



Easy-A High-Fidelity PCR Master Mix

Instruction Manual

Catalog #600640 (100 reactions), #600642 (400 reactions)

Revision C.0

For Research Use Only. Not for use in diagnostic procedures.

600640-12



Agilent Technologies

LIMITED PRODUCT WARRANTY

This warranty limits our liability to replacement of this product. No other warranties of any kind, express or implied, including without limitation, implied warranties of merchantability or fitness for a particular purpose, are provided by Agilent. Agilent shall have no liability for any direct, indirect, consequential, or incidental damages arising out of the use, the results of use, or the inability to use this product.

ORDERING INFORMATION AND TECHNICAL SERVICES

Email

techservices@agilent.com

World Wide Web

www.genomics.agilent.com

Telephone

Location	Telephone
United States and Canada	800 227 9770
Austria	01 25125 6800
Benelux	02 404 92 22
Denmark	45 70 13 00 30
Finland	010 802 220
France	0810 446 446
Germany	0800 603 1000
Italy	800 012575
Netherlands	020 547 2600
Spain	901 11 68 90
Sweden	08 506 4 8960
Switzerland	0848 8035 60
UK/Ireland	0845 712 5292
All Other Countries	Please visit www.agilent.com/genomics/contactus

Easy-A High-Fidelity PCR Master Mix

MATERIALS PROVIDED

Materials provided	Quantity	
	Catalog #600640 ^a	Catalog #600642 ^b
Easy-A 2 × master mix (0.1 U/μl) ^c	2.5 ml (250 U)	4 × 2.5 ml (1000 U)

^a Catalog #600640 provides enough PCR reagents for 100, 50μl PCR reactions.

^b Catalog #600642 provides enough PCR reagents for 400, 50μl PCR reactions.

^c The total Mg²⁺ concentration present in the final 1 × dilution of the 2 × Easy-A master mix is 2 mM.
The total dNTP concentration present in the final 1 × dilution is 800 μM (200 μM of each dNTP).

Storage: The Easy-A high-fidelity PCR master mix should be stored at –20°C upon receipt. Store the PCR master mix at 4°C after thawing. Once thawed, full activity is guaranteed for 3 months.

NOTICES TO PURCHASER

Use of this product is licensed under U.S. Patent Nos. 5,587,287 and corresponding foreign patents Europe, Hong Kong, and Japan

Limited Label License: The purchase of this product conveys to the purchaser only the limited, non-transferable right to use the purchased quantity of the product for the purchaser's own research by the purchaser only under patent claims outside the U.S. corresponding to the expired U.S. Patent No. 5,436,149 (claims 6-16). No rights are granted to the purchaser to sell, modify for resale or otherwise transfer this product, either alone or as a component of another product, to any third party. Takara Bio reserves all other rights, and this product may not be used in any manner other than as provided herein. For information on obtaining a license to use this product for purposes other than research, please contact Takara Bio Inc., Seta 3-4-1, Otsu, Shiga 520-2193, Japan (Fax +81-77-543-9254).

NOTICE TO PURCHASER: LIMITED LICENSE

Use of this product is covered by one or more of the following US patents and corresponding patent claims outside the US: 6,258,569, 6,171,785, 6,127,155, 6,030,787, 5,994,056, 5,876,930, 5,804,375, 5,789,224, 5,773,258 (claims 1 and 6 only), 5,723,591, 5,677,152 (claims 1 to 23 only), 5,618,711, 5,538,848, and claims outside the US corresponding to expired US Patent No. 5,079,352. The purchase of this product includes a limited, non-transferable immunity from suit under the foregoing patent claims for using only this amount of product for the purchaser's own internal research. No right under any other patent claim and no right to perform commercial services of any kind, including without limitation reporting the results of purchaser's activities for a fee or other commercial consideration, is conveyed expressly, by implication, or by estoppel. This product is for research use only. Diagnostic uses under Roche patents require a separate license from Roche. Further information on purchasing licenses may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA

INTRODUCTION

The Easy-A high-fidelity PCR master mix* is a 2× formulation of the Easy-A high-fidelity PCR cloning enzyme*, an optimized PCR reaction buffer, magnesium, and dNTPs. The Easy-A high-fidelity PCR cloning enzyme is a proprietary thermostable DNA polymerase preparation specifically designed for improved cloning with the StrataClone PCR Cloning Kit or with the TOPO TA Cloning® vector and other T-/U-vectors.[†] The Easy-A PCR cloning enzyme possesses both terminal transferase and exonuclease activities, generating PCR products containing 3'-A overhangs with a higher rate of accuracy. This enables 5-minute, high-fidelity cloning into T-vectors and U-/U*-vectors with an efficiency equivalent to that of *Taq* DNA polymerase. Additionally, the enzyme is provided in an antibody-based hotstart format for increased PCR sensitivity and yield from a variety of templates, while allowing room temperature setup. The Easy-A enzyme amplifies targets up to 10 kb from plasmid DNA and up to 5 kb from genomic DNA.

