

**The Validation of the OneStep ABACard[®] HemaTest[™]
for the Forensic Identification of Human Blood**

Sarah Mikolajczyk

The Northern Illinois Police Crime Laboratory, 1677 Old Deerfield Road, Highland Park, Illinois, 60034



The goal of this validation was to evaluate the use of the OneStep ABACard[®] HemaTest[™] for the qualitative detection of human blood within a forensics laboratory in lieu of the conventional cross-over electrophoretic technique. This rapid immunoassay detects human hemoglobin (hHb) which, in addition to the initial presumptive testing, can indicate the presence of human blood. Miscellaneous body fluid samples, species samples, various substrates with human blood, human blood samples processed with luminol, and non-probative casework samples were all evaluated with the OneStep ABACard[®] HemaTest[™]. The sensitivity of these kits were also established.