Bac-to-Bac[™] TOPO[™] Cloning Kit USER GUIDE

Five-minute cloning of blunt PCR products for expression in insect cells

Catalog Number A11098, A11099, A11100, A11101, A11338 Publication Number MAN0000698 Revision B.0



Manufacturer: Life Technologies Corporation | 5781 Van Allen Way | Carlsbad, CA 92008

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES 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	Date	Description
B.0	24 April 2018	Rebrand to Gibco. Add use of
		ExpiFectamine [™] Sf Transfection
		Reagent.
A.0	12 August 2015	Baseline for this revision.

Revision history: Pub. No. MAN0000698

Information for European customers using Mach1[™]-T1^R Cells

The Mach1[™]-T1^R *E. coli* strain is genetically modified to carry the *lac*Z∆M15 *hsd*R *lac*X74 *rec*A *end*A *ton*A genotype. As a condition of sale, this product must be in accordance with all applicable local legislation and guidelines including EC Directive 90/219/EEC on the contained use of genetically modified organisms.

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.

TRADEMARKS: All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. Triton is a trademark of Union Carbide Corporation. Falcon is a trademark of Corning, Inc.

©2018 Thermo Fisher Scientific Inc. All rights reserved.

Contents

Contents	3
Quick reference	4
Blunt-End TOPO [™] Cloning procedure for experienced users	4
Product information	5
Contents and storage	5
Description of the system	8
Experiment outline	11
Methods	
Generate blunt-end PCR products	12
Perform blunt-end TOPO [™] cloning reaction	19
Transform One Shot [™] Mach1 [™] T1 ^R Chemically Competent <i>E. coli</i>	21
Analyze positive clones	22
Next steps	23
Appendix A: Troubleshooting	25
Troubleshooting	25
Appendix B: Vectors	27
Map of pFastBac [™] /NT-TOPO [™]	27
Map of pFastBac [™] /CT-TOPO [™]	
Map of pFastBac [™] /HBM-TOPO [™]	29
Map of pFastBac [™] Gus control plasmid	
Appendix C: Support protocols	
Recipes	
One Shot [™] electroporation	
Perform the control reactions	
Control for protein expression in insect cells	
Appendix D: Ordering information	
Accessory products	
Appendix E: Safety	
Chemical safety	
Biological hazard safety	40
Documentation and support	41
Customer and technical support	41

Blunt-End TOPO[™] Cloning procedure for experienced users

Introduction	This o Cloni for th	quick reference sheet is ng procedure. If you ar e first time, follow the	provided for experienced us re performing the blunt-end detailed protocols provided i	ers of the blunt-end TOPO™ ГОРО™ Cloning procedure n the manual.	
Step			Action		
Generate PCR product	Ger you Ma liga	Generate PCR products using a thermostable proofreading DNA polymerase and your own protocol. End the PCR reaction with a final 7 to 30 minute extension step. Make sure that your PCR primers do not contain 5′ phosphates as they will inhibit ligation into your vector.			
Perform the Blunt-end TOPO [™] Cloning Reaction	1.	1. Set up one of the following blunt-end TOPO [™] Cloning reactions using the reagents in the order shown. For electroporation, dilute Salt Solution 4-fold to prepare a Dilute Salt Solution.			
		Reagent	Chemical transformation	Electroporation	
		Fresh PCR product	0.5 to 4 µL	0.5 to 4 μL	
		Salt Solution	1 µL	-	
		Dilute Salt Solution	-	1 μL	
		Sterile Water	to a final volume of 5 μ L	to a final volume of 5 μL	
		TOPO [™] Vector	1 µL	1 µL	
		Total volume	6 µL	6 µL	
		Note: The best insert:ve	ector ratio in a TOPO [™] Cloning r	reaction is 1:1 to 2:1.	
	2. 3.	Mix gently and incuba Place on ice and proce Competent <i>E. coli</i> , belo	te for 5 minutes at room tem ed to transform One Shot™ N ow.	perature. Iach1™ T1 ^ℝ Chemically	
Transform One Shot [™] Chemically Competent <i>E. coli</i>	1.	For each transformatio Competent <i>E. coli</i> on ic	n, thaw one vial of One Shot e.	[™] Mach1 [™] T1 ^R Chemically	
	2.	2. Add 2 μ L of the TOPO TM Cloning reaction into a vial of One Shot TM chemically competent <i>E. coli</i> and mix gently.			
	3.	3. Incubate the vial(s) on ice for 30 minutes.			
	4.	4. Heat-shock the cells for 30 seconds at 42°C without shaking.			
	5.	Immediately transfer t	he vial(s) to ice, and incubate	e on ice for 2 minutes.	
	6.	6. Add 250 μL of room temperature S.O.C. medium to each vial.			
	7.	7. Cap the vial(s) tightly and shake horizontally at 37°C for 1 hour (225 rpm).			
	8.	Spread 25–100 μL from containing 100 μg/mL	n each transformation on a pr ampicillin, and incubate ove	re-warmed LB agar plate ernight at 37°C.	
Control reaction	We re	ecommend using the Co	ontrol PCR Template and the	Control PCR Primers	

included with the kit to perform the control reaction. See the protocol on pages 34–35 for instructions.

Product information

Contents and storage

Types of products	This manual, supplied with the products listed below, only provides TOPO [™] cloning strategy for the direct insertion of blunt-end PCR products into pFastBac [™] TOPO [™] plasmid vectors.

For information on expressing your protein of interest from your TOPO[™] expression construct, refer to the Bac-to-Bac[™] TOPO[™] Expression System User Guide (Pub. No. MAN0000699) supplied with the Bac-to-Bac[™] N-His TOPO[™], Bac-to-Bac[™] C-His TOPO[™] Expression System, or Bac-to-Bac[™] HBM Secreted Expression System kits.

These manuals are also available at **thermofisher.com/manuals**.

Product	Quantity	Cat. No.
Bac-to-Bac™ N-His TOPO™ Cloning Kit	1 kit	A11099
Bac-to-Bac™ C-His TOPO™ Cloning Kit	1 kit	A11098
Bac-to-Bac™ HBM TOPO™ Cloning Kit	1 kit	A11338

Shipping and storage

Bac-to-Bac[™] TOPO[™] Cloning Kits are shipped on dry ice. In addition to the Bac-to-Bac[™] TOPO[™] Cloning Kit manual, each kit contains two boxes as described below. All reagents are guaranteed for six months if stored properly. Upon receipt, store boxes as detailed:

Box	Item	Storage
1	Bac-to-Bac™ N-His TOPO™, Bac-to-Bac™ C-His TOPO™,	-30°C to -10°C
	or Bac-to-Bac™ HBM TOPO™ Cloning Reagents	
2	One Shot [™] Mach1-T1 ^R Chemically Competent <i>E. coli</i>	–85°C to –68°C

Bac-to-	Bac™ TOPO™
Cloning	Kit reagents

The cloning reagents for the Bac-to-Bac[™] TOPO[™] Cloning Kits (Box 1) are listed below. **Store the contents of Box 1 at −30°C to −10°C.**

