

## Detection of Mutations in the *CYP21* Gene Using the DCode™ System

Gita Ohlsson and Marianne Schwartz, Department of Clinical Genetics, Juliane Marie Center, University Hospital, Rigshospitalet, Copenhagen, Denmark

### Introduction

Congenital adrenal hyperplasia (CAH) due to 21-hydroxylase deficiency is caused by mutations in the gene *CYP21* encoding the enzyme steroid 21-hydroxylase. In addition to deletions, approximately 20 different point mutations have been reported (Barbat et al. 1995, Ezquieta et al. 1995, Speiser et al. 1992, Wedell et al. 1994) and novel mutations are still detected (Kirby-Keyser et al. 1997, Lajic and Wedell 1996, Lajic et al. 1997, Levo and Partanen 1997). This makes genetic diagnosis as well as carrier detection of 21-hydroxylase deficiency a complicated matter. We have used the denaturing gradient gel electrophoresis method (DGGE) using the DCode universal mutation detection system to detect mutations in the coding sequence and intron-exon junctions of *CYP21* (Guldberg and Güttler 1993, Sheffield et al. 1989).

### Methods

#### Samples

DNA from healthy individuals and patients with 21-hydroxylase deficiency was isolated from peripheral blood lymphocytes using the salting-out method by Miller et al. (1988).

#### Polymerase Chain Reaction (PCR)

For amplification of *CYP21* and selection against *CYP21P*, PCR was performed according to Wedell and Luthman (1993). The resulting two PCR products were purified and subsequently used as template for the amplification of DNA fragments for DGGE. The PCR reactions were performed with GC-clamped primers. The GC density of almost all *CYP21* exons are very high, requiring long GC clamps to obtain a single melting domain of the sequence of interest.

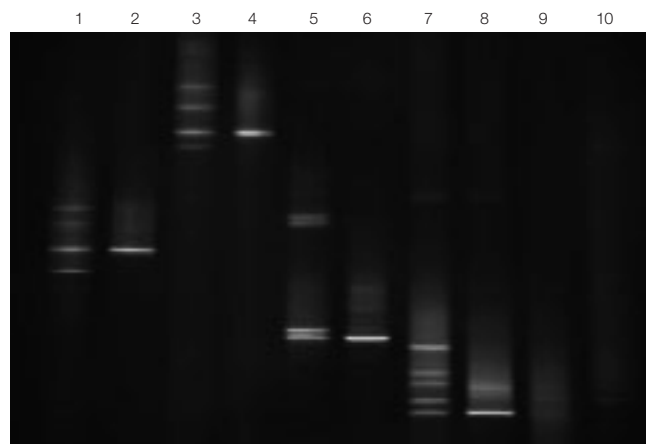
#### Site-Directed Mutagenesis

In order to obtain mutant control samples for all exons, it was necessary to introduce mutations in some DNA fragments. This was done by PCR-based site-directed mutagenesis (Kuipers et al. 1991, Landt et al. 1990). The mutant PCR product was

mixed with the corresponding wild-type PCR product; the mixture was placed at 96°C for 10 min to denature the PCR products and subsequently left at room temperature for gradual renaturation, thereby generating heteroduplex molecules.

#### DGGE

DGGE was carried out using the DCode system (Bio-Rad Laboratories, Inc.). PCR product (15 µl) was loaded on a gel containing a polyacrylamide gradient ranging from 6–12% and a gradient of urea and formamide. The gels were run in 1x TAE buffer, at 80 V overnight. After electrophoresis, the gels were stained in TAE buffer containing ethidium bromide, and subsequently, the resolved bands were visualized by ultraviolet (UV) transillumination.



