

Booster Mix

Reference: ME-0076-02

Licensed for PCR

Products and procedures described in this protocol are intended for research purposes only.

Batch details

Units per vial : ME-0076-02 2 x 1.2 ml

Description

- Compatible with all commercially available thermostable DNA polymerases
- Ideal for difficult templates

Booster Mix is a special 2x additive for use in reactions involving any thermostable DNA polymerase, and is designed to dramatically improve reaction specificity. Booster Mix provides an optimized composition of reagents, and is ideally suited to dirty/difficult templates with GC or AT rich DNA, repetitive sequences or sequences with a high level of secondary structure.

Booster Mix acts as a melting agent by allowing the DNA polymerase and oligonucleotides greater access to the template DNA. Booster Mix does not contain magnesium, dNTPs, or buffer components. In some cases it may be necessary to optimize the Magnesium concentration.

Package contents

Reagent	Volume	Description
2x Booster Mix	2x 1.2 ml	ME-0076-02

Shipping conditions

Shipping at ambient temperature has no detrimental effect on the performance of this enzyme (if lower than 35 °C).

Storage conditions

Storage at -20 °C is recommended.

Directions for use

- Compose the reaction mix, containing buffer, dNTPs, Mg²⁺, template DNA, primers, DNA polymerase.
- Add the Booster Mix at the volume of half of the final volume of the reaction (e.g. 25 µl per 50 µl final volume, etc).
- Add ddH₂O up to the final volume and mix with pipetting.

NB: Booster Mix should not be used in combination with any other additives for polymerase reactions.

Related products

Reagent	Package size	Reference
dNTP Mix 20 mM total	1X 20 µmol 5X 20 µmol 10X 20 µmol	NU-0010-10 NU-0010-50 NU-0010-100
dNTP Set 5 mM each NTP	4X 5 µmol 4X 25 µmol	NU-0020-10 NU-0020-50
Accustar™ DNA polymerase	250 U 500 U	ME-0067-02 ME-0067-05
Authentik™ DNA polymerase	250 U 500 U	ME-0077-02 ME-0077-05
HotGoldStar DNA Polymerase	500 Units	ME-0064-05
SilverStar DNA Polymerase	500 Units	ME-0074-05

For further information please contact our Customer Help Desk:

For Europe:

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

For USA:

E-mail: info.usa@eurogentec.com
Tel: +1 858 793 26 61



Eurogentec Headquarters

LIEGE SCIENCE PARK • 4102 SERAING • Belgium • Tel: +32 4 372 76 65
Fax: +32 4 372 75 00 • info@eurogentec.com • www.eurogentec.com

Eurogentec North America

3347 Industrial Ct • Suite A • San Diego, CA 92121 • USA • Tel: +1 858 793 26 61
Fax: +1 858 793 26 66 • info.usa@eurogentec.com • www.eurogentec.com

Ref. TDS-ME-0076-0408-V1

Booster Mix

Reference: ME-0076-02

Licensed for PCR

Products and procedures described in this protocol are intended for research purposes only.

Batch details

Units per vial : ME-0076-02 2 x 1.2 ml

Description

- Compatible with all commercially available thermostable DNA polymerases
- Ideal for difficult templates

Booster Mix is a special 2x additive for use in reactions involving any thermostable DNA polymerase, and is designed to dramatically improve reaction specificity. Booster Mix provides an optimized composition of reagents, and is ideally suited to dirty/difficult templates with GC or AT rich DNA, repetitive sequences or sequences with a high level of secondary structure.

Booster Mix acts as a melting agent by allowing the DNA polymerase and oligonucleotides greater access to the template DNA. Booster Mix does not contain magnesium, dNTPs, or buffer components. In some cases it may be necessary to optimize the Magnesium concentration.

Package contents

Reagent	Volume	Description
2x Booster Mix	2x 1.2 ml	ME-0076-02

Shipping conditions

Shipping at ambient temperature has no detrimental effect on the performance of this enzyme (if lower than 35 °C).