Item	Concentration	Amount
pFastBac™/NT-T0P0™ vector	20 mL at 10 ng/µL in	20 µL
(only with Cat. Nos. A11099)	50% glycerol	
or	50 mM Tris-HCl, pH 7.4 (at 25°C)	
pFastBac [™] /CT-TOPO [™] vector	1 mM EDTA	
(only with Cat. Nos. A11098)	2 mM DTT	
	0.1% Triton™ X-100	
(only with Cat. Nos. A11338)	100 μg/mL BSA	
	30 µM bromophenol blue	
10X PCR Buffer	100 mM Tris-HCl, pH 8.3 (at 42°C)	100 µL
	500 mM KCl	
	25 mM MgCl ₂	
	0.01% gelatin	
dNTP Mix	12.5 mM each dATP, dCTP, dGTP,	10 µL
	and dTTP; neutralized at pH 8.0 in	
	water	
Salt Solution	1.2 M NaCl	50 µL
	0.06 M MgCl ₂	
Sterile Water	-	1 mL
Control PCR template	50 ng/µL in TE buffer, pH 8.0*	10 µL
Control PCR primers	100 ng/µL each in TE buffer, pH	10 µL
	8.0	
Polyhedrin forward sequencing primer	100 ng/µL in TE buffer, pH 8.0	20 µL
SV40 polyA reverse sequencing	100 ng/µL in TE buffer, pH 8.0	20 µL
primer		•
pFastBac™ Gus control plasmid	0.2 ng/µL in TE buffer, pH 8.0	20 µL

*TE buffer, pH 8.0: 10 mM Tris-HCl, 1 mM EDTA, pH 8.0

Kit contents and storage, Continued

Mach1™T1^R Competent Cells

The following reagents are included in the One ShotTM Mach1TM T1^R Chemically Competent *E. coli* module (Box 2). Transformation efficiency of One ShotTM Mach1TM T1^R *E. coli* cells is $\geq 1 \times 10^9$ cfu/µg DNA. Store cells at -85°C to -68°C.

Reagent	Composition	Amount
One Shot [™] Mach1 [™] T1 ^R Chemically Competent <i>E. coli</i>	-	21 × 50 µL
S.O.C. Medium (may be stored at room temperature or 2°C to 8°C)	2% tryptone 0.5% yeast extract 10 mM NaCl 2.5 mM KCl 10 mM MgCl ₂ 10 mM MgSO4	6 mL
pUC19 Control DNA	20 mM glucose 10 pg/µL in 5 mM Tris-HCl, 0.5 mM EDTA, pH 8.0	50 µL

 $\begin{array}{lll} \mbox{Genotype of} & & \\ \mbox{Mach1}^{\tt m}\mbox{T1}^{\tt R} & & F^-\,\phi 80(\textit{lacZ})\Delta M15\;\Delta\textit{lacX74}\;\textit{hsdR}(r_{K}^-\,m_{K}^+)\;\Delta\textit{recA1398}\;\textit{endA1}\;\textit{tonA} \\ \end{array}$

IMPORTANT!The parental strain of Mach1[™]-T1^R E. coli is the non-K-12, wild-type W strain
(ATCC #9637, S. A. Waksman). Although the parental strain is generally
classified as Biosafety Level 1 (BL-1), we recommend that you consult the safety
department of your institution to verify the Biosafety Level.

Primers

Bac-to-Bac ${}^{\scriptscriptstyle\rm TM}$ TOPO ${}^{\scriptscriptstyle\rm TM}$ Cloning Kits contain the following primers to sequence your insert.

Primer	Sequence
Polyhedrin forward primer	5'-AAATGATAACCATCTCGC-3'
SV40 polyA reverse primer	5'-GGTATGGCTGATTATGATC-3'

Gus control
plasmidBac-to-Bac^T TOPO^T Cloning Kit includes the control expression plasmid
pFastBac^T Gus, which contains the Gus gene. When the recombinant baculovirus
produced from the control plasmid is used to infect insect cells, it allows the
expression of β -glucuronidase, which can be used in a rapid, qualitative assay for
expression. See page 36 for details.

Description of the system

Product featuresThe Bac-to-BacTM N-His TOPOTM, Bac-to-BacTM C-His TOPOTM, and Bac-to-BacTM HBM
TOPOTM Cloning Kits include the pFastBacTM/NT-TOPOTM, pFastBacTM/CT-TOPOTM,
and pFastBacTM/HBM-TOPOTM cloning vectors, respectively. These vectors provide
the means for the direct insertion of a blunt-end PCR product into a plasmid
vector in a highly efficient, 5-minute, one-step cloning reaction (TOPOTM Cloning).
The ability to clone blunt-end PCR products allows the use of proofreading
polymerases to amplify the gene of interest. No ligase, post-PCR procedures, or
PCR primers containing specific sequences are required.

How Topoisomerase I works Topoisomerase I from *Vaccinia* virus binds to duplex DNA at specific sites and cleaves the phosphodiester backbone after 5'-CCCTT in one strand (Shuman, 1991). The energy from the broken phosphodiester backbone is conserved by formation of a covalent bond between the 3' phosphate of the cleaved strand and a tyrosyl residue (Tyr-274) of topoisomerase I. The phospho-tyrosyl bond between the DNA and enzyme can subsequently be attacked by the 5' hydroxyl of the original cleaved strand, reversing the reaction and releasing topoisomerase (Shuman, 1994). TOPO[™] Cloning exploits this reaction to efficiently clone PCR products (see diagram below).



Continued on next page

Description of the system, Continued

TOPO [™] cloning	 The pFastBac[™]/NT-TOPO[™], pFastBac[™]/CT-TOPO[™], and pFastBac[™]/HBM-TOPO[™] cloning vectors are supplied linearized with <i>Vaccinia</i> virus DNA topoisomerase I covalently bound to the 3′ end of each DNA strand (referred to as "TOPO[™]-activated" vector). DNA topoisomerase I facilitates the cloning of bluntend PCR products encoding the gene of interest into the cloning vectors, which can then be transformed into chemically competent cells or electroporated directly into electrocompetent cells for analysis. Once the transformants are analyzed for correct orientation and reading frame, the recombinant pFastBac[™] TOPO[™] vector can be used for generating recombinant bacmid DNA to be transfected into insect cells to generate baculovirus for protein expression and further downstream applications. For more information on generating bacmid DNA and protein expression in insect cells, refer to the Bac-to-Bac[™] TOPO[™] Expression System User Guide (Pub. No. MAN0000699) available at thermofisher.com, or contact Technical Support.
Features of the pFastBac™/ NT-TOPO™ and	The pFastBac ^{m} /NT-TOPO ^{m} and pFastBac ^{m} /CT-TOPO ^{m} vectors contain the following elements. These features have been functionally tested and the vectors have been fully sequenced.
CT-TOPO [™] vectors	 Strong polyhedrin (P_H) promoter for high-level baculovirus-based protein expression in insect cells
	 TOPO[™] Cloning site for rapid and efficient cloning of blunt-end PCR products amplified with proofreading polymerases
	 N- or C-terminal polyhistidine tag (pFastBac[™]/NT-TOPO[™] and pFastBac[™]/CT-TOPO[™], respectively) for simple purification of recombinant proteins
	 TEV protease cleavage site for removal of the polyhistidine tag following protein purification using AcTEV[™] protease
	 SV40 polyadenylation signal for efficient transcription termination and polyadenylation of the recombinant transcript
	• Mini-Tn7 elements for site-specific transposition of your gene into the baculovirus shuttle vector (bacmid DNA) propagated in <i>E. coli</i>
	 Ampicillin (<i>bla</i>) resistance gene (β-lactamase) for selection of transformants in <i>E. coli</i>
	 pUC origin for high copy replication and maintenance of the plasmid in <i>E. coli</i>
	Gentamicin resistance gene for selection of transformants containing recombinant bacmid DNA
	For vector maps of pFastBac ^{TM} /NT-TOPO ^{TM} and pFastBac ^{TM} /CT-TOPO ^{TM} , see pages 27 and 28, respectively.
	Continued on next page

