

Application of the Experion™ Automated Electrophoresis System DNA Assays to VNTR Analysis, SNP Typing, and Bacterial 16S rRNA Gene Analysis

Michiei Oto¹ and Wataru Suda²

¹ Department of Biotechnology, Tokyo Technical College, 4-2-3 Higashinakano, Nakano-ku, Tokyo 164-8787, and ² Graduate School of Horticulture, Chiba University, 648 Matsudo, Matsudo-shi, Chiba 271-8510, Japan

Introduction

The Experion automated electrophoresis system is a microfluidics-based electrophoretic system that can be applied to high-speed analysis of DNA fragments. In this report we evaluate the performance of the Experion system in terms of reproducibility and detection limit for DNA fragments. We apply the Experion system to three types of commonly performed DNA analyses: variable number of tandem repeats (VNTR) analysis, SNP typing of human aldehyde dehydrogenase type 2 gene (*ALDH2*), and 16S rRNA gene analysis of a bacterial community by amplified rDNA restriction analysis (ARDRA).

Materials and Methods

DNA Samples

DNA molecular standards used in this work were *HincII* digested double-stranded DNA ϕ X174 (Toyobo Life Science) and the EZ Load™ DNA molecular mass ruler (Bio-Rad Laboratories, Inc). Human genomic DNA was extracted from cheek cells of volunteers using InstaGene™ matrix (Bio-Rad). DNA samples for ARDRA were extracted from isolated bacteria and an environmental bacterial community using conventional methods (Johnson 1991).

VNTR Analysis

VNTR analysis was performed on the D1S80 locus. A D1S80 DNA sequence was amplified using 0.50 μ M of specific primers (Kasai et al. 1990) (D1SF, forward: 5'-GAAACTGGCCTCCAAACACTGCCCGCCG-3' and D1SR, reverse: 5'-GTCTTGTTGGAGATGCACGTGCCCTTGC-3'), 200 μ M dNTP, and 0.5 U AmpliTaq Gold (Applied Biosystems

Inc.) in 50 mM Tris-HCl, pH 8.0, 2 mM MgCl₂. Forty cycles of 30 sec at 94°C, 30 sec at 65°C, and 4 min at 72°C were carried out in an iCycler® thermal cycler (Bio-Rad). The PCR products were analyzed with the Experion system using the Experion DNA 1K analysis kit.

SNP Analysis of Human *ALDH2* Gene

SNP typing was performed on the human *ALDH2* gene using restriction fragment length polymorphism (RFLP) (Nachamkin et al. 2001). The exon 12 sequence of the *ALDH2* gene was amplified using 0.2 μ M of specific primers (forward: 5'-CAAATTACAGGGTCAACTGCT-3' and reverse: 5'-CCACACTCACAGTTTTCTCTT-3'), 200 μ M dNTP, and 0.5 U Discoverase (Invitrogen Corporation) in 50 mM Tris-SO₄, pH 8.0, 18 mM (NH₄)₂SO₄, 2 mM MgSO₄. Forty cycles of 15 sec at 94°C, 15 sec at 58°C, and 60 sec at 72°C were carried out in an iCycler iQ® real-time PCR detection system. The PCR products were digested with the restriction endonuclease *MbolI* followed by analysis with the Experion system using the Experion DNA 1K analysis kit.

ARDRA of the 16S rRNA Gene

The nearly full-length 16S rRNA gene sequence was amplified using 0.2 μ M of specific primers (27f, forward: 5'-AGAGTTTGATCMTGGCTCAG-3' and 1525r, reverse: 5'-AAGGAGGTGWTCARCC-3'), 200 μ M dNTP, and 0.5 U Discoverase (Invitrogen) in 50 mM Tris-SO₄, pH 8.0, 18 mM (NH₄)₂SO₄, 2 mM MgSO₄. Forty cycles of 30 sec at 94°C, 30 sec at 58°C, and 60 sec at 72°C were carried out in an iCycler thermal cycler. The PCR products were digested with restriction endonucleases (*HaeIII*, *AluI*, and *Sau3AI*), followed by analysis with the Experion system using the Experion DNA 1K analysis kit. ARDRA is based on the RFLP technique (Nachamkin et al. 2001).

