American Society of Crime Laboratory Directors Laboratory Accreditation Board (ASCLD/LAB)



Promega 16th International Symposium on Human Identification

September 26-29, 2005 Grapevine, Texas

ASCLD/LAB Accreditation Programs Legacy and International

Presentation Focus:

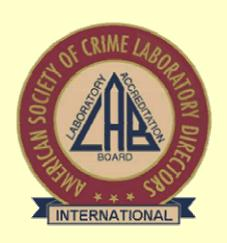
- Review Significant Changes from 2003 ASCLD/LAB Legacy Accreditation Manual to 2005 Accreditation Manual
- Review Timetable for Implementation of 2005 Legacy Manual
- Review Timetable for Updates to ASCLD/LAB-International Program Documents
- Additional Comments

Dual-Track Accreditation Program

- On April 1, 2004, ASCLD/LAB began offering accreditation under two separate programs Legacy and International.
- ASCLD/LAB continues to offer accreditation under the existing program Legacy Program. At this time the plan is to accept applications for accreditation under this program until April 1, 2014.
- ASCLD/LAB is also accepting applications for accreditation under the ASCLD/LAB-International Program which is an ISO 17025 based accreditation program.

ASCLD/LAB-International

- Based on ISO/IEC 17025
- Plus Supplemental Requirements based on Essential Requirements from ASCLD/LAB Legacy program and ILAC G 19 (Guidelines for Forensic Science Laboratories)
- Accreditation in the International program is for five years with annual Surveillance visits.
- Opens accrediting bodies up to 3rd party review.



301 Laboratories currently Accredited by ASCLD/LAB

as of 9-24-05

- 13 under ASCLD/LAB-International (significant number of applications)
- 289 under ASCLD/LAB Legacy Program
- 175 State
- 80 Local
- 23 Federal
- 14 Private
- 9 International

Significant Changes in the 2003 Manual

- Twenty-two revisions to the ASCLD/LAB Legacy Accreditation Manual were approved by the Delegate Assembly.
- The 2005 Manual will be published and distributed by **October 1, 2005**.
- ASCLD/LAB will accept applications under the 2005 manual immediately upon distribution.
- ASCLD/LAB will require all Legacy applications to be under the 2005 Manual on/or after **January** 1, 2006.

Summary of Relevant Revisions

- **Documentation of Compliance** must be maintained by the laboratory throughout the period of accreditation (five years).
- Standards for Individual Characteristic Databases (CODIS, NIBIN, AFIS) established.
- Defines Individual Characteristic

 Databases and Individual Characteristic

 Database Samples

- Annual Accreditation Audit Report due on anniversary of accreditation rather than April 1
- Reports of work must be generated.
- Reports must contain conclusions and opinions which address the purpose of the analysis.
- The significance of associations must be communicated clearly and properly qualified.

• Serology redefined as: A sub-discipline of biology, which is concerned with the identification of biological materials through the use of various tests.

• Evidence, physical defined as: Anything detectable by sensory, physical, chemical, optical or electronic means, including those things in a digital or multimedia form, that provides factual information about crime.

- A laboratory which must be accredited to satisfy legislation such as the "Justice for All Act" may seek accreditation in DNA only for a two-year period.
- The accreditation in a single discipline is for two years after which the laboratory must seek accreditation in all disciplines in which it provides services.

Updates to International Program Documents

- ISO/IEC 17025: 2005 was published by ISO in **May** 2005.
- The ASCLD/LAB-International Supplemental Requirements document will be revised to align it with the new ISO/IEC 17025 document and the 2005 Legacy Manual.
- Target publication date for the new ASCLD/LAB-International Supplemental Requirements document is January 1, 2006.
- Compliance with ISO/IEC 17025:2005 and the 2006 Supplemental Requirements will be **phased in during** 2006.

