



## Zero Blunt<sup>®</sup> PCR Cloning Kit

# A high efficiency system for cloning blunt-ended PCR products

Catalog numbers K2700-20, K2700-40, K2750-20, and K2750-40

Document Part Number 250162 Publication Number MAN0000040 Revision 5.00



For research use only. Not for use in diagnostic procedures.

### Contents

Contents	
Kit Contents and Storage	4
Introduction	7
Product Overview	7
Methods	8
Produce Blunt-End PCR Products	
Clone into pCR <sup>®</sup> -Blunt	
Transform Competent Cells	
Transform Competent Cells, Continued	
Analyze Transformants	
Blunt-Ending PCR Fragments	
Troubleshoot	
Appendix	
Perform the Control Reactions	
Recipes	
Zeocin <sup>™</sup> Selective Antibiotic	
Map and Features of pCR <sup>®</sup> -Blunt	
Accessory Products	
Technical Support	
Purchaser Notification	
References	

## Kit Contents and Storage

Type of kits	This manual is supplied with the following kits.		
	Kit	Quantity	Cat. no.
	Zero Blunt <sup>®</sup> PCR Cloning Kit	20 reactions	K2750-20
		40 reactions	K2750-40
	Zero Blunt <sup>®</sup> PCR Cloning Kit with One Shot <sup>®</sup>	20 reactions	K2700-20
	TOP10 Chemically Competent E. coli	40 reactions	K2700-40
Shipping and storage	The Zero Blunt <sup>®</sup> PCR Cloning Kits are shipped on a Store the Zero Blunt <sup>®</sup> PCR Cloning reagents at <b>–30</b> ° TOP10 Competent Cells (supplied with Cat. nos. Kito <b>–68°C</b> .	℃ to −10°C and th	
Product use	For research use only. Not for use in diagnostic p	rocedures.	
		Continu	ued on next page

#### Kit Contents and Storage, Continued

## Zero Blunt<sup>®</sup> PCR cloning reagents

Zero Blunt<sup>®</sup> PCR Cloning reagents (Box 1) are listed in the following table. Forty reaction kits are supplied as two 20-reaction kits.**Store reagents at –30°C to –10°C.** 

Item	Concentration	Amount
pCR <sup>®</sup> -Blunt vector, linearized with blunt ends	25 ng/μL in TE Buffer, pH 8	3 × 10 μL, 750 ng
ExpressLink™ T4 DNA Ligase⁺	5.0 Weiss units/µL	25 μL
5X Express Link™ T4 DNA Ligase Buffer**	$\begin{array}{l} 5X\ T4\ DNA\ Ligase\ Buffer\ (50\ mM\\ Tris-HCl,\ pH\ 7.6\ ,\ 50\ mM\ MgCl_2,\ 5\\ mM\ ATP,\ 5\ mM\ dithiothreitol,\ 25\\ \%\ (w/v)\ polyethylene\ glycol-8000) \end{array}$	200 µL
dNTP Mix	12.5 mM dATP 12.5 mM dCTP 12.5 mM dGTP 12.5 mM dTTP Neutralized to pH 8.0	10 μL
Control DNA Template	$0.1 \mu\text{g}/\mu\text{L}$ in TE Buffer, pH 8	10 µL
Sterile Water	Deionized, autoclaved	1 mL
M13 Forward (–20) Primer	0.1 μg/μL in TE Buffer, pH 8	20 µL
M13 Reverse Primer	$0.1 \mu\text{g}/\mu\text{L}$ in TE Buffer, pH 8	20 µL

## Sequence of primers

The following table lists the sequence of the M13 Sequencing Primers supplied with the kit

Primer	Sequence	pMoles Supplied
M13 Forward (-20)	5'-GTAAAACGACGGCCAG-3'	407
M13 Reverse	5´-CAGGAAACAGCTATGAC-3´	385

#### Kit Contents and Storage, Continued

One Shot<sup>®</sup> reagents

One Shot<sup>®</sup> TOP10 Chemically Competent *E. coli* are supplied with Cat. nos. K2700-20 and K2700-40 and contain the following reagents. Forty reaction kits are supplied as two 20-reaction kits.

The transformation efficiency for TOP10 cells is  $1 \times 10^9$  cfu/µg DNA.

Store competent cells at -85°C to -68°C.

Item	Composition	Amount
S.O.C. Medium	2% Tryptone	6 mL
(may be stored at 2°C to 8°C	0.5% Yeast Extract	
or 15°C to 30°C)	10 mM NaCl	
	2.5 mM KCl	
	10 mM MgCl2	
	10 mM MgSO4	
	20 mM glucose	
TOP10 cells		$21\times 50~\mu L$
pUC19 Control DNA	$10 \text{ pg/}\mu\text{L}$ in 5 mM Tris-HCl,	50 µL
-	0.5 mM EDTA, pH 8	

Genotype of TOP10

TOP10: F<sup>-</sup> mcrA  $\Delta$ (mrr-hsdRMS-mcrBC)  $\Phi$ 80lacZ $\Delta$ M15  $\Delta$ lacX74 recA1 araD139  $\Delta$ (ara-leu)7697 galU galK rpsL (Str<sup>R</sup>) endA1 nupG

