

## DGGE Analysis for Noninvasive Studies of Primate Diet: A Prototype for At-Distance Diet Analyses of Natural Populations

Jodi Irwin<sup>1</sup> and Cristián Orrego<sup>2</sup>, <sup>1</sup> Department of Cell and Molecular Biology, Northwestern University Medical School, Chicago, IL 60611-3008;

<sup>2</sup> Conservation Genetics Laboratory, Department of Biology, San Francisco State University, San Francisco, CA 94132

### Introduction

The applications of denaturing gradient gel electrophoresis (DGGE) to the investigation and diagnosis of human genetic diseases are well known (Fodde and Losekoot 1994). More recently, however, DGGE has also become a valuable tool for addressing questions pertaining to natural history. It has been applied to studies of population genetics (Lessa and Applebaum 1993) and phylogeography (Lessa 1993), and it has been particularly useful in analyses of microbial communities (Muyzer et al. 1993, Murray et al. 1996).

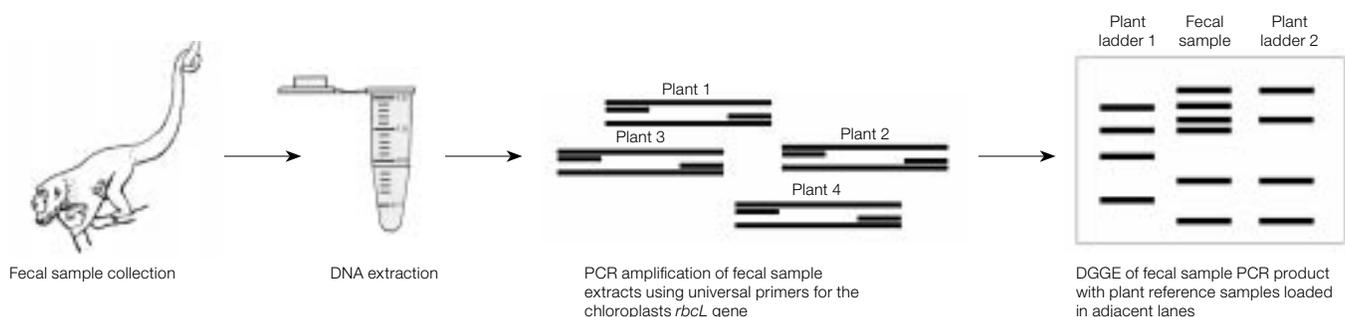
We demonstrate the use of DGGE as a molecular technique that can be employed in the absence of behavioral data for examining the diet of free-ranging howler monkeys (*Alouatta pigra*) (Figure 1). The difficulties associated with the collection of behavioral data for diet studies are well known. These include problems related to observing monkeys through thick foliage, the large distances separating observer from subject, and the difficulty of identifying tree species with very similar leaf morphologies. In order to circumvent these problems, we investigated the potential of a polymerase chain reaction (PCR)/DGGE method for augmenting knowledge of the howler diet. We hope that the development of a noninvasive technique for examining diet will facilitate studies of wild populations, particularly those for which behavioral data are not available.

### Methods

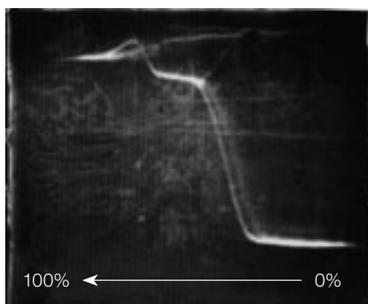
Genomic DNA was extracted from fecal samples and reference plants, and the chloroplast *rbcL* gene amplified via the PCR with one GC-clamped primer. PCR mixtures contained 1x PC2 PCR buffer (Ab Peptides), 0.1 units/ $\mu$ l KlenTaq DNA polymerase (Ab Peptides),  $MgCl_2$  (5.0 mM), 0.25  $\mu$ M each primer, and 1.5% dimethyl sulfoxide. Thermal cycling conditions were: an initial denaturation at 94°C for 4 min; 3 cycles of 94°C, 1 min; 62°C, 1 min; 80°C, 45 sec; 4 cycles of 94°C, 1 min; 64°C, 1 min; 80°C, 45 sec; 32 cycles of 94°C, 1 min; 68°C, 1 min; 80°C, 45 sec; and a 5 min final extension at 80°C. PCR products were subsequently subjected to DGGE, using the DCode™ universal mutation detection system from Bio-Rad, to resolve the various plant sequences produced by PCR amplification.

To determine the melting temperature of PCR products, they were first analyzed on a perpendicular denaturing gradient gel poured with Bio-Rad's Model 475 gradient former. The gel was 6.5% 37.5:1 acrylamide/bis-acrylamide, and contained a gradient of 1–100% denaturant. Approximately 100  $\mu$ l of PCR product were loaded on a gel, and electrophoresis performed at 130 V and 60°C for 3 hr (Figure 2).

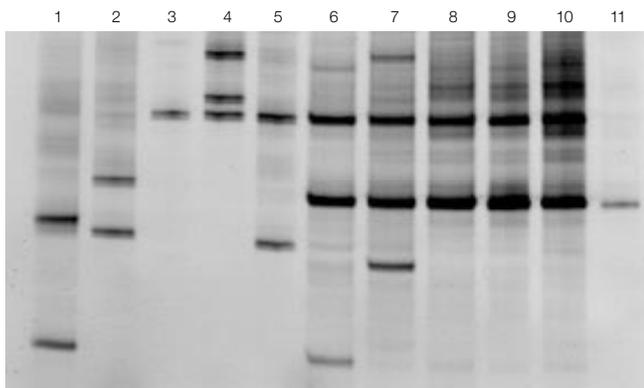
Optimal resolution of fragments on subsequent parallel gels was obtained with gradients ranging from 30–40% denaturant, with electrophoresis at 130 V and 60°C for approximately 4 hr (Figure 3). Gels were stained with ethidium bromide, and visualized by UV fluorescence.



