

Reverse Transcriptase Core kit Technical Data Sheet

Reference: RT-RTCK-05

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

Storage conditions

For long term storage the Reverse Transcriptase Core kit should be stored at -65 °C to -75 °C in a constant temperature freezer. When stored under these conditions the reagents are stable for 1 year.

For short term storage the Reverse Transcriptase Core kit can be stored at -15 °C to -25 °C for 6 months.

Kit contents

The Reverse Transcriptase Core kit contains enough reagents for up to 500 - 10 µl reactions.

Reagent	Volume	Description
10x reaction buffer (black cap)	2.5 ml	One bottle of RT reaction buffer KCl and Tris-HCl
EuroScript reverse transcriptase (white cap)	125 µl	One tube of Moloney Murine leukemia virus reverse transcriptase, 6250 U at 50 U/µl
RNase Inhibitor (purple cap)	200 µl	One tube of RNase inhibitor 4000U at 20 U/µl
2.5 mM dNTP Mix (green cap)	1.25 ml	One tube of dATP, dCTP, dGTP and dTTP in autoclaved, deionized water titrated with NaOH to pH 7.0
25 mM MgCl ₂ (orange cap)	1.5 ml	One tube of 25 mM MgCl ₂
Oligo d(T) ₁₅ VN (yellow cap)	250 µl	One tube containing 50 µM oligodeoxynucleotides of sequence d(T) ₁₅ VN in 10 mM Tris-HCL, pH 8.3
Random nonamers (pink cap)	250 µl	One tube containing 50 µM short oligonucleotides of random sequence (d(N) ₉) in 10 mM Tris-HCL pH 8.3
RNase free water (plain cap)	1.75 ml	One tube of DEPC water

Procedure for Two step RT qPCR reaction

1- Thaw all required reagents necessary for the RT step completely and put them on ice, except for the EuroScript, which should be kept in the freezer until required for use. Mix all reagents well by inversion and spin them down prior to pipeting.

2- Prepare the RT Reaction Mix (sufficient for 200 ng total RNA per 10 µl RT step)

Component	Volume (µl)	Final concentration
10x reaction buffer	1	1x
25 mM MgCl ₂	2	5 mM (or as required)
2.5 mM dNTP	2	500 µM each dNTP
Random nonamer*	0.5	2.5 µM
RNase Inhibitor	0.2	0.4 U/µl
EuroScript RT	0.25	1.25 U/µl
RNase free water	3.75	-
Template	1	10 pg - 200 ng Total RNA
Total Mix	10 µl	

*Note: random nonamers, oligo d(T)₁₅ VN or sequence specific primers can be used for primers. For nonamers and oligo d(T)₁₅ VN the final concentration should be 200 nM.

2.1- To correct for dispensing losses prepare an excess of reaction mix (for example 100 reactions reaction mix for 96 reactions). Add all components together, except for the template. Mix thoroughly by inversion. Spin down.

2.2- Add the reaction mix to the reaction vial. reaction set up should be done on ice

2.3- Add the template to individual reactions, gently mix by inversion. Spin down. A negative control containing no RNA template should always be included. Optionally a no RT-control should be set up in tubes / wells, which do not contain the EuroScript RT / RNase Inhibitor.

2.4- Program the Real-Time thermocycler using the following recommended parameters:

Initial step*	10 min 25 °C
Reverse Transcriptase step	30 min 48 °C
Inactivation of the RT enzyme	5 min 95 °C

* Only if random nonamers or oligo d(T)₁₅ VN are used are used.

The RT Core kit can be combined with any Eurogentec qPCR kits

3- Thaw all required reagents necessary for the PCR step completely and put them on ice. Mix all reagents well by inversion and spin them down prior to pipeting.

4- Prepare the PCR Reaction Mix

4.1- In case of a probe assay:

Component	Volume (µl)		Final concentration
Reaction buffer	5*	25**	1x
50 mM MgCl ₂	5	-	5 mM (or as required)
5 mM dNTP	2	-	200 µM each dNTP
Forward primer	5	5	-
Reverse primer	5	5	-
Probe	5	5	-
HotGoldStar	0.25	-	0.025 U / µl
RNase free water	7.75	-	-
Total Mix	40 µl		

* If using a qPCR Core kit

** If using a qPCR MasterMix

Add 40 µl of PCR reaction Mix to 10 µl or a dilution of the 1st strand reaction mix.

