

# Expand 20 kb<sup>PLUS</sup> PCR System

Deoxynucleoside-triphosphate: DNA deoxynucleotidyltransferase, *E.C. 2.7.7.7*

Cat. No. 11 811 002 001

200 U for approximately 40 reactions

 Version 15

Content version: October 2018

Store the kit at -15 to -25°C

## 1. What this Product Does

### Number of PCR Reactions

The kit is designed for approximately 40 reactions (with a final reaction volume of 50 µl each).

### Kit Contents

Vial	Label	Contents
1	Expand 20 kb <sup>PLUS</sup> Enzyme mix	40 µl Enzyme storage buffer: 20 mM Tris-HCl, pH 7.5 (+25°C), 100 mM KCl, 10 mM dithiothreitol (DTT), 0.1 mM EDTA, 0.5% Nonidet P40 (v/v), 0.5% Tween 20 (v/v), 50% glycerol (v/v)
2	Expand 20 kb <sup>PLUS</sup> reaction buffer	1 ml 10 x conc. with 275 mM MgCl <sub>2</sub>
3	MgCl <sub>2</sub> 25 mM Stock Solution	1 ml
4	Human Genomic DNA	12.5 µl (0.2 mg/ml) in 10 mM Tris-HCl, pH 7.5 and 0.1 mM EDTA Store at +2 to +8 °C
5	Human β-globin control primer forward (HβG forw.)	10 µl (158 ng/µl) in double-distilled water 5'-CAC AAG GGC TAC TGG TTG CCG ATT-3'
6	Human β-globin control primer reverse (HβG rev.)	10 µl (198 ng/µl) in double-distilled water 5'-AGC TTC CCA ACG TGA TCG CCT TTC TCC CAT-3'

### Storage and Stability

Store the kit components excluding the human genomic DNA at -15 to -25°C. When properly stored, the kit is stable until the expiration date printed on the label.

⚠ Always thaw and equilibrate all buffers at +37°C to +56°C before use. Vortex thoroughly. If crystals have formed, incubate at +37°C to +56°C until they are dissolved.

⚠ The supplied human control DNA must be stored at +2 to +8 °C since multiple freezing and thawing will degrade the DNA.

### Additional Equipment and Reagents Required

- Template DNA, gene-specific PCR primer pair, dNTP Mix\*
- Water, PCR grade\*
- Thermal block cycler (e.g., Applied Biosystems GeneAmp PCR System 9600)
- 0.2 ml thin-walled PCR tubes
- Sterile reaction tubes for preparing master mixes and dilutions

### Application

Polymerase chain reaction (PCR).

This PCR system is an improvement of the Barnes Technology (1,2,3) and shows good performance for the amplification of fragments longer than 20 kb.

The Expand 20 kb<sup>PLUS</sup> PCR System is composed of a special enzyme mix containing thermostable Taq DNA polymerase and Tgo DNA polymerase, a thermostable DNA polymerase with proofreading activity. This powerful polymerase mixture and associated buffer system is designed to give a high yield of PCR product when fragments longer than 20 kb need to be amplified.

The Expand 20 kb<sup>PLUS</sup> PCR System contains human control DNA and human control β-globin primers which allow the amplification of a 23 kb fragment.

These reagents may serve as a control reaction but can also be used to test the quality of human template DNA's and/or the respective primer pairs.

### Enzyme Properties

Volume activity	5 U/µl
Optimal enzyme concentration	varies from 2.5 to 7.5 U per 50 µl reaction
Standard enzyme concentration	5 U (1 µl) per 50 µl reaction
Optimal elongation temperature	+68°C
Standard Mg <sup>2+</sup> concentration	2.75 mM (as MgCl <sub>2</sub> ) when using 500 µM dNTP's each.
PCR product size	~ 28 kb
PCR Cloning	T/A cloning
Incorporation of dUTP	no
Repair of mismatched primers at 3' end	yes, due to the 3'-5' exonuclease activity of the proofreading polymerase

\* available from Roche Diagnostics

## 2. How To Use this Product

### 2.1 Before You Begin

#### General Considerations

The optimal conditions (incubation times and temperatures, concentrations of enzyme, template DNA, Mg<sup>2+</sup> concentration) depend on the system used and must be determined for each system. In particular, to ensure optimal reaction efficiency, you should titrate the Mg<sup>2+</sup> concentration and the amount of enzyme used per assay. In addition, increasing the cycle number may improve the yield of amplified DNA.

