pEF-DEST51 Gateway[™] Vector

A destination vector for cloning and expression of C-terminal fusion proteins in mammalian cells

Catalog number 12285011

Publication Number MAN0000240 Revision A.0



For Research Use Only. Not for use in diagnostic procedures.

Information in this document is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Revision history: MAN0000240

Revision	Date	Description
A.0	12 April 2017	Corrected pEF-DEST51 sequence and related maps, updated branding, legal/regulatory language
1.0	7 July 2010	Basis for this revision

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.



Manufacturer: Life Technologies Corporation | 5791 Van Allen Way | Carlsbad, CA 92008

Trademarks: All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

© 2017 Thermo Fisher Scientific Inc. All rights reserved.

Contents

Kit contents and storage	iv
Methods	1
Product overview	1
Using pEF-DEST51	
Transfection	6
Expression and analysis	8
Create stable cell lines	
Appendix	11
Recipes	
Blasticidin	
Human EF-1 α promoter	14
Map and features of pEF-DEST51	
Map of pEF/GW-51/lacZ	
Accessory products	
Documentation and support	20
Obtaining support	
References	21

Kit contents and storage

Shipping and storage	pEF-DEST51 and pEF/GW-51/ <i>lacZ</i> are shipped at room temperature. Upon receipt, store at –20°C. Products are guaranteed for six months from date of shipment when
	stored properly.

Contents

The pEF-DEST51 Gateway ${}^{\scriptscriptstyle\rm TM}$ Vector components are listed below.

Item	Concentration	Volume
pEF-DEST51 Vector	150 ng/µL in 10 mM Tris-HCl, 1 mM EDTA, pH 8.0	40 µL
pEF/GW-51/ <i>lacZ</i> Control Plasmid	0.5 μg/μL in 10 mM Tris-HCl, 1 mM EDTA, pH 8.0	20 µL

Methods

Product overview

Description	pEF-DEST51 is a 7.5 kb vector derived from pEF6/V5-His and adapted for use with the Gateway [™] Technology. It is designed to allow high-level, constitutive expression in a variety of mammalian hosts.
Features	pEF-DEST51 contains the following elements:Human elongation factor 1α-subunit promoter (hEF-1α)
	for high-level expression across a broad range of species and cell types (page 14)
	 Two recombination sites, <i>att</i>R1 and <i>att</i>R2, downstream of the EF-1α promoter 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 6× His tag for detection and purification (optional)
	• Bovine growth hormone (BGH) polyadenylation sequence for proper termination and processing of the recombinant transcript
	• f1 intergenic region for production of single-strand DNA in F plasmid-containing <i>E. coli</i>
	 SV40 early promoter and origin for expression of the blasticidin resistance gene and stable propagation of the plasmid in hosts expressing the SV40 large T antigen
	• EM7 promoter for expression of the blasticidin resistance gene in <i>E. coli</i>
	• Blasticidin resistance gene for selection of stable cell lines
	• The pUC origin for high copy replication and maintenance of the plasmid in <i>E. coli</i>
	• The ampicillin (<i>bla</i>) resistance gene for selection in <i>E. coli</i>
	For a map of pEF-DEST51, see page 15.

Product 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 [™] cloning technology, simply:
	 Clone your gene of interest into a Gateway[™] entry vector to create an entry clone.
	 Generate an expression clone by performing an LR recombination reaction between the entry clone and a Gateway[™] destination vector (<i>e.g.</i> pEF-DEST51).
	3. Transfect your expression clone into the cell line of choice for transient or constitutive expression of your gene of interest.
	For more information on the Gateway [™] System, refer to the Gateway [™] Technology User Guide. This user guide is available for download from www.thermofisher.com or by contacting Technical Support (page 20).

Using pEF-DEST51

Q Important	The pEF-DEST51 vector is supplied as a supercoiled plasmid. Although we have 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.
Propagate pEF-DEST51	If you wish to propagate and maintain pEF-DEST51, we recommend using Library Efficiency TM DB3.1 TM Competent Cells available for transformation (see page 18 for ordering). The DB3.1 TM <i>E. coli</i> strain is resistant to CcdB effects and can support the propagation of plasmids containing the <i>ccd</i> B gene.
	Note: DO NOT use general <i>E. coli</i> cloning strains including TOP10 or $DH5\alpha^{TM}$ for propagation and maintenance as these strains are sensitive to CcdB effects.
Entry clone	To recombine your gene of interest into pEF-DEST51, you should have an entry clone containing your gene of interest. For your convenience, we offer the pENTR [™] Directional TOPO [™] Cloning Kit (see page 11) for 5-minute cloning of your gene of interest into an entry vector. For more information on entry vectors available, go to www.thermofisher.com or contact Technical Support (page 20).
	For detailed information on constructing an entry clone, refer to the specific entry vector manual. For detailed information on performing the LR recombination reaction, refer to the Gateway [™] Technology user guide.

