



qPCR ProbesMaster lowROX

Master mix for real-time PCR with low ROX
2 x conc. master mix

Cat. No.	Amount
PCR-361S	2 x 1,25 ml (250 reactions x 20 µl)
PCR-361L	10 x 1,25 ml (1250 reactions x 20 µl)

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Additional Storage Conditions: avoid freeze/thaw cycles, store dark
Storage at 4 °C for up to 3 months possible.

Shelf Life: 12 months

Form: liquid

Concentration: 2x conc.

Description:

qPCR ProbesMaster lowROX is designed for quantitative real-time analysis of DNA samples using DNA probe based detection. The master mix is recommended for use with Dual Labeled Fluorescent Probes, e.g. TaqMan®, Molecular Beacons or FRET probes. It provides an easy-to-handle and powerful tool for quantification of sample DNA in a broad dynamic range of up to 6 orders of magnitude with exceptional sensitivity and precision.

The mix contains all reagents required for qPCR (except template, primer and labeled fluorescent probe) in a premixed 2x concentrated ready-to-use solution. The high specificity and sensitivity of the mix based on an optimized hot-start polymerase. Its activity is blocked by antibody at ambient temperature and switched on automatically at the onset of the initial denaturation. The thermal activation prevents the extension of nonspecifically annealed primers and primer-dimer formation at low temperatures during PCR setup.

The reaction chemistry of the kit is optimized for instruments that are compatible with the evaluation of a low ROX reference signal.

ROX reference dye:

The qPCR ProbesMaster with lowROX contains 50 nM ROX passive reference dye in the final assay. The dye does not take part in the PCR reaction but allows to normalize for non-PCR related signal variation and provides a baseline in multiplex reactions.

Dual-labeled DNA probes:

Real-time PCR technology based on dual-labeled DNA probes provides a high sensitive and high specific PCR system with multiplexing capability. It requires two standard PCR primers and the DNA probe that hybridizes to an internal part of the amplicon. The sequence of the dual-labeled DNA probe should avoid secondary structure and primer-dimer formation.

Content:

qPCR ProbesMaster lowROX (red cap)

Hot Start Polymerase Ab+, dATP, dCTP, dGTP, dTTP, KCl, (NH₄)₂SO₄, MgCl₂, additives and stabilizers, 100 nM ROX

PCR-grade water



qPCR ProbesMaster lowROX

Master mix for real-time PCR with low ROX
2 x conc. master mix

Preparation of the qPCR master mix:

The preparation of a master mix is crucial in quantitative PCR reactions to reduce pipetting errors. Prepare a master mix of all components except template as specified. A reaction volume of 20-50 µl is recommended for most real-time instruments. Prepare 13 volumes of master mix for 12 samples or a triple-set of 4 samples. Pipet with sterile filter tips and minimize the exposure of the labeled DNA probe to light. Perform the setup in an area separate from DNA preparation or analysis. No-template controls should be included in all amplifications.

component	20 µl assay	50 µl assay	final conc.
qPCR ProbesMaster with lowROX	10 µl	25 µl	1x
primer forward (10 µM) ¹⁾	0.6 µl	1.5 µl	300 nM
primer reverse (10 µM) ¹⁾	0.6 µl	1.5 µl	300 nM
dual-labeled probe (10 µM) ²⁾	0.4 µl	1 µl	200 nM
template DNA	x µl	x µl	<500 ng/assay
PCR-grade water	fill up to 20 µl	fill up to 50 µl	-

¹⁾ The optimal concentration of each primer may vary from 100 to 500 nM.

²⁾ Optimal results may require a titration of DNA probe concentration between 50 and 800 nM.

Dispensing the master mix:

Vortex the master mix thoroughly to assure homogeneity and dispense the mix into real-time PCR tubes or wells of the PCR plate.

Addition of template DNA:

Add the remaining x µl of sample/template DNA to each reaction vessel containing the master mix and cap or seal the tubes/plate. Do not exceed 500 ng DNA per reaction as final concentration. Tubes or plates should be centrifuged before cycling to remove possible bubbles.

Recommended cycling conditions:

Initial denaturation and polymerase activation	95 °C	2 min	1x
Denaturation	95 °C	15 sec	35-45x
Annealing and elongation	60-65 °C ⁴⁾	1 min ⁵⁾	35-45x

⁴⁾ The annealing temperature depends on the melting temperature of the primers and DNA probe used.

⁵⁾ The elongation time depends on the length of the amplicon. A time of 1 min for a fragment of up to 500 bp is recommended.

For optimal specificity and amplification an individual optimization of the recommended parameters, especially of the annealing temperature may be necessary for each new combination of template DNA, primer pair and DNA probe.

Related Products:

Dual-labeled DNA probes
Custom primers