As a starting point for developing your assays, use the following guidelines:

#### Amount of Enzyme

- Optimal enzyme concentration range from 2.5 to 7.5 U per assay.
- The recommended starting concentration is 5 U (1 μl).

#### dNTP/Mg<sup>2+</sup> Concentration

- The combination of 2.75 mM MgCl<sub>2</sub> (concentration of the supplied buffer) with 500 μM dNTPs (each) is recommended. Nevertheless in some cases titration of Mg ions (adding additional Mg ions) may be necessary to obtain optimal results.

ⓐ dNTP concentration: always use balanced solutions of all four dNTP's.

#### Dilution Buffer

- The optimal buffer for dilution of the template DNA is either double-distilled water or 5 to 10 mM Tris (pH 7 to 8).

ⓑ Avoid dissolving the template in TE buffer because EDTA chelates Mg<sup>2+</sup>.

#### Primers

- The potential for secondary structure and dimer formation should be minimized. Typical primers for long PCR amplifications have a length of 22 to 34 nucleotides with balanced melting temperatures > +60°C.
- Such primers permit the use of higher temperatures to enhance reactions specificity. This can be critical as the amplification of long targets will be compromised by preferential amplification of shorter non-specific fragments. The design of primers suitable for the amplification of > 20 kb fragments is very critical. The following forward primer TGC TGC TCT GTG CAT CCG AGT G can be used, in combination with the enclosed HbG reverse primer to amplify a 29.8 kb fragment from the human globin gene. The annealing temperature is +60°C.

#### Hot Start

- Do not use AmpliWax because of resulting difficulties with volume reduction.

#### Cloning

- The obtained PCR fragments have mainly a 3'-single A overhang.

#### Sample Material

Template DNA, *e.g.*, human genomic DNA\*

- The quality (length and purity) of the template influence dramatically the performance of PCR. Therefore, it is recommended to check the length of the DNA by agarose gel electrophoresis. DNA fragments should be longer than 50 kb. A recommended procedure to get high molecular weight DNA is described in ref. 6 and 7.
- Keep denaturation steps as short as possible and denaturation temperature as low as possible (*i.e.*, +92°C).

ⓐ If possible, linearize circular templates.

### 2.2 Preparation of the Reaction Mixes

For a larger number of reactions, we recommend that you prepare two reaction mixes. This circumvents the need of "Hot Start" and avoids that the enzyme interacts with primers and template in the absence of dNTPs which could lead to partial degradation of primer and template through the 3'-5' exonuclease activity of Tgo DNA Polymerase.

- ⓐ It is also recommended to prepare a Master Mix for setting up multiple reactions. The Master Mix typically contains all of the components needed for all PCR tests to be performed at a volume 10% greater than that required for the total number of PCR assays.

### 2.3 Procedure

Please refer to the following table.

The use of 50 μl reaction volumes is recommended. Smaller volumes, *i.e.*, 30 μl are possible. Avoid evaporation by using mineral oil or a thermal block cycler with heated lid.

- 1 Thaw the components and place them on ice. Mix briefly and centrifuge all reagents before starting.

- 2 • Prepare two mixes or reagents in sterile microfuge tubes (on ice):

- **Mix 1** (for one reaction):

Reagent	Vol.	Control	Final conc.
Water, PCR Grade or double-distilled water	add up to 25 μl	19.3 μl	
dNTP mix, 10 mM	2.5 μl	2.5 μl	500 μM (of each dNTP)
Upstream primer	variable	-	400 nM
Downstream primer	variable	-	400 nM
β-globin control forward primer	-	1 μl	
β-globin control reverse primer	-	1 μl	
Template DNA	variable	-	250 to 500 ng genomic DNA
Human genomic DNA	-	1.2 μl	250 ng
<b>Final volume</b>	<b>25 μl</b>	<b>25 μl</b>	

