



---

## RTS™ Linear Template Fab Kit Manual

---

For the generation of linear expression templates by PCR  
and easy cloning of PCR products into the pIX3.0 Vector

RTS Linear Template Kit Plus

RTS pIX3.0 Vector

RTS Linear Template Fab Kit, August, 2025

© 2025 biotechrabbit, all rights reserved.

This document and the product it describes are subject to change without prior notice. This document does not represent a commitment on the part of biotechrabbit GmbH or its distributors.

Trademarks: Coomassie® (ICI [Imperial Chemical Industries] Organics Inc.); Eppendorf®, Eppendorf-Netheler-Hinz GmbH; ProteoMaster™, RTS™, ExactiFi™, GenUP™, biotechrabbit GmbH.

For Research Purposes Only. Proteins expressed using the RTS, and data derived therefrom that would enable the expression of such proteins (collectively, "Expressed Proteins"), may be used only for the internal research of the purchaser of this system. Expressed Proteins may not be sold or transferred to any third party without the written consent of biotechrabbit GmbH.

The purchase price of this product includes a limited, non-exclusive, non-transferable license under U.S. patents 6.168.931 and 6.337.191 and their foreign counterparts, exclusively licensed by a member of the biotechrabbit GmbH.

The continuous-exchange cell-free (CECF) technology applied in the RTS 100 Wheat Germ CECF, RTS 500 Wheat Germ CECF, RTS 100 E. coli Disulfide, RTS 500 E. coli Disulfide, RTS 500 ProteoMaster E. coli HY and RTS 9000 E. coli HY products is based on patented technology (U.S. Patent 5,478,730). The purchase price of this product includes practicing a cell-free expression achieving continuous production of a polypeptide in the presence of a semi-permeable barrier and related processes described in said patents.

# Contents

<b>Contents</b> .....	<b>3</b>
<b>Materials supplied</b> .....	<b>4</b>
<b>Shipping and storage conditions</b> .....	<b>4</b>
<b>Product limitations</b> .....	<b>5</b>
<b>Technical Assistance</b> .....	<b>5</b>
<b>Quality assurance</b> .....	<b>5</b>
<b>Safety Information</b> .....	<b>6</b>
<b>Product warranty</b> .....	<b>6</b>
<b>Introduction</b> .....	<b>7</b>
The RTS System	8
RTS Platform - Complementary Prokaryotic and Eukaryotic Systems	8
Identification of optimal constructs in a single day	12
<b>Principle and Procedure</b> .....	<b>13</b>
Generating PCR Products for Use in <i>In Vitro</i> Translation Reactions	14
<b>Protocol: Two-Step PCR Procedure for Generating an Expression Template</b> .....	<b>18</b>
Important points before starting	18
Procedure: First PCR using gene-specific primers	19
Procedure: Second PCR using adapter primers	21
<b>Troubleshooting Guide</b> .....	<b>23</b>
<b>References</b> .....	<b>24</b>
<b>Appendix: Adapter Primer Sequences and Multiple Cloning Sites</b> .....	<b>25</b>
Cloning PCR Products Generated with the RTS Linear Template Fab Kit into the RTS pIX3.0 Vector	26
<b>Ordering information</b> .....	<b>28</b>

## Materials supplied

---

### RTS Linear Template Fab Kit

<b>Ordering number</b>	<b>BR1402201</b>
<b>Number of reactions</b>	<b>96 reactions</b>
No tag Sense Primer (yellow screw-cap)	200 µl
No tag Sense Primer Signal Peptide (blue screw-cap)	200 µl
No tag Antisense Primer (colorless screw-cap)	200 µl
His6 tag Antisense Primer (brown screw-cap)	200 µl
XE-Solution (green screw-cap)	400 µl

---

### RTS pIX3.0 Vector

<b>Ordering number</b>	<b>BR1402701</b>
pIX3.0 Vector (white screw-cap)	25 µg (0.5 µg/µl)

---

## Shipping and storage conditions

The RTS Linear Template Fab Kit is shipped on dry ice and should be stored immediately upon receipt at  $-20^{\circ}\text{C}$  in a constant-temperature freezer upon arrival.

When stored under the above conditions and handled correctly, all kits are stable until the expiration date printed on the label. Avoid repeated freezing and thawing.

---

## Product limitations

RTS Linear Template Fab Kit and RTS 100 E. coli HY Kits and RTS 100 insect cell-based kits are developed, designed and sold for research purposes only. They are not to be used for human diagnostic or drug purposes or to be administered to humans unless expressly cleared for that purpose by the Food and Drug Administration in the USA or the appropriate regulatory authorities in the country of use. All due care and attention should be exercised in the handling of many of the products and materials described in this text.

## Technical Assistance

Our Technical Service is staffed by experienced scientists with extensive practical and theoretical expertise in molecular biology. If you have any questions or experience any difficulties regarding products for cell-free protein synthesis in general, please do not hesitate to contact us.

We encourage you to contact us if you have any suggestions about product performance or new applications and techniques.

For technical assistance and more information please contact biotechrabbit Technical Service: [support@biotechrabbit.com](mailto:support@biotechrabbit.com)

## Quality assurance

biotechrabbit products are manufactured using quality chemicals and materials that meet our high standards. All product components are subjected to rigorous quality assurance testing process:

- Component testing: each component is tested to ensure the composition and quality meet stated specifications.
- Performance testing: each product is tested to ensure it meets the stated performance specification.

Additional quality information is available from [www.biotechrabbit.com](http://www.biotechrabbit.com). Certificate of analysis sheets for biotechrabbit products can be obtained on request.

---

## Safety Information

All due care and attention should be exercised in the handling of the products. We recommend all users of *E. coli* cell-free protein synthesis kit products to adhere to the NIH guidelines that have been developed for recombinant DNA experiments, or to other applicable guidelines. Specifically, always wear a suitable lab coat, disposable gloves, and protective goggles when working with chemicals.

## Product warranty

biotechrabbit is committed to providing products that improve the speed, ease-of-use and quality of enabling technologies.

biotechrabbit guarantees the performance of all products in the manner described in our product literature. The purchaser must determine the suitability of the product for its particular use.

This warranty is in place of any other warranty or guarantee, expressed or implied, instituted by law or otherwise. biotechrabbit provides no other warranties of any kind, expressed or implied, including warranties of merchantability and fitness for a particular purpose. Under no circumstance shall biotechrabbit be responsible for any direct, indirect, consequential or incidental damages or loss arising from the use, misuse, results of use or inability to use its products, even if the possibility of such loss, damage or expense was known by biotechrabbit.

