene Fusion Expression System That Circumvents Inclusion Body Formation in the *E. coli* Cytoplasm. Bio/Technology *11*, 187-193.

Lee, N. (1980) Molecular Aspects of *ara* Regulation. In *The Operon*, J. H. Miller and W. S. Reznikoff, eds. (Cold Spring Harbor, N.Y.: Cold Spring Harbor Laboratory), pp. 389-410.

Lee, N., Francklyn, C., and Hamilton, E. P. (1987). Arabinose-Induced Binding of AraC Protein to *araI*₂ Activates the *araBAD* Operon Promoter. Proc. Natl. Acad. Sci. USA *84*, 8814-8818.

Lindner, P., Bauer, K., Krebber, A., Nieba, L., Kremmer, E., Krebber, C., Honegger, A., Klinger, B., Mocikat, R., and Pluckthun, A. (1997). Specific Detection of His-tagged Proteins With Recombinant Anti-His Tag scFv-Phosphatase or scFv-Phage Fusions. BioTechniques 22, 140-149. Lu, Z., DiBlasio-Smith, E. A., Grant, K. L., Warne, N. W., LaVallie, E. R., Collins-Racie, L. A., Follettie, M. T., Williamson, M. J., and McCoy, J. M. (1996). Histidine Patch Thioredoxins. J. Biol. Chem. 271, 5059-5065.

Miyada, C. G., Stoltzfus, L., and Wilcox, G. (1984). Regulation of the *ara*C Gene of *Escherichia coli*: Catabolite Repression, Autoregulation, and Effect on *ara*BAD Expression. Proc. Natl. Acad. Sci. USA *81*, 4120-4124.

Ogden, S., Haggerty, D., Stoner, C. M., Kolodrubetz, D., and Schleif, R. (1980). The *Escherichia coli* L-Arabinose Operon: Binding Sites of the Regulatory Proteins and a Mechanism of Positive and Negative Regulation. Proc. Natl. Acad. Sci. USA 77, 3346-3350.

Russell, C. B., Stewart, R. C., and Dahlquist, F. W. (1989). Control of Transducer Methylation Levels in *Escherichia coli*: Investigation of Components Essential for Modulation of Methylation and Demethylation Reactions. J. Bacteriol. *171*, 3609-3618.

San Millan, J. L., Boyd, D., Dalbey, R., Wickner, W., and Beckwith, J. (1989). Use of *phoA* Fusions to Study the Topology of the *Escherichia coli* Inner Membrane Protein Leader Peptidase. J. Bacteriol. *171*, 5536-5541.

Schleif, R. S. (1992). DNA Looping. Ann. Rev. Biochem. 61, 199-223.

Southern, J. A., Young, D. F., Heaney, F., Baumgartner, W., and Randall, R. E. (1991). Identification of an Epitope on the P and V Proteins of Simian Virus 5 That Distinguishes Between Two Isolates with Different Biological Characteristics. J. Gen. Virol. 72, 1551-1557.

Wallace, B. J., and Kushner, S. R. (1984). Genetic and Physical Analysis of the Thioredoxin (*trxA*) Gene of *Escherichia coli* K-12. Gene 32, 399-408.

^{© 2002–2008, 2010} Invitrogen Corporation. All rights reserved.

invitrogen

Corporate Headquarters Invitrogen Corporation 5791 Van Allen Way Carlsbad, CA 92008 T: 1 760 603 7200 F: 1 760 602 6500 E: tech_support@invitrogen.com

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