- **Mix 2** (for one reaction):

Reagent	Volume	Control	Final conc.
Water, PCR Grade or double-distilled water	19 μl	19 μl	
Expand 20 kb <sup>plus</sup> reaction buffer, 10×	5 μl	5 μl	1×
Expand 20 kb <sup>plus</sup> enzyme mix	1 μl	1 μl	5 U
<b>Final volume</b>	<b>25 μl</b>	<b>25 μl</b>	

- 3 • Combine Mix 1 and Mix 2 in a thin-walled PCR tube (on ice).  
• Gently vortex the mixture to produce a homogeneous reaction, then centrifuge briefly to collect sample at the bottom of the tube.  
• Continue to thermal cycling immediately.

### 2.4 Thermal Cycling

1. Place samples in the thermal block cycler, and start cycling using one of the thermal profiles below.

	Temperature	Time	Cycle No.
Initial Denaturation	92°C	2 min	1
Denaturation	92°C	10 sec	10
Annealing	variable (62°C <sup>a</sup> for control reaction)	30 sec	
Elongation	68°C (for control reaction)	18 min <sup>b</sup>	
Denaturation	92°C	10 sec	20 <sup>c</sup>
Annealing	variable (62°C <sup>a</sup> for control reaction)	30 sec	
Elongation	68°C (for control reaction)	18 min <sup>b</sup> + 10 sec cycle elongation for each successive cycle <sup>c</sup>	
Final Elongation	68°C	7 min	1
Cooling	4°C	unlimited time	

a) Optimal annealing temperature depends on the melting temperature of the primers and the system used. Appropriate primers should have annealing temperatures > +60°C.

b) Elongation time depends on fragment length. we recommend the following times:

PCR fragment length (kb):	15	20	25	30	35	40	45
Elongation Time (min):	11	14	17	20	23	27	30

Do not forget to extend the elongation time for each new cycle.

c) The number of cycles depends on the amount of template (copies of target) DNA used. For human genomic DNA, we get good results with 250 ng of template using 30 cycles (in total). However, an increase of the cycle number up to 35 or 40 may increase the yield of the amplified DNA.

2. After cycling, the samples may be frozen for later use. Possible further procedures:

- Check the PCR product on an agarose gel for size and specificity using an appropriate size marker\*.
- Purify the PCR product with the High Pure PCR Product Purification Kit\*.

Ⓢ The obtained PCR fragments have mainly a 3'-single A overhang.

Ⓢ The thermal profiles were developed for the Applied Biosystems GeneAmp PCR System 9600. Other thermal block cyclers may require different profiles.

### 3. Troubleshooting

	Possible Cause	Recommendation
<b>Little or no PCR product</b>	DNA template problems	Check quality and concentration of template: <ul style="list-style-type: none"> <li>• Use highest purified template (phenolisation).</li> <li>• Analyze an aliquot on an agarose gel to check for possible degradation.</li> <li>• Make a control reaction on template with an established primer pair or PCR system.</li> <li>• Check or repeat purification of template.</li> <li>• Circular templates should be linearized if possible.</li> </ul>
	MgCl <sub>2</sub> concentration too low	Increase the MgCl <sub>2</sub> concentration in 0.25 mM steps (solution supplied).
	Cycle conditions not optimal	<ul style="list-style-type: none"> <li>• Check annealing temperature and denaturation temperatures. If necessary decrease annealing temperature.</li> <li>• Increase cycle number.</li> <li>• Make sure that the final elongation step was carried out.</li> </ul>
	Primer design not optimal	<ul style="list-style-type: none"> <li>• Design alternative primers.</li> <li>• Both primers should have nearly the same melting temperatures.</li> </ul>
	Primer concentration not optimal	<ul style="list-style-type: none"> <li>• Both primers must have the same concentration.</li> <li>• Titrate primer concentration.</li> </ul>
	Primer quality or storage problems	<ul style="list-style-type: none"> <li>• If you are using an established primer pair, check their performance under established PCR conditions (with a control template).</li> <li>• Make sure primers are not degraded.</li> <li>• Always store primers at -15 to -25°C and as a stock solutions.</li> </ul>

