## INITIAL INVESTIGATION OF THE ERASE SPERM ISOLATION KIT

Nicole Hinton, M.S., Stephanie Kuntz, M.S., Sarah Ozanick, Ph.D., and Gretchen DeGroot, M.S. Wisconsin State Crime Lab, DNA Analysis, 1578 S 11<sup>th</sup> St, Milwaukee, WI 53204, USA

The Erase Sperm Isolation Kit from PTC Laboratories is a differential extraction system that uses a nuclease to selectivity degrade epithelial cell DNA prior to sperm cell lysis. This method reportedly provides single source male autosomal profiles from samples that would be expected to yield mixtures due to the presence of high concentrations of epithelial cell DNA. We have processed duplicate samples using either the Wisconsin State Crime Laboratory protocol for a differential extraction or the Erase Sperm Isolation Kit protocol. Substantial differences in the amount of DNA recovered and the profiles generated by the two methods will be presented. **3**