

1-Step CHO High-Yield IVT Trial Kit

88893

2595.0

Number	Description
88893	1-Step CHO High-Yield IVT Trial Kit , contains sufficient reagents to perform 2 reactions (100µL each)

Kit Contents	Cap Color	88893X
CHO Lysate	Silver seal	1 lyophilized vial
Accessory Proteins	Green	25µL
5X Reaction Mix	Yellow	40µL
4X Dialysis Buffer	Clear	700µL
Positive Control DNA: pCFE-GFP (0.5µg/µL, 10µg)	Clear	20µL
pT7CFE1-NHis-GST-CHA Expression Vector (0.5µg/µL, 10µg)	Clear	20µL

Kit Contents	88893Y
Microdialysis Device	2 each
Nuclease-free Water	50mL

Note: Completely read the instructions before proceeding with the protocols.

Storage: Upon receipt store 88893X at -80°C and 88893Y at room temperature. 88893X is shipped with dry ice. 88893Y is shipped at ambient temperature.

Table of Contents

Introduction	2
Procedure Summary.....	2
Important Product Information	2
Additional Materials Required.....	2
Protocol for using the 1-Step CHO High-Yield IVT Trial Kit	3
A. Protein Expression.....	3
B. Determination of Protein Expression Level	4
C. Purification of IVT-expressed Proteins	4
Troubleshooting	5
Additional Information	6
A. pT7CFE1-NHis-GST-CHA Vector Cloning Sites and Sequence Features	6
B. Vector DNA Clean-up and Concentration Protocol	6
C. Expression-ready Clones for Use with the 1-Step CHO High-Yield IVT Kits	6
Related Thermo Scientific Products	7
General References	7

Introduction

The Thermo Scientific™ 1-Step CHO High-Yield IVT Trial Kit is a mammalian *in vitro* translation (IVT) system based on CHO cell lysates. The kit contains all of the cellular components required for protein synthesis, including ribosomes, initiation factors, elongation factors and tRNA. When supplemented with the included proprietary Accessory Proteins, Reaction Mix and a DNA template cloned into the Thermo Scientific™ pT7CFE1-NHis-GST-CHA Vector, this system can synthesize protein for up to 16 hours.

The benefits of *in vitro* protein expression over traditional *in vivo* systems include the ability to express toxic proteins, synthesize proteins faster and label protein with modified amino acids. The optimized kit contains a T7 promoter and an EMCV internal ribosome entry site (IRES) to facilitate high levels of *in vitro* protein expression in a cap-independent fashion. Using a vector containing the EMCV IRES element is critical for obtaining high expression levels in this CHO *in vitro* protein expression system.

Procedure Summary

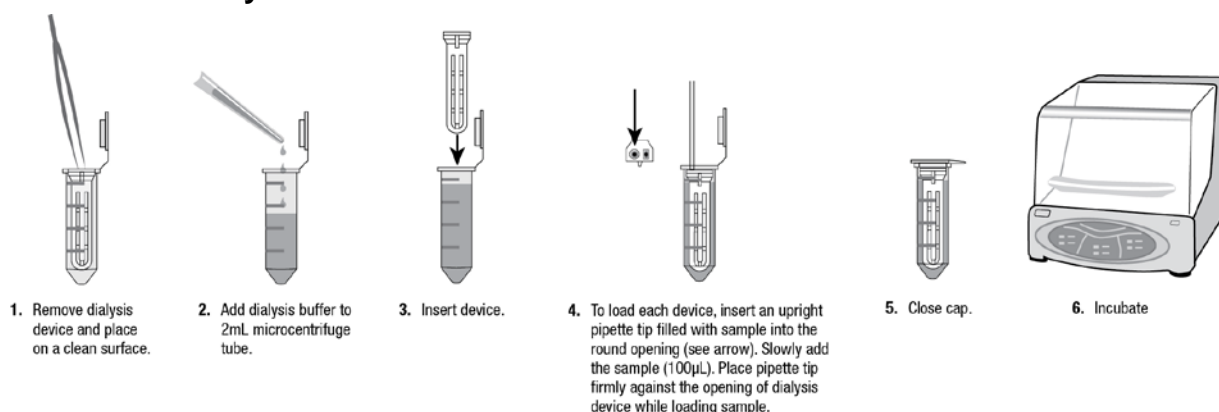


Figure 1. The Thermo Scientific 1-Step CHO High-Yield IVT Trial Kit protocol.