Using pEF-DEST51, continued

Points to consider before recombining	Your insert should contain a Kozak consensus sequence with an ATG initiation codon for proper initiation of translation (Kozak, 1987; Kozak, 1991; Kozak, 1990). An example of a Kozak consensus sequence is provided below. Other sequences are possible, but the G or A at position –3 and the G at position +4 are the most critical for function (shown in bold). The ATG initiation codon is shown underlined.
	(G/A)NNATGG
	If you wish to include the V5 epitope and 6x His tag, your gene in the entry clone should not contain a stop codon. In addition, the gene should be designed to be in frame with the C-terminal epitope tag after recombination. Refer to the Recombination Region on page 5.
	If you do NOT wish to include the V5 epitope and 6x His tag, your gene should contain a stop codon in the entry clone.
Recombine 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 ^{TM} LR Clonase ^{TM} enzyme mix (see page 20 for ordering). 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 chloramphenicol (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 TM Technology user guide to set up the LR Clonase TM reaction, transform <i>E. coli</i> , and select for the expression clone.
Note	The presence of the EM7 promoter and the blasticidin resistance gene allows for selection of <i>E. coli</i> transformants using blasticidin. For selection, use Low Salt LB agar plates containing 100 μ g/mL blasticidin (see page 11 for a recipe). For blasticidin to be active, the salt concentration of the

medium must remain low (< 90 mM) and the pH must be 7.0.

Using pEF-DEST51, continued

Confirm the expression clone			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 µg/mL chloramphenicol. A true expression clone will not grow in the presence of chloramphenicol.													
Reco regio	ombi on	nati	on	The recombination region of the expression clone resulting from pEF-DEST51 × entry clone is shown below.												
				Fea •	tures Shac trans recor pEF- The : corre	s of t led r sferre mbin DES nucle espoi	the re egion ed fro ation T51 v eotid nd to	ns con om th n. Nc vecto es or base	bina rresp ne en on-sh r. n eith es 172	tion ond try c aded er sid 25 an	regia to the lone l regi de of ad 34	on: iose l into ions a the 08, re	DNA pEF- are d shad	sequ DES erive ed re tivel	ience T51 k ed fro egion y, of	es by om the the
					рег-	DES	151 1	vecto	r seq	uenc	ce.					
1639	TCAG AGTC	GTGT GTGT CACA	CG T GC T	GAGO	GAATI CTTA <i>F</i>	TA GO	CTTGO	GTACI CATGA	AAT TTA	ACGA	ACTC TGAG	ACTA TGAI	TAGO ATCC	GA G CT C	GACCC CTGGG	AAGCT
							1	725								
1699	GGCI	AGGI TCC	AA G TT C	GAAC	GATCA CTAGI	A CA T GI	AGTI TCAF	<u>T</u> GTA ACAI	A CAA CAA	AAAA TTTT	AGCA CGT	GGCI	'N \N	G	ENE	NAC
			24	00					<i>att</i> B1							
	pro	ala	phe	leu	tyr	lys	val	val	asp	leu	glu	gly	pro	arg	phe	glu
3400	CCA	GCT	TT <u>C</u>	TTG	TAC	AAA	GTG	GTT	GAT	CTA	GAG	GGC	CCG	CGG	TTC	GAA
	GGT	CGA	AAG	attB2	ATG	111	CAC	CAA	CTA	GAT	CTC	CCG	GGC	GCC	AAG	CIT
	~1	1		+10			V5 e	pitope	1.011	~1	1.011			+ h m		+ h m
3448	GGT	AAG	CCT	ATC	CCT	AAC	CCT	CTC	CTC	GGT GGT	CTC	GAT	TCT	ACG	CGT	ACC
	CCA	TTC	GGA	TAG	GGA	TTG	GGA	GAG	GAG	CCA	GAG	CTA	AGA	TGC	GCA	TGG
				6xH	is tag											
	gly	his	his	his	his	his	his	***								
3496	GGT CCA	CAT GTA	CAT GTA	CAC GTG	CAT GTA	CAC GTG	CAT GTA	TGA ACT	GTT1 CAAZ	AAA(דידידי	CC C(36 G(GAC	ATCAC	G CCI	CGAC	CTGT GACT
	0.011	~ * * *	- + + I	010	- 111	010	- + + 1		~~ 14 14					. 501	-0010	