PCR PROTOCOL

1. Add the PCR reaction components in order while mixing gently. Table I provides an example reaction mixture for the amplification of a typical single-copy chromosomal target. The recipe listed is for one reaction and must be adjusted for multiple samples.

* U.S. Patent Nos. 6,734,293, 6,444,428, 6,183,997.

[†] Use of these cloning vector products may require licenses from third parties in certain countries.

TABLE I Reaction Mixture for a Typical PCR Amplification

Component	Amount per reaction
Distilled water (dH ₂ O)	22 µl
Primer #1 (100 ng/µl) ^a	1 µl
Primer #2 (100 ng/µl) ^a	1 µl
DNA template (100 ng/µl) ^b	1 µl
Easy-A 2× master mix	25 µl
Total reaction volume	50 µl

^a Primer concentrations between 0.2 and 0.5 µM are recommended (this corresponds to 100–250 ng for typical 18- to 25-mer oligonucleotide primers in a 50-µl reaction volume).

^b The amount of DNA template required varies depending on the type of DNA being amplified. Generally 50–100 ng of genomic DNA template is recommended. Less DNA template can be used for amplification of lambda (1–30 ng) and vector (0.1–10 ng) PCR targets or for amplification of multicopy chromosomal genes (10–100 ng).

- Aliquot 50 µl of the reaction mixture into the appropriate number of sterile thin-wall PCR tubes or standard 0.5-ml microcentrifuge tubes.
- Perform PCR using optimized cycling conditions. Suggested cycling parameters are indicated in Table II.
- Analyze the PCR amplification products on a 0.7–1.0% (w/v) agarose gel.
- Use 2 µl of PCR product for cloning into the StrataClone PCR Cloning Kit vector arms. Use 0.5–1.5 µl of PCR product for cloning into T-/U-vectors, following the manufacturer's recommendations.

TABLE II PCR Cycling Parameters for a Typical PCR Amplification^{a,b}

Segment	Number of Cycles	Temperature	Duration
1	1	95°C	2 minutes
2	30	95°C	40 seconds
		Primer T _m – 5°C ^c	30 seconds
		72°C	1 minute for targets ≤ 1 kb; 1 minute per kb for targets > 1 kb and ≤ 5 kb
3	1	72°C	7 minutes

^a Thin-wall PCR tubes are highly recommended.

^b Optimized cycling parameters are not necessarily transferable between thermal cyclers designed by different manufacturers; therefore, each manufacturer's recommendations for optimal cycling parameters should be consulted.

^c The annealing temperature may require optimization. Typically annealing temperatures will range between 55° and 72°C.¹

TROUBLESHOOTING

Observation	Solution(s)
No product or low yield	Increase extension time to 2 minutes per kb of PCR target.
	Use cosolvents such as DMSO in a 1–10% (v/v) final concentration for GC-rich templates.
	Lower the annealing temperature in 5°C increments.
	Denaturation times of 30–60 seconds at 94–95°C are usually sufficient while longer denaturation times may damage the DNA template; use the shortest denaturation time compatible with successful PCR on the thermal cycler.
	Remove extraneous salts from the PCR primers and DNA preparations.
	Use the recommended primer concentrations between 0.3 and 0.5 µM (corresponding to 100–200 ng for typical 18- to 25-mer oligonucleotide primers in a 50-µl reaction volume).
	Check the melting temperature, purity, GC content, and length of the primers.
Multiple bands	Increase the annealing temperature in 5°C increments.
Artificial smears	Reduce the extension time.

REFERENCES

- Innis, M. A., Gelfand, D. H., Sninsky, J. J. and White, T. J. (1990). *PCR Protocols: A Guide to Methods and Applications*. Academic Press, New York.

ENDNOTES

TOPO TA Cloning[®] is a registered trademark of Invitrogen Corp.

MSDS INFORMATION

Material Safety Data Sheets (MSDSs) are provided online at <http://www.genomics.agilent.com>. MSDS documents are not included with product shipments.