



---

## RTS™ 100 *E. coli* HY Kit Manual

---

For cell-free expression of functionally active protein from  
circular and linear templates

RTS 100 *E. coli* HY Kit Manual, 21.03.2025

© 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; GenUp™, ProteoMaster™, RTS™, 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, RTS 500 Wheat Germ, RTS 100 *E. coli* Disulfide, RTS 500 *E. coli* Disulfide, RTS 500 *E. coli* and RTS 9000 *E. coli* 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

Product description	4
Product limitations	4
Materials supplied	4
Additional materials	6
Shipping and storage conditions	6
Safety information	6
Quality assurance	7
Product warranty	7
Product principle	8
Protocol 1: Preparation of DNA for <i>in vitro</i> expression	11
Protocol 2: Protein synthesis reaction	14
Short protocol: Reconstitution of reaction components and running a reaction	21
Optimizing SDS-polyacrylamide gel electrophoresis (SDS-PAGE)	23
Typical results	23
Application: Radioactive labeling with L-[ <sup>35</sup> S] Methionine	24
References	25
Troubleshooting guide	28

## Product specifications

The RTS 100 *E. coli* HY Kit is designed for:

- rapid parallel protein synthesis reactions
- compatibility with PCR generated templates and plasmids
- rapid optimization of expression constructs
- rapid functional testing of PCR-generated mutations
- expression of toxic gene products
- synthesis of truncated gene products from PCR-generated or restriction enzyme digested DNA, for epitope or functional domain mapping
- successful synthesis of proteins in the molecular weight range from 10–120 kDa
- Incorporation of radiolabeled Methionine into the protein

## Product description

The RTS 100 *E. coli* HY Kit provides the components and procedures necessary for 6 (cat. no. BR1400106), 24 (cat. no. BR1400101) or 96 (cat. no. BR1400102) coupled transcription/translation reactions of 50  $\mu$ l.

## Product limitations

The RTS 100 *E. coli* HY Kit is developed, designed, and sold for research purposes only. It is 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 the materials described in this text.

## Materials supplied

Kit	Catalog no.	Size
RTS 100 <i>E. coli</i> HY Kit	BR1400106	6 reactions
RTS 100 <i>E. coli</i> HY Kit	BR1400101	24 reactions
RTS 100 <i>E. coli</i> HY Kit	BR1400102	96 reactions

RTS 100 *E. coli* HY Kit

Kit	Contents and function	No. included		
		6 reactions	24 reactions	96 reactions
Ordering number		BR1400106	BR1400101	BR1400102
<i>E. coli</i> Lysate; 100 <i>E. coli</i> (Bottle 1, red cap)	→ Lysate from <i>E. coli</i> ; stabilized and lyophilized  → Contains components for transcription and translation	1 vial	1 bottle	4 bottles
Reaction Mix; 100 <i>E. coli</i> (Bottle 2, green cap)	→ Substrate mix to prepare Reaction Solution; stabilized and lyophilized	1 vial	1 bottle	4 bottles
Amino Acids; 100 <i>E. coli</i> (Bottle 3, brown cap)	→ Mix of 19 amino acids without Methionine to prepare Reaction Solution; stabilized and lyophilized	1 vial	1 bottle	4 bottles
Methionine; 100 <i>E. coli</i> (Bottle 4, yellow cap)	→ Methionine to prepare Reaction Solution; stabilized and lyophilized	1 vial	1 bottle	1 bottle
Reconstitution Buffer; 100 <i>E. coli</i> (Bottle 5, white cap)	→ Buffer solution for the reconstitution of bottles 1, 2, 3, and 4	1 vial	1 bottle	4 bottles
Control Vector GFP; 100 <i>E. coli</i> (Vial 6, colorless cap)	→ Lyophilized plasmid GFP (green fluorescent protein) expression vector with C-terminal His <sub>6</sub> -tag for the control reaction	1 vial	1 vial	1 vial
Reaction Tubes + Caps or Microplate + Adhesive Films; 100 <i>E. coli</i>	→ Reaction tubes or microplate	8 x 200 µl reaction tubes	24 x 200 µl reaction tubes	Module with 12 strips (8 wells)

## Additional materials

To perform the protocols described in this manual, the following additional materials must be provided by the user:

- Eppendorf® ThermoMixer® C (or Eppendorf® Thermomixer Comfort). Alternatively, a water bath or heater adjustable to 30°C can be used. Optimal results are obtained if the reaction is shaken during incubation.
- Pipets 0–10 µl, 10–200 µl, 200–1,000 µl,
- Pipette tips autoclaved at 121°C for 20 min (RNase free)
- Besides the template vector coding for the protein of interest, pIVEX cloning vectors (cat. no. BR1400701) or the RTS Linear Template Kit Plus (cat. no. BR1402401) PCR reagents to create a linear DNA expression construct are needed. N- and C-terminal tags for affinity purification of synthesized protein can be added to the gene of interest by using the RTS Linear Template Kit Plus. The addition of the N-terminal tags often greatly enhances the yield of the synthesized protein.
- [<sup>35</sup>S]-Methionine (>1,000 Ci/mmol at 15 mCi/ml) is required for optional radioactive labeling
- Deionized DNase- and RNase-free water is required for reconstitution of the GFP Control Vector
- Acetone is required for precipitating samples prior to SDS-gel electrophoresis.

For convenience, additional materials to be supplied by the user are listed at the beginning of the protocol for which they are required.

## Shipping and storage conditions

The RTS 100 *E. coli* HY Kit is shipped on dry ice.

The RTS 100 *E. coli* HY Kit and components should be stored in the dark at -15 to -25°C and are stable until the expiration date printed on the label. Avoid repeated freezing and thawing.

## Safety information

All due care and attention should be exercised in the handling of this product. We recommend all users of biotechrabbit 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.

## 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 biotechrabbit ([www.biotechrabbit.com](http://www.biotechrabbit.com)). Certificates of Analysis are available on request.

## 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.

# Protocols

## Product principle

### Introduction

The Rapid Translation System (RTS) workflow (Figure 1) combines a series of new technologies for efficient and optimized protein expression. They overcome the limitations that often restrict the use of cell-free systems. These innovations include software-based template optimization, generation of stable expression templates without cloning, optimization of *in vitro* expression conditions and high yield *in vitro* expression using the continuous exchange cell free principle (CECF) and an optimized lysate biochemistry.

The RTS 100 *E. coli* HY Kit allows the expression of up to 20 µg functionally active protein in a 50 µl reaction from circular and linear templates within 6 hours. It is therefore suited for the rapid testing and optimization of expression constructs as well as for the screening of many expressions in parallel.

Note: This kit cannot introduce post-translational glycosylation, phosphorylation, disulfide bond formation or signal sequence cleavage.

