

## **UPDATE ON BIOLOGY/ DNA SCIENTIFIC AREA COMMITTEE ACTIVITIES**

George Herrin, Biology/DNA Scientific Area Committee

In 2014, the National Institute of Standards and Technology formed the Organization of Scientific Area Committees (OSAC) to produce consensus documentary standards and guidelines to improve quality and consistency of work in the forensic science community. The OSAC structure includes five Scientific Area Committees (SAC), one of which is the Biology/DNA SAC Scientific Area Committee. This Biology/DNA SAC includes three subcommittees: Biological Methods, Biological Data Interpretation and Reporting, and Wildlife Forensics. A short overview of the OSAC structure and goals, followed by a status update of the various action items from each of the three Biology/DNA subcommittees will be presented.

The presentation will be followed the next morning by a public 90 minute question and answer session in which members of the SAC will receive comments and/or respond to questions about the subcommittee initiatives.