

# Bac-to-Bac<sup>®</sup> HBM TOPO<sup>®</sup> Secreted Expression System

An efficient site-specific transposition system to generate recombinant baculovirus for high-level secreted protein expression

Catalog no. A11339

**Revision date: 15 July 2009** Manual part no. A11341

MAN0001704

**User Manual** 

### Contents

Kit Contents and Storage	iv
Introduction	1
Description of the System	1
Bac-to-Bac® HBM TOPO® Secreted Expression System Components	3
Experiment Outline	5
Methods	7
Culturing Insect Cells	7
General Guidelines	7
Generating the Recombinant pFastBac <sup>™</sup> HBM Vector	9
Generating the Recombinant Bacmid	10
Transforming DH10Bac <sup>™</sup> E. coli	
Isolating Recombinant Bacmid DNA	
Analyzing Recombinant Bacmid DNA by PCR	15
Producing Recombinant Baculovirus	18
Transfecting Insect Cells	
Isolating P1 Viral Stock	21
Amplifying Your Baculoviral Stock	23
Performing a Viral Plaque Assay	25
Expressing Your Recombinant Protein	
Analyzing Recombinant Protein	
Expected Results	40
Optimizing Protein Expression	41
Large-Scale Expression of Recombinant Protein	
Troubleshooting	44
Appendix	49
Recipes	
Map of pFastBac <sup>™</sup> /HBM-TOPO <sup>®</sup>	
Map of pFastBac <sup>™</sup> Gus Control Plasmid	
Accessory Products	
Technical Support	56
Purchaser Notification	
References	

#### Kit Contents and Storage

#### System Components

Each Bac-to-Bac<sup>®</sup> HBM TOPO<sup>®</sup> Secreted Expression System contains the components listed below. See the next page for a detailed description of other reagents supplied with each system.

Component	Amount
Bac-to-Bac® HBM TOPO® Cloning Kit	1 kit
One Shot <sup>®</sup> Mach1-T1 <sup>R</sup> Chemically Competent <i>E. coli</i>	1 kit
MAX Efficiency <sup>®</sup> DH10Bac <sup>™</sup> Competent E. coli	4 kits
Cellfectin <sup>®</sup> II Reagent	1 each
Bac-to-Bac® TOPO® Cloning Kit manual	1 each
Bac-to-Bac <sup>®</sup> HBM TOPO <sup>®</sup> Secreted Expression System manual	1 each

#### Shipping/Storage

The Bac-to-Bac<sup>®</sup> HBM TOPO<sup>®</sup> Secreted Expression System is shipped in four boxes as described below. Upon receipt, store each box as detailed below. All reagents are guaranteed for six months if stored properly.

Box	Item	Shipping	Storage
1	Bac-to-Bac <sup>®</sup> HBM TOPO <sup>®</sup> Cloning Kit	Dry ice	-20°C
2	One Shot <sup>®</sup> Mach1-T1 <sup>R</sup> Chemically Competent <i>E. coli</i>	Dry ice	–80°C
3	MAX Efficiency <sup>®</sup> DH10Bac <sup>™</sup> Competent E. coli	Dry ice	-80°C
4	Cellfectin <sup>®</sup> II Reagent	Gel ice	4°C

### Kit Contents and Storage, Continued

Bac-to-Bac<sup>®</sup> HBM TOPO<sup>®</sup> Cloning Kit The cloning reagents for the Bac-to-Bac<sup>®</sup> HBM TOPO<sup>®</sup> Cloning Kit are listed below. **Store the contents at –20°C.** 

Item	Concentration	Amount
pFastBac <sup>™</sup> /HBM-TOPO <sup>®</sup>	20 mL at 10 ng/µl in	20 µl
vector	50% glycerol	
	50 mM Tris-HCl, pH 7.4 (at 25°C)	
	1 mM EDTA	
	2 mM DTT	
	0.1% Triton <sup>®</sup> X-100	
	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 <sub>2</sub>	
	0.01% gelatin	
dNTP Mix	12.5 mM each dATP, dCTP, dGTP, and dTTP; neutralized at pH 8.0 in water	10 µL
Salt Solution	1.2 M NaCl	50 µL
	60 mM MgCl <sub>2</sub>	
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 8.0	10 µL
Polyhedrin forward sequencing primer	100 ng/μL in TE buffer, pH 8.0	20 µL
SV40 polyA reverse sequencing primer	100 ng/μL in TE buffer, pH 8.0	20 µL
pFastBac <sup>™</sup> 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

#### One Shot<sup>®</sup> Mach1<sup>™</sup> T1<sup>R</sup> Competent E. coli

The following reagents are included in the One Shot<sup>®</sup> Mach1<sup>™</sup> T1<sup>R</sup> Chemically Competent *E. coli* kit. Transformation efficiency of One Shot<sup>®</sup> Mach1<sup>™</sup> T1<sup>R</sup> *E. coli* cells is  $\ge 1 \times 10^9$  cfu/µg DNA. Store cells at -80°C.

Reagent	Composition	Amount
One Shot <sup>®</sup> Mach1 <sup>™</sup> T1 <sup>R</sup> Chemically Competent <i>E. coli</i>	-	$21\times 50~\mu L$
S.O.C. Medium	2% tryptone	6 mL
(may be stored at room	0.5% yeast extract	
temperature or 4°C)	10 mM NaCl	
	2.5 mM KCl	
	10 mM MgCl <sub>2</sub>	
	10 mM MgSO <sub>4</sub>	
	20 mM glucose	
pUC19 Control DNA	10 pg/μL in 5 mM Tris-HCl, 0.5 mM EDTA, pH 8.0	50 µL

## Genotype of

 $F^- \phi 80(lacZ) \Delta M15 \Delta lacX74 hsdR(r_K^-m_K^+) \Delta recA1398 endA1 tonA$ 

Mach1<sup>™</sup> T1<sup>R</sup>

#### MAX Efficiency<sup>®</sup> DH10Bac<sup>¬</sup> Competent E. coli

MAX Efficiency<sup>®</sup> DH10Bac<sup>™</sup> Competent E. coli (Box 3) have a transformation efficiency of  $1 \times 10^8$  cfu/µg DNA. Store at -80°C.

Item	Composition	Amount
MAX Efficiency <sup>®</sup> DH10Bac <sup>™</sup> Competent <i>E. coli</i>	-	4  kits (4 × 5 reactions)
pUC19 Control DNA	$10\ pg/\mu L$ in 5 mM Tris-HCl, $0.5\ mM$ EDTA, pH $8.0$	100 µL

 $F^-$  mcrA Δ(mrr-hsdRMS-mcrBC)  $\phi$ 80lacZΔM15 ΔlacX74 recA1 endA1 araD139 Δ(ara, Genotype of DH10Bac<sup>®</sup> *leu*)7697 galU galK λ<sup>-</sup> rpsL nupG/bMON14272/pMON7124

Cellfectin<sup>®</sup> II Cellfectin® II Reagent is a proprietary cationic lipid formulation that offers the highest transfection efficiencies and protein expression levels on the widest variety Transfection of adherent and suspension insect cell lines. Reagent

Amount supplied: 1 mL

**Composition:** 1 mg/mL transfection reagent in membrane-filtered water Storage conditions: 4°C (do not freeze)

### Introduction

## **Description of the System**

System Overview	• The Bac-to-Bac <sup>®</sup> HBM TOPO <sup>®</sup> Secreted Expression System provides a rapid, effective method of generating recombinant baculoviruses for secreted expression of your protein of interest. The Bac-to-Bac <sup>®</sup> HBM TOPO <sup>®</sup> Secreted Expression System combines the ease of <b>blunt-end</b> TOPO <sup>®</sup> cloning with the efficiency of site-specific transposition technology of the Bac-to-Bac <sup>®</sup> System.
Advantages of the Bac-to-Bac <sup>®</sup> HBM TOPO <sup>®</sup> Secreted	Using the Bac-to-Bac <sup>®</sup> HBM TOPO <sup>®</sup> Secreted Expression System to generate a recombinant baculovirus provides the following advantages over the traditional method using homologous recombination:
Expression System	• Enables the cloning of the gene of interest as a <b>blunt-end PCR product</b> in a highly efficient one-step reaction, thus allowing the use of proofreading polymerases in the PCR amplification step
	• Requires less than 2 weeks to identify and purify a recombinant baculovirus, compared to the 4–6 weeks required to generate a recombinant baculovirus using homologous recombination
	• Reduces the need for multiple rounds of plaque purification, because the recombinant virus DNA isolated from selected colonies is not mixed with parental, non-recombinant virus
	<ul> <li>Permits rapid and simultaneous isolation of multiple recombinant baculoviruses</li> <li>Allows secreted expression of protein variants for structure/function studies</li> </ul>
Purpose of This Manual	This manual provides an overview of the Bac-to-Bac <sup>®</sup> HBM TOPO <sup>®</sup> Secreted Expression System, and provides instructions and guidelines to:
	1. Transform the pFastBac <sup>™</sup> construct containing your gene of interest into MAX Efficiency <sup>®</sup> DH10Bac <sup>™</sup> competent <i>E. coli</i> to generate recombinant bacmid.
	2. Transfect the recombinant bacmid DNA into the insect cell line of choice to produce recombinant baculovirus particles.
	3. Amplify and titer the baculoviral stock, and use this stock to infect insect cells to express your recombinant protein.
	Detailed instructions for cloning your gene of interest into the pFastBac <sup>™</sup> /HBM-TOPO <sup>®</sup> vector are provided in the Bac-to-Bac <sup>®</sup> TOPO <sup>®</sup> Cloning Kit manual (part no. A10605) supplied with the Bac-to-Bac <sup>®</sup> HBM TOPO <sup>®</sup> Secreted Expression System.
	The Bac-to-Bac <sup>®</sup> TOPO <sup>®</sup> Cloning Kit manual is also available for downloading at www.invitrogen.com or from Technical Support (see page 56).

#### Description of the System, Continued



The Bac-to-Bac<sup>®</sup> HBM TOPO<sup>®</sup> Secreted Expression System is designed to help you create a recombinant baculovirus for high-level secreted expression of your gene of interest in insect cells. Although the system has been designed to help you to easily produce recombinant baculovirus and express your protein of interest, use of the system is geared towards those users who are familiar with baculovirus biology and insect cell culture. We highly recommend that users possess a working knowledge of viral and tissue culture techniques.

For more information about baculovirus biology, refer to published reference sources (King & Possee, 1992; Luckow, 1991; O'Reilly *et al.*, 1992). For more information about insect cell culture, refer to the *Guide to Baculovirus Expression Vector Systems (BEVS) and Insect Cell Culture Techniques* available from Invitrogen at www.invitrogen.com or by contacting Technical Support (see page 56).

# Bac-to-Bac<sup>®</sup> HBM TOPO<sup>®</sup> Secreted Expression System Components

Components of the Bac-to-Bac<sup>®</sup> HBM TOPO<sup>®</sup> Secreted Expression System The Bac-to-Bac<sup>®</sup> HBM TOPO<sup>®</sup> Secreted Expression System facilitates rapid and efficient generation of recombinant baculoviruses (Ciccarone *et al.*, 1997) by combining the ease of TOPO<sup>®</sup> cloning with the efficiency of the Bac-to-Bac<sup>®</sup> System. Based on a method developed by Luckow *et al.* (Luckow *et al.*, 1993), the Bac-to-Bac<sup>®</sup> HBM TOPO<sup>®</sup> Secreted Expression System takes advantage of the site-specific transposition properties of the Tn7 transposon to simplify and enhance the process of generating recombinant bacmid DNA. The following major system components are described in detail below:

- pFastBac<sup>™</sup>/HBM-TOPO<sup>®</sup> plasmid—allows the rapid generation of an expression construct containing the gene of interest under the control of a baculovirus-specific strong polyhedrin (P<sub>H</sub>) promoter and in frame with the Honey Bee Mellitin (HBM) secretion signal coding sequence. One Shot<sup>®</sup> Mach1<sup>™</sup> T1R Chemically Competent *E. coli* enable same-day isolation of recombinant pFastBac<sup>™</sup> expression construct from the transformation mix.
- An *E. coli* host strain, DH10Bac<sup>™</sup>— contains a baculovirus shuttle vector (bacmid) and a helper plasmid to facilitate the generation of a recombinant bacmid following transposition of the pFastBac<sup>™</sup> expression construct.
- **Cellfectin® II Reagent**—for fast, efficient transfection of insect cells to generate recombinant baculovirus particles.

**Note:** A control expression plasmid, pFastBac<sup>TM</sup> Gus allows production of a recombinant baculovirus which, when used to infect insect cells, constitutively expresses  $\beta$ -glucuronidase.

#### pFastBac<sup>™</sup>/HBM-TOPO<sup>®</sup> vector

The first major component of the System is the pFastBac<sup>™</sup>/HBM-TOPO<sup>®</sup> vector into which your gene of interest is cloned.

After you amplify your gene of interest using a proofreading polymerase and clone it into the pFastBac<sup>™</sup>/HBM-TOPO<sup>®</sup> vector as a blunt-end PCR product, you transform One Shot<sup>®</sup> Mach1<sup>™</sup> T1<sup>R</sup> Chemically Competent *E. coli*. You then select and analyze transformants for the correct insertion of your blunt-end PCR products, and use the recombinant vector as a "donor plasmid" to generate a recombinant baculovirus.

The expression of the gene of interest is controlled by the *Autographa californica* multiple nuclear polyhedrosis virus (AcMNPV) polyhedrin (P<sub>H</sub>) promoter for high-level expression in insect cells. This expression cassette is flanked by the left and right arms of Tn7. The cassette also contains a gentamicin resistance gene and an SV40 polyadenylation signal to form a mini Tn7.

The presence of the N-terminal Honey Bee Mellitin (HBM) secretion signal coding sequence on the plasmid facilitates the secretion of the cloned gene product into the extracellular medium; the C-terminal polyhistidine tag allows easy purification of the secreted protein.

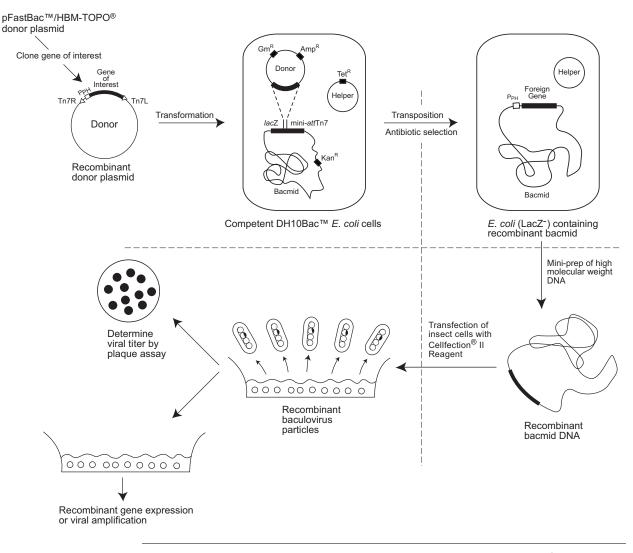
# Bac-to-Bac<sup>®</sup> HBM TOPO<sup>®</sup> Secreted Expression System Components, Continued

DH10Bac <sup>™</sup> <i>E. coli</i>	The second major component of the System is the DH10Bac <sup>TM</sup> <i>E. coli</i> strain that is used as the host for the pFastBac <sup>TM</sup> /HBM construct containing your gene of interest. DH10Bac <sup>TM</sup> cells contain a baculovirus shuttle vector (bacmid) with a mini- <i>att</i> Tn7 target site and a helper plasmid (see the next page for details).
	After the pFastBac <sup>™</sup> /HBM expression plasmid (the "donor plasmid") is transformed into DH10Bac <sup>™</sup> cells, transposition occurs between the mini-Tn7 element on the pFastBac <sup>™</sup> /HBM vector and the mini- <i>att</i> Tn7 target site on the bacmid to generate a recombinant bacmid. This transposition reaction occurs in the presence of transposition proteins supplied by the helper plasmid.
	After you have performed the transposition reaction, you isolate the high molecular weight recombinant bacmid DNA and transfect the bacmid DNA into insect cells using the Cellfectin <sup>®</sup> II reagent to generate a recombinant baculovirus that can be used for preliminary expression experiments. After the baculoviral stock is amplified and titered, this high-titer stock can be used to infect insect cells for large-scale expression of the recombinant protein of interest.
	For a schematic representation of the Bac-to-Bac <sup>®</sup> HBM TOPO <sup>®</sup> Secreted Expression System, see the diagram on page 5.
Baculovirus Shuttle Vector	The baculovirus shuttle vector (bacmid), bMON14272 (136 kb), present in DH10Bac <sup>™</sup> <i>E. coli</i> contains:
	• A low-copy number mini-F replicon.
	Kanamycin resistance marker.
	<ul> <li>A segment of DNA encoding the LacZα peptide from a pUC-based cloning vector into which the attachment site for the bacterial transposon, Tn7 (mini-<i>att</i>Tn7) has been inserted. Insertion of the mini-<i>att</i>Tn7 attachment site does not disrupt the reading frame of the LacZα peptide.</li> </ul>
	The bacmid propagates in <i>E. coli</i> DH10Bac <sup><math>\mathbb{M}</math></sup> as a large plasmid that confers resistance to kanamycin. This bacmid can complement a <i>lacZ</i> deletion present on the chromosome to form colonies that are blue (Lac <sup>+</sup> ) in the presence of a chromogenic substrate such as Bluo-gal or X-gal and the inducer, IPTG.
	Recombinant bacmids (composite bacmids) are generated by transposing a mini-Tn7 element from a pFastBac <sup>™</sup> donor plasmid to the mini- <i>att</i> Tn7 attachment site on the bacmid. The Tn7 transposition functions are provided by a helper plasmid (see below).
Helper Plasmid	DH10Bac <sup><math>TM</math></sup> <i>E. coli</i> also contain the helper plasmid, pMON7124 (13.2 kb), which encodes the transposase and confers resistance to tetracycline. The helper plasmid provides the Tn7 transposition function <i>in trans</i> (Barry, 1988).
Cellfectin <sup>®</sup> II Reagent	Cellfectin <sup>®</sup> II Reagent is a proprietary cationic lipid formulation that offers the highest transfection efficiencies and protein expression levels on the widest variety of adherent and suspension insect cell lines, including Sf9 and Sf21 cells.