#### Introduction

#### **Product Overview**

Description of the			
Description of the system	any blu pCR®-B LacZα of the <i>li</i> upon tr	To Blunt <sup>®</sup> PCR Cloning Kit is designed to clone blunt PCR fragme ont DNA fragment) with a low background of non-recombinants. For vector contains the lethal <i>E. coli ccdB</i> gene fused to the C-terr (Bernard <i>et al.,</i> 1994). Ligation of a blunt PCR fragment disrupts e $acZ\alpha$ - <i>ccdB</i> gene fusion permitting growth of only positive recomb ansformation. Cells that contain non-recombinant vector are killen asformation mixture is plated.	The ninus of xpression inants
<i>ccdB</i> gene	locus co amino a 1983). T segrega	B gene is found in the <i>ccd</i> ( <u>c</u> ontrol of <u>c</u> ell <u>d</u> eath) locus on the F pla ontains two genes, <i>ccd</i> A and <i>ccd</i> B, which encode proteins of 72 an acids respectively (Karoui <i>et al.</i> , 1983; Miki <i>et al.</i> , 1984; Ogura and <i>Che ccd</i> locus participates in stable maintenance of F plasmid by p tional killing of cells that do not contain the F plasmid (Jaffe <i>et al.</i> )	d 101 Hiraga, ost- , 1985).
	protein (topoise	dA protein acts as an inhibitor of CcdB protein. In the absence of , the CcdB protein becomes toxic as it interferes with bacterial DN omerase II), an essential enzyme that catalyzes the ATP-depender piling of DNA. Inhibiting DNA gyrase results in DNA breakage a	IA gyrase nt negative
Experimental outline			sformed either triction
	Chara		
	Step	Action	Page
	1	Action         Amplify your PCR product using a thermostable, proofreading polymerase and your own primers and parameters.	Page 8
		Amplify your PCR product using a thermostable, proofreading polymerase and your own primers and	
	1	Amplify your PCR product using a thermostable, proofreading polymerase and your own primers and parameters.	8

#### Methods

### **Produce Blunt-End PCR Products**

Guidelines for PCR	amplifying a pool of cDN abundance of the messag	NA is sufficient to use as a template for PCR. If A, the amount needed will depend on the relative e of interest in your mRNA population. For optimal ecommend using no more than 30 cycles of	
<b>Q</b> Important	blunt fragments. If you u	se to produce your PCR product as it will not produce se <i>Taq</i> polymerase or an enzyme mix containing <i>Taq</i> 15 for a protocol to blunt-end your PCR product before	
Required	Components required but n	ot supplied:	
materials	Thermostable proofr	ading polymerase	
	• 10X PCR buffer appr	opriate for your polymerase	
	Thermocycler		
	DNA template and primers for PCR product		
	Components supplied with the kit:		
	<ul> <li>dNTPs (adjusted to p</li> </ul>	H 8)	
Produce PCR	Perform the PCR in a 50	L volume containing:	
products	DNA Template	10–100 ng	
	10X PCR Buffer	5 µL	
	100 mM dNTPs	0.5 μL	
	Primers	1 μM each	
	Sterile water to	a total volume of 49 μL	
	Proofreading Polymerase	<u>(2 to 2.5 U/μL) 1 μL</u>	
	Total Volume	50 µL	
Gel purification	verify the quality and qu single, discrete band from proceeding. Take special long exposure to UV ligh eliminate multiple bands	The PCR reaction and use agarose gel electrophoresis to antity of your PCR product. If you do not obtain a a your PCR, you may gel-purify your fragment before care to avoid sources of nuclease contamination and t. Alternatively, you may optimize your PCR to and smearing (Innis <i>et al.</i> , 1990). The PCR Optimizer <sup>TM</sup> you optimize your PCR. Contact Technical Support for 7).	

## Clone into pCR<sup>®</sup>-Blunt

Introduction	After producing the desired PCR product, you are ready to clone it into pCR <sup>®</sup> -Blunt. Follow the following guidelines.
	If this is the first time you are using the Zero Blunt <sup>®</sup> PCR Cloning Kit, we recommend performing the control reactions to help you evaluate your results (see pages 18–20).
	We recommend using a 10:1 molar ratio of insert:vector as a starting point. You may need to increase this ratio to optimize cloning efficiency. If you are using the PCR product directly from the PCR reaction mixture, do not use more than 5 $\mu$ L of the reaction since salts in the PCR sample may inhibit the T4 DNA Ligase.
Calculating amount of PCR product to use	A range from 10:1 to 100:1 molar ratio of insert:vector is recommended to optimize blunt-ended PCR ligation efficiency. If necessary, the amount of vector in the ligation reaction can be reduced to as little as 5 ng. Use the following formula below to estimate the amount of PCR product needed for a 10:1 molar ratio of insert to 25 ng of pCR <sup>TM</sup> -Blunt vector.
	x ng insert = $\frac{(10) (y \text{ bp PCR product}) (25 \text{ ng linearized pCR}^{\otimes}-\text{Blunt})}{(25 \text{ ng linearized pCR}^{\otimes}-\text{Blunt})}$
	(3500 bp pCR <sup>®</sup> -Blunt)
	where $x$ ng is the amount of PCR product of $y$ base pairs to be ligated for a 10:1 insert:vector molar ratio.
Purify and concentrate PCR products	In most cases, the PCR reaction may be used directly in the ligation reaction; however, it may be necessary on occasion to purify and concentrate the PCR product. Follow the protocol below to purify your PCR product.
	1. Extract the PCR reaction with an equal volume of phenol/chloroform. Centrifuge 1 minute at $10,000 \times g$ and transfer the upper layer to a fresh, sterile tube.
	<ol> <li>Add 1/10 volume of 3 M sodium acetate, pH 5.6 and add 2 volumes of 100% ethanol at room temperature. You may also add 5 μg of carrier (e.g. glycogen) to help precipitate the DNA.</li> </ol>
	3. Flick the tube to mix and centrifuge at 10,000 × $g$ for 5 minutes.
	4. Remove the ethanol and rinse the pellet with 70% ethanol.
	5. Centrifuge and remove all the ethanol. Air-dry the pellet for ~5 minutes.
	6. Resuspend the pellet in $5-10\mu$ L of sterile water or TE buffer.
	7. Estimate the DNA concentration by electrophoresis or absorbance at 260 nm.
	Continued on next page

