



## **30–330-bp AFLP® DNA Ladder**

**Cat. No. 10832-012**

**Size: 300 applications**

**Store at -20°C**

### **Description**

The 30–330-bp AFLP® DNA ladder consists of 32 10-bp repeats. Additionally, there are diffuse bands at 10 bp and 20 bp, plus a fragment at 1668 bp. This ladder is specifically designed for sizing AFLP® DNA fragments and can be visualized on a 5–6% urea-polyacrylamide gel following electrophoresis of the end-labeled product.<sup>1,2</sup> Because both DNA strands are the same nucleotide composition, the denatured product produces a set of single-stranded oligonucleotides increasing in length by 10-bp increments. This product is easily radiolabeled using T4 polynucleotide kinase.

### **Storage Buffer**

10 mM Tris-HCl (pH 7.5)

1 mM EDTA

### **Recommended Protocol:**

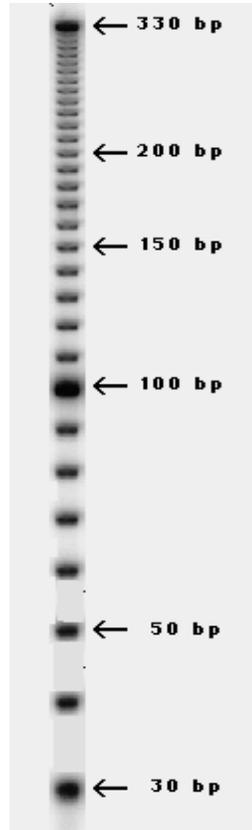
30–330-bp AFLP® DNA ladder is end-labeled with <sup>32</sup>P or <sup>33</sup>P with T4 polynucleotide kinase using the exchange reaction in the following protocol. Apply approximately 2 µl of ladder per lane on a 5–6% urea-polyacrylamide gel.

Part No. 10832012.pps

Rev. Date: 03/13/03

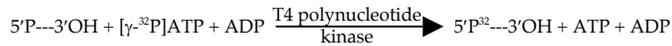
This product is distributed for laboratory research only. CAUTION: Not for diagnostic use. The safety and efficacy of this product in diagnostic or other clinical uses has not been established.

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

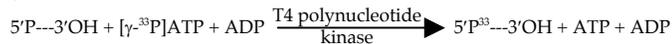


**Note:** The 100-bp band will appear two to three times brighter than other bands.

30–330-bp AFLP<sup>®</sup> DNA Ladder  
labeled with <sup>32</sup>P  
2 µl ladder / reaction  
6% polyacrylamide/urea gel

**Labeling of the 30–330-bp AFLP® DNA ladder with <sup>32</sup>P or <sup>33</sup>P**

or



1. Pipet into a 0.5-ml microcentrifuge tube on ice:
  - 2  $\mu$ l 30–330-bp AFLP® DNA ladder
  - 1  $\mu$ l 5X Exchange Reaction Buffer [250 mM imidazole (pH 6.4), 60 mM MgCl<sub>2</sub>, 5 mM 2-mercaptoethanol, 350  $\mu$ M ADP]
  - 1  $\mu$ l  $\gamma\text{--}^{32}P$  ATP (>3,000 Ci/mmole, 10  $\mu$ Ci/ $\mu$ l) or 1  $\mu$ l  $\gamma\text{--}^{33}P$  ATP (>3,000 Ci/mmole, 10  $\mu$ Ci/ $\mu$ l)
  - 1  $\mu$ l T4 polynucleotide kinase (10 U/ $\mu$ l)
2. Make sure all components are at the bottom of tube. Mix thoroughly but not vigorously. Centrifuge briefly.
3. Incubate for 10 min at 37°C.
4. Stop the reaction by heating the tube 15 min at 65°C.
5. This reaction will yield specific activities of approximately  $2 \times 10^6$  cpm/ $\mu$ l.
6. To 5  $\mu$ l of 30–330-bp AFLP® DNA ladder reaction mix, add an equal volume of TE buffer, pH7.5, and 25  $\mu$ l denaturing solution [98% (v/v) formamide, 10 mM EDTA (pH 8.0), 0.025% (w/v) bromophenol blue, 0.025% (w/v) xylene cyanol]. Total volume = 35  $\mu$ l.
7. Incubate at 70°C for 5 min.
8. Electrophorese 2  $\mu$ l of the sample in a denaturing 5–6% polyacrylamide/urea gel following the same protocol for AFLP® analysis. A set of single-stranded oligonucleotides (30–330 nucleotides) increasing in length of 10 nucleotide increments should be clearly visible after exposed in x-ray film.

**Note:** Diffuse bands at 10 bp and 20 bp are also visible.

**Quality Control**

Electrophoresis of the 30–330-bp AFLP® DNA ladder on a 4% Low Melting Point agarose gel shows that the bands between 30 bp to 330 bp are distinguishable. The 100-bp band is two to three times brighter than the rest of the bands.

**Additional Products**

<u>Product</u>	<u>Catalog no.</u>
T4 Polynucleotide Kinase	18004-010/18004-028
T4 Polynucleotide Kinase Exchange Buffer	10456-010
AFLP® Analysis System I	10544-013
AFLP® Analysis System II	10717-015

**References**

1. Vos, P., et. al. (1995) *Nucleic Acids Res.* 23, 4407.
2. Lin, J.-J., et. al. (1995) *FOCUS*® 17, 66.

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