

**Instruction Manual** 

# LUX<sup>™</sup> Fluorogenic Primers

#### For real-time PCR and RT-PCR

**Version F** 13 September 2004 25-0546

### **Table of Contents**

Introduction	1
Designing and Ordering Custom LUX <sup>™</sup> Primers	3
Storing and Reconstituting Primers	5
Certified LUX <sup>™</sup> Primer Sets for Housekeeping Genes	6
Real-Time qPCR	7
Multiplex Real-Time qPCR	11
Two-Step Real-Time qRT-PCR	12
One-Step Real-Time qRT-PCR	16
Troubleshooting	21
Accessory Products	23
Purchaser Notification	24
Technical Service	25
References	27

#### Introduction

#### Overview

LUX<sup>™</sup> (Light Upon eXtension) Primers are an easy to use, highly sensitive, and efficient method for performing real-time quantitative PCR (qPCR) and RT-PCR (qRT-PCR). Each primer pair in the LUX<sup>™</sup> system includes a fluorogenic primer with a fluorophore attached to its 3' end and a corresponding unlabeled primer. The fluorogenic primer has a short sequence tail of 4–6 nucleotides on the 5' end that is complementary to the 3' end of the primer. The resulting hairpin secondary structure provides optimal quenching of the fluorophore (see the figure below). When the primer is incorporated into double-stranded DNA during PCR, the fluorophore is dequenched and the signal increases by up to 10-fold.

LUX<sup>™</sup> Primers combine high specificity with simple design and streamlined protocols. LUX<sup>™</sup> Primers require no special probes or quenchers and are compatible with melting curve analysis of real-time qPCR products, which allows for the differentiation of amplicons and primer dimer artifacts by their melting temperatures. LUX<sup>™</sup> Primers are available with two different reporter dyes, which provides for multiplexing capability. You can custom-design LUX<sup>™</sup> Primers<sup>™</sup> from a target DNA sequence using Web- or desktop-based software, or order predesigned and validated Certified LUX<sup>™</sup> Primer Sets for Housekeeping Genes.



Labeling

Each fluorogenic LUX<sup>™</sup> primer is labeled with one of two reporter dyes—FAM (6-carboxy-fluorescein) or JOE (6-carboxy-4', 5'-dichloro-2', 7'-dimethoxy-fluorescein).

Applications	LUX <sup>™</sup> Primers can be used in real-time quantitative PCR and RT-PCR to quantify 100 or fewer copies of a target gene in as little as 1 pg of template DNA or RNA. They have a broad dynamic range of 7–8 orders. See the guidelines and sample protocols for qPCR on pages 7–10 and guidelines and sample protocols for qRT-PCR on pages 12–20. Multiplex applications use separate FAM and JOE-labeled primer sets to detect two different genes in the same sample. Typically, a custom-designed FAM-labeled primer set would be used to detect the gene of interest, and a JOE-labeled Certified LUX <sup>™</sup> Primer Set would be used to detect a housekeeping gene as an internal control. See the optimization guidelines for multiplex qPCR on page 11.
Instrument Compatibility	LUX <sup>™</sup> Primers are compatible with a wide variety of real-time qPCR instruments, including but not limited to the ABI PRISM <sup>®</sup> 7700, 7000, and 7900 and GeneAmp <sup>®</sup> 5700; the Bio-Rad iCycler <sup>™</sup> ; the Stratagene Mx4000 <sup>™</sup> and Mx3000 <sup>™</sup> ; the Cepheid Smart Cycler <sup>®</sup> ; the Corbett Research Rotor-Gene; and the Roche LightCycler <sup>®</sup> .
	At a minimum, the real-time qPCR instrument should:
	Detect fluorescence at each PCR cycle
	<ul> <li>Excite and detect FAM-labeled LUX<sup>™</sup> Primers near their excitation/emission wavelengths of 490/520 nm, and/or</li> </ul>
	<ul> <li>Excite and detect JOE-labeled LUX<sup>™</sup> Primers near their excitation/emission wavelengths of 520/550 nm</li> </ul>
	Refer to the specific instrument's user manual for operating instructions.
	ABI PRISM is a registered trademark of Applera Corporation. GeneAmp is a registered trademark of Roche Molecular Systems, Inc. LightCycler is a registered trademark of Idaho Technologies, Inc. iCycler, Mx4000, Mx3000, Rotor-Gene, and Smart Cycler are trademarks of their respective companies.

## Designing and Ordering Custom LUX<sup>™</sup> Primers

LUX <sup>™</sup> Designer Primer Design Software	To design and order custom LUX <sup>™</sup> Primers for your genes of interest, visit the Invitrogen LUX <sup>™</sup> Web site at <u>www.invitrogen.com/LUX</u> and follow the link to the LUX <sup>™</sup> Designer software. The software is available as either a Web-based application or a Microsoft <sup>®</sup> Windows <sup>®</sup> -compatible download. Follow the step-by-step instructions in the software to submit your target sequence and generate primer designs. LUX <sup>™</sup> Designer will automatically generate one or more primer designs based on each sequence you submit and the selected design parameters. The design software includes algorithms to minimize primer self-complementarity and interactions between primers. It also assigns rankings to the generated designs—based on primer melting temperature, hairpin structure, self-annealing properties, etc.—to aid in selection. When the designs have been generated, you can review them, select a design, select the fluorophore labels, and place your order.
Guidelines for Submitting a Target Sequence	<ul> <li>When you submit a target sequence containing your gene of interest, keep in mind the following design criteria:</li> <li>The optimal amplicon length for real-time qPCR ranges from 80 to 200 bases. You can specify a minimum, optimal, and maximum amplicon length when you submit the sequence.</li> <li>The target sequence should be at least 10 bases longer than the minimum amplicon size you select. The longer the sequence, the more likely that an optimal primer design can be developed.</li> <li>The sequence must contain only standard IUPAC (International Union of Pure and Applied Chemistry) letter abbreviations.</li> <li>When you select the design parameters, the default melting temperature range is 60–68°C. Do not change this default unless the design engine finds no primers in this range. For primers in this range, PCR annealing temperatures from 55° to 64°C are appropriate.</li> </ul>
_	When you first submit a sequence, the Disable Score-Based Rejection checkbox should <i>not</i> be checked; the resulting scores provide an important measure of primer suitability. <b>Scores in the range of 0.0–4.0 are acceptable.</b> If no primers with a score of 4.0 or lower can be generated from a sequence, you can disable score-based rejection and redesign the primers. See the LUX <sup>™</sup> Designer <b>Help</b> for additional guidance.
Selecting a Primer Design	After you submit your sequence, LUX <sup>™</sup> Designer will first generate one or more designs for the labeled primer. The labeled primer can be either the forward or the reverse primer. After you select a design for the labeled primer, you will be prompted to select a design for the corresponding unlabeled primer.

## **Designing and Ordering Custom LUX<sup>™</sup> Primers**, Continued

Selecting Labels	After you have selected a primer set (labeled and unlabeled) for a particular sequence, you can specify the particular label and synthesis scale. Custom LUX <sup>™</sup> Primers are provided in 50 nM or 200 nM synthesis scale. When selecting labels in a multiplex reaction, we recommend using the FAM label for your gene of interest and the JOE label for the housekeeping gene that you will use as the internal control. Certified LUX <sup>™</sup> Primer Sets for Housekeeping Genes are recommended for the JOE-labeled control gene.
Placing the Order	After you have selected the label and synthesis scale, you can submit your order to Invitrogen using the Web site or by e-mail or fax. Each primer order will be shipped directly from Invitrogen's Custom Primer Facilities. Labeled primers are supplied in an amber tube; unlabeled primers are supplied in a clear tube.
	Each primer ordered from Invitrogen's Custom Primer Facilities comes with a Certificate of Analysis (COA) verifying the amount and sequence.
Product Qualification	Custom LUX <sup>™</sup> Primers are tested post-synthesis by optical density (OD) ratio measurements and mass spectroscopy to ensure efficient dye labeling and correct molecular weight and composition.
	See the Certificate of Analysis shipped with each primer for more information.

