

Genetic Monitoring by Denaturant Gel Electrophoresis of *Piscirickettsia salmonis*, a Bacterial Disease of Farmed Salmonids

Sekou Heath, Sunny Pak, and Cristián Orrego (cob@sfsu.edu), Conservation Genetics Laboratory, Department of Biology, San Francisco State University, 1600 Holloway Avenue, San Francisco, CA 94132, USA; Sergio Marshall, Instituto de Biología, Universidad Católica de Valparaíso, Valparaíso, Chile

Introduction

Piscirickettsia salmonis is the agent of an aggressive transmissible bacterial disease affecting salmonid aquaculture in Chile (Almendras and Fuentealba et al. 1997, Fryer et al. 1992, Larenas et al. 1997). Commercial rearing of coho salmon (*Oncorhynchus kisutch*), rainbow trout (*Oncorhynchus mykiss*), and Atlantic salmon (*Salmo salar*) has been compromised in the aquafarms of southern Chile (Chiloé Island, Region Ten) since the emergence of this pathogen about 10 years ago. Efforts in various laboratories toward control of the disease include development of a vaccine. The success of this approach will depend in part on the degree of biological and genetic diversity within the *P. salmonis* lineage. For example, the effectiveness of a vaccine might be diminished by unpredictable temporal shifts of *P. salmonis* types invading salmon farms.

We have employed denaturing gradient gel electrophoresis (DGGE) and a simpler format, constant denaturing gel electrophoresis (CDGE), to monitor genetic diversity in *P. salmonis* obtained from salmon farms in southern Chile. Amplification and sequencing of a section of the internal transcribed spacer (ITS) between the 16S and 23S rRNA genes indicate that all reported strains of *P. salmonis* can be grouped into four sequence types. In this report, we demonstrate their distinction by CDGE. Initial surveys with this method show the presence of only one sequence type of *P. salmonis* in Chilean salmonid farms. The recurrent type groups with strain LF-89, the type strain (ATCC VR-1361) first described in the early 1990s (Fryer et al. 1990, Fryer et al. 1992), a few years after initial reports of the emergence of the salmonid rickettsial syndrome in Chile.

Methods

DNA was extracted by the Chelex® method from fish serum and tissue or infected cell lines (Marshall et al. 1998). Polymerase chain reaction (PCR) amplification was conducted as previously described with primers RTS1 and RTS4 positioned just inside the borders of the ITS with the 16S and 23S rRNA genes, to obtain a fragment 283 bp in length (Marshall et al. 1998). Amplification for subsequent electrophoresis on denaturant gels was done with the same primer pair, but with a 40-bp GC clamp (Murray et al. 1996) added at the 5'-end of primer RTS1, in order to produce a fragment suitable for detection of single-base changes (Sheffield et al. 1989). Sequences from fragments amplified without a GC clamp were obtained as previously described (Marshall et al. 1998).

The melting temperatures of representative PCR products were obtained on a perpendicular denaturing gradient gel cast with a Bio-Rad Model 475 gradient former in 16 cm long plates, which are part of the DCode™ universal mutation detection system (Bio-Rad Laboratories, Inc.). The gel was 6.0% 37.5:1 acrylamide/bis with a 10–40% urea-formamide gradient (Bio-Rad's Model 475 gradient delivery system instruction manual). Approximately 100 µl of the PCR product was distributed evenly along the top of the gel, and electrophoresis was conducted at 165 V and 56°C for 3.5 hr. Resolution of sequences was then assessed with parallel gels with a gradient from 22–25% denaturant. The CDGE format set at a denaturant level of 23.5% proved equally useful and was adopted as the method of choice for surveying natural isolates. Gels were stained with ethidium bromide, and separation patterns were revealed by ultraviolet (UV) fluorescence, digitized with the Gel Doc™ 1000 system (Bio-Rad), and analyzed with PC software provided with the system.

Results and Discussion

The migration behavior of amplified fragments from strains LF-89 and EM-90 during perpendicular DGGE at 56°C revealed two melting transitions (Figure 1), the first at inflection point 24% and a second at inflection point 27% denaturant (average of the two curves; Figure 1). The theoretical melting curves for the sequences in question obtained with WinMelt™ program (Bio-Rad) also displayed two melting domains, one evident at 62°C and a second between 62–63.5°C (Figure 2). The theoretical curve roughly corresponds to the 3% denaturant differential in the observed inflection points, should one adopt the relationship of 0.3°C/1% denaturant (Abrams and Stanton et al. 1992).

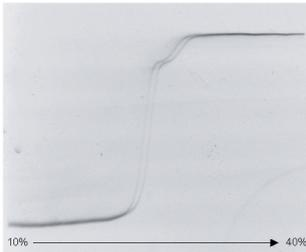


Fig. 1. Perpendicular denaturing gradient gel of amplified products obtained from strains LF-89 and EM-90. Melting theory predicts that LF-89 is the profile at left (see Figure 2).

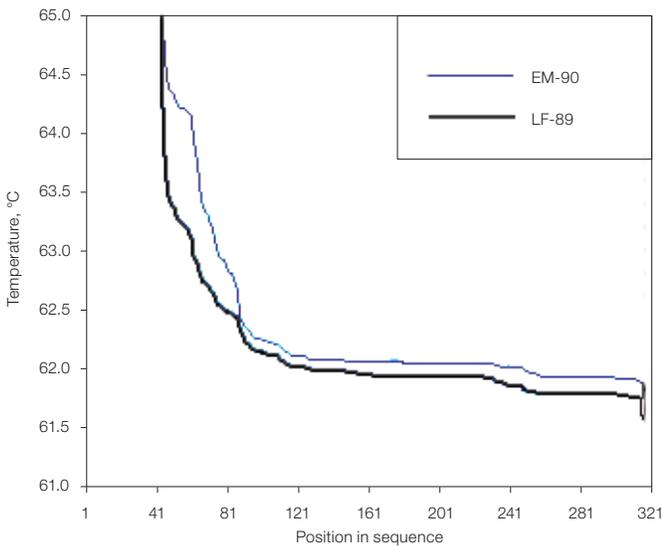


Fig. 2. Predicted melting profiles obtained with WinMelt software for strains LF-89 and EM-90 at 50% helicity with expanded y-axis. The x-axis covers the entire length of the amplified fragment, including the GC clamp at the left edge.