Results and Discussion

Performance of the Experion system

Resolution, Accuracy and Reproducibility — The fragment resolution, sizing accuracy, and reproducibility of the Experion system were evaluated using a *HincII* digest of double-stranded DNA ϕ X174.

The Experion DNA 1K assay sizing resolution range for DNA fragments is between 15 and 1500 bp, which is sufficient for RFLP analysis. Particularly for the ~200 bp DNA fragments, the Experion system exhibited sufficient resolution to detect size differences of less than 10 bp. Fragments of 291, 297, 335, 345, and 392 bp were clearly separated by electrophoresis on the Experion DNA 1K analysis kit while they appear as a single band on an agarose gel (Figure 1). The separation and analysis using the Experion system was performed in about 50 min, compared to agarose gel electrophoresis which takes about 120 min (Figure 2). The sizing accuracy, intra-chip deviation (%CV), and inter-chip deviation (%CV) were $\pm 1.03\%$, 0.62% , and 1.06% , respectively, indicating very good accuracy and reproducibility.

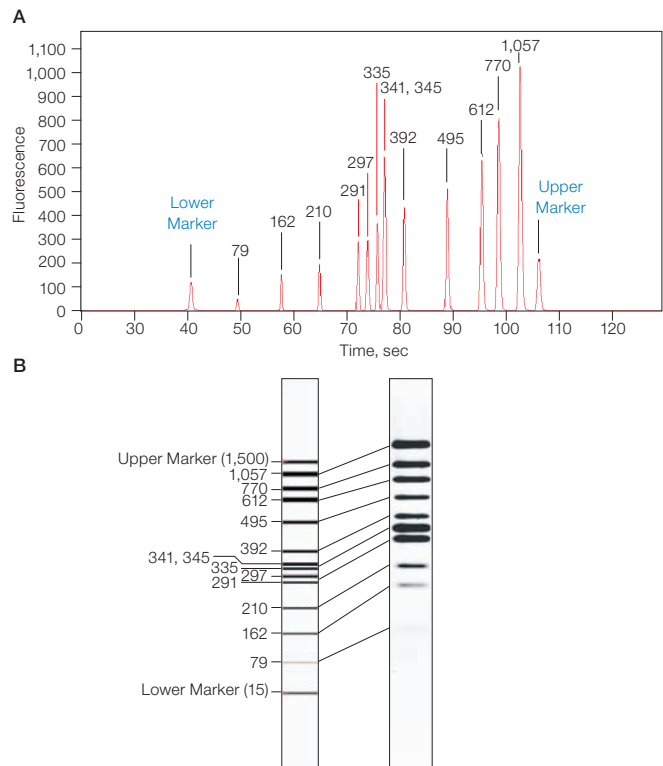


Fig. 1. Resolution of the Experion system and agarose gel electrophoresis. **A**, electropherogram of ϕ X174 *HincII* digested DNA; **B**, comparison of the Experion system gel image and 3% agarose gel image of ϕ X174 *HincII* digested DNA. The assay procedures are shown in Figure 2. The size of the bands is indicated in bp.

