

CASEWORK EXPERIENCE AND INTERESTING CASES WITH THE GLOBALFILER® CASEWORK AND EXPRESS KITS

Nicholas Homa, Kyra Groeblichhoff

St. Louis County Police Department Crime Laboratory

The GlobalFiler™ Casework and Express kits were brought online at SLCPD in June of 2014. Since then approximately 1000 cases have been analyzed with the kit. The re-formulated kit was brought online in May of 2015. This presentation will cover general trends observed as well as casework examples. Some interesting cases will also be presented. In addition, preliminary Quantifiler® Trio validation data will be included.