---

## Introduction

Proteins such as enzymes, antibodies, hormones, and structural elements play essential roles in nearly all biological processes. Therefore, great efforts have been made to develop technologies for the production of proteins using recombinant technology. Modern protein engineering methods, which include cloning of DNA sequences and the *in vivo* expression of genes, allow production of specific proteins in large amounts and also production of proteins with improved or altered biological activities.

Several factors must be carefully considered when producing recombinant proteins using *in vivo* expression methods. Cells must be transformed with an expression construct (e.g., plasmid DNA), and transformants containing the correct construct must be selected and cultivated. Overexpression of proteins that are toxic to the host cells can be difficult. Cell lysis and procedures used for purification of protein from whole cell lysates can be complicated; problems may arise because of aggregation or degradation of proteins within the cell.

In most cases these limitations can be overcome by the use of cell-free protein biosynthesis systems (also termed *in vitro* translation [IVT] systems), which are often seen as a very attractive alternative to classical *in vivo* expression systems. Cell-free expression generates proteins by coupled or successive transcription and translation in cell-free extracts of prokaryotic or eukaryotic cells. The advantages of cell-free expression systems include time savings, the possibility to produce proteins that are toxic or have modified or isotope-labeled amino acids, a high protein yield per unit volume, and the ability to adapt reaction conditions to the requirements of the synthesized protein (for example, the inclusion of cofactors). Compared with current cloning techniques, another important advantage offered by a cell-free expression system is the possibility of using PCR products as templates for protein synthesis. This greatly accelerates the protein production process, because no cloning steps are required. Moreover, there is no need for any specialized equipment: only an incubator, pipette, and reaction tubes are needed.

Proteins produced by cell-free expression can be used for the same wide variety of downstream applications as *in-vivo* produced proteins, including activity assays, structural and functional analyses, protein–protein interaction studies, and the expression and analysis of open reading frames.

## The RTS System

biotechrabbit's Rapid Translation System (RTS) is a scalable *in vitro* protein expression platform that produces large amounts of protein for characterization studies, functional assays, or structural analysis. Specific applications can easily be adapted with a truly "open" system based on bacterial as well as eukaryotic cell lysates. Expression of target proteins can be performed either using RTS expression vectors or using PCR-based generation of Linear Template in a screening manner with RTS Linear Template Kits thereby avoiding cloning procedures. PCR products generated with the RTS Linear Template Fab Kit (cat. no. BR1402201) and RTS Linear Template Kit Plus (cat. no. BR1402401) can be applied in the *E. coli*-based RTS 100 *E. coli* HY Kit (cat. no. BR1400101) and in the insect cell-based RTS 100 Insect Membrane Kits (cat. no. BR1401501).

With linear templates, synthesis can be scaled up to approximately 0.5 mg protein in 1 ml synthesis scale, while upscaling with Expression Vectors synthesis is possible up to 50 mg protein in 10 ml synthesis scale using our patented technology (continuous-exchange, cell-free) offering much higher levels of protein expression compared to other *in vitro* T7 RNA polymerase-based transcription/translation systems.

## RTS Platform - Complementary Prokaryotic and Eukaryotic Systems

The Rapid Translation System is designed to cover a broad range of individual proteins by a combined family of pro- and eukaryotic expression systems, eliminating expression limitations of previously available cell-free systems. Synthesis with outstanding high yields of up to 6 mg/ml is realized by RTS *E. coli*, which in most cases is the first choice for expression.

As an alternative, expression of proteins with complex requirements can be achieved using eukaryotic systems which benefit of chaperone systems suited for eukaryotic proteins. Another advantage of eukaryotic RTS is the absence of bacterial endotoxins avoiding laborious purification of proteins for cell-based assays.

Moreover, RTS Wheat Germ combines high synthesis yields up to 1 mg/ml and high success rate independent on codon usage and especially for large proteins which tend to fragmentation in other systems. RTS Insect benefits of endogenous ER based membrane vesicles suited for signal peptide directed translocation of glycosylated proteins and membrane proteins into natural membranes. This system yields protein synthesis up to 50 µg/ml.

In the Rapid Translation System a unique set of innovative technologies is combined to a powerful new protein expression approach. The entire workflow (Figures 1 and 2) has been analyzed to help reduce bottlenecks and accelerate the protein expression process. The different modules of the Rapid Translation System contribute to a new and revolutionary expression concept. A comparison of pro- and eukaryotic RTS Kit for screening and scaled-up protein synthesis is shown in Figure 3.



Figure 1. Workflow of RTS System starting with expression vectors.



Figure 2. Workflow of RTS System starting with linear templates.

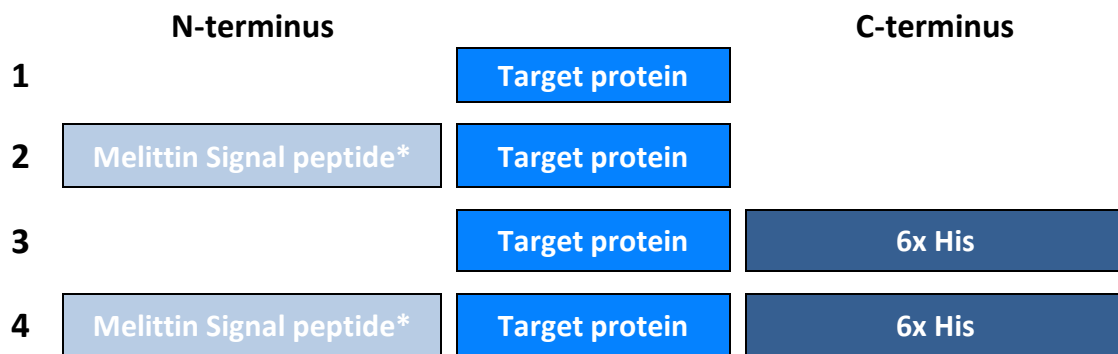
System	Application	Mode	Kit	Expected Yield/Reaction	Linear Templates	Available Reactions	CECF	Eppendorf Thermomixer?
<b>E. coli</b>	Expression and functionality optimization, disulfide bridges, supplements for enhanced solubility/functionality, membrane proteins in detergents, scale-up, purification, NMR, X-ray	Screening	RTS 100 <i>E. coli</i> HY	Up to 20 µg / 50 µl	Yes	24 / 96	No	Recommended
			RTS 100 <i>E. coli</i> Disulfide	Up to 80 µg / 50 µl	No	24	Yes	Required
		Scale-up	RTS 500 ProteoMaster <i>E. coli</i> HY	Up to 6 mg / 1 ml	No	5	Yes	Required
			RTS 500 <i>E. coli</i> Disulfide	Up to 2.5 mg / 1 ml	No	5	Yes	Required
			RTS 9000 <i>E. coli</i> HY	Up to 50 mg / 10 ml	No	1	Yes	Required
<b>Wheat Germ</b>	Screening, expression of non-optimized native eukaryotic sequences and large eukaryotic proteins, scale-up, purification, X-ray	Screening	RTS 100 Wheat Germ CECF	Up to 50 µg / 50 µl	Yes	24	Yes	Required
		Scale-up	RTS 500 Wheat Germ CECF	Up to 1 mg / 1 ml	No	5	Yes	Required
<b>Insect</b>	Screening, membrane proteins in natural membranes, glycosylation	Screening	RTS 100 insect Membrane	Up to 2 µg / 50 µl	Yes	5 / 20	No	Recommended

Figure 3. Comparison of pro- and eukaryotic RTS kits for screening and upscaling.