Advantages of the system include:

- fast
  - express 2-20 µg protein within 6 hours
- convenient
  - protein expression reaction preparation takes less than 30 minutes
  - no fermentation and therefore no safety regulations for recombinant organisms
- ability to express multiple proteins in parallel
- flexible
  - alter reaction conditions to overcome insolubility or folding problems
  - express toxic proteins
  - incorporate radiolabeled Methionine
- reliable
  - kits are tested to yield at least 300 µg/ml GFP
  - lyophilized bacterial lysate offers unprecedented stability
  - good reproducibility

## Description of procedure

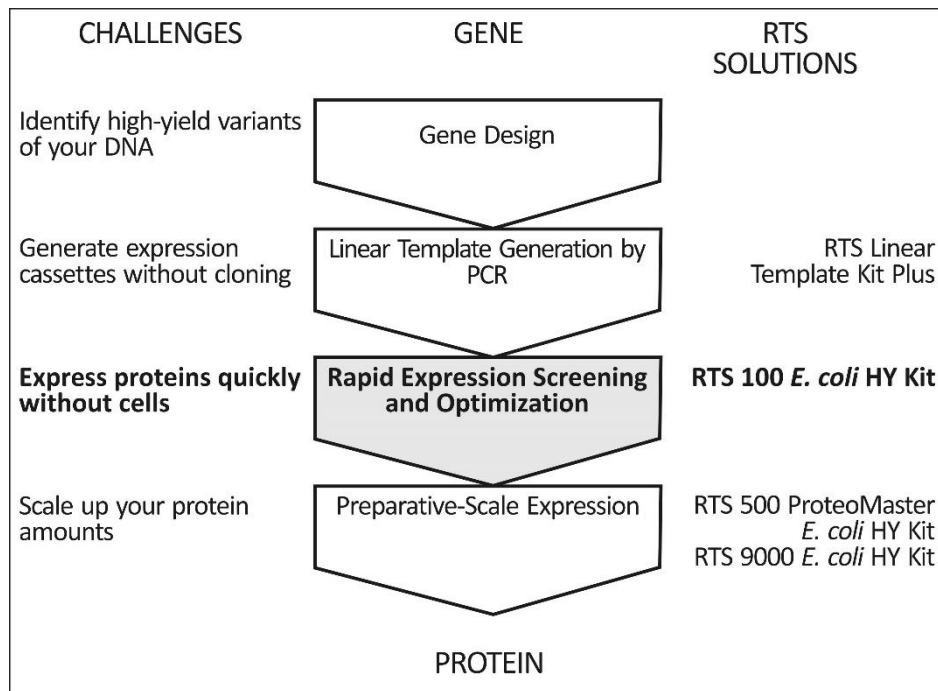


Figure 1. Integration of the RTS 100 *E. coli* HY Kit into the RTS workflow.

## Reaction principle

RTS uses a coupled transcription/translation reaction for *in vitro* protein synthesis. Transcription and translation take place simultaneously in the reaction: While the T7-RNA-polymerase transcribes the template gene, the ribosomes provided by the *E. coli* lysate start to translate the 5'-end of the nascent mRNA (Figure 2). This makes the system highly productive compared with the use of isolated mRNA (1). The *E. coli* lysate is prepared with some modifications according to the method of Zubay (2), resulting in the biochemically enhanced High-Yield (HY) *E. coli* lysate (3, 4). An *E. coli* strain with low exonuclease activity was selected and growth conditions were optimized to allow optimum protein expression from linear (PCR-generated) and plasmid templates.

Regulatory elements, such as T7 promoter, ribosomal binding site and T7 terminator are added to the gene of interest by PCR or cloning using a pIVEX vector. The resulting DNA template is then added to the Reaction Solution.

In a coupled *in vitro* reaction, the template DNA is first transcribed into mRNA by T7 RNA polymerase, followed by translation into protein by the ribosomal machinery present in the *E. coli* lysate. The expressed protein accumulates during the reaction and is harvested after 1-6 hours. The reaction will reach 90% completion after 4 hours.

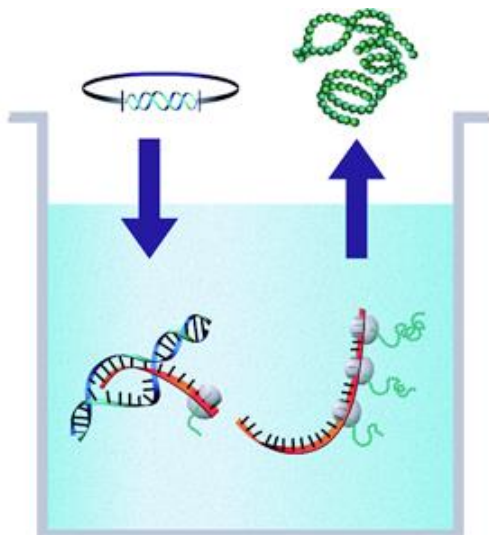


Figure 2. Schematic illustration of the coupled transcription/translation reaction. Firstly, an expression plasmid or linear template DNA carrying the gene of interest is added to the reaction compartment. In a coupled *in vitro* reaction, the DNA is first transcribed from the template vector into mRNA by T7 RNA polymerase, followed by translation by the ribosomal machinery present in the *E. coli* lysate. Expressed protein accumulates in the reaction compartment and is harvested after 1-6 hours.

## Template DNA

Linear and circular expression templates are suited. Use pIVEX vectors (cat. no. BR1400701), the RTS Linear Template Kit Plus (cat. no. BR1402401) or a simple one-step PCR to attach T7 transcription/ translation regulatory elements to the target gene. The PCR product can be easily cloned into the RTS pIX3.0 Vector (cat. no. BR1402701, see handbook of RTS Linear Template Kit Plus).

Other vectors, designed for prokaryotic *in vitro* protein expression and containing a T7 promoter as well as a ribosomal binding site may also be used. The requirements are described in Protocol 1, 'Use of expression vectors', page 13.

## Protocol 1: Preparation of DNA for *in vitro* expression

### Template optimization

The expression yield of a given gene is strongly influenced by its mRNA structure. Interactions between the coding sequence and the 5'-untranslated region can interfere with translation initiation and therefore have an impact on translation efficiency. In the case of low expression yields, biotechrabbit recommends that the sequence is optimized. For example, N-terminal tag sequences or silent mutations in the first few codons can be introduced using the RTS Linear Template Kit Plus and may help to increase expression and protein yield.

### Generation of expression templates by PCR: Principal applications

The use of PCR protocols instead of subcloning allows rapid expression from new or modified DNA-templates:

- generation of a DNA expression template by a two-step expression PCR protocol (by addition of a T7 promoter, a prokaryotic ribosomal binding site and other regulatory elements to the gene of interest)
- addition of Tag sequences for detection with antibodies and purification
- optimization of the gene sequence for expression (see Protocol 1, 'Use of expression vectors', page 13)
- introduction of mutations
- change of codon usage
- truncation of proteins
- domain screening

### Generation of a linear DNA expression template by expression PCR

The RTS Linear Template Kit Plus (cat. no. BR1402401) is recommended to generate a linear DNA template for the *in vitro* expression via a two-step expression PCR protocol. It provides a DNA for the introduction of the regulatory elements, as well as different N- or C-terminal tags.

#### Procedure

1. Select the terminus and kind of tag you wish to add to the gene of interest.
2. Design a gene-specific sense primer that contains 17–20 nucleotides homologous to the gene of interest and additional bases overlapping the chosen tag region, as indicated in the RTS Linear Template Kit Plus manual.
3. Design a gene-specific antisense primer that contains 17–20 nucleotides homologous to the gene of interest and additional bases overlapping the chosen tag region, as indicated in the RTS Linear Template Kit Plus. (For an example, see 'Typical results', page 23.)
4. Perform 30 cycles of a 25 µl PCR using these primers and the template DNA containing the gene of interest as indicated in the RTS Linear Template Kit Plus manual.

5. Perform 30 cycles of a second 25  $\mu$ l PCR using 100 ng (1-2  $\mu$ l) of the first PCR product as template together with the primers and DNA supplied with the RTS Linear Template Kit Plus containing regulatory elements for prokaryotic expression and C- or N-terminal tags.
6. Determine the concentration of the PCR product densitometrically on an agarose gel by comparison to DNA standards.
7. Use 500 ng PCR product for a 50  $\mu$ l *in vitro* protein synthesis reaction.

