

## **NON-INVASIVE PRENATAL PATERNITY TESTING USING NEXT GENERATION SEQUENCING AND BIOINFORMATICS BASED ANALYSIS**

Richard Chmelo<sup>1</sup>, Ph.D., Tudor Constantin<sup>2</sup>, Ph.D., Allison Ryan<sup>2</sup>, Ph.D., Dmitry Grudzinskiy<sup>2</sup>, Ph.D., Eser Kirkizlar<sup>2</sup>, Ph.D., Kacee Floyd<sup>1</sup>, MT, Jonathan Sheena<sup>2</sup>, Ph.D., Michael Baird<sup>1</sup>, Ph.D., and Matthew Rabinowitz<sup>2</sup>, Ph.D.

<sup>1</sup>DNA Diagnostics Center

<sup>2</sup>Natera

Prenatal paternity testing has been available for several decades using chorionic villus sampling or amniotic fluid collection by amniocentesis. Both of these procedures have a small risk of miscarriage to the developing fetus (1). The discovery that cell-free fetal DNA is present in the blood of pregnant women specific to the current pregnancy (2) offers the opportunity to develop a safer test for the determination of paternity prenatally. Here we describe the use of a set of SNPs interrogated by next generation sequencing combined with a version of the Parental Support algorithm develop by the Natera Corporation to determine paternity prenatally and outline the validation performed.

1. American College of Obstetricians and Gynecologists. Invasive prenatal testing for aneuploidy. ACOG Practice, Bulletin No. 88, December 2007. *Obstet Gynecol* 2007; 110:1459-1467.
2. Lo, YM, Corbetta, N, Chamberlain, PF, et al, Presence of fetal DNA in maternal plasma and serum. *Lancet* 1997; 350; 485-487