#### **Experiment Outline**

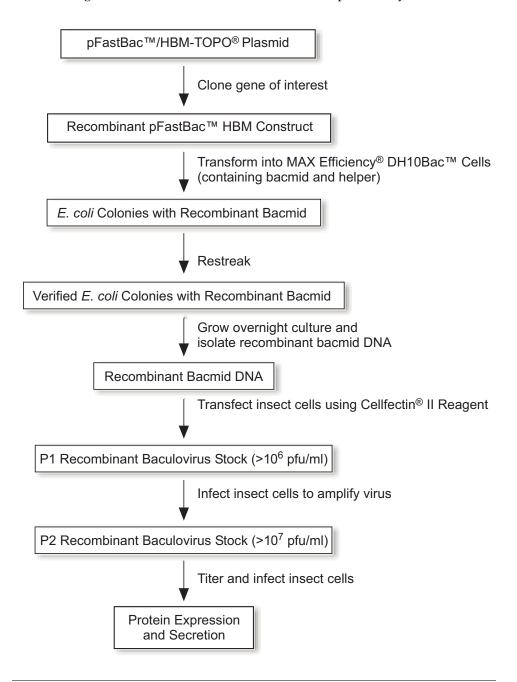
#### Diagram of the Bac-to-Bac<sup>®</sup> System

The figure below depicts the generation of recombinant baculovirus and the expression of your gene of interest using the Bac-to-Bac<sup>®</sup> HBM TOPO<sup>®</sup> Secreted Expression System.



#### Experiment Outline, Continued

**Flow Chart** The figure below illustrates the general steps required to express your gene of interest using the Bac-to-Bac<sup>®</sup> HBM TOPO<sup>®</sup> Secreted Expression System.



# Methods Culturing Insect Cells

### **General Guidelines**

Introduction	We recommend using <i>Spodoptera frugiperda</i> Sf9 or Sf21 insect cells as the host for your baculovirus transfer vector when you produce your recombinant bacmid.
	We recommend using High Five <sup>™</sup> cells for secreted expression of your protein from the recombinant bacmid, because they are particularly well suited for expression of secreted recombinant proteins.
	Before you start your transfection and expression experiments, be sure to have cultures of Sf9 or Sf21 and High Five <sup>™</sup> cells growing, and have frozen master stocks available. Sf9, Sf21, and High Five <sup>™</sup> cells and cell culture reagents are available separately from Invitrogen (see page 54 for ordering information).
	Note: High Five <sup><math>TM</math></sup> and Mimic <sup><math>TM</math></sup> Sf9 insect cells are suitable for use for expression only.
Using Serum-Free Medium	Insect cells may be cultured under serum-free conditions. We recommend using Sf-900 II SFM or Sf-900 <sup>™</sup> III SFM available from Invitrogen (see page 54) for culturing Sf9 and Sf21 cells. Sf-900 II SFM and Sf-900 <sup>™</sup> III SFM are protein-free media optimized for the growth and maintenance of Sf9 and Sf21 cells. For culturing High Five <sup>™</sup> cells under serum-free conditions, use Express Five <sup>®</sup> SFM (see page 54 for ordering). Express Five <sup>®</sup> SFM is optimized for the growth and maintenance of High Five <sup>™</sup> cells, as well as for the large-scale production and secretion of recombinant proteins expressed using the Bac-to-Bac <sup>®</sup> HBM TOPO <sup>®</sup> Secreted Expression System. For more information, refer to www.invitrogen.com or contact Technical Support (see page 56).
Insect Cell Culture Reference Guide	For guidelines and detailed information on insect cell culture, refer to the <i>Guide to Baculovirus Expression Vector Systems (BEVS) and Insect Cell Culture Techniques,</i> available for downloading at www.invitrogen.com or by contacting Technical Support (see page 56). This guide contains information on:
	Maintaining and passaging insect cells in adherent and suspension culture
	Freezing cells
	<ul> <li>Using serum-free medium (includes protocols to adapt cells to serum-free medium)</li> </ul>
	Scaling up cell culture
	Continued on next nage

## General Guidelines, Continued

General Guidelines	Insect cells are very sensitive to environmental factors. In addition to chemical and nutritional culture factors, physical factors can also affect insect cell growth; therefore optimization is required to maximize cell growth. Consider the following when culturing insect cells:
	• <b>Temperature:</b> The optimal range to grow and infect cultured insect cells is 27°C to 28°C.
	• <b>pH:</b> A range of 6.1 to 6.4 works well for most culture systems. Sf-900 II SFM will maintain a pH in this range under conditions of normal air and open-capped culture systems.
	• <b>Osmolality:</b> The optimal osmolality of medium for use with lepidopteran cell lines is 345 to 380 mOsm/kg.
	• Aeration: Insect cells require passive oxygen diffusion for optimal growth and recombinant protein expression. Active or controlled oxygenated systems require dissolved oxygen at 10% to 50% of air saturation.
	• Shear Fforces: Suspension culture generates mechanical shear forces. Growing insect cells in serum-containing media (10% to 20% FBS) generally provides adequate protection from cellular shear forces. If you are growing insect cells in serum-free conditions, supplementation with a shear force protectant such as PLURONIC <sup>®</sup> F-68 may be required.
	Note: Growing cells in Sf-900 II SFM or Sf-900 <sup>™</sup> III SFM does <b>not</b> require addition of shear force protectants.
Cells for Transfection	You need log-phase Sf9 or Sf21 cells with >95% viability to perform a successful transfection. Refer to page 19 to determine how many cells you will need for transfection.

# Generating the Recombinant pFastBac<sup>™</sup> HBM Vector

Introduction	To generate a recombinant plasmid containing your gene of interest for use in the Bac-to-Bac <sup>®</sup> HBM TOPO <sup>®</sup> Secreted Expression System, perform the following steps:
	<ol> <li>Generate a <b>blunt-end</b> PCR product containing your gene of interest with a thermostable <b>proofreading DNA polymerase</b> such as the Platinum<sup>®</sup> <i>Pfx</i> or the AccuPrime<sup>™</sup> <i>Pfx</i> DNA Polymerase.</li> </ol>
	<ol> <li>TOPO<sup>®</sup> Clone your blunt-end PCR product into the pFastBac<sup>™</sup>/HBM-TOPO<sup>®</sup> vector, and use the reaction to transform One Shot<sup>®</sup> Mach1<sup>™</sup> T1<sup>R</sup> Chemically Competent <i>E. coli</i>. Do not transform the ligation reaction into DH10Bac<sup>™</sup> cells.</li> </ol>
	<ol> <li>Pick colonies, isolate plasmid DNA (see below), and screen for insert directionality by sequencing expression clones with the primers provided in the kit.</li> </ol>
	For detailed instructions, refer to the Bac-to-Bac <sup>®</sup> TOPO <sup>®</sup> Cloning Kit manual (part no. A10605) supplied with this kit. The Bac-to-Bac <sup>®</sup> TOPO <sup>®</sup> Cloning Kit manual is also available at www.invitrogen.com or from Technical Support (see page 56).
Guidelines for Isolating DNA	1. Pick 10 overnight-grown colonies from the selective plates and culture them overnight in LB medium containing 100 $\mu$ g/mL ampicillin.
	<ol> <li>Isolate plasmid DNA using your method of choice. If you need ultra-pure plasmid DNA for automated or manual sequencing, we recommend using Invitrogen's PureLink<sup>™</sup> HiPure Mini Plasmid Purification Kit (see page 53 for ordering information).</li> </ol>
	You need 1 ng of purified recombinant plasmid (5 $\mu$ L at 200 pg/ $\mu$ L) to transform into DH10Bac <sup>TM</sup> <i>E. coli</i> for transposition into the bacmid (see page 11).
	<b>Note</b> : If you have used One Shot <sup>®</sup> Mach1 <sup>™</sup> T1 <sup>R</sup> 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.
Important	When generating the recombinant plasmid containing your gene of interest for use in the Bac-to-Bac <sup>®</sup> HBM TOPO <sup>®</sup> Secreted Expression System, you <b>must</b> follow certain design parameters for your PCR insert and the recommendations for the transformation procedure outlined in the Bac-to-Bac <sup>®</sup> TOPO <sup>®</sup> Cloning Kit manual.
	To ensure proper expression of your recombinant protein, it is imperative that you read the sections on generating the blunt-end PCR product, blunt-end TOPO <sup>®</sup> Cloning, transforming One Shot <sup>®</sup> Mach1 <sup>™</sup> T1 <sup>R</sup> Chemically Competent <i>E. coli</i> , and analyzing transformants in the Bac-to-Bac <sup>®</sup> TOPO <sup>®</sup> Cloning Kit manual before beginning.
pFastBac <sup>™</sup> /HBM- TOPO <sup>®</sup> vector	For a vector map of pFastBac <sup>™</sup> /HBM-TOPO <sup>®</sup> , see page 51. For more instructions on generating your recombinant plasmid containing your gene of interest, refer to the Bac-to-Bac <sup>®</sup> TOPO <sup>®</sup> Cloning Kit manual (part no. A10605) supplied with this kit, also available at www.invitrogen.com or by contacting Technical Support (see page 56).

#### **Generating the Recombinant Bacmid**

# Transforming DH10Bac<sup>™</sup> E. coli

transforming DH10Bac <sup>™</sup> cells are provided in this section.         Positive Control       The pFastBac <sup>™</sup> /HBM-TOPO <sup>®</sup> vector is supplied with the control plasmid pFastBac <sup>™</sup> Gus for use as a positive transfection and expression control. We recommend including the control plasmid in your DH10Bac <sup>™</sup> transformation experiments. For a map of the control plasmid, see page 52.         Materials Needed       Your purified pFastBac <sup>™</sup> /HBM construct (200 pg/mL in TE, pH 8.0; see page 9)         Positive expression control (i.e., pFastBac <sup>™</sup> Gus; use as a control for transposition)         MAX Efficiency <sup>®</sup> DH10Bac <sup>™</sup> chemically competent cells (supplied with the Bac-to-Bac <sup>®</sup> HBM TOPO <sup>®</sup> Secreted Expression System; use 1 tube of competer cells for every transformation)         pUC19 (supplied with the MAX Efficiency <sup>®</sup> DH10Bac <sup>™</sup> E. coli; use as a control for transformation, if desired)         LB agar plates containing kanamycin, gentamicin, tetracycline, Bluo-gal, and IPTG (3 freshly prepared plates for each transformation; see below)         LB agar plate containing 100 µg/mL ampicillin (for plating pUC19 transformation control)         S.O.C. Medium         42°C water bath         37°C shaking and non-shaking incubator	Introduction	After you have generated your pFastBac <sup><math>TM</math></sup> HBM construct containing your gene of interest, transform purified plasmid DNA into DH10Bac <sup><math>TM</math></sup> <i>E. coli</i> for transposition into the bacmid. Use blue/white selection to identify colonies containing the recombinant bacmid.
<ul> <li>pFastBac<sup>™</sup> Gus for use as a positive transfection and expression control. We recommend including the control plasmid in your DH10Bac<sup>™</sup> transformation experiments. For a map of the control plasmid, see page 52.</li> <li>Materials Needed <ul> <li>Your purified pFastBac<sup>™</sup>/HBM construct (200 pg/mL in TE, pH 8.0; see page 9)</li> <li>Positive expression control (i.e., pFastBac<sup>™</sup> Gus; use as a control for transposition)</li> <li>MAX Efficiency<sup>®</sup> DH10Bac<sup>™</sup> chemically competent cells (supplied with the Bac-to-Bac<sup>®</sup> HBM TOPO<sup>®</sup> Secreted Expression System; use 1 tube of competer cells for every transformation)</li> <li>pUC19 (supplied with the MAX Efficiency<sup>®</sup> DH10Bac<sup>™</sup> E. coli; use as a control for transformation, if desired)</li> <li>LB agar plates containing kanamycin, gentamicin, tetracycline, Bluo-gal, and IPTG (3 freshly prepared plates for each transformation; see below)</li> <li>LB agar plate containing 100 µg/mL ampicillin (for plating pUC19 transformation control)</li> <li>S.O.C. Medium</li> <li>42°C water bath</li> <li>37°C shaking and non-shaking incubator</li> </ul> </li> </ul>		Bac-to-Bac <sup>®</sup> HBM TOPO <sup>®</sup> Secreted Expression System. These cells are also available separately from Invitrogen (see page 53). Guidelines and instructions for
<ul> <li>page 9)</li> <li>Positive expression control (i.e., pFastBac<sup>™</sup> Gus; use as a control for transposition)</li> <li>MAX Efficiency<sup>®</sup> DH10Bac<sup>™</sup> chemically competent cells (supplied with the Bac-to-Bac<sup>®</sup> HBM TOPO<sup>®</sup> Secreted Expression System; use 1 tube of competer cells for every transformation)</li> <li>pUC19 (supplied with the MAX Efficiency<sup>®</sup> DH10Bac<sup>™</sup> <i>E. coli</i>; use as a control for transformation, if desired)</li> <li>LB agar plates containing kanamycin, gentamicin, tetracycline, Bluo-gal, and IPTG (3 freshly prepared plates for each transformation; see below)</li> <li>LB agar plate containing 100 µg/mL ampicillin (for plating pUC19 transformation control)</li> <li>S.O.C. Medium</li> <li>42°C water bath</li> <li>37°C shaking and non-shaking incubator</li> </ul>	Positive Control	pFastBac <sup>™</sup> Gus for use as a positive transfection and expression control. We recommend including the control plasmid in your DH10Bac <sup>™</sup> transformation
<ul> <li>transposition)</li> <li>MAX Efficiency<sup>®</sup> DH10Bac<sup>™</sup> chemically competent cells (supplied with the Bac-to-Bac<sup>®</sup> HBM TOPO<sup>®</sup> Secreted Expression System; use 1 tube of competer cells for every transformation)</li> <li>pUC19 (supplied with the MAX Efficiency<sup>®</sup> DH10Bac<sup>™</sup> E. coli; use as a control for transformation, if desired)</li> <li>LB agar plates containing kanamycin, gentamicin, tetracycline, Bluo-gal, and IPTG (3 freshly prepared plates for each transformation; see below)</li> <li>LB agar plate containing 100 µg/mL ampicillin (for plating pUC19 transformation control)</li> <li>S.O.C. Medium</li> <li>42°C water bath</li> <li>37°C shaking and non-shaking incubator</li> </ul>	Materials Needed	
<ul> <li>Bac-to-Bac<sup>®</sup> HBM TOPO<sup>®</sup> Secreted Expression System; use 1 tube of competer cells for every transformation)</li> <li>pUC19 (supplied with the MAX Efficiency<sup>®</sup> DH10Bac<sup>™</sup> E. coli; use as a control for transformation, if desired)</li> <li>LB agar plates containing kanamycin, gentamicin, tetracycline, Bluo-gal, and IPTG (3 freshly prepared plates for each transformation; see below)</li> <li>LB agar plate containing 100 µg/mL ampicillin (for plating pUC19 transformation control)</li> <li>S.O.C. Medium</li> <li>42°C water bath</li> <li>37°C shaking and non-shaking incubator</li> </ul>		
<ul> <li>for transformation, if desired)</li> <li>LB agar plates containing kanamycin, gentamicin, tetracycline, Bluo-gal, and IPTG (3 freshly prepared plates for each transformation; see below)</li> <li>LB agar plate containing 100 μg/mL ampicillin (for plating pUC19 transformation control)</li> <li>S.O.C. Medium</li> <li>42°C water bath</li> <li>37°C shaking and non-shaking incubator</li> </ul>		Bac-to-Bac® HBM TOPO® Secreted Expression System; use 1 tube of competent
<ul> <li>IPTG (3 freshly prepared plates for each transformation; see below)</li> <li>LB agar plate containing 100 μg/mL ampicillin (for plating pUC19 transformation control)</li> <li>S.O.C. Medium</li> <li>42°C water bath</li> <li>37°C shaking and non-shaking incubator</li> </ul>		I (- II , , ,
<ul> <li>transformation control)</li> <li>S.O.C. Medium</li> <li>42°C water bath</li> <li>37°C shaking and non-shaking incubator</li> </ul> Prepare LB agar plates containing 50 μg/mL kanamycin, 7 μg/mL gentamicin,		
<ul> <li>42°C water bath</li> <li>37°C shaking and non-shaking incubator</li> <li>Prepare LB agar plates containing 50 μg/mL kanamycin, 7 μg/mL gentamicin,</li> </ul>		
<ul> <li>37°C shaking and non-shaking incubator</li> <li>Prepare LB agar plates containing 50 μg/mL kanamycin, 7 μg/mL gentamicin,</li> </ul>		• S.O.C. Medium
Prepare LB agar plates containing 50 $\mu$ g/mL kanamycin, 7 $\mu$ g/mL gentamicin,		• 42°C water bath
Prepare LB agar plates containing 50 $\mu$ g/mL kanamycin, 7 $\mu$ g/mL gentamicin,		• 37°C shaking and non-shaking incubator
$-\frac{10 \mu\text{g}}{\text{mL}}$ tetracycline, 100 $\mu\text{g}$ /mL Bluo-gal, and 40 $\mu\text{g}$ /mL IPIG to select for transformants. See page 50 for instructions to prepare plates.	MMENO	10 $\mu$ g/mL tetracycline, 100 $\mu$ g/mL Bluo-gal, and 40 $\mu$ g/mL IPTG to select for

If you are preparing LB plates using a pre-mixed formulation, we use Luria Broth Base (see page 53) instead of Lennox L (LB). Using Lennox L plates reduces the color intensity and may lower the number of colonies obtained.