## Identification of optimal constructs in a single day

The PCR-based RTS Linear Template Fab Kit can be used to generate one or more transcription templates encoding target proteins with an optional C-terminal His6 affinity tag and additionally the option to encode a signal peptide (Figure 4).

The Melittin signal peptide originates from honey bee and is ideally suited to insert proteins into membraneous vesicles contained in the Insect lysate [2]; it may be used to overcome limited production of secreted-, posttranslationally modified-, membrane-, or disulfide-bonded proteins due to inefficient signal peptides. Please note, that the Melittin signal sequence can only be used in combination with the RTS 100 Insect Membrane Kit (cat. no. BR1401501). The PCR products are generated in a 2-step PCR procedure that uses target-specific primers (supplied by the user) and specially designed tag-encoding primers, which encode regulatory elements that optimize transcription and translation in cell-free expression reactions. Having identified the optimal expression template, optionally, the PCR product can be easily cloned into the RTS pIX3.0 Vector (cat. no. BR1402701).



\* only for use in combination with RTS 100 Insect Membrane Kit

Figure 4. The variety of constructs whose expression can be tested using the Linear Template and Protein Synthesis Kits.

The screening of such constructs in small-scale cell-free expression (IVT) reactions using *E. coli*-based protein synthesis kits (or the insect-cell based kits) saves a significant amount of time. The analysis of cell-free expressed proteins by western blotting with tag-specific antibodies enables the identification of the construct or reaction conditions that provide the highest ratio of soluble protein. This procedure also enables a number of different mutation or truncation forms to be tested to find the best-expressing construct. Expression efficiencies obtained using *in vitro* systems correlate well with those seen in *in vivo* systems, meaning that the best-expressing construct identified in the screening procedure will usually deliver optimal yields in scaled-up *in vivo* expression.

## Principle and Procedure

The RTS Linear Template Fab Kit uses a two-step PCR process to generate linear DNA templates for *in vitro* translation systems. Using specially designed primers, coding DNA sequence is both amplified and supplemented with regulatory elements required for optimal transcription and translation in cell-free expression systems. Specially designed 5' untranslated regions (UTRs) on the sense adapter primer sequences reduce the formation of secondary structure in the translation initiation region, one of the commonest causes of low expression rates. A His-tag can be added to the C-terminus, greatly simplifying protein purification and detection. The fast procedure enables researchers to discover the optimal template structure within a single working day (Figure 5).

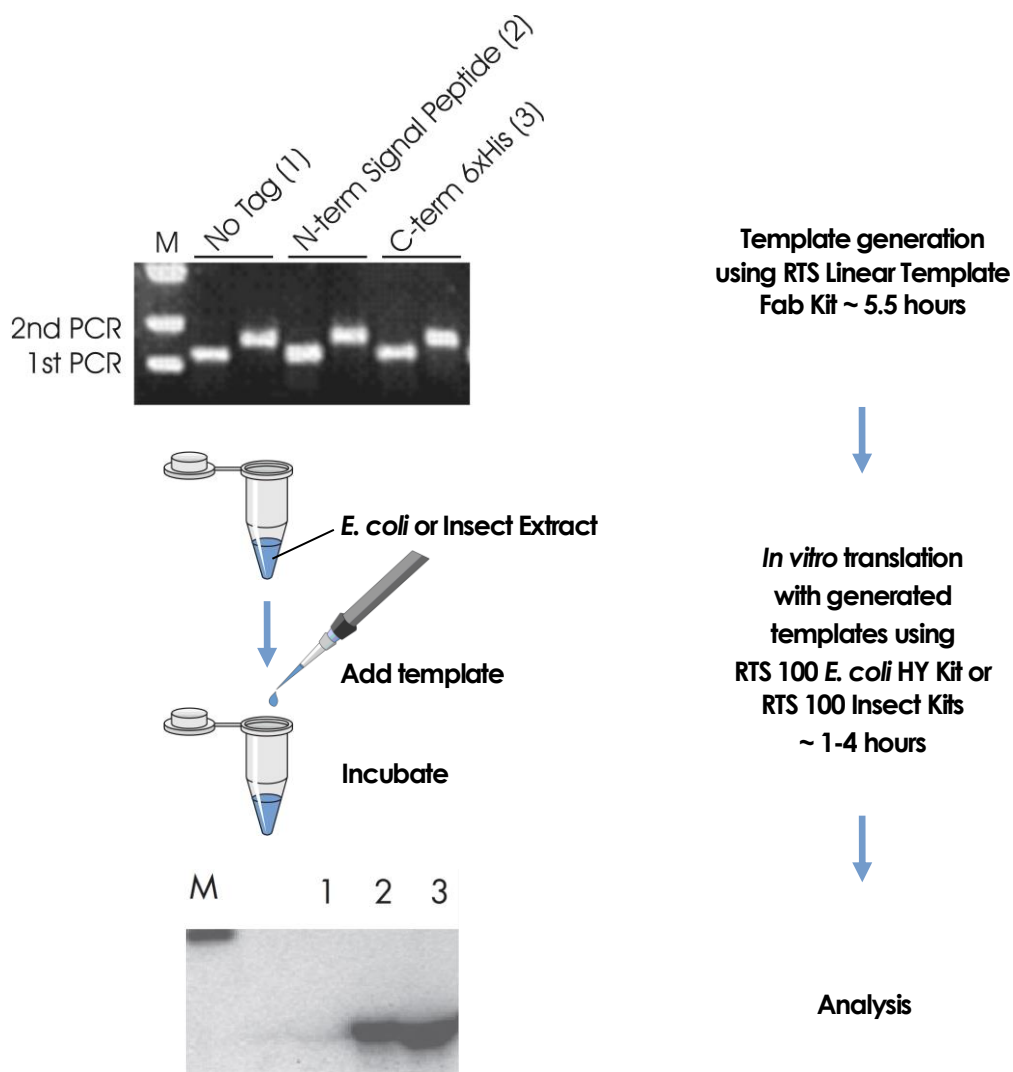


Figure 5. Identification of the optimal expression construct maximizes yields in large-scale *in vitro* reactions or *in vivo* expression. The western blot was probed with anti-His- and Strep-tag antibodies. Therefore, the untagged protein is not detected.