The predicted melting curves indicating higher melting values for EM-90 compared to LF-89 (Figure 2) were confirmed in the parallel gel format with a gradient ranging between 22–25% (results not shown) and also with the CDGE format at a set denaturant concentration of 23.5% (Figure 3). CDGE became the preferred format, given the reproducibility from run to run obtained by casting gels at a single urea-formamide concentration. CDGE revealed four migration classes: EM-90 as the sequence type having the highest melting value, and consequently the fastest migration; NOR-92; ATL-4-91; and the most slowly migrating class, including strains LF-89, C1-95, and SLGO-94.

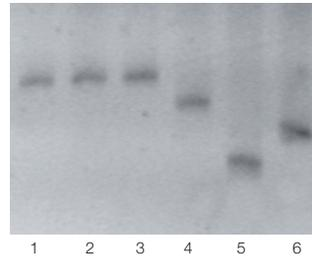


Fig. 3. CDGE of amplified fragments from *P. salmonis* strains. Lane 1, C1-95; lane 2, SLGO-94; lane 3, LF-89; lane 4, ATL-4-91; lane 5, EM-90; and lane 6, NOR-92.

Comparative analysis of sequences obtained from these strains confirmed the grouping of the fragments into four distinct melting classes on the denaturant gels. One sequence type was common to strains from farmed Chilean salmonids, including *P. salmonis* LF-89 (coho salmon), SLGO-94 (rainbow trout), and C1-95 (coho salmon), while strains EM-90 (isolated in Chile from Atlantic salmon), ATL-4-91 (Canada, Atlantic salmon), and NOR-92 (Norway, Atlantic salmon) each had unique sequences.

Surveys by CDGE of 44 samples (38 from coho salmon, 6 from rainbow trout) amplified from tissue biopsies obtained during 1997 and 1998 from five different salmon farms on Chiloé Island showed the presence only of *P. salmonis* that group with type strain LF-89. Earlier, Marshall et al. (1998) suggested, on the basis of DNA sequence data, that LF-89-like *P. salmonis* has been the prevalent type in the Chiloé region since the emergence of the disease during the 1980s.

The sequence data alluded to here differ at a few positions from those previously reported (GenBank accessions U36943-6 and U62103-4). This will be the subject of a future communication (Heath et al., unpublished). The congruence of melting behavior of the amplified fragments from the six *P. salmonis* strains and the corresponding DNA sequence data is consistent with our revisions.

The CDGE assay offers a rapid molecular approach to monitoring for bacterial types involved in the salmonid rickettsial syndrome. Conceivably, the appearance of a new *P. salmonis* type distinct from those already reported should be quickly evident from CDGE analysis. The CDGE protocol optimized here is sufficiently sensitive so that one base-pair substitution, as between NOR-92 and ATL-4-91 (Heath et al., unpublished), results in easily distinguishable differential mobility of the corresponding amplicons (Figure 3). The admittedly limited survey of salmon farms in Chile reported here is nevertheless reassuring in regard to the dependability of future vaccines against *P. salmonis*, given the observed constancy of type since the emergence of piscirickettsiosis about 10 years ago.

Acknowledgements

We are grateful to Ellen Prager (SFSU) for insightful editing, Vitalia Henríquez (UCV) for catalogued samples, Michael Mauel (University of Rhode Island) for purified DNA from *P. salmonis* strains, and Asmita Patel (Bio-Rad) for the WinMelt program. S.H., S.P., and C.O. were supported by Minority Biomedical Research Support grant GM52588 from the National Institutes of Health to C.O., and S.M. was supported by Fomento al Desarrollo Científico y Tecnológico (FONDEF)/Comisión Nacional de Investigación Científica y Tecnológica (CONICYT), project 1038. This work was in partial fulfillment of a Master's in science degree by S.H.

References

- Abrams ES and Stanton VP, Use of denaturing gradient gel electrophoresis to study conformational transitions in nucleic acids, *Methods Enzymol* 212, 71–104 (1992)
- Almendras FE and Fuentealba C, Salmonid rickettsial septicemia caused by *Piscirickettsia salmonis*: a review, *Diseases of Aquatic Organisms* 29, 137–144 (1997)
- Fryer JL et al., Isolation of a rickettsiales-like organism from diseased coho salmon (*Oncorhynchus kisutch*) in Chile, *Fish Pathol* 25, 107–114 (1990)
- Fryer JL et al., *Piscirickettsia salmonis* gen. nov., sp. nov., the causative agent of an epizootic disease in salmonid fishes, *Int J Syst Bacteriol* 42, 120–126 (1992)
- Larenas J et al., <http://aquatic.unizar.es/n1/art505/piscirick.html> (1997)
- Marshall S et al., Minimally invasive detection of *Piscirickettsia salmonis* in cultivated salmonids via the PCR, *Appl Environ Microbiol* 64, 3066–3069 (1998)
- Murray AE et al., 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)
- 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)

GenBank is a trademark of the US Dept of Health and Human Services. Practice of the polymerase chain reaction (PCR) may require a license. Information in this tech note was current as of the date of writing (2000) and not necessarily the date this version (rev C, 2006) 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