**Note:** Use Bluo-gal instead of X-gal for blue/white selection. Bluo-gal generally produces a darker blue color than X-gal.

# Transforming DH10Bac<sup>™</sup> *E. coli*, Continued

Preparing for Transformation	For each transformation, you will need one vial of competent cells and three selective plates.				
	• Equilibrate a water bath to 42°C.				
	• Warm selective plates at 37°C for 30 minutes.				
	• Warm the S.O.C. Medium to room temperature.				
Transformation Procedure	Follow the procedure below to transform MAX Efficiency <sup>®</sup> DH10Bac <sup>™</sup> chemically competent <i>E. coli</i> cells with your recombinant pFastBac <sup>™</sup> HBM construct. We recommend including positive controls for transposition (i.e., pFastBac <sup>™</sup> Gus) and transformation (i.e., pUC19) in your experiment to help you evaluate your results.				
	1. Thaw <b>on ice</b> one vial of MAX Efficiency <sup>®</sup> DH10Bac <sup>™</sup> competent <i>E. coli</i> cells for each transformation.				
	<ol> <li>For each transformation, add the appropriate amount of plasmid DNA to 100 µL of DH10Bac<sup>™</sup> cells and mix gently. Do not pipet up and down to mix.</li> </ol>				
	<ul> <li>Your recombinant pFastBac<sup>™</sup>/HBM construct: 1 ng (5 μL)</li> </ul>				
	<ul> <li>pFastBac<sup>™</sup> Gus control plasmid: 1 ng</li> </ul>				
	• pUC19 control: 50 pg (5 μL)				
	3. Incubate the cells on ice for 30 minutes.				
	4. Heat-shock the cells for 45 seconds at 42°C without shaking.				
	5. Immediately transfer the tubes to ice and chill for 2 minutes.				
	6. Add 900 μL of room temperature S.O.C. Medium.				
	7. <b>For pFastBac<sup>™</sup> transformations:</b> Shake the tubes at 37°C at 225 rpm for 4 hours.				
	For pUC19 transformation: Shake the tube at 37°C at 225 rpm for 1 hour.				
	<ol> <li>For each pFastBac<sup>™</sup> transformation: Prepare 10-fold serial dilutions of the cells (10<sup>-1</sup>, 10<sup>-2</sup>, 10<sup>-3</sup>) with S.O.C. Medium. Plate 100 µL of each dilution on an LB agar plate containing 50 µg/mL kanamycin, 7 µg/mL gentamicin, 10 µg/mL tetracycline, 100 µg/mL Bluo-gal, and 40 µg/mL IPTG.</li> </ol>				
	<b>For the pUC19 transformation:</b> Dilute the cells 1:100 with S. O.C. Medium. Plate 100 $\mu$ L of the dilution on an LB agar plate containing 100 $\mu$ g/mL ampicillin.				
	9. Incubate the plates for 48 hours at 37°C. Pick white colonies for analysis (see the next page for recommendations).				
	<b>Note:</b> We do <b>not</b> recommend picking colonies earlier than 48 hours. Incubating the plates for less than 48 hours may create difficulty distinguishing between white and blue colonies.				

## Transforming DH10Bac<sup>™</sup> E. coli, Continued

# 

Insertions of the mini-Tn7 into the mini-*att*Tn7 attachment site on the bacmid disrupt the expression of the LacZ $\alpha$  peptide, so colonies containing the recombinant bacmid are white in a background of blue colonies that harbor the unaltered bacmid. **Select white colonies for analysis.** True white colonies tend to be large. To avoid selecting false positives, choose the largest, most isolated white colonies. Avoid picking colonies that appear gray or are darker in the center, because they can contain a mixture of cells with empty bacmid and recombinant bacmid.

#### Verifying the Phenotype

- Pick 10 white colonies and restreak them on fresh LB agar plates containing 50 μg/mL kanamycin, 7 μg/mL gentamicin, 10 μg/mL tetracycline, 100 μg/mL Bluo-gal, and 40 μg/mL IPTG. Incubate the plates overnight at 37°C.
  - 2. From a single colony confirmed to have a white phenotype on restreaked plates containing Bluo-gal and IPTG, inoculate a liquid culture containing 50 μg/mL kanamycin, 7 μg/mL gentamicin, and 10 μg/mL tetracycline.
  - 3. Isolate recombinant bacmid DNA for analysis using the procedure provided on the next page.
  - 4. Analyze the recombinant bacmid DNA to verify successful transposition to the bacmid. We recommend using PCR to analyze your bacmid DNA (see **Analyzing Recombinant Bacmid DNA by PCR**, page 15).

**Note:** It is possible to verify successful transposition to the bacmid by using agarose gel electrophoresis to look for the presence of high molecular weight DNA. This method is less reliable than performing PCR analysis, because high molecular weight DNA can be difficult to visualize.



You may also use other methods to prepare purified recombinant bacmid DNA for analysis and transfection. However, bacmid DNA must be clean and free from phenol and sodium chloride because contaminants may kill the insect cells, and salt interferes with lipid complexing, decreasing the transfection efficiency.

The PureLink<sup>™</sup> HiPure Plasmid Prep Kits, available separately from Invitrogen, allow the purification of all types and sizes of plasmid DNA, including BAC, bacmids, and ssM13 DNAs. PureLink<sup>™</sup> HiPure Plasmid Prep Kits are ideally suited for bacmid purification (see page 53 for ordering information).

## Isolating Recombinant Bacmid DNA

Introduction	The PureLink <sup>™</sup> HiPure Plasmid DNA Miniprep Kit allows you to purify high quality bacmid DNA from DH10Bac <sup>™</sup> <i>E. coli</i> (see page 53 for ordering information). The isolated bacmid DNA is suitable for use in insect cell transfections. Note: We do not recommend the PureLink <sup>™</sup> HiPure Precipitator Module or the PureLink <sup>™</sup> HiPure Plasmid Filter Mini/Midi/Maxiprep Kits for isolating bacmid DNA.
Before Starting	<ul> <li>Inoculate a single white bacterial colony into 2 mL LB medium with 50 μg/mL kanamycin, 7 μg/mL gentamicin, and 10 μg/mL tetracycline. Incubate the culture at 37°C in a shaking water bath at 250 rpm overnight.</li> <li>Verify that RNase A is added to the Resuspension Buffer (R3) and that the Lysis Buffer (L7) contains no precipitates.</li> </ul>
Equilibrating the Column	Place the PureLink <sup>™</sup> HiPure Mini column on the PureLink <sup>™</sup> Nucleic Acid Purification Rack (see the manual supplied with the rack for more details). Apply 2 mL Equilibration Buffer (EQ1) to the column. Allow the solution in the column to drain by gravity flow.
Preparing the Cell Lysate	<ol> <li>Harvest 1.5 mL bacterial cells by centrifuging at 9,000 × g for 15 minutes. Remove all medium.</li> <li>Add 0.4 mL Resuspension Buffer (R3) containing RNase A to the pellet and resuspend the cells until homogeneous. Transfer the cell suspension to a centrifuge tube.</li> <li>Add 0.4 mL Lysis Buffer (L7). Mix gently by inverting the capped tube five times. Do not vortex. Incubate at room temperature for 5 minutes.</li> <li>Add 0.4 mL Precipitation Buffer (N3) and mix immediately by inverting the capped tube until the mixture is homogeneous. Do not vortex.</li> <li>Centrifuge the mixture at &gt;15,000 × g at room temperature for 10 minutes. Note: If the pellet does not adhere to the bottom of the tube, incubate the tube at room temperature for 5 minutes to allow the separation of the lysate and gelatinous pellet. Pipet the clear lysate into a sterile tube and centrifuge at &gt;15,000 × g for 5 minutes at room</li> </ol>
Binding and Washing the DNA	<ol> <li>Load the supernatant resulting from Step 5 (see above) onto the equilibrated column. Allow the solution in the column to drain by gravity flow.</li> <li>Wash the column twice with 2.5 mL Wash Buffer (W8). Allow the solution in the column to drain by gravity flow after each wash. Discard the flow-through.</li> </ol>

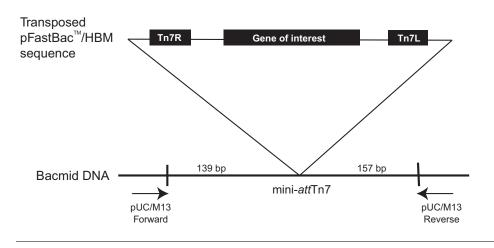
## Isolating Recombinant Bacmid DNA, Continued

Eluting and	1. Place a sterile centrifuge tube (elution tube) under the column.
Precipitating DNA	<ol> <li>Add 0.9 mL Elution Buffer (E4) to the column to elute DNA. Allow the solution to drain by gravity flow. Do not force out any remaining solution.</li> <li>The elution tube contains the purified DNA. Discard the column.</li> </ol>
	3. Add 0.63 mL isopropanol to the elution tube. Mix, then place the tube on ice for 10 minutes.
	4. Centrifuge the mixture at >15,000 $\times$ g at 4°C for 20 minutes. Carefully remove and discard the supernatant.
	5. Resuspend the DNA pellet in 1 mL 70% ethanol.
	6. Centrifuge at >15,000 × g at 4°C for 5 minutes. Carefully remove and discard the supernatant.
	7. Air-dry the pellet for 10 minutes.
	<ol> <li>Resuspend the DNA pellet in 40 μL TE Buffer (TE). Allow the pellet to dissolve for at least 10 minutes on ice. To avoid shearing the DNA, pipet only 1 or 2 times to resuspend.</li> </ol>
	9. Store the bacmid DNA at <b>4°C</b> .
Important	You may store your bacmid DNA at –20°C if you avoid frequent freeze/thaw cycles, which decrease the transfection efficiency. To store your purified bacmid DNA at –20°C, aliquot the bacmid DNA into separate tubes in TE Buffer, pH 8.0 to avoid more than one freeze/thaw cycle. Do <b>not</b> store the bacmid DNA in a frost-free freezer. You may also store the purified bacmid DNA for up to 2 weeks at 4°C in TE Buffer, pH 8.0.
	You may prepare glycerol stocks of DH10Bac <sup>™</sup> <i>E. coli</i> containing the bacmid DNA from mid-logarithmic phase culture grown from white colonies picked during the blue-white screening, and store the glycerol stocks at –80°C for future bacmid DNA isolation.

#### Analyzing Recombinant Bacmid DNA by PCR

#### Introduction

Recombinant bacmid DNA is greater than 135 kb in size. Since restriction analysis is difficult to perform with DNA of this size, we recommend using PCR analysis to verify the presence of your gene of interest in the recombinant bacmid. Use the pUC/M13 Forward and Reverse primers (sequences given below) that hybridize to sites flanking the mini-*att*Tn7 site (see figure below). This section provides guidelines and instructions for performing PCR using the pUC/M13 Forward and Reverse primers.



# PCR Analysis with pUC/M13 Primers

To verify the presence of your gene of interest in the recombinant bacmid using PCR, you may:

- Use the pUC/M13 Forward and Reverse primers (see sequences below).
- Use a combination of the pUC/M13 Forward or Reverse primer and a primer that hybridizes within your insert.

Invitrogen does not supply the pUC/M13 Forward and Reverse primers; you must have these primers custom synthesized.

Primer	Sequence
pUC/M13 Forward	5'-CCCAGTCACGACGTTGTAAAACG-3'
pUC/M13 Reverse	5'-AGCGGATAACAATTTCACACAGG-3'

**DNA Polymerase** You may use any DNA polymerase of your choice for PCR including Platinum<sup>®</sup> *Taq* DNA Polymerase. If the expected PCR product is > 4 kb, we recommend using a polymerase mixture such as Platinum<sup>®</sup> *Taq* DNA Polymerase High Fidelity for best results. See page 53 for ordering information.

### Analyzing Recombinant Bacmid DNA by PCR, Continued

Generating the PCR Product	Use the procedure below to amplify your recombinant bacmid DNA using the pUC/M13 Forward and Reverse primers and Platinum <sup>®</sup> <i>Taq</i> polymerase. If you are using a combination of the pUC/M13 Forward or Reverse primers and a primer specific for your gene, determine the amplification conditions to use. If you are using another polymerase, follow the manufacturer's recommendations for the polymerase you are using.		
	Note: Amplification conditions may need to be optimized	ed if your insert is > 4 kb.	
	1. For each sample, set up the following 50 µL PC microcentrifuge tube:	CR reaction in a 0.5 mL	
	Recombinant bacmid DNA (100 ng)	1 µL	
	10X PCR Buffer (appropriate for enzyme)	5 µL	
	10 mM dNTP Mix	1 µL	
	50 mM MgCl <sub>2</sub>	1.5 µL	
	PCR Primers (1.25 μL each 10 μM stock)	2.5 μL	
	Sterile Water	38.5 µL	
	<u>Platinum<sup>®</sup> <i>Taq</i> polymerase (5 units/µL)</u>	0.5 µL	
	Total Volume	50 mL	

2. Amplify using the following cycling parameters:

Step	Time	Temperature	Cycles
Initial Denaturation	3 minutes	94°C	1X
Denaturation	45 seconds	94°C	
Annealing	45 seconds	55°C	25–35X
Extension	5 minutes	72°C	
Final Extension	7 minutes	72°C	1X

3. Remove 5–10  $\mu$ L from the reaction and analyze by agarose gel electrophoresis.

#### Analyzing Recombinant Bacmid DNA by PCR, Continued

# What You ShouldIf transposition has occurred and you have used the pUC/M13 Forward and<br/>Reverse primers for amplification, you should see a PCR product of the following<br/>size on the agarose gel:

Bacmid transposed with	Size of PCR Product
Bacmid alone	~300 bp
pFastBac <sup>™</sup> /HBM-TOPO <sup>®</sup>	~2,500 bp + size of your insert
pFastBac™ Gus	~4,200 bp

If you have used a combination of the pUC/M13 Forward or Reverse primer and a gene-specific primer for amplification, determine the expected size of your PCR product. Refer to the diagram on page 15 to help you calculate the expected size of your PCR product.