Transfection

Introduction	This section provides general information for transfecting your expression clone into the mammalian cell line of choice. We recommend that you include the positive control vector pEF/GW-51/ <i>lacZ</i> and a mock transfection (negative control) in your experiments to evaluate your results.
Plasmid preparation	Once you have generated your expression clone, you must isolate plasmid DNA for transfection. Plasmid DNA for transfection into eukaryotic cells must be very clean and free from phenol and sodium chloride. Contaminants will kill the cells, and salt will interfere with lipid complexing, decreasing transfection efficiency. We recommend isolating plasmid DNA using the PureLink [™] HiPure MiniPrep Kit (up to 30 µg DNA), the PureLink [™] HiPure MidiPrep Kit (up to 150 µg DNA) (see page 18), or CsCl gradient centrifugation.
Methods of transfection	For established cell lines (<i>e.g.</i> HeLa), consult original references or the supplier of your cell line for the optimal method of transfection. We recommend that you follow exactly the protocol for your cell line. Pay particular attention to medium requirements, when to pass the cells, and at what dilution to split the cells. Further information is provided in <i>Current Protocols in Molecular Biology</i> (Ausubel <i>et al.</i> , 1994).
	Methods for transfection include calcium phosphate (Chen and Okayama, 1987; Wigler <i>et al.</i> , 1977), lipid-mediated (Felgner <i>et al.</i> , 1989; Felgner and Ringold, 1989) and electroporation (Chu <i>et al.</i> , 1987; Shigekawa and Dower, 1988) For cationic lipid-based transfection, we recommend using Lipofectamine [™] 2000 Reagent. For electroporation of primary, stem cell and difficult-to-transfect cells, we recommend the Neon [™] Transfection System (see page 18). For more information, go to www.thermofisher.com or contact Technical Support (page 20).
	Continued on next page

Transfection, continued

To propagate and maintain the plasmid:

- Use the supplied 0.5 µg/µL stock solution in TE buffer, pH 8.0 to transform a *recA*, *endA E*. *coli* strain like TOP10, DH5α[™], JM109, or equivalent.
- Select transformants on LB agar plates containing 50–100 µg/mL ampicillin.
- 3. Prepare a glycerol stock of a transformant containing plasmid for long-term storage.

Expression and analysis

Introduction	Expression of your gene of interest from the expression clone can be performed in either transiently transfected cells or stable cell lines (see page 10 for guidelines to create stable cell lines). A sample protocol is provided below. Other protocols are suitable.
Prepare cell lysates	 To lyse cells: Wash cell monolayers (~5 × 10⁵ to 1 × 10⁶ cells) once with phosphate-buffered saline (PBS, Cat. No. 10010023). Scrape cells into 1 mL PBS and pellet the cells at 1500 × <i>g</i> for 5 minutes. Resuspend in 50 μL Cell Lysis Buffer (see page 12 for a recipe). Other cell lysis buffers are suitable. Vortex. Incubate cell suspension at 37°C for 10 minutes to lyse the cells.
	 Note: You may prefer to lyse the cells at room temperature or on ice if degradation of your protein is a potential problem. 5. Centrifuge the cell lysate at 10,000 × <i>g</i> for 10 minutes at 4°C to pellet nuclei and transfer the supernatant to a fresh tube. Assay the lysate for protein concentration. Note: Do not use protein assays utilizing Coomassie Blue or other dyes. NP-40 interferes with the binding of the dye with the protein. 6. Add SDS-PAGE sample buffer (see page 12 for a recipe) to a final concentration of 1X and boil the sample for 5 minutes. 7. Load 20 µg of lysate onto an SDS-PAGE gel and electrophorese. Use the appropriate percentage of acrylamide to resolve your fusion protein.
Polyacrylamide gel electrophoresis	To facilitate separation and visualization of your recombinant fusion protein by polyacrylamide gel electrophoresis, a wide range of pre-cast NuPAGE [™] and Novex [™] Tris-Glycine polyacrylamide gels and electrophoresis apparatus are available for purchase. For more information, go to www.thermofisher.com or contact Technical Support (page 20)