Important Product Information

- Use the included Thermo Scientific pT7CFE1-NHis-GST-CHA Vector (Product No. 88871) for cloning and expressing the target gene. See the Additional Information Section for additional vector choices, cloning sites and expression-ready clones.
- CHO Lysate is provided as a lyophilized lysate. Reconstitute contents on ice, as recommended in the Protocol Section. Aliquot if necessary and quickly store at -80°C. All components of the kit are stable for up to five freeze-thaw cycles as long as the contents are stored at -80°C immediately after use.
- Undiluted lysate and reactions containing lysate will appear cloudy before and after incubation. Accessory Proteins and Reaction Mix may also appear clear to cloudy upon thawing; mix thoroughly but gently before and after adding each component to the IVT reaction. Undiluted 4X Dialysis Buffer may appear cloudy; mix well before and after dispensing.
- Avoid RNase contamination by wearing gloves; working in a clean, dust-free environment; and using RNase-free tips and microcentrifuge tubes.

Additional Materials Required

- DNA preparation kit (e.g., Thermo Scientific™ GeneJET™ Plasmid Maxi Prep Kit, Product No. K0503 or K0492)
- Western immunoblot accessories for detecting expressed protein
- FITC filter-containing device to observe the expression of GFP in positive-control reactions
- 1.5mL RNase-free microcentrifuge tubes for assembling reactions
- RNase-free pipette tips
- Optional: Shaker/incubator capable of maintaining temperature at 30°C

Protocol for Using the 1-Step CHO High-Yield IVT Trial Kit

A. Protein Expression

- With the exception of 4X Dialysis Buffer, thaw all other reagents in the kit contents of 88893X and maintain on ice. Thaw 4X Dialysis Buffer at 25-30°C for a maximum of 15 minutes and, after making a 1X mixture, maintain the diluted buffer at room temperature.

Note: Store any unused 88893X kit components at -80°C.

Note: The 4X Dialysis Buffer may appear cloudy. Mix or vortex gently. Do not centrifuge before use. Once diluted, the 1X Dialysis Buffer will become clear within minutes.

- Remove dialysis device as shown in Step 1 of Procedure Summary. Combine room temperature 4X Dialysis Buffer and Nuclease-free Water (volumes per Table 1) in the provided 2mL microcentrifuge tube.

Table 1. Reconstitution of the Dialysis Buffer.

Component	μL
4X Dialysis Buffer	350
Nuclease-free Water	1050
Total	1400

- Place the dialysis device inside the tube containing 1X Dialysis Buffer as shown in the Procedure Summary Section. Confirm dialysis membrane is completely immersed. Incubate at 30°C for 30 minutes.
- Reconstitution of the lyophilized CHO Lysate vial:** Add 110 μ L of Nuclease-free Water provided in the kit to the lyophilized vial. Let it stand for 2-3 minutes and then slowly pipette up and down the contents of the tube at least 5 times to completely reconstitute the lysate. Store at -80°C any unused portion of the CHO Lysate.
- Prepare IVT reactions using Table 2. Add the reagents in the order listed into a 1.5mL RNase/DNase-free tube at room temperature. Mix the CHO Lysate with Accessory Proteins and incubate for 5-10 minutes at room temperature before adding the remaining components. Gently mix the reaction after each reagent addition.

Table 2. Components of the IVT reaction.

Component	GFP Control (μL)	Target Protein (μL)
CHO Lysate	50	50
Accessory Proteins	10	10
5X Reaction Mix	20	20
pCFE-GFP DNA (0.5 μ g/ μ L)	8	—
Cloned DNA (0.5 μ g/ μ L)	—	8
Nuclease-free Water	12	12
Total	100	100

- Transfer the supernatant into the empty dialysis device as described in the Procedure Summary Section.
- Close the lid of the microcentrifuge tube over the dialysis device. Incubate the reaction for 6-16 hours at 30°C.