## Storing and Reconstituting Primers

Primer Storage and Stability	Store primers at −20°C in the dark. LUX <sup>™</sup> Primers are stable for:		
otability	• >12 months when stored at $-20^{\circ}$ C in lyophilized form.		
	• >6 months when stored at $-20^{\circ}$ C in solution.		
	Stability can be extended by storing at –70°C.		
	Be careful to minimize the exposure of labeled LUX <sup>™</sup> Primers to direct light, as this can reduce their fluorescent intensity.		
Reconstituting Primers	Custom LUX <sup>™</sup> Primers are provided lyophilized in 50-nmole or 200-nmole synthesis scale. To reconstitute primers, centrifuge the tube for a few seconds to collect the oligonucleotide in the bottom of the tube. Carefully open, add an appropriate volume of TE buffer or ultrapure water, close the tube, rehydrate for 5 minutes, and vortex for 15 seconds.		
	We recommend that you rehydrate primers at concentrations greater than 10 $\mu$ M. To prepare a 100 $\mu$ M primer stock solution, multiply the primer amount in nmoles by ten to determine the volume of diluent in $\mu$ l.		
	After reconstitution, store the primer stock at $-20^{\circ}$ C in the dark, where it will be stable for 6 months or more.		

## Certified LUX<sup>™</sup> Primer Sets for Housekeeping Genes

Certified LUX™Certified LUX™ Primer Sets for Housekeeping Genes are predesigned primerPrimer Sets forSets for genes that are commonly used as internal controls for normalizingHousekeepingreal-time qRT-PCR experiments. These primer sets have been optimized and<br/>functionally validated to provide accurate, reproducible results using standard<br/>LUX™ protocols. They are supplied ready to use in TE buffer.

Each Certified LUX<sup>TM</sup> Primer Set includes a FAM- or JOE-labeled LUX<sup>TM</sup> primer and a corresponding unlabeled primer. Each primer (labeled and unlabeled is supplied at 100  $\mu$ l and a concentration of 10  $\mu$ M. Available sets are listed below. For additional information, visit <u>www.invitrogen.com/LUX</u>.

Product	GenBank Accession no.	Forward/ Reverse Label	Cat. no. FAM label	Cat. no. JOE label	Relative expression	CDS Location	PCR Product Size Range
Human genes					•		<u> </u>
18S rRNA	X03205	Forward	115HM-01	115HM-02	++++	n/a	101-150 bp
hβ-ACTIN	NM_001101	Forward	101H-01	101H-02	+++	Exons 2/3	101-150 bp
hATPSase	NM_001686	Forward	108H-01	108H-02	+++	n/a	101-150 bp
hB2M	NM_004048	Forward	113H-01	113H-02	+++	Exons 1/2	101-150 bp
hGAPDH	NM_002046	Forward	100H-01	100H-02	+++	Exons 4/5	151–200 bp
hPGK1	NM_000291	Forward	109H-01	109H-02	+++	n/a	50-100 bp
hPPIA	NM_021130	Forward	106H-01	106H-02	+++	Exons 2/3	50-100 bp
hRPL4	NM_000968	Reverse	103H-01	103H-02	+++	Exons 8/9	101-150 bp
hEEF1G	NM_001404	Forward	107H-01	107H-02	++	n/a	50-100 bp
hHPRT1	NM_000194	Reverse	105H-01	105H-02	++	Exons 5/6	50-100 bp
hSDHA	NM_004168	Forward	102H-01	102H-02	++	Exons 12/13	50-100 bp
hTFRC	NM_003234	Forward	111H-01	111H-02	++	Exons 10/11	101-150 bp
hGUS	NM_000181	Forward	112H-01	112H-02	+	Exons 7/8	101-150 bp
hHMBS	NM_000190	Forward	110H-01	110H-02	+	Exons 2/3	50-100 bp
hTBP	NM_003194	Forward	104H-01	104H-02	+	Exons 3/4	101-150 bp
hUBC	NM_021009	Forward	114H-01	114H-02	+	n/a	50-100 bp
Mouse/rat gene	s	1			1		
18S rRNA	X03205	Forward	115HM-01	115HM-02	++++	n/a	101-150 bp
mβ-ACTIN	NM_007393	Forward	101M-01	101M-02	+++	Exons 2/3	101-150 bp
mB2M	X01838	Forward	113M-01	113M-02	+++	n/a	50-100 bp
mEEF1G	AF321126	Forward	107M-01	107M-02	+++	n/a	101-150 bp
mGAPDH	NM_008084	Forward	100M-01	100M-02	+++	Exons 4/5	151–200 bp
mPGK1	NM_008828	Forward	109M-01	109M-02	+++	Exons 1/2	101-150 bp
mPPIA	NM_008907	Reverse	106M-01	106M-02	+++	Exons 1/2	50-100 bp
mRPL4	NM_022510	Forward	103M-01	103M-02	+++	Exons 2/3	151-200 bp
mHPRT1	NM_013556	Forward	105M-01	105M-02	++	Exons 6/7	50-100 bp
mSDHA	AF095938	Forward	102M-01	102M-02	++	Exons 6/7	50-100 bp
mATPSase	NM_016774	Forward	108M-01	108M-02	+	n/a	50-100 bp
mGUS	NM_010368	Forward	112M-01	112M-02	+	Exons 7/8	50-100 bp
mHMBS	XM_129404	Reverse	110M-01	110M-02	+	Exons 4/5	50-100 bp
mTBP	NM_013684	Forward	104M-01	104M-02	+	Exons 3/4	101-150 bp
mTFRC	NM_011638	Forward	111M-01	111M-02	+	Exons 2/3	101-150 bp
Drosophila gen	es	Γ					
d18S rRNA	AY037174	Reverse	115D-01	115D-02	++++	n/a	101-150 bp
dActin	NM_079486	Forward	101D-01	101D-02	+++	Exons 2/3	50–100 bp

## Real-Time qPCR

Introduction	Real-time qPCR uses genomic DNA, cDNA, or plasmid DNA as starting material. This section provides guidelines and example protocols for performing real-time qPCR using LUX <sup>™</sup> Primers.
Template Specifications	The target template for real-time qPCR is linear single-stranded or double-stranded DNA, cDNA, or circular DNA (such as plasmids) that has been linearized. The amount of DNA typically ranges from $10^2$ to $10^7$ copies or 1 pg to 10 µg of template.
_	See page 12 for instructions on generating cDNA using reverse transcription as part of two-step real-time qRT-PCR.
Primer Concentration	For optimal PCR conditions, primer titrations of 50–500 nM per primer are recommended. The sample reactions on pages 9–10 use 200 nM of each primer, equivalent to 1 $\mu$ l of a 10- $\mu$ M primer solution in a 50- $\mu$ l reaction.
Magnesium Concentration	The optimal Mg <sup>++</sup> concentration for a given target/primer/polymerase combination can vary between 1 mM and 10 mM, but is usually in the range of 3 mM. See the sample reactions on pages 9–10.
dNTP Concentration	The optimal concentration of dATP, dCTP, dGTP, and dTTP is 200 $\mu$ M each. If dUTP is used in place of dTTP, its optimal concentration is 400 $\mu$ M.
DNA Polymerase	We recommend using a "hot-start" DNA polymerase, preferably one that has been optimized for real-time qPCR. Platinum <sup>®</sup> Quantitative PCR SuperMix UDG (Catalog no. 11730-017) is a 2X-concentrated, ready-to-use mixture containing all components except primers and template. It uses Platinum <sup>®</sup> <i>Taq</i> DNA polymerase and has been specifically formulated to provide optimal performance in real-time qPCR systems.
ROX Reference Dye	We recommend using ROX Reference Dye (Cat. no. 12223-023) to normalize the fluorescent reporter signal in real-time qPCR for instruments that are compatible with this option. ROX Reference Dye can be used to adjust for non- PCR-related fluctuations in fluorescence between reactions, and provides a stable baseline in multiplex reactions. It is composed of a glycine conjugate of 5-carboxy-X-rhodamine, succinimidyl ester (25 $\mu$ M) in 20 mM Tris-HCl (pH 8.4), 0.1 mM EDTA, and 0.01% Tween <sup>®</sup> 20.
	ROX is supplied at 50X concentration. Add 1 $\mu$ l of ROX for every 50 $\mu$ l of reaction volume. For convenience and to reduce pipetting errors, you can premix a solution of ROX and Platinum <sup>®</sup> Quantitative PCR SuperMix-UDG. Add 1 $\mu$ l of ROX for every 25 $\mu$ l of SuperMix-UDG. Store mixture at either -20°C or 4°C in the dark.

