

LI-COR AFLP® Systems

Cost Analysis of Samples

There are five basic steps in the AFLP® process:

- 1) Restriction digest of genomic DNA.
- 2) Ligation of oligonucleotide adapters.
- 3) Pre-selective amplification.
- 4) Selective amplification.
- 5) Electrophoresis of the fragments on a LI-COR DNA Analysis instrument.

1) - LI-COR Template Preparation Kit - Used for steps **1-3** of the AFLP® process and prepares 100 templates in large enough quantities to do approximately 5000 reactions each.

2) - LI-COR Selective Amplification Kit - Used for step number **4** of the AFLP® process. This kit selectively amplifies 25 different templates with 64 different primer combinations for a total of 1600 reactions.

	Cost (\$)	# of Templates	# of Reactions (Samples) per Template	# of Reactions (Samples)	Cost (\$) to prepare each Template	Cost (\$) per Reaction (Sample)
1. Template Preparation Kit	\$300	100	5000	500,000	\$3.00	NA
2. Selective Amplification Kit	\$450	25	64	1600	NA	\$0.28
Total	\$750					~ \$0.28

3) - LI-COR Startup Kit - Includes both a template preparation kit and a selective amplification kit at an economical price.

	Cost (\$)	# of Templates	# of Reactions	Cost (\$) per Template	Cost (\$) per Reaction
LI-COR Startup Kit	\$675	Same as above	Same as above	Same as above	\$0.25

LI-COR AFLP® Systems

Throughput of Samples

	# of Sample Lanes	# of Samples 700 Channel	# of Samples 800 Channel	Total # of Samples per Gel	# of Gels per 8 Hour Day	# of Runs per gel	Total # of runs per 8 Hour Day	# of Samples per 8 Hour Day
High Throughput	96	96	96	192	1	2	2	384
Medium Throughput	64	64	64	128	1	2	2	256
Low Throughput	48	48	48	96	1	2	2	192

The following is a list of some key reasons why researchers purchase LI-COR's system for their AFLP® research.

- 1) 110 nm separation between IRDyes™ eliminates spectral overlaps seen in four dye instruments
- 2) Direct access to our genomics technical support team via phone and e-mail
- 3) LI-COR is licensed for AFLP® technology development through Keygene, N.V.
- 4) AFLP® Kits are quality controlled for consistent, reliable performance on LI-COR instruments
- 5) LI-COR produces the only AFLP® Expression Analysis Kit on the market
- 6) To be able to immediately extract bands from the gel using our Odyssey Infrared Imaging System in either current research, or future research
- 7) Powerful Saga MX software is designed specifically for the AFLP® researcher, eliminating hours of analysis time as found in other systems

AFLP® is a registered trademark of KeyGene, N.V.
IRDyes is a registered trademark of Li-cor Biosciences