

***In Vitro* Director™ PCR System – 5' Anchor™**

 Product Number **ID0040**
 Storage Temperature –20 °C

Product Information
TECHNICAL BULLETIN
Product Description

Sigma's *In Vitro* Director™ PCR System – 5' Anchor™ kit provides all reagents needed for the rapid generation of high quality, expression-ready DNA templates that can be used with *in vitro* transcription/translation systems to produce native proteins. *In vitro* synthesized proteins are very useful for rapid analysis of gene functions, such as enzymatic activity, protein-protein interaction, function domain mapping, and structure analysis.

The PCR system uses the Director method, a novel method that facilitates the directional ligation of transcription/translation regulatory sequences (Anchors) to the PCR products. It differs from conventional PCR kits in that it uses a specially formulated dNTP mix containing dATP α S and dGTP α S. After PCR, the cohesive 5' termini of the amplicon are produced by Exonuclease III digestion instead of being generated by traditional restriction enzyme digests. Incorporation of dA/GTP α S into the amplicon protects the amplicon from overdigestion by Exonuclease III.¹

The kit contains paired 5' and 3' Anchors as well as universal primers complementary to the Anchor sequences. The 5' Anchor is a double-stranded DNA containing the T7 promoter, Kozak sequence, and a *Hind* III overhang. The 3' Anchor is also a double-stranded DNA containing a *Bgl* II overhang and terminator sequence.

The universal process of rapidly generating expression-ready DNA templates involves five steps: a first round PCR using gene-specific primers, Exonuclease III digestion, PCR clean-up, ligation of the PCR amplicon to the 5' and 3' Anchors, and a second round of PCR using Anchor primers.

Reagents

The kit contains sufficient reagents to perform 25 reactions.

Kit Components	Product Number	Quantity
JumpStart™ REDAccuTaq™ LA DNA Polymerase	D 1938	2 vials (125 μ l)
10x AccuTaq™ LA DNA Polymerase Buffer	B 0174	0.5 ml
ExoClone™ dNTP Mix (20x)	E 4280	1vial (62.5 μ l)
Control PCR Template	D 4314	10 μ l (1 ng/ μ l)
Control RDC Primer-F	P 4986	25 μ l (20 μ M)
Control RDC Primer-R	P 5111	25 μ l (20 μ M)
Exonuclease III	E 1131	25 μ l (100U/ μ l)
dNTP Mix, 10 mM	D 7295	0.2 ml
Water Molecular Biology Reagent	W 4502	1.5 ml
5' Anchor™	A 7103	50 μ l (1 ng/ μ l)
3' Anchor™	A 7228	50 μ l (1 ng/ μ l)
Anchor™ Primer- Forward	A 7478	25 μ l (20 μ M)
Anchor™ Primer-Reverse	A 7353	25 μ l (20 μ M)

Reagents and Equipment Required but not Provided

- Template DNA and 5' phosphorylated gene-specific primers with *Hind* III/*Bgl* II overhangs
- Pipettes and aerosol resistant pipette tips
- 0.5 or 0.2 ml thin-walled PCR micro-centrifuge tubes
- Thermal cycler
- GenElute™ PCR Clean-up Kit (Product Code NA1020) or equivalent
- T4 DNA Ligase with 10x Buffer (Product Code D 1567)
- *In vitro* TNT® Wheat Germ Extract System (Product Code TN0100) or TNT T7 Quick for PCR DNA rabbit reticulocyte system (Promega)

Precautions and Disclaimer

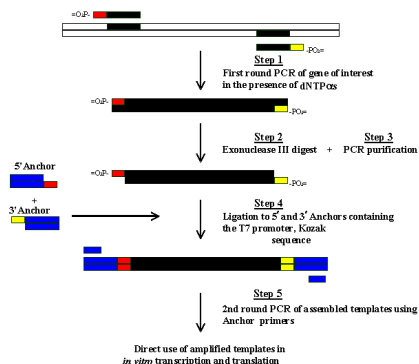
This kit is for laboratory use only, not for drug, household or other uses. Please consult the Material Safety Data Sheet for information regarding hazards and safe handling practices.

Storage/Stability

All components should be stored at -20°C . Avoid multiple freeze-thaw cycles.

Procedure

Outline



General Considerations

Template DNA

The kit is compatible with different sources of DNA template, including plasmid DNA, cDNA, and genomic DNA. However, PCR performance may be affected by the complexity and purity of the DNA template.