Expression and analysis, continued

Detect recombinant fusion proteins	To detect expression of your recombinant fusion protein by Western blot analysis, you can use the Anti-V5 antibodies or the Anti-His(C-term) antibodies (see page 18) or an antibody to your protein of interest. In addition, the Positope [™] Control Protein is available for use as a positive control for detection of fusion proteins containing a V5 epitope or a polyhistidine (6x His) tag.
	For more information, go to www.thermofisher.com or contact Technical Support (page 20).
Assay for β-galactosidase	If you use the pEF/GW-51/ <i>lac</i> Z plasmid as a positive control vector, you may assay for β -galactosidase expression by Western blot analysis or activity assay using cell lysates (Miller, 1972). β -Gal Antiserum, the β -Gal Assay Kit, and the β -Gal Staining Kit are available for fast and easy detection of β -galactosidase expression (see page 18).
Note	The C-terminal peptide containing the V5 epitope and the polyhistidine region will add approximately 5 kDa to your protein.
Purify recombinant fusion proteins	The presence of the C-terminal polyhistidine (6x His) tag in your recombinant fusion protein allows use of a metal- chelating resin such as ProBond [™] to purify your fusion protein. The ProBond [™] Purification System and bulk ProBond [™] resin are available (see page 19 for ordering information). Refer to the ProBond [™] Purification System manual for protocols to purify your fusion protein. The Ni-NTA Agarose is available for purification of proteins containing a polyhistidine (6x His) tag (see page 18).
	Note: Other metal-chelating resins and purification methods are suitable.

Create stable cell lines

Introduction

The pEF-DEST51 vector contains the blasticidin resistance gene to allow selection of stable cell lines. To create stable cell lines, transfect your construct into the mammalian cell line of choice and select for foci using blasticidin. General guidelines are provided below.



To obtain stable transfectants, we recommend that you linearize your pEF-DEST51 construct before transfection. While linearizing the vector may not improve the efficiency of transfection, it increases the chances that the vector does not integrate in a way that disrupts elements necessary for expression in mammalian cells. To linearize your construct, cut at a unique site that is neither located within a critical element nor within your gene of interest.

Blasticidin	Blasticidin S HCl is a nucleoside antibiotic isolated from <i>Streptomyces griseochromogenes</i> which inhibits protein synthesis in both prokaryotic and eukaryotic cells (Takeuchi <i>et al.</i> , 1958; Yamaguchi <i>et al.</i> , 1965). Resistance is conferred by expression of either one of two blasticidin S deaminase genes: <i>bsd</i> from <i>Aspergillus terreus</i> (Kimura <i>et al.</i> , 1994) or <i>bsr</i> from <i>Bacillus cereus</i> (Izumi <i>et al.</i> , 1991). These deaminases convert blasticidin S to a non-toxic deaminohydroxy derivative (Izumi <i>et al.</i> , 1991).
Disstisidin	Blacticidin is available separately (see page 18 for ordering)

Blasticidin selection quidelines Blasticidin is available separately (see page 18 for ordering). Use as follows:

- 1. Prepare blasticidin in sterile water and filter-sterilize the solution.
- 2. Use 2.5–10 μ g/mL of blasticidin in complete medium.
- 3. Test varying concentrations of blasticidin on your cell line to determine the concentration that kills your cells (kill curve). Cells differ in their susceptibility to blasticidin. Complete selection can take up to 10 days of growth in selective medium.

See page 13 for instructions on how to prepare and store blasticidin.

Appendix

Recipes

LB (Luria- Bertani) medium and plates	Composition: 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.				
	 Autoclave on liquid cycle for 20 minutes at 15 psi. Allow solution to cool to 55°C and add antibiotic (100 μg/mL ampicillin) if needed. 				
	4. Store at room temperature or at 4°C.				
	LB agar plates1. Prepare LB medium as above, but add 15 g/L agar before autoclaving.				
	2. Autoclave on liquid cycle for 20 minutes at 15 psi.				
	 After autoclaving, cool to ~55°C, add antibiotic (100 μg/mL of ampicillin), and pour into 10 cm plates. 				
	4. Let harden, then invert and store at 4°C.				
Low salt LB medium with Blasticidin	Low Salt LB Medium: 10 g Tryptone 5 g NaCl 5 g Yeast Extract				
	 Combine the dry reagents above and add deionized, distilled water to 950 mL. Adjust pH to 7.0 with 1 N NaOH. Bring the volume up to 1 liter. For plates, add 15 g/L agar before autoclaving. 				
	 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 blasticidin to $100 \ \mu g/mL$ final concentration.				
	 Store plates at 4°C in the dark. Plates containing blasticidin are stable for up to 2 weeks. 				