Note: Although protein expression is complete within 6 hours for most proteins tested, incubating up to 16 hours may increase expression of some proteins. Optimal time to express each protein must be determined empirically.

Note: Shaking the dialysis device containing the reaction mixture at ~600-900rpm during incubation may increase total protein yields up to 30%. Use a small table-top shaker/incubator such as an Eppendorf™ Thermomixer™ Shaker.

- Resulting reactions may be stored on ice for same-day use. For long-term storage, transfer the reaction contents from the dialysis device and store separately in a 1.5mL microcentrifuge tube at -20°C or colder. Occasionally, a small white precipitate may be visible, which can be removed by centrifuging the reactions at 10,000 \times g for 2 minutes.
- Proteins expressed using this kit may be purified using the purification guidelines provided in the Product Blog article “Choosing a vector and purification method for *in vitro* protein expression” on our website at: thermoscientific.com/pierce.

B. Determination of Protein Expression Level

Note: The GFP control protein is from the copepod *Pontellina plumata*. This GFP is not reactive to antibodies generated against *Aequorea victoria* GFP (i.e., EGFP or other EGFP mutants).

Note: Use polyclonal antibodies to TurboGFP (Product No. PA5-22688).

1. Visualize or quantitate the GFP control protein using one of the following methods:

Quick visual detection: Place the GFP reaction tubes directly under a microscope or imaging equipment containing a FITC filter (ex/em: 482/502nm); alternatively, spot a small volume (1-2 μ L) on a piece of plastic wrap or laboratory film and visualize with fluorescent imaging equipment.

Fluorescent plate reader: Place sample directly into a white or black 96- or 384-well plate. Evaluate signal using a fluorescent plate reader at ex/em: 482/502nm. To quantitate GFP, compare the fluorescence to a recombinant GFP standard curve.

2. Visualize or quantitate non-fluorescent protein expression using one of the following methods:

Fast Western immunoblot analysis: This is a quick protocol consisting of transfer and detection of proteins separated on SDS-PAGE using ultra-sensitive Thermo Scientific™ SuperSignal™ Substrate. A detailed protocol and reagents required for Western blot detection can be found at thermoscientific.com/pierce; search using “fast Western blot.”

SDS-PAGE analysis: Separate proteins by SDS-PAGE and stain using Thermo Scientific™ GelCode™ Blue Stain Reagent (Product No. 24590), Thermo Scientific™ Imperial™ Protein Stain (Product No. 24615) or Thermo Scientific™ PageBlue™ Protein Staining Solution (Product No. 24620) (Figure 2).

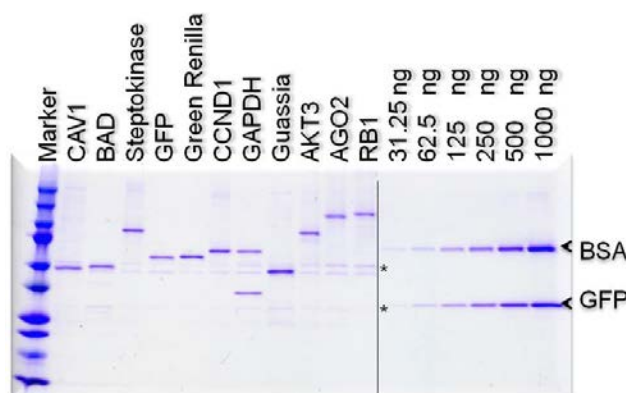


Figure 2. Purification of N-terminal GST fusion proteins with immobilized glutathione. Genes cloned into pT7CFE1-NHis-GST-CHA were used to express GST-fusion proteins for 6 hours following the procedure described above. Purification of GST-fused proteins was performed as described with 50mM glutathione using instructions provided in the Product Blog article “Choosing a vector and purification method for *in vitro* protein expression” on our website at thermoscientific.com/pierce. The additional bands denoted with a * and found with the purified proteins were previously identified as cellular proteins eEF1G and GSTM3 by mass spectrometry. These proteins are known to bind to the glutathione column and co-elute with GST-tagged proteins. Approximately 500ng of each of the purified proteins were separated by SDS-PAGE and stained using the Thermo Scientific™ Pierce™ Power Stainer (Product No. 22833).