Primer Design

The gene specific primers require 5' phosphorylation and *Hind* III/*Bgl* II overhangs. The 5' forward primer requires the addition of 5 basepairs corresponding to the *Hind* III overhang (5'-P-AGCTT - target sequence - 3'). The 3' Reverse primer requires the addition of 5 basepairs corresponding to the *Bgl* II overhang (5'-P-GATCT - target sequence - 3'). Primers should be ≥ 20 bp (specific target sequence plus 5 bases for overhangs) to achieve adequate specificity; however, the length can vary depending on the sequence content of the selected region. Phosphorylation is necessary for efficient ligation and can be done during primer synthesis. The T_m , GC content of primers, and complexity of the template DNA are important considerations in successful PCR. We recommend the

T_m of your primers be between $65-75^{\circ}\text{C}$ with the GC content between 40-60%. The T_m difference of forward and reverse primers should not exceed $\pm 5^{\circ}\text{C}$.

DNA Polymerases

The kit uses JumpStart REDAccuTaq LA DNA polymerase for PCR amplification. This special enzyme blend offers several advantages over conventional Taq polymerases: excellent amplification for long-distance PCR (up to 22 kb for complex genomic DNA and 40 kb for less-complex templates); high fidelity (6.5x high than standard Taq polymerase and comparable to Pfu polymerase); hot-start mechanism using JumpStart Taq antibody to prevent nonspecific product formation, improving specificity; and direct loading of PCR products onto an agarose gel without the addition of loading buffers, saving time in product characterization. Other DNA polymerases are also compatible with this kit; however, additional PCR optimization may be required to obtain comparable results.

PCR Conditions

PCR conditions need to be determined empirically for each set of specific primers and template. It is critical to optimize PCR conditions for high specificity, i.e., generating a single product band, to avoid the need for further gel purification. A "hot-start" mechanism using JumpStart Taq antibody, which significantly improves PCR specificity, has been incorporated in this kit. For most applications, you should only need to adjust the annealing temperature and extension time according to the T_m of the primers and the length of the template. However, other techniques such as using "touchdown" PCR may further improve the specificity of the amplification.³

Step 1. First Round PCR

Amplify gene of interest using phosphorylated gene specific primers (5' forward primer = 5'-P-AGCTT - target sequence, 3' Reverse primer = 5'-P-GATCT - target sequence) and PCR system. Mix the following components in a 0.2 ml or a 0.5 ml PCR tube on ice.

Control reaction:

5 μl	10x AccuTaq LA DNA Polymerase Buffer
2.5 μl	ExoClone dNTP mix (20x) (teal cap)
1 μl	Control RDC Primer-F (20 μM)
1 μl	Control RDC Primer-R (20 μM)
1 μl	Control PCR template (1 ng/ μl)
2.5 μl	JumpStart REDAccuTaq LA DNA Polymerase
37 μl	Water
50 μl	Total volume

Sample reaction:

5 μ l	10x AccuTaq LA DNA Polymerase Buffer
2.5 μ l	ExoClone dNTP mix (20x) (teal cap)
1 μ l	Gene specific primer-forward (20 μ M)
1 μ l	Gene specific primer-reverse (20 μ M)
x μ l	Target sequence template*
2.5 μ l	JumpStart REDAccuTaq LA DNA Polymerase
38-x μ l	Water
50 μ l	Total volume

*Note: if using a simple template such as plasmid, 1 ng target template should be sufficient; if using a complex template, such as a cDNA mixture, 50-500 ng template is needed.

Control PCR:

Initial denaturation:	95 °C, 5 minutes
30 cycles:	
denaturation	94 °C, 30 seconds
annealing	62 °C, 30 seconds
extension	68 °C, 2 minutes
Final extension:	68 °C, 7 minutes
Soak:	4 °C

Sample PCR:

The PCR amplification profile will vary depending on the T_m value of the specific primer set and the length of the template. We recommend setting the annealing temperature 3-5 °C below the T_m of the gene specific primers and the extension time at 1-1.5 minute per kb of target length. If the T_m of the gene specific primers is between 65-72 °C, the control conditions can be used. However, if the T_m of the gene specific primers is above 75 °C, use a two-step PCR profile (94 °C for denaturation and 68 °C or 72 °C for annealing/extension).

Step 2. Exonuclease III digestion

After PCR, add 1 μ l of Exonuclease III to each 50 μ l PCR (final concentration of 2 U/ μ l). Mix thoroughly and incubate at 37 °C for 10 minutes. After the digestion remove a 3-5 μ l aliquot of each reaction for agarose gel analysis.

