

# Taq DNA Polymerase, 5 U/ $\mu$ l

From *Thermus aquaticus* BM, recombinant (*E. coli*)  
Deoxynucleoside-triphosphate: DNA deoxynucleotidyltransferase, EC 2.7.7.7

<b>Cat. No. 11 146 165 001</b>	100 U
<b>Cat. No. 11 146 173 001</b>	500 U
<b>Cat. No. 11 418 432 001</b>	4 × 250 U
<b>Cat. No. 11 596 594 001</b>	10 × 250 U
<b>Cat. No. 11 435 094 001</b>	20 × 250 U

 **Version 24**  
Content version: April 2016

Store at –15 to –25°C

## 1. What this Product Does

### Number of Reactions

If 1.25 U are used per 50  $\mu$ l reaction, Taq DNA Polymerase is designed for approximately:

- 80 reactions (Cat. No. 11 146 165 001)
- 400 reactions (Cat. No. 11 146 173 001)
- 800 reactions (Cat. No. 11 418 432 001)
- 2,000 reactions (Cat. No. 11 596 594 001)
- 4,000 reactions (Cat. No. 11 435 094 001)

### Contents

Label	Contents
Taq DNA Polymerase (5 U/ $\mu$ l)	<ul style="list-style-type: none"> <li>• 20 <math>\mu</math>l (100 U pack size)</li> <li>• 2 × 50 <math>\mu</math>l (500 U pack size)</li> <li>• 4 × 50 <math>\mu</math>l (1,000 U pack size)</li> <li>• 10 × 50 <math>\mu</math>l (2,500 U pack size)</li> <li>• 20 × 50 <math>\mu</math>l (5,000 U pack size)</li> </ul> Enzyme storage buffer: 20 mM Tris-HCl, 1 mM DTT, 0.1 mM EDTA, 0.1 M KCl, 0.5% Nonidet P40 (v/v), 0.5% Tween 20 (v/v), 50% glycerol (v/v), pH 8.0 (+4°C)
PCR Reaction Buffer with MgCl <sub>2</sub> , 10 × conc.	<ul style="list-style-type: none"> <li>• 1 ml (100 U pack size)</li> <li>• 3 × 1 ml (500 U pack size)</li> <li>• 6 × 1 ml (1,000 U pack size)</li> <li>• 15 × 1 ml (2,500 U pack size)</li> <li>• 30 × 1 ml (5,000 U pack size)</li> </ul> Buffer composition: 100 mM Tris-HCl, 15 mM MgCl <sub>2</sub> , 500 mM KCl, pH 8.3 (+20°C)

### Storage and Stability

The undiluted solutions are stable when stored at –15 to –25°C until the expiration date printed on the label.

### Additional Equipment and Reagents Required

Additional equipment and reagents required to perform PCR reactions with Taq DNA Polymerase, but not provided, include:

- standard laboratory equipment
  - nuclease-free, aerosol-resistant pipette tips
  - pipettes with disposable, positive-displacement tips
  - sterile reaction tubes for preparing PCR mixes and dilutions
  - standard benchtop microcentrifuge
- for the PCR reaction
  - thermal block cycler (e.g., the Applied Biosystems GeneAmp PCR System 9600)
  - PCR primers
  - template DNA

- PCR reaction vessels (thin-walled PCR tubes or plates are recommended)
- PCR Nucleotide Mix, PCR Grade\*
- Water, PCR Grade\*

### Application

- Polymerase Chain Reaction (PCR): Taq DNA Polymerase activity is stable during prolonged incubation at high temperatures (+95°C) and can therefore be used to amplify DNA fragments by PCR.
- DNA labeling reactions (4, 5)
- Sequencing / cycle sequencing (4, 6)

