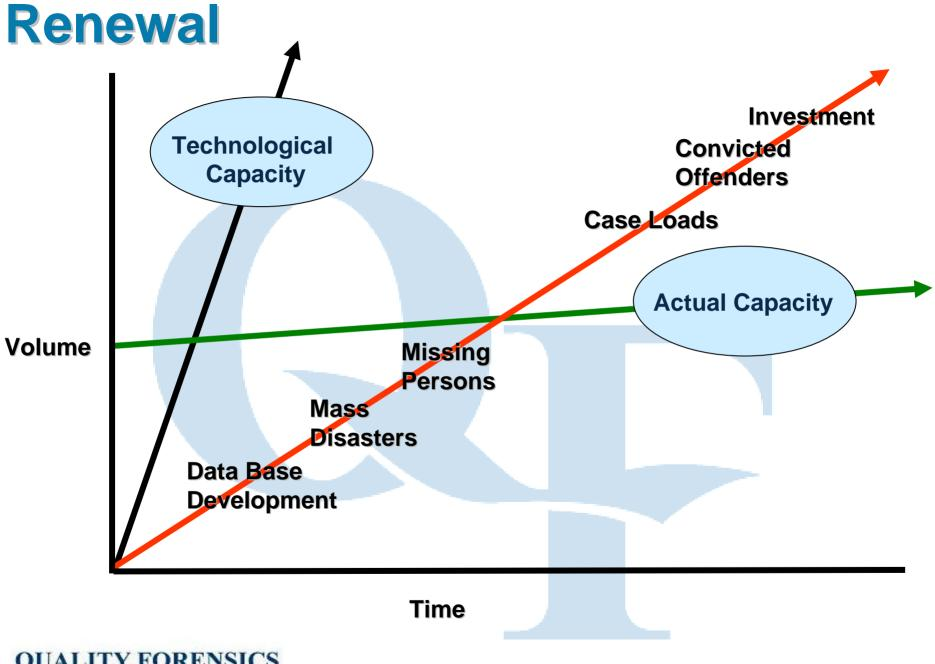
# Forensic Laboratory Renewal Process *FLRP*<sup>TM</sup>

Partnering Forensic Science with Management Science

Lucy Davis Houck

Quality Assurance Director









#### What's Different with FLRP?

In the past adopting new science, new technology or just making process changes was difficult if not impossible in forensic laboratories due to a number of systemic factors;

- Organizational resistance
- Lack of internal implementation expertise
- Lack of funding
- Use of expensive consulting services
- Operational workloads

FLRP is designed to address these and other issues by taking an approach that is bottom up and inclusive. The program uses modern facilitated processes that are designed to inexpensively and quickly assess, design, and transfer targeted professional implementation knowledge so that your organization can adapt and internalize the process renewals with minimal external assistance.





### Why the Approach Works

- Highly accelerated.
- Quantifiable results.
- Focus on high-leverage improvement areas.
- Driven by laboratory personnel from the ground up.
- High degree of implementation success.
- Creates and nourishes internal commitment.
- Focuses on the critical success factors.
- Respects the scientific process.
- Conforms to the Forensic Quality Assurance process.
- Maintains ongoing context that is tied to the actual process.







### Determine Where You Are



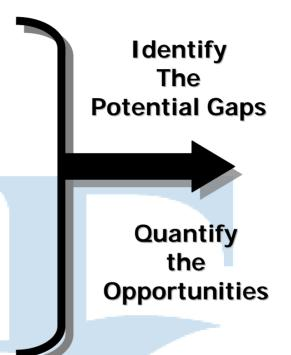
- Current State Process Review
- Current State Process Map
- Objectives
- Potential Quick Hits
- Enablers
- Inhibitors
- Information/Communications

**Technology Assessment** 

Recommendations and

rationale for Future State

**Design or Status Quo.** 



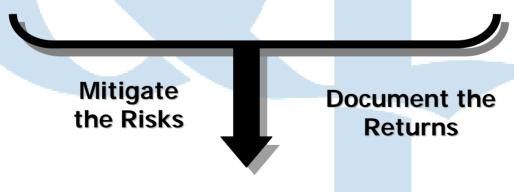






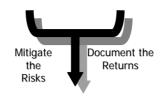
### **Determine Where You Want to Be**

- Validated Current State Process Map
- Future State Process Map
- Objectives
- Stretch Targets
- Assumptions
- Constraints
- Information/CommunicationsTechnology Requirements Definition
- Risk Management Plan









