

## Detection of Variation in Highly Polymorphic MHC Genes by Denaturing Gradient Gel Electrophoresis Using the DCode™ System

Kristi Miller, Tobi Ming, and Ruth Withler, Pacific Biological Station, Department of Fisheries and Oceans, Nanaimo, BC V9R 5K6, Canada

### Introduction

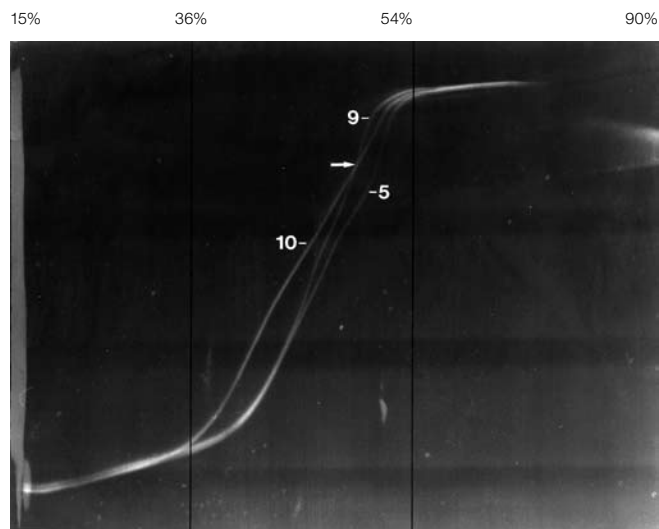
We describe a rapid and sensitive method for the detection of nucleotide sequence variation, which can be used for large-scale screening of population markers. Denaturing gradient gel electrophoresis (DGGE) detects sequence variants of amplified fragments by the differences in their melting behavior (Fischer and Lerman 1983). By adding a GC clamp to one of the primers, nearly all single-base substitutions can be detected (Sheffield et al. 1989). Although DGGE has primarily been utilized for the detection of single-base mutations in disease studies, it offers great potential for use in population analysis of genetic markers with higher levels of sequence variation. Herein, we describe our DGGE analysis of an MHC class I locus in salmon.

The methodology described was developed to identify the number and distribution of MHC class I A1 (exon 2) alleles among chinook salmon (*Oncorhynchus tshawytscha*) populations (Miller and Withler 1997, 1998; Miller et al. 1997). DGGE detects 25 of the 28 identified A1 sequences, with a nucleotide sequence variation between them of 1–16 nt plus a 2 codon indel. By creating a unique network of control alleles, 22 of these alleles can be rapidly and accurately resolved by a single-gel run condition. This technique has been used in surveys of over 20,000 salmon, representing two MHC markers (class I A1 and A2) in two salmon species (chinook and coho salmon). A single person in our laboratory now analyzes 160 salmon per day with DGGE, using two DCode universal mutation detection systems.

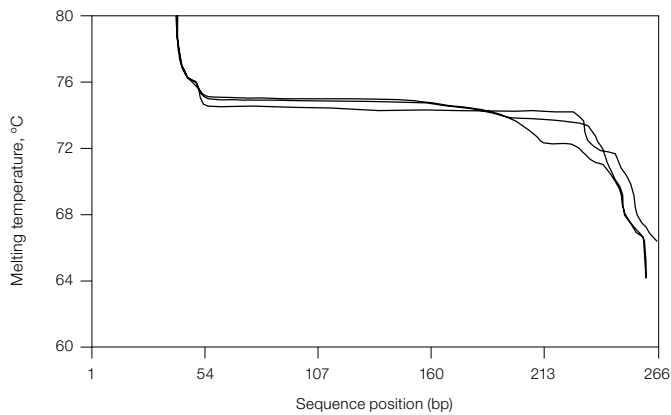
### Methods

Single locus MHC class I A1 primers were used to amplify fragments of 301–307 bp (including primers) (Miller et al. 1997). Because a large proportion of the sequence variation and the lower melting domain reside on the 3' end of the fragment, a GC clamp was added to the 5' primer.