**Fig. 1. DGGE analysis of *CYP21*.** Each exon is represented by a mutant and a normal control sample. Lanes 1–2, exon 2; lanes 3–4, exon 5; lanes 5–6, exon 6; lanes 7–8, exon 7. For exon 2 and 5, the mutant control samples were obtained by site-directed mutagenesis at the following nucleotide positions: exon 2: 2011 (A→C) and exon 5: 2851 (T→G). Nucleotide positions are given in accordance with *CYP21*, GenBank accession numbers: M12792; M23280. For exon 6, the mutant control sample represents the cluster-E6 mutation (lane 5) and for exon 7, the Val281Leu mutation (lane 7).

## Results

The results of the DGGE analysis of exons 2, 5, 6, and 7 are presented in Figure 1. The presented DGGE analysis was carried out using a denaturing gradient of 30–70%. To validate the method, one wild-type control sequence and one heterozygous mutant sequence was analyzed for each exon. For exons 2 and 5, the mutant controls were generated by site-directed mutagenesis at the following nucleotide positions: exon 2: 2011 (A→C) and exon 5: 2851 (T→G). Nucleotide positions are given in accordance with CYP21, GenBank: accession numbers M12792; M23280. The mutant control sample for exon 6 contains the cluster-E6 mutation and for exon 7 the Val281Leu mutation.

## Discussion

To perform complete genotyping of 21-hydroxylase disease alleles and reliable carrier diagnosis, it is essential to use a technique that not only detects the most frequent previously identified point mutations but also both undefined and rare point mutations. The DGGE analysis fulfills these criteria, and therefore constitutes a fast and reliable procedure for genetic analysis of 21-hydroxylase deficiency.

## References

- Barbat B et al., Screening of CYP21 gene mutations in 129 French patients affected by steroid 21-hydroxylase deficiency, *Hum Mutat* 5, 126–30 (1995)
- Ezquieta B et al., Analysis of steroid 21-hydroxylase gene mutations in the Spanish population, *Hum Genet* 96, 198–204 (1995)
- Guldberg P and Güttler F, A simple method for identification of point mutations using denaturing gradient gel electrophoresis, *Nucleic Acids Res* 21, 2261–2262 (1993)
- Kirby-Keyser L et al., E380D: a novel point mutation of CYP21 in an HLA-homozygous patient with salt-losing congenital adrenal hyperplasia due to 21-hydroxylase deficiency, *Hum Mutat* 9, 181–182 (1997)
- Kuipers OP et al., Improved site-directed mutagenesis method using PCR, *Nucleic Acids Res* 19, 4558 (1991)
- Lajic S and Wedell A, An intron 1 splice mutation and a nonsense mutation (W23X) in CYP21 causing severe congenital adrenal hyperplasia, *Hum Genet* 98, 182–184 (1996)
- Lajic S et al., A cluster of missense mutations at Arg356 of human steroid 21-hydroxylase may impair redox partner interaction, *Hum Genet* 99, 704–709 (1997)
- Landt O, Grunert HP, and Hahn U, A general method for rapid site-directed mutagenesis using the polymerase chain reaction, *Gene* 96, 125–128 (1990)
- Levo A and Partanen J, Novel nonsense mutation (W302X) in the steroid 21-hydroxylase gene of a Finnish patient with the salt-wasting form of congenital adrenal hyperplasia, *Hum Mutat* 9, 363–365 (1997)
- Miller SA et al., A simple salting out procedure for extracting DNA from human nucleated cells, *Nucleic Acids Res* 16, 1215 (1988)
- 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)
- Speiser PW et al., Disease expression and molecular genotype in congenital adrenal hyperplasia due to 21-hydroxylase deficiency, *J Clin Invest* 90, 584–595 (1992)
- Wedell A et al., Mutational spectrum of the steroid 21-hydroxylase gene in Sweden: implications for genetic diagnosis and association with disease manifestation, *J Clin Endocrinol Metab* 78, 1145–1152 (1994)
- Wedell A and Luthman H, Steroid 21-hydroxylase deficiency: two additional mutations in salt-wasting disease and rapid screening of disease-causing mutations, *Hum Mol Genet* 2, 499–504 (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](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