## Clone into pCR<sup>®</sup>-Blunt, Continued

Ligate into pCR <sup>®</sup> -Blunt		llow the following protocol to ligate your PC Determine the volume of PCR sample need PCR product (see page 9).	
	2.	Set up the following $10\mu L$ ligation reaction:	
		pCR <sup>®</sup> -Blunt (25 ng)	1 μL
		Blunt PCR product	1–5 μL
		5X ExpressLink <sup>™</sup> T4 DNA Ligase Buffer	2 µL
		Sterile water to a total volum	ne of 9 µL
	_	ExpressLink <sup>TM</sup> T4 DNA Ligase (5 U/ $\mu$ L)	1 μL
		Total Volume	10 µL
	1	Incubate the ligation reaction at room temp 5 minutes. Longer ligation times increase th Control Ligation Reaction section below. Pr <b>Cells</b> , page 11. Proceed to <b>Transform Comp</b> page.	e cloning efficiency, see the oceed to <b>Transform Competent</b>

Note: You may store your ligation reaction at  $-20^{\circ}$ C until you are ready for transformation

## **Transform Competent Cells**

Introduction	After ligating your insert into pCR <sup>®</sup> -Blunt, you are ready to transform the construct into competent <i>E. coli</i> . One Shot <sup>®</sup> TOP10 cells are provided with Cat. nos. K2700-20 and K2700-40 to facilitate transformation. A protocol to transform One Shot <sup>®</sup> cells is provided in this section. To transform another competent strain, refer to the manufacturer's instructions.
<i>E. coli</i> host strain	We recommend using One Shot <sup>®</sup> TOP10 <i>E. coli</i> for transformation, however other host strains are suitable (see the following <b>Important</b> note). TOP10 does <b>not</b> contain <i>lac</i> I <sup>q</sup> ; therefore, expression from the <i>lac</i> promoter is constitutive. <b>There is no need to use IPTG or X-gal to identify possible recombinants.</b> If you use a strain that contains <i>lac</i> I <sup>q</sup> , you will need to include IPTG to induce expression from the <i>lac</i> promoter since expression is required to produce the lethal phenotype.
<b>Q</b> Important	Do not use any <i>E. coli</i> strain that contains the complete Tn5 transposon (e.g. DH5 F'IQ). This transposon contains resistance markers to both Kanamycin and Zeocin <sup>TM</sup> selective antibiotic and will not allow selection of positive recombinants. Do not use any <i>E. coli</i> strain that contains the F' episome (e.g. TOP10F'). These strains contain the <i>ccd</i> A gene and will prevent negative selection with the <i>ccd</i> B gene.
Zeocin <sup>™</sup> selection	The Zeocin <sup>TM</sup> resistance gene in pCR <sup>®</sup> -Blunt allows you to select <i>E. coli</i> transformants using Zeocin <sup>TM</sup> antibiotic. For selection, use Low Salt LB agar plates containing 25 µg/mL Zeocin <sup>TM</sup> selective antibiotic (see page 21 for a recipe). Note that for Zeocin <sup>TM</sup> selective antibiotic to be active, the salt concentration of the bacterial medium must remain low (less than 90 mM) and the pH must be 7.5. We recommend growing positive recombinants in SOB medium containing 25 µg/mL Zeocin <sup>TM</sup> selective antibiotic for better growth and plasmid yields (see page 23 for a recipe). For more information on storing and handling Zeocin <sup>TM</sup> selective antibiotic, refer to page 23.

## Transform Competent Cells, Continued

Required	Components required but not supplied:
materials	• LB plates containing 50 µg/mL kanamycin or Low Salt LB plates containing 25 µg/mL Zeocin <sup>™</sup> selective antibiotic (see page 21 for a recipe) (two for each transformation)
	• 42°C water bath
	• 37°C shaking and non-shaking incubator
	General microbiological supplies (e.g. plates, spreaders)
	• SOB Medium, optional (warmed to room temperature; see page 22 for recipe) <i>Components supplied with the kit:</i>
	<ul> <li>S.O.C. medium (warmed to room temperature)</li> </ul>
	• Chemically competent <i>E. coli</i> suitable for transformation (One Shot <sup>®</sup> TOP10 Competent Cells (supplied with Cat. nos. K2700-20 and K2700-40)
	• Positive control, optional ( <i>e.g.</i> pUC19)
Prepare for	• Equilibrate a water bath to 42°C.
transformation	• Warm the vial of S.O.C. medium or SOB medium to room temperature.
	• Warm selective plates at 37°C for 30 minutes
One Shot <sup>®</sup> transformation	Follow the protocol below to transform One Shot <sup>®</sup> TOP10 Competent Cells. To transform another strain, refer to the manufacturer's instructions.
protocol	1. Thaw on ice one 50-µL vial of One Shot <sup>®</sup> TOP10 cells for each transformation.
	<ol> <li>Pipet 2 μL of each ligation reaction directly into the vial of competent cells and mix by stirring gently with a pipette tip. Do not mix cells by pipetting.</li> </ol>
	3. Incubate the vials on ice for 30 minutes. Store the remaining ligation mixtures at $-20^{\circ}$ C.
	4. Heat shock the cells for 45 seconds at 42°C without shaking. Immediately place the vials on ice for 2 minutes.
	5. Add 250 µL of room temperature S.O.C. medium to each vial.
	<ol> <li>Shake the vials horizontally at 37°C for 1 hour at 225 rpm in a shaking incubator.</li> </ol>
	7. Spread 10–100 µL from each transformation vial on LB plates containing 50 µg/mL kanamycin or Low Salt LB plates containing 25 µg/mL Zeocin <sup>™</sup> selective antibiotic. We recommend plating two different volumes to ensure well-spaced colonies. For plating smaller volumes, add 20 µL S.O.C. to ensure even spreading.
	8. Incubate plates overnight at 37°C. Proceed to <b>Analyze Transformants</b> on page 13
Expected results	For an insert size of 400–800 bp, you should obtain 50–2000 colonies per plate depending on the volume plated and of these, greater than 80% should contain inserts. Note that ligation and transformation efficiency depend on insert size. As insert size increases, the efficiency will decrease.

