

Detection of Apolipoprotein Gene Variants by Denaturing Gradient Gel Electrophoresis Using the DCode™ System

Clive Pullinger, Arghavan Shahidi, Andrea Verhagen, and John Kane^{1,2}
 Cardiovascular Research Institute, ¹ Department of Medicine,
² Department of Biochemistry and Biophysics, University of California,
 San Francisco, California

Introduction

In a number of fields, techniques are required for searching candidate genes for mutations that underlie hereditary diseases. We have used the denaturing gradient gel electrophoresis method (DGGE) (Myers et al. 1987) using the Bio-Rad DCode universal mutation detection system to identify a number of new, as well as known, mutations and polymorphisms in the genes for apolipoprotein A-I (apoA-I) and apolipoprotein B (apoB).

Methods

Genomic DNA was prepared from patients with dyslipidemia. Segments of the apoA-I and apoB genes were amplified from genomic DNA using a temperature cycler (Ericomp, Inc.). Both oligonucleotide primers had GC clamps (Myers et al. 1985). Following restriction endonuclease digestion, the samples were subjected to DGGE.

The amplification reactions were performed in 50 mM Tris-HCl, pH 9 (at 25°C), 20 mM (NH₄)₂SO₄, 1.5 mM MgCl₂, 200 μM of each dNTP, 1 unit of Hot Tub polymerase (Amersham Life Science, Inc.), 15 pmol of each primer, and a minimum of 200 ng of DNA in a total volume of 50 μl. After initial denaturation at 96°C for 1 min, PCR was carried out for a total of 33 cycles at an annealing temperature of 62°C (apoA-I) or 58°C (apoB), in each case for 30 sec. The denaturing and elongation steps were 96°C for 30 sec and 72°C for 60 sec.

A final elongation step of 8 min at 72°C was then carried out. The temperature cycler was programmed to denature the amplification products at 96°C for 8 min and to then perform slow cooling to 40°C. Without this last step, heteroduplex DNA may not be visible on the DGGE gels.

In the case of apoA-I, a 420 bp region at the 5' end of the gene, containing the core promoter and the first exon, was amplified using two GC-clamped primers, such that the final PCR product was 532 bp in length. Similarly, 465 bp of the low-density lipoprotein (LDL) receptor binding domain of apoB was amplified with a final size of 591 bp. In each case, 10 μl of the PCR product was digested in a volume of 30 μl with either Avall (apoA-I) or EcoRI (apoB). Figure 1 shows the melting profile graphs for the 5' and 3' fragments generated after digestion of the apoB PCR product (Lerman and Silverstein 1987). These graphs were produced using Bio-Rad MacMelt™ software and were used to predict the percentage range of the denaturing gradient to be used in the gels. Multiple aliquots of the digests were loaded onto 7.5% acrylamide (38.5:1) DGGE gels at hourly intervals. These travel gels were run in the DCode system apparatus at a constant temperature of 56°C and 200 V for a total of 5 hr to determine the optimal running time. This was found to be 3 hr for the apoA-I fragments and 4 hr for apoB. The gels were stained for 30 min in SYBR Green and photographed under epi-illumination at 254 nm (Karlsen et al. 1995). A 35–65% denaturing gradient was used for apoA-I and a 30–50% gradient for apoB.

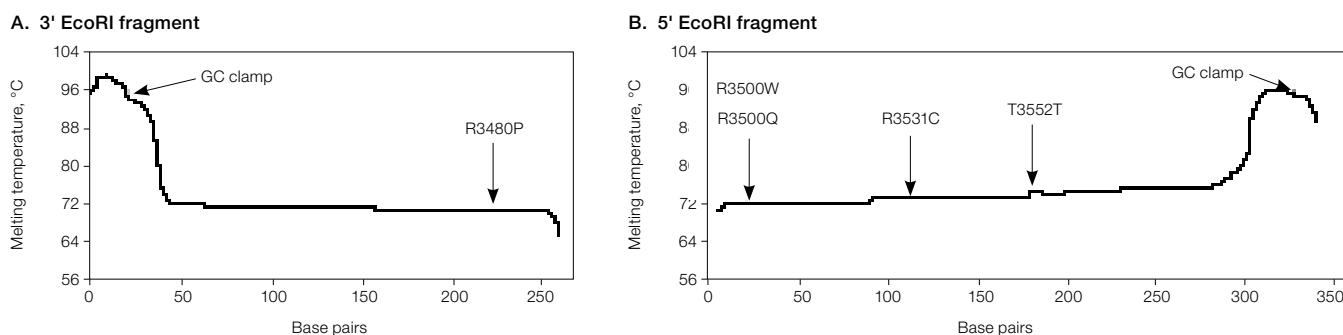


Fig. 1. Melting profiles of a portion of the receptor binding domain of the apoB gene amplified using GC-clamped primers. The graphs were generated using MacMelt software. A total of five single-base substitution mutations are shown in A and B.

Results

While no new mutations were found in the apoA-I promoter region, we were able to detect, using DGGE, a common polymorphism at nucleotide -75 (Figure 2), which had been previously shown to be associated with differences in plasma high-density lipoprotein (HDL) levels.