DGGE was performed on the DCode system. We determined initial DGGE conditions by electrophoresing three of the amplified chinook salmon A1 chinook alleles (5, 9, and 10), differing from one another by 7–12 nt and a 2 codon indel on a perpendicular gradient gel (Figure 1). Dissociation of A1 alleles occurred over a 36–54% gradient. The presence of three distinctive but overlapping curves indicated that all three A1 alleles melted at different concentrations of denaturant and could be differentiated on parallel DGGE gels, but might shift relative positions if run into different portions of the denaturant gradient. Further analysis of these alleles by WinMelt™ software again showed three distinctive curves, with two melting domains still apparent (despite the GC clamp), and elucidated the crossover point between alleles 9 and 10 at nt position 207 (Figure 2). This crossover point directly coincided with the location of the 2 codon insert, present in allele 10, but absent in allele 9. Thus, it appeared that the 2 codon insert and the presence of two melting domains may cause shifting of some alleles on parallel gels.



**Fig. 1. Perpendicular DGGE of A1 alleles 5, 9, and 10 run on a chemical gradient of 15–90%. Melting curves for A1 alleles ranged from 36–54%. Curves for alleles 9 and 10 cross at 51.5% denaturant.**



**Fig. 2. Predicted melting curves for alleles 5, 9, and 10 calculated using WinMelt software.** Alleles differ by 7 to 12 nt and a 2 codon indel.

Time-series analyses on 40–65% (7.5% acrylamide) parallel gradient gels were used to determine the run time that separated most of the 28 A1 alleles (homoduplexes only) and to assess changes in the relative positions of alleles at different denaturant concentrations. Virtually all of the alleles were differentiated by DGGE, with the exception of two sets of alleles with a single C-G transversion within 10 bp of the clamped primer and one set of alleles with a single A-T transversion at nt 237. The remaining 5 sets of alleles with single-base differences at positions ranging from nt 85 to 282 (24 to 243, not including 5'-clamped primer) were differentiated, as well as all alleles with a greater number of differences. Thus, the range of denaturant gradient used and the addition of the GC clamp facilitated sampling of sequence variation over virtually the entire A1 fragment.

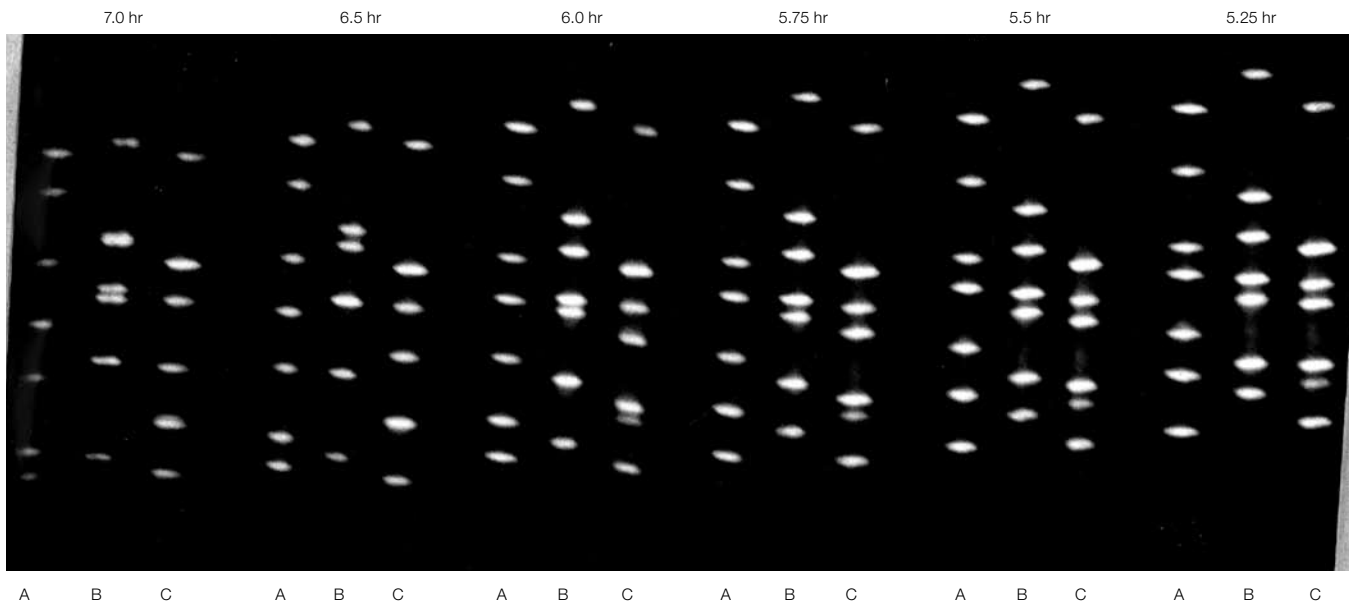
As expected from the presence of two melting domains within the A1 sequence, time-series analysis revealed a number of alleles which shifted relative positions on the gel when electrophoresed into different ranges of denaturant concentration. Moreover, some alleles that migrated differentially through the gel at low concentrations of denaturant became indistinguishable when electrophoresed in high denaturant concentrations (Figure 3). As a result, there was no single run time that consistently separated all of the alleles. In the time series shown in Figure 3, the best separation of alleles was obtained at a run time of 5.5 hr, which placed the DNA fragments between 48% and 53% denaturant, a range in which dissociation of the A1 low-melting domain was completed and the high-melting domain was initiated. At this run time, 18 alleles were distinctive, and 5 sets of alleles migrated to similar positions and were consequently binned into 5 allelic groups.