Description of the system, Continued

Features of the pFastBac™/	The pFastBac [™] /HBM-TOPO [™] vector contains the following elements. These features have been functionally tested and the vector has been fully sequenced.
HBM-TOP0 [™] vector	 Strong polyhedrin (P_H) promoter for high-level baculovirus-based protein expression in insect cells
	 TOPO[™] Cloning site for rapid and efficient cloning of blunt-end PCR products amplified with proofreading polymerases
	• N-terminal Honey Bee Melittin (HBM) secretion signal coding sequence for secretion of the cloned gene product into the extracellular medium
	• C-terminal polyhistidine tag for simple purification of recombinant proteins
	 TEV protease cleavage site for removal of the polyhistidine tag following protein purification using AcTEV[™] protease
	 SV40 polyadenylation signal for efficient transcription termination and polyadenylation of the recombinant transcript
	• Mini-Tn7 elements for site-specific transposition of your gene into the baculovirus shuttle vector (bacmid DNA) propagated in <i>E. coli</i>
	• Ampicillin (<i>bla</i>) resistance gene (β-lactamase) for selection of transformants in <i>E. coli</i>
	 pUC origin for high copy replication and maintenance of the plasmid in E. coli
	 Gentamicin resistance gene for selection of transformants containing recombinant bacmid DNA

For the vector map of pFastBac[™]/HBM-TOPO[™], see page 29.

Experiment outline

Experiment outline To TOPO[™] Clone your gene of interest into pFastBac[™] TOPO[™] vectors, perform the following steps:

- 1. Generate a PCR product containing your gene of interest with a thermostable proofreading DNA polymerase (e.g., AccuPrime[™] *Pfx* DNA Polymerase, Platinum[™] *Pfx* DNA Polymerase).
- 2. TOPO[™] Clone your blunt-end PCR product into the pFastBac[™]/NT-TOPO[™], pFastBac[™]/CT-TOPO[™], or pFastBac[™]/HBM-TOPO[™] vector, and use the reaction to transform One Shot[™] Mach1[™] T1[®] Chemically Competent *E. coli*.
- 3. Pick colonies, isolate plasmid DNA, and screen for insert directionality by sequencing expression clones with primers provided in the kit.



Methods

Generate blunt-end PCR products

Introduction	This kit is specifically designed to clone blunt-end PCR products generated by thermostable proofreading polymerases such as AccuPrime ^{TM} <i>Pfx</i> DNA Polymerase and Platinum ^{TM} <i>Pfx</i> DNA Polymerase. Follow the guidelines below to design your PCR primers and to produce your blunt-end PCR product.	
Note	Do not add 5´ phosphates to your primers for PCR. The PCR product synthesized will not ligate into pFastBac [™] TOPO [™] vectors.	
Considerations for pFastBac™/ NT-TOPO™	The cloning of a blunt-end PCR product into a pFastBac [™] /NT-TOPO [™] vector is a rapid and efficient process. However, to ensure proper expression of your recombinant protein, pay attention to the general considerations outlined below:	
	• The pFastBac [™] /NT-TOPO [™] vector contains the ATG start codon immediately upstream of the N-terminal polyhistidine tag; therefore, it is not necessary to include the initiation codon when designing your insert. However, your insert may include an ATG. Since the ribosome generally recognizes the first ATG, unless the internal ATG is in a particularly good context, the ATG in front of the polyhistidine will be used for initiation.	
	• It is best to eliminate the untranslated leader sequence from your insert and have the ATG start codon as close to the polyhedrin promoter as possible.	
	Your insert must contain a stop codon.	
	 Do not include the Kozak sequence in the insert cloned into the pFastBac[™]/NT-TOPO[™] vector, because this sequence is not required for translation initiation in insect cells. 	
	• If the gene of interest contains a polyadenylation signal, the first signal from the gene is recognized and the second signal in the vector is ignored. This does not affect expression.	
	• The cloning step presents the only limitation to the size of the insert. While the baculovirus genome can accommodate inserts of considerable size, large plasmids are more difficult to transform into <i>E. coli</i> .	
	Continued on next page	

Cloning site of
pFastBac™/The vector sequence is available at thermofisher.com or by contacting Technical
Support.NT-TOPO™ vectorSupport.



Considerations for pFastBac™/ CT-TOPO™	The cloning of a blunt-end PCR product into a pFastBac [™] /CT-TOPO [™] vector is a rapid and efficient process. However, to ensure proper expression of your recombinant protein, it is important to pay attention to the general considerations outlined below:		
	• When using the pFastBac [™] /CT-TOPO [™] vector, your insert must contain the ATG initiation codon.		
	• It is best to eliminate the untranslated leader sequence from your insert, and have the ATG start codon as close to the polyhedrin promoter as possible.		
	• It is not necessary to include the Kozak sequence in the insert cloned into the pFastBac [™] /CT-TOPO [™] vector; this sequence is not required for translation initiation in insect cells. However, in some cases, the Kozak sequence acts as an enhancer.		
	• Do not include a stop codon in the reverse primer for PCR if you want to use the C-terminal polyhistidine tag, because the pFastBac [™] /CT-TOPO [™] vector contains a stop codon immediately downstream of the C-terminal polyhistidine tag.		
	• If the gene of interest contains a polyadenylation signal, the first signal from the gene is recognized and the second signal in the vector is ignored. This does not affect expression.		
	• You can design your reverse primer to include a stop codon to omit the C-terminal tag encoded by the pFastBac [™] /CT-TOPO [™] vector, and express your protein in its native state.		
	• Similarly, you can design your PCR product to encode a different C-terminal tag followed by a stop codon to eliminate the pFastBac [™] /CT-TOPO [™] polyhistidine tag.		
	• The cloning step presents the only limitation to the size of the insert. While the baculovirus genome can accommodate inserts of considerable size, large plasmids are more difficult to transform into <i>E. coli</i> .		

Cloning site of
pFastBac™/The vector sequence is available at thermofisher.com or by contacting Technical
Support.CT-TOPO™ vector



Considerations for pFastBac™/ HBM-T0P0™	The cloning of a blunt-end PCR product into a pFastBac [™] /HBM-TOPO [™] vector is a rapid and efficient process. However, to ensure proper expression of your recombinant protein, it is important to pay attention to the general considerations outlined below:		
	• The pFastBac [™] /HBM-TOPO [™] vector contains the ATG start codon immediately upstream of the N-terminal HBM secretion signal sequence; therefore it is not necessary to include the initiation codon when designing your insert. However, your insert may include an ATG. Since the ribosome generally recognizes the first ATG, unless the internal ATG is in a particularly good context, the ATG in front of the polyhistidine will be used for initiation.		
	• Do not include the Kozak sequence in the insert cloned into the pFastBac [™] /HBM-TOPO [™] vector; this sequence is not required for translation initiation in insect cells.		
	• Do not include a stop codon in the reverse primer for PCR if you want to use the C-terminal polyhistidine tag, because the pFastBac [™] /HBM-TOPO [™] vector contains a stop codon immediately downstream of the C-terminal polyhistidine tag.		
	• If the gene of interest contains a polyadenylation signal, the first signal from the gene is recognized and the second signal in the vector is ignored. This does not affect expression.		
	• You can design your reverse primer to include a stop codon to omit the C- terminal tag encoded by the pFastBac [™] /HBM-TOPO [™] vector, and express your protein in its native state.		
	• Similarly, you can design your PCR product to encode a different C-terminal tag followed by a stop codon to eliminate the pFastBac [™] /HBM-TOPO [™] polyhistidine tag.		
	• The cloning step presents the only limitation to the size of the insert. While the baculovirus genome can accommodate inserts of considerable size, large plasmids are more difficult to transform into <i>E. coli</i> .		

