EUROPE

LIEGE SCIENCE PARK • 4102 Seraing • Belgium • Tel.: +32 4 372 74 00 Fax: +32 4 372 75 00 • info@eurogentec.com • www.eurogentec.com



ANASPEC - 34801 Campus Drive • Fremont, CA 94555 • USA Tel.: 510-791-9560 • toll-free: 800-452-5530 • Fax: 510-791-9572 info.anaspec@eurogentec.com • www.eurogentec.com



Red'y'Star Mix PK-0073-02R • PK-0073-02RSA

Eurogentec products are sold for research or laboratory use only and are not to be administrated to humans or used for medical diagnostics.

Description

The ready to use 2x Red'y'Star Mix contains HotGoldStar *Taq* DNA polymerase, dNTPs, $MgCl_2$ and buffer. To prepare amplification, only add your primers, template DNA and water to the Mix. The 2x Red'y'Star Mix also contains a red dye buffer to enable immediate loading on agarose or polyacrylamide gels following the DNA amplification.

Without activation, HotGoldStar *Taq* DNA polymerase completely lacks any activity below 74°C ("HotStart" activity avoids non specific priming at low temperature.)

Package contents

Reagent	Volume	Description
Red'y'Star Mix PK-0073-02R PK-0073-02RSA	5 x 1ml 1 ml	2 X PCR Mix: HotGoldstar, dNTPs, MgCl ₂ , buffer, red loading dye
MgCl ₂ 25 mM	1 ml	Additional 25 mM MgCl ₂ for optimization if needed (see table behind)

Shipping conditions

Shipped on dry ice.

Storage conditions and stability

Red'y'star Mix can be stored at -20 °C (in a constant temperature freezer) for 24 months or at 4 °C for 3 months. Do not repeat more than 10 freeze/thaw cycles.

Quality control

Each lot is tested for activity by PCR. Using λ DNA as template we guarantee an amplification of at least 10⁵ fold.

Unit Definition

One unit of enzyme is defined as the amount required for incorporation of 10 nmoles of dNTPs into acid - insoluble material after 30 minutes amplification at 72 °C under the standard reaction conditions.

Reaction conditions

25 μ l of the 2x Red'y'Star Mix diluted to a final volume of 50 μ l will give a reaction medium that contains 1 units of HotGoldStar DNA polymerase, 200 μ M dNTPs, 1.5 mM MgCl₂, 50 mM KCl, 15 mM Tris-HCl (pH 8.0 at 25 °C), 0.01% (v/v) stabilizer and red dye loading buffer.

Procedure

1. Thaw vial, mix and place on ice.

2. To 25 μl of Red'y'Star Mix, add 0.1 nmol of primers, <1 μg of template DNA and H_2O to bring the total reaction volume to 50 $\mu l.$

Standard cycling conditions

HotGoldStar Activation:	10 min at 95°C

35 cycles:		
Denaturation	ר	20 sec at 94°C
Annealing		20 sec at 62°C
suggested	Elongation	30 sec at 72°C

No amplification will be observed without the 10 minutes activation step at 95 $^{\circ}\mathrm{C}$

Time and temperature for denaturation and annealing steps depend on the type of machine and primers. We advise that you check primer design using primer design software.

EUROPE

LIEGE SCIENCE PARK • 4102 Seraing • Belgium • Tel.: +32 4 372 74 00 Fax: +32 4 372 75 00 • info@eurogentec.com • www.eurogentec.com



ANASPEC - 34801 Campus Drive • Fremont, CA 94555 • USA Tel.: 510-791-9560 • toll-free: 800-452-5530 • Fax: 510-791-9572 info.anaspec@eurogentec.com • www.eurogentec.com



MgCl, optimization

MgCl ₂ final concentration	2× PCR Mix	MgCl ₂ (25 mM)	Primers, template, H ₂ O
1.5 mM	25 µl	0 µl	25 µl
2.0 mM	25 µl	1 µl	24 µl
2.5 mM	25 µl	2 µl	23 µl
3.0 mM	25 µl	3 µl	22 µl
3.5 mM	25 µl	4 µl	21 µl
4.0 mM	25 µl	5 µl	20 µl
4.5 mM	25 µl	6 µl	19 µl
5.0 mM	25 µl	7 µl	18 µl

Related products

Reagent	Pack size	Reference
dNTP Mix 20 mM total	1 x 20µmol 5 x 20µmol	NU-0010-10 NU-0010-50
dNTP Set 100 mM total	4 x 25µmol	NU-0020-50
HotGoldStar Mix	5 x 1 ml	PK-0073-02
SmartLadder DNA ladder	1000 lanes	MW-1700-10
Molecular Biology Grade Agarose	500 g 1000 g	EP-0010-05 EP-0010-10
Mupid [®] -One electrophoresis system	1	MU-0041

For further information please contact our Customer Help Desk:

TDS-PK0073-02R-1111-V4

For Europe:

E-mail: info@eurogentec.com Tel: +32 4 372 76 65

For USA:

E-mail: info.anaspec@eurogentec.com Tel.: 510-791-9560 • toll-free: 800-452-5530

NOTICE TO PURCHASER : LIMITED LICENSE

Roche Enzyme Products Conveying Polymerase Patent Rights • Use of this product is covered by one or more of the following US patents and corresponding patent claims outside the US: 5,789,224, 5,618,771, 6,127,155 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, no right to perform any patented method, 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.

Chemically-Modified Hot-Start Polymerase Reagents and Kits • Use of this product is covered by one or more of the following US patents and corresponding patent claims outside the US: 6,127,155 5,677,152 (claims 1 to 23 only), 5,773,258 (claims 1 and 6 only), 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.

Mupid® is a registered trademark of Advance Co., Ltd.