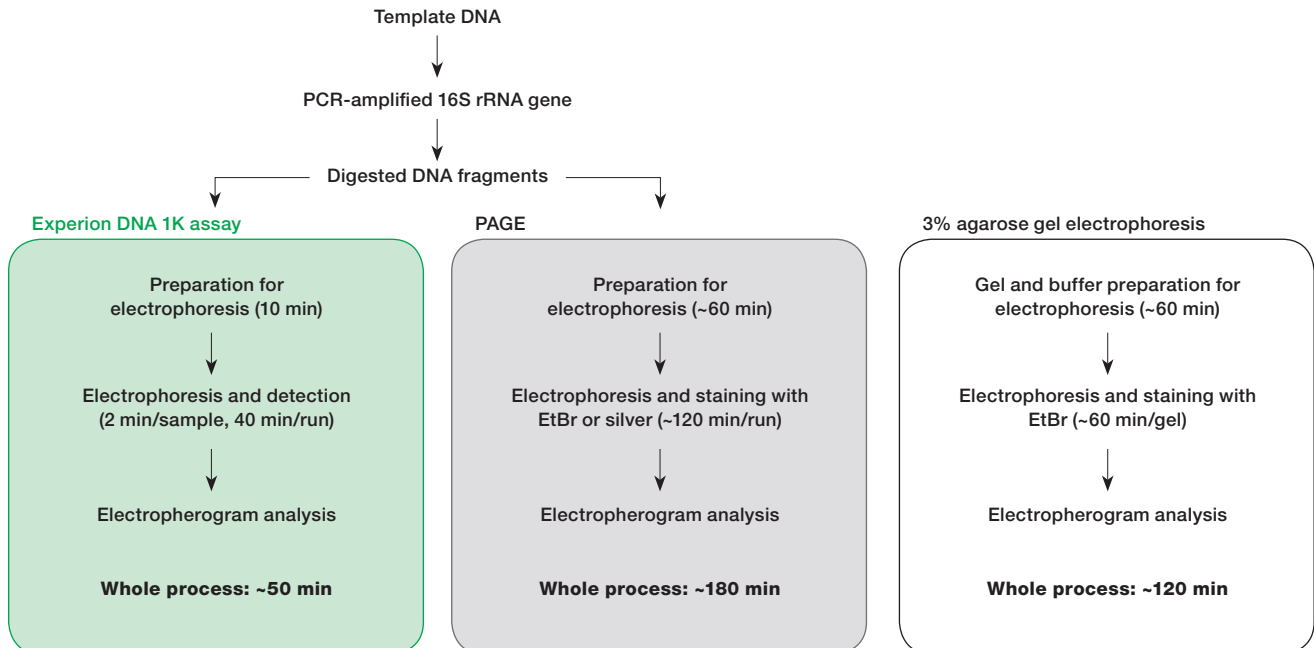


Fig. 2. Comparison of workflow and assay time. The workflow and process times for PCR-amplified DNA electrophoresis with the Experion system, polyacrylamide gel, and agarose gel are compared.

Limits of Detection — To compare the detection limits of the Experion system and agarose gel electrophoresis fivefold serial dilutions of the EZ Load DNA molecular mass ruler were analyzed with both systems. The results show that with the Experion system the limit of detection of a 100 bp fragment is 80 pg, while it is more than 400 pg on an agarose gel (Figure 3).

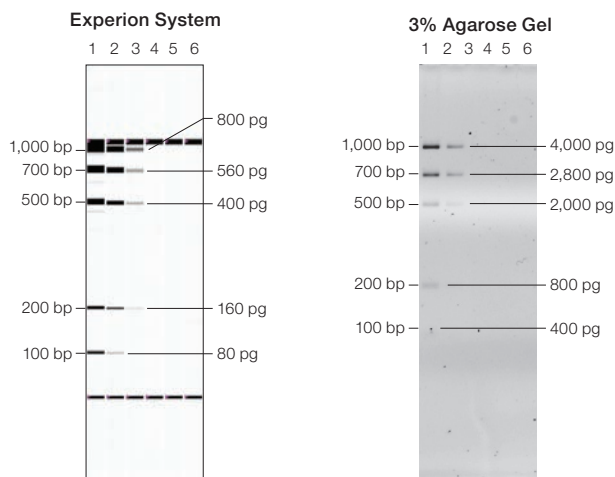


Fig. 3. Limit of detection of Experion system and agarose gel electrophoresis. Fivefold serial dilutions of EZ Load DNA molecular mass ruler run on the Experion system (left) and on a 3% ethidium bromide-stained agarose gel (right).

