

## iQ™ Supermix

100 x 50 µl reactions	170-8860
500 x 50 µl reactions	170-8862
1,000 x 50 µl reactions	170-8864

For research purposes only  
Store at -20°C

### Storage and Stability

Store the iQ Supermix at -20°C in a constant temperature freezer. Avoid repeated freeze/thaw cycles. When stored under these conditions the supermix is stable for one year after ship date. You may aliquot the supermix and store a portion at 4°C for ready use. At 4°C the supermix is stable for six months.

### Kit Contents

iQ Supermix is a 2X mix for real-time PCR applications, containing a proprietary buffer and stabilizers specifically optimized for qPCR using fluorescent-labeled probes for detection. This mix provides the highest level of specificity to reduce the occurrence or delay the detection of primer-dimer and other non-specific artifacts. The antibody-mediated hot-start iTaq™ DNA polymerase is activated after an initial three minute denaturation step at 95°C.

Reagent	Kit Size	Volume	Description
iQ SYBR Green (Violet Cap)	100 reactions	1.25 ml x 2	2X reaction buffer with dNTPs, iTaq DNA polymerase, 6 mM MgCl <sub>2</sub> , and stabilizers
	500 reactions	1.25 ml x 10	
	1,000 reactions	1.25 ml x 20	

iQ Supermix comes optimized for real-time PCR applications. If you would like to have extra MgCl<sub>2</sub>, a 50 mM MgCl<sub>2</sub> solution is available free of charge upon request. Please request catalog number 170-8872 for 1.25 ml of this solution.

### Quality Control

The iQ Supermix is free of contaminating DNase, RNase, exonuclease, and protease activities.

Functionally, the iQ supermix gives a dose response over four logs of input genomic DNA concentration (500 ng to 50 pg) in a real-time PCR assay using a single-copy gene, IL-1β.

## Reaction Set Up

Thaw all components at room temperature. Mix vigorously, and then centrifuge to collect contents to the bottom of the tube before using.

Component	Volume per reaction	Final concentration
iQ Supermix	25 $\mu$ l	1X
Primer 1	x $\mu$ l	100 nM–500 nM
Primer 2	x $\mu$ l	100 nM–500 nM
Probe	x $\mu$ l	100 nM–500 nM
Sterile water	x $\mu$ l	
DNA template	x $\mu$ l	
<hr/>		
Total Volume	50 $\mu$ l	

## Reagents and Materials Not Supplied

Microcentrifuge tubes, screw capped  
Pipette tips, preferably aerosol barrier tips  
Vortexer  
Microcentrifuge  
Optical plates/tubes  
Sterile water  
Reaction primers and probes  
Real-time PCR detection system

## Recommendations for Optimal Results

Due to the sensitivity of quantitative PCR, results can be easily affected by pipetting errors.

- Always prepare a master mix of iQ Supermix containing the primers and probe.
- Add the template DNA sample to aliquots of the master mix for optimal reproducibility of replicate samples.
- This allows you to pipet once into the sample well or tube.

Individual pipetting of replicate samples is not recommended.

To learn more about Bio-Rad's complete solution for amplification, visit our website:

**[www.bio-rad.com/amplification](http://www.bio-rad.com/amplification)**

### NOTICE TO PURCHASER: LIMITED LICENSE

A license to perform the patented 5' Nuclease Process for research is obtained by the purchase of (i) both Authorized 5' Nuclease Core Kit and Licensed Probe, (ii) a Licensed 5' Nuclease Kit, or (iii) license rights from Applied Biosystems.

This product is an Authorized 5' Nuclease Core Kit. 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, (Claims 1-23 only), 5,773,258, (claims 1 and 6 only), 5,407,800, 5,322,770, 5,310,652, 5,210,015, 5,487,972, 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. Separate purchase of a Licensed Probe would convey rights under the applicable claims on US Patents Nos. 5,538,848, 5,723,591, 5,876,930, 6,030,787, 6,258,569, 5,804,375 (claims 1-12 only), and 6,214,979, and corresponding claims outside the United States. No right under any other patent claim 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 from the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

iQ is a trademark of Bio-Rad Laboratories.