Cloning site of
pFastBac™/The vector sequence is available at thermofisher.com or by contacting Technical
Support.HBM-TOPO™ vector

	Polyhedrin forv primer binding	vard site Polyhedrin promoter Peu
001	ATCATGGAGA TAATTAAAAT GATAACCATC TAGTACCTCT ATTAATTTA CTATTGGTAG	C TCGCAAATAA ATAAGTATTT TACTGTTTTC GTAACAGTTT G AGCGTTTATT TATTCATAAA ATGACAAAAG CATTGTCAAA
	Polyhedrin p	promoter P _{PH}
071	TGTAATAAAA AAACCTATAA ATATTCCGGA ACATTATTTT TTTGGATATT TATAAGGCCT	A TTATTCATAC CGTCCCACCA TCGGGCGCGG ATCCACCGGT F AATAAGTATG GCAGGGTGGT AGCCCGCGCC TAGGTGGCCA
	HBM see	cretion signal sequence
141	ATGAAATTCT TAGTCAACGT TGCCCTTGTT TACTTTAAGA ATCAGTTGCA ACGGGAACAA	T TTTATGGTCG TATACATTTC TTACATCTAT GCGGATCGAT A AAATACCAGC ATATGTAAAG AATGTAGATA CGCCTAGCTA
	ATG Start Codon	Melittin cleavage site TEV recognition
211	Blunt PCR product	AAG GGC GAA AAC TTG TAC TTT CAA GGC CAT CAC TTC CCG CTT TTG AAC ATG AAA GTT CCG GTA GTG Lys Gly Glu Asn Leu Tyr Phe Gln Gly His His
	6x His	AcTEV™ cleavage site
243	CAT CAC CAT CAC TAG CTC GAG GCA GTA GTG GTA GTG ATC GAG CTC CGT His His His His *** Leu Glu Ala	TGCGGTACCA AGCTTGTCGA GAAGTACTAG AGGATCATAA ACGCCATGGT TCGAACAGCT CTTCATGATC TCCTAGTATT
		nylation signal ————
313	TCAGCCATAC CACATTTGTA GAGGTTTTAC AGTCGGTATG GTGTAAACAT CTCCAAAATG -SV40 reverse primer binding site	TTGCTTTAAA AAACCTCCCA CACCTCCCCC TGAACCTGAA AACGAAATTT TTTGGAGGGT GTGGAGGGGG ACTTGGACTT

PCR reaction	After you have designed primers to amplify your gene of interest, you are ready to generate your PCR product for TOPO TM Cloning into pFastBac TM TOPO TM vectors. Note: You must use a thermostable proofreading DNA polymerase such as Platinum TM <i>Pfx</i> DNA Polymerase or AccuPrime TM <i>Pfx</i> DNA Polymerase to produce your blunt-end PCR product. <i>Taq</i> Polymerase has a terminal transferase activity that adds a single 3'-A overhang to each end of the PCR product, thus rendering it unsuitable for blunt-end TOPO TM Cloning.	
Required materials	 Thermostable proofreading polymerase 10X PCR buffer appropriate for your polymerase Thermocycler DNA template and primers for your PCR product Note: dNTPs (adjusted to pH 8) are provided in the kit. 	
Generate blunt-end PCR products	 Set up a 25 µL or 50 µL PCR reaction using the guidelines: Follow the instructions and recommendations provided by the manufacturer of your thermostable proofreading polymerase to produce blunt-end PCR products. Use the cycling parameters suitable for your primers and template. Make sure to optimize PCR conditions to produce a single, discrete PCR product. Use a 7–30 minute final extension to ensure that all PCR products are completely extended. After completing the PCR reaction, place the tube on ice or store at -20°C for up to 2 weeks. 	
Check the PCR product	After you have produced your blunt-end PCR product, use agarose gel electrophoresis to verify the quality and quantity of your PCR product. Be sure you have a single, discrete band of the correct size. If you do not have a single, discrete band, follow the manufacturer's recommendations for optimizing your PCR with the polymerase of your choice. Alternatively, you may gel-purify the desired product using the PureLink [™] Quick Gel Extraction Kit, available separately. We also offer the E-Gel [™] CloneWell [™] SYBR [™] Safe gels, which allow the isolation of DNA bands without any additional gel purification steps. For ordering information, See page 37.	

Perform blunt-end TOPO[™] cloning reaction

Introduction	After you have produced the desired PCR product, you are ready to $TOPO^{\mathbb{M}}$ Clone your blunt-end insert into the pFastBac ^{\mathbb{M}} TOPO ^{\mathbb{M}} vector (pFastBac ^{\mathbb{M}} /NT- TOPO ^{\mathbb{M}} , CT-TOPO ^{\mathbb{M}} , or HBM-TOPO ^{\mathbb{M}}), and use the recombinant vector to transform competent <i>E. coli</i> . It is important to have everything you need to set up the reaction so that you can obtain the best results. We suggest that you read this entire section and the next section about transformation before beginning.
	Note: If this is the first time you are TOPO [™] Cloning, perform the control reactions detailed on pages 34 and 35 in parallel with your samples.
Salt in the TOPO™ Cloning reaction	Perform TOPO [™] Cloning in a reaction buffer containing salt (i.e., using the stock salt solution provided in the kit).
	Note: The amount of salt added to the TOPO [™] Cloning reaction varies depending on whether you plan to transform chemically competent cells (provided) or electrocompetent cells.
	• If you are transforming One Shot [™] Mach1 [™] T1 ^R Chemically Competent <i>E. coli</i> (included in the kit), use the stock Salt Solution as supplied, and set up the TOPO [™] Cloning reaction as directed on the next page.
	 If you are transforming electrocompetent <i>E. coli</i> (available separately; see page 37), the amount of salt in the TOPO[™] Cloning reaction must be reduced to 50 mM NaCl, 2.5 mM MgCl₂ to prevent arcing during electroporation. Dilute the stock Salt Solution 4-fold with water to prepare a 300 mM NaCl, 15 mM MgCl₂ Dilute Salt Solution. Use the Dilute Salt Solution to set up the TOPO[™] Cloning reaction as directed on the upcoming section.
Note	We have found that including salt (200 mM NaCl, 10 mM MgCl₂) in the TOPO [™] Cloning reaction increases the number of transformants 2- to 3-fold. In addition, incubating the reaction mixture for greater than 5 minutes in the presence of salt can also increase the number of transformants. In experiments without salt, the number of transformants decreases as the incubation time increases beyond 5 minutes.
	Including salt in the TOPO [™] Cloning reaction allows for longer incubation times because it prevents topoisomerase I from rebinding and potentially nicking the DNA after ligating the PCR product and dissociating from the DNA. The result is more intact molecules leading to higher transformation efficiencies.
	Continued on next page

Perform blunt-end TOPO[™] cloning reaction, Continued

Final volume

Required materials	Your PCR product (freshly prepared) pFastBac [™] /NT-TOPO [™] , pFastBac [™] /CT-TOPO [™] , or pFastBac [™] /HBM-TOPO [™] vector Salt Solution or Dilute Salt Solution Sterile Water		
Perform the TOPO™ Cloning reaction	 Set up your TOPO[™] Cloning reaction (6 µL) as described in the table below. The best insert:vector ratio in a TOPO[™] Cloning reaction is 1:1 to 2:1. Note: The blue color of the TOPO[™] vector solution is normal. 		
	Reagent	Chemically Competent <i>E. coli</i>	Electrocompetent <i>E. coli</i>
	PCR Product	0.5 to 4 µL	0.5 to 4 µL
	Salt Solution	1 µL	-
	Dilute Salt Solution	-	1μL
	Sterile Water	To a total volume of 5 μL	To a total volume of 5 μL
	pFastBac™ T0P0™ vector	1 µL	1μL

Note: Store all reagents at -20° C when finished. Salt solution and water can be stored at room temperature or 4° C.

