

Microbial Diversity in Ground and Surface Water Analyzed by Denaturing Gradient Gel Electrophoresis Using the DCode™ System

Bernd Eschweiler and Beate Kilb, Institute for Water Research, Schwerte, Germany

Introduction

Bacterial diversity in environmental samples is usually determined by a characterization of isolated strains. A problem for the analysis and characterization of microbial communities is the inability to culture most of the bacteria species present in the sample. Therefore, isolated bacteria may account for only a minor portion of the total bacterial diversity originally present in the sample. This problem is particularly severe in oligotrophic habitats like ground water, where approximately only 0.1-1% of the bacterial species are culturable (Amann et al. 1995). A new approach in microbial ecology is based on the analysis of bacterial genetic information without cultivation. This cultureindependent approach has greatly enhanced the ability to assess bacterial diversity in ecosystems such as ground and surface water environments. After isolation of total bacterial DNA, variable regions of the 16S rRNA gene are amplified by the polymerase chain reaction (PCR). The similar-sized PCR products are separated by subsequent denaturing gradient gel electrophoresis (DGGE), and the resulting diversity pattern are analyzed and compared (Kuhlmann et al. 1997, Eschweiler et al. 1998).

Methods

DNA from surface and ground water was isolated and purified as described earlier. Two universal bacterial 16S rDNA primers were used to amplify a 527 bp fragment from total genomic DNA. PCR was performed in a total volume of 100 µl containing 1x PCR buffer (10 mM Tris-HCl, pH 8.3, 50 mM KCl, 1.5 mM MgCl₂, 0.001% gelatin), 200 µM each dNTP, 0.5 µM each primer, 0.3 mg/ml bovine serum albumin (BSA), and 2.5 U AmpliTaq DNA polymerase (PerkinElmer, Inc.), 10 µl DNA solution. The temperature cycle for the PCR was 60 sec of denaturation at 94°C, 60 sec of annealing (see below), and 90 sec of primer extension at 72°C. During an initial touchdown cycle, the annealing temperature was lowered from 65°C to 55°C in intervals of 1°C per cycle. The additional annealing cycles were done at 55°C. Ten PCR cycles were performed for the touchdown procedure and then 20 additional cycles at the actual annealing temperature of 55°C. This touchdown procedure reduces the formation of spurious by-products

during the amplification process (Don et al. 1991, Muyzer et al. 1993). DGGE was performed with the DCode universal mutation detection system (Bio-Rad Laboratories, Inc.). The polyacrylamide gels (7.5%) contained a denaturing gradient from 40% to 70% (100% denaturant: 7 M urea and 40% formamide) and were run in 0.5x TAE buffer (40 mM Tris base, pH 7.4, 20 mM sodium acetate, 1 mM EDTA). Gels were poured with a stacker on top (7.5% polyacrylamide, no denaturant). The optimum denaturant gradient was determined by performing perpendicular DGGE according to the DCode instruction manual. Electrophoresis was performed at constant voltage (70 V) and temperature (57°C) for 16 hr. After electrophoresis, the gels were stained with silver nitrate (Heukeshoven and Dernick 1985), dried, and photographed.

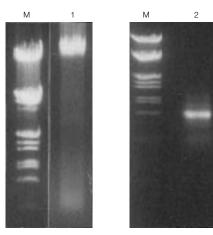


Fig. 1. Agarose gel electrophoresis from isolated and PCR-amplified DNA from surface water. Lane 1, isolated DNA from river Ruhr; lane 2, PCR-amplified DNA. M, size standard (λ DNA x HindIII x EcoRI).

Results and Discussion

Isolated and PCR-amplified DNA from surface water is shown in Figure 1. Separation of 16S rDNA amplified fragments by DGGE revealed the great diversity in ground and surface water samples. Complex banding patterns could be observed in samples with different hydrochemical conditions. These patterns were characteristic for each sample and showed the differences and common features in species composition.



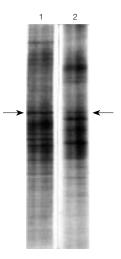


Fig. 2. DGGE patterns from ground water with different hydrochemical characteristics. Lane 1, anoxic ground water; lane 2, oxic ground water. Arrows indicate example of bands found in both samples.

The DGGE patterns from ground water with different hydrochemical characteristics are shown in Figure 2. Patterns from anoxic ground water are different from those obtained with oxic ground water. Some bands were found in both samples, indicating bacterial populations that could adapt to both habitats. In order to demonstrate temporal changes in microbial populations, DGGE patterns from surface water (river Ruhr) were recorded over a period of six months. Figure 3 shows changing patterns where certain species of bacteria are found over a limited period of time, while others are found in all samples.

The complex DGGE patterns are further analyzed and compared by digitized gel imaging using the Model GS-700 imaging densitometer and Multi-Analyst™ software (Bio-Rad). DGGE patterns from surface and ground water reveal a specific and unique bacterial population depending on the hydrochemical properties. In addition, DGGE patterns are used to demonstrate temporal and spatial variations in species composition. Sudden changes in DGGE patterns might indicate the introduction of harmful compounds (for example, pesticides and heavy metals) to surface and ground water. Analysis by DGGE is also suitable for a subsequent species identification by sequencing individual bands (Ferris et al. 1996).

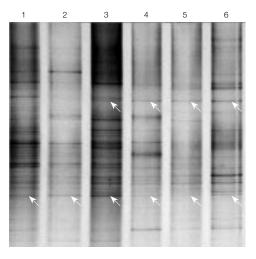


Fig. 3. DGGE patterns from surface water. Lanes 1–6, surface water (river Ruhr) from March through August, respectively. Arrows indicate species found over a limited period of time (top) and species found in all samples (bottom).

References

Amann RI et al., Phylogenetic identification and in situ detection of individual microbial cells without cultivation, Microbiol Rev 59, 143-169 (1995)

Don RH et al., 'Touchdown' PCR to circumvent spurious priming during gene amplification, Nucleic Acids Res 19, 4008 (1991)

Eschweiler B et al., pp 129-134 in Peters JH et al. (eds), Artificial recharge of groundwater: proceedings of the third International Symposium on Artificial Recharge of Groundwater: TISAR 98, AA Balkema, Rotterdam, Brookfield (1998)

Ferris MJ et al., Denaturing gradient gel electrophoresis profiles of 16S rRNAdefined populations inhabiting a hot spring microbial mat community, Appl Environ Microbiol 62, 340-346 (1996)

Heukeshoven J and Dernick R, Simplified method for silver staining of proteins in polyacrylamide gels and the mechanism of silver staining, Electrophoresis 6, 103-112 (1985)

Kuhlmann B et al., Molekularbiologische Methoden zur Bestimmung der mikrobiellen Diversität in Grundwassersytemen, Vom Wasser 89, 205-214 (1997)

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)

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 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

Bulletin 2366 Rev B US/EG 06-0373 0308 Sig 0308