

OSAC Biology/DNA Scientific Area Committee Status Report

George Herrin, Jr
Chair, Biology/DNA Scientific Area Committee

The Biology/DNA Scientific Area Committee (SAC) of the Organization of Scientific Area Committees (OSAC), is responsible for developing and reviewing standards and guidelines related to the biological analysis of forensic evidence. The Biology/DNA SAC includes subcommittees on Biological Methods, Biological Data Interpretation and Reporting, and Wildlife Forensics. Documents covering a number of topics including training and education of analysts, terminology, several validation processes, and general standards for wildlife forensic laboratories have been in development since the formation of the OSAC in 2014. Several documents drafted within the OSAC process have been submitted to a Standards Development Organization (SDO) since late 2016. This presentation will provide updated information to the forensic community on the status of the standards and guidelines documents developed by the three subcommittees.