C. Purification of IVT-expressed Proteins

Proteins expressed using this kit may be purified using the purification guidelines provided in the Product Blog article “Choosing a vector and purification method for *in vitro* protein expression” on our website at: thermoscientific.com/pierce.

Troubleshooting

Problem	Possible Cause	Solution
GFP not detected by fluorescence in positive control reaction	Incorrect filter set used	The excitation/emission wavelengths of GFP are 482/502nm
	Lysates became inactive	Store unused lysate in nuclease-free tubes at -80°C; do not exceed five cycles of freezing and thawing
No expression of target protein	Incorrect vector used	Use cloning vector pT7CFE1-NHis-GST-CHA provided in the kit to clone and express the gene of interest Note: The 1-Step CHO High-Yield IVT Kits are optimized using the pCFE1 vector and its derivatives; for a complete listing, please visit thermoscientific.com/pierce . See also Section C in Additional Information for the readily available clone collection
	CHO Lysate, Accessory Proteins and Reaction Mix were stored at a suboptimal temperature	Store unused CHO Lysate, Accessory Proteins and Reaction Mix in nuclease-free tubes at -80°C; do not exceed five cycles of freezing and thawing
	Poor-quality DNA	Ethanol precipitate the DNA to remove trace amounts of inhibitors or salts; see the Additional Information Section for the recommended protocol
	Degradation of mRNA in the translation reaction	Maintain an RNase-free environment by wearing gloves; working in a clean, dust-free environment; and using RNase-free tips and microcentrifuge tubes
	Protein was sensitive to proteases	Add Thermo Scientific™ Halt™ Protease Inhibitor Single-Use Cocktail, EDTA-free (100X) (Product No. 78425) at 0.5X to the reaction mix in Table 2
Low yield of target proteins	Incorrect incubation temperature	Perform reactions at 30°C
	Incorrect order of reagent addition	Incubate CHO Lysate with Accessory Proteins for 5-10 minutes before adding remaining components to improve target protein expression
Smaller band size than predicted	Stop codons were in genes of interest	Ensure the cloned genes do not have a stop codon in the open reading frame
Protein appears to be degraded	Proteins were susceptible to proteases	Add Halt Protease Inhibitor Single-Use Cocktail, EDTA-free (100X) (Product No. 78425) at 0.5X to the reaction mix in Table 2
Low protein yield after purification	Reaction scale was too small	Follow guidelines provided in the Product Blog article “Choosing a vector and purification method for <i>in vitro</i> protein expression” on our website at: thermoscientific.com/pierce
		Increase reaction size
Low protein yield after purification	Affinity tag was not accessible	Use different affinity purification for the tagged protein
		Purify protein under denaturing conditions (e.g., 8M urea) using the Thermo Scientific™ HisPur™ Cobalt Purification Kit (Product No. 90090)

Additional Information

A. pT7CFE1-NHis-GST-CHA Vector Cloning Sites and Sequence Features

The 1-Step CHO High-Yield IVT Kits have been optimized using the pT7CFE1-NHis-GST-CHA cloning vector, which is designed for high-level protein expression. In addition to multiple purification tags, it contains an HRV 3C cleavage site for tag removal. For a complete listing of pT7CFE1 expression vector derivatives, visit thermoscientific.com/pierce; search using “expression vectors.”