Note: If necessary, use a PCR product purification kit (e.g. GenUP™ PCR/Gel Cleanup Kit, cat. no. BR0700501 or cat. no. BR0700301) to remove potentially inhibitory primer-dimers. Do not purify the PCR product using agarose gels because this treatment inhibits *in vitro* protein synthesis.

### Addition of a C-terminal epitope tag sequence by a one-step PCR protocol

Standard PCR reactions may be used to introduce epitope tags or mutations into the already cloned gene of interest.

#### Procedure

1. Design a sense primer that is located upstream of a T7 promoter sequence in a prokaryotic expression plasmid with the gene of interest.  
This plasmid has to contain a T7 promoter sequence and the necessary regulatory elements as detailed in 'Template optimization', page 11.
2. Design an antisense primer that contains the T7 terminator sequence, a spacer region of at least 25 nucleotides distance to the stop codon followed by the tag sequence and a 15–20 base-pair sequence homologous to the gene of interest.
3. Perform a PCR using these primers and follow steps 4 and 5 of the procedure on this page.
4. Use 500 ng (~5  $\mu$ l) of the PCR product for a 50  $\mu$ l *in vitro* protein synthesis reaction.
5. Protocol for a 50  $\mu$ l PCR. We recommend using Hot Start Taq DNA Polymerase, (cat. no. BR0200102). Combine the following (example for Hot Start Taq DNA Polymerase):

31  $\mu$ l H<sub>2</sub>O

5  $\mu$ l 10 $\times$  Reaction Buffer

1.5  $\mu$ l 50 mM MgCl<sub>2</sub>

1  $\mu$ l 10 mM dNTP Mix

5  $\mu$ l sense primer (3 pmol/ $\mu$ l)

5  $\mu$ l antisense primer (3 pmol/ $\mu$ l)

1  $\mu$ l of template DNA (5 ng/ $\mu$ l)

0.5  $\mu$ l 5 U/ $\mu$ l Hot Start Taq DNA Polymerase

Run the following PCR cycles:

5 min / 94°C

+ 30x (30 sec / 94°C + 45 sec / 53°C + 1 min/kb / 72°C)

5 min / 72°C

+ 4°C

## Use of expression vectors

### Necessary vector elements

Any vector or linear DNA to be used in combination with RTS must include the following elements and structural features:

- target gene under control of T7 promoter located downstream of a RBS (ribosomal binding site) sequence
- distance between T7 promoter and start ATG should not exceed 100 base pairs
- distance between the RBS sequence and start ATG should not be more than 5–8 base pairs
- T7 terminator sequence at the 3'-end of the gene

### General recommendations

The pIVEX vector family has been developed and optimized for use in the RTS. biotechrabbit recommends cloning target genes into a pIVEX vector prior to expression (see 'Generation of expression templates', page 11).

Maps of some of the available pIVEX vectors are shown schematically in Figure 3. For more information, visit [www.biotechrabbit.com](http://www.biotechrabbit.com), cat. no. BR1400701

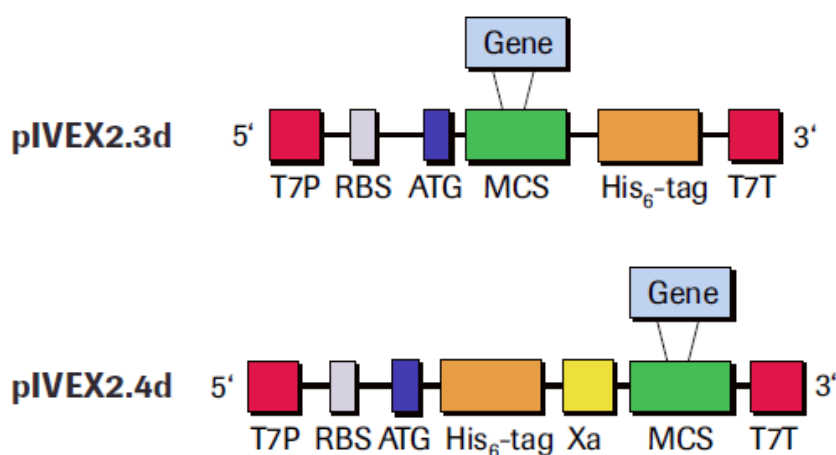


Figure 3. Functional elements of cloning vectors.

**T7P:** T7 Promoter; **RBS:** Ribosome binding site; **ATG:** Start codon; **Xa:** Factor Xa restriction protease cleavage site; **MCS:** Multiple cloning site for the insertion of the target gene; **T7T:** T7 Terminator.

## Purity of the plasmid preparation

Plasmids obtained from commercially available DNA preparation kits (e.g. GenUP™ Plasmid Kit, cat. no. BR0700201) are usually pure enough to be used as template in the RTS. If DNA is not pure enough ( $OD_{260/280} < 1.7$ ), use phenol extraction to remove traces of RNase from the preparation, which may enhance its performance in the expression reaction.

Note: do not purify DNA-fragments from agarose gels, as this treatment inhibits *in vitro* protein synthesis.

## Protocol 2: Protein synthesis reaction

### Equipment and reagents required

- DNA template: Prepare and purify the DNA template as described in Protocol 1, page 11.
- Use the Eppendorf® ThermoMixer® C (or Eppendorf® Thermomixer Comfort) or a water bath or an incubator, adjusted to 30°C. Shaking of the reaction solution during the reaction results in higher yields.
- Calibrated pipets
- UV lamp (360 nm) for the detection of GFP (control reaction)
- RNase-free plastic and glassware

### Reagent notes

- Do not combine reagents from different kit lots
- Reconstitute only the bottles needed for the experiment
- For reconstitution of bottles 1 to 4, use only the Reconstitution Buffer supplied with the kit (bottle 5). For reconstitution of the plasmid (bottle 6), use sterile DNase- and RNase-free water
- Store reconstituted solutions of bottles 1 to 4 at –15 to –25°C. The reagents can withstand two freeze–thaw cycles without significant decrease in activity. Lysate (bottle 1) should be frozen in liquid nitrogen and subsequently stored at –15 to –25°C
- Reconstitution Buffer can be thawed in a water bath at 25°C
- Store reconstituted plasmid at –15 to –25°C.
- Reconstitute the lyophilized reagents or thaw solutions immediately before use
- Keep reagents and working solutions on ice

## Procedure

1. Kit format for **6 reactions** (BR1400106): reconstitute the reaction components according to Table 1.

Kit format for **24 and 96 reactions** (BR1400101 and BR1400102): reconstitute the reaction components according to Table 2.

Table 1. Reaction components BR1400106 (6 reactions)

Solution	Contents	Reconstitution procedure
1	<i>E. coli</i> Lysate; 100 <i>E. coli</i> (6 rxns) (Vial 1, red cap)	Reconstitute the lyophilizate with 90 $\mu$ l Reconstitution Buffer. Do not vortex!
2	Reaction Mix; 100 <i>E. coli</i> (6 rxns) (Vial 2, green cap)	Reconstitute the lyophilizate with 75 $\mu$ l Reconstitution Buffer
3	Amino Acids w/o Methionine; 100 <i>E. coli</i> (6 rxns) (Vial 3, brown cap)	Reconstitute the lyophilizate with 90 $\mu$ l Reconstitution Buffer
4	Methionine; 100 <i>E. coli</i> (6 rxns) (Vial 4, yellow cap)	Reconstitute the lyophilizate with 82.5 $\mu$ l Reconstitution Buffer
5	Reconstitution Buffer; 100 <i>E. coli</i> (6 rxns) (Vial 5, white cap)	→ 0.4 ml → Ready-to-use solution → The solution is stable at 2–8°C, but can also be stored at –15 to –25°C
6	Control Vector GFP; 100 <i>E. coli</i> (6 rxns) (Vial 6, colorless cap)	→ Briefly centrifuge down the contents of the vial → Reconstitute the lyophilizate (15 $\mu$ g of Control Vector GFP) with 15 $\mu$ l of sterile DNase- and RNase-free water → The solution is stable at –15 to –25°C

All reconstituted solutions should be clear, with the exception of the *E. coli* lysate, which remains cloudy.