The RTS pIX3.0 Vector (see vector map on page 27) has been developed to enable easy cloning of PCR products generated using the RTS Linear Template Fab Kit and the RTS Linear Template Kit Plus. Its multiple cloning site (MCS) is compatible with restriction sites in the

sense and antisense adapter primers supplied with the RTS Linear Template Kits (see page 26). Once cloned into pIX3.0, expression constructs can be used to generate larger amounts of protein in large-scale protein synthesis reactions. Alternatively, the vector can be used to transform *E. coli* cells for conventional *in vivo* expression.

## Generating PCR Products for Use in *In Vitro* Translation Reactions

The RTS Linear Template Fab Kit uses a two-step procedure to generate PCR products suitable for *in vitro* translation in RTS 100 *E. coli* HY Kits and RTS 100 Insect Membrane Kits. In the first step, defined 5'-tails are added to PCR products using gene-specific primers. The 5'-tails serve as hybridization sites for primers used in a second PCR, in which DNA is amplified using adapter primers that code for regulatory elements required for optimal expression in prokaryotic-cell extracts. These elements include a T7 promoter, ribosomal binding site, and T7 terminator. The resulting PCR products contain multiple cloning sites that are compatible for cloning into the Vector pIX3.0 (Figure 6).

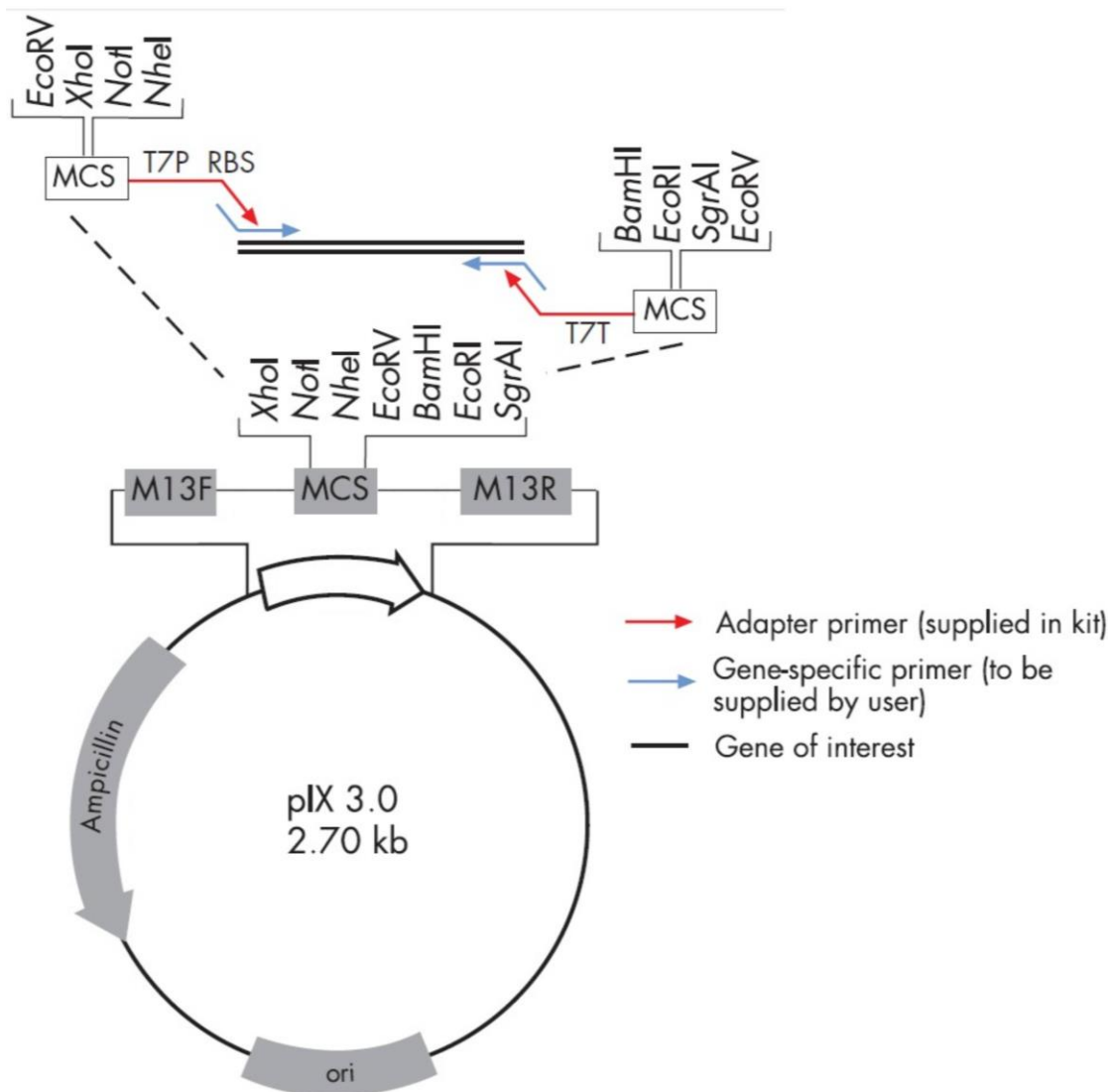


Figure 6. Straightforward cloning into vector RTS pIX3.0 for expression scale-up. T7P: T7promoter; RBS: Ribosome binding site; T7T: T7 terminator; MCS: multiple cloning site.