## Analyze Transformants

Analyze positive clones	<ol> <li>Pick at least 10 transformants for analysis.</li> <li>Grow colonies overnight in 2 mL LB or SOB medium containing 50 µg/mL kanamycin or Low Salt LB containing 25 µg/mL Zeocin<sup>™</sup> selective antibiotic.</li> <li>Isolate plasmid and analyze by restriction digestion or sequencing to verify the correct insert and orientation. We recommend using the Pure Link<sup>™</sup> HQ Mini Plasmid Purification Kit for purifying your plasmid DNA (see page 26 for ordering information).</li> </ol>		
Sequence	You may sequence your pCR <sup>®</sup> -Blunt construct to confirm the presence and correct orientation of the insert. The M13 Forward and Reverse sequencing primers are included in the kit to help you sequence the insert. Refer to the diagram on page 24 for the location of the primer binding sites.		
Analyze transformants by PCR	You may wish to use PCR to directly analyze positive transformants. For PCR primers, use either the M13 Forward (–20) or the M13 Reverse primer and a primer that hybridizes within your insert. If you are using this technique for the first time, we recommend performing restriction analysis in parallel. Artifacts may be obtained because of mispriming or contaminating template. The protocol is provided below for your convenience. Other protocols are suitable.		
	Materials Needed		
	PCR SuperMix High Fidelity (see page 26)		
	Appropriate forward and reverse PCR primers (20 µM each)		
	Procedure		
	<ol> <li>For each sample, aliquot 48 μL of PCR SuperMix High Fidelity into a 0.5-mL microcentrifuge tube. Add 1 μL each of the forward and reverse PCR primer.</li> </ol>		
	<ol> <li>Pick 10 colonies and resuspend them individually in 50 µL of the PCR cocktail from step 1 of this procedure. Don't forget to make a patch plate to preserve the colonies for further analysis.</li> </ol>		
	3. Incubate the reaction for 10 minutes at 94°C to lyse the cells and inactivate nucleases.		
	4. Amplify for 20–30 cycles.		
	5. For the final extension, incubate at 72°C for 10 minutes. Store at 4°C.		
	6. Visualize by agarose gel electrophoresis.		
<b>Q</b> Important	If you have problems obtaining transformants or the correct insert, perform the control reactions described on pages 18–20. These reactions will help you troubleshoot your experiment. Refer to <b>Troubleshoot</b> on page 16 for additional tips.		

## Analyze Transformants, Continued

Long-Term storage	After identifying the correct clone, be sure purify the colony and make a glycerol stock for long-term storage. We recommend that you store a stock of plasmid DNA at $-20^{\circ}$ C.		
	1.	Streak the original colony on LB plates containing 50µg/mL kanamycin.	
	2.	Isolate a single colony and inoculate into 1–2 mL of LB containing 100 μg/mL ampicillin or 50μg/mL kanamycin.	
	3.	Grow until culture reaches stationary phase.	
	4.	Mix 0.85 mL of culture with 0.15 mL of sterile glycerol and transfer to a cryovial.	
	5.	Store at -80°C.	

## **Blunt-Ending PCR Fragments**

Introduction	If you used <i>Taq</i> polymerase or an enzyme mix containing <i>Taq</i> polymerase to generate your PCR product, you will need to blunt-end the PCR product (i.e. remove the 3' A-overhangs) before cloning into pCR <sup>®</sup> -Blunt To clone PCR products generated using <i>Taq</i> polymerase, we recommend the TOPO TA Cloning <sup>®</sup> Kit (see page 26). The cloning vector supplied with this kit, pCR <sup>®</sup> -Blunt, is provided linearized with 3' T-overhangs which will pair with the 3' A-overhangs on <i>Taq</i> -generated PCR products.		
Before starting	<ol> <li>Prepare a 10-mM dNTP solution by diluting 1 μL of a 100-mM dNTP solution into 9 μL sterile water. Place on ice.</li> <li>Equilibrate at water bath or heat block to 72°C.</li> </ol>		
Procedure	<ol> <li>Purify and concentrate the PCR product as described on page 9. <i>Taq</i> polymerase must be removed to prevent 3' extension from competing with blunt-ending.</li> <li>Set up the following ~10 μL reaction:         <ul> <li>PCR product (or any DNA fragment)</li> <li>8 μL</li> <li>10 mM dNTPs</li> <li>1 μL</li> <li>10X PCR Buffer (user supplied)</li> <li>1 μL</li> <li>Thermostable, proofreading polymerase (0.5 unit)</li> <li>0.25 μL</li> </ul> </li> <li>Mix by flicking the tube and add 2 drops of mineral oil. Centrifuge briefly.</li> <li>Incubate at 72°C for 20 to 30 minutes.</li> </ol>		
	5. Use 1–5 $\mu$ L directly in the ligation reaction, page 10.		
Using other enzymes	If you do not have a thermostable, proofreading polymerase, you may try T4 DNA Polymerase to blunt end your fragment. Follow the manufacturer's instructions to blunt-end your PCR product, then proceed to the ligation reaction on page 10.		