Possible Cause	Recommendation
Annealing temperature too low	Increase annealing temperature, never exceed +68°C.
Multiple bands or background smear	<ul style="list-style-type: none"> <li>• Review primer design.</li> <li>• Titrate primer concentration (0.1 to 0.6 μM).</li> <li>• Both primers must be present in the reaction at the same concentration.</li> <li>• Perform nested PCR with nested primers.</li> </ul>
Enzyme concentration too high	Reduce amount of enzyme
DNA template problems	Use serial dilutions of template.
Cycle conditions not optimal	Reduce the number of cycles (not recommended for human or similar complex DNA).

### 4. Additional Information on this Product

#### References

- 1 Barnes, W.M. (1994) *Proc. Natl. Acad. Sci USA* **91**, 2216-2220
- 2 Cheng, S. *et al.* (1994) *Proc. Natl. Acad. Sci USA* **91**, 5694-5699
- 3 Frey, B. & Suppmann B. (1995) *Biochemica* **2**; 8-9
- 4 Lindahl, T. (1993) *Nature* **362**, 709-715
- 5 Lindahl, T. *et al.* (1972) *Biochemistry* **11**, 3611-3618
- 6 Cheng, S. *et al.* (1995) *PCR Methods and Applications* **4**, 294-298
- 7 Huder, J. B. *et al.* (2002); Identification and Characterization of Two Closely Related Unclassifiable Endogenous Retroviruses in Python (Python molurus and Python curtus) *J. Virol.* **76**, 7607-7615.
- 8 Simonic, T. *et al.* (2000) cDNA cloning of turtle prion protein. *FEBS Letters* **469**, 33-38.

#### Quality Control

Each lot of Expand 20 kb<sup>PLUS</sup> PCR System is function tested in PCR. Routinely, the Expand 20 kb<sup>PLUS</sup> PCR System is used in combination with human genomic DNA and specific human β-globin primers to amplify a 23 kb PCR fragment.

### 5. Supplementary Information

#### 5.1 Conventions

##### Text Conventions

To make information consistent and memorable, the following text conventions are used in this document:

Text Convention	Use
Numbered Instructions	Steps in a process that usually occur in the order listed labeled ①, ②, etc.
Numbered Instructions	Steps in a procedure that must be performed in the order listed labeled 1, 2, etc.
Asterisk *	Denotes a product available from Roche Diagnostics

#### Symbols

In this document, the following symbols are used to highlight important information:

Symbol	Description
Ⓢ	Information Note: Additional information about the current topic or procedure.
⚠	Important Note: Information critical to the success of the procedure or use of the product.

#### 5.2 Changes to Previous Version

- Editorial changes.