Recipes, continued

Cell lysis buffer	50 mM Tris, pH 7.8 150 mM NaCl 1% Nonidet P-40				
	1.	This solution can be prepared from the following common stock solutions. For 100 mL, combine			
		1 M Tris base5 mL5 M NaCl3 mLNonidet P-401 mL			
	2.	 Bring the volume up to 90 mL with deionized water and adjust the pH to 7.8 with HCl. 			
	3.	Bring the volume up to 100 mL. Store at room temperature.			
	Tc leı	prevent proteolysis, you may add 1 mM PMSF, 1 μM upeptin, or 0.1 μM aprotinin before use.			
4X SDS-PAGE sample buffer	1.	Combine the following reagents: 0.5 M Tris-HCl, pH 6.85 mLGlycerol (100%)4 mL β -mercaptoethanol0.8 mLBromophenol Blue0.04 gSDS0.8 g			
	2.	Bring the volume to 10 mL with sterile water.			
	3.	Aliquot and freeze at -20° C until needed.			

Blasticidin

Molecular weight, formula, and structure The formula for blasticidin S is $C_{17}H_{26}N_8O_5\text{-}HCl$, and the molecular weight is 458.9. The diagram below shows the structure of blasticidin.



How to handle Blasticidin	Always wear gloves, mask, goggles, and protective clothing (<i>e.g.</i> a laboratory coat) when handling blasticidin. Weigh out blasticidin and prepare solutions in a hood.
Prepare and store stock	Blasticidin may be obtained separately in 50 mg aliquots. Blasticidin is soluble in water. Sterile water is generally used to prepare stock solutions of 5 to 10 mg/mL.
Socutions	• Dissolve blasticidin in sterile water and filter-sterilize the solution.
	• Aliquot in small volumes suitable for one time use and freeze at -20°C for long-term storage or store at 4°C for short-term storage.
	• Aqueous stock solutions are stable for 1–2 weeks at 4°C and 6–8 weeks at –20°C.
	• pH of the aqueous solution should be 7.0 to prevent inactivation of blasticidin.
	• Do not subject stock solutions to freeze/thaw cycles (do not store in a frost-free freezer).
	• Upon thawing, use what you need and store the thawed stock solution at 4°C for up to 2 weeks.
	• Medium containing blasticidin may be stored at 4°C for up to 2 weeks.

Human EF-1 α promoter

Descripti	on Ti pi 19	he diagram romoter use 990; Uetsuki	below show d in pEF vec <i>et al.,</i> 1989).	s all the feat ctors (Mizus)	ures of the H hima and Na	EF-1α agata,
	_	— 5' end of hum	an EE-1a promo	for		
459	 AAGGAGTGGG	AATTGGCTCC	GGTGCCCGTC	AGTGGGCAGA	GCGCACATCG	CCCACAGTCC
519	CCGAGAAGTT	GGGGGGAGGG	GTCGGCAATT	GAACCGGTGC	CTAGAGAAGG	TGGCGCGGGG
579	TAAACTGGGA TATA box	AAGTGATGTC	GTGTACTGGC	TCCGCCTTTT Start of Tra		GGGGGAGAAC
639	CGTATATAAG	TGCAGTAGTC	GCCGTGAACG	I TTCTTTTTCG	CAACGGGTTT	GCCGCCAGAA
	5´	end of Intron 1				Exon
699	CACAGGTAAG	TGCCGTGTGT	GGTTCCCGCG	GGCCTGGCCT	CTTTACGGGT	TATGGCCCTT
759	GCGTGCCTTG	AATTACTTCC	ACCTGGCTGC	AGTACGTGAT	TCTTGATCCC	GAGCTTCGGG
819	TTGGAAGTGG	GTGGGAGAGT	TCGAGGCCTT	GCGCTTAAGG	AGCCCCTTCG	CCTCGTGCTT
879	GAGTTGAGGC	CTGGCCTGGG	CGCTGGGGGCC	GCCGCGTGCG	AATCTGGTGG	CACCTTCGCG
939	CCTGTCTCGC	TGCTTTCGAT	AAGTCTCTAG	CCATTTAAAA	TTTTTGATGA	CCTGCTGCGA
999	CGCTTTTTTT	CTGGCAAGAT	AGTCTTGTAA	ATGCGGGCCA	AGATCTGCAC	ACTGGTATTT
1059	CGGTTTTTGG	Gecc <u>(ceeec</u>	<u>GGCGA</u> CGGGG	CCCGTGCGTC	CCAGCGCACA	tgttcggd <u>GA</u>
1119	<u>Sp 1</u> <u>GGCGGGG</u> CCT	GCGAGCGCGG	CCACCGAGAA	TCGGACGGGG	GTAGTCTCAA	GCTGGCCGGC
1179	CTGCTCTGGT	GCCTGGCCTC	GCGCCGCCGT	GTATCGCCCC	GCCCTGGGCG	GCAAGGCTGG
1239	CCCGGTCGGC	ACCAGTTGCG	TGAGCGGAAA	GATGGCCGCT	TCCCGGCCCT	GCTGCAGGGA
1299	GCTCAAAATG	GAGGACGCGG	CGCTCGGGAG	Sp 1 AQCGGGCGGG	TGAGTCACCC	ACACAAAGGA
1359	AAAGGGCCTT	TCCGTCCTCA	GCCGTCGCTT	Ap 1 CATG <u>TGACTC</u>	CACGGAGTAC	CGGGCGCCGT
1419	CCAGGCACCT	CGATTAGTTC	TCGAGCTTTT	GGAGTACGTC	GTCTTTAGGT	TGGGGGGAGG
1479	GGTTTTATGC	GATGGAGTTT	CCCCACACTG	AGTGGGTGGA	GACTGAAGTT	AGGCCAGCTT
1539	GGCACTTGAT	GTAATTCTCC	TTGGAATTTG	CCCTTTTTGA	GTTTGGATCT	TGGTTCATTC
				s end of intro		
1599	TCAAGCCTCA	GACAGTGGTT	CAAAGTTTTT	TTCTTCCATT	TCAGGTGTCG	TGA
					5' end of	Exon 2