**Fig. 1. Schematic outline of the molecular method used to examine the howler monkey diet.** The technique produces PCR products from different plants that are identical in size, but may differ in sequence composition.



**Fig. 2. Perpendicular denaturing gradient gel of fecal sample PCR product.** The two inflection points visible in the gel represent two melting domains in the fragment. The existence of two melting domains was also evident from analysis with MacMelt™ software (Bio-Rad). The two bands visible in the first domain represent at least two plant DNAs amplified in the PCR.

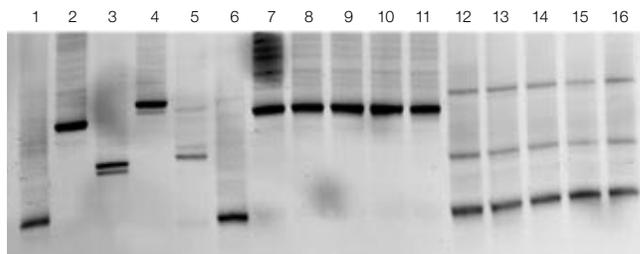


**Fig. 3. DGGE of fecal samples from captive animals and the plants they are known to have eaten.** The *rbcl* gene was amplified from individual plants, and the PCR products were combined as follows prior to loading: Lane 1, plants 10 and 11; lane 2, plants 12 and 13; lane 3, plant 14; lane 4, plants 15–17; lane 5, plants 18 and 19; lanes 6–10, correspond to PCR products from fecal samples 1–5 and from plant 20, respectively. The additional 7 plants eaten by the monkeys were not loaded on this gel, yet are distinguishable from each other and from plants 10–20. Note the identical migration of plants 14, 17, and 18.

## Results and Discussion

Our results indicate that DGGE is an effective technique for resolving unique sequences from mixed template PCR amplifications. Analyses of eighteen plant species consumed by captive animals showed that DGGE permitted resolution of 15 of the 18 species, or over 80% (Figure 3). The three unresolved species possessed identical sequences (unpublished) in the melting domain under examination, and thus differentiation could not be expected.

Additional results from samples obtained from captive animals whose diets were known, indicate that, on average, only about 30% of all plants ingested by the monkeys were detected by PCR/DGGE. Further experiments on constructed plant mixtures confirmed that a moderate amount of amplification bias occurs in PCR amplifications of mixed templates (Figure 4). Experiments to assess the preferential amplification observed in both the control fecal samples and the constructed plant mixtures suggest multiple factors at play, including template concentration and differential primer fidelity to various templates.



**Fig. 4. PCR amplification of the *rbcl* gene from individual plants and from plant mixtures.** Lanes 1–6, individual plants 1–6, respectively; lanes 7–11, replicates of mixture of plants 1–6; lanes 12–16, replicates of mixture of plants 1,3,4,5, and 6.

These data suggest that the PCR/DGGE approach tested in this investigation should be useful in situations where a universal detection system is not required, or where nothing is known about the diet of an animal. Further optimization of the technique should permit greater fragment resolution and enhanced sensitivity. Yet, any future investigations utilizing these methods should carefully consider the effects of PCR bias (Reysenbach et al. 1992, Ferris et al. 1996, Suzuki and Giovannoni 1996).

## References

- Ferris MJ et al., Denaturing gradient gel electrophoresis profiles of 16S rRNA-defined populations inhabiting a hot spring microbial mat community, *Appl Environ Microbiol* 62, 340–346 (1996)
- Fodde R and Losekoot M, Mutation detection by denaturing gradient gel electrophoresis (DGGE), *Hum Mutat* 3, 83–94 (1994)
- Lessa EP, Analysis of DNA sequence variation at population level by polymerase chain reaction and denaturing gradient gel electrophoresis, *Methods Enzymol* 224, 419–428 (1993)
- Lessa EP and Applebaum G, Screening techniques for detecting allelic variation in DNA sequences, *Mol Ecol* 2, 119–129 (1993)
- Murray AE, Hollibaugh JT and Orrego C, Phylogenetic compositions of bacterioplankton from two California estuaries compared by denaturing gradient gel electrophoresis of 16S rDNA fragments, *Appl Environ Microbiol* 62, 2676–2680 (1996)
- Muyzer G et al., Profiling of complex microbial populations by denaturing gradient gel electrophoresis analysis of polymerase chain reaction-amplified genes coding for 16S rRNA, *Appl Environ Microbiol* 59, 695–700 (1993)
- Reysenbach AL et al., Differential amplification of rRNA genes by polymerase chain reaction, *Appl Environ Microbiol* 58, 3417–3418 (1992)
- Suzuki MT and Giovannoni SJ, Bias caused by template annealing in the amplification of mixtures of 16S rRNA genes by PCR, *Appl Environ Microbiol* 62, 625–630 (1996)
- We are grateful to Hal Markowitz (San Francisco State University) for the opportunity to conduct this work. This project was supported by a U.S. Department of Education GAANN Predoctoral Fellowship, an SFSU mini-grant, and Oceanic Society Expeditions. This work was in partial fulfillment of a master's of science degree by Jodi Irwin.

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