Adapter primers that encode an optional C-terminal His6 affinity tag and an optional N-terminal signal peptide are contained in the kit. Addition of affinity tags to constructs greatly facilitates purification and detection of expressed proteins. Addition of affinity tags to constructs greatly facilitates purification and detection of expressed proteins. The kit contains enough reagents for 96 two-step PCRs, with each reaction yielding enough expression template for 3–4 *in vitro* translation reactions. The final PCR product can be added to the *in vitro* translation reaction without any further purification steps. The user must provide a DNA template encoding the protein of interest and two gene-specific PCR primers

Plasmid DNA, genomic DNA mixtures, or cDNA mixtures can be used as a template for the first PCR. Alternatively, cDNA can be generated by reverse transcription PCR (RT-PCR) using a gene-specific antisense primer (with defined 5' tail sequence, see Table 1), total RNA, and a reverse transcriptase (e.g. MMuLV Reverse Transcriptase, cat. no. BR0400201). Rules for the design of the gene-specific PCR primers containing the relevant 5' tails are given below. For optimal expression using PCR products in *in vitro* translation reactions, XE-Solution is provided. XE-Solution is added to *in vitro* translation reactions where it protects linear DNA from degradation by exonucleolytic nucleases.

### Strategy for designing gene-specific primers

Prerequisites for successful PCR include the design of optimal gene-specific primer pairs, the use of appropriate primer concentrations, and the correct storage of primer solutions. Primers, for best results, should be HPLC-purified.

The final PCR product added to the *in vitro* translation reaction is generated by a two-step PCR procedure. In the first PCR, primers must be designed that are not only specific for the protein of interest, but also provide 5' tails that will act as hybridization sites for adapter primers used in the second PCR (see Figure 6 on page 14 and Tables 1 and 2). The first step in designing primers is to decide whether a His6 affinity tag should be attached to the C-terminus. Use the 5'-end sequence information in Table 1 and the information below to design forward (sense) and reverse (antisense) primers for protein constructs with or without a C-terminal His6-tag and with or without a N-terminal signal peptide.

### Length

The length of the gene-specific sequence should be 17–20 bases (see Table 1). This may be reduced or increased in some cases to give primers suitable for the annealing temperature of 50°C.

### Melting temperature ( $T_m$ )

The optimal melting temperature ( $T_m$ ) for primers used with the RTS Linear Template Fab Kit is 55°C. The optimal annealing temperature is 5°C below  $T_m$ .

Simplified formula for estimating melting temperature ( $T_m$ ):

$$T_m = 2^\circ\text{C} \times (\text{A}+\text{T}) + 4^\circ\text{C} \times (\text{G}+\text{C})$$

Whenever possible, design primer pairs with similar  $T_m$  values.

**Table 1. Sense and Antisense Primer Pairs Required for Second-Round PCR.**

<b>Desired feature(s)</b>	<b>Gene-specific sense primer</b>	<b>Gene-specific antisense primer</b>
<b>No tag</b>	5'-AGA AGG AGA TAA ACA + <b>ATG</b> + 17 nt target sequence (ATG = start codon)	5'-C TTG GTT AGT TAG TTA + <b>TTA</b> + 20 nt target sequence (TTA = stop codon)
<b>C-terminal 6xHis tag</b>	5'-AGA AGG AGA TAA ACA + <b>ATG</b> + 17 nt target sequence (ATG = start codon)	5'-TG GTG ATG GTG ACC CCA + 20 nt target sequence (ensure that a stop codon from target sequence does not prevent tag expression)
<b>Melittin signal peptide without C-terminal tag</b>	5'-GTA TAC ATT TCT TAC ATC TAT GCG GAC + 20 nt sense strand of target sequence	5'-C TTG GTT AGT TAG TTA + <b>TTA</b> + 20 nt target sequence (TTA = stop codon)
<b>Melittin signal peptide and C-terminal 6xHis tag</b>	5'-GTA TAC ATT TCT TAC ATC TAT GCG GAC + 20 nt sense strand of target sequence	5'-TG GTG ATG GTG ACC CCA + 20 nt target sequence (ensure that a stop codon from target sequence does not prevent tag expression)

**Table 2. Sense and Antisense Primer Pairs Required for Second-Round PCR.**

<b>Desired feature(s)</b>	<b>N-terminus sense primer* (yellow screw-cap)</b>	<b>C-terminus antisense primer* (brown screw-cap)</b>
<b>No tag</b>	No tag Sense Primer	No tag Antisense Primer
<b>C-terminal 6xHis tag</b>	No tag Sense Primer	6xHis tag Antisense Primer
<b>Melittin signal peptide without C-terminal tag</b>	N-term. Signal Sense Primer	No tag Antisense Primer
<b>Melittin signal peptide and C-terminal 6xHis tag</b>	N-term. Signal Sense Primer	6xHis tag Antisense Primer

\* Nucleotide sequences of primers can be found in the Appendix on page 25.

# Protocol: Two-Step PCR Procedure for Generating an Expression Template

This protocol is made up of two separate PCR procedures. In the first PCR, protein-specific sequence is used as a template. The primers used in this first PCR add sequences that will serve as hybridization sites in a second round of PCR (see Figure 6). In the second PCR, adapter primers (see Table 2) are used to add sequence that encodes regulatory elements required for efficient expression and optional affinity tags.

## Important points before starting

- Decide which DNA polymerase you want to use. The kit is adapted to HotStar HiFidelity DNA Polymerase (Qiagen, cat. no. 202602). However, other DNA polymerases may also give satisfactory results (e.g. Hot Start *Taq* DNA Polymerase, cat. no. BR0200101). If a non-hot-start DNA polymerase is used, preheat the thermocycler to the polymerases denaturing temperature. Subsequently put the PCR tubes in the thermocycler and start the temperature program immediately.
- Wear gloves for all working steps in order to protect the reaction components from contaminating DNA and nucleases.
- Use DNase- and RNase-free filter pipet tips.
- Avoid using DNA templates that already contain promoter and terminator elements of the phage T7 gene 10. To avoid contamination of the second PCR with these elements, remove them from the protein-coding sequence using restriction enzymes. Separate the reaction products by agarose gel electrophoresis and purify the target sequence band from the agarose gel using a gel extraction kit (e.g. GenUP PCR/Gel Cleanup Kit, cat. no. BR0700501). Use the purified target sequence DNA for the first PCR.
- When using HotStar HiFidelity DNA Polymerase, to amplify PCR products <2 kb, use values marked with a ● and for PCR products >2 kb, use the values marked with a ▲.
- The optimal Mg<sup>2+</sup> concentration should be determined empirically, but in most cases the standard concentration provided in the PCR buffer of the purchased enzyme will produce satisfactory results.
- Lyophilized primers should be dissolved in a small volume of TE buffer (10 mM Tris·HCl; 1 mM EDTA, pH 8.0) to make a concentrated stock solution.
- Prepare small aliquots of primer working solutions (10 μM) to avoid repeated thawing and freezing. Store all primer solutions at –20°C.
- It is recommended to use a thermocycler with heated lid.
- All experimental procedures – if not stated otherwise – should be performed on ice.

