

## Detection of *p53* Gene in Breast Cancer by Denaturing Gradient Gel Electrophoresis and the DCode™ System

Akira Ichikawa and Seiji Igarashi, Tochigi Cancer Center, Tochigi, Japan

### Introduction

Denaturing gradient gel electrophoresis (DGGE) is one of the most consistently used mutation scanning methods. It has evolved enormously over the last decade to be widely used, particularly in gene diagnostic laboratories. A practical method on this topic has been reported recently (Fodde and Losekoot 1994). The more frequently used mutation screening methods are DGGE and single-strand conformational polymorphism (SSCP). However, the sensitivity of DGGE is far higher (~95%) than that of SSCP, making it an attractive technique for screening unknown mutations.

The principle of DGGE is based on the fact that DNA duplexes melt at high temperature in the presence of a gradient of chemical denaturant. DNA melting is sequence specific and occurs in discrete segments called melting domains. Since melting involves breaking the hydrogen bonds that hold the base pair (bp) together, a G-C rich region melts at a higher temperature than an A-T rich region.

A polymerase chain reaction (PCR) product begins to melt from the region which has the lowest melting point (lowest  $T_m$ ) first when it is exposed to chemical denaturant and high temperature. If there are point mutations in the PCR product, the melting point will change, and thus its mobility on a polyacrylamide gel will be different from the mobility of the wild-type DNA. In other words, the mobility of the DNA fragment-containing mutation in the gel will be different from that of the wild type. In this paper we describe the use of DGGE to screen *p53* mutations in breast cancer.

### Methods

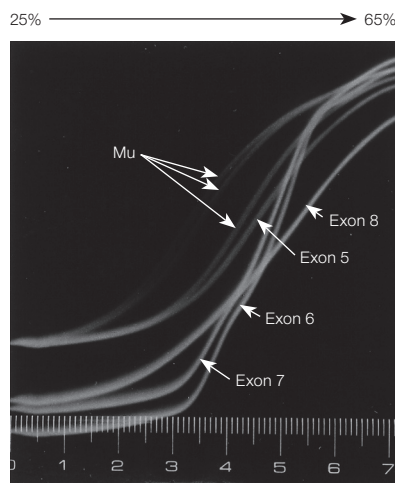
One hundred and two genomic DNA samples obtained from breast cancers were examined from exon 5, 6, 7, and 8 of the *p53* gene. They were amplified by PCR with a pair of primers which contains a 40 bp GC clamp in one of them (Andersen and Borresen 1995). All products were analyzed both in perpendicular gradient gels and parallel gradient gels.

### Perpendicular Gradient Gel Analysis

A 7.5 x 10 cm, 1 mm thick 25–65% denatured gradient gel was made using 6% acylamide/bis (37.5:1) in 1x TAE buffer (50 mM Tris, 25 mM acetic acid, 1.25 mM EDTA). Four different PCR products each from exon 5, 6, 7, and 8 were mixed in a 25  $\mu$ l volume, and the total sample volume made up to 100  $\mu$ l. We added 100  $\mu$ l 2x gel loading dye (70% glycerol, 0.05% Bromophenol Blue, 0.05% Xylene Cyanole, 2 mM EDTA) to these samples, and electrophoresed them on the DCode universal mutation detection system at 130 V for 2.5 hr at 56°C. After electrophoresis, the gels were stained in a 1:10,000 dilution of SYBR Green I (Molecular Probes, Inc.) in 1x TAE buffer for 45 min and destained in 1x TAE buffer for 45 min. The gels were imaged under ultraviolet (UV) transillumination.

### Parallel Gradient Gel Analysis

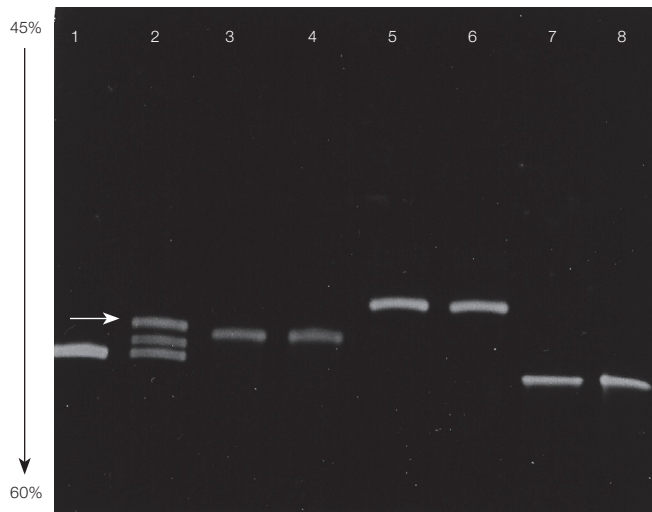
A 16 x 16 cm, 1 mm thick 45–60% denatured gradient gel was made using 6% acylamide/bis (37.5:1) in 1x TAE buffer. Five  $\mu$ l of PCR products (~200–300 ng) was mixed with 5  $\mu$ l 2x gel loading dye and electrophoresed on the DCode system at 150 V for 3.5 hr at 60°C. The post-run analysis was the same as perpendicular gradient gel analysis.



**Fig. 1. The typical case of perpendicular analysis using the DCode system.** Each wild-type exon is shown by an arrow, and mu of exon 5 is a mutant DNA. At this point in our investigations, the mutant DNA can be detected, but exon cannot be specified.

## Results and Discussion

The optimum concentration of denaturant to be used for parallel gradient gels was determined from the perpendicular gradient gel analysis. All exons were separated and had a melting transition using a 25–65% gradient (Figure 1). Therefore a gradient of 45–60% was used for parallel gel analysis (Figure 2).



**Fig. 2. Parallel gradient gel analysis using the DCode system.** Lane 1, wild-type DNA of exon 5; lane 2, sample DNA of exon 5; lane 3, wild-type DNA of exon 6; lane 4, sample DNA of exon 6; lane 5, wild-type DNA of exon 7; lane 6, sample DNA of exon 7; lane 7, wild-type DNA of exon 8; lane 8, sample DNA of exon 8. The arrow in lane 2 shows a mutant of exon 5, which is the same as the one showed in Figure 1 (mu).

DNA from 34 cases was found to have a mutation on perpendicular gels and in 33 cases on parallel gels. The sample missed showed a mutation of exon 7 on perpendicular gels but not on parallel gels. A possible explanation for this could be that the parallel gel gradient of 45–60% was too wide for that particular sample. A narrow gradient may be required to identify the mutation of exon 7 on parallel gels.

We chose to use parallel gradient method because it provides high-throughput analysis. The detection of mutation was unclear in some cases where the rate of mutated allele was low or the genomic DNA was extracted from paraffin-embedded materials. In such cases, we used SYBR Green I instead of ethidium bromide to stain the gels. This improved sensitivity and enabled us to see the bands clearly.

These experiments indicate that DGGE is a powerful technique to screen mutations. The DCode system is easy to use and reliable, two important factors when scanning numerous mutations.

## References

Andersen TI and Borresen AL, Alterations of the *TP53* gene as a potential prognostic marker in breast carcinomas. Advantages of using constant denaturant gel electrophoresis in mutation detection, *Diagn Mol Pathol* 4, 203–211 (1995)

Fodde R and Losekoot M, Mutation detection by denaturing gradient gel electrophoresis (DGGE), *Hum Mutat* 3, 83–94 (1994)

SYBR is a trademark of Molecular Probes, Inc.

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

Information in this tech note was current as of the date of writing (1998) and not necessarily the date this version (rev B, 2007) was published.

**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Web site [www.bio-rad.com](http://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