

AFLP® Core Reagent Kit

Cat. No.: 10482-016

Size: 50 templates

Store at -20°C.

Description:

Amplified Restriction Fragment Polymorphism (AFLP®) is a powerful technique used for DNA fingerprinting (1,2). The AFLP® Analysis Systems contain both the AFLP® Core Reagent Kit and an AFLP® Primer Kit designed for a specific application. The AFLP® Core Reagent Kit contains reagents for template preparation, including restriction enzymes, adapters, ligase and appropriate buffers. Primers, a manual and other reagents required for the AFLP® amplification reactions are contained in the AFLP® Primer Kits.

Components	Part No.	Amount
<i>EcoR</i> I/ <i>Mse</i> I [1.25 units/μl each in 10 mM Tris-HCl (pH 7.4), 50 mM NaCl, 0.1 mM EDTA, 1 mM DTT, 0.1 mg/ml BSA, 50% glycerol (v/v), 0.1% Triton® X-100]	51114	100 μl
5X Reaction Buffer [50 mM Tris-HCl (pH 7.5), 50 mM Mg-acetate, 250 mM K-acetate]	51115	250 μl
Distilled Water	50837	1.25 ml
Adapter/Ligation Solution [<i>EcoR</i> I/ <i>Mse</i> I adapters, 0.4 mM ATP, 10 mM Tris-HCl (pH 7.5), 10 mM Mg-acetate, 50 mM K-acetate]	51116	1.2 ml
T4 DNA Ligase [1 unit/μl in 10 mM Tris-HCl (pH 7.5), 1 mM DTT, 50 mM KCl, 50% (v/v) glycerol]	Y01301	50 μl
TE Buffer [10 mM Tris-HCl (pH 8.0), 0.1 mM EDTA]	50282	4.5 ml
Tomato DNA (100 ng/μl)	51117	10 μl
<i>Arabidopsis</i> DNA (100 ng/μl)	51118	10 μl

Quality Control:

All components were tested in the following protocol using the control DNA. Template prepared in this protocol was amplified using the primers and reagents provided in AFLP® Primer Kits and displayed the specified DNA fingerprint.

Instructions for Use:(For more detailed information, please refer to the AFLP® Analysis System manual)

1. Add 100-500 ng of your sample DNA (volume of sample DNA must be 18 μl or less) to a 1.5-ml microcentrifuge tube. Add 5 μl 5X Reaction Buffer, 2 μl *EcoR* I/*Mse* I and Distilled Water to a final volume of 25 μl.
2. For control reaction: Combine 2.5 μl control DNA, 5 μl 5X Reaction Buffer, 2 μl *EcoR* I/*Mse* I and 15.5 μl Distilled Water to a 1.5-ml microcentrifuge tube.
3. Mix gently and collect reaction by brief centrifugation. Incubate the mixture 2 h at 37°C.
4. Incubate the mixture for 15 min at 70°C to inactivate the restriction endonucleases. Place the tube on ice and collect contents by brief centrifugation.

Doc. Rev. 102601

This product is distributed for laboratory research use 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 800 955 6288

5. To a tube of digested DNA from Step 3, add 24 µl Adapter/Ligation Solution and 1 µl of T4 DNA Ligase.
6. Mix gently at room temperature, centrifuge briefly to collect contents, and incubate at 20°C for 2 h.
7. Perform a 1:10 dilution of the ligation mixture as follows:
Take 10 µl of the reaction mixture and transfer to a 1.5-ml microcentrifuge tube.
Add 90 µl TE buffer and mix well.
8. Diluted ligation mixture is then used for pre-amplification using AFLP® primers and other reagents provided in AFLP® Primer Kits.
9. The unused portion of the reaction mixture may be stored at -20°C.

References:

1. Zabeau, M. and Vos, P. (1993) *European Patent Application*, publication number EP 0534858.
2. Lin, J.J. and Kuo, J. (1995) *Focus*® 17, 66.

AFLP® is a registered trademark of Keygene n.v.

Triton® is a registered trademark of Rohm & Haas, Co.