Evaluation of VNTR Analysis, SNP Typing, and Bacterial 16S rRNA Gene Analysis

VNTR Analysis — Typical electrophoretic profiles of amplified D1S80 DNA fragments are shown in Figure 4. The profiles were reproducible (data not shown) and coincided with those obtained on conventional PAGE analysis, indicating that data of the same quality can be obtained with the Experion assay.

SNP Genotyping — Codon 487 of the *ALDH2* gene has three kinds of SNPs, which are categorized as G/G homozygote, G/A heterozygote, and A/A homozygote. These polymorphisms are found mainly in Asian people including Japanese (Shibuya and Yoshida 1988). To assess the utility of the Experion system in SNP genotyping we performed a genotyping test on the DNA of 10 individuals. A 135 bp fragment containing codon 487 was amplified, digested with *Mbol*I, and analyzed using an Experion DNA 1K analysis kit. Since G and A peaks were completely separated, analysis with the Experion system enabled typing of these three kinds of SNPs (Figure 5) within 2 min (per sample), showing that quality data, suitable for practical high-speed analysis of SNPs, can be obtained. For DNA samples derived from 10 individuals, the SNP typing data with the Experion system correlated well with those obtained on DNA sequencing (data not shown).

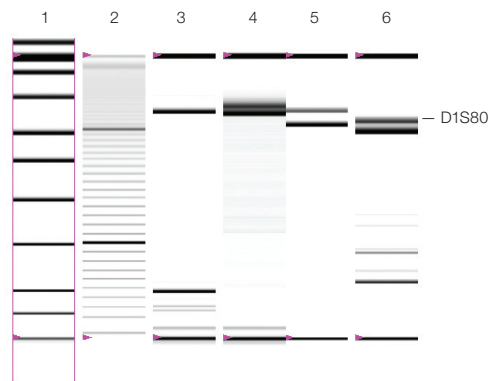


Fig. 4. D1S80 analysis using the Experion system. Lanes 1 and 2, 1K ladder and 20 bp ladder, respectively. Lanes 3–6, amplified D1S80 DNA from different individuals.

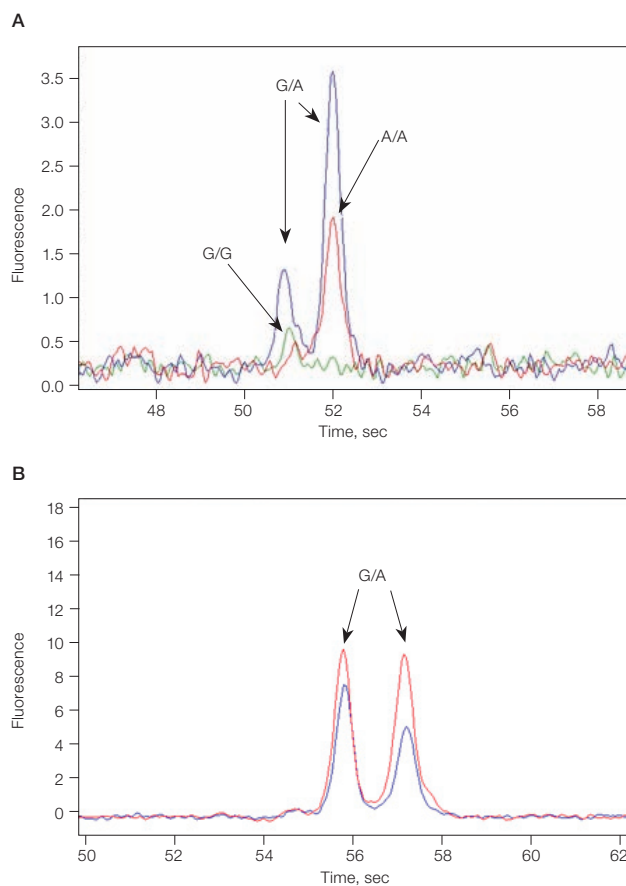
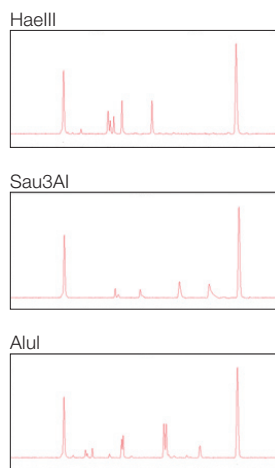


