



## AFLP<sup>®</sup> Primer E-AGG

**Cat. No.: 10760-015**

**Size: 46 µl**

**Conc: 27.8 ng/µl**

**Store at -20°C.**

Description:

AFLP<sup>®</sup> *Eco*R I Primer E-AGG is provided as an adjunct product to the AFLP<sup>®</sup> Analysis System I for plant genomic DNA fingerprinting using the Amplified Restriction Fragment Polymorphism (AFLP<sup>®</sup>) technology. In AFLP<sup>®</sup>, selective nucleotides are added to the 3'-end of the PCR primers such that only a subset of preamplified restriction fragments are recognized and amplified. These amplified fragments are then analyzed by denaturing polyacrylamide gel electrophoresis to generate the DNA fingerprint. AFLP<sup>®</sup> Primer E-AGG is designed for the identification of polymorphisms in plants having genome size ranges of  $5 \times 10^8$  to  $6 \times 10^9$  bp when paired with AFLP<sup>®</sup> *Mse* I Primers containing 3 selective nucleotides.

Components:

50137 AFLP<sup>®</sup> Primer E-AGG

Quality Control:

The primer was demonstrated to be functional in a standard AFLP<sup>®</sup> assay.

Protocol:

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

Limited Label License No. 17:

The AFLP<sup>®</sup> 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.: 102601

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