### Enzyme Properties

Volume Activity	5 U/ $\mu$ l
Optimal Enzyme Concentration	Varies between 0.5 and 2.5 U per 50 $\mu$ l reaction
Standard Enzyme Concentration	1.25 U per 50 $\mu$ l reaction
Optimal pH	Around 9 (adjusted at +20°C)
Optimal Elongation Temp.	Around +72°C
Optimal Mg <sup>2+</sup> Concentration	Varies between 1.5 and 5 mM (as MgCl <sub>2</sub> )
Standard Mg <sup>2+</sup> Concentration	1.5 mM (as MgCl <sub>2</sub> ) when used with 200 $\mu$ M of each dNTP
Size of PCR Products	Enzyme optimally amplifies up to 3 kb products. (PCR is possible up to 10 kb, but yield diminishes as DNA fragment length increases.)
PCR Cloning	T/A-cloning (Enzyme adds a single, overhanging A.)
Incorporation of Modified Nucleotides	Enzyme accepts modified nucleotides like radiolabeled nucleotides, DIG-dUTP, biotin-dUTP.
Thermostability	Enzyme retains over 80% activity after 30 cycles (1 min +95°C, 1 min +37°C, 3 min +72°C).

## 2. How To Use this Product

### 2.1 Before You Begin

#### General considerations

The optimal conditions (incubation times and temperatures, concentration of enzyme, template DNA, Mg<sup>2+</sup>) vary from system to system and must be determined for each individual experimental system (7). At the very least, you should titrate the Mg<sup>2+</sup> concentration and the amount of enzyme used per assay to ensure optimal efficiency of DNA synthesis.

As a starting point, use the following guidelines:

- Optimal enzyme concentration: 0.5 to 2.5 U/50  $\mu$ l. A concentration of 1.25 U/50  $\mu$ l will usually produce satisfactory results.
- Optimal  $Mg^{2+}$  concentration can vary between 1.5 mM and 5 mM. In most cases a  $Mg^{2+}$  concentration of 1.5 mM will produce satisfactory results (2, 3) if you use 200  $\mu$ M of each dNTP.
- dNTP concentration: Always use equal concentrations of all four dNTPs. The final concentration of each dNTP should be between 50 and 500  $\mu$ M; the most commonly used concentration is 200  $\mu$ M. If you increase the dNTP concentration, you must also increase the  $Mg^{2+}$  concentration.

### Sample Material

- Use any template DNA (*e.g.*, genomic or plasmid DNA, cDNA) suitable for PCR in terms of purity, concentration, and absence of inhibitors. For reproducible isolation of nucleic acids use:
  - either the MagNA Pure 96 Instrument, the MagNA Pure LC Instrument, or the MagNA Pure Compact Instrument together with a dedicated nucleic acid isolation kit (for automated isolation)
  - or a High Pure nucleic acid isolation kit (for manual isolation)
- Ⓢ For details see the Roche Applied Science Biochemicals catalog or home page, [www.roche-applied-science.com](http://www.roche-applied-science.com).
- Use 10 to 500 ng complex genomic DNA or 0.1 to 10 ng plasmid DNA/cDNA. Recommended starting concentrations are 250 ng genomic DNA or 1 ng plasmid DNA.
- ⚠ Store the template DNA either in Water, PCR Grade or 5 to 10 mM Tris-HCl (pH 7 to 8). Avoid dissolving the template in TE buffer because EDTA chelates  $Mg^{2+}$ .

## 2.2 Procedure

### Preparation of PCR Master Mixes

Prepare two PCR master mixes: the first one contains enzyme and reaction buffer, the second one contains all other reaction components. This circumvents the need for hot start and avoids that the enzyme interacts with primers or template during the reaction setup. If you are setting up multiple reactions, the volume of each master mix should be 110% of the volume needed for all your samples (*e.g.*, to prepare Master Mix 2 below for 10 reactions, make 275  $\mu$ l of the mix.) The extra volume allows for losses during pipetting.

