

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

Ubiquitin C Control kit Technical Data Sheet

Reference: RT-CKYD-UBC

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

Storage conditions

The Ubiquitin C 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 Ubiquitin C Control kit contains enough reagents for up to $250 - 50 \ \mu l$ reactions.

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

The Ubiquitin C 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 UBC Primer Mix (plain cap) and 1 µl of UBC Probe (brown cap).

If required a mix containing the appropriate amount of both primer and probe can be prepared. Dispense 4 μ l of this mix per 50 μ l 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 Ubiquitin C Control Kit has been developed to detect the human UBC gene (cDNA/ gDNA). The obtained amplicon is 76 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 incativation	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 developped control kits that can be used for either cDNA or cDNA and gDNA.

18S rRNA Control Kit

human, rat, mouse, rabbit cDNA an gDNA very high expression level

RT-CKYD-18S or RT-CKFT-18S

23S rRNA Control kit

F.coli

very high expression level

RT-CKFT-23S

28S rRNA Control Kit

human, mouse cDNA and gDNA very high expression level

RT-CKYD-28S

ß-actin Control Kit

human cDNA and gDNA high expression level RT-CKYD-ACTB

GAPDH Control kit

human cDNA

high expression level

RT-CKYD-GAPD

ß-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 gPCR and RTgPCR (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 LISE ONLY

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 Luf (Roche*), has an up-front fee component and art anning-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 (PCRT) 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 must be obtained by the rod with user in order to have a complete license. These rights under the up-front fee component must be obtained by the up-front fee component must b

The use of TagMan fluorogenic probes in 5' nuclease assays is covered by US patents 5.210,015 and 5.487,972, owned by Molecular Systems, inc., and by US patent 5.538,848, owned by The Perkin-Elmer Corporation. Purchase of this product do provide a license to use this patented technology. A license to practise this technology must be obtained from PE Biosystems Lincoh Centre Drive, Foster Chy, CA 94404. or from Roche Molecular Systems inc., 1145 Atalantic Avenue, Almeda, 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.