FBI Quality Assurance Standards

- In 2000, ASCLD/LAB signed an agreement with the FBI that for all ASCLD/LAB inspections involving DNA analysis, the FBI's DNA Audit Document will be used as a supplement to ASCLD/LAB standards for both the Legacy and International Programs.
- ISO/IEC 17025 ASCLD/LAB Supplement to 4.1.2 Note regarding FBI Audit Document use when DNA is part of laboratories scope.

Criteria File

Useful in preparing for and conducting the on-site.

<u>Legacy Program:</u>

- Must be completed by the laboratory and made available prior to on-site inspection
- A file (electronic or hard copy) which documents compliance with each of the criteria of the accreditation program
- Available in interactive format on the ASCLD/LAB website

Conformance File

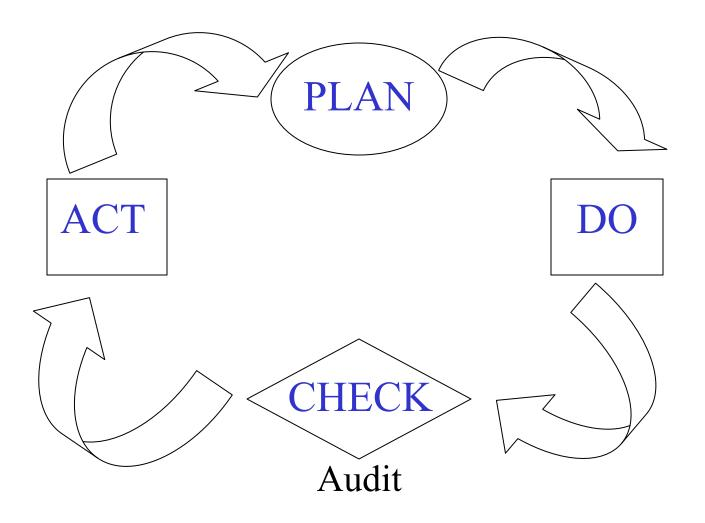
International Program:

Includes all applicable documentation of conformance with each requirement or the location of the policies and procedures that reflect conformance.

- Allows feedback to laboratory prior to setting visit
- Must be created by the laboratory and made available prior to on-site assessment.
- Part of the laboratories internal audit or GAP analysis which assists in identifying missing elements and readiness for the on-site assessment.
- Optional pre-assessment visit by lead assessor (not required) to conduct a gap-analysis.

Process Managing (Deming)

Laboratory Inspection/Assessment



ASCLD/LAB Audit Process

- Internal audits to assure a consistent high quality product.
- **Legacy** Draft inspection reports are audited post inspection.
- International Assessment reports are audited prior to Summation Conference with the laboratory.
- FBI DNA Quality Assurance Audit Document is audited by an FBI Trained Auditor. Newly added audit process.

Audit-Draft Inspection Report (Legacy)

- At the completion of the inspection, a **Draft inspection** report is formulated.
- The Draft report is <u>pre-decisional</u>, pending consideration of the Board of Directors.
- No report is left.
- The "draft" report to the ASCLD/LAB Executive Director.
- The report is forwarded to staff inspectors and Board Coordinator who will participate in an audit, for comment and questions.
- All comments received are compiled and discussed during an Audit Committee conference usually within a week.
- Audited Report sent to the laboratory.

Audited Summary Assessment Report (International)

- A Quality Review Panel will conduct an audit of the Summary Assessment Report, consisting of the CAR's for each non-conformity (level 1 and 2).
- Prior to the summation conference and presentation of the Report to the laboratory director.
- A copy is left with the laboratory.
- Full assessment report will be completed by the Lead Assessor within 15 days of the Summation Conference.

Sampling Plan

- ISO 17025 Clause 5.7.1
- "The laboratory shall have a sampling plan and procedures for sampling....."
- Based on appropriate statistical methods
- ASCLD/LAB Supplemental Note 3.

Further Information



- Rodney H. Andrus,
 Staff Inspector
- randrus@ascld-lab.org
- Website: ascld-lab.org