Fig. 5. *ALDH2* SNP analysis using the Experion system. **A**, electrophoretic profile of a *Mbol*I digested PCR fragment containing codon 487 of *ALDH2*. — G/G homozygous DNA, — G/A heterozygous DNA, and — A/A homozygous DNA; **B**, G/A heterozygous DNA from different samples.

16S rDNA Analysis — ARDRA with the Experion system was optimized and then used to identify isolated bacteria and to model the identification of a bacterium in an environmental sample. The optimized analytical procedure was performed within 2 min per sample in a total of 50 min (Figure 2), indicating this is a much quicker analytical technique than conventional PAGE (180 min). ARDRA electropherograms from the Experion system and conventional PAGE are shown in Figure 6. These data indicated that the resolution with the Experion system was the same or better than for PAGE.

PCR-amplified 16sRNA genes from *E. coli*, *B. subtilis*, and *S. aureus* were digested with HaeIII, Sau3AI, and AluI and analyzed by electrophoresis on the Experion system and by 10% PAGE. The sizes of the fragments estimated from their electrophoretic migration pattern were compared with the sizes of the fragments deduced from the gene sequence. Results shown in Table 1 indicate that size estimation with the Experion system is very accurate.

A. Experion System



B. PAGE

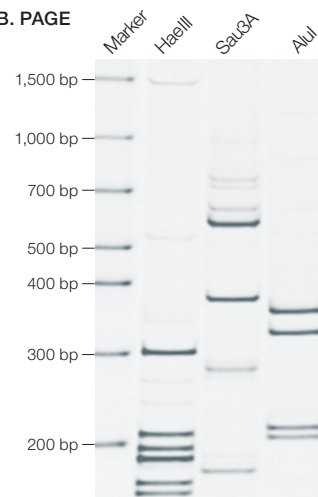


Fig. 6. Typical ARDRA profile. 16S rRNA gene from *E. coli* K12 was amplified and digested with HaeIII, Sau3AI, and AluI, followed by electrophoretic analyses with the Experion system and 10% PAGE. **A**, electropherograms obtained with the Experion system; **B**, electrophoretic profile on PAGE.

Table 1. DNA fragment size estimation. The size of DNA fragments (in base pair number) determined experimentally by electrophoresis on the Experion system (Experion) or on a 10% polyacrylamide gel (PAGE) is compared to the predicted size based on sequence data (In silico).

			Fragment sizes (bp)									
			34	39	67	154	161	167	180	205	210	317
HaeIII	<i>E. coli</i>	In silico	34	39	67	154	161	167	180	205	210	317
	Experion	—	—	67	157	165	177	180	207	207	317	
	PAGE	—	—	112	130	165	180	205	305	—	—	
HaeIII	<i>B. subtilis</i>	In silico	22	155	309	457	599	—	—	—	—	—
	Experion	—	155	314	447	610	—	—	—	—	—	
	PAGE	—	120	305	375	560	—	—	—	—	—	
HaeIII	<i>S. aureus</i>	In silico	310	1,234	—	—	—	—	—	—	—	—
	Experion	312	1,392	—	—	—	—	—	—	—	—	
	PAGE	305	1,200	—	—	—	—	—	—	—	—	
Sau3AI	<i>E. coli</i>	In silico	7	12	174	267	424	651	—	—	—	—
	Experion	—	—	178	266	434	652	—	—	—	—	
	PAGE	—	—	140	280	335	560	—	—	—	—	
	<i>B. subtilis</i>	In silico	7	12	51	119	175	187	296	695	—	—
	Experion	—	—	54	117	176	189	300	689	—	—	
	PAGE	—	—	—	—	140	175	295	630	—	—	
	<i>S. aureus</i>	In silico	7	12	74	119	175	223	352	582	—	—
	Experion	—	—	75	116	175	225	365	583	—	—	
	PAGE	—	—	—	140	170	235	350	530	—	—	
AluI	<i>E. coli</i>	In silico	75	86	99	122	207	211	361	374	—	—
	Experion	79	85	100	—	203	209	363	373	586	—	
	PAGE	—	—	—	—	210	215	325	355	—	—	
	<i>B. subtilis</i>	In silico	73	173	186	207	208	265	430	—	—	
	Experion	78	174	184	205	210	261	435	—	—		
	PAGE	—	155	195	—	245	—	385	—	—		
	<i>S. aureus</i>	In silico	51	74	84	86	87	124	209	214	615	
	Experion	—	—	—	86	86	122	201	212	646	—	
	PAGE	—	—	—	—	—	135	210	215	620	—	