#### Troubleshoot

Introduction	If you do not obtain the results you expect, use the following table to troubleshoot your experiment. We recommend performing the control reactions (pages 18–20) to help you evaluate your results.			
Observation	Cause	Solution		
No PCR product	Either the thermostable polymerase is inactive or the conditions for your PCR are not optimal.	Perform the control reactions on pages 18–20 to test the activity of the polymerase. If the polymerase is active, you may need to optimize the conditions of your PCR reaction.		
Few or no transformants obtained	Loss of DNA during precipitation	Be careful not to lose the DNA pellet when precipitating and washing. You may wish to add carrier to help precipitate the DNA (e.g. glycogen).		
	Molar ratio of insert to vector is incorrect	Determine the concentration of insert and use a molar ratio of at least 10:1 (insert:vector). Increase the molar ratio up to 100:1 if necessary.		
	Low transformation efficiency of <i>E. coli</i> strain	Check transformation efficiency with the pUC19 control vector included with the One Shot <sup>®</sup> Kit. One Shot <sup>®</sup> TOP10 cells should yield $\sim 1 \times 10^9$ transformants/µg DNA.		
	A polymerase that adds 3' A-overhangs was used (e.g. <i>Taq</i> polymerase)	If you used an enzyme mix containing <i>Taq</i> polymerase, you will need to blunt-end your PCR product (see page 15).		
	Too much of the amplification reaction was added to the ligation.	The high salt content of PCR reactions can inhibit ligation. Do not use more than 5 $\mu$ L of the PCR mixture in the ligation reaction.		
High background of transformants that do not contain inserts	Nuclease contamination in reagents	Autoclave all reagents used for cloning (especially water).		
	Absence of insert allows vector to religate	Increase the molar ratio of insert to vector (see page 9).		
	Insert too large for efficient transformation	Try electroporation with electrocompetent cells. Do not use One Shot <sup>®</sup> Chemically Competent Cells for electroporation.		
	Difficult insert to clone by the usual methods	Some DNA sequences may be difficult to replicate and propagate (i.e. inverted repeats, AT- or GC- rich).		

## Troubleshoot, Continued

Observation	Cause	Solution	
Thin "lawn" of cells on plate	Insufficient amount of antibiotic in plate medium	Use LB plates containing 50 µg/mL kanamycin or Low Salt LB plates containing 25 µg/mL Zeocin <sup>™</sup> selective antibiotic. Let medium cool sufficiently (~55°C) before adding antibiotics.	
	High salt in LB medium when selecting on Zeocin <sup>™</sup> selective antibiotic	Make sure you make LB medium with 5 g/liter NaCl (see page 21 for a recipe).	
	Too many cells plated	Plate a smaller volume of the transformation mixture or dilute 10-fold with S.O.C. before plating.	
DNA bands run at a larger molecular weight than expected and seem slightly smeared	Protein bound to DNA	Perform phenol/chloroform extraction during plasmid preparation of after restriction digest. You may also use the PureLink <sup>™</sup> HQ Mini Plasmid Purification Kit (see page 26).	
Low plasmid yield	Medium is not sufficiently rich enough for optimal growth.	Use SOB medium containing antibiotic (see page 22 for recipe).	

#### Appendix

#### **Perform the Control Reactions**

Introduction	hel pro	e recommend performin p you evaluate your res oducing a control PCR p e control PCR product in	ults. Performi roduct using	ng the control re the reagents incl	eactions inv	olve	
Produce the	1.	Set up a 50µL PCR as follows:					
control PCR		Control DNA Template (100 ng)				1 μL	
product		10X PCR Buffer				5 μL	
		100 mM dNTPs			0.	5 µL	
		M13 Reverse Primer (2		2 μL			
		M13 Forward (–20) Primer (200 ng)				2 µL	
		Sterile Water			38.5 μL		
		Proofreading polymerase (2.5 units)				<u>1 µL</u>	
		Total Volume			5	0 μL	
	2.	Amplify using the follo					
		Step	Time	Temperature	Cycles		
		Initial Denaturation	2 minutes	94°C	1	-	
		Denature	1 minute	94°C		-	
		Anneal	1 minute	55°C	25		
		Extend	1 minute	72°C			
		Final Extension	7 minutes	72°C	1	-	
	3.	Remove 10 µL from the A discrete 800-bp band			ose gel elec	ctrophoresis.	

4. Estimate the amount of DNA by measuring against a known standard run on the same gel. You should get a concentration of about 20–30 ng/µL for your Control PCR Product. Proceed to the **Control Ligation Reaction** on page 19.

## Perform the Control Reactions, Continued

Control ligation reaction	Using the control PCR product produced on the previous page, set up the following ligation reaction.		
	. Set up the 10 μL Control Ligation Reaction as follows. Note that the insert:vector molar ratio is 10:1.		
	5X ExpressLink <sup>TM</sup> T4 DNA Ligase Buffer $2 \mu L$		
	pCR <sup>®</sup> -Blunt vector (25 ng/ $\mu$ L) 1 $\mu$ L		
	Control PCR Product (~60 ng) 3 µL		
	Sterile water 3 µL		
	ExpressLink <sup>TM</sup> T4 DNA Ligase (5 U/ $\mu$ L) 1 $\mu$ L		
	Total Volume 10 µL		
	<ol> <li>Incubate the Control Ligation Reaction at room temperature for 1 hour. (See Note about incubation times on next page)</li> </ol>		
	Transform 1 µL of the Control Ligation Reaction into One Shot <sup>®</sup> TOP10 Competent Cells or into another suitable competent <i>E. coli</i> strain.		
	Plate 10–50 μL of each transformation mix on LB agar plates containing 50 μg/mL kanamycin or Low Salt LB plates containing 25 μg/mL Zeocin <sup>™</sup> selective antibiotic.		
	. Incubate plates overnight at 37°C.		
Transformation control	Zero Blunt <sup>®</sup> PCR Cloning Kits supplied with One Shot <sup>®</sup> TOP10 Competent Cells will also be supplied with pUC19 plasmid for use as a transformation control. Transform one vial of One Shot <sup>®</sup> TOP10 cells with 10 pg of pUC 19 using the protocol on page 12. Plate 10–50 $\mu$ L of the transformation mixture on LB plates containing 100 $\mu$ g/mL ampicillin. Transformation efficiency should be greater than 1 × 10 <sup>9</sup> cfu/ $\mu$ g DNA.		
	Continued on the next pa	ıge	

#### Perform the Control Reactions, Continued

Note about incubation times

The total number of transformants was reduced with shorter ligation incubation times. In the tables below, ligation reactions were performed using either Accuprime<sup>®</sup> PFx-generated blunt insert **(Table 1)** or Platinum<sup>®</sup> PFx-generated blunt insert **(Table 2)**.

