

## **DEVELOPMENT OF LIMS FOR FORENSIC LABORATORIES IN JAPAN**

Kanako Kitamori, Nippon Control System Corporation

We make the SimpDoc: a Laboratory Information Management System (LIMS), and we have been work with forensic laboratories of police in all over Japan. LIMS is an indispensable tool to batch analysis work in recent years. The analysis data is a large amount, and you are not allowed to make a mistake. However, the workflow is vary depending on the size and role of the laboratory. And we think there are some cultural differences between Asia and Western countries. For example, a way of thinking about permission of each workers, necessary paper documents, and others. Especially in Japan, it is important to fit to user's workflow when make a system like LIMS. In this way, we are studying to make our LIMS to a more efficient and flexible software.