To evaluate the sensitivity of 16S rDNA detection, an *E. coli* 16S rDNA-HaeIII digest was mixed with an *E. coli* 16S rDNA-AluI digest in different ratios and analyzed by electrophoresis with the Experion system. The *E. coli* 16S rDNA-AluI digest could be detected even when representing as little as 11% of the DNA present in the mixture (green line in Figure 7), indicating that specific DNA fragments might be detected in a DNA sample from an environmental bacterial community if the bacterium represents 10% of the bacterial community.

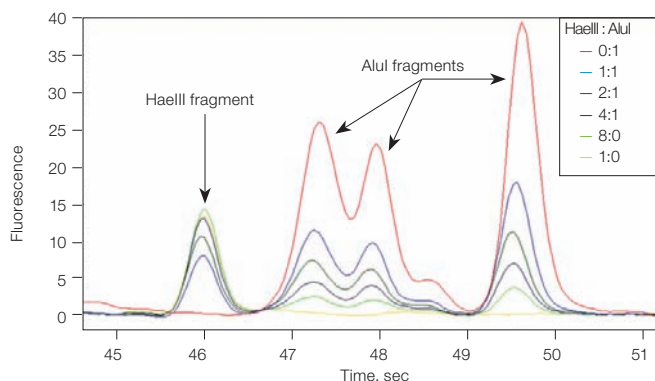


Fig. 7. Limit of detection in the ARDRA using the Experion system. An *E. coli* 16S rDNA HaeIII digest was mixed with an *E. coli* 16S rDNA AluI digest in different ratios (inset) and electrophoresed with the Experion system.

To assess the use of the Experion system to identify a bacterium in an environmental bacterial community, various amounts (0.4, 10, and 50 ng) of *E. coli* 16S rDNA-Sau3AI digest were spiked into 20 ng of a 16S rDNA-Sau3AI digest from a microbial community prior to analysis with the Experion system (Figure 8). The spiked DNA was detectable as a peak, indicating the possibility of identifying a specific bacterium in an environmental bacterial community by using the Experion system as well as conventional 16S rDNA analytical techniques including denaturing gradient gel electrophoresis (DGGE) (Muyzer et al. 1993), and single-strand conformational polymorphism (SSCP) (Oto et al. 2006).

