



# SCRIPT RT-qPCR SybrMaster

RT-real-time-PCR mix with SYBR® Green fluorescent DNA stain

Cat. No.	Amount
PCR-520S	2 x 1,25 ml (2x conc.)
PCR-520L	10 x 1,25 ml (2x conc.)

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Additional Storage Conditions: avoid freeze/thaw cycles, store dark

stable at 4 °C for up to 4 weeks

Shelf Life: 12 months

Form: liquid

Concentration: 2x conc.

Spectroscopic Properties:  $\lambda_{exc}$  494 nm (bound to DNA),  $\lambda_{em}$  521 nm (bound to DNA)

### **Description:**

SCRIPT RT-qPCR SybrMaster is designed for quantitative real-time analyses of RNA templates using the fluorescent DNA stain SYBR® Green. The ready-to-use mix is based on a genetically engineered reverse transcriptase with enhanced thermal stability providing increased specificity, high cDNA yield and improved efficiency for highly structured and long cDNA fragments.

The 2x conc. mix contains all reagents required for RT-qPCR (except template and primers) allow fast and easy preparation with a minimum of pipetting steps. The premium quality enzymes and the optimized reaction buffer ensure superior real time PCR results.

RT-qPCR is used to amplify double-stranded DNA from singlestranded RNA templates to allow a rapid real-time quantification of RNA targets. In the reverse transcription step the reverse transcriptase synthesizes single-stranded DNA molecules (cDNA) complementary to the RNA template. In the first cycle of the PCR step the hot-start DNA polymerase synthesizes DNA molecules complementary to the cDNA, thus generating a double-stranded DNA template. The hot-start polymerase activity is blocked at ambient temperature and switched on automatically at the onset of the initial denaturation. The thermal activation prevents the extension of non-specifically annealed primers and primer-dimer formations at low temperatures during PCR setup.

One-step RT-qPCR offers tremendous convenience when applied to analysis of targets from multiple samples of RNA and minimizes the risk of contaminations.

# Content:

### SCRIPT RT-qPCR SybrMaster

Ready-to-use mix of SCRIPT Reverse Transcriptase, Hot Start Polymerase, RNase Inhibitor, dNTPs, reaction buffer, SybrGreen DNA intercalotor dye and stabilizers.

### **PCR-grade Water**

## SYBR® Green fluorescent DNA stain:

SYBR® Green fluorescent DNA stain is a superior DNA intercalator dye specially developed for DNA analysis applications like real-time PCR (qPCR). Upon binding to DNA, the non-fluorescent dye becomes highly fluorescent while showing only lowest inhibition to the PCR process. The dye is stable both thermally and hydrolytically, providing convenience during routine handling.

#### Sensitivity:

Targets can generally be detected from <1 pg to 20 ng poly(A) RNA (mRNA) or 10 pg to 1  $\mu$ g total RNA. Even lower amounts of RNA may be successfully amplified by using highly expressed transcripts.

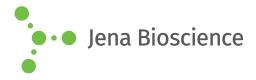
#### Preparation of the RT-qPCR assay

Add the following components to a nuclease-free microtube and mix the components by pipetting gently up and down. In general, water, RNA and primers should be mixed together before adding the master mix.



Jena Bioscience GmbH Löbstedter Str. 71 | 07749 Jena, Germany | Tel.:+49-3641-6285 000 https://www.jenabioscience.com







# SCRIPT RT-qPCR SybrMaster

RT-real-time-PCR mix with SYBR® Green fluorescent DNA stain

com- ponent	stock conc.	final conc.	20 µl assay	50 µl assay
PCR- grade water	-	-	fill up to 20 µl	fill up to 50 µl
RNA template <sup>1)</sup>	-	<100 ng	xμl	xμl
forward Primer <sup>2)</sup>	10 µM	400 nM	0.8 µl	2 µl
reverse Primer <sup>2)</sup>	10 µM	400 nM	0.8 µl	2 µl
SCRIPT RT-qPCR SybrMaster	2x 4)	1x	10 µl	25 µl

<sup>1)</sup> up to 100 ng polyA RNA or total RNA

<sup>2)</sup>The optimal concentration for primers and probe may vary from100 to 500 nm.

<sup>4)</sup> The Mix already contains RNase inhibitor that may be essential when working with low amounts of starting RNA.

Continue with reverse transcription and thermal cycling as recommended.

# Reverse transcription and thermal cycling:

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

reverse transcription <sup>5)</sup>	50-55 °C	10-15 min	1x
initial denaturation <sup>6)</sup>	95°C	5 min	1x
denaturation	95°C	15 sec	35-45x
annealing <sup>7)</sup>	55-65°C	20 sec	35-45x
elongation <sup>8)</sup>	72°C	30 sec	35-45x

<sup>5)</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>6)</sup> An initial denaturation time of 5 min is recommended to inactivate the reverse transcriptase

<sup>7)</sup> The annealing temperature depends on the melting temperature of the primers.

<sup>8)</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.

SYBR® is a registered trademark of Invitrogen Corporation, Carlsbad, California, USA

