

## **IDENTIFICATION OF ALLEGED FATHER (AF) USING INCEST CHECKER SOFTWARE IN A SEXUAL ASSAULT CASE: A CASE REPORT**

Gang-nam Jin<sup>1</sup>, Dong-sub Lee<sup>2</sup>, Eun-jung Lee<sup>1</sup>, Man-il Kim<sup>1</sup>, Jung-yoon Lee<sup>1</sup>

<sup>1</sup>Forensic DNA Division, Gwangju institute, National Forensic Service

<sup>2</sup>Forensic DNA Division, Seoul institute, National Forensic Service

The female victim, who was 14 years old, was pregnant by being raped from someone. The female victim's father was suspected as the biological father of the victim's fetus, but the victim denied that suspicion. The Prosecutor cautiously asked a genetic analysis on the abortive fetus, on the victim and on her father and mother, but the exclusion at five loci was found. So, we have developed user friendly software "Incest Checker" to solve this problem. Finally, we have proposed that the most probable situation would be the existence of victim's brother. The victim's brother admitted the crime later.

Keywords: Incest; Alleged father; genetic analysis software