

## Denaturing Gradient Gel Electrophoresis of Natural Virus Samples on the DCode™ System

Steven Short and Curtis Suttle, Department of Botany, The University of British Columbia, Vancouver, British Columbia

### Introduction

Denaturing gradient gel electrophoresis (DGGE) is a method of separating DNA fragments of the same size that differ in sequence. In DGGE, polyacrylamide gels are cast with linear gradients of denaturants. A 100% denaturing gel solution is defined as having 40% (v/v) formamide and 7 M urea. DNA fragments of the same size that differ in sequence will not resolve in agarose gels, but will migrate to different positions in a DGGE gel as the DNA denatures.

DGGE makes use of the fact that the melting behavior of double-stranded DNA (dsDNA) is affected by temperature, concentration of denaturant (e.g., urea and/or formamide), and interactions of adjacent base pairs stacked in a helix (Myers et al. 1987). Therefore, discrete segments or domains within a dsDNA molecule denature under different conditions, according to their base composition. As a result of the formation of branched DNA molecules, partial denaturation reduces electrophoretic mobility. Resolution of dsDNA samples of the same length, but differing in sequence, occurs by running the DNA in a polyacrylamide gel that has a continuously increasing concentration of denaturant. Because DNA strands that differ in sequence partially melt at different concentrations of denaturant, their migration slows at different positions in the gel.

In this experiment, we show that parallel DGGE analysis on the DCode universal mutation detection system can be used to analyze sequence differences among virus genes amplified from related viruses and from natural virus communities. Furthermore, we found that a GC clamp on polymerase chain reaction (PCR) primers used for amplification was not required for DGGE with these products. This approach can be used to genetically fingerprint natural virus communities.

### Methods

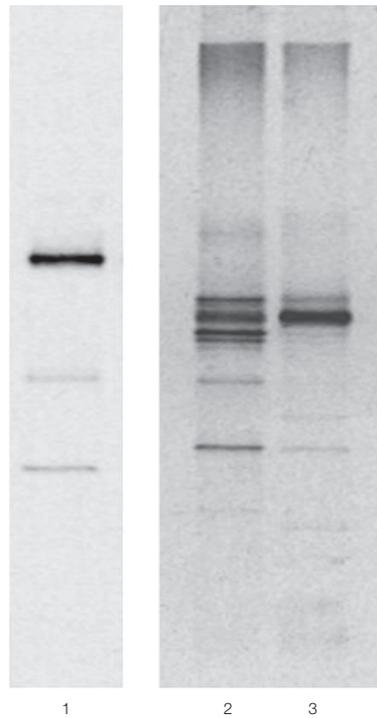
DNA polymerase gene fragments were amplified from a combined template consisting of DNA from cultured viruses (MpV-SP1, MpV-PB8, and CVA-1) belonging to the family Phycodnaviridae. Template DNA was also obtained from natural marine virus samples collected from two inlets on the west coast of Vancouver Island, Canada. Target sequences from all templates were amplified by PCR using degenerate primers that are specific for members of the virus family Phycodnaviridae (Chen and Suttle 1995). We did not need GC clamps to amplify these longer PCR products. Reactions conducted on the combined template and the two natural virus samples all produced an end product of approximately 700 base pairs.

DGGE was conducted using a 6% acrylamide/bis (37.5:1) gel with a parallel denaturing gradient range of 0–60% in 1x TAE (40 mM Tris, 20 mM acetic acid, 1 mM EDTA). The gradient gel was cast using the Bio-Rad Model 475 gradient delivery system. Approximately 300 ng of the cultured virus product was mixed with 12 µl of 2x loading dye (70% glycerol, 0.05% Bromophenol Blue, 0.05% Xylene Cyanol, 2 mM EDTA), while 1 µg of PCR products from the natural communities was combined with 20 µl of 2x loading dye. Electrophoresis was conducted using the DCode system at 135 V for 6 hr at a buffer temperature of 58°C. After electrophoresis, the gel was stained in a 1:10,000 dilution of SYBR Green I (Molecular Probes, Inc.) in 1x TAE buffer overnight. The gel was imaged under ultraviolet (UV) transillumination.

## Results and Discussion

Figure 1 shows virus DNA polymerase gene fragments run on a parallel DGGE gel. Lane 1 reveals that PCR amplification from combined templates produced three fragments that correspond to the individual viruses. It has been previously demonstrated that when amplified individually, each of the three viruses produces a unique DGGE band (Short and Suttle 2000). Lanes 2 and 3 reveal unique banding patterns for each virus sample, suggesting that virus community composition is spatially variable. In a related study, DGGE of natural samples revealed both spatial and temporal differences in the genetic composition of natural virus samples (Short and Suttle 1999). Notably, when parallel DGGE-separated mixtures of similar PCR products (Figure 1) were run on a standard agarose gel, they migrated as a single band. Thus, the ability of parallel DGGE to resolve similar-sized fragments produced in a single PCR reaction makes it possible to detect unique sequences in samples of varying genetic diversity. Our data indicate that DGGE can be used to distinguish DNA polymerase gene fragments of the same size from related yet genetically distinct viruses. Furthermore, DGGE distinguishes natural virus samples obtained from different marine locations.

DGGE is a reliable technique to screen the genetic diversity of aquatic virus samples using the DCode system. Reproducible gradients are easily cast with the Bio-Rad gradient delivery system, and buffer temperature can be held constant by the DCode system, ensuring reproducible electrophoresis conditions.



**Fig. 1. Parallel DGGE separation of virus DNA polymerase gene fragments run on the DCode system.** Lane 1, PCR products amplified from the combined templates of MpV-SP1, MpV-PB8, and CVA-1 DNA; lane 2, PCR products amplified from a virus sample from Grappler Inlet, British Columbia, Canada; lane 3, PCR products amplified from a virus sample from Bamfield Inlet, British Columbia, Canada.

## References

- Chen F and Suttle CA, Amplification of DNA polymerase gene fragments from viruses infecting microalgae, *Appl Environ Microbiol* 61, 1274–1278 (1995)
- Myers RM et al., Detection and localization of single base changes by denaturing gradient gel electrophoresis, *Methods Enzymol* 155, 501–527 (1987)
- Short SM and Suttle CA, Denaturing gradient gel electrophoresis resolves virus sequences amplified with degenerate primers, *Biotechniques* 28, 20–26 (2000)
- Short SM and Suttle CA, Use of the polymerase chain reaction and denaturing gradient gel electrophoresis to study diversity in natural virus communities, *Hydrobiologia* 401, 19–33 (1999)

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 (1999) and not necessarily the date this version (rev B, 2006) 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