Table 2. Reaction components BR1400101 (24 reactions) and BR1400102 (96 reactions)

Solution	Contents	Reconstitution procedure	For use in
1	<i>E. coli</i> Lysate; 100 <i>E. coli</i> (Bottle 1, red cap)	Reconstitute the lyophilizate with 0.36 ml of Reconstitution Buffer (bottle 5), mix carefully by rolling or gentle shaking. Do not vortex!	Step 2 Solution 7
2	Reaction Mix; 100 <i>E. coli</i> (Bottle 2, green cap)	Reconstitute the lyophilizate with 0.30 ml of Reconstitution Buffer (bottle 5), mix by rolling or shaking.	Step 2 Solution 7
3	Amino Acids w/o Methionine; 100 <i>E. coli</i> (Bottle 3, brown cap)	Reconstitute the lyophilizate with 0.36 ml of Reconstitution Buffer (bottle 5), mix by rolling or shaking.	Step 2 Solution 7
4	Methionine; 100 <i>E. coli</i> (Bottle 4, yellow cap)	Reconstitute the lyophilizate with 0.33 ml of Reconstitution Buffer (bottle 5), mix by rolling or shaking.	Step 2 Solution 7
5	Reconstitution Buffer; 100 <i>E. coli</i> (Bottle 5, white cap)	→ 1.6 ml → Ready-to-use solution → The solution is stable at 2–8°C, but can also be stored at –15 to –25°C	Solutions 1, 2, 3, and 4
6	Control Vector GFP; 100 <i>E. coli</i> (Vial 6, colorless cap)	→ Briefly centrifuge down the contents of the vial → Reconstitute the lyophilizate (50 µg of Control Vector GFP) with 50 µl of sterile DNase- and RNase-free water → The solution is stable at -15 to -25°C	GFP control reaction, see page 18

All reconstituted solutions should be clear, with the exception of the *E. coli* lysate, which remains cloudy.

## 2. Prepare the working solution according to Table 3.

Table 3. Working solution

Solution	Contents	Preparation of working solution for one 50 $\mu$ l reaction	For use in
7	Reaction Solution	<p>Into one of the reaction tubes supplied, pipet the following components:</p> <ul style="list-style-type: none"> <li>→ 12 <math>\mu</math>l <i>E. coli</i> Lysate</li> <li>→ 10 <math>\mu</math>l Reaction Mix</li> <li>→ 12 <math>\mu</math>l Amino Acids</li> <li>→ 1 <math>\mu</math>l Methionine</li> <li>→ 5 <math>\mu</math>l Reconstitution Buffer</li> <li>→ 10 <math>\mu</math>l DNA in water or TE buffer</li> </ul> <p>Template amount:</p> <ul style="list-style-type: none"> <li>→ 0.5 <math>\mu</math>g circular template</li> <li>→ 0.5 <math>\mu</math>g linear template derived from one-step or two-step PCR</li> <li>→ 0.5 <math>\mu</math>g linear template derived from two-step PCR</li> <li>→ A premix of solutions 1–5 without DNA is recommended for multiple parallel reactions</li> <li>→ Mix carefully by rolling or gentle shaking; do not vortex!</li> </ul>	Running an experiment, see page 18

## Running an experiment, standard reaction

1. Close the reaction tubes, or the modules of the microplate, with the supplied caps or adhesive film, respectively.
2. Place the loaded reaction tubes into the Eppendorf® ThermoMixer® C (or Eppendorf® Thermomixer Comfort), an incubator or a water bath at 30°C.  
Note: If using modules, place the loaded modules within the frame of the microplate into the Eppendorf® ThermoMixer® C (or Eppendorf® Thermomixer Comfort), or an incubator at 30°C. A water bath is unsuitable for modules.
3. After 4–6 hours remove the reactions from the Eppendorf® ThermoMixer® C (or Eppendorf® Thermomixer Comfort), incubator or water bath.
4. Store Reaction Solution at –15°C to –25°C or at 0–4°C until purification or further processing.
5. See ‘Supporting information’, page 23 for sample preparation for SDS-PAGE.

### Points to consider

Temperature: The optimal temperature for most proteins is 30°C. However, lower temperatures may be used for proteins that tend to aggregate.

Time: Protein synthesis continues for up to 6 hours. However, the reaction is 90% complete after 4 hours.

## GFP control reaction

1. Reconstitute bottles 1 to 4 and bottle 6 according to Table 1 (kit format 6 reactions), or Table 2 (kit format 24 or 96 reactions) pages 15 and 16.
2. Prepare Reaction Solution in one of the reaction tubes according to Table 3, page 17.
3. Add 1 µg (1 µl) of reconstituted Control Vector GFP (vial 6) and 9 µl water or TE-buffer.
4. Start the reaction: Temperature: 30°C; Time: 6 hours.
5. After the run, store the Reaction Solution in a 2 ml vial for 24 hours at 2–8°C for maturation of GFP.
6. The fluorescence of GFP (excitation wavelength 395 nm, emission wavelength 504 nm) can be observed using an UV lamp (360 nm).
7. Apply 2–5 µl of the reaction onto SDS-polyacrylamide gels.
8. Run the gel, and then stain with Coomassie® Blue.  
Note: The GFP protein can also be detected on Western blots by using an anti-His<sub>6</sub> antibody.

### Oxidation of GFP

GFP needs molecular oxygen to form the fluorophore post-translationally (5). The yield of properly-folded fluorescent GFP is further increased by storing the Reaction Solution after the expression for 24 hours at 2–8°C.

After 24 hours of storing at 2–8°C the maturation is almost completed.

## Solubility of GFP

A remarkable fraction of the synthesized GFP will be in the pellet fraction. This is caused by the nature of the GFP gene used in the control reaction. GFP expression and detection via UV are for qualitative purpose only and should not be used to quantify the expression yield.

## Radioactive labeling

1. Reconstitute bottles 1 to 4 and vial 6 according to Table 1 (kit format 6 reactions), or Table 2 (kit format 24 or 96 reactions), pages 15 and 16.
2. Dilute 10  $\mu\text{l}$  of the reconstituted Methionine solution (Table 1 or Table 2) with 990  $\mu\text{l}$  of nuclease-free water to yield a 1 mM Methionine solution.
3. For one radioactive reaction prepare the following Reaction Solution in one of the reaction tubes:
  - 6  $\mu\text{l}$  *E. coli* Lysate
  - 5  $\mu\text{l}$  Reaction Mix
  - 6  $\mu\text{l}$  Amino Acids
  - 1.25  $\mu\text{l}$  1 mM Methionine solution (see Note)
  - 2  $\mu\text{l}$  of a L- $^{35}\text{S}$ Methionine 15 mCi/ml
  - 2.5  $\mu\text{l}$  Reconstitution Buffer
  - 0.25  $\mu\text{g}$  of plasmid DNA or 0.25  $\mu\text{g}$  of linear template in 2.25  $\mu\text{l}$  of water or TE buffer

Note: Addition of unlabeled Methionine to the labeling reaction is required to prevent premature termination for larger proteins or proteins with many Methionine residues.
4. Start the reaction: Temperature: 30°C; Time: Stop the reaction after 60 min.

