

Gene Copy Number analysis using semi-quantitative multiplex PCR-based assay on capillary electrophoresis systems

Stéphane Jankowski, Lichen Xu, Erica Currie-Fraser
Applied Biosystems, 850 Lincoln Centre Drive, Foster City, CA, 94404 USA.



INTRODUCTION

Large genomic rearrangements such as duplications and deletions have been recognized as pathogenic mutations for many diseases. These types of mutations are thought to represent 5.5% of reported mutations⁽¹⁾. However, given that mutation scans have not included searches for deletions and duplications, it seems likely that these figures are an underestimate of the actual number⁽¹⁾. Detection of genomic rearrangements is technically challenging and is typically done using techniques such as Southern blot analysis or Fluorescent In Situ Hybridization (FISH). These techniques often require high quantities of DNA or can be time-consuming and laborious, therefore limiting the efficiency of molecular screening. To better facilitate the detection of such Copy Number Variations (CNVs), researchers have developed simple, semi-quantitative methods using a multiplex PCR-based assay of short fluorescent fragments^(1,2) on Applied Biosystems capillary electrophoresis (CE) platforms. In this poster, we highlight this technique on the Applied Biosystems 3730/3730xl DNA Analyzers.

MATERIALS AND METHODS

CNV assays consist of the simultaneous amplification and fluorescent labeling of short, specific DNA loci, using a limited number of cycles to allow accurate quantitation in the exponential amplification range. Each multiplex PCR will yield a pattern composed of fluorescent peaks, with each peak corresponding to a specific DNA locus. The comparison of fluorescence is done between the same peaks generated from different samples, controls and suspect samples (Figure 1).

Targeted loci for BRCA1, BRCA2, MLH1/MSH2 genes and 9p21 region were amplified using dye-labelled primers from DNA that had been isolated from blood. Amplified samples were then run on an Applied Biosystems capillary electrophoresis platform, and the data was analyzed using GeneMapper® Software version 4.0.

GeneMapper® Software version 4.0 provides new features especially useful for Copy Number analysis:

- 1. AFLP® Analysis Method** - for sample normalization by sum of signal (Figure 2)
- 2. Report Manager** - for sample to sample ratio calculation, and to append a "normal", "deletion" or "duplication" result. Vertical calculation divides the peak height of the sample amplicon by that of the same amplicon in the control. An example is shown in Figure 3. Note that the number of rows will depend on the number of samples that are being assayed. Also, the row location of the control sample needs to be specified as indicated in Figure 3. In the example shown, the control sample was the first sample in a set of 8. Set up an analysis, as shown in Figure 4, by setting an appropriate threshold to identify candidate samples with deletions or duplications in a specific genomic region.
- 3. Dye Scale** - for graphic normalization, preferred for low-throughput analysis or for some specific signal adjustment.

Figure 1. Example of deletions after control amplicon normalization. Reduction of the peak heights corresponds to the deleted regions, while an increase of the peak heights suggests a partial duplication of a given region. The peaks are being displayed using the custom plot colors feature in GeneMapper® Software.