6 µL

6μL

2. Mix the reaction gently, and incubate it for 5 minutes at room temperature (22°C–23°C).

Note: For most applications, 5 minutes yields a sufficient number of colonies for analysis. The length of the TOPO[™] Cloning reaction can be increased from 30 seconds to several hours. For routine subcloning of PCR products, 30 seconds may be sufficient. For larger PCR products (>3 kb), increasing the reaction time, may yield more colonies.

3. Place the reaction on ice and proceed to transformation.

Note: You may store the TOPOTM Cloning reaction overnight at -20° C.

Transform One Shot[™] Mach1[™] T1^R Chemically Competent *E. coli*

Introduction	Or to Ch als de Aj Pro sec	Ince you have performed the Blunt-End TOPO TM Cloning reaction, you are ready of use your construct to transform competent <i>E. coli</i> . One Shot TM Mach1 TM T1 ^R Themically Competent <i>E. coli</i> cells are included with the kit (Box 2). You may also transform electrocompetent cells (for ordering information, see page 37) if esired following the One ShotTM electroporation protocol provided in the ppendix (see page 32). rotocols for transforming chemically competent <i>E. coli</i> are provided in this ection. For instructions on performing control reactions, see page 34. ote: Do not use One Shot TM Mach1 TM T1 ^R Chemically Competent <i>E. coli</i> for electroporation.			
IMPORTANT!	If y ess gro	rou are transforming One Shot [™] Mach1 [™] -T1 ^R Chemically Competent <i>E. coli</i> , it is ential that selective plates are pre-warmed to 37° prior to spreading for optimal owth of cells.			
Required materials	•	TOPO [™] Cloning reaction from Perform the TOPO[™] Cloning reaction , Step 2 (page 20)			
	•	S.O.C. medium at room temperature			
	•	42°C water bath			
	•	LB plates containing 100 μ g/mL ampicillin, pre-warmed to 37°C			
	٠	37°C shaking and non-shaking incubator			
	•	Competent cells (thawed on ice)			
One Shot [™] Mach1 [™] T1 [®] Chemical	1.	Thaw on ice one vial of One Shot ^{M} Mach1 ^{M} T1 ^{R} Chemically Competent <i>E. coli</i> for each transformation.			
Transformation	2.	Add 2 µL of the TOPO [™] Cloning reaction (Step 2, page 20) into a vial of One Shot [™] Mach1 [™] T1 ^ℝ Chemically Competent <i>E. coli</i> and mix gently. IMPORTANT! Do not mix by pipetting up and down.			
	3.	Incubate the vial(s) on ice for 30 minutes.			
	4.	Heat-shock the cells for 30 seconds at 42°C without shaking.			
	5.	Immediately transfer the vial(s) to ice, and incubate them on ice for 2 minutes.			
	6.	Add 250 µL of room temperature S.O.C. medium to each vial.			
	7.	Cap the vial(s) tightly and shake them horizontally at 37°C for 1 hour (225 rpm).			
	8.	Spread 25–100 μ L from each transformation on a pre-warmed selective plate. We recommend plating two different volumes to ensure that at least one plate will have well-spaced colonies.			
	9.	Invert the plate(s) and incubate at 37°C. With ampicillin selection, visible colonies should appear within 8 hours.			
	10.	Pick ~10 colonies for analysis (see Analyze positive clones , page 22). An efficient TOPO [™] Cloning reaction produces several hundred colonies.			

Analyze positive clones

Introduction	After transforming your pFastBac ^{TM} /NT-TOPO ^{TM} , pFastBac ^{TM} /CT-TOPO ^{TM} , or pFastBac ^{TM} /HBM-TOPO ^{TM} construct into <i>E. coli</i> , select and sequence several colonies using the specific primers included in the kit to determine the orientation of the insert.		
Analyze positive clones	 Pick 10 overnight-grown colonies from the selective plates and culture them overnight in LB medium containing 100 μg/mL ampicillin. Isolate plasmid DNA using your method of choice. If you need ultra-pure plasmid DNA for automated or manual sequencing, we recommend using 		
	 the PureLink[™] HiPure Mini Plasmid Purification Kit (see page 37 for ordering information). Analyze plasmid DNA by sequencing. 		
Note	If you have used One Shot [™] Mach1 [™] T1 ^R Chemically Competent <i>E. coli</i> for your transformation, you can prepare plasmid DNA 4 hours after inoculating a single, overnight-grown colony in the selective media of choice. Note that this feature is not limited to ampicillin selection.		
Sequencing	To confirm that your gene of interest is in the correct orientation, you may sequence your expression construct using the Polyhedrin forward and SV40 polyA reverse primers included with the kit. Refer to page 7 for the sequences of the primers and pages 13, 15, and 17 for the location of the primer binding sites of pFastBac [™] /NT-TOPO [™] , pFastBac [™] /CT-TOPO [™] , and pFastBac [™] /HBM-TOPO [™] , respectively.		
PCR analyze transformants	You may also determine the orientation of your insert by PCR amplification using a pair of primers, where one primer binds outside the TOPO [™] Cloning site (e.g., polyhedrin forward primer), while the other is internal to your blunt-end PCR insert. The PCR product, the size of which will depend on the orientation of the insert, can then easily be visualized on an agarose gel.		
Long-term storage	 Once you have identified the correct clone, be sure to purify the colony and make a glycerol stock for long-term storage. Streak the original colony out for single colonies on an LB plate containing 100 μg/mL ampicillin. 		
	2. Isolate a single colony and inoculate into 1–2 mL of LB containing $100 \ \mu g/mL$ ampicillin.		
	3. Grow at 37°C with shaking until culture reaches stationary phase.		
	4. Mix 0.85 mL of culture with 0.15 mL of sterile glycerol.		
	5. Vortex and transfer to a labeled cryovial.		
	6. Freeze the tube in liquid nitrogen or dry ice/ethanol bath and store at -80° C.		
	We also recommend that you store a stock of plasmid DNA at -20° C.		

