



# iProof™ High-Fidelity Master Mix

Catalog #	Kit	Kit Contents	Kit Size
172-5310	iProof HF Master Mix	Master mix — 2.5 ml (2 x 1.25 ml vials)	100 x 50 µl reactions
172-5320	iProof GC Master Mix	DMSO (100%) — (1 x 0.5 ml vial)	

For research purposes only.

## Storage and Stability

Guaranteed for six months in a constant temperature freezer at  $-20^{\circ}\text{C}$ . For convenience, the mix can be refrozen or stored at  $4^{\circ}\text{C}$  for three months.

## Kit Contents

iProof high-fidelity PCR master mixes are 2x concentrated, ready-to-use supermixes containing all components (0.04 U/µl iProof, 400 µM dNTPs, and 3 mM  $\text{MgCl}_2$ ), except primers and template, for high-fidelity, long, or fast PCR for amplicons up to 37 kb.

The iProof high-fidelity DNA polymerase in these master mixes incorporates a patented fusion protein technology for speed and accuracy, making this an ideal choice for cloning. The error rate of iProof polymerase is determined to be 50-fold lower than that of *Thermus aquaticus*, and six-fold lower than that of *Pyrococcus furiosus*. iProof polymerase produces blunt-end DNA products.

## Reaction Setup

1. Thaw supermix and other frozen reaction components at room temperature. Mix thoroughly, centrifuge briefly to collect solution at the bottom of tubes, and store on ice.
2. Prepare reaction mix on ice using PCR tubes according to the recommendations in Table 1.

Table 1. Reaction setup.		
Component	Volume per 50 µl Reaction*	Final Concentration
iProof HF or GC master mix (2x)	25 µl	1x
Forward and reverse primers	Variable	0.5 µM each primer (0.2 to 1.0 µM)
DNA template	Variable	Plasmid DNA, lambda, or BAC DNA: 10 ng–1 µg Genomic DNA: 500–50 ng
H <sub>2</sub> O	Variable	—
Total reaction mix volume	50 µl	—

Scale all components proportionally according to sample number and reaction volumes.

- Seal tubes and vortex 30 sec or more to ensure thorough mixing of the reaction components. Spin the tubes to remove any air bubbles and collect the reaction mixture in the vessel bottom.
- Program the PCR cycling protocol on the PCR instrument according to Table 2.

Step	Temperature	Time	Cycle
Initial denaturation	98 °C	30 sec–3 min	1
Denaturation	98 °C	5–10 sec	25–35
Annealing	45–72 °C (optimized temperature)	10–30 sec	
Extension	72 °C	15–30 sec/kb	
Final extension	72 °C	5–10 min	1

### Notes about Cycling Conditions and Reaction Components

- For difficult DNA templates, a 3 min initial denaturation time is recommended
- When using iProof master mix, a general rule is to anneal primers (>20 nt) for 10–30 sec at 3 °C above the primer with the lowest melting temperature ( $T_m$ ). Primer  $T_m$  should be calculated using the nearest-neighbor method, as results can vary significantly depending on the method used. For primers  $\leq 20$  nt, use an annealing temperature equal to the primer with the lowest  $T_m$
- Template extension temperature must be at 72 °C and extension time depends on amplicon length (15–30 sec/kb) and complexity. Do not exceed 1 min per kb for amplicons that are >5 kb
- The error rate in HF buffer is  $4.4 \times 10^{-7}$ , which is lower than the error rate in GC buffer ( $9.5 \times 10^{-7}$ )
- Adding 3% DMSO aids in template denaturation for GC-rich amplicons. Further optimization of DMSO should be made in 2% increments. Ten-percent DMSO will require lowering the annealing temperature by 5.5–6.0 °C

### Quality Control

iProof HF and GC master mixes demonstrate robust amplification of long amplicons. Stringent specifications are maintained to ensure lot-to-lot consistency.

To learn more about Bio-Rad's complete list of solutions for amplification, visit [www.bio-rad.com/amplification](http://www.bio-rad.com/amplification).

### NOTICE TO PURCHASER: LIMITED LICENSE

Use of this product is covered by one or more of the following U.S. patents and corresponding patent claims outside the U.S.: 5,789,224; 5,618,711; and 6,127,155, and claims outside the U.S. corresponding to U.S. Patent number 5,079,352. 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, no right to perform any patented method, 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.