Step 3. Purification of PCR products

If PCR results reveal a single amplification band with a reasonable yield, proceed with a post-reaction clean-up using GenElute PCR Clean-up Kit) or equivalent system.

If PCR results reveal more than one band, gel extraction and purification should be performed in order to obtain a specific PCR product. The GenElute Gel Extraction kit or equivalent system may be used.

If no PCR product is observed, check the PCR conditions. Refer to the troubleshooting guide in this technical bulletin for assistance with optimizing PCR conditions.

Step 4. Ligation

Ligate Anchors to the amplified gene. Typically, 1-2 μ l of amplified gene product is all that is necessary in the ligation reaction. However, longer genes (>1 kb) may require increasing the amount added to the ligation reaction to ensure adequate ligation efficiency.

Ligation reaction:

14-x μ l	Water
x μ l	First round PCR product (20-100 ng)
2 μ l	5' Anchor (1 ng/ μ l)
2 μ l	3' Anchor (1 ng/ μ l)
2 μ l	10x T4 DNA Ligase Buffer
1 μ l	T4 DNA Ligase
21 μ l	Total volume

Mix thoroughly and incubate at room temperature for 30 minutes.

Step 5. Second Round PCR

Perform second round amplification of assembled template using Anchor primers. Mix the following components in a 0.2 ml or a 0.5 ml PCR tube on ice.

Control and sample reaction:

2 μ l	ligation reaction
1 μ l	Anchor primer-Forward (20 μ M)
1 μ l	Anchor primer-Reverse (20 μ M)
2.5 μ l	JumpStart REDAccuTaq LA DNA Polymerase
5 μ l	AccuTaq LA DNA Polymerase Buffer
2.5 μ l	dNTP mix, 10 mM (yellow cap)
36 μ l	Water
50 μ l	Total volume

Control and sample PCR:

Initial denaturation:	95 °C, 5 minutes
35 cycles:	
denaturation	94 °C, 30 seconds
annealing	60 °C, 30 seconds
extension	68 °C, 2 minutes 30 seconds
Final extension:	68 °C, 10 minutes
Soak:	4 °C

Note: If the assembled template is >2 kb, it may be necessary to modify the PCR profile with a longer extension time.

Remove a 3-5 μ l aliquot of each reaction to check on an agarose gel. Refer to the troubleshooting guide in this technical bulletin if a single band is not visible.

Step 6. Purification (optional) and Use of Expression-Ready Template

Samples may be purified before use in *in vitro* transcription/translation reactions using the GenElute PCR Clean-up Kit (Product Code NA 1020) or equivalent system. This step is optional for some transcription/translation kits. TNT[®] T7 Quick for PCR DNA (Promega L5540) does not require purification. However, purification is recommended for use in the TNTT7 Coupled Wheat Germ Extract System (Product Code TN0100).

Templates are ready for use in any *in vitro* T7 transcription system or any coupled transcription/translation system for eukaryotic expression. Please refer to the manufacturer's specific instructions regarding use of PCR products in the particular transcription/translation system.

References

1. Putney, S. D., et al., A DNA fragment with an alpha-phosphorothioate nucleotide at one end is asymmetrically blocked from digestion by exonuclease III and can be replicated *in vivo*. Proc. Natl. Acad. Sci. USA, **78**, 7350-7354 (1981).
2. Olsen, D. B. and Eckstein, F., Incomplete primer extension during *in vitro* DNA amplification catalyzed by *Taq* polymerase; exploitation for DNA sequencing. Nucleic Acids Res., **17**, 9613-9620 (1989).
3. Don, R. H. et al., "Touchdown" PCR to circumvent spurious priming during gene amplification. Nucleic Acids Res., **19**, 4008 (1991).