Map and features of pEF-DEST51

Map of pEF-DEST51 The map below shows the elements of pEF-DEST51. DNA from the entry clone replaces the region between bases 1725 and 3408. The nucleotide sequence of pEF-DEST51 is available from **www.thermofisher.com** or by contacting Technical Support (page 20).



(c) = complementary strand

Map and features, continued

Features of
pEF-DEST51pEF-DEST51 (7450 bp) contains the following elements. All
features have been functionally tested.

Feature	Benefit
Human elongation factor 1 \Box (hEF-1 α) promoter	Allows expression of recombinant proteins in a broad range of mammalian cell types (Goldman <i>et al.,</i> 1996; Mizushima and Nagata, 1990)
T7 promoter	Allows <i>in vitro</i> transcription in the sense orientation
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 recombinant fusion proteins by the Anti-V5 antibodies (Southern <i>et al.</i> , 1991)
C-terminal polyhistidine tag	Allows purification of recombinant proteins 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)
Bovine growth hormone (BGH) polyadenylation signal	Allows efficient transcription termination and polyadenylation of mRNA (Goodwin and Rottman, 1992)
f1 origin	Allows rescue of single-stranded DNA
SV40 early promoter and origin	Allows high-level expression of the blasticidin resistance gene and episomal replication in cells expressing the SV40 large T antigen
EM7 promoter	Allows expression of the blasticidin resistance gene in <i>E. coli</i>
Blasticidin resistance gene	Allows selection of stable transfectants in mammalian cells (Kimura <i>et al.</i> , 1994)
SV40 early polyadenylation signal	Allows efficient transcription termination and polyadenylation of mRNA
pUC origin	Allows high-copy number replication and growth in <i>E. coli</i>
Ampicillin resistance gene	Allows selection of transformants in <i>E. coli</i>

Map of pEF/GW-51/lacZ

Description

pEF/GW-51/*lacZ* is an 8904 bp control vector containing the gene for β -galactosidase. pEF/GW-51/*lacZ* was constructed using the GatewayTM LR recombination reaction between an entry clone containing the *lacZ* gene and pEF-DEST51. β -galactosidase is expressed as a fusion to the C-terminal tag. The fusion protein is approximately 121 kDa in size.

