

# ABI PRISM<sup>®</sup> BigDye<sup>®</sup> Terminator v3.1 Cycle Sequencing Kits



Product P/N 4336923, 4336921, 4336919, 4336917, 4336915  
Insert P/N 4336875B Printed in U.S.A.

**For Research Use Only.**  
**Not for use in diagnostic procedures.**

## Kit Contents:

ABI PRISM<sup>®</sup> BigDye<sup>®</sup> Terminator v3.1 Ready Reaction mix  
pGEM<sup>®</sup> - 3Zf(+) Control Template, 0.2 µg/µL  
21 M13 Control Primer (forward), 0.8 pmol/µL

For Research Use Only. Not for use in diagnostic procedures.

## NOTICE TO PURCHASER: LIMITED LICENSE

A license under the process claims of U.S. patents or their foreign counterpart claims, has an up-front fee component and a running-royalty component. The purchase price of a ABI PRISM<sup>®</sup> BigDye<sup>®</sup> Terminator v3.1 Cycle Sequencing Kit includes limited, non-transferable rights under the running-royalty component to use only this amount of the product to practice the DNA sequence and fragment analysis processes described in said patents when this product is used in conjunction with an Authorized DNA sequence analysis instrument whose use is covered under the up-front fee component of these patents. No other rights are hereby granted expressly, by implication, or by estoppel, or under any other patent rights owned or licensable by Applied Biosystems. Further information relating to the purchase of licenses for DNA sequence and fragment analysis and other applications may be obtained by contacting the Director of Licensing at Applied Biosystems, 850 Lincoln Centre Drive, Foster City, CA 94404, USA.

## NOTICE TO PURCHASER: LIMITED LICENSE

The purchase of the ABI PRISM<sup>®</sup> BigDye<sup>®</sup> Terminator v3.1 Cycle Sequencing Kit includes a limited, nontransferable, non-exclusive license (without the right to resell, repackage, or sublicense) under one or more US patents and corresponding foreign patent claims and patent applications, to use this product solely with an Applied Biosystems commercial automated DNA sequencing machine or other authorized automated DNA sequencing machine that has been authorized under these patents by Applied Biosystems. No license is hereby granted for the use of this kit or the reagents therein, in any other automated sequencing machine. Such license is granted solely for research and other uses that are not unlawful. No other license is granted expressly, impliedly, or by estoppel. 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 About Limited License

This kit (reagent) is sold pursuant to a limited sublicense from Amersham International plc under one or more U.S. Patent Nos. 5,498,523 and 5,614,365, and corresponding foreign patents and patent applications. The purchase of this kit (reagent) includes a limited non-exclusive sublicense (without the right to resell, repackage or further sublicense) under such patent rights to use this reagent for DNA sequencing or fragment length analysis solely with an Applied Biosystems commercial automated sequencing machine or other authorized DNA sequencing machines that have been authorized for such use by Applied Biosystems, or for manual DNA sequencing. No license is hereby granted for use of this kit, or the reagents therein, in any other automated sequencing machine. Such sublicense is granted solely for research or other uses that are not unlawful. No other license is granted expressly, impliedly, or by estoppel. For information concerning the availability of additional license to practice the patented methodologies, contact: Amersham Life Science, Inc., Vice President, Regulatory Affairs, P.O. Box 22400, Cleveland, Ohio 44122. Patents are pending in countries outside the United States.

## NOTICE TO PURCHASER: LIMITED LICENSE

Use of the ABI PRISM<sup>®</sup> BigDye<sup>®</sup> Terminator v3.1 Cycle Sequencing Kit is covered by US patent claims and patent claims outside the US. 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 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.

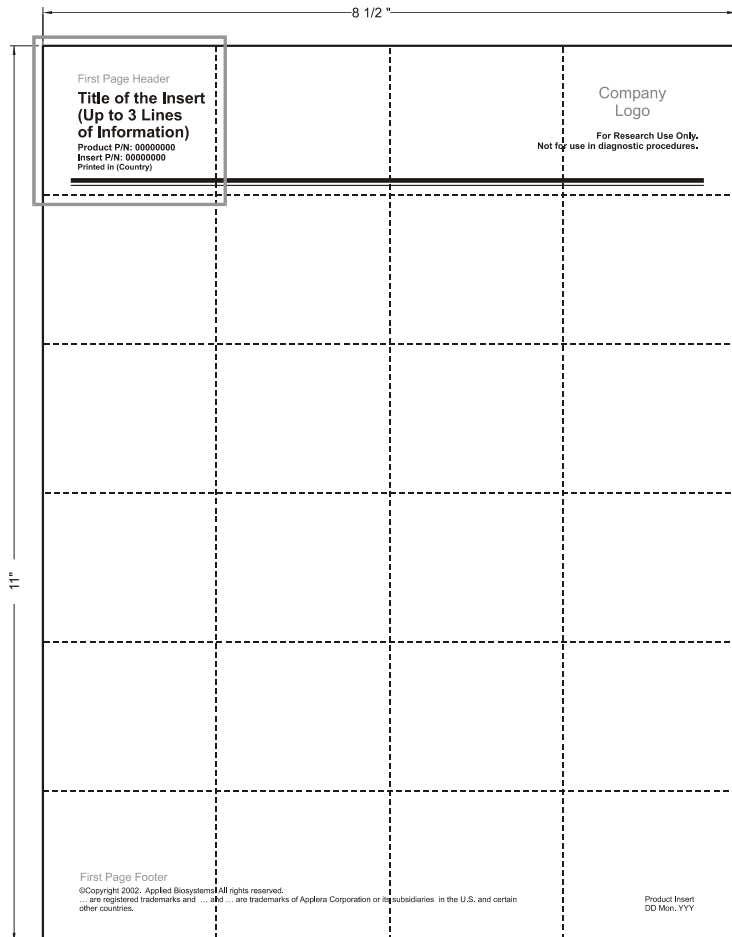
# GUIDELINE, INSERT FOLD INSTRUCTIONS



Notes :

1. Specification is intended for full and half sheet inserts, two sheet max.
2. Material: 50# Offset (White)
3. Dimensions: Full Sheet (8 1/2" x 11") folded size = 1.833" x 2.125" +/- .125"
4. Dimensions: Half Sheet (5 1/2" x 8 1/2") folded size = 1.833" x 2.125" +/- .125"
5. Two sheet inserts to have pages 2 and 3 facing each other.
6. After final folding, upper left-hand corner of page 1 is to face out, with title and part number displayed.
7. Folded inserts to be packaged as follows:
  - Bundled in groups of 10ea.
  - Shrink-wrap 10ea groups for 100 inserts total.
  - Label shrink-wrapped bundles [Part# & Rev (font large enough to be legible at 8 feet), Qty, Manufacturer, and Barcode (#128), if appropriate].
  - Seal shrink-wrapped bundles in an appropriately sized corrugated shipper (not to exceed 25#, or have outside dimensions greater than 12" x 13 1/2" x 4").
  - Appropriately label shipper [Part# & Rev (font large enough to be legible at 8 feet), Description, P.O.#, Qty, Mfg Date, Manufacturer, Country of Origin, and Barcode (#128) if appropriate] on at least one end panel (12" x 13 1/2" side).

Folding instructions ARE NOT and MUST NOT BE PRINTED as part of this insert.



THIS DOCUMENT CONTAINS PROPRIETARY INFORMATION OF APPLIED BIOSYSTEMS AND IS TO BE RETAINED IN CONFIDENCE, NOT TO BE REPRODUCED, COPIED, OR INCORPORATED IN ANY MANNER.  
© APPLIED BIOSYSTEMS 2002, 2010