## Producing Recombinant Baculovirus

## **Transfecting Insect Cells**

Introduction	After you have confirmed that your recombinant bacmid contains the gene of interest, you are ready to transfect insect cells to produce recombinant baculovirus. This section provides guidelines and instructions for transfecting insect cells.	
Cellfectin <sup>®</sup> II Reagent	We recommend using a cationic lipid such as Cellfectin <sup>®</sup> II Reagent for transfection. Cellfectin <sup>®</sup> II Reagent is a proprietary cationic lipid formulation that offers the highest transfection efficiencies and protein expression levels on the widest variety of adherent and suspension insect cell lines, including Sf9 and Sf21 cells.	
	Cellfectin <sup>®</sup> II Reagent is supplied with the Bac-to-Bac <sup>®</sup> HBM TOPO <sup>®</sup> Secreted Expression System, and is also available separately from Invitrogen. See page 53 for ordering information.	
Insect Cell Lines for Transfection	We recommend using Sf9 or Sf21 cells for transfection and identification of recombinant plaques. High Five <sup>™</sup> and Mimic <sup>™</sup> Sf9 cells are not recommended because they generally transfect less efficiently. However, once you have generated your baculovirus stock, we recommend using High Five <sup>™</sup> cells for secreted expression studies (see Secretion in High Five <sup>™</sup> Cells, page 32).	
Media for Transfection	For the highest transfection efficiency, we recommend performing the transfection in Grace's Insect Cell Culture Medium, Unsupplemented (see page 54 for ordering information). Note that the Grace's Insect Cell Culture Medium <b>should not</b> contain supplements or fetal bovine serum (FBS), because the supplements and the proteins in the FBS interferes with the Cellfectin <sup>®</sup> II Reagent, inhibiting the transfection.	
	Note: If you are culturing Sf9 or Sf21 cells in Sf-900 II SFM or Sf-900 <sup>™</sup> III SFM, you can perform the transfection in unsupplemented Grace's Medium, and then easily switch back to Sf-900 II SFM or Sf-900 <sup>™</sup> III SFM after transfection.	
Positive Control	If you have generated a recombinant bacmid from the pFastBac <sup>TM</sup> Gus control plasmid, we recommend including this positive control in your transfection and expression experiments to help you evaluate your results. In this bacmid, the gene encoding $\beta$ -glucuronidase is expressed under the control of the strong polyhedrin (P <sub>H</sub> ) promoter. After transfection, you may assay expression of $\beta$ -glucuronidase as appropriate (see page 39).	

#### Transfecting Insect Cells, Continued

	(500 ng/ $\mu$ L in TE Buffer, pH 8.0)			
	<ul> <li>Purified recombinant bacmid DNA construct (if desired, 500 ng/µL in 7</li> </ul>	1		
	• Sf9 or Sf21 cells cultured in the appr	ropriate medium		
	• Cellfectin <sup>®</sup> II Reagent (store at 4°C u	intil use)		
	• Grace's Insect Cell Medium, Unsupplemented (see page 54), media should not contain supplements, FBS, or antibiotics			
	• 6-well tissue culture plates and othe	er tissue culture supplies		
	• 1.5 mL sterile microcentrifuge tubes	3		
	<ul> <li>Complete growth medium for culturing insect cells (<i>e.g.</i>, Sf-900 II SFM, Sf-900<sup>™</sup> III SFM, TNM-FH, Grace's Supplemented Insect Cell Culture Medium, or other suitable medium)</li> </ul>			
~MENO	Calculate the number of Sf9 or Sf21 cells that you need for your transfection experiment and expand cells accordingly. Make sure your cells are healthy with greater than 95% viability and are growing in the logarithmic phase with a density of $1.5 \times 10^6$ – $2.5 \times 10^6$ cells/mL before proceeding to transfection.			
	experiment and expand cells accordingl greater than 95% viability and are grow	y. Make sure your cells are healthy with ing in the logarithmic phase with a		
Transfection Conditions	experiment and expand cells accordingl greater than 95% viability and are grow	y. Make sure your cells are healthy with ing in the logarithmic phase with a efore proceeding to transfection. s in Sf9 or Sf21 cells using the following tions as a starting point for your fection efficiency and low non-specific by varying DNA and Cellfectin <sup>®</sup> II		
	experiment and expand cells accordingl greater than 95% viability and are grow density of 1.5 × 10 <sup>6</sup> –2.5 × 10 <sup>6</sup> cells/mL be We generally produce baculoviral stock transfection conditions. Use these condi transfection. To obtain the highest trans effects, optimize transfection conditions	y. Make sure your cells are healthy with ing in the logarithmic phase with a efore proceeding to transfection. s in Sf9 or Sf21 cells using the following tions as a starting point for your fection efficiency and low non-specific by varying DNA and Cellfectin <sup>®</sup> II		
	experiment and expand cells accordingly greater than 95% viability and are grow density of $1.5 \times 10^6$ – $2.5 \times 10^6$ cells/mL be We generally produce baculoviral stock transfection conditions. Use these condi- transfection. To obtain the highest trans- effects, optimize transfection conditions Reagent concentrations, and cell density	y. Make sure your cells are healthy with ing in the logarithmic phase with a efore proceeding to transfection. s in Sf9 or Sf21 cells using the following tions as a starting point for your fection efficiency and low non-specific by varying DNA and Cellfectin <sup>®</sup> II 7.		
	experiment and expand cells accordingl greater than 95% viability and are grow density of 1.5 × 10 <sup>6</sup> –2.5 × 10 <sup>6</sup> cells/mL be We generally produce baculoviral stock transfection conditions. Use these condi transfection. To obtain the highest trans effects, optimize transfection conditions Reagent concentrations, and cell density <b>Condition</b>	y. Make sure your cells are healthy with ing in the logarithmic phase with a efore proceeding to transfection. s in Sf9 or Sf21 cells using the following tions as a starting point for your fection efficiency and low non-specific by varying DNA and Cellfectin <sup>®</sup> II 7. Amount		

Amount of Cellfectin® II Reagent8 μL (can vary from 1.5 to 9 μL)Note: This procedure is for insect cells in a 6-well format. All amounts and volumes are<br/>given on a per well basis.

#### Important Guidelines for Transfection

- Use Grace's Insect Cell Culture Medium, Unsupplemented to seed all cells in plate for Sf9 and Sf21 cells grown in Grace's Insect Cell Culture Medium, Supplemented (with 10% FBS).
- With Cellfectin<sup>®</sup> II, you do not have to remove the medium from cells and wash cells prior to adding the DNA-lipid complex to cells.
- The DNA-lipid complex formation time is shorter (~15–30 minutes) when using Cellfectin<sup>®</sup> II as compared to Cellfectin<sup>®</sup> reagent.
- **Do not** add antibiotics during transfection. Adding antibiotics during transfection causes cell death.

## Transfecting Insect Cells, Continued

Transfection Procedure	For Sf9 or Sf21 insect cells cultured in Supplemented Grace's Insect Medium containing 10%FBS, use the following protocol to prepare your cells for transfection in a 6-well format. All amounts and volumes are given on a per well basis. To transfect cells in other tissue culture formats, first determine the optimal conditions to use.		
	1.	Verify that the Sf9 or Sf21 cells are in the log phase $(1.5 \times 10^6 - 2.5 \times 10^6 \text{ cells/mL})$ with greater than 95% viability.	
	2.	Prepare transfection samples:	
		• If the cell density is in range of $1.5 \times 10^{6}$ – $2.5 \times 10^{6}$ cells/mL and the culture is without antibiotics:	
		a. Add 2 mL of Grace's Insect Medium, Unsupplemented (without antibiotics and serum) in each well.	
		<ul> <li>b. Seed 8 × 10<sup>5</sup> Sf9 or Sf21 cells from Step 1 per well. Do not change medium or wash the cells. The medium carried over will enhance the transfection efficiency.</li> </ul>	
		c. Allow cells to attach for 15 minutes at room temperature in the hood. Proceed to step 3.	
		• If the cell density is <b>not</b> in this range or the cell culture contains antibiotics:	
		<ul> <li>a. Prepare 10 mL of plating medium by mixing 1.5 mL Supplemented Grace's Insect Medium containing 10% FBS (without antibiotics) and 8.5 mL Grace's Insect Medium, Unsupplemented (without FBS and antibiotics).</li> </ul>	
		b. Plate $8 \times 10^5$ Sf9 or Sf21 cells from Step 1 per well.	
		c. Allow cells to attach for 15 minutes at room temperature in the hood.	
		d. Remove the medium. Add 2.5 mL plating medium from step 2a per well. Proceed to step 3.	
	3.	For each transfection sample, prepare complexes as follows:	
		a. Mix Cellfectin <sup>®</sup> II before use, and dilute 8 μL in 100 mL Grace's Medium, Unsupplemented (without antibiotics and serum). Vortex briefly to mix. <b>Note</b> : You may leave this mixture at room temperature for up to 30 minutes.	
		<ul> <li>b. Dilute 1 µg baculovirus DNA in 100 µL Grace's Medium, Unsupplemented (without antibiotics and serum). Mix gently.</li> </ul>	
		<ul> <li>Combine the diluted DNA with diluted Cellfectin<sup>®</sup> II (total volume ~210 μL). Mix gently and incubate for 15–30 minutes at room temperature.</li> </ul>	
	4.	Add ~210 µL DNA-lipid mixture or transfection mixture (Step 3c) <b>dropwise</b> onto the cells from Step 2a or 2c. Incubate cells at 27°C for 3–5 hours.	
	5.	Remove the transfection mixture and replace it with 2 mL of complete growth medium (e.g., Grace's Insect Medium, Supplemented and 10% FBS). Using antibiotics is optional.	
	6.	Incubate the cells at 27°C for 72 hours or until you see signs of viral infection.	

#### **Isolating P1 Viral Stock**

Introduction	Budded virus should be released into the medium 72 hours after transfection. However, if your transfection efficiency was not optimal, cells may not show all of the signs of viral infection until 4 or 5 days post-transfection. Beginning at 72 hours after transfection, visually inspect the cells daily for signs of infection (see below).
Characteristics of	Insect cells infected with baculovirus typically display the following characteristics

**Characteristics of Infected Cells** Insect cells infected with baculovirus typically display the following characteristics when visually observed using an inverted phase microscope at 250–400X magnification. The time points provided below assume that the transfection was successful (i.e., transfection efficiency was high).

Signs of Infection	Phenotype	Description
Early (first 24 hours)	Increased cell diameter	A 25–50% increase in cell diameter may be seen.
	Increased size of cell nuclei	Nuclei may appear to "fill" the cells.
Late (24–72 hours)	Cessation of cell growth	Cells appear to stop growing when compared to a cell-only control.
	Granular appearance	Signs of viral budding; vesicular appearance to cells.
	Detachment	Cells release from the plate or flask.
Very Late (>72 hours)	Cell lysis	Cells appear lysed, and show signs of clearing in the monolayer.

#### Preparing the P1 Viral Stock

1. When the transfected cells (from Step 6, previous page) demonstrate signs of late stage infection (e.g., 72 hours post-transfection), collect the medium containing the virus from each well (~2 mL) and transfer to sterile 15 mL snap-cap tubes. Centrifuge the tubes at 500 × g for 5 minutes to remove cells and large debris.

2. Transfer the clarified supernatant to fresh 15 mL snap-cap tubes. **This is the P1 viral stock**. Store at 4°C, protected from light. See the next page for additional storage information.

**Note:** To concentrate your viral stock to obtain a higher titer, filter your viral supernatant through a 0.2  $\mu$ m, low protein binding filter after the low-speed centrifugation step, if desired.

### Isolating P1 Viral Stock, Continued

Storing Viral Stocks	• Store viral stock at 4°C, protected from light.
	• If medium is serum-free (e.g., Sf-900 II SFM, Sf-900 <sup>™</sup> III SFM), add fetal bovine serum to a final concentration of 2%. Serum proteins act as substrates for proteases.
	• For long-term storage, store an aliquot of the viral stock at –80°C for later reamplification.
	• Do <b>not</b> store routinely used viral stocks at temperatures below 4°C. Repeated freeze/thaw cycles can result in a 10- to 100-fold decrease in virus titer.
The Next Step	Once you have obtained your clarified P1 baculoviral stock, you may:
	• Amplify the viral stock (see the next section for details). This procedure is recommended to obtain the highest viral titers and optimal results in your expression studies.
	• Determine the titer of your viral stock (see <b>Performing a Viral Plaque Assay</b> , page 25).
	• Plaque purify your recombinant baculovirus, if desired (see <b>Performing a Viral Plaque Assay</b> , page 25).
	• Use the P1 viral stock to infect your Sf9 or Sf21 cells for preliminary expression experiments (see below).
Note	To perform small-scale or preliminary expression experiments, it is possible to proceed directly to expression studies by using the P1 viral stock to infect your cells. Note that the MOI is unknown if viral titer is not determined, and the

conditions may not be reproducible.

amount of viral stock is limited without viral amplification; therefore, expression

# Amplifying Your Baculoviral Stock

Introduction	The P1 viral stock is a small-scale, low-titer stock. You may use this stock to infect cells to generate a high-titer P2 stock. The titer of the initial viral stock obtained from transfecting Sf9 or Sf21 cells generally ranges from $1 \times 10^6$ to $1 \times 10^7$ plaque forming units (pfu)/mL. Because amplification allows production of a P2 viral stock with a titer ranging from $1 \times 10^7$ to $1 \times 10^8$ pfu/mL, we generally recommend it. This section provides guidelines and protocols for amplifying the recombinant baculovirus to prepare a P2 viral stock.
Materials Needed	Sf9 or Sf21 cells cultured in the appropriate growth medium
	P1 baculoviral stock
	• Any appropriate tissue culture vessel (see <b>Important Note</b> below)
	<ul> <li>Tissue culture reagents</li> </ul>
	<ul> <li>27°C humidified incubator</li> </ul>
Important	To amplify your P1 viral stock, you may infect Sf9 or Sf21 cells growing in suspension or monolayer culture. Depending on your needs, you may amplify your P1 viral stock at any scale, but remember that you may be limited by the amount of P1 viral stock available.
	We generally amplify our P1 viral stock in a 10 mL suspension culture at $2 \times 10^6$ cells/mL or in 6-well tissue culture plates at $2 \times 10^6$ cells/well. Calculate the number of Sf9 or Sf21 cells that you need for infection and expand cells accordingly. Make sure that the cells are healthy, of low passage (5–20), and have >95% viability before proceeding to infection.
Multiplicity of Infection (MOI)	To amplify your viral stock, infect cells at a multiplicity of infection (MOI) ranging from 0.05 to 0.1. MOI is defined as the number of virus particles per cell. Use the following formula to calculate how much viral stock to add to obtain a specific MOI:
	Inoculum required (mL) = $\frac{MOI (pfu/cell) \times number of cells}{titer of viral stock (pfu/mL)}$
	<b>Note:</b> If you have not determined the titer of your P1 viral stock, you may assume that the titer ranges from $1 \times 10^6$ to $1 \times 10^7$ pfu/mL.
Example	To infect a 10 mL culture at $2 \times 10^6$ cells/mL with an MOI of 0.1 using a P1 viral stock at $5 \times 10^6$ pfu/mL:
	Inoculum required (mL) = $\frac{(0.1 \text{pfu/cell}) \times (2 \times 10^7 \text{cells})}{5 \times 10^6 \text{pfu/mL}}$
	Inoculum required (mL) = 0.4 mL

# Amplifying Your Baculoviral Stock, Continued

Important considerations	For successful amplification of your baculovirus, pay attention to several key points:
	<ul> <li>Use Sf9 or Sf21 cells that are in excellent health, low passage (5–20), log-phase growth, and have &gt;95% viability.</li> </ul>
	• Use sterile P1 baculoviral stock that is free of contaminants.
	• Use a low MOI between 0.05–0.1. Higher MOI reduces baculovirus quality.
	• Harvest the virus when 70–80% of cells are dead.
	• You <b>cannot</b> amplify the baculovirus indefinitely, because the baculovirus acquires deleterious mutations with each passage. Usually, P3 is highest usable passage.
Amplification	Follow the guidelines below to amplify your P1 viral stock in a 6-well plate.
Procedure	1. On the day of infection, prepare your Sf9 or Sf21 cell suspension and plate cells at $2 \times 10^6$ cells/well. Incubate cells at room temperature for 1 hour to allow attachment.
	2. After 1 hour, inspect cells under an inverted microscope to verify attachment.
	3. Add the appropriate amount of P1 viral stock to each well.
	4. Incubate the cells for 48 hours in a 27°C humidified incubator.
	5. 48 hours post-infection, collect 2 mL of medium containing virus from each well and transfer the virus to sterile 15 mL snap-cap tubes. Centrifuge the tubes at $500 \times g$ for 5 minutes to remove cells and large debris and to obtain clarified baculoviral stock.
	<b>Note:</b> It is possible to harvest virus at later times after infection (e.g., 72 hours). Because optimal harvest times can vary, determine them for each baculoviral construct. Remember that culture viability decreases over time as cells lyse.
	6. Transfer the supernatant to fresh 15 mL snap-cap tubes. This is the <b>P2 viral stock.</b> Store at 4°C, protected from light. For long-term storage, you may store an aliquot of the P2 stock at –80°C, protected from light. See page 22 for storage guidelines.
	7. Proceed to the next section to determine the titer of your P2 viral stock.
Scaling Up the Amplification Procedure	After you have generated a high-titer P2 baculoviral stock, you may scale-up the amplification procedure to any volume of your choice. To produce this high-titer P3 stock, scale up the amount of cells and volume of virus used appropriately, and follow the guidelines and procedure outlined in this section.
Generating High- Titer Stocks From Frozen Master Stock	If you have stored your viral master stock at $-80^{\circ}$ C, we recommend amplifying this stock to generate another high-titer stock for use in expression experiments. Viral titers generally decrease over time when virus is stored at $-80^{\circ}$ C. Follow the guidelines and amplification procedure detailed in this section.