### Real-Time qPCR, Continued

Bovine Serum Albumin (BSA)	Bovine serum albumin (BSA) is required in qPCR reactions on the Roche LightCycler <sup>®</sup> because the glass capillaries in the LightCycler <sup>®</sup> have a high surface-to-volume ratio and the glass surface binds molecules like <i>Taq</i> DNA polymerase, effectively removing them from the reaction. The addition of BSA blocks this surface binding.			
	Nonacetylated BSA is strongly recommended because acetylated BSA will inhibit PCR at the concentrations required in LightCycler <sup>®</sup> reactions. This inhibition is most likely due to the transfer of acetyl groups to essential components of the PCR, like the <i>Taq</i> DNA polymerase.			
_	To ensure that the BSA does not contain RNA or DNA, we recommend using ultrapure, molecular biology-grade nonacetylated BSA from Panvera/Invitrogen (Cat. nos. P2489 and P2046).			
Melting Curve Analysis	Melting curve analysis is strongly recommended during qPCR to identify the presence of primer dimers and analyze the specificity of the reaction. Program your real-time instrument to perform this analysis after thermocycling.			
	Melting curve analysis identifies the change in fluorescent signal that occurs as double-stranded DNA (dsDNA) dissociates or "melts" into single-stranded DNA. By identifying the temperature at which the dsDNA dissociates, you can distinguish smaller artifacts like primer dimers with a lower annealing temperature from larger amplicons with a higher annealing temperature.			
	The presence of primer dimers in samples containing template decreases PCR efficiency and obscures analysis and determination of cycle thresholds. The formation of primer dimers most often occurs in no-template controls, where the polymerase enzyme is essentially idle, and in this case the quantitative analysis of the template samples is not affected. Melting curve analysis of no-template controls can discriminate between primer dimers and spurious amplification due to contaminating nucleic acids in reagent components.			
Instrument Settings	Follow the manufacturer's instructions for configuring your real-time qPCR instrument for use with LUX <sup>™</sup> Primers. Note the following general settings:			
	• Program your instrument to perform melting curve analysis at the end of thermocycling, if this option is available.			
	• The quencher setting on the instrument should reflect the fact that LUX <sup>™</sup> Primers do not contain a quencher.			
	• We recommend using ROX Reference Dye, if your instrument is compatible with this option. Adjust the instrument settings accordingly.			
	Additional guidelines and settings for specific instruments are available at <u>www.invitrogen.com/lux</u> ; click on <b>Instrument Protocols</b> .			
=				

### Real-Time qPCR, Continued

Protocol for Instruments Using PCR Tubes or Plates	The following protocol uses Platinum <sup>®</sup> Quar ROX reference reagent. It has been optimized instruments that use PCR tubes or plates. A LightCycler <sup>®</sup> is provided on the following p <b>Note:</b> The following protocol uses a 50-µl re may be used, depending on the requirement proceeding, see the real-time qPCR guideling multiplex reactions, see the guidelines on particular 1. To reduce well-to-well variation, preparticular ingredients except template. The follow volumes for one reaction and 50 reaction	ntitative PCR SuperMix-UDG with ed for use with real-time qPCR protocol for the Roche age. action volume; smaller volumes ts of your instrument. Before uses on the previous pages. For age 11. re a Master Mix of all the reaction <i>r</i> ing table provides Master Mix ns (scale up or down as needed):
	Component Platinum <sup>®</sup> Quantitative PCR SuperMix-UDC ROX Reference Dye (optional) Labeled LUX <sup>™</sup> Primer (10 µM) Unlabeled primer (10 µM) Sterile distilled water <sup>2</sup> <sup>1</sup> Final concentration: 0.03 U/µl Platinum <sup>®</sup> Taq (pH 8.4), 50 mM KCl, 3 mM MgCl <sub>2</sub> , 200 µM dC 400 µM dUTP, 1 U UDG <sup>2</sup> or use DNase/RNase Free Distilled Water (C	Vol/1 rxn         Vol/50 rxns           G <sup>1</sup> 25 μl         1250 μl           1 μl         50 μl           to 40 μl         to 2000 μl           DNA polymerase, 20 mM Tris-HCl           GTP, 200 μM dATP, 200 μM dCTP,           rat. No. 10977-015).
	<ol> <li>Program the real-time qPCR instrument</li> </ol>	t as follows:
	<b>3-Step Cycling (recommended)</b> $50^{\circ}$ C, 2 min hold (UDG treatment) $95^{\circ}$ C, 2 min hold 45 cycles of: $95^{\circ}$ C, 15 s $55^{\circ}$ C, 30 s $72^{\circ}$ C, 30 s	<b>2-Step Cycling (optional)</b> 50°C, 2 min hold (UDG treatment) 95°C, 2 min hold 45 cycles of: 95°C, 15 s 60-65°C, 30-45 s
	Melting Curve Analysis Refer to instrument docu	(recommended)
	<ol> <li>Add 40 μl of the Master Mix to an optica 96-well PCR plate.</li> </ol>	al PCR tube or each well of a
	4. Add 10 $\mu$ l of template in TE or sterile dE seal the tube/plate.	$H_2O$ to each reaction vessel. Cap or
	5. Gently mix and make sure that all comp tube/plate wells. Centrifuge briefly if n	ponents are at the bottom of the needed.
	6. Place reaction in the real-time qPCR inst Collect and analyze results.	trument and run the program.
		Continued on next page

### Real-Time qPCR, Continued

Protocol for the Roche LightCycler <sup>®</sup>	The following protocol uses Platinum <sup>®</sup> Quantitative PCR SuperMix-UDG and has been optimized for the Roche LightCycler <sup>®</sup> . Consult the LightCycler <sup>®</sup> documentation for detailed instructions on preparing the capillary tubes and operating the instrument. FAM-labeled LUX <sup>™</sup> Primers are also compatible with Roche enzyme mixes.				
	No of use gu	<b>ote:</b> JOE-labeled LUX <sup>™</sup> Primers are no the LightCycler <sup>®</sup> ; use FAM-labeled pr es a 20-μl reaction volume. Before pro idelines on the previous pages.	t compatil imers onl ceeding, s	ble with the c y. The follow see the real-tin	urrent version ing protocol me qPCR
	1.	To reduce well-to-well variation, pro- ingredients except template. The fol reaction and 34 reactions (scale as ne	epare a M lowing tal eeded):	aster Mix of a ble provides	all the reaction volumes for one
	Co	omponent		<u>Vol/1 rxn</u>	<u>Vol/34 rxns</u>
	Pla	atinum <sup>®</sup> Quantitative PCR SuperMix-	$UDG^1$	10 µl	340 µl
	FA	M-labeled LUX <sup>™</sup> Primer (10 µM)		1 µl	34 µl
	Ur	nlabeled primer (10 μM)		1 µl	34 µl
	Во	wine serum albumin $(5 \text{ mg/ml})^2$		1 µl	34 µl
	Pla	atinum <sup>®</sup> Taq DNA Polymerase <sup>3</sup>		0.12 μl	4 µl
	Ste	erile distilled water <sup>4</sup>		to 18 μl	to 612 μl
	<sup>1</sup> Final concentration: 0.03 U/ $\mu$ l Platinum <sup>®</sup> <i>Taq</i> DNA polymerase, 20 mM Tris-HCl (pH 8.4), 50 mM KCl, 3 mM MgCl <sub>2</sub> , 200 $\mu$ M dGTP, 200 $\mu$ M dATP, 200 $\mu$ M dCTP, 400 $\mu$ M dUTP, 1 U UDG				
		nos. P2489 and P2046).	e BSA (10%)	solution) from	Panvera (Cat.
		<sup>3</sup> Total units of Platinum <sup>®</sup> <i>Taq</i> DNA Polym from Platinum <sup>®</sup> Quantitative PCR SuperM	nerase in th ⁄lix-UDG)	e reaction is 1.2	2 (including 0.6 U
		<sup>4</sup> or use DNase/RNase Free Distilled Wate	er (Cat. No	. 10977-015).	
	2.	Set the fluorescence on the Roche Li	ghtCycler	® to the F1 ch	annel.
	3.	Program the instrument as follows:			
		Thermal Cycling	Melting	Curve Analy	ysis (optional)
		Program choice: Amplification Analysis mode: Quantification Cycling: 50°C, 2 min hold (UDG treatment) 95°C, 2 min hold 45 cycles of: 94°C, 5 s 55°C, 10 s (single acquire) 72°C, 10 s	<i>Program</i> <i>Analysis</i> Cycling: 95°C, 0 s 55°C, 15 95°C, 0 40°C, 0 s	<i>choice:</i> Meltin <i>mode:</i> Meltin sec (increase 0. continuous	ng curve g curves 1°C/s with acquisition)
	4.	Add 18 µl of Master Mix to each cap	illary tub	e of the Light	Cycler <sup>®</sup> .
	5.	Add 2 $\mu$ l of template to each tube, an	nd cap the	e tube.	
	6. Centrifuge the tubes at $700 \times g$ for 5 seconds.				
	7.	Place the reaction tubes in the rotor	of the Lig	htCvcler <sup>®</sup> and	d run the

 Place the reaction tubes in the rotor of the LightCycler<sup>®</sup> and rur program. Collect and analyze results.