## Procedure: First PCR using gene-specific primers

Please note, that the protocol is an example for **HotStar HiFidelity DNA Polymerase**. When using another enzyme, please follow the instructions of the corresponding DNA polymerase.

1. **Thaw 5x HotStar HiFidelity PCR Buffer (contains dNTPs)<sup>†</sup>, primer solutions, and, if required, 25 mM MgSO<sub>4</sub>. Mix the solutions thoroughly before use.**
2. **Prepare a master mix according to Table 3.**
3. **Mix the master mix thoroughly, and dispense appropriate volumes into PCR tubes.**

**Table 3. Reaction composition** (Example for HotStar HiFidelity DNA Polymerase)

Component	Volume/reaction	Final concentration
<b>Master Mix</b>		
5x HotStar HiFidelity PCR Buffer (contains dNTPs) <sup>†</sup>	5 $\mu$ l	1x
Sense primer	0.75 – 2.5 $\mu$ l	0.3 – 1 $\mu$ M
Antisense primer	0.75 – 2.5 $\mu$ l	0.3 – 1 $\mu$ M
HotStar HiFidelity DNA Polymerase (2.5 units/ $\mu$ l) <sup>‡</sup>	● 0.5 $\mu$ l ▲ 1 $\mu$ l	● 1.25 U ▲ 2.5 U
RNase-free water	Variable	—
<b>Template DNA</b>		
Template DNA (added in step 4)	Variable	50–500 ng genomic DNA 1–100 ng cDNA 0.5–5.0 ng plasmid DNA
<b>Total volume</b>	<b>25 <math>\mu</math>l</b>	

<sup>†</sup> Contains optimized concentration of dNTPs, 7.5 mM MgSO<sub>4</sub>, and Factor SB.

<sup>‡</sup> Dependent on expected PCR product length. In general, use ● 0.5  $\mu$ l HotStar HiFidelity DNA Polymerase when amplifying PCR products <2 kb and ▲ 1  $\mu$ l enzyme when amplifying PCR products >2 kb.

4. **Add template DNA to the individual tubes containing the master mix.**
5. **When using a thermal cycler with a heated lid, do not use mineral oil. Proceed directly to step 6. Otherwise, overlay with approximately 25  $\mu$ l mineral oil.**
6. **Program the thermal cycler according to the manufacturer's instructions.**  
Each PCR program must start with an initial heat-activation step at 95°C for 5 min.

**Table 4. Cycling protocol 1<sup>st</sup> PCR** (Example for HotStar HiFidelity DNA Polymerase)

	<b>Time</b>	<b>Temp.</b>	<b>Comments</b>
<b>Initial activation step</b>	5 min	95°C	HotStar HiFidelity DNA Polymerase is activated by this heating step.
<b>3-step cycling</b>			
Denaturation	15 s	94°C	
Annealing	1 min	50°C	Approximately 5°C below T <sub>m</sub> of primers.
Extension	1 min/kb (PCR products of 1–2 kb)	72°C	
	2 min/kb (PCR products >2kb)	68°C	
<b>Number of cycles</b>	30–45 cycles		The number of cycles is dependent on the origin of the template DNA (see footnote)
<b>Final extension</b>	10 min	72°C	

\* Use 30 cycles if amplifying PCR products from plasmid DNA and 40–45 cycles when amplifying PCR products from genomic DNA or cDNA.

**7. Place the PCR tubes in the thermal cycler and start the cycling program.**

After amplification, samples can be stored overnight at 2 to 8°C or at -20°C for longer storage.

**8. Analyze 1 µl PCR product on a 0.8–1.5% agarose gel.**

The product of the first PCR should be the dominant band. Use 1 µl (approximately 100 ng DNA) of the first PCR product as template for the second PCR.

## Procedure: Second PCR using adapter primers

Please note, that the protocol is an example for **HotStar HiFidelity DNA Polymerase**. When using another enzyme, please follow the instructions of the corresponding DNA polymerase.

1. **Thaw 5x HotStar HiFidelity PCR Buffer (contains dNTPs)<sup>†</sup>, primer solutions, and, if required, 25 mM MgSO<sub>4</sub>. Mix the solutions thoroughly before use.**
2. **Prepare a master mix according to Table 5.**
3. **Mix the master mix thoroughly, and dispense appropriate volumes into PCR tubes.**

**Table 5. Reaction composition** (Example for HotStar HiFidelity DNA Polymerase)

Component	Volume/reaction	Final concentration
<b>Master Mix</b>		
5x HotStar HiFidelity PCR Buffer (contains dNTPs) <sup>†</sup>	5 $\mu$ l	1x
Sense Adapter Primer*	2 $\mu$ l	
Antisense Adapter Primer*	2 $\mu$ l	
HotStar HiFidelity DNA Polymerase (2.5 units/ $\mu$ l) <sup>‡</sup>	● 0.5 $\mu$ l ▲ 1 $\mu$ l	● 1.25 U ▲ 2.5 U
RNase-free water	Variable	—
<b>Template DNA</b>		
Template DNA (added in step 4)	1–2 $\mu$ l product from first PCR (~100 ng)	
<b>Total volume</b>	<b>25 <math>\mu</math>l</b>	—

<sup>†</sup> Contains optimized concentration of dNTPs, 7.5 mM MgSO<sub>4</sub>, and Factor SB.

<sup>‡</sup> Dependent on expected PCR product length. In general, use ● 0.5  $\mu$ l HotStar HiFidelity DNA Polymerase when amplifying PCR products <2 kb and ▲ 1  $\mu$ l enzyme when amplifying PCR products >2 kb.

\* For possible adapter primer combinations see Table 2, page 17.

4. **Add template DNA (1–2  $\mu$ l of the first PCR, ~100 ng) to the individual tubes containing the master mix.**
5. **When using a thermal cycler with a heated lid, do not use mineral oil. Proceed directly to step 6. Otherwise, overlay with approximately 25  $\mu$ l mineral oil.**
6. **Program the thermal cycler according to the manufacturer's instructions.**

Each PCR program must start with an initial heat-activation step at 95°C for 5 min.

**Table 6. Cycling protocol 2<sup>nd</sup> PCR** (Example for HotStar HiFidelity DNA Polymerase)

	Time	Temp.	Comments
<b>Initial activation step</b>	5 min	95°C	HotStar HiFidelity DNA Polymerase is activated by this heating step.
<b>3-step cycling</b>			
Denaturation	15 s	94°C	
Annealing	1 min	50°C	Approximately 5°C below T <sub>m</sub> of primers.
Extension	1 min/kb (PCR products of 1–2 kb)	72°C	
	2 min/kb (PCR products >2kb)	68°C	
<b>Number of cycles</b>	30 cycles		
<b>Final extension</b>	10 min	72°C	

**7. Place the PCR tubes in the thermal cycler and start the cycling program.**

After amplification, samples can be stored overnight at 2 to 8°C or at –20°C for longer storage.