4.2- In case of a SYBR® green I assay

Component	Volume (µl)		Final concentration
Reaction buffer	5*	25**	1x
50 mM MgCl ₂	5	-	5 mM (or as required)
5 mM dNTP	2	-	200 µM each dNTP
Forward primer	5	5	-
Reverse primer	5	5	-
diluted SYBR green I	1.5	-	-
HotGoldStar	0.25	-	0.025 U / µl
RNase free water	7.75	-	-
Total Mix	40 µl		

* If using a qPCR Core kit for SYBR® green I

** If using a qPCR MasterMix Plus for SYBR® green I

Add 40 µl of PCR reaction Mix to 10 µl or a dilution of the 1st strand reaction mix.

4.3- To correct for dispensing losses prepare an excess of reaction mix (for example 100 reactions reaction mix for 96 reactions). Add all components together, except for the template. Mix thoroughly by inversion. Spin down.

4.4- Add 10 µl or a dilution of the first strand reaction mix, 5 µl of the template control (plus 5 µl of water or buffer) for your positive control and 10 µl of water or buffer for your negative control in to your PCR tubes / 96-well plate / 384-well plate.

4.5- Add 40 µl of the reaction mix to the reaction vial, close the vial and mix gently on a stirrer or spin down. Ensure that no bubbles are present in the reaction vial. Reaction set up can be done at room temperature.

4.6- Program the Real-Time thermocycler using the following recommended parameters:

UNG step	2 min. 50 °C
HotGoldStar activation/UNG inactivation	10 min. 95 °C
40 Cycles	15 sec. 95 °C 1 min. 60 °C
Hold	50 °C forever

For any further informations concerning the qPCR step, in terms of primer and probe design, primer and probe concentrations, MgCl₂ concentration or concerning optimization of the reaction, please refer to the instruction of the Eurogentec qPCR Core kits or qPCR MasterMixes.

Further information available through Eurogentec web site, www.eurogentec.com.

- Troubleshooting Guide for qPCR and RTqPCR (under the "Technical Resources / Troubleshooting Guide" section).

- Primers and probe design (please refer to our Troubleshooting Guide).

- "Your One-stop-shop Real-Time qPCR supplier" handbook (under the "Technical Resources / Documentation" section).

- MSDSs, (under the "Technical Resources / MSDS" section)

- Certificates of Analysis (please contact us).

For any further information required 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 66

FOR RESEARCH USE ONLY

NOTICE TO PURCHASER: LIMITED LICENSE

A license under U.S. Patents 4,683,202, 4,683,195 and 4,965,188 or their foreign counterparts, owned by Roche Molecular Systems, Inc. and F. Hoffmann-La Roche Ltd (Roche®), has an up-front fee component and a running-royalty component. The purchase price of this product includes limited, non-transferrable rights under the running-royalty component to use only this amount of the product to practice the Polymerase Chain Reaction ("PCR") and related processes described in said patents solely for the research and development activities of the purchaser when this product is used in conjunction with a thermal cycler whose use is covered by the up-front fee component. Rights to the up-front fee component must be obtained by the end user in order to have a complete license. These rights under the up-front fee component may be purchased from Applied Biosystems or obtained by purchasing an Authorized Thermal Cycler. No right to perform or offer commercial services of any kind using PCR, including without limitation reporting the results of purchaser's activities for a fee or other commercial consideration, is hereby granted by implication or estoppel. Further information on purchasing licenses to practice the PCR Process may be obtained by contacting the Director of Licensing at Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404 or the Licensing Department at Roche Molecular Systems, Inc., 1145 Atlantic Avenue, Alameda, California 94501.

The use of certain fluorogenic probes in 5' nuclease assays may be covered by U.S. Patents 5,210,015 and 5,487,972, owned by Roche Molecular Systems, Inc. and by U.S. Patent 5,538,848, owned by The Perkin-Elmer Corporation. The purchase of Eurogentec's product does not provide a license to use this patented technology. A license must be obtained by contacting the Director of Licensing Applied Biosystems, 850 Lincoln Centre Drive, Foster City, CA 94404 or the Licensing Department at Roche Molecular Systems Inc., 1145 Atlantic Avenue, Alameda, CA 94501.

Use of UDG employs U.S. Patents 5,035,996, 5,945,313, 5,683,896 and their foreign counterparts licensed to Eurogentec, S.A. from Invitrogen Corporation.

