VALIDATION AND APPLICATION OF COMMERCIAL MINISTR TYPING KIT FOR KOREAN NATIONAL MISSING PEOPLE IDENTIFICATION PROJECT

K.S. Kim¹, Y.J. Kim¹, S.H. Kim¹, S.B. Hong¹, J.H. Kim¹, N.Y. Kim¹, K.W. Park¹, J.J. Kim¹, W. Kim², M.S. Han¹

National Institute of Scientific investigation is an official DNA identification center for missing people and unidentified human remains in South Korea. We conducted the DNA analysis over 17,000 missing people and 1,500 human remains for last 5 years. But we have met the problematic cases which conventional STR analysis systems could not detect STR genotype of highly degraded DNA specimens such as human skeletal remains.

Recently developed ABI AmpF{STR®MiniFiler™ kit is a good alternative for these highly degraded DNA samples because we use 17 STR loci (13 CODIS, D21338, D19S433, Penta E, and Penta D) for missing people database.

We participated in the validation project for the recently developed MiniFiler™ kit as South Korea partner, and validated the commercial MiniSTR typing kit for its application to forensic casework samples. Here we report on the usefulness of the MiniFiler™ kit in STR genotyping of highly degraded samples.

¹Korean National Institute of Scientific Investigation, Seoul, Republic of Korea,

²Dankook University, Cheonan, Republic of Korea