## Multiplex Real-Time qPCR

Multiplex Real-Time qPCR	In multiplex real-time qPCR, different sets of primers with different fluorogenic labels are used to amplify separate genes in the template DNA. Multiplexing with LUX <sup>™</sup> Primers offers simplified PCR kinetics and increased reaction efficiency when compared with probe-based technologies, because only two oligos are used per target.				
LUX <sup>™</sup> Primers have been tested in multiplex reactions using a FAM- primer set for the gene of interest and a JOE-labeled set for a houseke gene used as an internal control to normalize between different react recommend using Certified LUX <sup>™</sup> Primer Sets for Housekeeping Gen internal control (see page 6).					
	Note: We recommend selecting a housekeeping gene expression level of your gene of interest. The relative predesigned, certified LUX <sup>™</sup> Primer Sets are shown o	that matches the relative expression levels of n page 6.			
	In a standard multiplex reaction, you can include the additional primers at the same volumes and concentration as the primers in a singleplex reaction, as shown in the example mixture below:				
	ComponentPlatinum® Quantitative PCR SuperMix-UDG (2X)ROX Reference Dye (50X)TemplateForward primer 1 (FAM label) (10 μM)Reverse primer 1 (10 μM)Forward primer 2 (JOE label) (10 μM)Reverse primer 2 (10 μM)Sterile distilled waterAll other reaction volumes remain the same.Follow the thermal cycling guidelines provided in PrUsing PCR Tubes or Plates on page 9. If you have di	Volume $25 \mu$ l $1 \mu$ l $10 \mu$ l $1 \mu$ lto 50 μl			
	multiplex reaction using these guidelines, see the opt	imization hints below.			
Optimizing Multiplex	If you notice a decline in PCR efficiency in your multiplex real-time qPCR, you can optimize the reaction by performing the steps listed below.				
Conditions	<b>Note:</b> We recommend that you perform one optimization step and then repeat the reaction to test for efficiency before moving on to the next step:				
	1. Reduce the primer concentration of the gene with the highest expression levels (typically the housekeeping gene) to $1/2$ the primer concentration of the other gene. For example, in a standard 50-µl reaction, you would add the primers for the less abundant gene at 1 µl each, and add the primers for the more abundant gene at 0.5 µl each.				
	. Increase the $MgCl_2$ in the reaction from 3 mM to 6 mM.				
	<ol> <li>Double the amount of polymerase enzyme (to 0.06 U/μl of reaction volume). If you are using Platinum<sup>®</sup> Quantitative PCR SuperMix-UDG, add Platinum<sup>®</sup> <i>Taq</i> DNA polymerase stand-alone enzyme (Catalog no. 10966-018) to double the amount of enzyme.</li> </ol>				
	4. Increase the dNTP concentrations in the reaction	to 400 μM each.			

### Two-Step Real-Time qRT-PCR

Introduction	Real-time qRT-PCR uses RNA as starting material in a reverse transcription reaction to generate first-strand cDNA. The cDNA is then quantified in a separate real-time qPCR reaction.		
	In two-step qRT-PCR, first-strand synthesis is performed, and then the reaction is transferred to a separate tube for the qPCR reaction. This section provides guidelines for two-step qRT-PCR and an optimized protocol using the SuperScript <sup>™</sup> III Platinum <sup>®</sup> Two-Step qRT-PCR Kit.		
Note	For the real-time qPCR portion of the two-step protocol, see additional guidelines on primers, magnesium, dNTPs, ROX Reference Dye, BSA, melting curve analysis, and instrument settings on pages 7–8.		
Template Specifications	The target template for two-step qRT-PCR is total RNA or mRNA. High- quality, intact RNA is essential for full-length, high-quality cDNA synthesis and accurate quantification. Starting material can range from 10 pg to 1 $\mu$ g total RNA. Then use 1 pg to 10 $\mu$ g of the cDNA from the first-strand reaction in the qPCR step.		
	The purity and integrity of the starting RNA have a direct impact on results. RNase and genomic DNA contamination are the most common problems, and purification methods should include RNase inhibitors and DNase digestion to minimize these.		
	We recommend using the Micro-to-Midi Total RNA Purification System (Catalog no. 12183-018) or TRIzol <sup>®</sup> reagent (Catalog no. 15596-026) to isolate total RNA. High-quality total RNA can be purified from as little as 100 cells up to $10^7$ cells or 200 mg of tissue.		
	Isolation of mRNA is typically not required, but can be performed using the FastTrack <sup>®</sup> 2.0 mRNA Isolation Kit (Catalog no. K1593-02).		
Enzyme Specifications	For two-step qRT-PCR, we recommend using a high-specificity, high-yield reverse transcriptase such as SuperScript <sup>™</sup> III Reverse Transcriptase and a "hot-start" DNA polymerase such as Platinum <sup>®</sup> <i>Taq</i> DNA Polymerase. The SuperScript <sup>™</sup> III Platinum <sup>®</sup> Two-Step qRT-PCR Kit (Catalog nos. 11734-050 and 11734-068) includes SuperScript <sup>™</sup> III RT, Platinum <sup>®</sup> <i>Taq</i> DNA Polymerase, and all the other necessary components for two-step qRT-PCR except the RNA. See the example protocol on page 13.		
	For first-strand cDNA synthesis alone, we recommend the SuperScript <sup>™</sup> III First-Strand Synthesis System for RT-PCR (Catalog no. 18080-051).		

Removing Genomic DNA from RNA Samples	We RN (Ca for acc	strongly recommend that you decrease th A sample by performing a digest with DN atalog no. 18068-015), as described below. up to 1 μg of RNA; for larger amounts of ordingly.	ne genomic D Vase I, Amplif The DNase I c RNA, increase	NA content in the fication Grade digest is designed e volumes
	Co	mbine the following in a tube on ice:		
	C RI D D	omponent NA template Nase reaction buffer Nase I, Amplification Grade EPC-treated ddH20	<b>Conc.</b> — 10Χ 1 U/μl	<b>Volume</b> x μl 1 μl 1 μl to 10 μl
	1.	Incubate at room temperature for 15 min		
	2.	Add 1 µl of 25-mM EDTA solution to the 65°C for 10 min to inactivate the DNase I	reaction mixtu	are and incubate at
Reverse Transcription Protocol	The fro: 117 RN hex <b>No</b> tem Thi	e following protocol for generating first-st m the SuperScript <sup>™</sup> III Platinum <sup>®</sup> Two-Ste 34-050 and 11734-068). The RT Enzyme Mix aseOUT <sup>™</sup> . The 2X RT Reaction Mix contai camers (2.5 ng/µl), 10 mM MgCl <sub>2</sub> , and dN te: The <i>E. coli</i> RNase H digestion step is in plate from the cDNA:RNA hybrid molec is has been shown to increase PCR sensitiv	rand cDNA u p qRT-PCR K contains Supe ns oligo(dT) <sub>20</sub> TPs. ncluded to rem ule after first-s vity.	ses components it (Catalog nos. erScript <sup>™</sup> III RT and (2.5 μM), random nove the RNA strand synthesis.
	1.	Combine the following kit components in reactions, a master mix without RNA ma 2X RT Reaction Mix	n a tube on ice ny be preparec 10 μl	e. For multiple 1:
		RNA (10 pg to 1 µg) DEPC-treated water to 2	2 μ1 x μl 20 μl	
	2.	Gently mix tube contents and incubate at	t 25°C for 10 n	nin.
	3. Incubate tube at 42°C for 50 min.			
	4. Terminate the reaction at 85°C at 5 min, and then chill on ice.			
	5. Add 1 μl (2 U) of <i>E. coli</i> RNase H and incubate at 37°C for 20 min.			
	Sto the	re the reaction at -20°C until use. Proceed following page.	to the real-tin	ne qPCR protocol on
-				

Real-Time qPCR Protocol for Instruments Using PCR Tubes or Plates The following real-time qPCR protocol uses components from the

SuperScript<sup>™</sup> III Platinum<sup>®</sup> Two-Step qRT-PCR Kit (Catalog nos. 11734-050 and 11734-068). It has been optimized for use with real-time qPCR instruments that use tubes or plates. See the guidelines on pages 7–8. A protocol for the Roche LightCycler<sup>®</sup> is provided on the following page.