Storage conditions

Storage at -20 °C is recommended.

Directions for use

- Compose the reaction mix, containing buffer, dNTPs, Mg²⁺, template DNA, primers, DNA polymerase.
- Add the Booster Mix at the volume of half of the final volume of the reaction (e.g. 25 µl per 50 µl final volume, etc).
- Add ddH₂O up to the final volume and mix with pipetting.

NB: Booster Mix should not be used in combination with any other additives for polymerase reactions.

Related products

Reagent	Package size	Reference
dNTP Mix 20 mM total	1X 20 µmol 5X 20 µmol 10X 20 µmol	NU-0010-10 NU-0010-50 NU-0010-100
dNTP Set 5 mM each NTP	4X 5 µmol 4X 25 µmol	NU-0020-10 NU-0020-50
Accustar™ DNA polymerase	250 U 500 U	ME-0067-02 ME-0067-05
Authentik™ DNA polymerase	250 U 500 U	ME-0077-02 ME-0077-05
HotGoldStar DNA Polymerase	500 Units	ME-0064-05
SilverStar DNA Polymerase	500 Units	ME-0074-05

For further information please contact our Customer Help Desk:

For Europe:

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

For USA:

E-mail: info.usa@eurogentec.com
Tel: +1 858 793 26 61



Eurogentec Headquarters

LIEGE SCIENCE PARK • 4102 SERAING • Belgium • Tel: +32 4 372 76 65
Fax: +32 4 372 75 00 • info@eurogentec.com • www.eurogentec.com

Eurogentec North America

3347 Industrial Ct • Suite A • San Diego, CA 92121 • USA • Tel: +1 858 793 26 61
Fax: +1 858 793 26 66 • info.usa@eurogentec.com • www.eurogentec.com

Ref. TDS-ME-0076-0408-V1

Belgium	United Kingdom	France	Germany • Austria	The Netherlands	Switzerland	Eurogentec North America Inc.
Eurogentec Headquarters +32 4 372 74 00 +32 4 372 75 00 info@eurogentec.com	Eurogentec Ltd. +44 (0) 1794 511 411 +44 (0) 1794 522 417 info.uk@eurogentec.com	Eurogentec France s.a.s.u. +33 2 41 73 33 73 +33 2 41 73 10 26 info@eurogentec.com	Eurogentec Deutschland GmbH +49 221 258 94 55 +49 221 258 94 54 info@eurogentec.com	Eurogentec Nederland b.v. +31 43 352 06 98 +31 43 354 19 65 info@eurogentec.com	Eurogentec s.a. succursale de Genève +32 4 372 74 00 +49 221 258 94 55 +32 4 372 75 00 +49 221 258 94 54	3347 Industrial Court - Suite A San Diego, CA 92121, USA 877 387 6436 858 793 2666 info.usa@eurogentec.com

Belgium	United Kingdom	France	Germany • Austria	The Netherlands	Switzerland	Eurogentec North America Inc.
Eurogentec Headquarters +32 4 372 74 00 +32 4 372 75 00 info@eurogentec.com	Eurogentec Ltd. +44 (0) 1794 511 411 +44 (0) 1794 522 417 info.uk@eurogentec.com	Eurogentec France s.a.s.u. +33 2 41 73 33 73 +33 2 41 73 10 26 info@eurogentec.com	Eurogentec Deutschland GmbH +49 221 258 94 55 +49 221 258 94 54 info@eurogentec.com	Eurogentec Nederland b.v. +31 43 352 06 98 +31 43 354 19 65 info@eurogentec.com	Eurogentec s.a. succursale de Genève +32 4 372 74 00 +49 221 258 94 55 +32 4 372 75 00 +49 221 258 94 54	3347 Industrial Court - Suite A San Diego, CA 92121, USA 877 387 6436 858 793 2666 info.usa@eurogentec.com