Note: The Methionine added in this reaction is used up after 1 hour. For most applications the yield achieved with this protocol is sufficient. If more radioactive protein is required, please adapt conditions.
5. Apply 2–5  $\mu\text{l}$  of the reaction samples onto SDS-polyacrylamide gels.

Note: For optimum results precipitate the proteins before applying onto SDS-polyacrylamide gels (see 'Supporting information', page 23).
6. After the separation, dry the gel and apply for autoradiography (3–20 hours exposure).

### Protocol 3: Addition of supplements

Depending on the nature of the target protein, it may be necessary to add chemicals (e.g. detergents) or supplements (e.g. chaperones) in order to produce functional protein (examples in Table 4.). The design of this kit allows the addition of up to 10 µl of supplement solution to the reaction solution.

For efficient optimization of reaction conditions biotechrabbit recommends to study the effects (yield, solubility) of additional chemicals or supplements in the RTS 100 *E. coli* HY Kit. If a positive effect is observed in the 50 µl reaction volume of the RTS 100 *E. coli* HY Kit, the conditions can be applied in the 1 ml reaction volume of the RTS 500 ProteoMaster *E. coli* HY Kit.

Please also note that alternatively, aggregated proteins can often be reconstituted using reagents exemplarily shown in Table 4, because they are salted out in native conformation in contrast to the formation of inclusion bodies containing denatured protein after overexpression *in vivo*. In these cases, after cell-free protein synthesis precipitated protein should be pelleted by centrifugation (e.g. for 10 min at 16,000x g). This will also partially purify the target protein. The protein pellet should then be reconstituted in the reagent by strong shaking for several hours, while avoiding the generation of foam.

Table 4. Examples of chemicals and supplements to improve solubility and yield

Chemical/Supplement	Suggested concentration in reaction and feeding solution	Used for
Brij35	0.05 – 0.2%	membrane proteins, hydrophobic and other proteins
Brij78	0.05 - 0.2%	
DDM (n-Dodecyl β-D-maltoside)	0.05 - 0.2%	
CHAPS	10 mM	
Sorbitol	0.5 M	non-membrane proteins
DnaK supplement (cat. no. BR1401601)	please refer to RTS DnaK Supplement Short Instruction	
GroE supplement (cat. no. BR1401701)	please refer to RTS GroE Supplement Short Instruction	

## Supporting information

### Short protocol: Reconstitution of reaction components and running a reaction

Table 6. Short protocol steps, kit format for **6 reactions** (BR1400106)

Step	Contents	Reconstitution procedure
1	<i>E. coli</i> Lysate; 100 <i>E. coli</i> (6 rxns) (Vial 1, red cap)	Reconstitute the lyophilizate with 90 $\mu$ l Reconstitution Buffer. Do not vortex!
2	Reaction Mix; 100 <i>E. coli</i> (6 rxns) (Vial 2, green cap)	Reconstitute the lyophilizate with 75 $\mu$ l Reconstitution Buffer
3	Amino Acids w/o Meth.; 100 <i>E. coli</i> (6 rxns) (Vial 3, brown cap)	Reconstitute the lyophilizate with 90 $\mu$ l Reconstitution Buffer
4	Methionine; 100 <i>E. coli</i> (6 rxns) (Vial 4, yellow cap)	Reconstitute the lyophilizate with 82.5 $\mu$ l Reconstitution Buffer
5	Control Vector GFP; 100 <i>E. coli</i> (6 rxns) (Vial 6, colorless cap)	→ Briefly centrifuge down the contents of the vial → Reconstitute the lyophilizate with 15 $\mu$ l of sterile DNase- and RNase-free water
6	Reaction Solution	Mix the following components: → 12 $\mu$ l <i>E. coli</i> Lysate → 10 $\mu$ l Reaction Mix → 12 $\mu$ l Amino Acids → 1 $\mu$ l Methionine → 5 $\mu$ l Reconstitution Buffer → 10 $\mu$ l DNA in water or TE-buffer (0.5 $\mu$ g circular or linear template)
7	Start the reaction	Incubate for 4–6 hours at 30°C

Table 6. Short protocol steps, kit formats for **24 and 96 reactions** (BR1400101 and BR1400102)

Step	Contents	Reconstitution procedure
1	<i>E. coli</i> Lysate; 100 <i>E. coli</i> (Bottle 1, red cap)	Reconstitute the lyophilizate with 0.36 ml Reconstitution Buffer. Do not vortex!
2	Reaction Mix; 100 <i>E. coli</i> (Bottle 2, green cap)	Reconstitute the lyophilizate with 0.30 ml Reconstitution Buffer
3	Amino Acids; 100 <i>E. coli</i> (Bottle 3, brown cap)	Reconstitute the lyophilizate with 0.36 ml Reconstitution Buffer
4	Methionine; 100 <i>E. coli</i> (Bottle 4, yellow cap)	Reconstitute the lyophilizate with 0.33 ml Reconstitution Buffer
5	Control Vector GFP; 100 <i>E. coli</i> (Vial 6, colorless cap)	<ul style="list-style-type: none"> <li>→ Briefly centrifuge down the contents of the bottle</li> <li>→ Reconstitute the lyophilizate with 50 µl of sterile DNase- and RNase- free water</li> </ul>
6	Reaction Solution	Mix the following components: <ul style="list-style-type: none"> <li>→ 12 µl <i>E. coli</i> Lysate</li> <li>→ 10 µl Reaction Mix</li> <li>→ 12 µl Amino Acids</li> <li>→ 1 µl Methionine</li> <li>→ 5 µl Reconstitution Buffer</li> <li>→ 10 µl DNA in water or TE-buffer (0.5 µg circular or linear template)</li> </ul>
7	Start the reaction	Incubate for 4–6 hours at 30°C

## Optimizing SDS-polyacrylamide gel electrophoresis (SDS-PAGE)

### Precipitation of proteins prior to SDS-PAGE

The Reaction Solution contains polymers that interfere with the separation of proteins with an apparent molecular weight between 20 and 30 kDa. biotechrabbit recommends that samples are precipitated with acetone prior to the addition of SDS-PAGE sample buffer.

#### Procedure

1. To a 5  $\mu$ l sample of Reaction Solution add 50  $\mu$ l of  $-20^{\circ}\text{C}$  cold acetone, mix and incubate on ice for 5 min.
2. Centrifuge for 5 min at 10,000 rpm.
3. Discard the supernatant and air dry the pellet for 10 min (Speedvac may be used).
4. Dissolve the pellet in 20  $\mu$ l of SDS-PAGE sample buffer, heat for 5 min at  $95^{\circ}\text{C}$  and apply 5–20  $\mu$ l onto a SDS gel.

To distinguish between the soluble and pellet fractions of the protein of interest, add the following steps:

1. Centrifuge a 5  $\mu$ l sample of the Reaction Solution for 5 min at 10,000 rpm.
2. Pipet the supernatant into a separate tube.
3. Add 50  $\mu$ l ice-cold acetone to each tube and proceed as described above.

