

Authorization Statement

for Thermal Cyclers: MJ Mini™, MyCycler™, and T100™ Cyclers

Purchase of this instrument, Serial No. _____, conveys a limited non-transferable immunity from suit for the purchaser's own internal research and development and for use in applied fields under U.S. Patent No. 5,475,610 (claims 1, 44, 158, 160–163, and 167 only), or corresponding claims in its non-U.S. counterpart, owned by Applied Biosystems. No right is conveyed expressly, by implication, or by estoppel under any other patent claim, such as claims to apparatus, reagents, kits, or methods such as 5' nuclease methods. 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.

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