1. To reduce well-to-well variation, prepare a Master Mix of all the reaction ingredients except template:

<u>Component</u>	<u>Vol/1 rxn</u>	<u>Vol/50 rxns</u>
Platinum <sup>®</sup> Quantitative PCR SuperMix-UDG <sup>1</sup>	25 µl	1250 µl
ROX Reference Dye (optional)	1 µl	50 µl
Labeled LUX <sup>™</sup> Primer (10 µM)	1 µl	50 µl
Unlabeled primer (10 µM)	1 µl	50 µl
Sterile distilled water	to 45 µl	to 2250 µl

<sup>1</sup>Final concentration: 0.06 U/μl Platinum<sup>®</sup> *Taq* DNA polymerase, 20 mM Tris-HCl (pH 8.4), 50 mM KCl, 3 mM MgCl<sub>2</sub>, 200 μM dGTP, 200 μM dATP, 200 μM dCTP, 400 μM dUTP, 0.04 U/μl UDG

- 2. Program the real-time qPCR instrument as follows:
  - **Thermal Cycling** 50°C, 2 min hold (UDG treatment) 95°C, 2 min hold 45 cycles of: 95°C, 15 s 55°C, 30 s 72°C, 30 s **Melting Curve Analysis**

Refer to instrument documentation

- 3. Add 45  $\mu$ l of the Master Mix to an optical PCR tube or each well of a 96-well PCR plate.
- 4. Add ~5  $\mu$ l (10<sup>2</sup> to 10<sup>7</sup> copies or 1 pg to 10  $\mu$ g) of the cDNA from the firststrand synthesis reaction (step 5, page 13) to each reaction vessel. Cap or seal the tube/plate.
- 5. Gently mix and make sure that all components are at the bottom of the tube/plate wells. Centrifuge briefly if needed.
- 6. Place reaction in the real-time qPCR instrument and run the program. Collect and analyze results.

Protocol for the Roche LightCycler <sup>®</sup>		The following protocol uses components from the SuperScript <sup>™</sup> III Platinum <sup>®</sup> Two-Step qRT-PCR Kit (Catalog nos. 11734-050 and 11734-068) and has been optimized for the Roche LightCycler <sup>®</sup> . Consult the LightCycler <sup>®</sup> documentation for detailed instructions on preparing the capillary tubes and operating the instrument. FAM-labeled LUX <sup>™</sup> Primers are also compatible with Roche enzyme mixes.			
		Note: JOE-labeled LUX <sup>™</sup> Primers are not compatible with the current version of the LightCycler <sup>®</sup> ; use FAM-labeled primers only. The following protocol uses a 20-µl reaction volume. Before proceeding, see the real-time qPCR guidelines on pages 7–8.			
	1. To reduce well-to-well variation, prepare a Master Mix of all the reaction ingredients except template. The following table provides volumes for one reaction and 34 reactions (scale as needed):				
	Co Pla FA Un BS Pla Ste	mponent tinum <sup>®</sup> Quantitative PCR SuperMix-U M-labeled LUX <sup>™</sup> Primer (10 μM) labeled primer (10 μM) A, UltraPure (5 mg/ml) tinum <sup>®</sup> Taq DNA Polymerase <sup>2</sup> rile distilled water	JDG <sup>1</sup>	Vol/1 rxn 10 μ1 1 μ1 1 μ1 1 μ1 0.12 μ1 to 18 μ1	<u>Vol/34 rxns</u> 340 μl 34 μl 34 μl 34 μl 4 μl to 612 μl
		<sup>1</sup> Final concentration: 0.06 U/μl Platinum <sup>e</sup> (pH 8.4), 50 mM KCl, 3 mM MgCl <sub>2</sub> , 200 μN 400 μM dUTP, 0.04 U/μl UDG <sup>2</sup> Total units of Platinum <sup>®</sup> Taq DNA Polym	M dGTP, 20 erase in the	polymerase, 2 0 μM dATP, 2 e reaction is 1.:	0 mM Tris-HCl 200 μM dCTP, 2 (including 0.6 U
	2.	Set the fluorescence on the Roche Lig	ahtCycler <sup>®</sup>	<sup>®</sup> to the F1 ch	annel.
	3.	Program the instrument as follows:			
		Thermal Cycling	Melting	Curve Analy	ysis (optional)
	4	Program choice: Amplification Analysis mode: Quantification Cycling: 50°C, 2 min hold (UDG treatment) 95°C, 2 min hold 45 cycles of: 94°C, 5 s 55°C, 10 s (single acquire) 72°C, 10 s	Program of Analysis of Cycling: 95°C, 0 s 55°C, 15 s 95°C, 0 40°C, 0 s	<i>choice:</i> Meltin <i>mode:</i> Meltin sec (increase 0. continuous	ng curve g curves 1°C/s with acquisition)
	+. -	A LL 2 L $(10^2 + 10^7)$			
	5.	Add 2 $\mu$ I (10 <sup>2</sup> to 10 <sup>7</sup> copies or 1 pg to strand synthesis reaction (step 5, page	10 μg) of t ge 13) to ea	the cDNA from the cDNA from the cDNA from the cDNA from the contract of the contract of the contract of the contract of the cDNA from the contract of the cDNA from the contract of the contra	om the first- l cap the tube.
	6.	Centrifuge the tubes at $700 \times g$ for 5 s	seconds.		
	7.	Place the reaction tubes in the rotor of program. Collect and analyze results	of the Ligh 5.	ntCycler <sup>®</sup> and	l run the
-					

### One-Step Real-Time qRT-PCR

Introduction	In one-step real-time qRT-PCR, cDNA synthesis and PCR are performed in a single reaction tube using total RNA or mRNA as starting material. The optimized enzyme mixture includes both a reverse transcriptase and a DNA polymerase. This section provides guidelines and an example protocol for performing one-step real-time qRT-PCR using LUX <sup>™</sup> Primers. The example protocol uses the SuperScript <sup>™</sup> III Platinum <sup>®</sup> One-Step Quantitative RT-PCR System for superior specificity and sensitivity with LUX <sup>™</sup> Primers.
Primer Concentration	For optimal one-step qRT-PCR, primer titrations of 50–500 nM per primer are recommended. The 50- $\mu$ l sample reaction on page 19 uses 200 nM of each primer, equivalent to 1 $\mu$ l of a 10 $\mu$ M primer solution. Also see the note below.
	In one-step qRT-PCR, the reverse primer drives the reverse transcription reaction. We have found that doubling the concentration of the reverse primer from 200 nM to 400 nM can in some cases decrease the cycle threshold for detecting a given target concentration, and thus increase sensitivity. See pages 3–4 for guidance on primer design.
Template Specifications	The target template for one-step real-time qRT-PCR is RNA—usually total cellular RNA or mRNA. The amount of template typically ranges from 1 pg to 100 ng per assay. The purity and integrity of the RNA have a direct impact on results. RNase and genomic DNA contamination are the most common problems, and purification methods should be designed to avoid these. We recommend using the Micro-to-Midi Total RNA Purification System (Catalog no 12183-018) or TRIzol® reagent (Catalog no. 15596-026) to isolate total RNA. High-quality total RNA can be purified from as little as 100 cells up to 10 <sup>7</sup> cells or 200 mg of tissue. To isolate mRNA, we recommend using the FastTrack® 2.0 mRNA Isolation
Enzyme Specifications	Kit (Catalog no. K1593-02). The one-step qRT-PCR enzyme mix should contain an optimized mixture of a high-specificity, high-yield reverse transcriptase and a hot-start DNA polymerase. We recommend using the SuperScript <sup>™</sup> III Platinum <sup>®</sup> One-Step Quantitative RT-PCR System (Catalog nos. 11732-020 and -088), which uses a SuperScript <sup>™</sup> III RT/Platinum <sup>®</sup> <i>Taq</i> enzyme mix. See the sample reactions on pages 19–20.
Magnesium Concentration	The optimal MgCl <sub>2</sub> concentration for a given target/primer/polymerase combination can vary between 1 mM and 10 mM, but is usually in the range of 3 mM (see the sample reaction on page 19).
dNTP Concentration	The optimal concentration of dATP, dCTP, dGTP, and dTTP is 200 $\mu$ M each. If dUTP is used in place of dTTP, its optimal concentration is 400 $\mu$ M.

