



## SCRIPT RT-qPCR ProbesMaster Lyophilisate

lyophilised RT-real-time-PCR Master Mix for dual labeled probes

Cat. No.	Amount
PCR-159S	192 reactions x 20 µl
PCR-159L	960 reactions x 20 µl

### For general laboratory use.

**Shipping:** shipped at ambient temperature

**Storage Conditions:** store at ambient temperature

**Additional Storage Conditions:** Store in an aluminium-coated bag or on a dry place.

Lyophilisates may hydrate at humidity levels >70 % when sealing is opened.

**Shelf Life:** 6 months in sealed package

### Description:

SCRIPT RT-qPCR ProbesMaster Lyophilisate is designed for performing maximum sensitive and specific RT-qPCRs convenient in single tubes. The enzyme mix is based on a genetically engineered RT polymerase with enhanced thermal stability resulting in an increased specificity, higher cDNA yield and an improved efficiency for highly structured and long cDNA fragments.

The lyophilisate contains all reagents required for RTPCR (except template and primer) in one tube to ensure fast and easy preparation with a minimum of pipetting steps. The premium quality enzyme mix, ultrapure dNTPs and the optimized complete reaction buffer ensure superior amplification results.

SCRIPT RT-qPCR ProbesMaster is used to amplify double-stranded DNA from single-stranded RNA templates. In the RT step the reverse transcriptase synthesizes single-stranded DNA molecules (cDNA) complementary to the RNA template. In the first cycle of the PCR step Taq DNA polymerase synthesizes DNA molecules complementary to the cDNA, thus generating a double-stranded DNA template. During subsequent rounds of cycling the DNA polymerase exponentially amplifies this double-stranded DNA template.

In one-step RT-qPCR all components of RT and PCR are mixed in one tube prior to starting the reaction and thus carried out sequentially without opening the tube.

This offers tremendous convenience when applied to analysis of single targets from multiple samples of RNA and minimizes the risk of contaminations.

### Content:

SCRIPT RT-qPCR ProbesMaster Lyophilisate

Preloaded lyophilisates containing SCRIPT Reverse Transcriptase, Hot Start Polymerase Ab+, dNTPs, Reaction Buffer, MgCl<sub>2</sub> and stabilizers.

### PCR grade water

### Handling

SCRIPT RT-qPCR ProbesMaster Lyophilisate is delivered in PCR reaction tube strips or 96-well plates preloaded with a complete qPCR master mix in a dry, room temperature stable format. The lyophilisate combines highest performance with convenience of use and stability. There is no need for freezing, thawing or pipetting on ice. The few remaining pipetting steps minimize the risk of errors or contaminations.

Each vial contains all components (except primers and template) required for a 20 µl RT-qPCR assay.

To perform the assay, only fill up the vials with a mix of primers and RNA template.

### Sensitivity

Targets can generally be detected from 10 pg to 500 ng polyA RNA or 10 pg to 1 µg total RNA. Even lower amounts of RNA may be successfully amplified by using highly expressed transcripts.



## SCRIPT RT-qPCR ProbesMaster Lyophilisate

lyophilised RT-real-time-PCR Master Mix for dual labeled probes

### RT-qPCR assay preparation

#### 1. Preparation of the RNA/Primer Mix

Add the following components to a nuclease-free microtube and mix by pipetting gently up and down:

Component	stock conc.	final conc.	1 assay
RNA tem-plate		10 pg - 1 µg	
forward Primer	10 µM	400 - 600 nM	0.8 - 1.2 µl
reverse Primer	10 µM	400 - 600 nM	0.8 - 1.2 µl
RNase-free water			fill up to 20 µl

#### 2. Denaturation and primer annealing (optional)

*Please note: Sample denaturation is particularly recommended for RNA targets that exhibit a high degree of secondary structure, for self- or cross-complementary primers and for initial experiments with new targets. For many standard combinations of RNA and primers heat treatment may be omitted with no negative effect on results.*

Incubate the mixture at 70°C for 5 min and place it at room temperature for 5 min.

#### 3. Dispensing the master mix

Dispense 20 µl of the RNA/Primer Mix to each lyophilisate containing tube or well of the plate.

### Reverse transcription and thermal cycling:

Place the vials in a PCR cycler and start the following program.

reverse transcription <sup>1)</sup>	50-55 °C	10-15 min	1x
initial denaturation <sup>2)</sup>	95°C	5 min	1x
denaturation	95°C	15 sec	35-45x
annealing and elongation	60-65 °C <sup>3)</sup>	1 min <sup>4)</sup>	35-45x

<sup>1)</sup> A reverse transcription time of 10 min is recommended for optimal amplicon lengths between 100 and 200 bp. Longer amplicons up to 500 bp may require a prolonged incubation of 15 min. Add 3 min for each additional 100 bp. The optimal temperature depends on the structural features of the RNA. Increase the temperature to 55°C for difficult templates with high secondary structure. Note that optimal reaction time and temperature should be adjusted for each particular RNA.

<sup>2)</sup> An initial denaturation time of 5 min is recommended to inactivate the reverse transcriptase

<sup>3)</sup> The annealing temperature depends on the melting temperature of

the primers and DNA probe used.

<sup>4)</sup> The elongation time depends on the length of the amplicon. A time of 1 min for a fragment of 1,000 bp is recommended.

For optimal specificity and amplification an individual optimization of the recommended parameters may be necessary. Note that optimal reaction times and temperatures should be adjusted for each particular RNA / primer pair.