CAP/JWM 04/04

Troubleshooting guide

Problem	Possible cause	Solution
A first round PCR product is not observed	A PCR component may be missing or degraded	A positive control should always be run to ensure components are functioning.
	Too few PCR cycles performed	Increase the number of cycles in 3-5 cycle increments
	Annealing temperature may be too high	Decrease the annealing temperature in 2-4 °C increments
	Primer design may not be optimal	Confirm the accuracy of the sequence information. If the primer is less than 15 nucleotides long, try to lengthen the primer to 20 nucleotides (in addition to the 5 base pairs added to the 5' end)
	Not enough DNA template	If increasing the number of cycles has no effect, repeat the reaction with a higher concentration of template
	Poor quality DNA template	Evaluate the template integrity by agarose gel electrophoresis. It may be necessary to repurify the template using methods that minimize shearing and nicking
	Denaturation temperature may be too high or too low	Optimize the denaturation temperature by raising or lowering it in 1 °C increments
	Extension time may be too short	Increase the extension time in 1 minute increments, especially for long templates
	Mg ⁺⁺ levels may be too low	Optimize the magnesium concentration by using a PCR optimization kit (Product Code M 1932)
Multiple PCR products are observed	Too many PCR cycles performed	By reducing the number of cycles, nonspecific bands may be eliminated
	Primer design may not be optimal	Confirm the accuracy of the sequence information. If the primer is less than 15 nucleotides long, try to lengthen the primer to 20 nucleotides (in addition to the 5 base pairs added to the 5' end)
	Annealing temperature may be too low	Increase the annealing/extension temperature in 2-3 °C increments
A Second Round PCR product is not observed	A ligation reaction component may be missing or degraded	A positive control should be run to ensure components are functional
	The PCR product may contain impurities	Repurify the PCR product
	Exonuclease III incubation time may be too short	Increase the exonuclease III digestion to 30 minutes
	Ligation time may be too short	Increase the ligation time up to 4 hours

Patent pending.

LICENSE AGREEMENTS

Purchase of this product is accompanied by a limited license for use in the Polymerase Chain Reaction (PCR) process for research purposes only and in conjunction with a thermal cycler whose use in the automated performance of the PCR process is covered by an up-front license fee, either by payment to Applied Biosystems or as purchased, i.e., and authorized thermal cycler.

Antibody licensed for *in vitro* research use under U.S. Patent No. 5,338,671 and 5,587,287, and corresponding patents in other countries.

Licensed under U.S. Patent Number 5,436,149 owned by Takara Shuzo Co., Ltd.

The enclosed DNA expression vector and/or antibody are specifically adapted for a method of producing selected protein molecules covered by one or more of the following patents owned by Sigma-Aldrich Co.: U.S. Patent Nos. 5,011,912, 4,703,004, 4,782,137 and 4,851,341; EP Patent No. 150,126 (Austria, Belgium, Switzerland, France, United Kingdom, Italy, Netherlands and Sweden); EP Patent No. 335,899 (Belgium, Switzerland, Germany, France, United Kingdom, Italy, Luxembourg and Sweden); German Patent No. P3584260.1; Canadian Patent No. 1,307,752; and Japanese Patent Nos. 1,983,150 and 2,665,359. Your payment includes a limited license under these patents to make only the following uses of these products:

A. Vector License: You may use the enclosed vector to transform cells to produce proteins containing the amino acid sequence DYKDDDDK for research purposes provided, however, such research purposes do not include binding an unlicensed antibody to any portion of this amino acid sequence nor using such proteins for the preparation of antibodies having an affinity for any portion of this amino acid sequence.

B. Antibody License: You may only use the enclosed antibody for research purposes to perform a method of producing a protein in which the protein is expressed in a host cell and purified by use of the antibody in accordance with a claim in one of the above patents in force in a country where the use actually occurs so long as: (1) you perform such method with a DNA expression vector licensed from Sigma-Aldrich Co.; and (2) you do not bind (or allow others to bind) an unlicensed antibody to any DYKDDDDK epitope of any fusion protein that is produced by use of the method.

This license does not include any rights under any other patents. You are not licensed to use the vector and/or antibody in any manner or for any purpose not recited above. As used above, the term "unlicensed antibody" means any antibody which Sigma-Aldrich Co. has not expressly licensed pursuant to Paragraph B, above. Sigma-Aldrich Co. hereby expressly retains all rights in the above listed patents not expressly licensed hereunder.

If the terms and conditions of this License Agreement are acceptable to you, then you may open the vessel(s) containing the vector and/or antibody and, through such act of opening a vessel, will have shown your acceptance to these terms and conditions.

If the terms and conditions of this License Agreement are not acceptable to you, then please return the vessel(s) unopened to Sigma-Aldrich Co. for a complete refund of your payment.

For additional licensing information or to receive a copy of any of the above patents, please contact the Sigma-Aldrich Co. licensing department at telephone number 314-771-5765.

FLAG and ANTI-FLAG are registered trademarks of Sigma-Aldrich Co. JumpStart™, Accutaq™, REDAccuTaq™, ExoClone™, Director™, Anchor™, and GenElute™ are trademarks of Sigma-Aldrich Co. TNT® is a registered trademark of Promega Corp.

Sigma brand products are sold through Sigma-Aldrich, Inc.

Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see reverse side of the invoice or packing slip.