### Typical results

#### Yield from various template generation methods

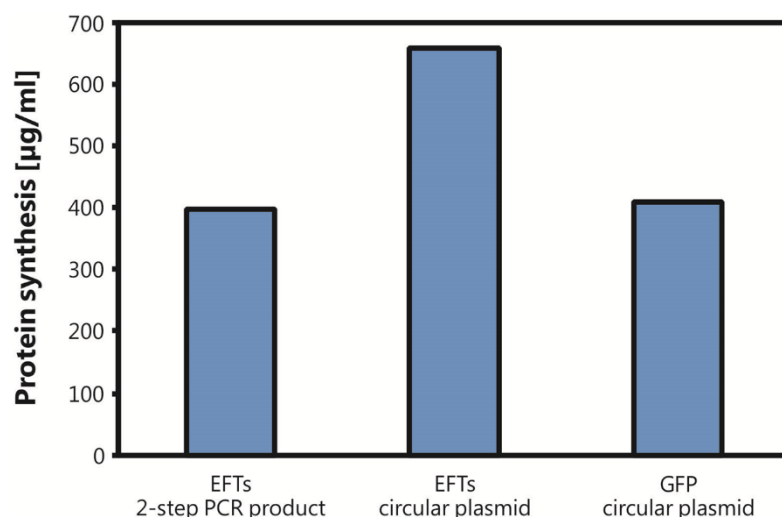


Figure 4. An EFTs coding DNA template (generated by 2-step PCR or by cloning into the pIVEX 2.3 expression plasmid) and a GFP coding vector (pIVEX 2.3) were used for the *in vitro* protein synthesis reactions. The 2-step PCR product was generated using the RTS Linear Template Kit Plus with C-terminal His<sub>6</sub>-Tag primers for the second PCR and the following primers for the first PCR:

Sense primer: 5' GAAGGAGATAAACAATGGCTGAAATTACCGCATC 3'

Antisense primer: 5' TGGTGATGGTGGTGACCCCAAGACTGCTTGGACATCGCA 3'

Up to 400  $\mu\text{g}/\text{ml}$  of protein from 2-step PCR, and up to 650  $\mu\text{g}/\text{ml}$  from plasmid template were obtained in a 4-hour reaction.

## Application: Optimization of protein expression using the RTS workflow

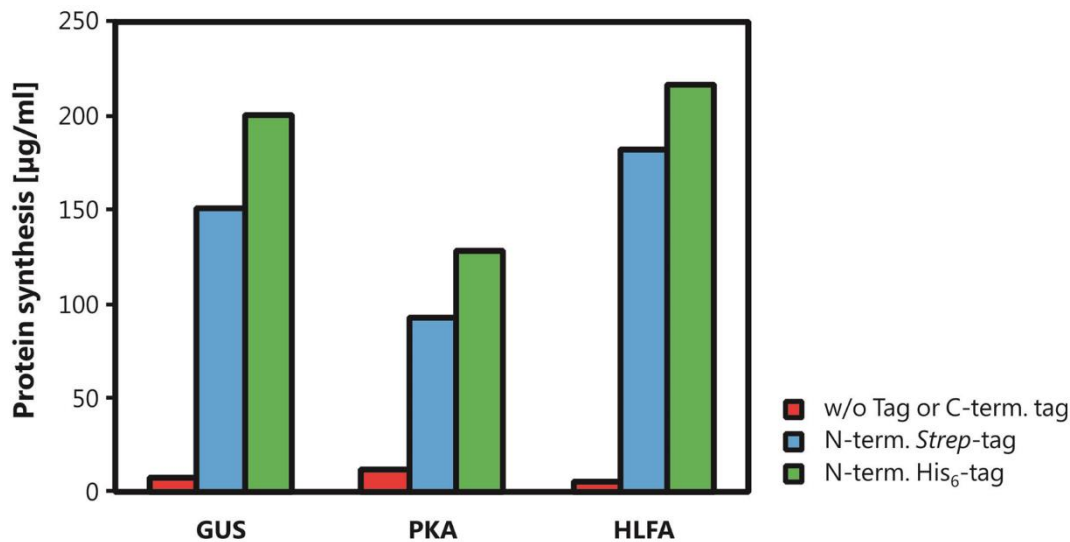


Figure 5. Optimization of the expression of human proteins  $\beta$ -Glucuronidase, Protein kinase A and liver Fatty acid binding protein.

Linear templates were generated with the RTS Linear Template Kit Plus. The proteins were expressed with the RTS 100 *E. coli* HY Kit. The figure shows protein yields of wild type and optimized gene variants. Both, N-terminal *Strep*-tag and N-terminal His<sub>6</sub>-tag overcome limited protein synthesis because of mRNA secondary structures, which impede translation initiation of the wild type genes.

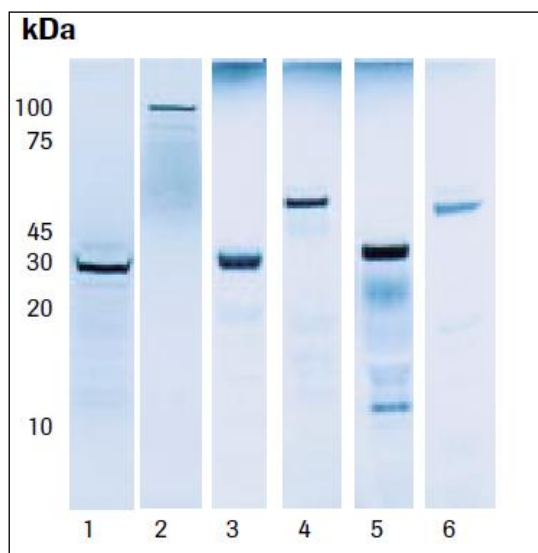
Application: Radioactive labeling with L-[<sup>35</sup>S] Methionine

Figure 6. Autoradiography of *in vitro* synthesized proteins from different species labeled with L-[<sup>35</sup>S] Methionine.

**Lane 1:** GFP (*A. victoria*); **Lane 2:**  $\beta$ -Galactosidase (*E. coli*); **Lane 3:** Fc gamma receptor (human); **Lane 4:** Rec. Plasminogen activator (human); **Lane 5:** Single chain antibody fragment (mouse); **Lane 6:** Rhodanese (bovine).

The indicated genes were cloned into pIVEX2.3 vectors and used for *in vitro* expression without further purification. After 30 minutes at 30°C, 2  $\mu$ l samples were separated on 10% SDS polyacrylamide gels. After drying, the gels were applied to Kodak X-OMAT AR film for autoradiography (3-hour exposure time).