# Performing a Viral Plaque Assay

Introduction	We recommend you perform a plaque assay to determine the titer of your viral stock. You may also perform a plaque assay to purify a single viral clone, if desired. In this procedure, you infect cells with dilutions of your viral stock and identify focal points of infection (plaques) on an agarose overlay. You may also titer your viral stock by the end-point dilution method described in O'Reilly <i>et. al.</i> , 1992
Experimental	To determine the titer of a baculoviral stock:
Outline	1. Plate Sf9 or Sf21 cells in 6-well plates.
	2. Prepare 10-fold serial dilutions of your baculoviral stock.
	3. Add the different dilutions of baculovirus to Sf9 cells and infect cells for 1 hour.
	4. Remove the virus and overlay the cell monolayer with Plaquing Medium.
	5. Incubate the cells for 7–10 days, stain (if desired), and count the number of plaques in each dilution.
Factors Affecting	A number of factors can influence viral titers including:
Viral Titer	• <b>The size of your gene of interest:</b> Titers generally decrease as the size of the insert increases.
	• The transfection efficiency: For the highest transfection efficiency, we recommend transfecting Sf9 or Sf21 cells using Cellfectin <sup>®</sup> II Reagent. Prepare DNA: lipid complexes in Grace's Insect Medium, Unsupplemented (see pages 18–20 for details).
	• The age of your baculoviral stock: Viral titers may decrease with long-term storage at 4°C or -80°C. If your baculoviral stock has been stored for 6 months to 1 year, we recommend titering or re-titering your baculoviral stock prior to use in an expression experiment.
	• <b>The number of freeze/thaw cycles:</b> If you are storing your viral stock at – 80°C, viral titers can decrease as much as 10% with each freeze/thaw cycle.
	• <b>Improper storage of your baculoviral stock:</b> For routine use, baculoviral stocks should be aliquoted and stored at 4°C, protected from light.
	Continued on next page

Materials Needed	• Your clarified baculoviral stock (store at 4°C until use)
	• Sf9 or Sf21 cells cultured in the appropriate medium (30 mL of log-phase cells at $5 \times 10^5$ cells/mL for each baculoviral stock to be titered)
	• Sf-900 II SFM, Sf-900 <sup>™</sup> III SFM or other appropriate complete growth medium (see <b>Note</b> below)
	• Sf-900 Medium (1.3X) (100 mL) or other appropriate plaquing medium (see <b>Note</b> below)
	• 4% Agarose Gel (specifically formulated for optimal insect cell growth)
	• Sterile, cell-culture grade, distilled water
	• 100 mL sterile, glass bottle
	• 6-well tissue-culture plates (2 plates for each viral stock to be titered)
	Sterile hood
	• Waters baths at 40°C and 70°C
	Microwave oven (optional)
	• 27°C humidified incubator
	• Neutral Red (Sigma, Cat. no. N7005)
	See page 54 for ordering information.
Note	If you are culturing your Sf9 or Sf21 cells in serum-supplemented media (i.e., complete TNM-FH), you should have the following reagents on hand:
	Grace's Insect Cell Culture Medium, Supplemented
	Grace's Insect Cell Culture Medium (2X)
	• Fetal Bovine Serum (FBS), Qualified, Heat-Inactivated

See page 54 for ordering information.

Preparing the Plaquing Medium	Plaquing medium consists of a mixture of culture medium and agarose. Plaquing medium is used to immobilize the infected cells for the plaque assay. Prepare plaquing medium immediately before use, following the procedure below. If you are culturing the Sf9 cells in Sf-900 II SFM or Sf-900 <sup>™</sup> III SFM, prepare Sf-900 Plaquing Medium. If you are culturing cells in TNM-FH, prepare Grace's Plaquing Medium. Note: Other Plaquing Media are suitable.		
	<ol> <li>Melt the 4% Agarose Gel by placing the bottle in a 70°C water bath for 20 to 30 minutes or heating the agarose in a microwave oven. While the 4% agarose gel is melting, place the following in the 40°C water bath:</li> </ol>		
	• Empty, sterile 100 mL bottle		
	• Sf-900 Medium (1.3X) or Grace's Insect Cell Culture Medium (2X), as appropriate		
	2. After the 4% agarose gel has liquefied, move the agarose gel, medium, and empty 100 mL bottle to a sterile hood.		
	3. Working quickly, prepare the plaquing medium as follows:		
	<b>Sf-900 Plaquing Medium:</b> Combine 30 mL of Sf-900 Medium (1.3X) and 10 mL of the melted 4% Agarose Gel in the empty 100 mL bottle and mix gently.		
	<b>Grace's Plaquing Medium:</b> Add 20 mL of heat-inactivated FBS to the 100 mL bottle of Grace's Insect Medium (2X) and mix. Combine 25 mL of the Grace's Insect Medium (2X) containing serum with 12.5 mL of cell-culture grade, sterile, distilled water and 12.5 mL of the melted 4% Agarose Gel in the empty 100 mL bottle and mix gently.		
	4. Return the bottle of plaquing medium to the 40°C water bath until use.		
	Continued on next page		

Plaque Assay Procedure	Use the procedure below to perform a plaque assay in 6-well plate format to determine the titer of your pFastBac <sup>™</sup> /HBM baculoviral stock. If you have generated a baculoviral stock of the expression control pFastBac <sup>™</sup> Gus, we recommend titering this stock as well. Remember to include a negative control (no virus) in your experiment. <b>Note:</b> The amounts provided in this procedure are suitable to titer one baculoviral stock (two 6-well plates per viral stock). To titer more than one baculoviral stock, scale up the reagent quantities accordingly.		
	1.	On the day of infection, harvest Sf9 or Sf21 cells and prepare a 30 mL cell suspension at $5 \times 10^5$ cells/mL in Sf-900 II SFM (or other complete growth medium). Aliquot 2 mL of cell suspension into each well of two 6-well plates. If you are including a negative control, you need another 6-well plate.	
	2.	Allow the cells to settle to the bottom of the plate and incubate, covered, at room temperature for 1 hour.	
	3.	Following the 1 hour incubation, observe the cell monolayers using an inverted microscope. Sf9 cells should be attached and at 50% confluence.	
	4.	Prepare an 8-log serial dilution (10 <sup>-1</sup> to 10 <sup>-8</sup> ) of the clarified baculoviral stock in Sf-900 II SFM <b>or</b> Grace's Insect Cell Culture Medium, Supplemented, without FBS, as appropriate.	
		Sequentially dilute 0.5 mL of the baculoviral stock or previous dilution in 4.5 mL of medium in 12 mL disposable tubes, finishing with 8 tubes of diluted viral stock (i.e., $10^{-1}$ , $10^{-2}$ , $10^{-3}$ , $10^{-4}$ , $10^{-5}$ , $10^{-6}$ , $10^{-7}$ , $10^{-8}$ ). Use the dilutions $10^{-4}$ to $10^{-8}$ in your assay.	
	5.	Move the 6-well plates containing Sf9 cells and the tubes of diluted virus to the sterile hood. Label the plates, in columns of 2 (1 sample well plus 1 duplicate) as follows: no virus (negative control), 10 <sup>-4</sup> , 10 <sup>-5</sup> , 10 <sup>-6</sup> , 10 <sup>-7</sup> , 10 <sup>-8</sup> .	
	6.	Remove the medium from each well, discard it, and immediately replace it with 1 mL of the appropriate virus dilution. As a negative control, add the appropriate medium without virus.	
	7.	Incubate the cells with the virus for 1 hour at room temperature.	
	8.	Move the cells and the bottle of plaquing medium from the 40°C water bath (Step 4, previous page) to a sterile hood.	
	9.	Sequentially starting from the highest dilution $(10^{-8})$ to the lowest dilution $(10^{-4})$ , remove the medium containing virus from the wells and replace it with 2 mL of plaquing medium. Work quickly to avoid desiccation of the cell monolayer.	
	10.	Allow the agarose overlay to harden for 1 hour at room temperature before moving the plates.	
	11.	Incubate the cells in a 27°C humidified incubator for 7–10 days until plaques are visible and ready to count. To stain plaques to facilitate counting, see the next page. To calculate the titer, see page 30.	
		Continued on next page	



To improve the visualization of plaques, stain the plates using Neutral Red. **Crystalline Blue and other plaque staining dyes containing organic solvents are not recommended because they kill the host cells**. To stain plaques, you may do one of the following:

- Prepare an agarose solution containing Neutral Red and overlay this solution on the plates 4 days post-infection. Count plaques 7–10 days post-infection.
   or
- Prepare a Neutral Red solution and add it to plates for 1–2 hours just prior to counting plaques (7–10 days post-infection).

**Important:** If you plan to plaque purify your baculovirus, **do not** stain plaques, because Neutral Red is a known mutagen that can alter your recombinant virus.

Neutral Red Staining Procedure

#### Preparing a Neutral Red Agarose Overlay (for use on Day 4)

- 1. Prepare a 1 mg/mL Neutral Red solution in Sf-900 II SFM (or other appropriate complete growth medium). Filter-sterilize.
- 2. Combine the reagents below in a 50 mL tube and place in a 40°C water bath.

1 mg/mL Neutral Red solution	1.5 mL
Sf-900 II SFM	16.5 mL

- 3. Microwave 4% Agarose Gel until melted, then place it in a 40°C water bath for 5 minutes.
- 4. Move the 50 mL tube of Neutral Red solution and the 4% agarose gel to a sterile hood. Add 6 mL of 4% agarose gel to the Neutral Red solution.
- 5. Add 1 mL of the Neutral Red overlay to each well containing plaquing overlay. Once the agarose has hardened, return plates to a 27°C humidified incubator until plaques are ready to count. Plaques appear as clear spots on a red monolayer.

#### Preparing a Neutral Red Stain (for use on Day 7–10 prior to counting plaques)

- 1. Prepare a 1 mg/mL Neutral Red solution in cell-culture grade, distilled water.
- 2. Add 0.5 mL of Neutral Red solution to each well containing plaquing overlay. Incubate for 1 to 2 hours at room temperature.
- 3. Gently remove excess stain with a pipette or blotter and count the plaques. Plaques will appear as clear spots in a nearly clear gel against a red background.

Calculating the Titer	Count the number of plaques present in each dilution, then use the following formula to calculate the titer (plaque forming units (pfu)/mL) of your viral stock. Note that the optimal range to count is 3 to 20 plaques per well of a 6-well plate.
	titer (pfu/mL) = number of plaques × dilution factor × $\frac{1}{mL \text{ of inoculum/well}}$
Example	If you add 1 mL of inoculum and observe 20 plaques in the well containing the $10^{-6}$ viral dilution, the titer of the viral stock is:
	titer (pfu/mL) = 20 plaques × $10^6$ × $\frac{1}{1 \text{ mL of inoculum/well}}$
	titer (pfu/mL) = $2 \times 10^7$ pfu/mL
What You Should See	When titering pFastBac <sup>™</sup> /HBM baculoviral stocks, we generally obtain titers ranging from:
	• $1 \times 10^6$ to $1 \times 10^7$ pfu/mL for P1 viral stocks
	• $1 \times 10^7$ to $1 \times 10^8$ pfu/mL for P2 viral stocks
	<b>Note:</b> If the titer of your baculoviral stock is less than $1 \times 10^6$ pfu/mL or $1 \times 10^7$ pfu/mL for a P1 or P2 viral stock, respectively, we recommend producing a new baculoviral stock.
	For tips and guidelines to optimize your viral yield, see <b>Factors Affecting Viral Titer</b> , page 25, and the <b>Troubleshooting</b> section, page 47.
	Continued on next page

Plaque Purification	You may generate a viral stock from a single viral clone by plaque purifying your baculovirus, if desired. Use a protocol of your choice or the procedure below.		
	Materials Needed		
	• Plate containing well-spaced viral plaques (from <b>Plaque Assay Procedure</b> , Step 11, page 28; <b>do not</b> stain plates with Neutral Red)		
	• Log phase Sf9 or Sf21 cells at greater than 95% viability		
	Sterile Pasteur pipette and bulb		
	Procedure		
	1. Follow Steps 1–3 in the <b>Plaque Assay Procedure</b> , page 28, to seed Sf9 or Sf21 cells.		
	2. Using a sterile Pasteur pipette and bulb, carefully pick a clear plaque and transfer the agarose plug (containing virus) to a 1.5 mL microcentrifuge tube containing 500 μL of complete growth medium. Mix well by vortexing.		
	3. Add 100 $\mu$ L of the agarose plug solution to each well.		
	4. Incubate the cells in a 27°C humidified incubator for 72 hours.		
	5. Collect the medium containing virus from each well (~2 mL) and transfer to sterile 15 mL snap-cap tubes. Centrifuge the tubes at 500 × g for 5 minutes to remove cells and large debris.		
	6. Transfer the clarified supernatant to fresh 15 mL snap-cap tubes. This is your plaque-purified viral stock.		
	7. Proceed to <b>Amplifying Your Baculoviral Stock</b> , page 23.		

## Expressing Your Recombinant Protein

Introduction	Once you have generated a pFastBac <sup><math>M</math></sup> /HBM baculoviral stock with a suitable titer (e.g., $1 \times 10^8$ pfu/mL), you are ready to use the baculoviral stock to infect High Five <sup><math>M</math></sup> insect cells and assay for expression of your recombinant protein. The following guidelines and recommendations are provided for your convenience. If you need more details about the techniques discussed, refer to <i>Current Protocols in Molecular Biology</i> , Unit 16.9-16.11 (Ausubel, <i>et al.</i> , 1994), <i>The Baculovirus Expression System: A Laboratory Guide</i> (King and Possee, 1992), or <i>Baculovirus Expression Vectors: A Laboratory Manual</i> (O'Reilly, <i>et al.</i> , 1992).
	<ul> <li>We recommend that you:</li> <li>Use High Five<sup>™</sup> cells adapted to suspension culture in serum-free medium for expression of secreted proteins</li> <li>Perform a time course of expression to determine the maximum point of expression</li> <li>Have a detection method for your protein</li> </ul>
High Five <sup>™</sup> Cells	<ul> <li>High Five<sup>™</sup> cells (see page 54 for ordering information) are particularly well suited for expression of secreted recombinant proteins. This cell line (BT1-TN-5B1-4) was originally developed by the Boyce Thompson Institute, Ithaca, NY and originated from the egg cells of the cabbage looper, <i>Trichoplusia ni</i>, the native host of AcMNPV (Davis <i>et al.</i>, 1992). This cell line has the following characteristics:</li> <li>Grows well in monolayer and doubles in less than 24 hours for ease of use</li> <li>Adaptable to suspension culture and serum-free medium for high-level protein expression and purification</li> <li>Provides 5–10 fold higher secreted expression than Sf9 cells (Davis <i>et al.</i>, 1993)</li> <li>For more information about High Five<sup>™</sup> cells or a protocol for adaptation to suspension culture, refer to the Growth and Maintenance of Insect Cell Lines manual (part no. 25-0127) available at www.invitrogen.com or from Technical Support (see page 56).</li> </ul>
Secretion in High Five <sup>™</sup> Cells	Expression of secreted alkaline phosphatase (SEAP) using its native secretion signal was evaluated in eight different cell lines including High Five <sup>™</sup> , Sf9, and Sf21. On a per cell basis, High Five <sup>™</sup> cells produced 20-fold more protein than Sf21 cells and 23-fold more protein than Sf9. Since High Five <sup>™</sup> cells are larger than Sf9 or Sf21 cells and the assays were performed on adherent cells, the amount of SEAP was also determined per milliliter of culture medium. In this case, High Five <sup>™</sup> cells produced 5-fold more SEAP than Sf9 cells and 8-fold more than Sf21 (Davis <i>et al.</i> , 1993).

