

Authorization Statement

for Real-Time Detection Modules: Chromo4™ Detection Module and iCycler iQ®, iQ™5, MyiQ™, MyiQ™2, DNA Engine Opticon®, DNA Engine Opticon 2, MiniOpticon™, CFX96™, and CFX384™ Real-Time Upgrades

NOTICE TO PURCHASER

This Chromo4, DNA Engine Opticon, or DNA Engine Opticon 2 detection module, when combined with a DNA Engine® thermal cycler for which the applicable real-time thermal cycler royalty fee has been paid, constitutes a real-time thermal cycler licensed under U.S. Patent No. 6,814,934 and corresponding claims in any Canadian counterpart patent thereof owned by Applied Biosystems Corporation, for use solely in research and all applied fields except veterinary in vitro diagnostics. These license rights are effective only if this detection module is combined with a Bio-Rad thermal cycler for which the applicable real-time thermal cycler royalty fee has been paid and not with any other thermal cycler. 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

This iCycler iQ, iQ5, MyiQ, or MyiQ2 detection module, when combined with an iCycler® thermal cycler for which the applicable real-time thermal cycler royalty fee has been paid, constitutes a real-time thermal cycler licensed under U.S. Patent No. 6,814,934 and corresponding claims in any Canadian counterpart patent thereof owned by Applied Biosystems Corporation, for use solely in research and all applied fields except veterinary in vitro diagnostics. These license rights are effective only if this detection module is combined with a Bio-Rad thermal cycler for which the applicable real-time thermal cycler royalty fee has been paid and not with any other thermal cycler. 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

This MiniOpticon detection module, when combined with an MJ Mini™ thermal cycler for which the applicable real-time thermal cycler royalty fee has been paid, constitutes a real-time thermal cycler licensed under U.S. Patent No. 6,814,934 and corresponding claims in any Canadian counterpart patent thereof owned by Applied Biosystems Corporation, for use solely in research and all applied fields except veterinary in vitro diagnostics. These license rights are effective only if this detection module is combined with a Bio-Rad thermal cycler for which the applicable real-time thermal cycler royalty fee has been paid and not with any other thermal cycler. 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

This CFX96 or CFX384 detection module, when combined with a C1000™ or C1000 Touch™ thermal cycler for which the applicable real-time thermal cycler royalty fee has been paid, constitutes a real-time thermal cycler licensed under U.S. Patent No. 6,814,934 and corresponding claims in any Canadian counterpart patent thereof owned by Applied Biosystems Corporation, for use solely in research and all applied fields except veterinary in vitro diagnostics. These license rights are effective only if this detection module is combined with a Bio-Rad thermal cycler for which the applicable real-time thermal cycler royalty fee has been paid and not with any other thermal cycler. 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