

Detection of Mutant *K-ras* in a Kindred With Hereditary Pancreatic Cancer by Denaturing Gradient Gel Electrophoresis

David Crispin,¹ Ru Chen,¹ Michael Kimmey,² and Teresa Brentnall²
Departments of ¹ Pathology and ² Medicine, Division of Gastroenterology,
University of Washington, Seattle, Washington 98195 USA

Introduction

Pancreatic cancer is the fourth leading cause of cancer death in the US, and its frequency is rising. At the time of diagnosis, 96–99% of patients are incurable and will shortly die. Current methods to evaluate patients for pancreatic adenocarcinoma include endoscopic retrograde cholangiopancreatography (ERCP), abdominal CT, ultrasound, and serum markers. These methods of diagnosis are often insensitive or equivocal in early disease. Tumorigenesis is believed to involve the *K-ras* oncogene and the *DCC*, *p16*, *APC*, *bcl-2*, and *p53* tumor suppressor genes; screening for mutations or loss of heterozygosity in these genes may provide better diagnostic tests that are more sensitive and specific (Rozenblum et al. 1997). The study of families in which cancer is inherited in an autosomal dominant fashion has provided considerable insight into understanding the molecular basis for pancreatic cancer. We have previously reported an extensive kindred in which pancreatic cancer is inherited in an autosomal dominant fashion and is associated with development of pancreatic insufficiency prior to the diagnosis of cancer (Evans et al. 1995). We have utilized the DCode™ universal mutation detection system for denaturing gradient gel electrophoresis (DGGE) to identify patient samples with *K-ras* mutations.

Methods

Genomic DNA was isolated from tissue or from fluid obtained from patients at ERCP. *K-ras* exon 1 was amplified from genomic DNA using a thermal cycler and primers as described by Imai et al. (1991). Amplifications were carried out using 300 ng of DNA template in a buffer containing 10 mM Tris-HCl, 1.5 mM MgCl₂, 50 mM KCl, pH 8.3, with 2 U of Taq DNA polymerase (Boehringer Mannheim), 200 M dNTPs, and 15 pmol of each primer in a total volume of 50 L. Reactions were denatured for 3 min at 95°C, which was followed by 37 cycles with the following profile: 95°C (20 sec), 55°C (60 sec), 72°C (40 sec). The 110 base pair (bp) product was run on a 3% agarose gel and visualized using ethidium bromide staining.

Successful samples were run on a 10% acrylamide 0–80% perpendicular DGGE gel at 150 V for 2 hr at 56°C, then stained with ethidium bromide to find the optimal conditions for parallel DGGE (Figure 1). This was determined to be 30–60%. Samples were then run on a parallel DGGE gel at 56°C for 4–5 hr at 150 V, and stained with ethidium bromide. Positive samples were sequenced using dye terminator chemistry and run on an ABI PRISM instrument (PerkinElmer, Inc.).

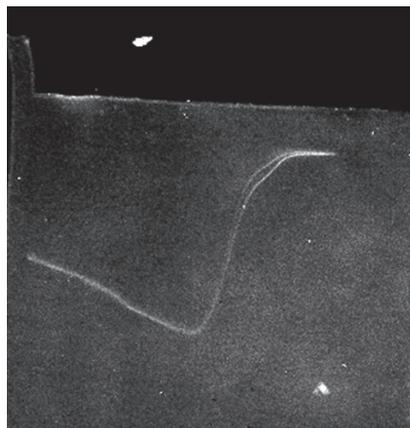


Fig. 1. 0–80% perpendicular DGGE with wild-type and mutant *K-ras* products. The fragment melts at a denaturant concentration of 48%.

Table 1. *K-ras* mutations in family X.

| III.1 PreOp Juice* | III.2 PreOp Juice* | IV.19 Metaplasia | III.15 Dysplasia | III.16 Cancer In Situ | III.6 Cancer | III.17 Cancer | III.19 Metastasis |
|-----------------------|-----------------------|---------------------|---------------------|--------------------------|---------------------|---------------------|----------------------|
| Codon 12 GGC→GAT | Codon 12 GGT→GAT | Codon 13 GGC→GAC | None | Codon 12 GGC→XXX | Codon 13 GGC→GAC | Codon 13 GGC→AGC | None |

*For those patients who had not undergone pancreatectomy yet, pancreatic juice was obtained at ERCP and tested for *K-ras* mutation.

Results

Exon 1, containing codons 12 and 13, of the *K-ras* gene was examined for the presence of *K-ras* mutations by DGGE. Constitutional and pancreatic tissues or pancreatic juice in eight of the family members were evaluated. All samples that were positive by DGGE (Figure 2) were confirmed by DNA sequencing. Three of these individuals had *K-ras* mutations in codon 13 present in pancreatic cancer or precancerous tissue, and three had mutations in codon 12. There was no evidence of *K-ras* mutation in the metastatic pancreatic cancer tissue from individual III.19, even when DNA was subcloned, or from the ERCP fluid or dysplastic tissue from III.15 (Table 1).

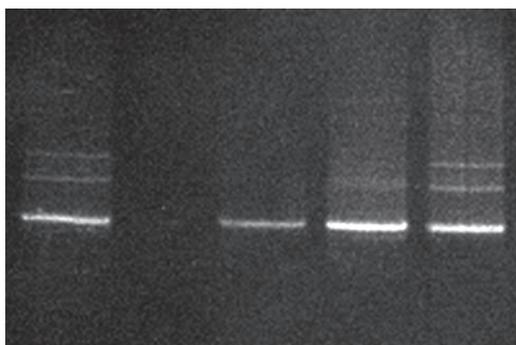


Fig. 2. Parallel DGGE of *K-ras* exon 1. Mutant and wild-type samples were run with a 30–60% denaturing gradient. Lane 1, positive control; lane 2, negative control; lane 3, ERCP fluid from patient III.1; lane 4, ERCP fluid from patient III.16.

Discussion

K-ras mutations are a common event in pancreatic adenocarcinoma. Screening for *K-ras* mutations along with a panel of other markers may prove useful in early diagnosis of this disease. DGGE is a practical tool in screening samples for the presence of such mutations.

References

- Evans JP et al., Familial pancreatic adenocarcinoma: association with diabetes and early molecular diagnosis, *J Med Genet* 32, 330–335 (1995)
- Imai M et al., *K-ras* codon 12 mutations in biliary tract tumors detected by polymerase chain reaction denaturing gradient gel electrophoresis, *Cancer* 73, 2727–2733 (1991)
- Rozenblum E et al., Tumor-suppressive pathways in pancreatic carcinoma, *Cancer Res* 57, 1731–1734 (1997)

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, 2007) was published.

BIO-RAD

**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