For the homoduplex A1 allele fragments, increasing migration distance through the denaturing gradient was strongly correlated with increasing overall GC content. However, the presence/absence of the 2-codon insert, and the GC content of the last 70 nt (distal to the GC clamp) also had apparent influence on allelic migration. Prior to any allele dissociation, alleles separated according to size, with the alleles containing the two-codon insert migrating slower than those without the insert. However, once alleles began to dissociate, the insert was apparently less influential on allele migration than was the GC content, and some of the alleles that contained the insert and a high overall GC content migrated past alleles that did not contain the insert. In addition, because much of the gradient sampled variation in the low-melt domain, GC content distal to the GC clamp was more influential on the overall migration distances of the homoduplexes than GC content close to the clamp.

Alleles that shifted (exchanged relative positions in a ladder of migration distance) often did not share the presence/absence of the 2-codon insert and/or differed in their relative GC content in the low- and the high-melt domains. The most notoriously shifting allele (9; fourth band from the bottom of lane A in Figure 3) shifts position with four other alleles so readily it can actually shift across a single gel if the denaturing gradient is not completely even. Careful monitoring of the standard allele sets is necessary to accurately score this allele, and it is often used as a marker for gel-to-gel run uniformity.

To facilitate rapid and accurate scoring of alleles on population gels, including those that shift relative positions during different run times, subsets of the “known” alleles were combined into three standard sets to use as controls in the following way: The relative positions of alleles run under optimal conditions were determined from the time-series analysis, and the fastest (latest melting) to slowest (earliest melting) alleles were alternately placed into each one of the 3 standard sets. Alleles that were known to shift relative positions at different denaturing concentrations (as identified in the time-series analysis) were placed in different standard sets. The A, B, and C lanes in the time series are the actual standard sets used in the population gels.

On the population gels, allele standard sets were loaded in three adjacent lanes in the center and on each side of the gel, with 20 individuals from one or two populations run in the remaining lanes. Parallel DGGE gels were photographed with a charge-coupled device (CCD) camera utilizing a 1024 x 1024 cm pixel array format and analyzed with standard gel documentation software. Standard allele sets were used to confirm that the electrophoretic conditions were sufficient to resolve all of the standard alleles, to monitor the relative positions of shifting alleles on the gel, and to form a network



**Fig. 3. Time-series analysis of chinook A1 alleles run on a 45–60% denaturant gradient for 5–7 hours at 120 V.** Lanes A, B, and C contain subsets of 21 of the 22 alleles distinguished by DGGE (allele 24 is not shown). Only a single allele from comigrating pairs of alleles is shown. The maximum resolution of alleles occurred at 5.5 hours, which was the run time used for further population analyses.

on which to score individual alleles. Alleles within the standard lanes were assigned “DGGE allele” numbers ranging from 50 to 400, approximately corresponding with their (logarithmic) positions on the gel. New alleles identified in the population survey with denaturing points that did not coincide with any of the known DGGE alleles were added to standard lanes as they were found.