### Preparation of Master Mix 1

Step	Action																					
1	<ul style="list-style-type: none"> <li>• Thaw the reagents and store on ice.</li> <li>• Briefly vortex and centrifuge all reagents before setting up the reactions.</li> </ul>																					
2	Prepare a 10 $\times$ conc. solution of each respective primer. <ul style="list-style-type: none"> <li>Ⓢ If you are using, <i>e.g.</i>, the final concentration of 0.5 <math>\mu</math>M for each primer, the 10<math>\times</math> conc. solution would contain a 5 <math>\mu</math>M concentration of the respective primer.</li> </ul>																					
3	To a sterile reaction tube on ice, add the components in the order listed below: (For each 50 $\mu$ l reaction)																					
	<table border="1"> <thead> <tr> <th>Component</th> <th>Volume</th> <th>Final conc.</th> </tr> </thead> <tbody> <tr> <td>Water, PCR Grade</td> <td>to make a final vol. of 25 <math>\mu</math>l</td> <td></td> </tr> <tr> <td>PCR Grade Nucleotide Mix (10 mM of each dNTP)</td> <td>1 <math>\mu</math>l</td> <td>200 <math>\mu</math>M (of each dNTP)</td> </tr> <tr> <td>Downstream primer</td> <td>5 <math>\mu</math>l</td> <td>0.1 to 0.6 <math>\mu</math>M</td> </tr> <tr> <td>Upstream Primer</td> <td>5 <math>\mu</math>l</td> <td>0.1 to 0.6 <math>\mu</math>M</td> </tr> <tr> <td>Template DNA</td> <td>variable</td> <td>0.1 to 250 ng</td> </tr> <tr> <td><b>Final volume</b></td> <td><b>25 <math>\mu</math>l</b></td> <td></td> </tr> </tbody> </table>	Component	Volume	Final conc.	Water, PCR Grade	to make a final vol. of 25 $\mu$ l		PCR Grade Nucleotide Mix (10 mM of each dNTP)	1 $\mu$ l	200 $\mu$ M (of each dNTP)	Downstream primer	5 $\mu$ l	0.1 to 0.6 $\mu$ M	Upstream Primer	5 $\mu$ l	0.1 to 0.6 $\mu$ M	Template DNA	variable	0.1 to 250 ng	<b>Final volume</b>	<b>25 <math>\mu</math>l</b>	
Component	Volume	Final conc.																				
Water, PCR Grade	to make a final vol. of 25 $\mu$ l																					
PCR Grade Nucleotide Mix (10 mM of each dNTP)	1 $\mu$ l	200 $\mu$ M (of each dNTP)																				
Downstream primer	5 $\mu$ l	0.1 to 0.6 $\mu$ M																				
Upstream Primer	5 $\mu$ l	0.1 to 0.6 $\mu$ M																				
Template DNA	variable	0.1 to 250 ng																				
<b>Final volume</b>	<b>25 <math>\mu</math>l</b>																					
4	Mix and centrifuge briefly.																					

### Preparation of Master Mix 2

Step	Action
1	<ul style="list-style-type: none"> <li>• Thaw the reagents and store on ice.</li> <li>• Briefly vortex and centrifuge all reagents before setting up the reactions.</li> </ul>

Step	Action															
2	To a sterile reaction tube on ice, add the components in the order listed below: (For each 50 $\mu$ l reaction)															
	<table border="1"> <thead> <tr> <th>Component</th> <th>Volume</th> <th>Final conc.</th> </tr> </thead> <tbody> <tr> <td>Water, PCR Grade</td> <td>19.75 <math>\mu</math>l</td> <td></td> </tr> <tr> <td>PCR Reaction Buffer, 10<math>\times</math></td> <td>5 <math>\mu</math>l</td> <td>1<math>\times</math> (1.5 mM <math>MgCl_2</math>)</td> </tr> <tr> <td>Taq DNA Polymerase (5 U/<math>\mu</math>l)</td> <td>0.25 <math>\mu</math>l</td> <td>1.25 U/reaction</td> </tr> <tr> <td><b>Final volume</b></td> <td><b>25 <math>\mu</math>l</b></td> <td></td> </tr> </tbody> </table>	Component	Volume	Final conc.	Water, PCR Grade	19.75 $\mu$ l		PCR Reaction Buffer, 10 $\times$	5 $\mu$ l	1 $\times$ (1.5 mM $MgCl_2$ )	Taq DNA Polymerase (5 U/ $\mu$ l)	0.25 $\mu$ l	1.25 U/reaction	<b>Final volume</b>	<b>25 <math>\mu</math>l</b>	
Component	Volume	Final conc.														
Water, PCR Grade	19.75 $\mu$ l															
PCR Reaction Buffer, 10 $\times$	5 $\mu$ l	1 $\times$ (1.5 mM $MgCl_2$ )														
Taq DNA Polymerase (5 U/ $\mu$ l)	0.25 $\mu$ l	1.25 U/reaction														
<b>Final volume</b>	<b>25 <math>\mu</math>l</b>															
3	Mix and centrifuge briefly.															