Features:

- 10 unique restriction sites are provided in the multiple cloning site for cloning genes of interest (Figure 1)
- 5' UTR consisting of EMCV internal ribosome entry site (IRES) required for high-level protein expression
- Poly A sequence in the 3' region promotes mRNA stabilization and protection from nucleases
- T7 terminator ensures synthesis of accurate-sized mRNA transcripts

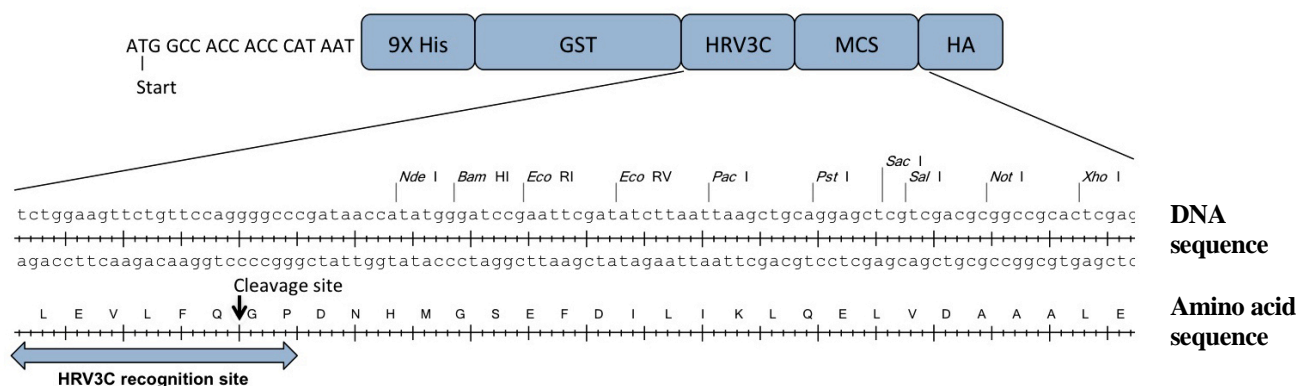


Figure 1. The Thermo Scientific pT7CFE1-NHis-GST-CHA Vector multiple cloning site, with the exception of Msc I, is common to all of the expression vectors used in the Thermo Scientific 1-Step High-Yield CHO IVT Kits. The translational start site is the ATG found upstream of the His tag region.

B. Vector DNA Clean-up and Concentration Protocol

Prepare DNA using a standard maxi- or mini-prep protocol. To avoid compromising protein expression yield, completely remove contaminating proteins and eliminate the RNase A used in many mini-prep protocols. Perform the following steps to precipitate and, subsequently, concentrate the DNA.

1. Add 1/10 volume of 3M sodium acetate, pH 5.5 and two volumes of ethanol. Thoroughly mix the reaction and incubate at -20°C for 15 minutes.
2. Centrifuge the mixture at 14,000 × g for 15 minutes. Remove the supernatant and wash the pellet once with 70% ethanol.
3. Centrifuge at 14,000 × g for 5 minutes. Using a fine tip, remove all of the supernatant, including the residual. Air-dry the pellet for 5 minutes at room temperature.
4. Resuspend the pellet in nuclease-free water before measuring the DNA concentration. DNA templates may be stored in a Tris-based buffer. It is not necessary to linearize the plasmid DNA before use.

C. Expression-ready Clones for Use with the 1-Step CHO High-Yield IVT Kits

- Custom cloning service; please visit thermoscientific.com/pierce and search for “cloning service.”
- The pANT7 vector library from the ASU Biodesign Institute DNASU Plasmid Repository is compatible with our 1-Step CHO High-Yield IVT Kit. Visit <http://dnasu.asu.edu/DNASU/Home.jsp> for information and ordering. Under advanced search options choose “pANT7” for vector selection.
- PCR templates: see Tech Tip #72: PCR protocol for generating optimized templates for Pierce Human *In Vitro* Expression Kits on our website.