Following the control ligation procedure using a 1:10 vector to insert ratio. Ligation reactions were incubated at room temperature for 5 minutes, 10 minutes, 15 minutes, 30 minutes and 1 hour, transformed into One Shot<sup>®</sup> Top10 cells and 50ul plated for colony formation. The data at each time point shows the total colony number for 3 replicates.

Table 1 - Accuprime <sup>®</sup> PFx		Table 2 - Platinum <sup>®</sup> PFx	
Time	Colonies	Time	Colonies
5 min	$40 \pm 12$	5 min	$30 \pm 19$
10 min	77 ± 17	10 min	$73 \pm 14$
15 min	$85 \pm 14$	15 min	$120 \pm 60$
30 min	$148 \pm 50$	30 min	$192 \pm 49$
60 min	$166 \pm 63$	60 min	$339 \pm 44$

# **Expected results** Pick 10 colonies and isolate plasmid DNA. Analyze the plasmids for the presence of insert by digesting the DNA with *Eco*R I to release the insert. Greater than 95% of the colonies should contain plasmid with the 800 bp insert.

## Recipes

LB (Luria-Bertani) medium and plates	Composition: 1.0% Tryptone 0.5% Yeast Extract 1.0% NaCl pH 7.0
	<ol> <li>For 1 liter, dissolve 10 g tryptone, 5 g yeast extract, and 10 g NaCl in 950 mL deionized water.</li> </ol>
	2. Adjust the pH of the solution to 7.0 with NaOH and bring the volume up to 1 liter.
	3. Autoclave on liquid cycle for 20 minutes at 15 psi. Allow the solution to cool to 55°C and add antibiotic if needed.
	4. Store at room temperature or at 4°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 (50 $\mu$ g/mL kanamycin or 50 $\mu$ g/mL ampicillin), and pour into 10-cm plates.
	4. Let the plates harden, then invert and store at 4°C in the dark.
Low salt LB medium	<ul> <li>Composition: <ol> <li>0.5% Tryptone</li> <li>0.5% Yeast Extract</li> </ol> </li> <li>0.5% NaCl pH 7.5 <ol> <li>For 1 liter, dissolve 10 g tryptone, 5 g yeast extract, and 5 g NaCl in 950 mL deionized water. For plates, add 15 g/L agar.</li> <li>Adjust the pH of the solution to 7.5 with NaOH and bring the volume up to 1 liter.</li> <li>Autoclave on liquid cycle for 20 minutes at 15 psi. Allow solution to cool to 55°C and add Zeocin<sup>™</sup> selective antibiotic to a final concentration of 25 µg/mL.</li> </ol> </li> <li>Store at room temperature or at 4°C.</li> </ul>
	Continued on next page

## Recipes, Continued

SOB medium with antibiotic	Composition: SOB (per liter)2%Tryptone0.5%Yeast Extract0.05%NaCl2.5 mMKCl10 mMMgCl2
	1. Dissolve 20 g tryptone, 5 g yeast extract, and 0.5 g NaCl in 950 mL deionized water.
	2. Prepare 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 of this procedure.
	3. Adjust pH to 7.5 with 5 M NaOH and add deionized water to 1 liter.
	4. Autoclave this solution, cool to ~55°C, and add 10 mL of sterile 1 M MgCl <sub>2</sub> .
	5. If desired, add kanamycin to a final concentration of 50 $\mu$ g/mL or Zeocin <sup>TM</sup> selective antibiotic to a concentration of 25 $\mu$ g/mL.
	6. Store at room temperature or 4°C.

#### Zeocin<sup>™</sup> Selective Antibiotic

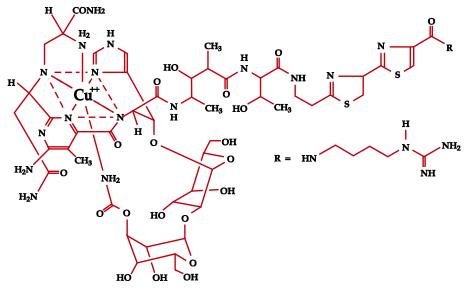
#### Introduction

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

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

#### Molecular weight, formula, and structure

The formula for Zeocin<sup> $^{\text{M}}$ </sup> selective antibiotic is  $C_{60}H_{89}N_{21}O_{21}S_3$  and the molecular weight is 1527.5. The structure of Zeocin<sup> $^{\text{M}}$ </sup> selective antibiotic is shown below.