## 2.3 PCR

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

Step	Action
1	<ul style="list-style-type: none"> <li>• For each reaction, combine 25 <math>\mu</math>l Master Mix 1 and 25 <math>\mu</math>l Master Mix 2 in a thin-walled PCR tube on ice.</li> <li>• Gently vortex the mixture to produce a homogeneous reaction, then centrifuge briefly to collect the solution at the bottom of the tube.</li> <li>⚠ Start thermal cycling immediately. Do not store complete reaction mixes on ice.</li> </ul>

- 2 Place your samples in a thermal block cycler and use either of the thermal profiles below to perform PCR.

• Thermal Profile A: fixed extension time			
	Cycles	Time	Temp
Initial Denaturation	1	2 min	94°C
Denaturation	25 to 30	15 to 30 sec	94°C
Annealing		30 to 60 sec	55 to 65°C
Elongation		45 sec to 3 min	72 or 68°C
Final Elongation	1	7 min	72 or 68°C
Cooling		indefinitely	4°C

• Thermal Profile B: gradually increasing extension time (This procedure ensures a higher yield of amplification products.)			
	Cycles	Time	Temp
Initial Denaturation	1	2 min	94°C
Denaturation	10	15 to 30 sec	94°C
Annealing		30 to 60 sec	55 to 65°C
Elongation		45 sec to 3 min	72 or 68°C
Denaturation	15 to 20	15 to 30 sec	94°C
Annealing		30 sec	55 to 65°C
Elongation		45 sec to 3 min	72 or 68°C
		+ 5 sec cycle elongation for each succ. cycle <sup>a</sup>	
Final Elongation	1	7 min	72 or 68°C
Cooling		indefinitely	4°C

<sup>a</sup> For example, cycle no. 11 is 5 sec longer than cycle 10, cycle no. 12 is 10 sec longer than cycle 10, cycle no. 13 is 15 sec longer than cycle 10, etc.

- Ⓢ The denaturation temperature can vary between +92°C and +95°C. The standard denaturation temperature is +94°C.

Optimal annealing temperature depends on the melting temperature of the primers and on the experimental system. For PCR products up to 1 kb, elongation temperature should be around +72°C; for PCR products larger than 1 kb, elongation temperature should be around +68°C.

- 3 After cycling, if the samples are not used immediately, store them frozen for later use.

- Ⓢ For best results, do the following:

- Check the PCR product on an agarose gel for size and specificity. Use an appropriate size marker.
- Purify the PCR product with the High Pure PCR Product Purification Kit\* (*e.g.*, before performing nested PCR).

## 2.4 DIG DNA Labeling

Digoxigenin 11-dUTP\* is incorporated into DNA by Taq DNA Polymerase. Please refer to Roche Applied Science DIG Kits, DIG Product Selection Guide or DIG Manuals for detailed protocols. For direct access please visit <http://www.roche-applied-science.com/DIG>.

## 3. Troubleshooting

	Possible Cause	Recommendation
<b>Little or no PCR product</b>	Difficult template <i>e.g.</i> , GC-rich templates	<ul style="list-style-type: none"> <li>Perform PCR with GC-RICH PCR System*.</li> <li>Add DMSO (final concentration, 8%) and reduce enzyme concentration (<i>e.g.</i>, use as little as 0.5 U per reaction).</li> </ul>
	DNA template problems	<ul style="list-style-type: none"> <li>Check quality and concentration of template:</li> <li>Analyze an aliquot on an agarose gel to check for possible degradation.</li> <li>Test the template with an established primer pair or PCR system.</li> <li>Check or repeat template purification.</li> </ul>
	Enzyme concentration too low	<ul style="list-style-type: none"> <li>Increase enzyme concentration to 2 U Taq DNA Polymerase per 50 µl reaction.</li> <li>If necessary, increase the amount of polymerase in 0.5 U steps.</li> </ul>
	MgCl <sub>2</sub> concentration too low	Increase the MgCl <sub>2</sub> concentration in 0.25 mM steps; the minimal acceptable concentration is 1.5 mM MgCl <sub>2</sub> .
	Cycle conditions not optimal	<ul style="list-style-type: none"> <li>Decrease annealing temperature.</li> <li>Increase cycle number.</li> <li>Make sure that the final elongation step is included in the program.</li> </ul>
Primer design not optimal	Design alternative primers.	
Primer concentration not optimal	<ul style="list-style-type: none"> <li>Both primers must have the same concentration.</li> <li>Titrate primer concentration (0.1 to 0.6 µM).</li> </ul>	
Primer quality or storage problems	<ul style="list-style-type: none"> <li>If you use an established primer pair, check performance in an established PCR system (<i>e.g.</i>, with a control template).</li> <li>Make sure that the primers are not degraded.</li> <li>Always store primers at -15 to -25°C.</li> </ul>	
Formation of primer dimers	<ul style="list-style-type: none"> <li>Use two Master Mixes, as directed in the protocol above.</li> <li>Use FastStart Taq DNA Polymerase* instead of Taq DNA Polymerase.</li> </ul>	