Several factors affect our ability to produce DGGE gels of sufficient quality to score reliably over time. The purity and age of the acrylamide and acrylamide denaturant solutions, the maximum voltage, and the overall run time all strongly affect the resolution and migration of fragments. Optimally, for our analysis, DGGE was the most reliable when run slowly (60–80 V) overnight (<17 hours). However, to maximize the throughput of the DCode system, we conduct the A1 analysis on daytime runs (at 120 V max) and analyze other loci on overnight runs. We use a single brand of acrylamide for analysis of each gene locus and run a time series on new batches of acrylamide for slight run time adjustments. Acrylamide over 6 months old and acrylamide denaturant solutions over 7 days old can be problematic and are avoided. In general, run times for A1 are decreased by 5 minutes per day after the acrylamide solution is made. Furthermore, in order to accurately reproduce the run-time conditions from the time series to the population gels, population gels were prerun for 1.5 hr prior to loading. A more refined time series (from 5 to 6 hr) would decrease the necessary prerun time.

### Results and Discussion

DGGE has been utilized for population analysis in only a limited number of studies (Brown et al. 1997; Norman et al. 1994), most merely to identify haplotypes among a few individuals for sequencing. Our study offers the first example of the potential use of DGGE for large scale population surveys of highly polymorphic DNA loci. In our laboratory, each technician runs 160 fish per day on DGGE (8 gels) using two DCode systems (80 fish each for daytime (A1 locus) and nighttime (A2 locus) runs). Ambiguous scoring of shifting alleles necessitates the rerunning of less than 2% of all fish analyzed. This rapid rate of analysis of nucleotide sequence variation makes it possible to quickly accumulate large databases and to use DGGE for applied, real-time fishery issues, such as stock identification in mixed stock fisheries, escapement enumeration within river systems, and the identification of escaped farmed fish, to name a few.

## Conclusions

The DGGE methodology described herein can be used on any region of DNA with nucleotide sequence variation. Best results are obtained with fragments of 300–500 bp, but larger fragments can be analyzed by first digesting them with restriction endonucleases (Sheffield et al. 1989). In DGGE, the detection of length variation is often superseded by nucleotide variation, but DGGE could be an invaluable tool for detecting homoplasmy in microsatellites with similar migration on acrylamide gels. The use of DGGE virtually eliminates the need to sequence large numbers of clones to examine sequence variability. Direct sequencing of fragments cut out of gradient gels can further expedite the process. Finally, by careful use of standard networks of known alleles, DGGE is a highly sensitive and expeditious technique for detecting nucleotide sequence variation in large-scale population surveys.

## References

- Brown JM et al., Phylogeography and host association in a pollinating seed parasite *Greya politella* (Lepidoptera: Prodoxidae), *Mol Ecol*, 6, 215–224 (1997)
- Fischer SG and Lerman LS, DNA fragments differing by single base-pair substitutions are separated in denaturing gradient gels: correspondence with melting theory, *Proc Natl Acad Sci USA* 80, 1579–1583 (1983)
- Miller KM and Withler RE, MHC diversity in Pacific salmon: population structure and trans-species allelism, *Hereditas*, 127, 83–95 (1997)
- Miller KM and Withler RE, The salmonid class I MHC: limited diversity in a primitive teleost, *Immunological Rev.*, 166, 279–293 (1998)
- Miller KM et al., T. D., Molecular evolution at MHC genes in two populations of chinook salmon *Oncorhynchus tshawytscha*, *Mol Ecol*, 6, 937–954 (1997)
- Norman JA et al., C. J., Mitochondrial DNA control region polymorphisms: genetic markers for ecological studies of marine turtles, *Mol. Ecol.*, 3, 363–373 (1994)
- Sheffield VC et al., Attachment of a 40-base-pair G + C-rich sequence (GC-clamp) to genomic DNA fragments by the polymerase chain reaction results in improved detection of single-base changes, *Proc. Natl. Acad. Sci. USA*, 86, 232–236 (1989)

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, 2008) 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