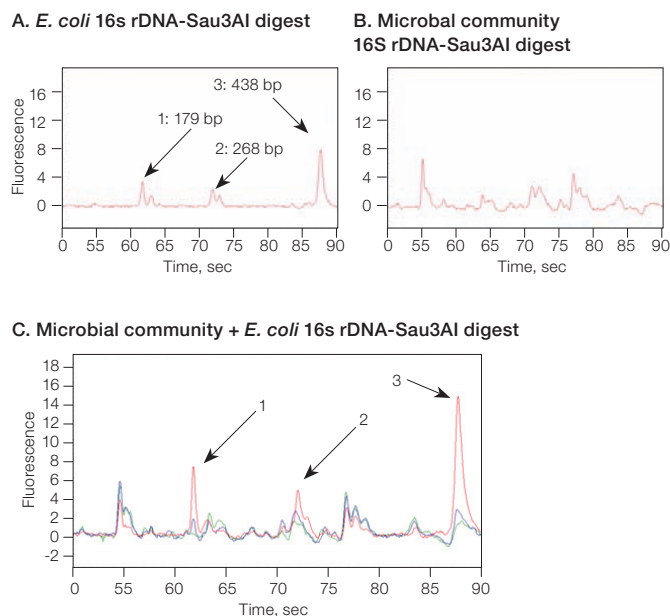


Fig. 8. Detection of target bacteria in a microbial community. Amplified 16S rDNA from *E. coli* digested with Sau3AI was spiked into 20 ng of microbial community DNA in different amounts (— 0.4, — 10, and — 50 ng). **A**, electropherogram of *E. coli* 16S rDNA; **B**, electropherogram of microbial community 16S rDNA; **C**, electropherogram of mix of *E. coli* and bacterial community 16S rDNA. **1**, **2**, and **3** indicate different peaks derived from the amplified *E. coli* 16S rDNA.

Conclusion

With the Experion system, it is possible to perform high-speed VNTR analysis, SNP typing, and bacterial 16S rRNA gene analysis. In particular, high-speed ARDRA with the Experion system may be a powerful tool for analyzing not only isolated bacteria but also microbial communities. Additionally, the Experion system is preferable for DNA length-based analyses, including sizing of PCR products and RFLP. Although the analytical formats such as the matrix concentration, buffer conditions, and running time are fixed, the Experion system is superior to gel electrophoresis in speed, resolution, detection limit, and simplicity.

References

Johnson JL (1991). Isolation and purification of nucleic acids. In *Nucleic Acid Techniques in Bacterial Systematics*, ed. (West Sussex, England: John Wiley & Son), pp. 1–19.

Kasai K et al. (1990). Amplification of a variable number of tandem repeats (VNTR) locus (PMCT118) by the polymerase chain reaction (PCR) and its application to forensic science. *J Forensic Sci* 35, 1196–1200.

Muyzer G et al. (1993). 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.

Nachamkin I et al. (2001). Agilent 2100 Bioanalyzer for restriction fragment length polymorphism analysis of the *Campylobacter jejuni* flagellin gene. *J Clin Microbiol* 39, 754–757.

Oto M et al. (2006). 16S rRNA gene-based analysis of microbial community by whole-genome amplification and minigel-single-strand conformation polymorphism technique. *Journal of Bioscience and Bioengineering* 102, 482–484.

Shibuya A and Yoshida A (1988). Frequency of the atypical aldehyde dehydrogenase-2 gene (ALDH2(2)) in Japanese and Caucasians. *Am J Hum Genet* 43, 741–743.

AmpliQ Gold is a registered trademark of Applied Biosystems. Discoverase is a trademark of Invitrogen Corporation.



LabChip and the LabChip logo are trademarks of Caliper Life Sciences, Inc. Bio-Rad Laboratories, Inc. is licensed by Caliper Life Sciences, Inc. to sell products using the LabChip technology for research use only. These products are licensed under U.S. Patent

Nos. 5,863,753, 5,658,751, 5,436,134, and 5,582,977, and pending patent applications, and related foreign patents, for internal research and development use only in detecting, quantitating, and sizing macromolecules, in combination with microfluidics, where internal research and development use expressly excludes the use of this product for providing medical, diagnostic, or any other testing, analysis, or screening services, or providing clinical information or clinical analysis, in any event in return for compensation by an unrelated party.

Information in this tech note was current as of the date of writing (2010) and not necessarily the date this version (rev A, 2010) was published.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 31 3689 6600
Canada 905 364 3435 **China** 86 20 8732 2339 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00 **France** 01 47 95 69 65
Germany 089 31 884 0 **Greece** 30 210 777 4396 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **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