Three mutations in the LDL receptor binding domain of the apoB gene that cause the disorder familial ligand-defective apoB (FDB) (Soria et al. 1989, Pullinger et al. 1995, Gaffney et al. 1995) were detected by DGGE (Figure 3, lanes 3, 4, and 5). In addition, a rare missense mutation at codon 3480 (Figure 3, lane 1) and a new silent mutation (Figure 3, lane 2) were detected.

Discussion

Detection of naturally mutant alleles of candidate genes and a study of their consequent effects can be expected to provide important clues to the precise biochemical roles of these gene products. For example, new causes of atherogenic dyslipidemia will be revealed in the search for mutations in proteins involved in lipid metabolism. We have shown that the DGGE approach is an excellent method to detect these mutations.

Acknowledgements

This work was supported by National Institutes of Health Grants HL14237 (Arteriosclerosis Specialized Center of Research), HL50782, and HL50779, by a gift from Donald and Susan Schleicher, and a grant from the Joseph Drown Foundation.

References

- Gaffney D et al., Independent mutations at codon 3500 of the apolipoprotein B gene are associated with hyperlipidemia, *Arterioscler Thromb Vasc Biol* 15, 1025-1029 (1995)
- Karlisen F et al, SYBR green I DNA staining increases the detection sensitivity of viruses by polymerase chain reaction, *J Virol Methods* 55, 153-156 (1995)
- Lerman LS and Silverstein K, Computational simulation of DNA melting and its application to denaturing gradient gel electrophoresis, *Methods Enzymol* 155, 482-501 (1987)
- Myers RM et al., Nearly all single base substitutions in DNA fragments joined to a GC-clamp can be detected by denaturing gradient gel electrophoresis, *Nucleic Acids Res* 13, 3131-3145 (1985)
- Myers RM et al., Detection and localization of single base changes by denaturing gradient gel electrophoresis, *Methods Enzymol* 155, 501-527 (1987)
- Pullinger CR et al., Familial ligand-defective apolipoprotein B. Identification of a new mutation that decreases LDL receptor binding affinity, *J Clin Invest* 95, 1225-1234 (1995)
- Soria LF et al., Association between a specific apolipoprotein B mutation and familial defective apolipoprotein B-100, *Proc Natl Acad Sci USA* 86, 587-591 (1989)

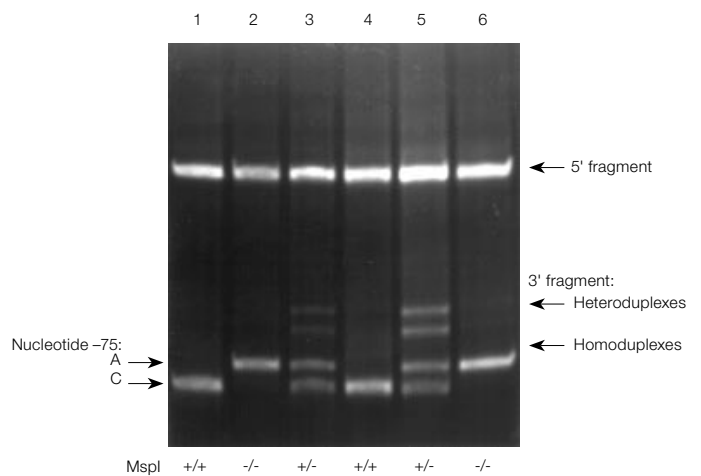


Fig. 2. DGGE gel of the apoA-I gene promoter showing an MspI polymorphism. Lanes 1 and 4 are homozygous for a C residue, and lanes 2 and 6 are homozygous for an A residue, at position -75. Lanes 3 and 5 are heterozygous subjects with the low-melting heteroduplexes clearly visible.

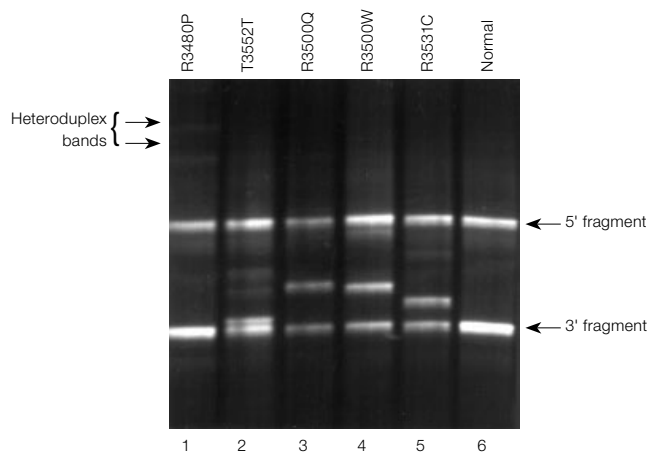


Fig. 3. DGGE gel of part of the receptor binding domain of the apoB gene showing five separate single-base substitutions. Four of the mutations, lanes 2 through 5, are on the 5' fragment (following EcoRI digestion) and, in each case, this band shows splitting. The fifth mutation (R3480P, lane 1) is on the 3' fragment and is only detectable because of the presence of heteroduplex bands.

Hot Tub is a trademark of GE Healthcare group companies.

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

Information in this tech note was current as of the date of writing (1997) and not necessarily the date this version (rev B, 2008) was published.

BIO-RAD

**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