Guidelines for Expression	<ul> <li>General guidelines are provided below to infect insect cells with the recombinant baculovirus to express your protein of interest. As with any expression system, optimal expression conditions depend on the characteristics of the protein being expressed.</li> <li>Cell line: We recommend using High Five<sup>™</sup> cells (see page 54 for ordering information) for expression of secreted recombinant proteins. Note: You may also use other cell lines such as Sf9, Sf21, or Mimic<sup>™</sup> Sf9, but your secreted expression levels will be lower.</li> <li>Culture Conditions: We generally culture High Five<sup>™</sup> cells in serum-free conditions using in Express Five<sup>®</sup> SFM (see page 54 for ordering information). You may grow your cells either in adherent or suspension culture using your culture vessel of choice.</li> <li>Depending on your application and the protein of interest, note that it may be necessary to supplement the culture post-infection with 0.1% to 0.5% FBS or BSA to protect the recombinant protein from proteolysis. Protein-based</li> </ul>
	<ul> <li>protease inhibitors are generally less expensive and more effective than many synthetic protease inhibitors.</li> <li>Infection Conditions: We recommend infecting cultures while cells are in the mid-logarithmic phase of growth at a density of 1 × 10<sup>6</sup> to 2 × 10<sup>6</sup> cells/mL. Make sure that the culture is not rate-limited by nutritional (<i>i.e.</i>, amino acid or carbohydrate utilization) or environmental factors (<i>i.e.</i>, pH, dissolved O<sub>2</sub>, or temperature) during infection.</li> <li>MOI: Optimal MOI will vary between cell lines, and the relative infection kinetics of the virus isolate or clone used. Establish a dose for each virus, medium, reactor, and cell line employed to determine the optimal infection parameters to use for protein expression. As a starting point, infect cells using an MOI of 5 to 10.</li> <li>Note: This MOI recommendation is unlike the generation of a high-titer stock, where a low MOI of 0.5–1.0 is recommended.</li> </ul>
	<ul> <li>Time course: We recommend performing a time course to determine the expression kinetics for your recombinant protein as many proteins may be degraded by cellular proteases released in cell culture. Maximum expression of secreted proteins is generally observed between 30 and 72 hours.</li> </ul>
Positive Control	If you have generated a high-titer viral stock from pFastBac <sup>™</sup> Gus control plasmid, include this recombinant baculovirus in your experiments for use as an expression control. After you have infected cells with the control virus, the gene encoding β-glucuronidase is constitutively expressed. The molecular weight of β-glucuronidase is 68.5 kDa. For a rapid but qualitative assay for β-glucuronidase expression, see page 39.

Seeding Densities	The table below gives approximate seeding densities and volumes for typical
and Volumes for	vessel sizes. Infection at these densities in the minimal volumes listed yields
Infections	optimal infection.
	<b>Minimal Volumes:</b> The total volumes used are lower than those used in general

al cell culture and maintenance, so that the virus added is concentrated and can infect cells more readily.

Cell Density: Cell density in adherent culture is approximately 50% confluent to allow maximal cell surface area for contact with virus and subsequent infection. However, to maximize the infection efficiency, we recommend that you determine the optimal cell density for the specific cell type you are using.

**MOI**: Use an MOI of 5–10 for a time course of protein expression or a large-scale protein preparation.

Type of Vessel	Cell Density	Final Volume (culture medium + added virus)
96-well plate	$3.0 \times 10^4$ cells/well	100 μL
24-well plate	$2.0 \times 10^5$ cells/well	500 μL
12-well plate	$4.0 \times 10^5$ cells/well	750 μL
6-well plate	$1.0 \times 10^6$ cells/well	1 mL
60 mm <sup>2</sup> plate	$2.5 \times 10^6$ cells/plate	3 mL
25 cm <sup>2</sup> flask	$2.0 \times 10^6$ cells/flask	5 mL
75 cm² flask	$6.0 \times 10^6$ cells/flask	10 mL
150 cm <sup>2</sup> flask	$1.2 \times 10^7$ cells/flask	15–20 mL
spinners (all)	$2.0 \times 10^6$ – $2.5 \times 10^6$ cells/mL	no more than half of the total volume of the flask

Amount of Virus to Add: The amount of virus to add depends on MOI.

**Calculating Virus** 

To calculate the volume of viral stock needed to achieve a given MOI;

#### Volumes

(MOI desired) (Total number of cells) Volume of virus = Titer of viral stock

For example, to infect a spinner with 50 mL of culture at a cell density of  $2 \times 10^6$  cells/mL (i.e., a total of  $1 \times 10^8$  cells) at an MOI of 5 using a high-titer virus stock at  $1 \times 10^8$  pfu (virions)/mL, you need:

- = 5 mL of your viral stock

(5 virions/cell)  $(1 \times 10^8 \text{ cells})$ 

 $1 \times 10^8$  virions/mL)

Determining Time Points for Optimal Secreted Protein Expression	When expressing a secreted protein, analyze the <b>supernatant</b> for the presence of <b>secreted</b> , <b>recombinant protein</b> and the lysates from the <b>cell pellet</b> for the presence of <b>unprocessed recombinant protein</b> . Compare the supernatant sample and the lysate sample to determine if recombinant protein is being secreted, how much protein has been secreted, and how much protein remains intracellular at different times during secretion. Using this data, optimize your time points for maximal secreted protein expression levels (see <b>Optimizing Protein Expression</b> , page 41). <b>Note</b> : β-glucuronidase expression from the pFastBac <sup>™</sup> Gus positive control baculovirus is intracellular (i.e., not secreted), because the gene product lacks the HBM secretion signal sequence.	
Protocol for Cells in Suspension Culture	The following procedure is designed to allow expression analysis from 50 mL of High Five <sup>™</sup> cells at a density of 2 × 10 <sup>6</sup> cells/mL cultured in a 100 mL spinner flask. <b>Note:</b> Use cells with a doubling time of 18–24 hours and a viability of 95%. Cells should be at a passage number less than 30, and they should not have been in spinner culture for more than 2 months.	
	Materials needed:	
	• High-titer pFastBac <sup><math>m</math></sup> /HBM baculoviral stock of known titer ( $\geq 10^8$ pfu/mL)	
	<ul> <li>High Five<sup>™</sup> cells</li> </ul>	
	Express Five <sup>®</sup> SFM	
	<ul> <li>One 100 mL spinner flask per sample</li> </ul>	
	<ul> <li>Microcentrifuge tubes</li> </ul>	
	Protocol:	
	<ol> <li>Add the pFastBac<sup>™</sup>/HBM baculoviral stock to the spinner flask at the desired MOI. Include the appropriate controls [mock-infected (uninfected) cells, pFastBac<sup>™</sup> Gus positive control baculovirus, previously characterized recombinant baculoviruses]. To calculate the amount of virus to add, see Calculating Virus Volumes, page 34.</li> </ol>	
	2. Incubate spinners at 27°C with a spin rate of 80 to 90 rpm.	
	3. Remove 1 mL aliquots of cells at designated time point(s) (see <b>Determining</b> <b>Time Points for Optimal Secreted Protein Expression</b> , above), and transfer each sample to a microcentrifuge tube.	
	4. Pellet cells at 800 × g for 10 minutes at 4°C. Keep samples at 4°C or on ice to prevent proteolysis.	
	5. Transfer supernatant to a new tube. Do not discard the cell pellet. Label each tube containing the supernatant and cell pellet.	
	<ol> <li>Proceed to Analyzing Recombinant Protein, page 37, or store the cell pellet and the supernatant at -80°C for analysis at a later time.</li> <li>Note: Storage at -80°C will reduce proteolysis of the recombinant protein. If you are performing a time course, lyse the cell pellet later when you have collected all time points.</li> </ol>	

#### The following procedure, adapted from Luckow and Summers, is designed to **Protocol for Adherent Cells** allow expression analysis in a 24-well format from recombinant baculovirus infected adherent cells harvested 24 to 96 hours post-infection. For other plate and/or flask sizes, adjust the cell seeding densities and volumes. Other protocols are also suitable. Materials needed: High-titer pFastBac<sup>M</sup>/HBM baculoviral stock of known titer ( $\geq 10^8$ pfu/mL) High Five<sup>™</sup> cells Express Five® SFM One 24-well plate per sample Microcentrifuge tubes **Protocol:** 1. Seed $2 \times 10^5$ High Five<sup>TM</sup> cells per well in a 24-well plate. Let cells attach for at least 30 minutes. 2. Remove the media and rinse the cells once with fresh growth media. Replace with 300 mL of fresh media. 3. Add the pFastBac<sup>M</sup>/HBM baculoviral stock to each well at the desired MOI. Include the appropriate controls [mock-infected (uninfected) cells, pFastBac<sup>™</sup> Gus positive control baculovirus, previously characterized recombinant baculoviruses]. To calculate the amount of virus to add, see Calculating Virus Volumes, page 34. 4. Incubate cells in a 27°C humidified incubator. 5. Harvest each well at the designated time point. Scrape the cells from each well and transfer the entire solution from each well to a microcentrifuge tube. 6. Pellet the cells at $800 \times g$ for 10 minutes at 4°C. Keep the samples at 4°C or on ice to prevent proteolysis. Transfer supernatant to a new tube. Label each tube. 7. 8. Proceed to Analyzing Recombinant Protein, page 37, or store the cell pellet and the supernatant at -80°C for analysis at a later time.

#### **Analyzing Recombinant Protein**

Introduction	The next step after harvesting baculovirus infected insect cells is to analyze the secreted expression of your protein by SDS-PAGE or western blot. In addition to analyzing the supernatant, we recommend that you analyze the cell lysate to determine if all of your protein is being secreted. Analyzing cell lysate can assist you in optimizing your MOI and time course for expression (see <b>Optimizing Protein Expression</b> , page 41). After you determine the optimal experimental parameters for secreted expression, you may proceed with <b>Large-Scale Expression</b> (page 42).			
Protease Inhibitors	We recommend that you add one or more protease inhibitors to each of the lysis buffers that are described in the protocol below. The following table summarizes recommended protease inhibitors, their method of action, and working concentrations.ProteaseMethod of ActionStock SolutionWorking			
	Inhibitor	Method of Action	Stock Solution	Concentration
	PMSF	Serine protease inhibitor	10 mg/mL in isopropanol	100 µg/mL
	Leupeptin	Serine and thiol protease inhibitor	50 μg/mL in deionized water	0.5 μg/mL
	Aprotinin	Serine protease inhibitor	50 μg/mL in deionized water	0.5 μg/mL
	Pepstatin A	Acid protease inhibitor	100 μg/mL in methanol	1μg/mL

You can store all of the above protease inhibitor solutions at -20°C except for PMSF. Store PMSF at room temperature in isopropanol. PMSF is not stable in aqueous solution; add it to the lysis buffer just before use.



**PMSF** (phenylmethylsulfonylfluoride) is very harmful if inhaled, swallowed, or contacted by the skin. Wear protective clothing and gloves when handling.

#### Preparing Cell Lysates

You may use any method to prepare cell lysates for analysis, including detergent lysis, sonication, or freeze-thaw lysis. The protocol on the next page, **Detergent Lysis**, provides a quick procedure for preparing lysates suitable for analyzing the secretion level of your recombinant protein. If you do not want to use detergent to lyse your cell samples (e.g., your protein is sensitive to detergent), lyse your cells by sonication and/or freeze-thawing.

## Analyzing Recombinant Protein, Continued

Detergent Lysis	Detergent lysis is a quick and efficient way to lyse cells and extract intracellular protein. The protocol below uses Triton <sup>®</sup> X-100, but you may also use Nonidet P-40 (NP-40).				
	1. Place all cell pellets from the time course on ice. Be sure to include the control sample.				
	2. Make up 2–5 mL of lysis buffer (0.1% Triton <sup>®</sup> X-100 in PBS or TBS). Use 100 $\mu$ L of lysis buffer for each 10 <sup>6</sup> cells.				
	3. Add each of the protease inhibitors (Leupeptin, Aprotinin and Pepstatin A) at the working concentrations described on the previous page. Perform this step on ice. Add PMSF just after adding the lysis buffer to the cell pellet (Step 5).				
	4. Add 100 $\mu$ L of lysis buffer for each 10 <sup>6</sup> cells in the pellet.				
	5. Add PMSF to each sample to a final concentration of $100 \mu\text{g/mL}$ .				
	6. Vortex each cell sample to break up the cell pellet and begin lysis.				
	7. <b>Lysis/Incubation:</b> Incubate all samples on ice for 30–45 minutes, then vortex them at 10 minute intervals to assist lysis.				
	<ol> <li>After all samples have been lysed, pellet cellular debris at 1,000 × g for 10 minutes at 4°C.</li> </ol>				
	<ol> <li>Check for lysis efficiency: To check for cell lysis, take a 10 μL sample, add 1 μL of Trypan Blue and load onto a hemacytometer. See Growth and Maintenance of Insect Cell Lines manual for protocols. All cells stain blue when lysis is complete.</li> </ol>				
	10. Transfer supernatant (lysate) to a new tube. Keep on ice. Proceed to <b>Detecting Recombinant Protein</b> , below.				
	<b>Note:</b> We recommend that you save the pellet from the lysate (the insoluble portion), because it may be useful for analysis if you cannot detect proteins in the lysate or the supernatant.				
Detecting Recombinant Protein	You may use any method of choice to detect your recombinant protein of interest including functional analysis or western blot. If you perform western blot analysis, you will need to have an antibody to your protein of interest.				
	The pFastBac <sup>™</sup> /HBM-TOPO <sup>®</sup> vector allows the expression of your recombinant protein of interest as a C-terminal 6×His fusion. You can use the antibodies listed on page 55 to detect your recombinant protein.				
Note	If you are using polyacrylamide gel electrophoresis to detect your recombinant protein, you should note that the presence of the C-terminal 6×His tag, and the Tobacco Etch Virus (TEV) recognition site will increase the size of your protein by at least 3 kDa.				
	The HBM secretion signal coding sequence is cleaved upon secretion, and thus does not change the size of your protein.				

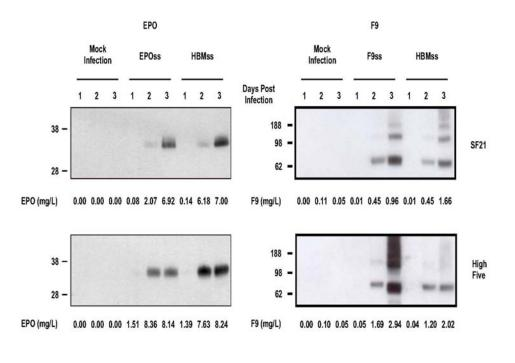
## Analyzing Recombinant Protein, Continued

Glycosylation	When expressing and purifying a glycosylated protein in a heterologous expression system, quickly determine whether the protein is glycosylated properly. Refer to published protocols for carbohydrate analysis of proteins to characterize glycosylated proteins of interest (Ausubel <i>et al.</i> , 1994). Further information about glycosylation in eukaryotes is also available in published literature (Varki & Freeze, 1994).
Purifying Recombinant Protein	The presence of the C-terminal 6×His tag in the pFastBac <sup>™</sup> /HBM-TOPO <sup>®</sup> vector allows the purification of your recombinant protein with a metal-chelating resin such as ProBond <sup>™</sup> and Ni-NTA available from Invitrogen (see page 55 for ordering information). Refer to the manual included with each product for guidelines to purify your fusion protein. These manuals are available for downloading at www.invitrogen.com or by contacting Technical Support (page 56).
Removing the 6×His Tag Using AcTEV <sup>™</sup> Protease	The pFastBac <sup>™</sup> /HBM-TOPO <sup>®</sup> vector contains a Tobacco Etch Virus (TEV) recognition site that allows the removal of the 6×His tag from your recombinant fusion protein using AcTEV <sup>™</sup> Protease available separately from Invitrogen (see page 53 for ordering information). Instructions for digestion are included with the product. For more information, refer to www.invitrogen.com or contact Technical Support (see page 56).
Assay for β-glucuronidase	<ul> <li>If you include the pFastBac<sup>™</sup> 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 development of blue color.</li> <li>Mix 5 µL of 20 mg/mL X-glucuronide solution (in DMSO or dimethylformamide) with 50 µL of cell-free medium.</li> <li>Monitor for development of blue color within 2 hours.</li> <li>Note: Other methods are also suitable.</li> </ul>
	Tore, o aler metrious are also butable.

#### **Expected Results**

#### Introduction

In the following experiments (see figures below), the relative secretion efficiency of proteins fused to the HBM signal sequence was compared to two ordinarily secreted proteins, human coagulation factor IX (F9) and erythropoietin (EPO). To asses the relative secretion efficiency, SF21 and High Five<sup>™</sup> cells were infected with recombinant viruses coding for EPO and F9 fused to their own or to HBM signal sequences as indicated. The culture media were collected at the indicated time points. Proteins were detected by Western blots using anti-his antibody. Protein yield (mg/L) was estimated by enzyme-linked immunosorbent assays using VisuLize<sup>™</sup> Factor IX Antigen Kit (Affinity Biologicals Inc, Ancaster, Ontario, Canada) and Quantikine<sup>®</sup> IVD<sup>®</sup> (R&D Systems, Minneapolis, MN).