	Possible Cause	Recommendation
<b>Multiple bands or background smear</b>	Annealing temperature too low	Increase annealing temperature (Longer primers have higher annealing temperatures).
	Primer design or concentration not optimal	<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 have the same concentration.</li> <li>Perform nested PCR with nested primers.</li> </ul>
	Difficult template ( <i>e.g.</i> , GC-rich template)	Perform PCR with GC-RICH PCR System*.
<b>PCR products in negative control experiments</b>	DNA template problems	Use serial dilution of template.
	Carryover contamination	<ul style="list-style-type: none"> <li>Replace all reagents, especially water.</li> <li>Use aerosol-resistant pipette tips.</li> <li>Set up PCR reactions in an area separate from that used for PCR product analysis.</li> <li>To eliminate carryover contaminants: Use dUTP* (600 µM) instead of dTTP (200 µM) and thermolabile UNG* (1 U/50 µl reaction); also, increase Mg<sup>2+</sup> concentration (to a maximum of 4 mM) to compensate for higher dNTP conc.</li> </ul>
<b>Problems specific to RT-PCR</b>	No product, additional bands, background smear	<ul style="list-style-type: none"> <li>The volume of cDNA template (RT-reaction) should not exceed 10% of the final volume of the PCR reaction.</li> <li>Follow troubleshooting tips above.</li> <li>Increase MgCl<sub>2</sub> in 0.25 mM steps.</li> </ul>

## 4. Additional Information on this Product

### How this Product Works

Taq DNA Polymerase (1, 2) is a highly processive 5'→3' DNA polymerase that lacks 3'→5' exonuclease activity (3). It is a single polypeptide chain with a molecular weight of approximately 95 kDa.

Taq DNA Polymerase was originally isolated from the thermophilic eubacterium *Thermus aquaticus* BM, a strain lacking Taq I restriction endonuclease. The enzyme was cloned in *E. coli*.

### Unit Definition

One unit Taq DNA Polymerase is defined as the amount of enzyme that incorporates 10 nmol of total deoxyribonucleosidetriphosphates into acid precipitable DNA within 60 min at +65°C under the assay conditions given above.

### Unit Assay

**Incubation buffer:** 67 mM Tris/HCl; pH 8.3 (+25°C), 5 mM MgCl<sub>2</sub>, 10 mM 2-mercaptoethanol, 0.2% polydocanol, 0.2 mg/ml gelatine, 0.2 mM each dATP, dGTP, dTTP and 0.1 mM dCTP.

**Incubation procedure:** M13mp9ss, M13 primer (17mer) and 1 µCi [ $\alpha$ -<sup>32</sup>P] dCTP are incubated with suitable dilutions of Taq DNA Polymerase in 50 µl incubation buffer for 60 min at +65°C. The amount of incorporated dNTPs is determined by trichloroacetic acid precipitation.

