# Transitioning from the National QA Standards to ISO Accreditation

Kathleen A. Savage, Ph.D. Forensic Quality Services

## Forensic Quality Services

- Not-for-profit corporation
- Separated from NFSTC January 2004
- FQS
  - Training and education
  - Needs analysis
  - Laboratory Services
- FQS-I
  - ISO Accreditation

#### Board of Directors and Members

#### Board of Directors

- Ben Perillo, (Chairman), West Palm Beach Sheriff's Laboratory
- Bill Tilstone, (President)
- Arthur Eisenberg, University of North Texas
- Kathy Lee, Georgia Bureau of Investigation
- Sudhir Sinha, Reliagene
- Manuel Valadez, Texas DPS

#### Members

- Barry A J Fisher, Los Angeles County Sheriff's Laboratory
- Clifton Vander Ark, Arizona DPS
- All organizations accredited by FQS-I

#### ISO Accreditation?

- International Organization for Standardization
- Standards are documented agreements of technical specifications or other criteria to ensure that materials, products, processes and services are fit for their purpose
- Credit, phone and 'smart' cards derived from an ISO standard

#### FQS-I

- ISO 17025 Accreditation to forensic labs including specialist areas
  - Crime Labs
  - DNA Labs
  - Environmental
  - Racing Labs
- Auditor Selection
  - QAS and NDIS Requirements
- Accreditation Decision Making Process

## NACLA Recognition

- National Cooperation for Laboratory Accreditation
- Nonprofit corporation
  - Coordinate laboratory accreditation activities within US
  - Serve as US link to worldwide lab accreditation system
- Evaluate ABs and grant recognition to those found to meet ISO/IEC guide 58 (soon to be ISO 17011)
- ABs judged competent then invited to sign NACLA Mutual Recognition Agreement (consider each other's accreditation as technically equivalent)

## NACLA Recognition

- FQS-I recognized by NACLA for ISO 17025 accreditation within the scope of forensic science
- One of only 7 recognized ABs and the only one with scope of forensic science
- Accreditations provided through FQS-I since NACLA requires independence of operations from the parent company

### NDIS Procedures Board Recognition

FQS-I recognized by the NDIS Procedures Board as meeting the requirements to accredit laboratories participating in NDIS or seeking to participate in NDIS.

## NIJ Grant Funding

"Are accredited by a nonprofit professional association actively involved in forensic science that is nationally recognized within the forensic science community, such as the American Society of Crime Laboratory Directors/Laboratory Accreditation Board (ASCLD/LAB) or Forensic Quality Services (FQS)"

Forensic Casework DNA Backlog Reduction Grant

## FQS-I Program Changes

- Legacy 'DNA certification of compliance' no longer being offered
- ISO accreditation only
- Other labs transitioned with no significant difficulties
- Laboratories can choose whether to apply for accreditation in all areas, or just a few, or even start with one so that they can introduce the benefit of accreditation in a step-wise manner to the laboratory.

## Gaps between ISO and the DNA Standards

- Document control ISO 4.3
- Review of requests ISO 4.4
- Some aspects of purchasing ISO 4.6
- Service to the customer ISO 4.7
- Complaints ISO 4.8
- Control of non-conforming testing ISO 4.9
- Preventive action ISO 4.11
- Management review ISO 4.14
- Safety DNA Standard 16

#### Conclusions

 FQS-I has successfully transitioned several laboratories from DNA certifications

Not difficult

Not something to be afraid of!

#### For Further Information

William J. Tilstone, Ph.D.PresidentForensic Quality Services

wjt@forquality.org

727.541.2228

Kathleen A. Savage, Ph.D.
Business Manager
Forensic Quality Services

kas@forquality.org

727.541.2228

www.forquality.org