# **Optimizing Protein Expression**

Introduction	A number of factors can influence determination of optimal expression conditions including the cell line, MOI, your application of interest, and the nature of your gene of interest. Use the following guidelines to determine the optimal conditions for expressing your recombinant protein of interest.
Time Course	Infect cells at a constant MOI and assay for recombinant protein expression at different time points post-infection (e.g., 24, 48, 72, 96 hours). We recommend that you assay for protein expression at 24-hour intervals initially, to get a general idea of when the protein is being expressed.
	Once you have determined a time frame where optimal protein expression occurs (e.g., between 48 and 72 hours), perform a second time course with selected intermediate time points (e.g., 52, 60, and 68 hours) to further optimize your expression levels.
Synchronous Infection	Synchronous infection is defined as the infection of all cells in a culture at the same time point. Therefore, a true time = zero is established. Achieving synchronous infection results in the maximum amount of protein being harvested at a given time point post-infection, because all cells in the culture are expressing protein at the same time. You must determine the maximum time point for each protein and for each cell line used. You may test different MOIs after the initial time course to achieve synchronous infection.
Multiplicity of Infection (MOI)	Infect a population of cells at varying MOIs ( <i>e.g.</i> , 1, 2, 5, 10, 20) and assay for protein expression. Use the MOI that provides the optimal level of recombinant protein expression.
	For example, if an MOI of 5 gives you protein over a wide range of times, but an MOI of 10 lyses all infected cells before sufficient protein can accumulate, try an MOI of 6 and/or an MOI of 8. The objective in trying different MOIs is to find the MOI that yields the highest protein levels and the least loss due to lysis and proteolysis.
Using Suspension Culture vs. Adherent Culture	The use of suspension culture (spinner or shake flask) versus adherent culture can increase the cell density per mL of culture, and therefore can potentially increase the relative yield of protein per mL of culture.

## Large-Scale Expression of Recombinant Protein

After successfully optimizing expression and secretion levels, you may proceed to large-scale expression of your recombinant protein. You may move up to larger vessels (1 liter or more) or go larger still and use airlift bioreactors and/or fermenters. This section summarizes the requirements and options that exist for large-scale secreted expression of protein using the baculovirus expression system.	
If you are scaling up your suspension culture up to 1 liter spinner flasks, see the sections on Suspension Cell Culture in the <b>Growth and Maintenance of Insect</b> <b>Cell Lines</b> manual. This manual provides information on how to adapt Sf9, Sf21, or High Five <sup>™</sup> insect cell lines to suspension culture, and protocols for maintaining and scaling up suspension cultures. It is available for downloading at www.invitrogen.com or by contacting Technical Support (see page 56). To scale up your culture to 1 liter spinner flasks (500 mL total culture volume), we	
recommend the following:	
1. Generate a large-scale, high-titer stock of the desired recombinant virus (see page 23). This stock will allow you to infect many large-scale cultures and ensure consistency in protein expression.	
2. Start with 100 or 250 mL spinner flasks (50–125 mL of insect cell culture) and scale up to 1 liter spinners with 500 mL of cell suspension.	
3. Seed cultures at $1 \times 10^6$ cells/mL and subculture cell suspension when the density reaches $2.0 \times 10^6$ - $2.5 \times 10^6$ cells/mL.	
4. Check cell viability daily to ensure the culture is >95% viable.	
5. Add Pluronic <sup>®</sup> F-68 to a final concentration of 0.1% in your spinner culture. Adding Pluronic <sup>®</sup> F-68 protects the cells from shearing forces, and it allows you to increase the impellar speed to 120 rpm for larger cultures. Increasing the impeller speed increases aeration of the culture for better growth.	

### Large-Scale Expression of Recombinant Protein, Continued

Large-ScaleThe following table summarizes other methods, requirements, benefits and<br/>references for scale-up production of recombinant protein using the baculovirus<br/>expression system.

Method	Requirements	Benefits	References
Stirred Bioreactor	<ul> <li>For a 5-liter Bioreactor:</li> <li>5-liter bioreactor.</li> <li>Sterilized tubing.</li> <li>Microbial air filters.</li> <li>High purity nitrogen, oxygen, and air.</li> <li>pH, dissolved oxygen and temperature probes.</li> <li>External dissolved oxygen controller.</li> <li>External pH controller.</li> <li>Peristaltic pump for acid/base lines.</li> <li>Linear recorder to monitor dissolved oxygen and pH control.</li> <li>Laminar flow hood in close proximity to the bioreactor.</li> </ul>	<ul> <li>Addresses increased oxygen needs of large-scale culture.</li> <li>Controlled growth and optimization of variables in the culture.</li> <li>Increased cell densities.</li> <li>Elevated protein production.</li> <li>Reproducible results for batch production of protein.</li> </ul>	(Tom <i>et al.</i> , 1995) (Murhammer & Goochee, 1988) (Maiorella <i>et al.</i> , 1988) (O'Reilly <i>et al.</i> , 1992)
Airlift Fermentor	<ul> <li>For a 5-liter Airlift Fermentor:</li> <li>5-liter airlift fermentor system.</li> <li>Dissolved oxygen control module.</li> <li>Dissolved oxygen electrode.</li> <li>Microbial air filters.</li> <li>1/4" stainless-steel tubing.</li> <li>Silicone tubing.</li> <li>Circulating water bath.</li> </ul>	<ul> <li>Addresses increased oxygen needs of large-scale culture.</li> <li>Control over environmental variables in the culture.</li> <li>Increased cell densities.</li> <li>Elevated protein production.</li> </ul>	(Maiorella <i>et al.,</i> 1988) (Murhammer & Goochee, 1988) (Onken & Weiland, 1983) (O'Reilly <i>et al.,</i> 1992)
Insect Larvae	Trichoplusia ni larvae and facilities to grow larvae.	<ul> <li>More accurate posttranslational modification of recombinant protein – does not rely on one cell type only.</li> <li>Higher levels than in cell culture.</li> <li>Inexpensive.</li> </ul>	(Medin <i>et al.,</i> 1990) (Wood <i>et al.,</i> 1993)

### Troubleshooting

Cloning into the pFastBac <sup>™</sup> /HBM- TOPO <sup>®</sup> Vector	For troubleshooting any problems you may encounter when generating your pFastBac <sup>™</sup> /HBM-TOPO <sup>®</sup> construct, refer to the Bac-to-Bac <sup>®</sup> TOPO <sup>®</sup> Cloning Kit manual (part no. A10605) supplied with this kit. The Bac-to-Bac <sup>®</sup> TOPO <sup>®</sup> Cloning Kit manual is also available at www.invitrogen.com or by contacting Technical Support (see page 56).

Generating Recombinant Bacmid DNA The table below lists some potential problems that you may encounter when generating the recombinant bacmid following transformation into DH10Bac<sup>TM</sup> *E. coli*.

Problem	Reason	Solution
No blue colonies (non- recombinant) obtained	Insufficient time for color development.	Wait at least 48 hours before identifying colony phenotypes.
(i.e., all colonies are white) <b>Note:</b> Although you will	Used X-gal instead of Bluo-gal in agar plates.	Use Bluo-gal in selective plates to increase the contrast between blue and white colonies.
pick white colonies, you should expect to see some blue colonies. Blue colonies contain non-	Insufficient growth after transposition.	Grow transformed cells in S.O.C. Medium for a minimum of 4 hours before plating.
recombinant bacmids.	Bluo-gal and IPTG omitted from plates.	Prepare fresh selective plates containing 50 µg/mL kanamycin, 7 µg/mL gentamicin, 10 µg/mL tetracycline, 100 µg/mL Bluo-gal, and 40 µg/mL IPTG.
	Too many colonies on the plate.	• Serially dilute the transformation mixture and plate to give well-separated colonies.
		• Adjust the serial dilutions of cells (10 <sup>-2</sup> to 10 <sup>-4</sup> ) to obtain well-spaced colonies.
	Plates too old or stored in light.	• Do not use plates that are more than 4 weeks old.
		• Store plates protected from light.
	Incubation period too short or temperature too low.	Wait at least 48 hours before picking colonies. Incubate plates at 37°C.

Problem	Reason	Solution	
All colonies are blue	DNA from your pFastBac™/HBM-TOPO®	• Use purified plasmid DNA for transformation.	
	construct used for transformation was of poor quality.	• Check the quality of your plasmid DNA; make sure that the DNA is not degraded.	
	Gentamicin omitted from plates.	Prepare fresh selective plates containing 50 µg/mL kanamycin, 7 µg/mL gentamicin, 10 µg/mL tetracycline, 100 µg/mL Bluo-gal, and 40 µg/mL IPTG.	
Few colonies obtained	Used LB medium for recovery/expression period.	Use S.O.C. Medium for the 4 hours growth time.	
	Recovery/expression time too short.	Increase the recovery time to > 4 hours at 37°C or 6 hours at 30°C.	
Poor blue/white colony	Agar not at correct pH.	Adjust pH of LB agar to 7.0.	
differentiation	Intensity of the blue color too weak.	• Use Bluo-gal, not X-gal.	
		<ul> <li>Increase the concentration of Bluo-gal to 300 μg/mL.</li> </ul>	
		• Use dark and light backgrounds to view plates.	
	Too many or too few colonies on plate.	Adjust the serial dilutions of cells to obtain an optimal number of colonies.	
	Incubation period too short or temperature too low.	• Do not pick colonies until 48 hours after plating.	
		• Incubate plates at 37°C.	
	IPTG concentration not optimal.	Optimize the IPTG concentration. A range of 20–60 $\mu$ g/mL IPTG generally gives optimal color development.	

#### Generating Recombinant Bacmid DNA, continued

Isolating Bacmid	The table below lists some potential problems and possible solutions to help you
DNA	troubleshoot recombinant bacmid DNA isolation.

Problem	Reason	Solution
Bacmid DNA is degraded	DNA stored improperly.	• Store purified bacmid DNA in aliquots at 4°C for no more than 2 weeks.
		• Do not freeze/thaw the bacmid DNA.
		<ul> <li>For long term storage of bacmid DNA, prepare glycerol stocks of DH10Bac<sup>™</sup> E. coli containing the verified bacmid DNA.</li> </ul>
	High molecular weight bacmid DNA handled improperly.	• When isolating bacmid DNA, do not vortex the DNA solution.
		• Do not resuspend DNA pellets mechanically; allow the solution to sit in the tube with occasional gentle tapping of the bottom of the tube.
Poor yield	Used incorrect antibiotic concentrations.	Grow transformed DH10Bac <sup>™</sup> cells in LB medium containing 50 µg/mL kanamycin, 7 µg/mL gentamicin, and 10 µg/mL tetracycline.
Bacmid DNA contains a mixture of recombinant bacmid and empty bacmid	Picked a colony that was gray or dark in the center.	Analyze more white DH10Bac <sup>™</sup> transformants and choose one that contains recombinant bacmid DNA only.

Transfecting	The table below lists some potential problems and possible solutions that may
Insect Cells	help you troubleshoot insect cell transfection.

Problem	Reason	Solution
Low yield of virus	Low transfection efficiency.	Use Invitrogen's Cellfectin <sup>®</sup> II     Reagent for transfection.
		• Perform transfection in unsupplemented Grace's Medium; make sure that no supplements, FBS, or antibiotics are present during transfection.
		<ul> <li>Harvest viral supernatant when signs of infection are visible (i.e., &gt;72 hours post-transfection).</li> </ul>
	Cells plated too sparsely.	Plate insect cells at the recommended cell density.
	Used too much or too little Cellfectin <sup>®</sup> II or other lipid reagent.	Optimize the amount of Cellfectin <sup>®</sup> II or other lipid reagent used.
	Time of incubation with DNA:lipid complexes too short or too long.	Optimize the incubation time (e.g., 3 to 8 hours).
	Recombinant bacmid DNA is degraded.	Check the quality of your recombinant DNA by agarose gel electrophoresis prior to transfection.
		<ul> <li>Prepare bacmid DNA using Invitrogen's PureLink<sup>™</sup> HiPure Miniprep or Maxiprep Kit (see page 53 for ordering information).</li> </ul>
		• Store purified bacmid at 4°C; <b>do</b> <b>not</b> freeze, because freezing the baculovirus decreases transfection efficiency.
	Bacmid DNA is not pure (i.e., contains recombinant bacmid and empty bacmid).	• Screen other DH10Bac <sup>™</sup> transformants and choose one that contains only recombinant bacmid.
		• Perform plaque purification to isolate recombinant baculovirus.

Expressing Your	The table below lists some potential problems and possible solutions that may
Protein	help you troubleshoot your expression experiments.

Problem	Reason	Solution
Low protein expression	Viral stock contains a mixture of recombinant and non- recombinant baculovirus.	Perform plaque purification to isolate recombinant baculovirus.
	Baculovirus not recombinant.	• Verify transposition of bacmid DNA by PCR analysis, using the pUC/M13 Forward and Reverse primers.
		• Re-transfect insect cells with new recombinant bacmid DNA.
	Used too low or too high viral titer.	Optimize infection conditions by varying the MOI.
	Time of cell harvest not optimal.	Perform a time course of expression to determine the optimal time to obtain maximal protein expression.
	Cell growth conditions and medium not optimal.	• Optimize culture conditions based on the size of your culture vessel and expression conditions.
		• Culture High Five <sup>™</sup> cells in Express Five <sup>®</sup> SFM for optimal cell growth and protein expression.
	Cell line not optimal.	Use High Five <sup>™</sup> Cells for highest level of secreted expression.
	Protein expression is not optimal.	Optimize protein expression by varying such parameters as incubation temperature and oxygenation.

## Appendix

Recipes			
Antibiotic Stor Solutions	premixed s	can be ordered in either dry powdered solution. Store these solutions according dations. For the antibiotics below, prep s directed:	g to the manufacturer's
	Antibiotic	Stock Solution Concentration	Storage
	Ampicillin	50 mg/mL in water; filter-sterilize	–20°C, protected from light
	Kanamycin	10 mg/mL in water; filter-sterilize	–20°C, protected from light
	Tetracycline*	10 mg/mL in 100% ethanol; filter- sterilize	–20°C, protected from light
	Gentamicin	7 mg/mL in water; filter-sterilize	–20°C, protected from light
IPTG	<ol> <li>Dissolv</li> <li>Adjust</li> <li>Filter-s</li> <li>Disper</li> <li>Store a</li> </ol>	procedure below to prepare a 200 mg/ ve 2 g of IPTG in 8 mL of sterile water. the volume of the solution to 10 mL w sterilize through a 0.22 micron filter. use the stock solution into 1 mL aliquots t $-20^{\circ}$ C.	ith sterile water. s.
Bluo-gal	<ol> <li>Dissolv to mak</li> <li>Import Disper</li> <li>Do not</li> </ol>	guidelines below to prepare a 20 mg/n we the Bluo-gal in dimethylformamide the a 20 mg/mL stock solution. Use a gla tant: Exercise caution when working w use solutions in a vented chemical hood filter the stock solution. t –20°C protected from light.	or dimethylsulfoxide (DMSO) ass or polypropylene tube. ith dimethylformamide.
			Continued on next page

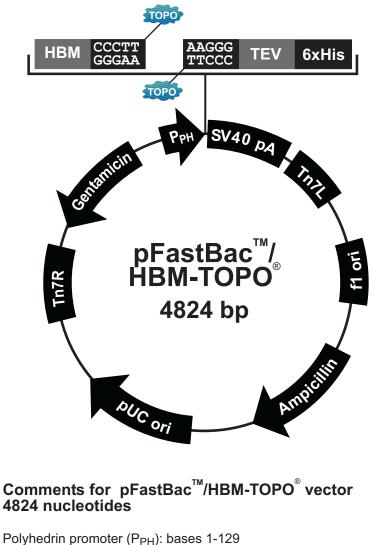
### Recipes, Continued

LB (Luria-Bertani) Medium	<ul> <li>Composition:</li> <li>1.0% Tryptone (casein peptone)</li> <li>0.5% Yeast Extract</li> <li>1.0% NaCl</li> <li>pH 7.0</li> <li>1. For 1 liter, dissolve 10 g tryptone, 5 g yeast extract, and 10 g NaCl in 950 mL deionized water.</li> <li>2. Adjust the pH of the solution to 7.0 with NaOH and bring the volume up to 1 liter.</li> <li>3. Autoclave on liquid cycle for 20 minutes. Allow solution to cool to ~55°C and add antibiotic if needed.</li> <li>4. Store at room temperature or at 4°C.</li> </ul>
LB (Luria-Bertani) Plates	<ul> <li>Follow the procedure below to prepare LB agar plates.</li> <li>Prepare LB medium as above, but add 15 g/L agar before autoclaving.</li> <li>Autoclave on liquid cycle for 20 minutes.</li> <li>After autoclaving, cool to ~55°C, add antibiotic(s) and pour into 10 cm plates.</li> <li>Let harden, then invert and store at 4°C, in the dark. Plates containing antibiotics are stable for up to 4 weeks.</li> <li>LB agar selective plates for DH10Bac<sup>™</sup> transformation</li> <li>Follow Steps 1–2 in the procedure above.</li> <li>After autoclaving, cool to ~55°C, and add the following: <ul> <li>50 µg/mL kanamycin</li> <li>7 µg/mL gentamicin</li> <li>10 µg/mL tetracycline</li> <li>100 µg/mL Bluo-gal</li> <li>40 µg/mL IPTG</li> </ul> </li> <li>Let harden, then invert and store at 4°C, in the dark. Tetracycline and Bluo-gal are light sensitive, so make sure that plates are stored protected from light.</li> </ul>

## Map of pFastBac<sup>™</sup>/HBM-TOPO<sup>®</sup>

Description

The map below shows the elements of pFastBac<sup>™</sup>/HBM-TOPO<sup>®</sup> vector. **The vector sequence is available for downloading from www.invitrogen.com or by contacting Technical Support (page 56).** 



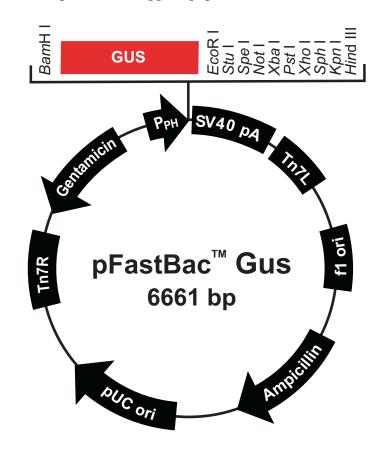
Honey Bee Mellitin (HBM) secretion signal: 141-210 TOPO cloning site: bases 215-216 TEV recognition site: bases 222-242 6xHis tag: bases 243-260 SV40 polyadenylation signal: bases 305-545 Tn7L: bases 574-739 f1 origin: bases 923-1377 Ampicillin resistance gene: bases 1508-2368 pUC origin: bases 2513-3186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

### Map of pFastBac<sup>™</sup> Gus Control Plasmid

#### Description

pFastBac<sup>T</sup> Gus is a 6,661 bp control vector that contains the *Arabidopsis thaliana* gene for  $\beta$ -glucuronidase (Gus) (Kertbundit *et al.*, 1991). The molecular weight of  $\beta$ -glucuronidase is 68.5 kDa.