ROX Reference Dye	We recommend using ROX Reference Dye (Cat. no. 12223-023) to normalize the fluorescent reporter signal for instruments that are compatible with this option. ROX Reference Dye can be used to adjust for non-PCR-related fluctuations in fluorescence between reactions, and provides a stable baseline in multiplex reactions. It is composed of a glycine conjugate of 5-carboxy-X- rhodamine, succinimidyl ester (25 $\mu$ M) in 20 mM Tris-HCl (pH 8.4), 0.1 mM EDTA, and 0.01% Tween <sup>®</sup> 20.
-	reaction volume.
Bovine Serum Albumin (BSA)	Bovine serum albumin (BSA) is required in qPCR reactions on the Roche LightCycler <sup>®</sup> because the glass capillaries in the LightCycler <sup>®</sup> have a high surface-to-volume ratio and the glass surface binds molecules like <i>Taq</i> DNA polymerase, effectively removing them from the reaction. The addition of BSA blocks this surface binding.
	Nonacetylated BSA is strongly recommended because acetylated BSA will inhibit PCR at the concentrations required in LightCycler <sup>®</sup> reactions. This inhibition is most likely due to the transfer of acetyl groups to essential components of the PCR, like the <i>Taq</i> DNA polymerase.
_	To ensure that the BSA does not contain RNA or DNA, we recommend using ultrapure, molecular biology-grade nonacetylated BSA from Panvera/Invitrogen (Cat. nos. P2489 and P2046).
Melting Curve Analysis	Melting curve analysis is strongly recommended during one-step qRT-PCR to identify the presence of primer dimers and analyze the specificity of the reaction. Program your real-time instrument to perform this analysis after thermocycling.
	Melting curve analysis identifies the change in fluorescent signal that occurs as double-stranded DNA (dsDNA) dissociates or "melts" into single-stranded DNA. By identifying the temperature at which the dsDNA dissociates, you can distinguish smaller artifacts like primer dimers with a lower annealing temperature from larger amplicons with a higher annealing temperature.
	The presence of primer dimers in samples containing template decreases reaction efficiency and obscures analysis and determination of cycle thresholds. The formation of primer dimers most often occurs in no-template controls, where the polymerase enzyme is essentially idle, and in this case the quantitative analysis of the template samples is not affected. Melting curve analysis of no-template controls can discriminate between primer dimers and spurious amplification due to contaminating nucleic acids in reagent components.

Instrument Settings	<ul> <li>Follow the manufacturer's instructions for construment for use with LUX<sup>™</sup> Primers. Note</li> <li>Program your instrument to perform methermocycling, if this option is available.</li> <li>The quencher setting on the instrument a Primers do not contain a quencher.</li> <li>We recommend using ROX Reference Dy compatible with this option. Adjust the instrument and settings for specific www.invitrogen.com/lux; click on Instrument.</li> </ul>	onfiguring your the following g elting curve analy should reflect the ye, if your instru instrument settir ic instruments a <b>ent Protocols</b> .	real-time qPCR eneral settings: ysis at the end of e fact that LUX <sup>™</sup> ument is ngs accordingly. re available at
Removing Genomic DNA from RNA Samples	We recommend that you decrease the genom sample by performing a digest with DNase I no. 18068-015), as described below. The DNa 1 µg of RNA; for larger amounts of RNA, inc Combine the following in a tube on ice:	nic DNA content , Amplification ( ise I digest is des crease volumes a	t in the RNA Grade (Catalog signed for up to accordingly.
	<ul> <li>Component RNA template DNase reaction buffer DNase I, Amplification Grade DEPC-treated ddH<sub>2</sub>0</li> <li>1. Incubate at room temperature for 15 min</li> <li>2. Add 1 µl of 25-mM EDTA solution to the 65°C for 10 min to inactivate the DNase</li> <li>To verify the absence of genomic DNA in the reaction identical to the reactions on pages 1° DNA polymerase (Catalog no. 10966-018) in RT/Platinum<sup>®</sup> Taq Mix.</li> </ul>	Conc. — 10X 1 U/μl n. e reaction mixtur I. e RNA sample, p 9–20, using 2 U o place of the Sup	Volume $x \mu l$ $1 \mu l$ $1 \mu l$ to $10 \mu l$ be and incubate at prepare a control of Platinum <sup>®</sup> Taq prescript <sup>™</sup> III

Protocol for Instruments Using PCR Tubes or Plates	The following protocol using the SuperScript <sup>™</sup> III Platinum <sup>®</sup> One-Step Quantitative RT-PCR System has been optimized for LUX <sup>™</sup> Primers. Further optimization may be required.				
	<b>Note:</b> Keep all components, reaction mixes and samples on ice. After assembly, transfer the reaction to a thermal cycler preheated to the cDNA synthesis temperature and immediately begin RT-PCR. We recommend performing the cDNA synthesis reaction at 50°C, but higher temperatures (up to 60°C) may be required for high GC content templates.				
	RNase inhibitor proteins, such as RNaseOUT™ (Catalog no. 10777-019), may be added to the reaction to safeguard against degradation of RNA.				
	1.	The following table provides Master Mix volume reaction size. Note that preparation of a master m quantitative applications to reduce pipetting error	es for a stand nix is <b>crucial</b> ors.	ard 50-µl in	
		Component         SuperScript™ III RT/Platinum® Taq Mix         2X Reaction Mix <sup>1</sup> ROX Reference Dye (optional)         Labeled LUX™ Primer (10 µM)         Unlabeled primer (10 µM) <sup>2</sup> RNaseOUT™ (optional)         Sterile distilled water <sup>1</sup> Supplied at 2X concentration: includes 0.4 mM of each dNTP <sup>2</sup> See the Important note on primer concentration on page 16.	<u>Vol/1 rxn</u> 1 μl 25 μl 1 μl 1 μl 1 μl 1 μl to 40 μl 2 and 6 mM Mg	<u>Vol/100 rxns</u> 100 μl 2500 μl 100 μl 100 μl 100 μl 100 μl 504	
	2.	Program the instrument with the following therm cDNA synthesis, use a 15-min incubation at 50°C	nal cycling p as a starting	rotocol (for 5 point):	
		cDNA synthesis: $50^{\circ}$ C for 15 min hold PCR: $95^{\circ}$ C for 2 min hold 40-50 cycles of: $95^{\circ}$ C, 15 s $60^{\circ}$ C, 30 s			
		<b>Melting Curve Analysis (optional)</b> Program according to instrument instructions			
	3.	For each reaction, add 40 $\mu$ l of the master mix to a tube or each well of a 96-well PCR plate on ice.	a 0.2-ml mic	rocentrifuge	
	4.	Add 10 $\mu$ l of sample RNA (1 pg to 1 $\mu$ g total RNA and cap or seal.	A) to each tu	pe/plate well,	
	5.	Gently mix and make sure that all components ar tube/plate wells. Centrifuge briefly if needed.	re at the bott	om of the	
	6.	Place reactions in a preheated thermal cycler prog above. Collect data and analyze results.	grammed as	described	
-			Contin	ued on next page	

Protocol for the Roche LightCycler <sup>®</sup>	Th Qu Ro LU No Lig Af cD per	e following protocol using the SuperS antitative RT-PCR System has been of che LightCycler <sup>®</sup> . Further optimizatio IX <sup>™</sup> Primers are also compatible with I ote: JOE-labeled primers are not compa- ghtCycler <sup>®</sup> ; use FAM-labeled primers of ter assembly, transfer the reaction to a NA synthesis temperature and immed rforming the cDNA synthesis reaction	cript <sup>™</sup> III P ptimized fo n may be r Roche enzy atible with only. thermal cy diately beg at 50°C, bu	'latinum <sup>®</sup> One- or LUX <sup>™</sup> Prime equired. FAM- me mixes. the current ve ycler preheated in RT-PCR. We ut higher temp	Step ers and the -labeled rsion of the d to the e recommend eratures (up
	to 60°C) may be required for high GC content templates. RNase inhibitor proteins, such as RNaseOUT <sup>™</sup> (Catalog no. 10777-019), may be added to the reaction to safeguard against degradation of RNA.				
	1.	The following table provides Master reaction size. Note that preparation quantitative applications to reduce p	Mix volun of a master vipetting er	nes for a stand mix is <b>crucial</b> rors.	ard 20-µl in
		Component         SuperScript <sup>™</sup> III RT/Platinum <sup>®</sup> Taq N         2X Reaction Mix <sup>1</sup> FAM-labeled LUX <sup>™</sup> Primer (10 µM) <sup>2</sup> Unlabeled primer (10 µM) <sup>3</sup> Bovine serum albumin (5 mg/ml) <sup>4</sup> Sterile distilled water <sup>1</sup> Includes 0.4 mM of each dNTP and 6 m <sup>2</sup> In the LightCycler <sup>®</sup> reaction, the LUX <sup>™</sup> N <sup>3</sup> See the Important note on primer conce <sup>4</sup> Validated with non-acetylated Ultrapur         nos. P2489 and P2046)	Mix M MgSO <sub>4</sub> Fluorogenic ntration on e BSA (10%	<u>Vol/1 rxn</u> 0.8 μl 10 μl 1 μl 1 μl 1 μl to 18 μl Primer must be page 16. solution) from F	<u>Vol/34 rxns</u> 27.2 μl 340 μl 34 μl 34 μl 34 μl to 612 μl FAM labeled. Panvera (Cat.
	2.	Set the fluorescence on the Roche Lig	ghtCycler®	to the F1 chan	nel.
	3.	Program the instrument as follows:			
		Thermal Cycling Program choice: Amplification Analysis mode: Quantification Cycling: 45°C, 30 min hold (cDNA synthesis) 95°C, 2 min hold 50 cycles of: 95°C, 5 s 55°C, 10 s (single acquire) 72°C, 10 s	Melting C Program c Analysis m Cycling: $95^{\circ}$ C, 0 s $55^{\circ}$ C, 15 s $95^{\circ}$ C, 0 $40^{\circ}$ C, 0 s	Curve Analysi hoice: Melting c hode: Melting c ec (increase 0.1% continuous ac	<b>s (optional)</b> curve urves C/s with quisition)
	4.	Add 18 $\mu l$ of Master Mix to each cap	illary tube	of the LightCy	cler <sup>®</sup> on ice.
	5.	Add 2 $\mu l$ of sample RNA (1 pg to 1 $\mu$ and cap the tube.	g total RN.	A) to each capi	illary tube
	6.	Centrifuge the tubes at $700 \times g$ for 5 s	seconds.		
	7.	Place the reaction tubes in the rotor of program. Collect and analyze results	of the Light s.	tCycler <sup>®</sup> and r	un the

