

Heteroduplex Analysis of Cystic Fibrosis Samples on the DCode™ System

Paul Zoller and Theresa Redila-Flores, Bio-Rad Laboratories, Inc., Hercules, CA 94547 USA

Introduction

Heteroduplex analysis (HA) is a commonly used mutation screening method due to its simplicity. The technique is based on conformational differences in double-stranded DNA caused by the formation of heteroduplex molecules (Nagamine et al. 1989). Heteroduplex molecules have a mismatch in the double strand, causing a distortion in its usual conformation. This distortion or altered conformation can be detected on polyacrylamide gels due to slower migration than the corresponding homoduplex molecules. Heteroduplexes are generated during the polymerase chain reaction (PCR) of a heterozygous individual or by adding mutant and wild-type DNA in the same PCR reaction or by denaturing and renaturing a mixture of mutant and wild-type DNA. Both mutant and wild-type samples are run on the same gel to analyze differences in mobility. Heteroduplex molecules with as little as one mismatch can show a different mobility in a gel from homoduplex molecules. Polyacrylamide analogs have been developed, such as DEM™ (detection-enhancing matrix) from Bio-Rad, which enhance the ability to detect mutations in heteroduplex samples when compared to conventional polyacrylamide gels (White et al. 1992).

In this experiment, we show that HA on the DCode universal mutation detection system can be used to analyze mutations in the cystic fibrosis gene.

Methods

The test samples consist of wild-type and mutant DNA samples from the cystic fibrosis gene exons 7 and 10. The heterozygous mutations from exon 7 were 1154insTC, two base insertion (+TC) and $\Delta F311$, three base deletion (-TTC). The mutations from exon 10 were a heterozygous $\Delta F508$, three base deletion (-CTT), a heterozygous compound sample Q493 (C to T) and $\Delta F508$ (-CTT), a homozygous $\Delta F508$ (-CTT), and a heterozygous $\Delta I507$, three base deletion (-ATC). Samples were provided by Dr L Silverman, University of North Carolina School of Medicine (Chapel Hill, NC). Genomic DNA from both wild-type and mutant samples were amplified by PCR to create end products of 289 bp for exon 7 and 369 bp for exon 10.

A 16 cm x 20 cm, 0.75 mm thick 1x DEM gel made up in 0.6x TBE buffer (54 mM Tris, 54 mM boric acid, 1.2 mM EDTA) was used. A 1.5 L volume of 0.6x TBE buffer was added to the lower electrophoresis tank, and 350 ml of 0.6x TBE buffer was added to the upper chamber in the DCode system. Five μ l (200–300 ng) of each amplified sample was mixed with 5 μ l 2x gel loading dye (70% glycerol, 0.05% Bromophenol Blue, 0.05% Xylene Cyanol, 2 mM EDTA) and electrophoresed on the DCode system at 100 V for 16 hr at room temperature. After electrophoresis, the gel was stained in 50 μ g/ml ethidium bromide in 0.6x TBE buffer for 5 min and destained in buffer for 10 min. The gel was imaged under ultraviolet (UV) transillumination.

Results and Discussion

Figure 1 shows the mutant and wild-type samples from the cystic fibrosis gene run on the DCode system. The mutant samples from exon 7 (Figure 1A) in lanes 2 and 3 resolve into two heteroduplex bands (upper bands) and a homoduplex band (lower band). The two heteroduplex bands in the mutant samples from exon 10 (Figure 1B) in lanes 2, 3, and 5 are resolved from the homoduplex bands. The $\Delta F508$ homozygous mutant in lane 4 has no heteroduplex bands, and the homoduplex bands migrate the same as the wild-type. Under these conditions, it was not possible to resolve the two homoduplex bands of the $\Delta F508$ mutant. This mutation may be detected if heteroduplexes are formed by mixing the mutant and wild-type samples together, then denaturing and renaturing the DNA.

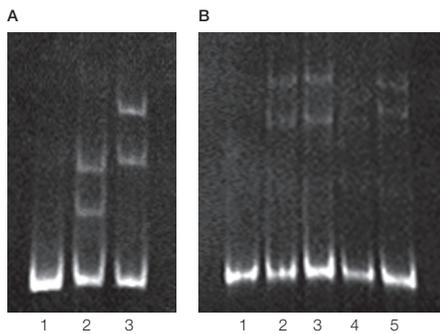


Fig. 1. Cystic fibrosis samples run on 1x DEM gels on the DCode system.
A, Exon 7 samples. Lane 1, wild-type DNA; lane 2, mutant sample 1154insTC; lane 3, mutant sample $\Delta F311$. **B,** Exon 10 samples. Lane 1, wild-type DNA; lane 2, mutant sample $\Delta F508$; lane 3, compound mutant sample Q493/ $\Delta F508$; lane 4, mutant sample $\Delta F508$; lane 5, mutant sample $\Delta I507$.

The ability to resolve the heteroduplex bands in the mutant samples makes it possible to distinguish between the mutant and wild-type samples. Heteroduplex analysis can be used as a method for screening heterozygous mutations in the cystic fibrosis gene using the DCode system.

References

- Nagamine CM et al., A PCR artifact: generation of heteroduplexes, *Hum Genet* 45, 337–339 (1989)
- White MB et al., Detecting single base substitutions as heteroduplex polymorphisms, *Genomics* 12, 301–306 (1992)

Practice of the polymerase chain reaction (PCR) may require a license.

Information in this tech note was current as of the date of writing (1999) and not necessarily the date this version (rev C, 2007) was published.



Bio-Rad
Laboratories, Inc.

Life Science
Group

Web site www.bio-rad.com USA 800 4BIORAD Australia 61 02 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 21 3237 9400
Canada 905 364 3435 China 86 21 6426 0808 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65
Germany 089 318 84 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 455 8800 India 91 124 4029300 Israel 03 963 6050
Italy 39 02 216091 Japan 03 6361 7000 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 0318 540666 New Zealand 0508 805 500
Norway 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700 Russia 7 495 721 14 04 Singapore 65 6415 3188 South Africa 27 861 246 723
Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 061 717 95 55 Taiwan 886 2 2578 7189 United Kingdom 020 8328 2000