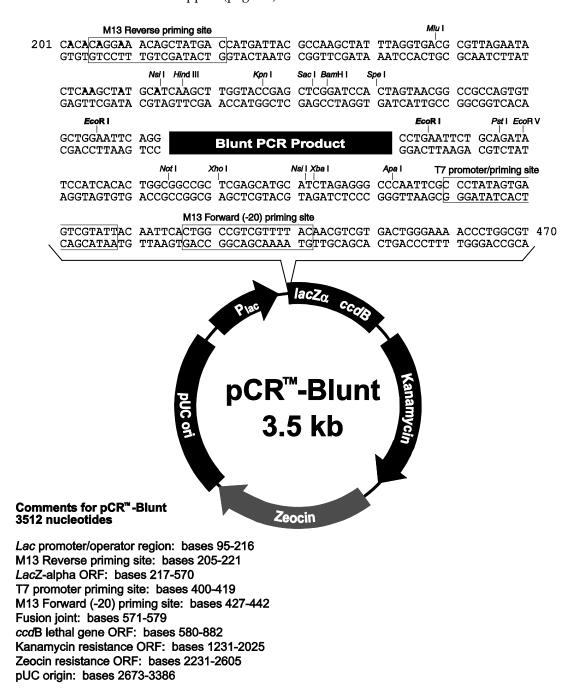


## Handling Zeocin<sup>™</sup> selective antibiotic

- High ionic strength and acidity or basicity inhibit the activity of Zeocin<sup>™</sup> selective antibiotic. Therefore, we recommend that you reduce the salt in bacterial medium and adjust the pH to 7.5 to keep the drug active (see page 21 for a recipe).
- Store Zeocin<sup>™</sup> selective antibiotic at −20°C and thaw on ice before use.
- Zeocin<sup>™</sup> selective antibiotic is light sensitive. Store the drug and plates or medium containing the drug in the dark.
- Wear gloves, a laboratory coat, and safety glasses when handling solutions that contain Zeocin<sup>™</sup>.
- Do not ingest or inhale solutions containing Zeocin<sup>™</sup>.

#### Map and Features of pCR<sup>®</sup>-Blunt

**pCR<sup>™</sup>-Blunt map** The following map summarizes the features of the pCR<sup>®</sup>-Blunt vector. Restriction sites that are only found in the polylinker are shown. **The complete sequence of the vector** is available from **www.lifetechnologies.com/support** or by contacting Technical Support (page 27).



## Map and Features of pCR®-Blunt, Continued

## Features of pCR<sup>®</sup>-Blunt

The table below describes the features of  $pCR^{\circledast}$ -Blunt. All features have been functionally tested.

Features	Benefit
lac promoter	Allows expression of $lacZ\alpha$ -ccdB gene fusion.
<i>lacZα-ccd</i> B fusion gene	Provides positive selection against non-recombinant vector. ccdB function is disrupted (inactivated) by an insert. Includes the M13 forward and reverse priming sites for sequencing of the insert.
Multiple Cloning Site	Allows insertion of PCR fragment to disrupt expression of the <i>ccd</i> B gene and allows convenient screening, restriction mapping, and excision of cloned insert.
T7 promoter	Allows <i>in vivo</i> or <i>in vitro</i> transcription and translation.
T7, M13 Forward (–20), M13 Reverse priming sites	Allows sequencing of the insert.
Kanamycin resistance gene	Allow selection and maintenance in <i>E. coli</i> .
<i>Sh ble</i> (Zeocin <sup>™</sup> resistance gene)	Allow selection and maintenance in <i>E coli</i> . Expression is driven using the <i>bla</i> ( $\beta$ -lactamase) promoter.
pUC origin	Allows high copy replication and maintenance of the plasmid in <i>E. coli</i> .

#### **Accessory Products**

## Additional products

The following table lists additional products that may be used with TOPO<sup>®</sup> TA Cloning Kits. For more information, visit **www.lifetechnologies.com/support** or contact Technical Support (page 27).

Item	Quantity	Cat. no.
TA Cloning Dual Promoter Kit	40 reactions	K460040
ExpressLink <sup>™</sup> T4 DNA Ligase	30 reactions	A13726
Platinum <sup>®</sup> Pfx DNA Polymerase	100 units	11708-013
	250 units	11708-021
	500 units	11708-039
AccuPrime <sup>™</sup> <i>Pfx</i> DNA Polymerase	200 reactions	12344-024
	1000 reactions	12344-032
AccuPrime <sup>™</sup> <i>Pfx</i> SuperMix	200 reactions	12344-040
PCR SuperMix High Fidelity	100 reactions	10790-020
The PCR Optimizer <sup>™</sup> Kit	100 reactions	K1220-01
PureLink <sup>®</sup> HQ Mini Plasmid Purification Kit	100 reactions	K2100-01
One Shot <sup>®</sup> TOP10 Chemically Competent	10 reactions	C4040-10
E. coli	20 reactions	C4040-03
	40 reactions	C4040-06
One Shot <sup>®</sup> TOP10 Electrocompetent <i>E. coli</i>	10 reactions	C4040-50
1	20 reactions	C4040-52
One Shot <sup>®</sup> Mach1 <sup>™</sup> -T1 <sup>R</sup> Chemically Competent <i>E. coli</i>	20 reactions	C8620-03
TOPO <sup>®</sup> TA Cloning <sup>®</sup> Kit	20 reactions	K4500-01
Kanamycin	5 g	11815-024
-	25 g	11815-032
	100 ml (10 mg/ml)	15160-054
Zeocin <sup>™</sup> Selection Reagent	8 x 1.25 ml	R250-01
~	50 ml	R250-05

