

PRODUCT NAME
One-Step RT-PCR Mastermix (ROX)

For research purposes only. Not for use in diagnostic procedures for clinical purposes. Please refer to the Certificate of Analysis for the expiration date and the lot number
FOR IN VITRO USE ONLY

Cat. No. GMO-RTWR-600

Product Description

The One-Step RT-PCR Mastermix is a HotStart Mastermix designed to simplify the preparation of a single tube reverse transcriptase real-time PCR assay. The Mastermix contains Hotstart Taq DNA polymerase, optimised reaction buffer, 5mM (final concentration) MgCl₂ and nucleotides (including dUTP), a tube of Reverse transcriptase combined with RNase Inhibitor and tube of a passive reference.

Storage and stability

The One-Step RT-PCR Mastermix reagents are stable for a minimum of twelve months when stored at -70°C in a constant temperature freezer. The Mastermix can be stored at -20°C but we recommend storage of the RT enzyme tube at -70°C.

Kit Contents

The Mastermix contains enough PCR reagents for up to 600 reactions (25µl).

Reagent	Volume
2X Reaction Buffer	7.5ml
50mM MgCl ₂	1.5ml
Reverse transcriptase combined with RNase Inhibitor	75µl
Passive reference (pink cap)*	0.5ml

* Passive reference stock solution for systems which require passive reference.

Materials Required But Not Supplied

The following items will be required in addition to the reagents supplied in the qPCR™ Mastermix.

Reagents :

Deionised or distilled water
 Unlabelled primers

Materials :

Optical 96-Well Reaction Plate
 Optical Tubes and Caps as needed
 Microcentrifuge
 Polypropylene tubes
 Pipettes, preferably positive displacement
 Pipette tips, with filter plugs
 Vortex
 Disposable Gloves

Usage

A typical 25µl PCR Mix would contain the following:

Component	Volume (µl)	Final concentration in 25µl
One-Step RT-PCR Mastermix	12.5	1x
Forward primer (1µM)	2.5	0.1 µM
Reverse primer (1µM)	2.5	0.1 µM
Probe	2.5	0.3 µM
Template	z	-
H2O	4.625- z	-
RT Enzyme/RNase Inhibitor	0.125	
Passive reference ((ROX)	0.25	3 µM
Total	25µl	

Typical cycling conditions for this kit are

1st strand reaction	30 min. 45°C
Hot Taq activation	10 min. 95°C
40 Cycles	15 sec. 95°C 60 sec. 60°C
Hold	25°C

Quality Control

The One-Step RT-PCR Mastermix has been tested for performance on an authorised real-time sequence detector.

Notice to purchaser : limited license

The purchase price of this product includes a limited, non-transferable license under US Patents 4.683.202, 4.683.195, 4.965.188 or their foreign counterparts, owned by Hoffmann-La Roche Inc. and F. Hoffmann-La Roche Ltd («Roche»), to use only this amount of the product to practice the Polymerase Chain Reaction («PCR») 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 an authorized thermal cycler. No right to perform or offer commercial services of any kind using PCR, including without limitation reporting the results of purchaser's activities for a fee or other commercial consideration, is hereby granted by implication or estoppel. Further information on purchasing licenses to practice the PCR process may be obtained by contacting the Director of Licensing at The Perkin-Elmer Corporation, 850 Lincoln Center Drive, Foster City, California 94404 or at Roche Molecular Systems Inc., 1145 Atlantic Avenue, Alameda, California 94501.

No rights are granted in the purchase of this product to perform 5' nuclease assays as covered under US patent or their foreign counterparts, owned by Hoffmann-La Roche Inc. and F. Hoffmann-La Roche Ltd.