Introduction	 After you obtain the correct pFastBac[™]/NT-TOPO[™], pFastBac[™]/CT-TOPO[™], or pFastBac[™]/HBM-TOPO[™] recombinant plasmid construct, create a recombinant bacmid to transfect into your insect cell line of choice to create a recombinant baculovirus. After amplifying and titering the baculovirus stock, you will be ready to use this stock to infect insect cells to express your protein of interest. For more information on generating bacmid DNA and intracellular protein expression in insect cells, refer to the Bac-to-Bac[™] TOPO[™] Expression System User Guide (Pub. No. MAN0000699). These manuals are available at thermofisher.com/manuals.
Bac-to-Bac™ TOPO™ Expression System	 The Bac-to-Bac[™] TOPO[™] Expression System provides a rapid and efficient method of generating recombinant baculoviruses, based on site-specific transposition of an expression cassette into a baculovirus shuttle vector (bacmid) propagated in <i>E. coli</i>. The major components of the Bac-to-Bac[™] TOPO[™] Expression System include: pFastBac[™]/NT-TOPO[™] or pFastBac[™]/CT-TOPO[™] plasmid that allows generation of an expression construct containing the gene of interest. A competent <i>E. coli</i> host strain, MAX Efficiency[™] DH10Bac[™], that contains a baculovirus shuttle vector (bacmid) and a helper plasmid, and that allows generation of a recombinant bacmid following transposition of your pFastBac[™] TOPO[™] expression construct.
	 pFastBac[™] Gus control expression plasmid that allows production of a recombinant baculovirus which, when used to infect insect cells, expresses β-glucuronidase.
	• A next-generation transfection reagent, ExpiFectamine [™] Sf, which provides high efficiency gene expression and baculovirus production with minimal cytotoxicity across adherent and suspension insect cell cultures. This versatile reagent enables the use of two flexible transfection protocols to simplify your workflow (see thermofisher.com/ExpiSf for more information).

Bac-to-Bac[™] HBM TOPO[™] Secreted Expression System The Bac-to-Bac[™] HBM TOPO[™] Secreted Expression System allows the rapid generation of recombinant baculoviruses for secreted protein expression. It contains the following major components:

- pFastBac[™]/HBM-TOPO[™] plasmid that allows generation of an expression construct containing the gene of interest in frame with the Honey Bee Melittin (HBM) secretion signal coding sequence for secretion of the cloned gene product into the extracellular medium
- A competent *E. coli* host strain, MAX Efficiency[™] DH10Bac[™], that contains a baculovirus shuttle vector (bacmid) and a helper plasmid, and that allows generation of a recombinant bacmid following transposition of your pFastBac[™]/HBM-TOPO[™] expression construct.
- pFastBac[™] Gus control expression plasmid that allows production of a recombinant baculovirus which, when used to infect insect cells, expresses β-glucuronidase.
- A next-generation transfection reagent, ExpiFectamine[™] Sf, which provides high efficiency gene expression and baculovirus production with minimal cytotoxicity across adherent and suspension insect cell cultures. This versatile reagent enables the use of two flexible transfection protocols to simply your workflow (see **thermofisher.com/ExpiSf** for more information).

Appendix A: Troubleshooting

Troubleshooting

Introduction

This table lists some potential problems solutions that may help you troubleshoot your TOPO[™] Cloning and expression of your gene of interest.

Observation	Possible cause	Action
Few or no colonies obtained from sample reaction, but transformation control yielded colonies	Incomplete extension during PCR	Include a final extension step of 7–30 minutes during PCR. Longer PCR products need a longer extension time.
	Excess or dilute PCR product used in the TOPO™ Cloning reaction	Reduce or concentrate the amount of PCR product.
	PCR primers contain 5' phosphates	Do not add 5' phosphates to your PCR primers.
	Large PCR product	 Increase the amount of PCR product used in the TOPO™ Cloning reaction. Increase the incubation time of TOPO™ Cloning reaction from 5 minutes to 30, 60, or 120 minutes. Or incubate overnight. Gel-purify the PCR product to remove primer-dimers or other artifacts.
	PCR reaction contains artifacts (i.e., not a single band on an agarose gel)	 Optimize your PCR conditions. Gel-purify your PCR product.
	PCR product contains 3' A-overhangs because you used <i>Taq</i> polymerase	Use a thermostable proofreading DNA polymerase such as Platinum [™] <i>Pfx</i> DNA Polymerase or AccuPrime [™] <i>Pfx</i> DNA Polymerase to produce your blunt-end PCR product.

Troubleshooting, Continued

Observation	Possible cause	Action
Large number of incorrect inserts cloned	PCR cloning artifacts	 Gel-purify your PCR product to remove primer-dimers and other artifacts. Optimize your PCR conditions.
		 Include a final extension step of 7–30 minutes during PCR.
High background and large number of satellite colonies after transformation.	Recovery period after transformation too long.	Reduce incubation period after transformation from 1 hour to 5–10 minutes at 37°C (225 rpm).
Few or no colonies obtained from sample reaction and the	One Shot [™] competent <i>E. coli</i> stored incorrectly	 Store One Shot[™] competent <i>E. coli</i> at -80°C.
transformation control gave no colonies		 If you are using another <i>E. coli</i> strain, follow the manufacturer's instructions.
	Insufficient amount of <i>E. coli</i> plated	Increase the amount of <i>E. coli</i> plated.
	Transformants plated on selective plates with the wrong antibiotic	Use LB ampicillin plates for selection.

Map of pFastBac[™]/NT-TOPO[™]

Description

The map of the elements of the pFastBac[™]/NT-TOPO[™] vector. The vector sequence is available from **thermofisher.com** or by contacting Technical Support.



Comments for pFastBac/NT-TOPO vector 4762 nucleotides

Polyhedrin promoter (P_{PH}): bases 1-129 Initiation ATG: bases 147-149 6xHis tag: bases 159-179 TEV recognition site: bases 189-209 TOPO cloning site: bases 212-213 SV40 polyadenylation signal: bases 262-502 Tn7L: bases 531-696 f1 origin: bases 880-1334 Ampicillin resistance gene: bases 1465-2325 pUC origin: bases 2470-3143 Tn7R: bases 3389-3613 Gentamicin resistance gene: bases 3680-4208 (complementary strand)

Map of pFastBac[™]/CT-TOPO[™]

Description

The map of the elements of pFastBac[™]/CT-TOPO[™] vector. The vector sequence is available from **thermofisher.com** or by contacting Technical Support.



Comments for pFastBac™/CT-TOPO™ vector 4754 nucleotides

Polyhedrin promoter (P_{PH}): bases 1-129 TOPO cloning site: bases 145-146 TEV recognition site: bases 152-169 6xHis tag: bases 173-190 SV40 polyadenylation signal: bases 235-475 Tn7L: bases 504-669 f1 origin: bases 853-1307 Ampicillin resistance gene: bases 1438-2298 pUC origin: bases 2443-3116 Tn7R: bases 3362-3586 Gentamicin resistance gene: bases 3653-4186 (complementary strand)

Map of pFastBac[™]/HBM-T0P0[™]

Description The map of the elements of $pFastBac^{T}/HBM$ -TOPOTM vector. The vector sequence is available from **thermofisher.com** or by contacting Technical Support.



Comments for pFastBac[™]/HBM-TOPO[™] vector 4824 nucleotides

Polyhedrin promoter (P_{PH}): bases 1-129 Honey Bee Mellitin (HBM) secretion signal: 141-210 TOPO cloning site: bases 215-216 TEV recognition site: bases 222-242 6xHis tag: bases 243-260 SV40 polyadenylation signal: bases 305-545 Tn7L: bases 574-739 f1 origin: bases 923-1377 Ampicillin resistance gene: bases 1508-2368 pUC origin: bases 2513-3186 Tn7R: bases 3432-3656 Gentamicin resistance gene: bases 3723-4256 (complementary strand)

Map of pFastBac[™] Gus control plasmid

Description

The pFastBacTM Gus plasmid is a 6,661 bp control vector that contains the *Arabidopsis thaliana* gene for β -glucuronidase (Gus). The molecular weight of β -glucuronidase is 68.5 kDa.

