

## **COMPARISON OF DNA QUANTIFICATION KITS THAT MONITOR THE INTEGRITY OF FORENSIC DNA SAMPLES**

Tetsushi Kitayama, National Research Institute of Police Science

For the purpose of assessing the quality and quantity of DNA, we had previously developed two quantitative PCR assays determining degree of DNA degradation using primers targeting short and long amplicons in human specific multi-copy regions. We have improved previous assays and developed a multiplexed system to monitor the integrity of DNA more precisely.

Recently, some DNA quantification kits which can monitor the degree of DNA degradation were made available. We compare sensitivity, precision, and ability to predict the integrity of forensic DNA samples between our quantification system and other quantification kits.