

Product P/N 4315974
Insert P/N 4363141 REV A
Printed in USA

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

The Matrix Standard Set DS-01 is used to generate the "multi-component matrix" required when analyzing dROX-, dTAMRA-, dR6G-, and dR110-labeled DNA fragments on the Applied Biosystems 3130 and 3100 Series Systems. The Data Collection Software uses the multi-component matrix to automatically analyze the four different colored fluorescent dye-labeled samples in a single capillary.

This kit contains one tube of Matrix Standard, which is sufficient for eight 16-capillary array runs. The Matrix Standard contains four sizes of DNA fragments labeled with four different colors. These standards are diluted in 1X TE buffer and are stable for one year when stored at 2°C to 8°C. (Do not freeze.)

Preparing the Matrix Standard Set DS-01 for the Applied Biosystems 3130 and 3100 Series Systems.

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.

1. Thoroughly mix the contents of the Matrix Standard tube and spin briefly in a microcentrifuge.
2. Dilute 5 µL of the tube contents with 195 µL of Hi-Di™ Formamide (P/N 4311320) in a 1.5 mL microcentrifuge tube.
3. Mix thoroughly and spin briefly in a microcentrifuge.
4. Denature and dispense reagents into a 96-well microtiter plate:

For convenience, we recommend dispensing the contents of the tube into a 96-well microtiter plate first, and then using a GeneAmp® PCR System 9600 thermal cycler for denaturation.

If a GeneAmp 9600/9700 thermal cycler is available for denaturation, follow steps A and B below.

- A) Dispense 10 µL of the Matrix Standard / Hi-Di™ formamide mixture into two columns (16 wells) of a 96-well microtiter plate.
- B) Cover the plate and denature at 95°C for 5 minutes. Immediately place on ice.

If a GeneAmp 9600/9700 thermal cycler is not available, follow steps C and D below.

- C) Heat the mixture at 95°C for 5 minutes to denature, and immediately place on ice.
- D) Dispense 10 µL of the denatured mixture into two columns (16 wells) of a 96-well microtiter plate.

5. Place the 96-well microtiter plate on the plate deck of the instrument.
6. For specifics on setting up a run, refer to your User's Manual or Getting Started Guide.

Ancillary Products Disclaimer

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.

