

MESA GREEN qPCR MasterMix Plus for SYBR® Assay Low ROX

Technical Data Sheet

Reference: RT-SY2X-06+WOU LR

Products and procedures described in this protocol are intended for research purposes only.

Storage conditions

For long term storage the MESA GREEN qPCR MasterMix Plus for SYBR® Assay Low ROX should be stored at -15 °C to -25 °C in a constant temperature freezer. When stored under these conditions the reagents are stable for 1 year.

For short term storage the MESA GREEN qPCR MasterMix Plus for SYBR® Assay Low ROX can be stored at 4 °C to 6 °C for 1 month.

The MESA GREEN qPCR MasterMix Plus for SYBR® Assay Low ROX reaction buffer should be protected from light whenever possible.

Kit contents

The MESA GREEN qPCR MasterMix Plus for SYBR® Assay Low ROX contains enough reagents for up to 600 - 50 µl reactions using the hotstart enzyme, Meteor Taq.

Reagent	Volume	Description
2x reaction buffer (Red cap)* (amber bottle)	15 ml	Two bottles (7,5 ml each) of reaction buffer, dNTPs (including dUTP), Meteor Taq DNA polymerase, MgCl ₂ (4mM final concentration), SYBR® Green I, stabilizers and passive reference
50 mM MgCl ₂ (clear cap)	3 ml	One bottle of 50 mM MgCl ₂

*The SYBR® Green I included in the 2x reaction buffer is light sensitive and should be kept away from light as much as possible.

Procedure

1- Thaw all required reagents completely and put them on ice. Mix all reagents well by inversion and spin them down prior to pipeting.

2- Prepare the reaction mix

Component	Volume (µl)	Final Concentration
2x reaction buffer	25	1x
Forward primer	5	(initially 100 up to 300 nM)*
Reverse primer	5	(initially 100 up to 300 nM)*
Template	5	-
Water	10	(volume is 50 µl minus all other components)
Total Mix	50 µl	

*Note 1: the primer concentrations are recommended as starting concentrations; always start at the lower end. These concentrations will be correct for many assays, but additional optimization may be required to obtain the best results with your primer set.

Note 2: Uracil-N-Glycosylase (reference RT-0610-03) can be added to a final concentration of 0.01 U/µl if required (0.5 µl of 1 U/µl UNG per 50 µl reaction). If the UNG is required please add the following step at point 6 before the Meteor Taq activation step: 2 min. at 50 °C.

- 3- To correct for dispensing losses prepare an excess of reaction mix (for example 100 reactions reaction mix for 96 reactions). Add all components together, except for the template. Mix thoroughly by inversion. Spin down.
- 4- Pipette 5 µl of the template DNA for your samples, 5 µl of the control DNA for your positive control and 5 µl of water or buffer for your negative control in to your PCR tubes / 96-well plate / 384-well plate.
- 5- Add 45 µl of the reaction mix to the reaction vial, close the vial and mix gently on a stirrer or spin down. Ensure that no bubbles are present in the reaction vial. Reaction set up can be done at room temperature.
- 6a- If the Real Thermocycler used has a “normal” block, program it using the following recommended parameters:

Meteor Taq activation	5 min. 95 °C
40 Cycles	15 sec. 95 °C 1 min. 60 °C
Hold at 50 °C forever or perform a meltcurve	

- 6b- If the Real Thermocycler used has a “fast” block, then it is possible to program it using the following recommended parameters:

Meteor Taq activation	5 min. 95 °C
40/45 Cycles	3 sec. 95 °C 45-60 sec. 60 °C
Hold at 50 °C forever or perform a meltcurve	

Technical information

Primer design guidelines

- GC content should be between 30 % and 80 % (ideally 40-60 %)
- avoid runs of identical nucleotides, especially of 3 or more Gs or Cs at the 3' end
- using the Primer Express® software the T_m should be 58 °C to 60 °C

Custom assay design

Commonly used concentrations are 100 nM for primers. Optimal results may require titration of primers. The purpose of such a process is to determine the minimum amount of primers required to obtain the most sensitive results with your assay.

