

Introduction of the Korean DNA Identification in National Forensic Services

Chong Min Choung, Yang Han Lee, Jong Jin Kim, Myun Soo Han
Forensic DNA Center, National Forensic Service, Seoul, Korea

Forensic DNA Center of the National Forensic Service in Seoul has been established in 1991 and internationally approved by Korea Laboratory Accreditation Schemer (KOLAS) in August 2004. Forensic DNA Center has 3 divisions-DNA analysis division, DNA index division, DNA information and ID division. These three divisions are organized 9 parts -Forensic DNA analysis section 1, Forensic DNA analysis section II, Forensic DNA analysis section III, DNA index section I, DNA index section II, Missing persons identification section, DNA paternity section, and DNA database management section. And we have four regional branches.

Our center performs DNA identification for all autopsy samples even there is identification or some evident to identify the body and search in the database first prior to record it into the DNA database. Forensic DNA center had completed more than 100,000 crime cases in 2011. And total of two hundred seventy seven unidentified persons were recorded into the DNA database in 2011. We would like to introduce Korean DNA identification system and current issues.