### 5.3 Ordering Information

	Product	Pack Size	Cat. No.	
Standard PCR	Taq DNA Polymerase	100 U	11 146 165 001	
		500 U	11 146 173 001	
		4 × 250 U	11 418 432 001	
		10 × 250 U	11 596 594 001	
		20 × 250 U	11 435 094 001	
	PCR Core Kit <sup>PLUS</sup>	1 kit	11 585 541 001	
	PCR Core Kit	1 kit	11 578 553 001	
	PCR Master	1 kit	11 636 103 001	
	Expand High Fidelity <sup>PLUS</sup> PCR System	125 U	03 300 242 001	
		2 × 250 U	03 300 226 001	
		10 × 250 U	03 300 234 001	
	Expand High Fidelity PCR System	100 U	11 732 641 001	
		2 × 250 U	11 732 650 001	
	10 × 250 U	11 759 078 001		
High Fidelity PCR Master	1 kit	12 140 314 001		
Maximum specificity	FastStart Taq DNA Polymerase, 5 U/μl	100 U	12 032 902 001	
		2 × 250 U	12 032 929 001	
		4 × 250 U	12 032 937 001	
		10 × 250 U	12 032 945 001	
		20 × 250 U	12 032 953 001	
	FastStart High Fidelity PCR System	125 U	03 553 426 001	
		2 × 250 U	03 553 400 001	
		10 × 250 U	03 553 361 001	
	High fidelity PCR	Pwo SuperYield DNA Polymerase	100 U	04 340 868 001
			2 × 250 U	04 340 850 001
Pwo SuperYield DNA Polymerase, dNTPack		100 U	04 743 750 001	
Pwo Master		10 × 250 μl	03 789 403 001	
Pwo DNA Polymerase		100 U	11 644 947 001	
		2 × 250 U	11 644 955 001	
Expand High Fidelity PCR System		100 U	11 732 641 001	
		2 × 250 U	11 732 650 001	
		10 × 250 U	11 759 078 001	
High Fidelity PCR Master		1 kit	12 140 314 001	
FastStart High Fidelity PCR System		125 U	03 553 426 001	
		2 × 250 U	03 553 400 001	
		10 × 250 U	03 553 361 001	
Expand High Fidelity <sup>PLUS</sup> PCR System	125 U	03 300 242 001		
	2 × 250 U	03 300 226 001		
	10 × 250 U	03 300 234 001		
Long template PCR	Expand Long Template PCR System	150 U	11 681 834 001	
		2 × 360 U	11 681 842 001	
		10 × 360 U	11 759 060 001	
Expand 20 kb <sup>PLUS</sup> PCR System	200 U	11 811 002 001		
Difficult templates & challenging assays	FastStart Taq DNA Polymerase, 5 U/μl	100 U	12 032 902 001	
		2 × 250 U	12 032 929 001	
		4 × 250 U	12 032 937 001	
		10 × 250 U	12 032 945 001	
		20 × 250 U	12 032 953 001	
	FastStart High Fidelity PCR System	125 U	03 553 426 001	
		2 × 250 U	03 553 400 001	
		10 × 250 U	03 553 361 001	
	GC-RICH PCR System	100 U	12 140 306 001	
	Ready-to-use mixes of all 4 nucleotides	PCR Nucleotide Mix	200 μl	11 581 295 001
		2,000 μl	11 814 362 001	
Ready-to-use mixes of all 4 nucleotides, buffer and polymerases	Pwo Master	10 × 250 μl	03 789 403 001	
	High Fidelity PCR Master	10 × 500 μl	12 140 314 001	
	PCR Master	10 × 500 μl	11 636 103 001	
Kits containing nucleotides, buffers and polymerases	PCR Core Kit	1 kit	11 578 553 001	
	PCR Core Kit <sup>PLUS</sup>	1 kit	11 585 541 001	

	Product	Pack Size	Cat. No.
DNA purification	High Pure PCR Template Preparation Kit	100 purifications	11 796 828 001
	High Pure PCR Product Purification Kit	50 purifications 250 purifications	11 732 668 001 11 732 676 001
Additional reagents	DNA MWM XV (Expand Marker)	50 μg (1 A <sub>260</sub> unit)	11 721 615 001
	Water, PCR Grade	25 ml (25 × 1 ml)	03 315 932 001
		25 ml (1 × 25 ml)	03 315 959 001
		100 ml (4 × 25 ml)	03 315 843 001
		Biotin-16-dUTP	50 μl (50 nmol)
	Fluorescein-12-dUTP	25 μl (25 nmol)	11 373 242 910
	Human Genomic DNA	100 μg	11 691 112 001
Bovine Serum Albumin	1 ml (20 mg)	10 711 454 001	

#### Disclaimer of License

For patent license limitations for individual products please refer to: [List of biochemical reagent products](#)

#### Trademarks

EXPAND, FASTSTART and HIGH PURE are trademarks of Roche. All third party product names and trademarks are the property of their respective owners.

#### Regulatory Disclaimer

For life science research only. Not for use in diagnostic procedures.

#### Contact and Support

To ask questions, solve problems, suggest enhancements and report new applications, please visit our [Online Technical Support Site](#).

To call, write, fax, or email us, visit [sigma-aldrich.com](http://sigma-aldrich.com), and select your home country. Country-specific contact information will be displayed.



Roche Diagnostics GmbH  
Sandhofer Strasse 116  
68305 Mannheim  
Germany