

THE LATEST TREND IN PROPERTY CRIME DNA ANALYSIS IN METROPOLITAN REGION OF KOREA AND INTRODUCTION OF EXTRAORDINARY CASEWORKS WHICH INCLUDES THE MIXTURE ANALYSIS.

NaYeon Kim, DNA Analysis Division, Seoul Institute, National Forensic Service

The forensic DNA analysis of criminal samples is very important not only in the severe case such as murder, robbery and sexual crimes but also in the ordinary life property crimes such as larceny, destruction and damage of property, fraud and so on. The proportion of property crimes in forensic DNA analysis demand has been continuously increased during last 10 years especially after the forensic DNA database launching and expansion in Republic of Korea.

The improvement of LCN DNA analysis technology and of database matching rate result in the dramatic uprising request for property crime DNA analysis, and now approximately 50% of DNA request from the crime scene is related with the property crimes in metropolitan region in Korea.

In this report, we summarized the latest trends in various property crime DNA analysis and database searching results during last two years in the metropolitan area in Republic of Korea, and introduced the special caseworks which include the serial offenders and the mixture interpretation.