The map of the elements of pFastBac[™] Gus control plasmid. The vector sequence is available from **thermofisher.com** or by contacting Technical Support.



Comments for pFastBac[™] Gus vector 6661 nucleotides

f1 origin: bases 2-457 Ampicillin resistance gene: bases 589-1449 pUC origin: bases 1594-2267 Tn7R: bases 2511-2735 Gentamicin resistance gene: bases 2802-3335 (complementary strand) Polyhedrin promoter (P_{PH}): bases 3904-4032 GUS ORF: bases 4081-5892 SV40 polyadenylation signal: bases 6047-6287 Tn7L: bases 6315-6480

Appendix C: Support protocols

Recipes				
Pre-mixed media	We carry pre-mixed growth media, such as imMedia ^{M} , in convenient pouches of in bulk. imMedia ^{M} is pre-mixed and pre-sterilized for convenient preparation or liquid medium or agar plates for <i>E. coli</i> growth, and it is available with or without IPTG and X-gal and a choice of three antibiotics: ampicillin, kanamycir or Zeocin ^{M} selection agent. For ordering information, see page 37.			
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.			
	3. Autoclave on liquid cycle for 20 minutes at 15 psi. Allow solution to cool to 55°C and add antibiotic if needed.			
	4. Store at room temperature or at 4°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, and pour into 10 cm plates.			
	4. Let harden, then invert and store at 4°C.			

One Shot[™] electroporation

Introduction	Although chemical transformation of <i>E. coli</i> is the most convenient method, electroporation is more efficient, and it is the method of choice for large plasmids. This section provides an optional protocol if you prefer to use electroporation. We offer a variety of high transformation-efficiency electrocompetent cells that are suitable for use with your recombinant pFastBac TM TOPO TM vector. For more information on electrocompetent <i>E. coli</i> cells, refer to thermofisher.com or contact Technical Support (page 37)			
	This section provides a protocol for transforming One Shot ^{TM} Electrocomp ^{TM} <i>E. coli</i> with your recombinant pFastBac ^{TM} TOPO ^{TM} vector.			
IMPORTANT!	Do not use One Shot TM Mach1 TM T1 ^R Chemically Competent <i>E. coli</i> for electroporation.			
Required materials	In addition to general microbiological supplies (e.g., plates, spreaders), you will need the following reagents and equipment.			
	• TOPO [™] Cloning reaction from Perform the TOPO[™] Cloning reaction , Step 2 (page 20)			
	Electroporator			
	• S.O.C. medium at room temperature			
	• LB plates containing 100 μg/mL ampicillin, pre-warmed to 37°C.			
	• Electrocompetent cells (thawed on ice)			
	Continued on next page			

One Shot[™] **Electroporation,** Continued

One Shot™ electroporation	1.	Add 2 μ L of the TOPO TM Cloning reaction to a vial (50 μ L) of One Shot TM Electrocomp TM <i>E. coli</i> and mix gently.			
protocol		Note: Do not mix by pipetting up and down. Avoid formation of bubbles.			
	2.	Carefully transfer cells and DNA to a chilled 0.1 cm cuvette.			
	3.	Electroporate your samples using your own protocol and electroporator.			
		Note: If you have problems with arcing, see the next page.			
	4.	Immediately add 250 µL of room temperature S.O.C. medium to the cuvette.			
	5.	Transfer the solution to a 15 mL snap-cap tube (e.g., Falcon [™]) and shake for at least 1 hour at 37°C to allow expression of the antibiotic resistance genes.			
	6.	Spread 10–50 μ L from each transformation on a pre-warmed selective plate and incubate overnight at 37°C. To ensure even spreading of small volumes add 20 μ L of S.O.C. medium. We recommend that you plate two different volumes to ensure that at least one plate will have well-spaced colonies. Incubate plates overnight at 37°C.			
	7.	Pick ~10 colonies for analysis (see Analyze positive clones , page 22). An efficient TOPO [™] Cloning reaction produces several hundred colonies.			
Note	Ad fin and ele cuv	Addition of the Dilute Salt Solution in the TOPO TM Cloning reaction brings the final concentration of NaCl and MgCl ₂ in the TOPO TM Cloning reaction to 50 mM and 2.5 mM, respectively. To prevent arcing of your samples during electroporation, the volume of cells should be between 50 µL and 80 µL (0.1 cm cuvettes) or 100 µL to 200 µL (0.2 cm cuvettes).			
	If y sug	If you experience arcing during transformation, try one of the following suggestions:			
	•	Reduce the voltage normally used to charge your electroporator by 10%			
	٠	Reduce the pulse length by lowering the load resistance to 100 ohms			
	•	Precipitate the TOPO [™] Cloning reaction and resuspend in water prior to electroporation			

Perform the control reactions

Introduction

We recommend performing the following control TOPO[™] Cloning reactions the first time you use the kit to help you evaluate your results. Performing the control reactions involves generating a 750 bp control PCR product and using the PCR product directly in a TOPO[™] Cloning reaction.

Produce the control 1. Set up the following reaction in a 50 μ L volume.

PCR product

Reagent	Amount
Control PCR Template	1 µL
10X PCR Buffer	5 µL
50 mM dNTPs	0.5 µL
Control PCR Primers (0.1 µg/µL each)	1 µL each
Sterile Water	40.5 μL
Thermostable Proofreading DNA Polymerase	1 µL
(e.g., Platinum™ <i>Pfx</i> DNA Polymerase)	
Total Volume	50 µL

2. Amplify the control PCR product using the following cycling parameters:

Step	Time	Temp.	Cycles
Initial Denaturation	2 min	94°C	1X
Denaturation	1 min	94°C	
Annealing	1 min	55°C	25X
Extension	1 min	72°C	
Final Extension	7 min	72°C	1X

3. Remove 10 µL from the reaction and analyze by gel electrophoresis on a 0.8% agarose gel. A discrete 750 bp band should be visible.

Perform the control reactions, Continued

Control TOPO™ Cloning reactions	1.	 Using the control PCR product generated in the steps above and the controvector, set up two 6 µL TOPO[™] Cloning reactions as described: 				
		Reagent	Vector only	Vector + PCR Insert		
		Control PCR Product	-	1 µL		
		Sterile Water	4 µL	3 µL		
		Salt Solution or Dilute Salt Solution	1 µL	1 µL		
		pFastBac [™] /NT-TOPO [™] , pFastBac [™] /CT-TOPO [™] , or pFastBac [™] /HBM-TOPO [™]	1 µL	1 µL		
	2.	2. Incubate the reactions at room temperature for 5 minutes, and place them on ice.				
	3.	 Use 2 µL of the reaction to transform two separate vials of One Shot[™] competent cells using the procedure on page 21. 				
	4. Spread 10–50 μ L of each transformation mix onto LB plates containing 100 μ g/mL ampicillin. When plating small volumes, add 20 μ L of S.O.C. medium to ensure even spreading. Be sure to plate two different volumes to ensure that at least one plate has well-spaced colonies.					
	5. Incubate plates overnight at 37°C.					
What you should see	The vector + PCR insert reaction should produce hundreds of colonies. 95% of these colonies should contain the 750 bp insert when analyzed by restriction digestion and agarose gel electrophoresis.					
	The vector-only reaction should yield very few colonies (<15% of the vector + PCR insert plate).					
Transformation control	The pUC19 plasmid is included to check the transformation efficiency of the One Shot ^{TM} Mach1 ^{TM} T1 ^R chemically competent cells.					
	1.	using the protocol on page 21.				
	2.	Plate 10 μ L of the transformatic containing 100 μ g/mL ampicil	on reaction plus 20 μL lin.	of S.O.C. on LB plates		
	The	e transformation efficiency shou	lld be 1×10^9 cfu/µg D	NA.		