### Troubleshooting

Problem	Cause	Solution
Signals are present in no- template controls, and/or multiple peaks are present in the melting curve graph	Template or reagents are contaminated by nucleic acids (DNA, cDNA)	Use melting curve analysis and/or run the PCR products on a 4% agarose gel in an area separate from the reaction assembly area to identify contaminants.
		To reduce the risk of contamination, take standard precautions when preparing your PCR reactions. Ideally, amplification reactions should be assembled in a DNA-free environment. We recommend using aerosol-resistant barrier tips.
	Amplification of PCR carryover products	Analyze the PCR product on a 4% agarose gel in an area separate from the reaction assembly area to identify contaminants.
		We recommend using a UDG-based carryover prevention system such as Platinum <sup>®</sup> Quantitative PCR SuperMix-UDG or the SuperScript <sup>™</sup> III Platinum <sup>®</sup> Two-Step qRT-PCR Kit (protocols provided on pages 9–10 and pages 12–15, respectively). Since dUTP is substituted for dTTP in the reaction cocktail, any amplified DNA will contain uracil. UDG prevents reamplification of PCR carryover products by removing uracil residues from single or double stranded DNA. dU-containing DNA that has been digested with UDG is unable to serve as template in future PCRs. UDG is inactivated at high temperature during PCR thermal cycling, thereby allowing amplification of genuine target sequence(s).
	Primer dimers or other primer artifacts are present	Use melting curve analysis of the PCR product to identify primer dimers by their lower melting point temperature. Confirm that your primer designs have low scores (0.0-4.0) when generated by the LUX <sup>™</sup> Designer to minimize self-annealing. Redesign primers if necessary.
		If you are redesigning primers, you can first try redesigning only the unlabeled primer to save the cost of the LUX <sup>™</sup> primer.
		Primer contamination or truncated or degraded primers can also lead to artifacts. Check the purity of your primers by gel electrophoresis. If agarose gels are used, we recommend cooling the gels before visualization with intercalating dyes.

## Troubleshooting, Continued

Problem	Cause	Solution
No amplification curve appears on the qPCR graph	There is no PCR product	Run the PCR product on a gel to determine whether PCR worked. Then proceed to the troubleshooting steps below.
No PCR product is evident, either in the qPCR graph or on a gel	The protocol was not followed correctly	Verify that all steps have been followed and the correct reagents, dilutions, volumes, and cycling parameters have been used.
	Template contains inhibitors, nucleases, or proteases, or has otherwise been degraded.	Purify or re-purify your template.
	Primer designs are not optimal	Confirm that you are using the correct primers for your sequence, the primer design scores are within the 0.0-4.0 range in the LUX <sup>™</sup> Designer, and the optimal melting temperatures have been specified. Redesign primers if necessary. When redesigning primers, note that you can first try redesigning only the unlabeled primer to save the cost of the LUX <sup>™</sup> primer.
PCR product is evident in the gel, but not on the qPCR graph	qPCR instrument settings are incorrect	Confirm that you are using the correct instrument settings (dye selection, reference dye, filters, acquisition points, etc.).
	Problems with your specific qPCR instrument	For instrument-specific tips and troubleshooting using LUX <sup>™</sup> Primers, see the instrument protocols at <u>www.invitrogen.com/lux</u> .
PCR efficiency is above 110%	Template contains inhibitors, nucleases, or proteases, or has otherwise been degraded.	Purify or re-purify your template. Inhibitors in the template may result in changes in PCR efficiency between dilutions
PCR efficiency is below 90%	The PCR conditions are suboptimal	Verify that the amount of primers you are using is correct and that the labeled primer has not been exposed to direct light. Verify that the reagents you are using have not been fragge theused multiple times and have not
		sat at room temperature for too long.
	Reagent concentration in a multiplex reaction may be limiting the rate of the reaction.	Perform a single-plex reaction using the same primers and template to check efficiency. Then determine which one of your primer sets should be in limiting concentration. See the multiplex guidelines on page 11.

#### **Accessory Products**

#### Products

The following products are available for use with  $LUX^{\mbox{\tiny M}}$  Primers in real-time qPCR and qRT-PCR protocols:

Product	Amount	Catalog no.
Platinum <sup>®</sup> Quantitative PCR SuperMix-UDG	100 rxns	11730-017
	500 rxns	11730-025
SuperScript <sup>™</sup> III Platinum <sup>®</sup> Two-Step qRT-PCR Kit	100 PCRs	11734-050
	500 PCRs	11734-068
SuperScript <sup>™</sup> III Platinum <sup>®</sup> One-Step Quantitative RT-	100 rxns	11732-020
PCR System	500 rxns	11732-088
Platinum <sup>®</sup> Taq DNA Polymerase	100 rxns	10966-018
	250 rxns	10966-026
	500 rxns	10966-034
	5,000 rxns	10966-083
Micro-to-Midi Total RNA Purification System	50 rxns	12183-018
TRIzol <sup>®</sup> Reagent	100 ml	15596-026
	200 ml	15596-018
Micro-FastTrack <sup>™</sup> 2.0 mRNA Isolation Kit	20 rxns	K1520-02
ROX Reference Dye	500 µl	12223-023
DNase I, Amplification Grade (1 U/µl)	100 U	18068-015
RNaseOUT <sup>™</sup> Recombinant Ribonuclease Inhibitor	5,000 U	10777-019
(40 U/µl)		
10 mM dNTP Mix	100 µl	18427-013
DEPC-treated water	4 x 1.25 ml	10813-012

#### **Purchaser Notification**

Limited Use Label License No. 114: LUX<sup>™</sup> Fluorogenic Primer The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for commercial purposes. The buyer may transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for the commercial purposes of the buyer, and that such collaborator agrees in writing (a) to not transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for commercial purposes. Commercial purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. Invitrogen Corporation will not assert a claim against the buyer of infringement of patents owned by Invitrogen based upon the manufacture, use or sale of a therapeutic, clinical diagnostic, vaccine or prophylactic product developed in research by the buyer in which this product or its components was employed, provided that neither this product nor any of its components was used in the manufacture of such product. If the purchaser is not willing to accept the limitations of this limited use statement, Invitrogen is willing to accept return of the products with a full refund. For information on purchasing a license to this product for purposes other than research, contact Licensing Department, 1600 Faraday Avenue, Carlsbad, California 92008. Phone (760) 603-7200. Fax (760) 602-6500.

Limited Use Label License No. 4: Products for PCR which do not include any rights to perform PCR This product is optimized for use in the Polymerase Chain Reaction (PCR) covered by patents owned by Roche Molecular Systems, Inc. and F. Hoffmann-La Roche, Ltd. ("Roche"). No license under these patents to use the PCR process is conveyed expressly or by implication to the purchaser by the purchase of this product. A license to use the PCR process for certain research and development activities accompanies the purchase of certain reagents from licensed suppliers such as Invitrogen, when used in conjunction with an Authorized Thermal Cycler, or is available from Applied Biosystems. 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 at Roche Molecular Systems, Inc., 1145 Atlantic Avenue, Alameda, California 94501.

#### **Technical Service**

#### World Wide Web

Visit the Invitrogen Web Resource using your World Wide Web browser. At the site, you can:

- Get the scoop on our hot new products and special product offers
- View and download vector maps and sequences
- Download manuals in Adobe® Acrobat® (PDF) format
- Explore our catalog with full color graphics
- Obtain citations for Invitrogen products
- Request catalog and product literature

Once connected to the Internet, launch your Web browser (Internet Explorer 5.0 or newer or Netscape 4.0 or newer), then enter the following location (or URL):

#### http://www.invitrogen.com

...and the program will connect directly. Click on underlined text or outlined graphics to explore. Don't forget to put a bookmark at our site for easy reference!