Primer Express® is a registered trademark of Applied Corporation.

qPCR Core kit Technical Data Sheet

Reference: RT-QP73-05

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

Storage conditions

For long term storage the qPCR Core kit should be stored at -15°C to -25 °C in a constant temperature freezer. When stored under these conditions the reagents are stable for 2 years.

For short term storage the qPCR Core kit 10x reaction buffer and MgCl₂ can be stored at 4 °C to 6 °C for 1 month. The Taq polymerase should be stored at -15 °C to -25 °C.

Kit contents

The qPCR Core kit contains enough reagents for up to 500 - 50 µl reactions using the hotstart enzyme, HotGoldStar.

Reagent	Volume	Description
10x Reaction buffer (red cap)	2.8 ml	Two tubes (1.4 ml) of reaction buffer, containing KCl and Tris-HCl and passive reference
50 mM MgCl ₂ (plain cap)	3 ml	Two tubes (1.5 ml) of 50 mM MgCl ₂
5 mM dNTP Mix (green cap)	1.25 ml	One tube of dATP, dCTP, dGTP, dTTP and dUTP in autoclaved, deionized water titrated with NaOH to pH 7.0
PCR enzyme (yellow cap)	125 µl	One tube containing 5 U/µl HotGoldStar

Procedure

- 1- Thaw all required reagents completely and put them on ice. Mix all reagents well by inversion and spin them down prior to pipeting.

2- Prepare the reaction mix

Component	Volume (µl)	Final Concentration
10x reaction buffer	5	1x
50 MgCl ₂	5	5 mM (or as required)
5 mM dNTP Mix	2	200 µM each dNTP
Forward primer	5	(starting with 300 nM)*
Reverse primer	5	(starting with 300 nM)*
Probe	5	(starting with 100 nM)*
HotGoldStar	0.25	0.025 U/µl
Template	5	-
Water	19.25	(volume is 50 µl minus all other components)
Total Mix	50 µl	

*Note 1: the primer and probe concentrations are recommended as starting concentrations. These concentrations will be correct for many assays, but additional optimization of the primer concentrations and primer-probe ratios may be required to obtain the best results with your primer-probe set.

Note 2: Uracil-N-Glycosylase (reference RT-0610-03) can be added to a final concentration of 0.01 u/µl (0.5 µl of 1 u/µl UNG per 50 µl reaction). If UNG is required, please add the following step in point 6 before the HotGoldStar activation step: 2 min at 50 °C.

- 3- To correct for dispensing losses prepare an excess of reaction mix (for example 100 reactions reaction mix for 96 reactions). Add all components together, except for the template. Mix thoroughly by inversion. Spin down.
- 4- Pipette 5 µl of the template DNA for your samples, 5 µl of the control DNA for your positive control and 5 µl of water or buffer for your negative control in to your PCR tubes / 96-well plate / 384-well plate.
- 5- Add 45 µl of the reaction mix to the reaction vial, close the vial and mix gently on a stirrer or spin down. Ensure that no bubbles are present in the reaction vial. Reaction set up can be done at room temperature.
- 6- Program the Real-Time thermocycler using the following recommended parameters:

HotGoldStar activation	10 min. 95 °C
40 Cycles	15 sec. 95 °C 1 min. 60 °C
Hold	50 °C forever

Technical information

Primer and probe design guidelines

Probes:

- avoid runs of identical nucleotides, especially of 4 or more Gs
- when using Primer Express® software, the T_m should be 65 °C to 67 °C
- avoid 5' end G (quenches the fluorophore)
- for genotyping, position of the polymorphism should be in the center of the probe, and the probe length should be adjusted such that both probes have the same T_m

Primers:

- GC content should be between 30 % and 80 % (ideally 40-60 %)
- avoid runs of identical nucleotides, especially of 3 or more Gs or Cs at the 3' end
- using the Primer Express® software the T_m should be 58 °C to 60 °C
- the primer should be placed as close as possible to the probe

Custom assay design

Commonly used concentrations are 300 nM for primers and 100 nM for probes. Optimal results may require titration of primers and probes or adjustment of the primer / probe ratio. The purpose of such a process is to determine the minimum amount of primers and probe required to obtain the most sensitive results with your assay.

Primer titration matrix

Titrate according to the Table 1, perform qPCR and select the concentration, which gives the lowest Ct value.

By doing this type of titration it is also possible to compensate for differences up to 2 °C in melt temperature of the primers.