The map below shows the elements of pFastBac<sup>™</sup> Gus control plasmid. **The vector sequence is available for downloading from www.invitrogen.com or by contacting Technical Support (page 56).** 



# Comments for pFastBac<sup>™</sup> 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<sub>PH</sub>): bases 3904-4032 GUS ORF: bases 4081-5892 SV40 polyadenylation signal: bases 6047-6287 Tn7L: bases 6315-6480

#### **Accessory Products**

#### Additional Products

All of the reagents supplied in the Bac-to-Bac® HBM TOPO® Secreted Expression System and other products suitable for use with the Bac-to-Bac® HBM TOPO® Secreted Expression System are available separately from Invitrogen. Ordering information for these reagents is provided below. For more information, refer to our website at www.invitrogen.com or contact Technical Support (see page 56).

Item	Quantity	Cat. no.
Bac-to-Bac® HBM TOPO® Cloning Kit	1 kit	A11338
MAX Efficiency <sup>®</sup> DH10Bac <sup>™</sup> Competent <i>E. coli</i>	$5 \times 100 \ \mu L$	10361-012
One Shot <sup>®</sup> Mach1 <sup><math>m</math></sup> -T1 <sup><math>R</math></sup> Chemically Competent <i>E. coli</i>	$21 \times 50 \ \mu L$	C8620-03
Cellfectin <sup>®</sup> II Reagent	1 mL	10362-100
Platinum <sup>®</sup> Pfx DNA Polymerase	100 units	11708-013
AccuPrime <sup>™</sup> <i>Pfx</i> DNA Polymerase	200 reactions	12344-024
<i>Pfx50</i> <sup>™</sup> DNA Polymerase	100 reactions	12355-012
Platinum <sup>®</sup> Taq DNA Polymerase High Fidelity	100 reactions	11304-011
PureLink <sup>™</sup> PCR Purification Kit	50 preps	K3100-01
PureLink <sup>™</sup> Quick Gel Extraction System	1 kit	K2100-12
PureLink™ HiPure Plasmid Miniprep Kit	25 preps 100 preps	K2100-02 K2100-03
PureLink™ HiPure Plasmid Midiprep Kit	25 preps 50 preps	K2100-04 K2100-05
PureLink <sup>™</sup> HiPure Plasmid Maxiprep Kit	10 preps 25 preps	K2100-06 K2100-07
Ampicillin Sodium Salt, irradiated	200 mg	11593-027
Kanamycin Sulfate	5 g 25 g	11815-024 11815-032
Kanamycin Sulfate (100X), liquid	100 mL	15160-054
Gentamicin Reagent Solution, liquid (50 mg/mL)	10 mL 10 × 10 mL	15750-060 15750-078
Bluo-gal	1 g	15519-028
Isopropylthio-β-galactoside (IPTG)	1 g	15529-019
S.O.C. Medium	10 × 10 mL	15544-034
(Miller's LB Broth Base) <sup>®</sup> Luria Broth Base, powder	500 g	12795-027
Water, distilled (cell-culture grade)	500 mL	15230-162
4% Agarose gel (optimal for insect cell growth)	40 mL	18300-012
Fetal Bovine Serum (FBS), Qualified, Heat Activated	100 mL	16140-063

#### Insect Cell Culture Products

A variety of insect cell lines and GIBCO<sup>®</sup> cell culture products are available from Invitrogen to facilitate baculovirus-mediated expression of your recombinant protein in insect cells. For more information about the insect cell lines and GIBCO<sup>™</sup> cell culture products, refer to www.invitrogen.com or contact Technical Support (see page 56).

Quantity	Cat. no.
$3 \times 10^{6}$ cells	B855-02
$1.5 \times 10^7$ cells	11496-015
$1.5 \times 10^7$ cells	11497-013
$1 \times 10^7$ cells	12552-014
500 mL	10902-096
500 mL	12658-019
100 mL	10967-032
1 liter	10486-025
500 mL	11595-030
500 mL	11605-094
100 mL	11667-037
100 mL	15070-063
100 mL	24040-032
	$\begin{array}{c c} 3 \times 10^{6} \text{ cells} \\ 1.5 \times 10^{7} \text{ cells} \\ 1.5 \times 10^{7} \text{ cells} \\ 1.5 \times 10^{7} \text{ cells} \\ 1 \times 10^{7} \text{ cells} \\ 500 \text{ mL} \\ 100 \text{ mL} \\ 10 \text{ m} \\ 100 \text{ m} \\ 10 $

Note: Reagents are also available in other sizes.

 $\ensuremath{\mathsf{PLURONIC}}\xspace^{\ensuremath{\mathbb{R}}}$  is a registered trademark of BASF Corporation.

#### Accessory Products, Continued

#### Detecting Recombinant Fusion Protein

If you have cloned your gene of interest in frame with C-terminal polyhistidine tag of the pFastBac<sup>™</sup>/HBM-TOPO<sup>®</sup> vector, you may detect expression of your recombinant fusion protein using the following antibodies. The amount of antibody supplied is sufficient for 25 western blots.

Product	Epitope	Cat. no.
Anti-His (C-term) Antibody	Detects the C-terminal	R930-25
Anti-His (C-term)-HRP Antibody	polyhistidine (6×His) tag:	R931-25
Anti-His (C-term)-AP Antibody	HHHHHH-COOH (requires the free carboxyl group for detection (Lindner <i>et al.</i> , 1997)	R932-25
Penta-His <sup>™</sup> mouse IgG₁ monoclonal Antibody	Detects both N- and C-terminal polyhistidine (6×His) tag	P21315

#### Purifying Recombinant Fusion Proteins

If you express your gene of interest as a fusion with the polyhistidine tag from the pFastBac<sup>™</sup>/HBM-TOPO<sup>®</sup> vector, you may use ProBond<sup>™</sup> or Ni-NTA resins to purify your recombinant fusion protein. See the table below for ordering information.

Item	Quantity	Cat. no.
ProBond <sup>™</sup> Nickel-chelating Resin	50 mL 150 mL	R801-01 R801-15
ProBond <sup>™</sup> Purification System	6 purifications	K850-01
Ni-NTA Agarose	10 mL 25 mL 100 mL	R901-01 R901-15 R901-10
Ni-NTA Purification System	6 purifications	K950-01
Purification Columns (10 mL polypropylene columns)	50 columns	R640-50
AcTEV <sup>™</sup> Protease	1,000 Units 10,000 Units	12575-015 12575-023

# **Technical Support**

Web Resources	Visit the Invitrogen website at <b>www.invitrogen.com</b> for:			
	<ul> <li>Technical resources, including manuals, vector maps and sequences, application notes, MSDSs, FAQs, formulations, citations, handbooks, etc.</li> </ul>			
	Complete	technical support contact informa	tion	
T	Access to	the Invitrogen Online Catalog		
	Additional product information and special offers			
Contact Us	For more information or technical assistance, call, write, fax, or email. Additional international offices are listed on our website (www.invitrogen.com).			
<b>Corporate Headquarters:</b> 5791 Van Allen Way Carlsbad, CA 92008 USA Tel: 1 760 603 7200 Tel (Toll Free): 1 800 955 6288 Fax: 1 760 602 6500 E-mail: techsupport@invitrogen.com		Japanese Headquarters: LOOP-X Bldg. 6F 3-9-15, Kaigan Minato-ku, Tokyo 108-0022 Tel: 81 3 5730 6509 Fax: 81 3 5730 6519 E-mail: jpinfo@invitrogen.com	<b>European Headquarters:</b> Inchinnan Business Park 3 Fountain Drive Paisley PA4 9RF, UK Tel: 44 (0) 141 814 6100 Tech Fax: 44 (0) 141 814 6117 E-mail: eurotech@invitrogen.com	
MSDS	Material Safety Data Sheets (MSDSs) are available on our website at www.invitrogen.com/msds.			
Certificate of Analysis	The Certificate of Analysis (CofA) provides detailed quality control information for each product and is searchable by product lot number, which is printed on each box. CofAs are .available on our website at www.invitrogen.com/support.			
Limited Warranty	<ul> <li>Invitrogen (a part of 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 an Invitrogen product or service, contact our Technical Support Representatives.</li> <li>All Invitrogen 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. This warranty limits the Company's liability to only the price of the product. No warranty is granted for products beyond their listed expiration date. No warranty is applicable unless all product components are stored in accordance with instructions. 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.</li> <li>Invitrogen 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, please report it to our Technical Support Representatives.</li> <li>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.</li> </ul>			

### **Purchaser Notification**

Introduction	Use of the Bac-to-Bac <sup>®</sup> HBM TOPO <sup>®</sup> Expression System and the pFastBac <sup>™</sup> /HBM-TOPO <sup>®</sup> vector is covered under the licenses detailed below.
Limited Use Label License No. 5: Invitrogen Technology	The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. The buyer may transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for any Commercial Purpose, and that such collaborator agrees in writing (a) not to transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes. Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. For products that are subject to multiple limited use label licenses, the terms of the most restrictive limited use label license shall control. Life Technologies Corporation will not assert a claim against the buyer of infringement of patents owned or controlled by Life Technologies Corporation which cover this product to ased upon the manufacture, use or sale of a therapeutic, clinical diagnostic, vaccine or prophylactic product developed in research by the buyer in which this product or its components was used in the manufacture of such product. If the purchaser is not willing to accept the limitations of this limited use statement, Life Techno
Limited Use Label License No. 21: Bac-to-Bac <sup>®</sup> and Bac-to-Bac <sup>®</sup> HT	This product is the subject of U.S. Patent No. 5,348,886. This product is sold under patent license from Monsanto for research purposes only and no license for commercial use is included. Requests for licenses for commercial manufacture or use should be directed to Director, Monsanto Corporate Research, 800 N. Lindbergh, St. Louis, Missouri 63167.
Limited Use Label License No. 22: Vectors and Clones Encoding Histidine Hexamer	This product is licensed under U.S. Patent Nos. 5,284,933 and 5,310,663 and foreign equivalents from Hoffmann-LaRoche, Inc., Nutley, NJ and/or Hoffmann-LaRoche Ltd., Basel, Switzerland and is provided only for use in research. Information about licenses for commercial use is available from QIAGEN GmbH, Max-Volmer-Str. 4, D-40724 Hilden, Germany.

## Purchaser Notification, Continued

Information for European Customers Using Mach1 <sup>™</sup> -T1 <sup>R</sup> Cells	The Mach1 <sup>™</sup> -T1 <sup>ℝ</sup> <i>E. coli</i> strain is genetically modified to carry the <i>lac</i> ZΔM15 <i>hsd</i> R <i>lac</i> X74 <i>rec</i> A <i>end</i> A <i>ton</i> 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.
Information for All Non-U.S. Customers Using Mach1 <sup>™</sup> -T1 <sup>R</sup> Cells	The parental strain of Mach1 <sup>™</sup> -T1 <sup>ℝ</sup> <i>E. coli</i> 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.
Information for European Customers Using DH10Bac <sup>™</sup> cells	The DH10Bac <sup><math>m</math></sup> strain is genetically modified and carries the pBR322-derived plasmid, pMON7124 ( <i>bom</i> <sup>+</sup> , <i>tra</i> <sup>-</sup> , <i>mob</i> <sup>-</sup> ). 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.

- 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*, Greene Publishing Associates and Wiley-Interscience, New York
- Barry, G. F. (1988) A Broad Host-Range Shuttle System for Gene Insertion into the Chromosomes of Gram-negative Bacteria. Gene 71, 75-84
- Ciccarone, V. C., Polayes, D., and Luckow, V. A. (1997) *Generation of Recombinant Baculovirus DNA in E. coli Using Baculovirus Shuttle Vector*. Methods in Molecular Medicine (Reischt, U., Ed.), 13, Humana Press Inc., Totowa, NJ
- Davis, T. R., Trotter, K. M., Granados, R. R., and Wood, H. A. (1992) Baculovirus Expression of Alkaline Phosphatase as a Reporter Gene for Evaluation of Production, Glycosylation, and Secretion. Bio/Technology 10, 1148-1150
- Davis, T. R., Wickham, T. J., McKenna, K. A., Granados, R. R., Shuler, M. L., and Wood, H. A. (1993) Comparative Recombinant Protein Production of Eight Insect Cell Lines. In Vitro Cell. Dev. Biol. 29A, 388-390
- Kertbundit, S., Greve, H. d., Deboeck, F., Montagu, M. V., and Hernalsteens, J. P. (1991) *In vivo* Random bglucuronidase Gene Fusions in *Arabidopsis thaliana*. Proc. Natl. Acad. Sci. USA *88*, 5212-5216
- King, L. A., and Possee, R. D. (1992) *The Baculovirus Expression System: A Laboratory Guide*, Chapman and Hall, New York, NY
- Lindner, P., Bauer, K., Krebber, A., Nieba, L., Kremmer, E., Krebber, C., Honegger, A., Klinger, B., Mocikat, R., and Pluckthun, A. (1997) Specific Detection of His-tagged Proteins With Recombinant Anti-His Tag scFv-Phosphatase or scFv-Phage Fusions. BioTechniques 22, 140-149
- Luckow, V. A. (1991) in *Recombinant DNA Technology and Applications* (Prokop, A., Bajpai, R. K., and Ho, C., eds), McGraw-Hill, New York
- Luckow, V. A., Lee, C. S., Barry, G. F., and Olins, P. O. (1993) Efficient Generation of Infectious Recombinant Baculoviruses by Site-Specific Transposon-Mediated Insertion of Foreign Genes into a Baculovirus Genome Propagated in *Escherichia coli*. J. Virol. 67, 4566-4579
- Luckow, V. A., and Summers, M. D. (1988) Signals Important for High-Level Expression of Foreign Genes in *Autographa californica* Nuclear Polyhedrosis Virus Expression Vectors. Virology 167, 56-71
- Maiorella, B., Inlow, D., Shauger, A., and Harano, D. (1988) Large-Scale Insect Cell Culture for Recombinant Protein Production. Bio/Technology 6, 1406-1410
- Medin, S. A., Hunt, L., Gathy, K., Evans, R. K., and Coleman, M. S. (1990) Efficient, Low-cost Protein Factories: Expression of Human Adenosine Deaminase in Baculovirus-infected Insect Larvae. Proc. Natl. Acad. Sci. USA 87, 2760-2764
- Murhammer, D., and Goochee, C. F. (1988) Scale-up of Insect Cell Cutures: Protective Effects of Pluronic F-68. Bio/Technology *6*, 1411-1415

- O'Reilly, D. R., Miller, L. K., and Luckow, V. A. (1992) *Baculovirus Expression Vectors: A Laboratory Manual*, W. H. Freeman and Company, New York, N. Y.
- Onken, U., and Weiland, P. (1983) Airlift Fermentors: Construction, Behaviour and Uses. Biotechnol. Proc. 1, 67-95
- Tom, R. L., Caron, A. W., Massie, B., and Kamen, A. (1995) Scale-up of Recombinant Virus and Protein Production in Stirred-tank Reactors. Methods in Molecular Biology (Richardson, C. D., Ed.), 39, Humana Press, Totowa, NJ
- Varki, A., and Freeze, H. H. (1994) The Major Glycosylation Pathways of Mammalian Membranes. A Summary. Subcell. Biochem. 22, 71-100
- Wood, H. A., Trotter, K. M., Davis, T. R., and R., H. P. (1993) Per Os Infectivity of Preoccluded Virions from Polyhedrin-Minus Recombinant Baculoviruses. J. Invert. Path. 62, 64-67

©2009 Life Technologies Corporation. All rights reserved. For research use only. Not intended for any animal or human therapeutic or diagnostic use.

# invitrogen®

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

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