Map of pEF/ GW-51/*lac*Z

The map below shows the elements of pEF/GW-51/*lacZ*. The nucleotide sequence of pEF/GW-51/*lacZ* is available from **www.thermofisher.com** or by contacting Technical Support (page 20).

lacZ V5 6xHis Stop BGH DA pEF/GW-51/*lac*Z 8904 bp Features of pEF/GW-51/lacZ 8904 nucleotides EF-1a promoter: bases 470-1653 T7 promoter: bases 1670-1689 attB1 recombination site: bases 1720-1744 PUC SV40 pA lacZ ORF: bases 1764-4823 attB2 recombination site: bases 4840-4864 V5 epitope: bases 4890-4931 6xHis tag: bases 4941-4958 BGH polyadenylation region: bases 4984-5211 f1 origin: bases 5257-5685 SV40 early promoter and origin: bases 5690-6034 EM7 promoter: bases 6069-6124 Blasticidin resistance gene: bases 6143-6541 SV40 early polyadenylation region: bases 6699-6829 pUC origin: bases 7212-7885 Ampicillin resistance gene (bla): bases 8030-8890 (c) bla promoter: bases 8891-85 (c) (c) = complementary strand

Accessory products

Additional
productsReagents suitable for use with the vector are available
separately. Ordering information for these reagents is
provided below. For more information, go to
www.thermofisher.com or call Technical Support (page 20).

Product	Amount	Catalog No.
pENTR [™] Directional TOPO [™] Cloning Kit	20 reactions	K240020
Gateway [™] LR Clonase [™] Enzyme Mix	20 reactions	11791019
One Shot [™] TOP10 Chemically Competent	10 reactions	C404010
Cells	20 reactions	C404003
One Shot [™] TOP10 Electrocompetent Cells	10 reactions	C404050
	20 reactions	C404052
Ling fastering™ 2000 Descent	1.5 mL	11668019
Liporectamine 2000 Reagent	0.75 mL	11668027
Neon [™] Transfection System	1 each	MPK5000
Blasticidin	50 mg	R21001
β-Gal Antiserum*	50 µL	R90125
β-Gal Assay Kit	100 reactions	K145501
β-Gal Staining Kit	1 kit	K146501
Positope [™] Control Protein	5 µg	R90050
Ni-NTA Agarose	10 mL	R90101
PureLink™ HiPure MiniPrep Kit	25 preps	K210002
PureLink™ HiPure MidiPrep Kit	25 preps	K210004

*Amount supplied is sufficient for 25 Westerns using 10 mL working solution per reaction.

Accessory products, continued

Detection of	Expression of your recombinant fusion protein can be
recombinant	detected using an antibody to the appropriate epitope. The
proteins	amount of antibody supplied is sufficient for 25 westerns.

Product	Epitope	Catalog No.
Anti-V5 Antibody	Detects 14 amino acid epitope derived from the P and V proteins of the paramyxovirus, SV5 (Southern <i>et al.</i> , 1991).	R96025
Anti-V5-HRP Antibody		R96125
Anti-V5-AP Antibody		R96225
	GKPIPNPLLGLDST	
Anti-His (C-term) Antibody	Detects the C-terminal polyhistidine (6x His) tag (requires the free carboxyl group for detection (Lindner <i>et al.,</i> 1997) HHHHHH-COOH	R93025
Anti-His(C-term)-HRP Antibody		R93125
Anti-His(C-term)-AP Antibody		R93225

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 (6x His) tag, you may use Immobilized Metal Affinity Chromatography (IMAC) to purify your recombinant fusion protein. The ProBond[™] Purification System or bulk ProBond[™] resin is available separately. See the table below for ordering information.

Product	Quantity	Catalog No.
ProBond [™] Nickel-chelating Resin	50 mL	R80101
	150 mL	R80115
ProBond [™] Purification System	6 purifications	K85001
ProBond [™] Purification System with Anti-His(C-term)-HRP Antibody	1 kit	K85301
ProBond [™] Purification System with Anti-V5-HRP Antibody	1 kit	K85401
Purification Columns (10 mL polypropylene columns)	50 columns	R64050

Documentation and support

Obtaining support

Technical support	For the latest services and support information for all locations, visit www.thermofisher.com . At the website, you can:				
	• Access worldwide telephone and fax numbers to contact Technical Support and Sales facilities				
	• Search through frequently asked questions (FAQs)				
	• Submit a question directly to Technical Support (thermofisher.com/support)				
	• Search for user documents, SDSs, vector maps and sequences, application notes, formulations, handbooks, certificates of analysis, citations, and other product support documents				
	Obtain information about customer training				
	• Download software updates and patches				
Safety Data Sheets (SDS)	Safety Data Sheets (SDSs) are available at thermofisher.com/support.				
	IMPORTANT! For the SDSs of chemicals not distributed by Thermo Fisher Scientific contact the chemical manufacturer.				
Limited Product Warranty	Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.thermofisher.com/us/en/home/global/terms-and- conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.				