## References

1. **DeVries, J.K. and G. Zubay.** 1967. Proc Natl Acad Sci U S A. 57, 1010.
2. **Zubay, G.** 1973. Annu. Rev. Genet. 7:267.
3. **Kim, D.M. and J.R. Swartz.** 2000. Biotechnol Lett 22:1537.
4. **Kim, D.M. and J.R. Swartz.** 2001. Biotechnol Bioeng 74:309.
5. **Coxon, A. and T.H. Bestor.** 1995. Chem Biol 2:119.
6. **Ahmed, A.K., et al.** 1975. J Biol Chem 250:8477.
7. **Odorzinsky, T.W. and A.Light.** 1979. J Biol Chem 254:4291.
8. **Rudolph, R. et al.** 1997. In: "Protein Function – A Practical Approach" Creighton, T.E. ed. Oxford University Press Inc. New York, pp 57.
9. **Sambrook J., et al.** 1989. "Molecular Cloning – A Laboratory Manual" Second Edition, Cold Spring Harbor Laboratory Press, New York.
10. **Ausubel, U.K., et al.** 1993. "Current Protocols In Molecular Biology" John Wiley & Sons Inc., New York.
11. **Abe, R., K. Shiraga, S. Ebisu, H. Takagi, and T. Hohsaka.** 2010. Incorporation of fluorescent non-natural amino acids into N-terminal tag of proteins in cell-free translation and its dependence on position and neighboring codons. J Biosci Bioeng 110:32-8.
12. **Babel, I., R. Barderas, A. Pelaez-Garcia, and J. I. Casal.** 2011. Antibodies on demand: a fast method for the production of human scFvs with minimal amounts of antigen. BMC Biotechnol 11:61.
13. **Beare, P. A., C. Chen, T. Bouman, J. Pablo, B. Unal, D. C. Cockrell, W. C. Brown, K. D. Barbian, S. F. Porcella, J. E. Samuel, P. L. Felgner, and R. A. Heinzen.** 2008. Candidate antigens for Q fever serodiagnosis revealed by immunoscreening of a Coxiella burnetii protein microarray. Clin Vaccine Immunol 15:1771-9.
14. **Berrier, C., I. Guilvout, N. Bayan, K. H. Park, A. Mesneau, M. Chami, A. P. Pugsley, and A. Ghazi.** 2011. Coupled cell-free synthesis and lipid vesicle insertion of a functional oligomeric channel MscL MscL does not need the insertase YidC for insertion in vitro. Biochim Biophys Acta 1808:41-6.
15. **Bertschinger, J., D. Grabulovski, and D. Neri.** 2007. Selection of single domain binding proteins by covalent DNA display. Protein Eng Des Sel 20:57-68.
16. **Blesneac, I., S. Ravaud, C. Juillan-Binard, L. A. Barret, M. Zoonens, A. Polidori, B. Miroux, B. Pucci, and E. Pebay-Peyroula.** 2012. Production of UCP1 a membrane protein from the inner mitochondrial membrane using the cell free expression system in the presence of a fluorinated surfactant. Biochim Biophys Acta 1818:798-805.
17. **Deniaud, A., L. Liguori, I. Blesneac, J. L. Lenormand, and E. Pebay-Peyroula.** 2010. Crystallization of the membrane protein hVDAC1 produced in cell-free system. Biochim Biophys Acta 1798:1540-6.
18. **Devenish, S. R. A., M. Kaltenbach, M. Fischlechner, and F. Hollfelde.** 2012. Droplets as Reaction Compartments for Protein Nanotechnology. Methods in Molecular Biology, Protein Nanotechnology: Protocols, Instrumentation, and Applications, Second Edition 996:269-286.
19. **Dortay, H., S. M. Schmockel, J. Fettke, and B. Mueller-Roeber.** 2011. Expression of human c-reactive protein in different systems and its purification from Leishmania tarentolae. Protein Expr Purif 78:55-60.
20. **Driguez, P., D. L. Doolan, A. Loukas, P. L. Felgner, and D. P. McManus.** 2010. Schistosomiasis vaccine discovery using immunomics. Parasites & Vectors 3:1-5.
21. **Elbaz, Y., T. Danieli, B. I. Kanner, and S. Schuldiner.** 2010. Expression of neurotransmitter transporters for structural and biochemical studies. Protein Expr Purif 73:152-60.
22. **Fujita, E., H. Dai, Y. Tanabe, Y. Zhiling, T. Yamagata, T. Miyakawa, M. Tanokura, M. Y. Momoi, and T. Momoi.** 2010. Autism spectrum disorder is related to endoplasmic reticulum stress induced by mutations in the synaptic cell adhesion molecule, CADM1. Cell Death Dis 1:e47.
23. **Goda, K., H. Ito, T. Kondo, and T. Oyama.** 2012. Fluorescence correlation spectroscopy to monitor Kai protein-based circadian oscillations in real time. J Biol Chem 287:3241-8.

24. **Helmy, O. M., M. M. M. Hussein, F. E. Murad, and H. A. Shoeb.** 2010. The preparation of 6x His-tagged granulocyte colony stimulating factor using an improved in vitro expression. *African Journal of Biotechnology* **9**:8566-8577.
25. **Ishido, T., N. Yamazaki, M. Ishikawa, and K. Hirano.** 2011. Characterization of DNA polymerase beta from *Danio rerio* by overexpression in *E. coli* using the in vivo/in vitro compatible pIVEX plasmid. *Microb Cell Fact* **10**:84.
26. **Jha, S. S., and A. A. Komar.** 2012. Using SecM arrest sequence as a tool to isolate ribosome bound polypeptides. *J Vis Exp* **19**.
27. **Kalantari-Dehaghi, M., G. J. Anhalt, M. J. Camilleri, A. I. Chernyavsky, S. Chun, P. L. Felgner, A. Jasinskas, K. M. Leiferman, L. Liang, S. Marchenko, R. Nakajima-Sasaki, M. R. Pittelkow, J. J. Zone, and S. A. Grando.** 2013. Pemphigus vulgaris autoantibody profiling by proteomic technique. *PLoS One* **8**:e57587.
28. **Karcher, D., and R. Bock.** 2009. Identification of the chloroplast adenosine-to-inosine tRNA editing enzyme. *Rna* **15**:1251-7.
29. **Kimelman, A., A. Levy, H. Sberro, S. Kidron, A. Leavitt, G. Amitai, D. R. Yoder-Himes, O. Wurtzel, Y. Zhu, E. M. Rubin, and R. Sorek.** 2013. A vast collection of microbial genes that are toxic to bacteria. *Genome Res* **22**:802-9.
30. **Krajicek, B. J., T. J. Kottom, L. Villegas, and A. H. Limper.** 2010. Characterization of the PcCdc42 small G protein from *Pneumocystis carinii*, which interacts with the PcSte20 life cycle regulatory kinase. *Am J Physiol Lung Cell Mol Physiol* **298**:L252-60.
31. **Kumar, G., and G. Chernaya.** 2009. Cell-free protein synthesis using multiply-primed rolling circle amplification products. *Biotechniques* **47**:637-9.
32. **Langlais, C., B. Guillaume, N. Wermke, T. Scheuermann, L. Ebert, J. LaBaer, and B. Korn.** 2007. A systematic approach for testing expression of human full-length proteins in cell-free expression systems. *BMC Biotechnol* **7**:64.
33. **Li, L., S. Munir, J. P. Bannantine, S. Sreevatsan, S. Kanjilal, and V. Kapur.** 2007. Rapid expression of *Mycobacterium avium* subsp. *paratuberculosis* recombinant proteins for antigen discovery. *Clin Vaccine Immunol* **14**:102-5.
34. **Llorens, J. V., J. B. Clark, I. Martinez-Garay, S. Soriano, R. de Frutos, and M. J. Martinez-Sebastian.** 2008. Gypsy endogenous retrovirus maintains potential infectivity in several species of *Drosophilids*. *BMC Evol Biol* **8**:302.
35. **Miyoshi, S., Y. Abe, M. Senoh, T. Mizuno, Y. Maehara, and H. Nakao.** 2011. Inactivation of *Vibrio vulnificus* hemolysin through mutation of the N- or C-terminus of the lectin-like domain. *Toxicon* **57**:904-8.
36. **Park, K. H., E. Billon-Denis, T. Dahmane, F. Lebaupain, B. Pucci, C. Breyton, and F. Zito.** 2011. In the cauldron of cell-free synthesis of membrane proteins: playing with new surfactants. *N Biotechnol* **28**:255-61.
37. **Rosenblum, G., C. Chen, J. Kaur, X. Cui, Y. E. Goldman, and B. S. Cooperman.** 2012. Real-time assay for testing components of protein synthesis. *Nucleic Acids Res* **40**:e88.
38. **Schaerli, Y., V. Stein, M. M. Spiering, S. J. Benkovic, C. Abell, and F. Hollfelder.** 2010. Isothermal DNA amplification using the T4 replisome: circular nicking endonuclease-dependent amplification and primase-based whole-genome amplification. *Nucleic Acids Res* **38**:e201.
39. **Singh, D., S. J. Chang, P. H. Lin, O. V. Averina, V. R. Kaberdin, and S. Lin-Chao.** 2009. Regulation of ribonuclease E activity by the L4 ribosomal protein of *Escherichia coli*. *Proc Natl Acad Sci U S A* **106**:864-9.
40. **Stein, V., and F. Hollfelder.** 2009. An efficient method to assemble linear DNA templates for in vitro screening and selection systems. *Nucleic Acids Res* **37**:e122.
41. **Stoevesandt, O., M. Vetter, D. Kastelic, E. A. Palmer, M. He, and M. J. Taussig.** 2011. Cell free expression put on the spot: advances in repeatable protein arraying from DNA (DAPA). *N Biotechnol* **28**:282-90.
42. **Suga, K., T. Tanabe, H. Tomita, T. Shimanouchi, and H. Umakoshi.** 2011. Conformational change of single-stranded RNAs induced by liposome binding. *Nucleic Acids Res* **39**:8891-900.

