



## AFLP® Pre-amp Primer Mix I

Cat. No.: 10792-018

Size: 1 ml

Store at -20°C.

### Description:

AFLP® Pre-amp Primer Mix I is provided as an adjunct product to the AFLP® Analysis System I (Cat. No. 10544-013) and AFLP® Starter Primer Kit (Cat. No. 10483-014) for plant genomic DNA fingerprinting using the Amplified Restriction Fragment Polymorphism (AFLP®) technology. In AFLP®, Pre-amp Primer Mix I is used in the preamplification reaction to amplify genomic DNA's following restriction digestion. It consists of a mixture of two AFLP® primers, each having one selective nucleotide (M-C; E-A) and dNTPs. The PCR product from preamplification is used as a template for a subsequent selective amplification reaction. AFLP® Pre-amp Primer Mix I is designed for use with plants having genome size ranges of either  $1$  to  $5 \times 10^8$  bp or  $5 \times 10^8$  to  $6 \times 10^9$  bp including tomato, corn, soybean, lettuce, barley, cotton, oilseed rape, potato, sunflower, pepper and brassica.

### Component:

50142 AFLP® Pre-amp Primer Mix I

### Quality Control:

This product was demonstrated to be functional in standard amplification reactions.

### Protocol:

Refer to the manual provided with the AFLP® Analysis System I and the AFLP® Starter Primer Kit for detailed protocols including guidelines for primer pair selection and a troubleshooting guide.

### Limited Label License No. 17:

The AFLP® technique is covered by patents or patented applications owned by Keygene n.v. This product is sold under license from Keygene n.v. This kit may be used for research purposes only. For use of this kit in plant breeding, contact the Director of Licensing, Invitrogen Corporation, 1600 Faraday Avenue, Carlsbad, California 92008. Phone (760) 603-7200. Fax (760) 602-6500. The use of this kit for any other purpose, whether commercial or noncommercial, including but not limited to the use for clinical, diagnostic, and/or therapeutic purposes; or for providing services to third parties, requires a license from Keygene n.v., P.O. Box 216, 6700 AE Wageningen, The Netherlands.

Doc. Rev: 08/20/01

This product is distributed for laboratory research only. CAUTION: Not for diagnostic use. The safety and efficacy of this product in diagnostic or other clinical uses has not been established.

For technical questions about this product, call the Invitrogen Tech-LineSM U.S.A. 800 955 6288