

β-2 microglobulin Control kit Technical Data Sheet

Reference: RT-CKYD-B2M

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

Storage conditions

The β-2 microglobulin Control 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 1 year.

Kit contents

The β-2 microglobulin Control kit contains enough reagents for up to 250 - 50 µl reactions.

Reagent	Volume	Description
B2M primer mix (plain cap)	750 µl	One tube of primer mix 10 µM
B2M probe (brown cap)	250 µl	One tube of 12.5 µM Yakima Yellow®-Eclipse® DarkQuencher probe

The β-2 microglobulin Control kit Yakima Yellow®-Eclipse® DarkQuencher should be combined with Eurogentec Reverse Transcriptase core kit (first strand reaction kit) reference RT-RTCK-05 followed by Eurogentec qPCR MasterMixes, reference RT-QP2X-03 or qPCR Core kits, reference RT-QP73-05. Alternatively a One step RTqPCR MasterMix Plus, reference RT-QPRT-032X can also be combined with the Control kit. Alternatively a qPCR MasterMix (RT-QP2X-03) can be combined with the control kit for quantitation of gDNA.

The kit is available with a Yakima Yellow®-Eclipse® DarkQuencher labelled probe detected in channel 2 for multiplex qPCR.

Procedure

We strongly recommend preparation of a reaction mix

Each 50 µl reaction requires the addition of 3 µl of B2M Primer Mix (plain cap) and 1 µl of B2M Probe (brown cap).

If required a mix containing the appropriate amount of both primer and probe can be prepared. Dispense 4 µl of this mix per 50 µl reaction.

Final concentration of primers is 600 nM.

Final concentration of probe is 250 nM.

Prepare an appropriate dilution series of chromosomal DNA in TE or PCR grade water and a no template control.

The β-2 microglobulin Control Kit has been developed to detect the human B2M gene (cDNA/ gDNA). The obtained amplicon is 112 bp in length, to guarantee an efficient PCR.

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 at 25 °C forever	

If UNG is used, program the Real-Time thermocycler using the following recommended parameters:

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

Other Eurogentec Control Kits available:

Eurogentec has developed control kits that can be used for either cDNA or cDNA and gDNA.

18S rRNA Control Kit
human, rat, mouse, rabbit cDNA and gDNA
very high expression level
RT-CKYD-18S or RT-CKFT-18S

23S rRNA Control kit
E. coli
very high expression level
RT-CKFT-23S

28S rRNA Control Kit
human, mouse cDNA and gDNA
very high expression level
RT-CKYD-28S

β -actin Control Kit
human cDNA and gDNA
high expression level
RT-CKYD-ACTB

GAPDH Control kit
human cDNA
high expression level
RT-CKYD-GAPD

β -2-microglobulin Control Kit
human cDNA and gDNA
high expression level
RT-CKYD-B2M

Ubiquitin C Control kit
human cDNA and gDNA
high expression level
RT-CKYD-UBC

Phospholipase A2 Control Kit
human cDNA
medium expression level
RT-CKYD-YWHAZ

Hypoxanthine phosphoribosyltransferase Control Kit
human cDNA
medium expression level
RT-CKYD-HPRT1

Ribosomal protein L13a Control Kit
human cDNA
medium expression level
RT-CKYD-RPL13A

Succinate dehydrogenase complex Control Kit
Human cDNA
medium expression level
RT-CKYD-SDHA

TATA box binding protein Control kit
human cDNA
low expression level
RT-CKYD-TBP

Hydroxymethyl-bilane synthase Control Kit
human cDNA
low expression level
RT-CKYD-HMBS

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 TaqMan fluorogenic probes in 5' nuclease assays is covered by US patents 5,210,015 and 5,487,972, owned by Roche Molecular Systems, Inc. and by US patent 5,538,848, owned by The Perkin-Elmer Corporation. Purchase of this product does not provide a license to use this patented technology. A license to practice this technology must be obtained from PE Biosystems, 850 Lincoln Centre Drive, Foster City, CA 94404, or from Roche Molecular Systems Inc., 1145 Atlantic Avenue, Alameda, CA 94501.

Use of UDG employs U.S. patents 5,035,996, 5,946,313, 5,683,896 and their foreign counterparts licensed to Eurogentec, S.A. from Invitrogen Corporation.

Eclipse® is a registered trademark of Epoch

Yakima Yellow® is a registered trademark of Epoch Biosciences, Inc.