Table 1: Primer titration matrix

Reverse	Forward		
	50 nM	300 nM	900 nM
50 nM	50 / 50	300 / 50	900 / 50
300 nM	50 / 300	300 / 300	900 / 300
900 nM	50 / 900	300 / 900	900 / 900

Primer-probe ratio matrix

Select optimal primer concentration as described in Table 1 and test with all probe concentrations described in Table 2. Select the concentration, which gives the lowest Ct value.

Table 2: Primer-probe ratio matrix

Opt. primers	Probe		
	50 nM	100 nM	250 nM
	50 / opt	100 / opt	250 / opt

MgCl₂ adjustment matrix

Standard MgCl₂ concentration is 3.5 mM but optimal MgCl₂ concentration can vary between assay. Always prefer optimizing the primer and probe concentrations before the MgCl₂ concentration.

Adjust the amount of water if MgCl₂ is added to the reaction.

Final MgCl ₂ concentration (mM)	MgCl ₂ to add (µl/50 µl)	10x reaction buffer (µl)
3	3	5
3.5	3.5	5
4	4	5
4.5	4.5	5
5	5	5
5.5	5.5	5
6	6	5
6.5	6.5	5
7	7	5
7.5	7.5	5
8	8	5

3-step protocol instead of 2-step protocol

Increasing extension time or performing a 3-step protocol can increase the ΔR_n and / or decrease the C_t of an assay, particularly when the PCR product is longer than 100 bp.

The protocol will be as follows:

HotGoldStar activation		10 min. 95 °C
40 Cycles	denaturation	15 sec. 95 °C
	annealing	20 sec. 60 °C
	extension	40 sec. 72 °C
Increase extension time with 10-second steps, if required.		

Further information available through Eurogentec web site, www.eurogentec.com.

- Manual for qPCR Core kit, reference RT-0000-04 (under the "Technical Resources / Manual" section).
- Troubleshooting Guide for qPCR and RTqPCR (under the "Technical Resources / Troubleshooting Guide" section).
- Primers and probe design (please refer to our Troubleshooting Guide).
- "Your One-stop-shop Real-Time qPCR supplier" handbook (under the "Technical Resources / Documentation" section).
- MSDSs, (under the "Technical Ressources / MSDS" section)
- Certificates of Analysis (please contact us).

For any further information required 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 66

FOR RESEARCH USE ONLY

NOTICE TO PURCHASER: LIMITED LICENSE

A license under U.S. Patents 4,683,202, 4,683,195 and 4,965,188 or their foreign counterparts, owned by Roche Molecular Systems, Inc. and F. Hoffmann-La Roche Ltd (Roche®), has an up-front fee component and a running-royalty component. The purchase price of this product includes limited, non-transferable rights under the running-royalty component to use only this amount of the product to practice the Polymerase Chain Reaction ("PCR") and related processes described in said patents solely for the research and development activities of the purchaser when this product is used in conjunction with a thermal cycler whose use is covered by the up-front fee component. Rights to the up-front fee component must be obtained by the end user in order to have a complete license. These rights under the up-front fee component may be purchased from Applied Biosystems or obtained by purchasing an Authorized Thermal Cycler. No right to perform or offer commercial services of any kind using PCR, including without limitation reporting the results of purchaser's activities for a fee or other commercial consideration, is hereby granted by implication or estoppel. Further information on purchasing licenses to practice the PCR Process may be obtained by contacting the Director of Licensing at Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404 or the Licensing Department at Roche Molecular Systems, Inc., 1145 Atlantic Avenue, Alameda, California 94501.

The use of certain fluorogenic probes in 5' nuclease assays may be covered by U.S. Patents 5,210,015 and 5,487,972, owned by Roche Molecular Systems, Inc. and by U.S. Patent 5,538,848, owned by The Perkin-Elmer Corporation. The purchase of Eurogentec's product does not provide a license to use this patented technology. A license must be obtained by contacting the Director of Licensing Applied Biosystems, 850 Lincoln Centre Drive, Foster City, CA 94404 or the Licensing Department at Roche Molecular Systems Inc., 1145 Atlantic Avenue, Alameda, CA 94501.

Use of UDG employs U.S. Patents 5,035,996, 5,945,313, 5,683,896 and their foreign counterparts licensed to Eurogentec, S.A. from Invitrogen Corporation.

Primer Express® is a registered trademark of Applied Biosystems Corporation.