**8. Analyze 1 µl of the first and second PCR on a 0.8–1.5% agarose gel.**

The introduction of regulatory elements and affinity-tag sequences in the second PCR adds approximately 160–200 bp to the first PCR product.

**9. Determine the yield of the second PCR by comparing the product band to the molecular weight marker bands.**

0.7 µg DNA (~7 µl of the second PCR) is required for a 50 µl *in vitro* translation reaction using RTS 100 *E. coli* HY Kits. 0.5 µg (~5 µl of the second PCR) is required for a 50 µl *in vitro* translation reaction using RTS 100 Insect Membrane Kits.

# Troubleshooting Guide

This troubleshooting guide may be helpful in solving any problems that may arise. The scientists in biotechrabbit Technical Service are always happy to answer any questions you may have about either the information or protocol(s) in this handbook. Many of the PCR-related problems in this guide are comprehensively covered in the instruction manuals of the corresponding DNA polymerase (e.g. HotStar HiFidelity DNA Polymerase PCR Handbook).

---

## Comments and suggestions

---

### ***First and second PCR***

#### **Little or no product**

a) DNA polymerase not activated	Check if PCR was started with an initial incubation step (typically some minutes at > 90°C).
b) Pipetting error or missing reagent	Repeat the PCR. Check the concentrations and storage conditions of reagents, including primers and dNTP mix.
c) Insufficient starting template	Increase amount of starting template used in PCR.
d) Primer concentration not optimal	Recalculate primer concentration.
e) Extension time too short	Increase extension time in increments of 1 min. Ensure that the extension time for large DNA fragments fits to the applied DNA polymerase.
f) Enzyme concentration too low	Some DNA polymerases require increased concentrations for larger DNA fragments.
g) Insufficient number of cycles	Increase the number of cycles in steps of 5 cycles.
h) Problems with starting template	Check the concentration, storage conditions, and quality of starting template. If necessary, make new serial dilutions of template nucleic acid from stock solutions. Repeat PCR using the new stock solutions.
i) Mg <sup>2+</sup> concentration not optimal	Perform PCR with different final concentrations of Mg <sup>2+</sup> from 1.5 to 5 mM.
j) Primer design not optimal	Review primer design, pages 15-16.
k) Incorrect dNTP concentration	Ensure that the dNTP concentration fits to the applied DNA polymerase. Increase dNTP concentration in increments of 50 µM of each dNTP. Do not exceed a concentration of 500 µM of each dNTP since this might lower PCR fidelity.

---

**Product is multi-banded**

a) PCR cycling conditions not optimal	Review primer design, see pages 15-16.
b) Enzyme concentration too low	Some DNA polymerases require increased concentrations for larger DNA fragments.
c) Extension time too short	Increase extension time in increments of 1 min. Ensure that the extension time for large DNA fragments fits to the applied DNA polymerase.
d) Mg <sup>2+</sup> concentration not optimal	Perform PCR with different final concentrations of Mg <sup>2+</sup> from 1.5 to 5 mM.
e) Primer concentration not optimal or primers degraded	Recalculate primer concentration. Particularly when performing highly sensitive PCR, check for possible degradation of the primers on a denaturing polyacrylamide gel.
f) Primer design not optimal	Review primer design, see pages 15-16.
g) Incorrect dNTP concentration	Ensure that the dNTP concentration fits to the applied DNA polymerase. Increase dNTP concentration in increments of 50 µM of each dNTP. Do not exceed a concentration of 500 µM of each dNTP since this might lower PCR fidelity.

---

**Product is smeared**

a) Insufficient starting template	Increase amount of starting template used in PCR.
b) Extension time too short	Increase extension time in increments of 1 min. Ensure that the extension time for large DNA fragments fits to the applied DNA polymerase.
c) Mg <sup>2+</sup> concentration not optimal	Perform PCR with different final concentrations of Mg <sup>2+</sup> from 1.5 to 5 mM.
d) dNTP concentration not optimal	Ensure that the dNTP concentration fits to the applied DNA polymerase. Increase dNTP concentration in increments of 50 µM of each dNTP. Do not exceed a concentration of 500 µM of each dNTP since this might lower PCR fidelity.
e) Primer design not optimal	Review primer design, see pages 15-16.

## References

1. Sambrook, J., Fritsch, E., and Maniatis, T. (1989) *Molecular Cloning — A laboratory Manual*. 2<sup>nd</sup> Ed. Cold Spring Harbor. Cold Spring Harbor Laboratory Press.
2. Merk, H., Gless, C., Maertens, B., Gerrits, M., and Stiege, W. (2012). Cell-free synthesis of functional and endotoxin-free antibody Fab fragments by translocation into microsomes. *Biotechniques* 53,153-160

# Appendix: Adapter Primer Sequences and Multiple Cloning Sites

## No tag Sense Primer

5'-ATGATATCTCGAGCGGCCGCTAGCTAATACGACTCACTATAGGGAGACCACAACGGTTTCCCTCTAGAA  
ATAATTTTGTTTAACTTTAAGAAGGAGATAAACA-3'

H<sub>2</sub>N — Protein

## No tag Antisense Primer

5'-ATGATATCACCGGTGAATTCGGATCCAAAAACCCCTCAAGACCCGTTTAGAGGCCCAAGGGGTACA  
GATCTTGGTTAGTTAGTTATTA-3'

Protein — Stop — COOH

## 6xHis tag Antisense Primer

5'-ATGATATCACCGGTGAATTCGGATCCAAAAACCCCTCAAGACCCGTTTAGAGGCCCAAGGGGTACA  
GATCTTGGTTAGTTAGTTATTAATGATGGTGATGGTGGTGACCCCA-3'

Protein — WGHHHHHH — Stop — COOH

## No tag Sense Primer Signal Peptide

5'-ATGATATCTCGAGCGGCCGCTAGCTAATACGACTCACTATAGGGAGACCACAACGGTTTCCCTCTAGAA  
ATAATTTTGTTTAACTTTAAGAAGGAGATAAACAATGAAATTCTTAGTCAACGTTGCCCTGTTTTTATGGT  
CGTATACATTTCTTACATCTATGCGGAC-3'

H<sub>2</sub>N — MKFLVNVALVFMVVYISYIYA \*D — Protein

\* signal peptide cleavage site

Underlined sequences hybridizes to 5' tails of gene-specific primers.