Primer titration matrix

Titrate according to the Table 1, perform qPCR and select the concentration, which gives the lowest Ct value. By doing this type of titration it is also possible to compensate for differences up to 2 °C in melt temperature of the primers.

Table 1: Primer titration matrix

Reverse	Forward		
	50 nM	100 nM	300 nM
50 nM	50 / 50	100 / 50	300 / 50
100 nM	50 / 100	100 / 100	300 / 100
300 nM	50 / 300	100 / 300	300 / 300

MgCl₂ adjustment matrix

Standard MgCl₂ concentration is 4 mM but optimal MgCl₂ concentration can vary between assay. If necessary use the 50 mM MgCl₂ tube. Always prefer optimizing the primer concentrations before the MgCl₂ concentration. Adjust the amount of water if MgCl₂ is added to the reaction.

Final MgCl ₂ concentration (mM)	MgCl ₂ to add (µl/50 µl)	2x reaction buffer (µl)
4	0	25
4.5	0.5	25
5	1	25
5.5	1.5	25
6	2	25

3-step protocol instead of 2-step protocol

Increasing extension time or performing a 3-step protocol can increase the ΔR_n and / or decrease the Ct of an assay, particularly when the PCR product is longer than 100 bp.

The protocol will be as follows:

Meteor Taq activation		5 min. 95 °C
40 Cycles	denaturation	15 sec. 95 °C
	annealing	20 sec. 60 °C
	extension	40 sec. 72 °C
Increase extension time with 10-second steps, if required.		

Further information available through Eurogentec web site, www.eurogentec.com.

- Troubleshooting Guide for qPCR and RTqPCR (under the “Technical Resources / Troubleshooting Guide” section).
- Primers and probe design (please refer to our Troubleshooting Guide).
- “Your One-stop-shop Real-Time qPCR supplier” handbook (under the “Technical Resources / Documentation” section).
- MSDS, (under the “Technical Resources / MSDS” section)
- Certificates of Analysis (please contact us).

For any further information required please contact our Customer Help Desk:

For Europe:

E-mail: info@eurogentec.com

Tel : +32 4 372 76 65

For USA:

E-mail: info.usa@eurogentec.com

Tel: +1 858 793 26 61

FOR RESEARCH USE ONLY

NOTICE TO PURCHASER: LIMITED LICENSE

Licensed dsDNA-Binding Dye Research Kits

Use of this product is covered by one or more of the following US patents and corresponding patent claims outside the US:

5,079,352, 5,789,224, 5,618,711, 6,127,155, 5,677,152, 5,773,258, 5,407,800, 5,322,770, 5,310,652, 5,994,056, 6,171,785, and claims outside the US corresponding to US Patent No. 4,889,818. The purchase of this product includes a limited, non-transferable immunity from suit under the foregoing patent claims for using only this amount of product for the purchaser's own internal research. No right under any other patent claim (such as apparatus or system claims in US Patent No. 6,814,934) and no right to perform commercial services of any kind, including without limitation reporting the results of purchaser's activities for a fee or other commercial consideration, is conveyed expressly, by implication, or by estoppel. This product is for research use only. Diagnostic uses under Roche patents require a separate license from Roche. Further information on purchasing licenses may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

Cyclic-substituted unsymmetrical cyanine dyes are covered by U.S. Patents 5,436,134 and 5,658,751 and licensed to Eurogentec S.A by Molecular Probes, Inc in the direct research field.

Use of UDG employs U.S. Patents 5,035,936, 5,945,313, 5,683,896 and their foreign counterparts licensed to Eurogentec, S.A from Invitrogen Corporation.

Primer Express® is a registered trademark of Applied Biosystems Corporation.