## **Technical Support**

Obtaining support	<ul> <li>For the latest services and support information for all locations, go to www.lifetechnologies.com/support.</li> <li>At the website, you can: <ul> <li>Access worldwide telephone and fax numbers to contact Technical Support and Sales facilities</li> <li>Search through frequently asked questions (FAQs)</li> <li>Submit a question directly to Technical Support (techsupport@lifetech.com)</li> <li>Search for user documents, SDSs, vector maps and sequences, application notes, formulations, handbooks, certificates of analysis, citations, and other product support documents</li> <li>Obtain information about customer training</li> <li>Download software updates and patches</li> </ul> </li> </ul>
Safety Data Sheets (SDS)	Safety Data Sheets (SDSs) are available at <b>www.lifetechnologies.com/support</b> .
Certificate of analysis	The Certificate of Analysis provides detailed quality control and product qualification information for each product. Certificates of Analysis are available on our website. Go to <b>www.lifetechnologies.com/support</b> and search for the Certificate of Analysis by product lot number, which is printed on the box.
Limited warranty	Life Technologies Corporation is committed to providing our customers with high- quality goods and services. Our goal is to ensure that every customer is 100% satisfied with our products and our service. If you should have any questions or concerns about a Life Technologies product or service, contact our Technical Support Representatives. All Life Technologies products are warranted to perform according to specifications stated on the certificate of analysis. The Company will replace, free of charge, any product that does not meet those specifications. <u>This warranty limits</u> <u>the Company's liability to only the price of the product</u> . No warranty is granted for products beyond their listed expiration date. No warranty is applicable unless all product components are stored in accordance with instructions. The Company reserves the right to select the method(s) used to analyze a product unless the Company agrees to a specified method in writing prior to acceptance of the order. Life Technologies makes every effort to ensure the accuracy of its publications, but realizes that the occasional typographical or other error is inevitable. Therefore the Company makes no warranty of any kind regarding the contents of any publications or documentation. If you discover an error in any of our publications, report it to our Technical Support Representatives. Life Technologies Corporation shall have no responsibility or liability for any special, incidental, indirect or consequential loss or damage whatsoever. The above limited warranty is sole and exclusive. No other warranty is made, whether expressed or implied, including any warranty of merchantability or fitness for a particular purpose.

#### **Purchaser Notification**

Limited Use Label License No. 54: <i>ccd</i> B-Fusion Vectors	ccdB selection technology is described in Bernard et al., "Positive Selection Vectors Using the F Plasmid ccdB Killer Gene" Gene 148 (1994) 71-74. 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). For licensing information for use in other than research, please contact: Out Licensing, Life Technologies Corporation, 5791 Van Allen Way, Carlsbad, California 92008; Phone (760) 603-7200 or e-mail <b>outlicensing@lifetech.com</b> .
Limited Use Label License: Research Use Only	The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial services of any kind, including, without limitation, reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact <b>outlicensing@lifetech.com</b> or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.

#### References

- Ausubel, F. M., Brent, R., Kingston, R. E., Moore, D. D., Seidman, J. G., Smith, J. A., and Struhl, K. (1994). Current Protocols in Molecular Biology (New York: Greene Publishing Associates and Wiley-Interscience).
- Berdy, J. (1980) Bleomycin-Type Antibiotics. In *Amino Acid and Peptide Antibiotics*, J. Berdy, ed. (Boca Raton, FL: CRC Press), pp. 459-497.
- Bernard, P., and Couturier, M. (1992). Cell Killing by the F Plasmid CcdB Protein Involves Poisoning of DNA-Topoisomerase II Complexes. J. Mol. Biol. 226, 735-745.
- Bernard, P., Gabant, P., Bahassi, E. M., and Couturier, M. (1994). Positive Selection Vectors Using the F Plasmid *ccd*B Killer Gene. Gene *148*, 71-74.
- Bernard, P., Kezdy, K. E., Melderen, L. V., Steyaert, J., Wyns, L., Pato, M. L., Higgins, P. N., and Couturier, M. (1993). The F Plasmid CcdB Protein Induces Efficient ATP-dependent DNA Cleavage by Gyrase. J. Mol. Biol. 234, 534-541.
- Drocourt, D., Calmels, T. P. G., Reynes, J. P., Baron, M., and Tiraby, G. (1990). Cassettes of the *Streptoalloteichus hindustanus ble* Gene for Transformation of Lower and Higher Eukaryotes to Phleomycin Resistance. Nucleic Acids Res. *18*, 4009.
- Gatignol, A., Durand, H., and Tiraby, G. (1988). Bleomycin Resistance Conferred by a Drug-binding Protein. FEB Letters 230, 171-175.
- Innis, M. A., Gelfand, D. H., Sninsky, J. J., and White, T. S. (1990) PCR Protocols: A Guide to Methods and Applications. Academic Press, San Diego, CA.
- Jaffe, A., Ogura, T., and Hiraga, S. (1985). Effects of the *ccd* Function of the F Plasmid on Bacterial Growth. J. Bacteriol. *163*, 841-849.
- Karoui, H., Bex, F., Greze, P., and Couturier, M. (1983). *Ham22*, a Mini-F Mutation Which is Lethal to Host Cell and Promotes *rec*A-Dependent Induction of Lambdoid Prophage. EMBO J. 2, 1863-1868.
- Miki, T., Yoshioka, K., and Horiuchi, T. (1984). Control of Cell Division by Sex Factor F in *E. coli*. I. The 42.84-43.5 F Segment Couples Cell Division of the Host Bacteria with Replication of Plasmid DNA. J. Mol. Biol. *174*, 605-625.
- Ogura, T., and Hiraga, S. (1983). Mini-F Plasmid Gene That Couples Host Cell Division to Plasmid Proliferation. Proc. Natl. Acad. Sci. USA *80*, 4784-4788.
- Sambrook, J., Fritsch, E. F., and Maniatis, T. (1989). Molecular Cloning: A Laboratory Manual, Second Edition (Plainview, New York: Cold Spring Harbor Laboratory Press).
- van Melderen, L., Bernard, P., and Couturier, M. (1994). Lon-Dependent Proteolysis of CcdA is the Key Control for Activation of CcdB in Plasmid-Free Segregant Bacteria. Mol. Microbiol. *11*, 1151-1157.
- White, B. A. (1993) PCR Protocols: Current Methods and Applications. In *Methods in Molecular Biology*, Vol. 15. (J. M. Walker, ed. Humana Press, Totowa, N. J.

©2013 Life Technologies Corporation. All rights reserved.

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

#### Notes





15 March 2013