Related Thermo Scientific Products

88894	1-Step CHO High-Yield Maxi IVT Kit
88859-71	pT7CFE1-based Expression Vectors
88899	Recombinant GFP Protein
88881-2	1-Step Human Coupled IVT Kits — DNA
88890-2	1-Step Human High Yield IVT Kits
MA4004	Mouse anti-Glutathione S-transferase Monoclonal Antibody (8-326)
MA121315	Mouse anti-6x-His Epitope Tag Monoclonal Antibody (HIS.H8)
26183	Mouse anti-HA Monoclonal Antibody (2-2.2.14)
MA4004	Mouse anti-Glutathione S-transferase Monoclonal Antibody (8-326)
PA5-22688	Anti-TurboGFP Polyclonal Antibody
88221	HisPur™ Ni-NTA Resin, see our website for all related products
89964	HisPur™ Cobalt Resin, see our website for all related products
16100	Pierce™ Glutathione Agarose, see our website for all related products
26182	Pierce™ Anti-HA Agarose, see our website for all related products
88836-7	Pierce™ Anti-HA Magnetic Beads
88821-2	Pierce™ Glutathione Magnetic Beads
88831-2	HisPur™ Ni-NTA Magnetic Beads
90410-1	Low Protein Binding Microcentrifuge Tubes, 2.0mL

General References

- Imataka, H., *et al.* (2009). Advantages of CHO cell-derived, cell-free protein synthesis systems (Japanese). *Seikagaku* **81**(4):303-7.
- Brödel, A.K., Sonnabend, A., Kubick, S. (2014) Cell-free protein expression based on extracts from CHO cells. *Biotechnol Bioeng* **111**(1):25-36.
- Kozak, M. (1983). Comparison of initiation of protein synthesis in prokaryotes, eukaryotes and organelles. *Microbiol Rev* **47**(1):1-45.
- Kozak, M. (2005). Regulation of translation via mRNA structure in prokaryotes and eukaryotes. *Gene* **361**:13-37.

Products are warranted to operate or perform substantially in conformance with published Product specifications in effect at the time of sale, as set forth in the Product documentation, specifications and/or accompanying package inserts ("Documentation"). No claim of suitability for use in applications regulated by FDA is made. The warranty provided herein is valid only when used by properly trained individuals. Unless otherwise stated in the Documentation, this warranty is limited to one year from date of shipment when the Product is subjected to normal, proper and intended usage. This warranty does not extend to anyone other than Buyer. Any model or sample furnished to Buyer is merely illustrative of the general type and quality of goods and does not represent that any Product will conform to such model or sample.

NO OTHER WARRANTIES, EXPRESS OR IMPLIED, ARE GRANTED, INCLUDING WITHOUT LIMITATION, IMPLIED WARRANTIES OF MERCHANTABILITY, FITNESS FOR ANY PARTICULAR PURPOSE, OR NON INFRINGEMENT. BUYER'S EXCLUSIVE REMEDY FOR NON-CONFORMING PRODUCTS DURING THE WARRANTY PERIOD IS LIMITED TO REPAIR, REPLACEMENT OF OR REFUND FOR THE NON-CONFORMING PRODUCT(S) AT SELLER'S SOLE OPTION. THERE IS NO OBLIGATION TO REPAIR, REPLACE OR REFUND FOR PRODUCTS AS THE RESULT OF (I) ACCIDENT, DISASTER OR EVENT OF FORCE MAJEURE, (II) MISUSE, FAULT OR NEGLIGENCE OF OR BY BUYER, (III) USE OF THE PRODUCTS IN A MANNER FOR WHICH THEY WERE NOT DESIGNED, OR (IV) IMPROPER STORAGE AND HANDLING OF THE PRODUCTS.

Unless otherwise expressly stated on the Product or in the documentation accompanying the Product, the Product is intended for research only and is not to be used for any other purpose, including without limitation, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses, or any type of consumption by or application to humans or animals.

Current product instructions are available at www.thermoscientific.com/pierce. For a faxed copy, call 800-874-3723 or contact your local distributor.

© 2014 Thermo Fisher Scientific Inc. All rights reserved. Eppendorf and Thermomixer are trademarks of Eppendorf AG. Unless otherwise indicated, all trademarks are property of Thermo Fisher Scientific Inc. and its subsidiaries. Printed in the USA.