43. **Tian, J., H. Gong, N. Sheng, X. Zhou, E. Gulari, X. Gao, and G. Church.** 2004. Accurate multiplex gene synthesis from programmable DNA microchips. *Nature* **432**:1050-4.
44. **Umakoshi, H., K. Suga, H. T. Bui, M. Nishida, T. Shimanouchi, and R. Kuboi.** 2009. Charged liposome affects the translation and folding steps of in vitro expression of green fluorescent protein. *J Biosci Bioeng* **108**:450-4.
45. **Umakoshi, H., T. Tanabe, K. Suga, H. T. Bui, T. Shimanouchi, and R. Kuboi.** 2011. Oxidative stress can affect the gene silencing effect of DOTAP liposome in an in vitro translation system. *Int J Biol Sci* **7**:253-60.
46. **Wang, Y., and Y. H. Zhang.** 2009. Cell-free protein synthesis energized by slowly-metabolized maltodextrin. *BMC Biotechnol* **9**:58.
47. **Wu, N., J. G. Oakeshott, C. J. Easton, T. S. Peat, R. Surjadi, and Y. Zhu.** 2011. A double-emulsion microfluidic platform for in vitro green fluorescent protein expression. *J Micromech Microeng* **21**:1-7.
48. **Yang, Z., R. Ma, L. Huang, X. Zhu, J. Sheng, J. Cai, W. Hu, and Z. Xu.** 2012. High-level production of soluble adenine nucleotide translocator in *E. coli* cell-free system. *Process Biochem* **47**:395-400.
49. **Zarate, X., D. C. Henderson, K. C. Phillips, A. D. Lake, and D. W. Galbraith.** 2010. Development of high-yield autofluorescent protein microarrays using hybrid cell-free expression with combined *Escherichia coli* S30 and wheat germ extracts. *Proteome Sci* **8**:32.
50. **Zhang, D., G. Liu, J. Xue, J. Lou, K. H. Nierhaus, W. Gong, and Y. Qin.** 2012. Common chaperone activity in the G-domain of trGTPase protects L11-L12 interaction on the ribosome. *Nucleic Acids Res* **40**:10851-65.
51. **Zichel, R., A. Mimran, A. Keren, A. Barnea, I. Steinberger-Levy, D. Marcus, A. Turgeman, and S. Reuveny.** 2010. Efficacy of a potential trivalent vaccine based on Hc fragments of botulinum toxins A, B, and E produced in a cell-free expression system. *Clin Vaccine Immunol* **17**:784-92.

## Troubleshooting guide

The following troubleshooting recommendations are designed to address unexpected or undesired results. To ensure optimal use, follow the guidelines and recommendations in the manual.

### The control reaction

Observation	No control protein visible
Possible cause	Kit expired
Resolving	Order a new kit.
Possible cause	Kit has not been stored at $-15$ to $-25^{\circ}\text{C}$
Resolving	Order a new kit.
Possible cause	Contamination with RNases
Resolving	Repeat experiment and ensure to work RNase-free at every step.

### Expression of the target protein

Observation	No protein band of the target protein, but normal expression of control protein
Possible cause	Protein concentration too low
Resolving	Load the maximum amount of sample on the gel.
Possible cause	Protein is insoluble
Resolving	Analyze the supernatant and pellet.

Possible cause	Tag is hidden
Resolving	Try different tag position or protein-specific antibody. If the protein is His <sub>6</sub> -tagged, try a different anti-histidine antibody.
Resolving	Check protein on a Coomassie-stained gel in addition to the western blot.
Possible cause	Cloning error
Resolving	Check the sequence.
Possible cause	Low purity of plasmid DNA
Resolving	Ensure that the absorbance ratio 260 nm/280 nm is at least 1.7.
Resolving	Perform a phenol extraction.
Resolving	Make a new plasmid preparation.
Possible cause	Contamination with RNases
Resolving	Repeat experiment and ensure to work RNase-free at every step.
Possible cause	No initiation of translation due to strong secondary structures of the mRNA
Resolving	Introduce different N- or C-terminal tags using the RTS Linear Template Kit Plus.
Possible cause	Expressed protein interferes with the translation or transcription process
Resolving	Express the protein of interest together with control protein. If control protein expression is inhibited, the active protein cannot be expressed using RTS.

---

Observation	Low expression yield
Possible cause	Expression time too short
Resolving	Extend expression time.
Possible cause	The tag has a negative influence on the folding of the protein
Resolving	Introduce different N- or C-terminal tags using the RTS Linear Template Kit Plus.
Possible cause	Amount of template DNA not optimal
Resolving	Vary (increase) DNA concentration to get optimum results.
Observation	Sufficient protein expression, but low yield of active protein
Possible cause	Incorrect folding of the protein due to: <ul style="list-style-type: none"> <li>→ dependence on co-factors</li> <li>→ necessity of disulfide bonds</li> <li>→ dependence on post-translational modifications</li> <li>→ dependence on chaperones</li> </ul>
Resolving	Add necessary co-factors.
Resolving	Proteins with more than a few disulfide bonds (it could be shown, that three disulfide bonds can be formed) may be not expressed in a functional form in the RTS 100 <i>E. coli</i> HY system. Use RTS 100 <i>E. coli</i> Disulfide Kit.
Resolving	Add chaperones (8).
Resolving	For proteins with up to 3 disulfide bonds, allow oxidation after the reaction for the formation of disulfide bonds (2, 3).
Resolving	The <i>E. coli</i> lysate cannot introduce post-translational modifications such as glycosylation, phosphorylation, or signal sequence cleavage.
Resolving	Use RTS GroE Supplement for proteins in the range between 20 and 30 kDa.

Observation	Product appears in the pellet after centrifugation
Possible cause	Aggregation
Resolving	Add/adjust chaperones or cofactors.
Resolving	Change experimental conditions (time, temperature, e.g. lower in steps of 4–14°C).
Resolving	Add mild detergents (e.g. up to 0.2% Brij 35, or 10 mM CHAPS for membrane proteins, see also Table 4 on page 20 and reference 9).
Observation	Several product bands on SDS-PAGE or product smaller than expected
Possible cause	Proteolytic degradation
Resolving	Dissolve 1 tablet of cOmplete mini EDTA-free (Roche) in 0.5 ml nuclease-free water and add 2 µl for a 50 µl reaction.
Possible cause	Internal initiation site
Resolving	Check sequence for internal Methionine codons (ATG), look for RBS-like sequence (variants of `AAGGAG`) 5 to 12 nucleotides upstream of the ATG. Mutate the RBS-like sequence, if possible replace the corresponding Methionine by point mutation.
Possible cause	Premature termination of the translation
Resolving	Check the sequence of the target gene regarding reading frame and mutations which might yield a stop codon.
Resolving	Search for strong secondary structures of the mRNA and eliminate them by conservative mutations.

## 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