Control for protein expression in insect cells

Assay for B-glucuronidase	If you include the pFastBac TM Gus baculoviral control construct in your expression experiment, you may assay for β -glucuronidase expression. To assess β -glucuronidase expression in a rapid manner, mix a small amount of media from the infected cells with the chromogenic indicator X-glucuronide, and observe the development of blue color.		
	1.	Mix 5 μ L of 20 mg/mL X-glucuronide solution (in DMSO or dimethylformamide) with 50 μ L of cell-free medium.	
	2.	Monitor for development of blue color within 2 hours.	

Note: Other methods are also suitable.

Accessory products

Additional products

The table below lists additional products that may be used with Bac-to-Bac[™] TOPO[™] Cloning Kits. For more information, refer to **thermofisher.com** or contact Technical Support.

Product	Amount	Cat. no.
Platinum™ <i>Pfx</i> DNA Polymerase	100 units 250 units 500 units	11708-013 11708-021 11708-039
AccuPrime [™] <i>Pfx</i> DNA Polymerase	200 reactions 1000 reactions	12344-024 12344-032
<i>Pfx50</i> [™] DNA Polymerase	100 reactions 500 reactions	12355-012 12355-036
One Shot™ TOP10 Chemically Competent <i>E. coli</i>	10 reactions 20 reactions	C4040-10 C4040-03
One Shot [™] Mach1 [™] -T1 ^R Chemically Competent <i>E. coli</i>	20 reactions	C8620-03
One Shot™ TOP10 Electrocompetent <i>E. coli</i>	10 reactions 20 reactions	C4040-50 C4040-52
PureLink [™] PCR Purification Kit	50 preps	K3100-01
PureLink™ Quick Gel Extraction System	1 kit	K2100-12
PureLink™ HiPure Plasmid Miniprep Kit	25 preps 100 preps	K2100-02 K2100-03
E-Gel™ CloneWell™ 0.8% SYBR Safe™, E-Gel™ iBase™ & E-Gel™ Safe Imager™ Starter Kit	1 kit	G6500ST
E-Gel™ CloneWell™ 0.8% SYBR™ Safe gels, 18-Pak	18 gels	G6618-08
E-Gel™ 1.2% Starter Pak (6 gels + Powerbase™)	1 kit	G6000-01
E-Gel [™] 1.2% 18 Pak	18 gels	G5018-01
PCR 0ptimizer™ Kit	1 kit	K1220-01
AcTEV [™] Protease	1000 Units 10,000 Units	12575-015 12575-023
imMedia [™] Amp Liquid	20 pouches (200 mL medium)	Q600-20
imMedia™ Amp Agar	20 pouches (8–10 plates)	Q601-20
LB Broth (1X), liquid	500 mL	10855-021
S.O.C. Medium	10 × 10 mL	15544-034
Ampicillin Sodium Salt, irradiated	200 mg	11593-027
Gentamicin, liquid	10 mL (50 mg/mL)	15750-060

WARNING!

WARNING! GENERAL SAFETY. Using this product in a manner not specified in the user documentation may result in personal injury or damage to the instrument or device. Ensure that anyone using this product has received instructions in general safety practices for laboratories and the safety information provided in this document.

- Before using an instrument or device, read and understand the safety information provided in the user documentation provided by the manufacturer of the instrument or device.
- Before handling chemicals, read and understand all applicable Safety Data Sheets (SDSs) and use appropriate personal protective equipment (gloves, gowns, eye protection, etc). To obtain SDSs, see the "Documentation and Support" section in this document.

WARNING!

GENERAL CHEMICAL HANDLING. To minimize hazards, ensure laboratory personnel read and practice the general safety guidelines for chemical usage, storage, and waste provided below, and consult the relevant SDS for specific precautions and instructions:

- Read and understand the Safety Data Sheets (SDSs) provided by the chemical manufacturer before you store, handle, or work with any chemicals or hazardous materials. To obtain SDSs, see the "Documentation and Support" section in this document.
- Minimize contact with chemicals. Wear appropriate personal protective equipment when handling chemicals (for example, safety glasses, gloves, or protective clothing).
- Minimize the inhalation of chemicals. Do not leave chemical containers open. Use only with adequate ventilation (for example, fume hood).
- Check regularly for chemical leaks or spills. If a leak or spill occurs, follow the manufacturer's cleanup procedures as recommended in the SDS.
- Handle chemical wastes in a fume hood.
- Ensure use of primary and secondary waste containers. (A primary waste container holds the immediate waste. A secondary container contains spills or leaks from the primary container. Both containers must be compatible with the waste material and meet federal, state, and local requirements for container storage.)
- After emptying a waste container, seal it with the cap provided.
- Characterize (by analysis if necessary) the waste generated by the particular applications, reagents, and substrates used in your laboratory.
- Ensure that the waste is stored, transferred, transported, and disposed of according to all local, state/provincial, and/or national regulations.
- **IMPORTANT!** Radioactive or biohazardous materials may require special handling, and disposal limitations may apply.

Biological hazard safety

WARNING!

WARNING! BIOHAZARD. Biological samples such as tissues, body fluids, infectious agents, and blood of humans and other animals have the potential to transmit infectious diseases. Conduct all work in properly equipped facilities with the appropriate safety equipment (for example, physical containment devices). Safety equipment can also include items for personal protection, such as gloves, coats, gowns, shoe covers, boots, respirators, face shields, safety glasses, or goggles. Individuals should be trained according to applicable regulatory and company/ institution requirements before working with potentially biohazardous materials. Follow all applicable local, state/provincial, and/or national regulations. The following references provide general guidelines when handling biological samples in laboratory environment.

- U.S. Department of Health and Human Services, Biosafety in Microbiological and Biomedical Laboratories (BMBL), 5th Edition, HHS Publication No. (CDC) 21-1112, Revised December 2009; found at: www.cdc.gov/biosafety/publications/bmbl5/BMBL.pdf
- World Health Organization, Laboratory Biosafety Manual, 3rd Edition, WHO/CDS/CSR/LYO/2004.11; found at: www.who.int/csr/resources/publications/biosafety/Biosafety7.pdf

Customer and technical support

Visit **thermofisher.com/support** for the latest in services and support, including:

- Worldwide contact telephone numbers
- Product support, including:
 - -Product FAQs
 - Software, patches, and updates
 - Training for many applications and instruments
- Order and web support
- Product documentation, including:
 - User guides, manuals, and protocols
 - Certificates of Analysis
 - Safety Data Sheets (SDSs; also known as MSDSs)
 - **Note:** For SDSs for reagents and chemicals from other manufacturers, contact the manufacturer.

Limited product warranty

Life Technologies and/or its affiliate(s) warrant their products as set forth in the Life Technologies General Terms and Conditions of Sale found on the Life Technologies web site 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**.