## Sequencing primers

M13 forward (-20)                      GTAAAACGACGGCCAGT

M13 reverse (-21)                      CAGGAAACAGCTATGAC

## Cloning PCR Products Generated with the RTS Linear Template Fab Kit into the RTS pIX3.0 Vector

The pIX3.0 vector is designed for cloning of PCR products generated by the RTS Linear Template Fab Kit and the RTS Linear Template Kit Plus. The PCR products contain multiple cloning sites that are compatible with the multiple cloning site of the RTS pIX3.0 Vector.

### RTS Sense Primer Multiple Cloning Site

*EcoRV XhoI NotI NheI*

5'-ATGATATCTCGAGCGGCCGCTAGCT-3'  
3'-TACTATAGAGCTCGCCGCGATCGA-5'

### RTS Antisense Primer Multiple Cloning Site

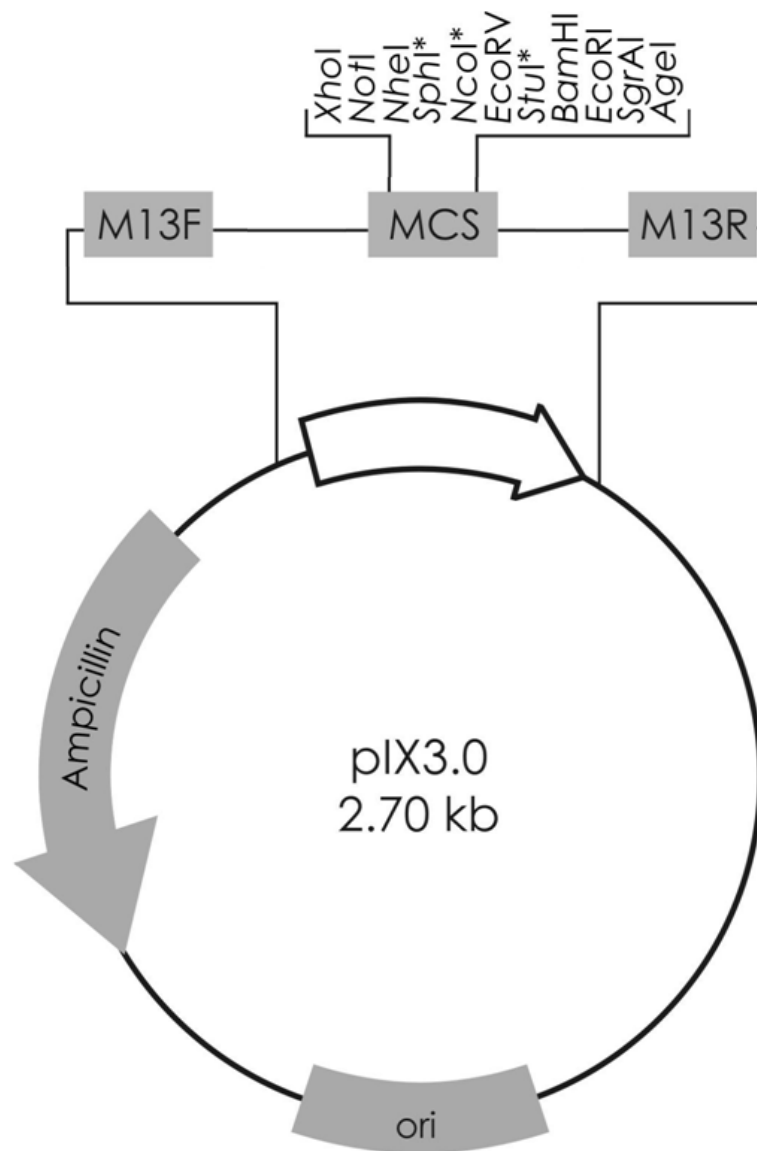
*BamHI EcoRI SgrAI/AgeI EcoRV*

5'-TGGATCCGAATTCTCCGGTGATATCAT-3'  
3'-ACCTAGGCTTAAGTGCCACTATAGTA-5'

### RTS pIX3.0 Vector Multiple Cloning Site

*XhoI NotI NheI EcoRV BamHI EcoRI SgrAI/AgeI*

5'-CTCGAGCGGCCGCTAGCATGCCATGGATATCAGGCCTGGATCCGAATTCACCGGTG-3'  
3'-GAGCTCGCCGCGATCGTACGGTACCTATAGTCCGGACCTAGGCTTAAGTGCCAC-5'



\* As long as they do not have a restriction site within the PCR insert, these restriction enzymes can be added directly to the vector–insert ligation reaction to minimize the number of false positive clones (religations).

Figure 7. The RTS pIX3.0 Vector.

## Ordering information

Product	Size	Order no.
RTS Linear Template Kit Plus	20 reactions	BR1402401
RTS pIX3.0 Vector	1 vector, 25 µg	BR1402701
RTS 100 <i>E. coli</i> HY Kit	24 reactions	BR1400101
RTS 100 <i>E. coli</i> HY Kit	96 reactions	BR1400102
RTS 500 ProteoMaster <i>E. coli</i> HY Kit	5 reactions	BR1400201
RTS 9000 <i>E. coli</i> HY Kit	1 reaction	BR1400301
RTS 100 <i>E. coli</i> Disulfide Kit	24 reactions	BR1400401
RTS 500 <i>E. coli</i> Disulfide Kit	5 reactions	BR1400501
RTS pIVEX <i>E. coli</i> His-tag, 2nd Gen. Vector Set	2 vectors, 10 µg each	BR1400701
RTS Wheat Germ LinTempGenSet, His6-tag	96 reactions	BR1401201
RTS pIVEX Wheat Germ His6-tag Vector Set	2 vectors, 10 µg each	BR1401301
RTS 100 Wheat Germ Kit	24 reactions	BR1402501
RTS 100 Wheat Germ CECF Kit	24 reactions	BR1401001
RTS 500 Wheat Germ CECF Kit	5 reactions	BR1401101
RTS 500 Adapter	1 adapter	BR1401901
RTS GroE Supplement	For 5 reactions of 1 ml	BR1401701
RTS DnaK Supplement	For 5 reactions of 1 ml	BR1401601
RTS Amino Acid Sampler	1 set	BR1401801
RTS 100 Insect Membrane Kit	5 reactions	BR1401501
RTS 100 Insect Membrane Kit	20 reactions	BR1401502
RTS Linear Template Fab Kit	96 reactions	BR1402201
RTS pIX4.0 Insect Vector	1 vector, 25 µg	BR1400901