#### 4.1 References

- Chien, A., Edgar, D. B. & Trela, J. M. (1976) Deoxyribonucleic acid polymerase from the extreme thermophile *Thermus aquaticus*. *J. Bacteriol.* **127**, 1550-1557.
- Lawyer, F. C. *et al.* (1989) Isolation, characterization and expression in *Escherichia coli* of the DNA polymerase gene from the extreme thermophile *Thermus aquaticus*. *J. Biol. Chem.* **264**, 6427-6437.
- Tindall, K. R. & Kunkel, T. A. (1988) Fidelity of DNA synthesis by the *Thermus aquaticus* DNA polymerase. *Biochemistry* **27**, 6008-6013.
- Innis, M. A., *et al.* (1988) DNA sequencing with *Thermus aquaticus* DNA polymerase and direct sequencing of polymerase chain reaction-amplified DNA. *Proc. Natl. Acad. Sci. USA* **85**, 9436-9440.
- Lo, Y.-M. D., Mehal, W. Z. & Fleming, K. A. (1988) Rapid production of vector-free biotinylated probes using the polymerase chain reaction. *Nucleic Acids Res.* **16**, 8719.
- Taq polymerase: increased enzyme versatility in DNA sequencing (1988) Applied Biosystems.
- Erllich, H. A. (ed.) (1989) PCR Technology: Principles and Application for DNA Amplification, Stockton Press, New York.
- Mesquita, P. (2003) Human MUC2 mucin gene is transcriptionally regulated by cdx homeodomain proteins in gastrointestinal carcinoma cell lines. *J. Biol. Chem.* **278**: 51549-51556.
- Zhu, Y. (2002) Hemin induces neuroglobin expression in neural cells. *Blood* **100**: 2494-2498.

#### 4.2 Quality Control

Each lot of Taq DNA Polymerase is tested for contaminating activities as described in the following:

##### Test Buffer

10 mM Tris-HCl, 1.5 mM MgCl<sub>2</sub>, 50 mM KCl, pH 8.3 (+20°C).

##### Absence of Endonucleases

Lambda DNA (1 µg) is incubated with Taq DNA Polymerase in 50 µl test buffer for 16 hours at +37°C. The amount of enzyme that shows no degradation of the lambda DNA is > 30 U.

##### Absence of Nicking Activity

Supercoiled pBR322 DNA (1 µg) is incubated with Taq DNA Polymerase in 50 µl test buffer for 4 hours at +37°C. The amount of enzyme that shows no relaxation of the supercoiled DNA is > 30 U.

##### Absence of Exonucleases

Different amounts of Taq DNA Polymerase are incubated in 100 µl test buffer containing [<sup>3</sup>H]-labeled DNA, overlaid with paraffin oil, for 4 hours at +65°C. The amount of enzyme that shows no exonuclease activity is > 15 U.

### 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 labeled ①, ②, etc.	Steps in a process that usually occur in the order listed
Numbered Instructions labeled ❶, ❷, etc.	Steps in a procedure that must be performed in the order listed
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.

#### Changes to Previous Version

- Editorial Changes

#### 5.2 Ordering Information

	Product	Pack Size	Cat. No.
<b>Nucleotides</b>	PCR Nucleotide Mix	200 µl	11 581 295 001
		2,000 µl	11 814 362 001
<b>DNA Purification</b>	High Pure PCR Template Purification 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
<b>Additional Reagents</b>	Taq DNA Polymerase, dNTPack	100 U	04 728 866 001
		500 U	04 728 874 001
		4 × 250 U	04 728 882 001
		10 × 250 U	04 728 904 001
		20 × 250 U	04 728 858 001
	Digoxigenin-11-dUTP (alkali-labile)	25 nmol (25 µl) 125 nmol (125 µl)	11 573 152 910 11 573 179 910
	Digoxigenin-11-dUTP (alkali-stable)	25 nmol (25 µl)	11 093 088 910
	Biotin-16-dUTP	50 nmol (50 µl)	11 093 070 910
	Fluorescein-12-dUTP	25 nmol (25 µl)	11 373 242 910
	Water, PCR Grade	25 ml (25 x 1 ml)	03 315 932 001
		25 ml (1 x 25 ml)	03 315 959 001
		100 ml (4 x 25 ml)	03 315 843 001

##### Regulatory Disclaimer

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

##### Trademarks

FASTSTART, MAGNA PURE, MAGNA LYSER, LIGHTCYCLER, LC and HIGH PURE are Trademarks of Roche. All other product names and trademarks are the property of their respective owners.

##### License Disclaimer

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

#### 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