Contact Us	For more information or technical assistance, please call, write, fax, or email.
	Additional international offices are listed on our Web page
	(www.invitrogen.com).

Corporate Headquarters:	Japanese Headquarters:	European Headquarters:
Invitrogen Corporation	Invitrogen Japan K.K.	Invitrogen Ltd
1600 Faraday Avenue	Nihonbashi Hama-Cho Park	3 Fountain Drive
Carlsbad, CA 92008	Bldg. 4F	Inchinnan Business Park
USA	2-35-4, Hama-Cho, Nihonbashi	Paisley PA4 9RF, UK
Tel: 1 760 603 7200	Tel: 81 3 3663 7972	Tel: +44 (0) 141 814 6100
Tel (Toll Free): 1 800 955 6288	Fax: 81 3 3663 8242	Tel (Toll Free in UK):
Fax: 1 760 602 6500	E-mail: j <u>pinfo@invitrogen.com</u>	0800 5345 5345
E-mail:		Fax: +44 (0) 141 814 6287
tech_service@invitrogen.com		E-mail: eurotech@invitrogen.com

MSDS Requests	To the	o request an MSDS, please visit our Web site ( <u>www.invitrogen.com</u> ) and follow the instructions below.	
	1.	On the home page, go to the left-hand column under 'Technical Resources' and select 'MSDS Requests'.	
	2.	Follow instructions on the page and fill out all the required fields.	
	3.	To request additional MSDSs, click the 'Add Another' button.	
	4.	All requests will be faxed unless another method is selected.	
	5.	When you are finished entering information, click the 'Submit' button. Your MSDS will be sent within 24 hours.	



#### Technical Service, Continued

#### **Limited Warranty**

Invitrogen is committed to providing our customers with high-quality goods and services. Our goal is to ensure that every customer is 100% satisfied with our products and our service. If you should have any questions or concerns about an Invitrogen product or service, please contact our Technical Service Representatives.

Invitrogen warrants that all of its products will perform according to the specifications stated on the certificate of analysis. The company will replace, free of charge, any product that does not meet those specifications. <u>This warranty</u> <u>limits Invitrogen Corporation's liability only to the cost of the product</u>. No warranty is granted for products beyond their listed expiration date. No warranty is applicable unless all product components are stored in accordance with instructions. Invitrogen reserves the right to select the method(s) used to analyze a product unless Invitrogen agrees to a specified method in writing prior to acceptance of the order.

Invitrogen makes every effort to ensure the accuracy of its publications, but realizes that the occasional typographical or other error is inevitable. Therefore Invitrogen makes no warranty of any kind regarding the contents of any publications or documentation. If you discover an error in any of our publications, please report it to our Technical Service Representatives.

Invitrogen assumes no responsibility or liability for any special, incidental, indirect or consequential loss or damage whatsoever. The above limited warranty is sole and exclusive. No other warranty is made, whether expressed or implied, including any warranty of merchantability or fitness for a particular purpose.

- Ailenberg, M., and Silverman, M. (2000) Controlled hot start and improved specificity in carrying out PCR utilizing touch-up and loop incorporated primers (TULIPS). *BioTechniques 29*, 1018–1024.
- Bustin, S. A. (2000) Absolute quantification of mRNA using real-time reverse transcription polymerase chain reaction assays. *J. Mol. Endocrinol.* 25, 169–193.
- Cardullo, R. A., Agrawal, S., Flores, C., Zamecnik, P. C., and Wolf, D. E. (1988) Detection of nucleic acid hybridization by nonradiative fluorescence resonance energy transfer. *Proc. Natl. Acad. Sci. USA 85*, 8790–8794.
- Crockett, A.O., and Wittwer, C.T. (2001) Fluorescein-labeled oligonucleotides for real-time pcr: using the inherent quenching of deoxyguanosine nucleotides. *Anal. Biochem.* 290, 89–97.
- Higuchi, R., Fockler, C., Walsh, P.S., and Griffith, R. (1992) Simultaneous amplification and detection of specific DNA sequences. *Biotechnology* 10, 413–417.
- Higuchi, R., Fockler, C., Dollinger, G., and Watson, R. (1993) Kinetic PCR analysis: real-time monitoring of DNA amplification reactions. *Biotechnology* 11, 1026–1030.
- Holland et al. (1991) Detection of specific polymerase chain reaction product by utilizing the 5'-3' exonuclease activity of Thermus aquaticus DNA polymerase. *Proc. Natl. Acad. Sci. USA 88*, 7276–7280.
- Kaboev, O. K., Luchkina, L. A., Tret'iakov, A. N., and Bahrmand, A.R. (2000) PCR hot start using primers with the structure of molecular beacons (hairpin-like structure). *Nucleic Acids Res.* 28, e94.
- Knemeyer, J.P., Marme, N., and Sauer, M. (2000) Probes for detection of specific DNA sequences at the single-molecule level. *Anal. Chem.* 72, 3717–3724.
- Murchie, A. I. H., Clegg, R. M., von Kitzing, E., Duckkett, D. R., Diekmann, S., and Lilley D. M. J. (1989) Fluorescence energy transfer shows that the four-way DNA junction is a right-handed cross of antiparallel molecules. *Nature* 341, 763–766.
- Myakishev, M.V., Khripin, Y., Hu, S., and Hamer, D. H. (2001) High-throughput SNP genotyping by allele-specific PCR with universal energy-transfer-labeled primers. *Genome Res.* 11, 163–169.
- Nazarenko, I., Lowe, B., Darfler, M., Ikonomi, P., Schuster, D., and Rashtchian, A. (2002) Multiplex quantitative PCR using self-quenched primers labeled with a single fluorophore. *Nucl. Acids Res.* 30, e37
- Nazarenko, I., Pires, R., Lowe, B., Obaidy, M., and Rashtchian, A. (2002) Effect of primary and secondary structure of oligodeoxyribonucleotides on the fluorescent properties of conjugated dyes. *Nucl. Acids Res.* 30, 2089–2095
- Nazarenko, I.A., Bhatnagar, S.K., and Hohman, R.J. (1997) A closed tube format for amplification and detection of DNA based on energy transfer. *Nucleic Acids Res.* 25, 2516–2521.
- Nuovo, G. J., Hohman, R. J., Nardone, G. A., and Nazarenko I. (1999) In situ amplification using universal energy transfer-labeled primers. *J. Histochem. Cytochem.* 47, 273–279.
- Todd, A. V., Fuery, C. J., Impey, H. L., Applegate, T. L. and Haughton, M.A. (2000) DzyNA-PCR: use of DNAzymes to detect and quantify nucleic acid sequences in a real-time fluorescent format. *Clin. Chem.* 46, 625–630.
- Tyagi, S., and Kramer, F.R. (1996) Molecular beacons: probes that fluoresce upon hybridization. *Nature Biotechnol.* 14, 303–308.
- Wittwer, C.T., Herrmann, M.G., Moss, A.A., Rasmussen, R.P. (1997) Continuous fluorescence monitoring of rapid cycle DNA amplification. *BioTechniques* 22, 130–138.



#### **United States Headquarters:**

Invitrogen Corporation 1600 Faraday Avenue Carlsbad, California 92008 Tel: 1 760 603 7200 Tel (Toll Free): 1 800 955 6288 Fax: 1 760 603 7229 Email: tech\_service@invitrogen.com

#### **European Headquarters:**

Invitrogen Ltd 3 Fountain Drive Inchinnan Business Park Paisley PA4 9RF, UK Tel (Free Phone Orders): 0800 269 210 Tel (General Enquiries): 0800 5345 5345 Fax: +44 (0) 141 814 6287 Email: eurotech@invitrogen.com

#### **International Offices:**

Argentina 5411 4556 0844 Australia 1 800 331 627 Austria 0800 20 1087 Belgium 0800 14894 Brazil 0800 11 0575 Canada 800 263 6236 China 10 6849 2578 Denmark 80 30 17 40

France 0800 23 20 79 Germany 0800 083 0902 Hong Kong 2407 8450 India 11 577 3282 Italy 02 98 22 201 Japan 03 3663 7974 The Netherlands 0800 099 3310 New Zealand 0800 600 200 Norway 00800 5456 5456

Spain & Portugal 900 181 461 Sweden 020 26 34 52 Switzerland 0800 848 800 Taiwan 2 2651 6156 UK 0800 838 380 For other countries see our website

www.invitrogen.com