

GeneScan™ 500 ROX SIZE STANDARD

Product P/N 4310361, 4310360
Insert P/N 4363111 REV A
Printed in USA



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

GeneScan™ -500 ROX® is an internal lane size standard developed for use with the Applied Biosystems fluorescence-based DNA electrophoresis systems. The use of an internal lane size standard enables automated data analysis and is also essential for achieving high run to run precision in sizing DNA fragments by electrophoresis. GeneScan™ -500 ROX® Size Standard is designed for sizing DNA fragments in the 35-500 bp range and provides 16 single-stranded labeled fragments of 35, 50, 75, 100, 139, 150, 160, 200, 250, 300, 340, 350, 400, 450, 490, and 500 bases. Each of the DNA fragments is labeled with a proprietary fluorophore, which results in a single peak when run under denaturing or native conditions.

Each kit contains two-200 µL tubes of the size standard and 1 tube of loading buffer. This is sufficient for 800 analyses when using the recommended loading amount of 0.5 µL. The kit may be stored up to six months at 2°C to 4°C. (Do not freeze)

Directions for Use

WARNING! CHEMICAL HAZARD. Hi-Di™ Formamide. Exposure causes eye, skin, and respiratory tract irritation. It is a possible developmental and birth defect hazard. Read the MSDS, and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

Sample Preparation

- 1) Before use, mix the contents of each tube thoroughly and centrifuge briefly to collect the liquid at the bottom of the tube. Typical loading cocktails are described below.

	377	310	3130 and 3100 Series Systems	3700	3730 Series Systems
Sample	0.4	0.5	0.5	0.5	0.5
Size Standard	0.4	0.5	0.25	0.5	0.5
Hi-Di™ Formamide	2.1	9.0	9.25	9.0	9.0
Loading Buffer	2.0	na	na	na	na
Loading volume	1.5	na	na	na	na

Note: Applied Biosystems recommends using the above ratios of DNA sample (PCR product) and size standard as a starting point only. Optimize these ratios if necessary, based on your experimental results.

The Hi-Di™ Formamide (4311320) is not included in the kit.

Heat the loading cocktail for 3 minutes at 95°C. Immediately chill on ice for a few minutes and load samples.

*The 250 bp peak is sensitive to small temperature variations on capillary electrophoresis instruments. The 250 bp fragment should not be used when defining the size standard in GeneScan™ or GeneMapper® Software.

Also on capillary electrophoresis instruments, the 340 bp peak is subject to large temperature variations. Fragment analysis primer peaks can often interfere with the detection of the 35 bp peak. Optimize your analysis based on this information.

Notice to Purchaser: Disclaimer of License

This product is optimized for use in the DNA sequencing or fragment analysis methods covered by patents owned or licensable by Applied Biosystems. No license under these patents to use the DNA sequencing or fragment analysis methods is conveyed expressly or by implication to the purchaser by the purchase of this product. A license to use the DNA sequencing or fragment analysis methods for certain research and development activities accompanies the purchase of certain Applied Biosystems reagents when used in conjunction with an authorized DNA sequencing machine, or is available from Applied Biosystems. Further information on purchasing licenses to practice the DNA sequencing or fragment analysis methods may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.



©Copyright 2004. Applied Biosystems All rights reserved.
Applied Biosystems and GeneMapper are registered trademarks and AB (Design), Applera, Hi-Di, GeneScan, and ROX are trademarks of Applera Corporation or its subsidiaries in the U.S. and/or certain other countries.

Product Insert
30/Nov/2004

VENDOR INFORMATION,
PRODUCT INSERT
FULL SHEET

Insert P/N 4363111 REV A
Printed in USA



GUIDELINE, INSERT FOLD INSTRUCTIONS

- 1 Specification is intended for Full Sheet, two sheets max.
- 2 Material: 50# Offset (White)
- 3 Dimensions: Full Sheet (8 1/2" x 11") folded size = 1.833" x 2.125" +/- .125"
- 4 Two sheet inserts to have pages 2 and 3 facing each other.
- 5 After final folding, upper left-hand corner of page 1 is to face out, with title and part number displayed.
- 6 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), Qty, Manufacturer, and Barcode Part #128C].
 - 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 24").
 - 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 Part # in 128C] on at least one end panel (12" x 13 1/2" side).

NOTICE TO PRINTER

Print page(s) prior
to this page **ONLY**.

This page is
folding instructions
which is NOT and
MUST NOT BE PRINTED
as part of this insert.

8 1/2"			
First Page Header Title of the Insert (Up to 3 Lines of Information) Product P/N: 00000000 Insert P/N: 00000000 Printed in (Country)		Company Logo For Research Use Only. Not for use in diagnostic procedures.	
First Page Footer ©Copyright 2002. Applied Biosystems. All rights reserved. ... are registered trademarks and ... are trademarks of Applied Biosystems Corporation or its subsidiaries in the U.S. and certain other countries.			Product Insert DD Mon. YYY
11"			

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 2004