

PRODUCT NAME
Universal Mastermix 7,5ml
Please refer to the Certificate of Analysis for the expiration date and the lot number <b>FOR IN VITRO USE ONLY</b>
<b>Cat. No. DMMLD2D600</b>

**Product description**

The Universal Mastermix is a Hotstart Mastermix designed to simplify the preparation of a real-time PCR assay. The Universal Mastermix contains a Hotstart Taq NA polymerase, optimised reaction buffer, 4mM (final concentration) MgCl<sub>2</sub> and nucleotides (dNTPs). In separate tubes, passive reference (ROX), SYBR® Green, and FITC.

**Storage and stability**

Store the Universal Mastermix at -15 to -25°C in a constant temperature freezer. When stored under these conditions the reagents are stable for a minimum of 12 months.

**Kit Contents**

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

Reagent	Volume
2X Reaction Buffer dNTPs (including dUTP), Hotstart DNA polymerase, MgCl <sub>2</sub> (4mM final concentration)	7.5ml
Passive reference Pink Cap, 500µl stock solution	0.5ml
SYBR® Green Brown tube	0.5ml
1mM FITC solution Green Cap, 1000µl stock solution	1 ml

Depending the applications and the system prepare a **stock solution** following the instructions in the table enclosed. (Universal Mastermix). This stock solution will be used to prepare the PCR Mastermix (see next page).  
Passive reference stock solution for systems which require passive reference.  
FITC for Biorad users working with SYBR® Green.

**Materials Required But Not Supplied**

The following items will be required in addition to the reagents supplied in the Universal 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

**FOR SYBR GREEN**

**Prepare a stock solution by adding**

System	Master Mix 2X	Rox	SYBR Green	FITC	Denaturation time	PCR cycles	Normal Cycles	Fast Cycles
7000-7300-7900 (ABI)	1000µl	40 µl	60 µl	0 µl	10 minutes at 95C	40	15' at 95C; 60' at 60C	5' at 95C; 60' at 60C (7900 ONLY)
7500 (ABI)	1000µl	2.8 µl	60 µl	0 µl	10 minutes at 95C	40	15' at 95C; 60' at 60C	5' at 95C; 60' at 60C
LC480 (Roche)	1000µl	0 µl	60 µl	0 µl	10 minutes at 95C	40		5' at 95C; 60' at 60C
Icycler (Biorad)	1000µl	0 µl	60 µl	10µl of a 1000 dilution of the FITC solution	10 minutes at 95C	40	15' at 95C; 60' at 60C	
Rotorgene 6000 (Corbett)	1000µl	0 µl	60 µl	0 µl	10 minutes at 95C	40	15' at 95C; 60' at 60C	

Use 12.5 µl of this stock solution and add primers and DNA to reach 25µl

**FOR DOUBLE DYE PROBE TAQMAN™**

**Prepare a stock solution by adding**

System	Master Mix 2X	Rox	SYBR Green	FITC	Denaturation time	PCR cycles	Normal Cycles	Fast Cycles
7000-7300-7900 (ABI)	1000µl	40 µl	0 µl	0 µl	10 minutes at 95C	40	15' at 95C; 60' at 60C	5' at 95C; 60' at 60C (7900 ONLY)
7500 (ABI)	1000µl	2.8 µl	0 µl	0 µl	10 minutes at 95C	40	15' at 95C; 60' at 60C	5' at 95C; 60' at 60C
LC480 (Roche)	1000µl	0 µl	0 µl	0 µl	10 minutes at 95C	40		5' at 95C; 60' at 60C
Icycler (Biorad)	1000µl	0 µl	0 µl	0 µl	10 minutes at 95C	40	15' at 95C; 60' at 60C	
Rotorgene 6000 (Corbett)	1000µl	0 µl	0 µl	0 µl	10 minutes at 95C	40	15' at 95C; 60' at 60C	
Stratagene Mx3000P/3005P	1000µl	0 µl	0 µl	0 µl	10 minutes at 95C	40	15' at 95C; 60' at 60C	

Use 12.5 µl of this stock solution and add primers and DNA to reach 25µl

### PCR Reaction Mix

A typical 25 µl PCR Reaction Mix would contain the following:

Component	Volume (µl)	Final concentration (in 25µl)
<b>2x Stock solution</b>	12.5	1x
Forward primer	2.5	0.1µM
Reverse primer	2.5	0.1µM
Probe when required	2.5	0.3µM
Water	x	-
Template	5	-
<b>Total Mix</b>	<b>25µl</b>	

### PCR Temperature Cycling Conditions

ATTENTION 10min initial deprotection time:

Normal cycles	
Hotstart enzyme Activation	10min. 95°C
45 Cycles	15sec. 95°C 60sec. 60°C
Hold	25°C forever
Fast cycles	
Hotstart enzyme Activation	10min. 95°C
40 Cycles	5sec. 95°C 60sec. 60°C
Hold	25°C forever

### Quality Control

The Universal Mastermix has been tested for performance on an authorised real-time sequence detector. Products and procedures described in this protocol are intended for research purposes only.

### 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 cyler. 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.

#### Diagenode sa. BELGIUM | EUROPE

LIEGE SCIENCE PARK  
Rue Bois Saint-Jean, 3  
4102 Seraing (Ougrée) - Belgium  
Tel: +32 4 364 20 50  
Fax: +32 4 364 20 51  
orders@diagenode.com  
info@diagenode.com

#### Diagenode Inc. USA | NORTH AMERICA

400 Morris Avenue, Suite 101  
Denville, NJ 07834 - USA  
Tel: +1 862 209-4680  
Fax: +1 862 209-4681  
orders.na@diagenode.com  
info.na@diagenode.com

Last update: 2009