

---

# Authorization Statement

for Real-Time Thermal Cyclers: iQ™5, MyiQ™, MyiQ™2, Chromo4™, DNA Engine Opticon®, DNA Engine Opticon 2, MiniOpticon™, CFX96™, CFX96 Touch™, CFX384™, CFX384 Touch™, and CFX Connect™ Systems

---

## NOTICE TO PURCHASER

This real-time thermal cycler is licensed under U.S. Patent No. 6,814,934 and corresponding claims in any counterpart Canadian patent thereof owned by Applied Biosystems, for use solely in research and all other applied fields except veterinary in vitro diagnostics. No rights are conveyed expressly, by implication or estoppel to any patents on real-time methods, including but not limited to 5' nuclease assays, or to any patent claiming a reagent or kit. For further information on purchasing license rights, contact the Director of Licensing at Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

## NOTICE TO PURCHASER

These products are covered by one or more of the following U.S. patents or their foreign counterparts owned by Eppendorf AG: U.S. Patent Nos. 6,767,512 and 7,074,367.

**BIO-RAD**

Effective June 10, 2011