- Ausubel, F. M., Brent, R., Kingston, R. E., Moore, D. D., Seidman, J. G., Smith, J. A., and Struhl, K. (1994). Current Protocols in Molecular Biology (New York: Greene Publishing Associates and Wiley-Interscience).
- Chen, C., and Okayama, H. (1987). High-Efficiency Transformation of Mammalian Cells by Plasmid DNA. Molec. Cell. Biol. 7, 2745-2752.
- Chu, G., Hayakawa, H., and Berg, P. (1987). Electroporation for the Efficient Transfection of Mammalian Cells with DNA. Nucleic Acids Res. 15, 1311-1326.
- Felgner, P. L., Holm, M., and Chan, H. (1989). Cationic Liposome Mediated Transfection. Proc. West. Pharmacol. Soc. *32*, 115-121.
- Felgner, P. L. a., and Ringold, G. M. (1989). Cationic Liposome-Mediated Transfection. Nature 337, 387-388.
- Goodwin, E. C., and Rottman, F. M. (1992). The 3'-Flanking Sequence of the Bovine Growth Hormone Gene Contains Novel Elements Required for Efficient and Accurate Polyadenylation. J. Biol. Chem. 267, 16330-16334.
- Izumi, M., Miyazawa, H., Kamakura, T., Yamaguchi, I., Endo, T., and Hanaoka, F. (1991). Blasticidin S-Resistance Gene (*bsr*): A Novel Selectable Marker for Mammalian Cells. Exp. Cell Res. 197, 229-233.
- Kimura, M., Takatsuki, A., and Yamaguchi, I. (1994). Blasticidin S Deaminase Gene from Aspergillus terreus (BSD): A New Drug Resistance Gene for Transfection of Mammalian Cells. Biochim. Biophys. ACTA 1219, 653-659.
- Kozak, M. (1987). An Analysis of 5'-Noncoding Sequences from 699 Vertebrate Messenger RNAs. Nucleic Acids Res. 15, 8125-8148.
- Kozak, M. (1991). An Analysis of Vertebrate mRNA Sequences: Intimations of Translational Control. J. Cell Biology 115, 887-903.
- Kozak, M. (1990). Downstream Secondary Structure Facilitates Recognition of Initiator Codons by Eukaryotic Ribosomes. Proc. Natl. Acad. Sci. USA 87, 8301-8305.
- Landy, A. (1989). Dynamic, Structural, and Regulatory Aspects of Lambda Sitespecific Recombination. Annu. Rev. Biochem. *58*, 913-949.
- 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 Histagged Proteins With Recombinant Anti-His Tag scFv-Phosphatase or scFv-Phage Fusions. BioTechniques 22, 140-149.
- Miller, J. H. (1972). Experiments in Molecular Genetics (Cold Spring Harbor, New York: Cold Spring Harbor Laboratory).
- Mizushima, S., and Nagata, S. (1990). pEF-BOS, a Powerful Mammalian Expression Vector. Nucleic Acids Res. *18*, 5322.

- Shigekawa, K., and Dower, W. J. (1988). Electroporation of Eukaryotes and Prokaryotes: A General Approach to the Introduction of Macromolecules into Cells. BioTechniques *6*, 742-751.
- 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.
- Takeuchi, S., Hirayama, K., Ueda, K., Sakai, H., and Yonehara, H. (1958). Blasticidin S, A New Antibiotic. The Journal of Antibiotics, Series A 11, 1-5.
- Uetsuki, T., Naito, A., Nagata, S., and Kaziro, Y. (1989). Isolation and Characterization of the Human Chromosomal Gene for Polypeptide Chain Elongation Factor-1α. J. Biol. Chem. *264*, 5791-5798.
- Wigler, M., Silverstein, S., Lee, L.-S., Pellicer, A., Cheng, Y.-C., and Axel, R. (1977). Transfer of Purified Herpes Virus Thymidine Kinase Gene to Cultured Mouse Cells. Cell 11, 223-232.
- Yamaguchi, H., Yamamoto, C., and Tanaka, N. (1965). Inhibition of Protein Synthesis by Blasticidin S. I. Studies with Cell-free Systems from Bacterial and Mammalian Cells. J. Biochem (Tokyo) 57, 667-677.

For support visit **thermofisher.com/support** or email **techsupport@lifetech.com**

thermofisher.com

12 April 2017