## Plan How Your Going to Get to Where You Want to Be

Create Partnerships

Create Consequences

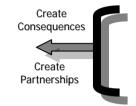
- Project Charter
- Objectives
- Assumptions
- Constraints
- Timeline
- Budget
- Accountabilities
- Project Schedule
- Project Budget
- Accountability Statements
- Status Accounting Process
- Management process
- Stakeholder Communication
   Plan
- Risk Management Plan





Manage the Change





## **Implement Your Plan**

- Status Reporting
  - Cost
  - Schedule
- Quality Review
- Best Practices Reviews
- Performance Measurement
- Change Management Reviews
- Risk Management Reviews
- Corrective Actioning



#### Determine Where You Are

#### Determine Where You Want to Be

- Current State Process Review
- Current State Process Map
- Objectives
- Potential Quick Hits
- Enablers
- Inhibitors
- Information/Communications Technology Assessment
- Recommendations and rationale for **Future State Design or Status Quo.**

Identify The **Potential Gaps** 



Quantify the Opportunities

- Validated Current State Process Map
- Future State Process Map
- Objectives
- Stretch Targets
- Assumptions
- Constraints
- Information/Communications **Technology Requirements Definition**
- Risk Management Plan

Measure the Results



Manage the Change

Implement Your Plan

- Status Reporting
  - •Cost
  - Schedule
  - Quality Review
  - Best Practices Reviews
  - Performance Measurement
  - Change Management Reviews
  - Risk Management Reviews
  - Corrective Actioning

**Document** the Returns



Mitigate the **Risks** 

Plan How Your Going to Get to Where You Want to Be

Create **Partnerships** 



Create Consequences

- Project Charter
  - Objectives, Assumptions, Constraints, Timeline, Budget
- Accountabilities
- Project Schedule
- Project Budget
- Accountability Statements
- Status Accounting Process
- Management process
- Stakeholder Communication Plan
- Risk Management Plan

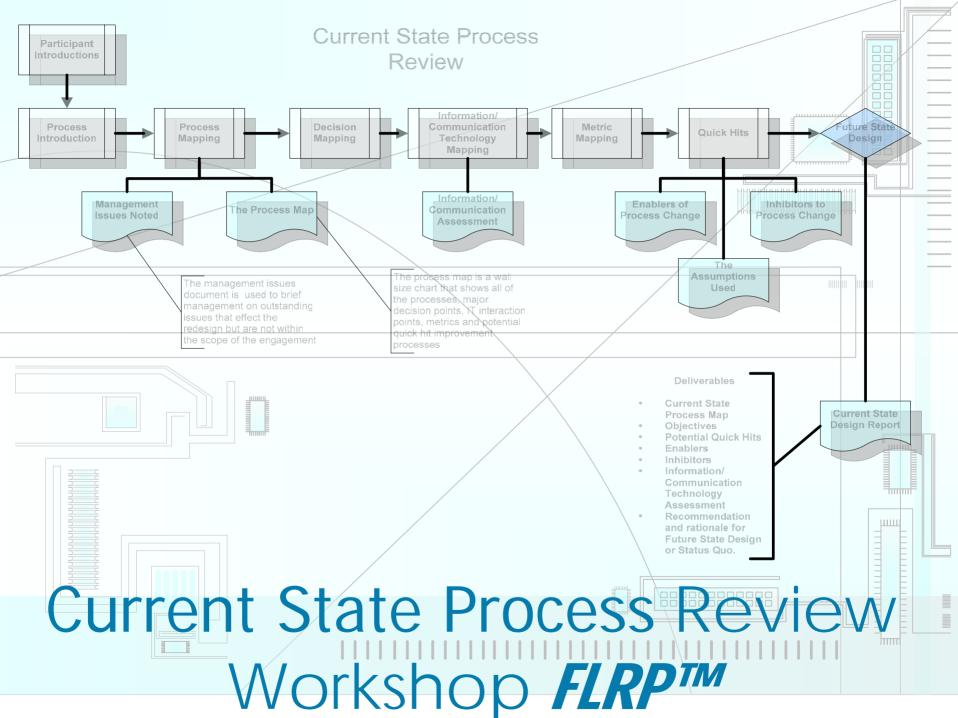
Proficiency Testing Services

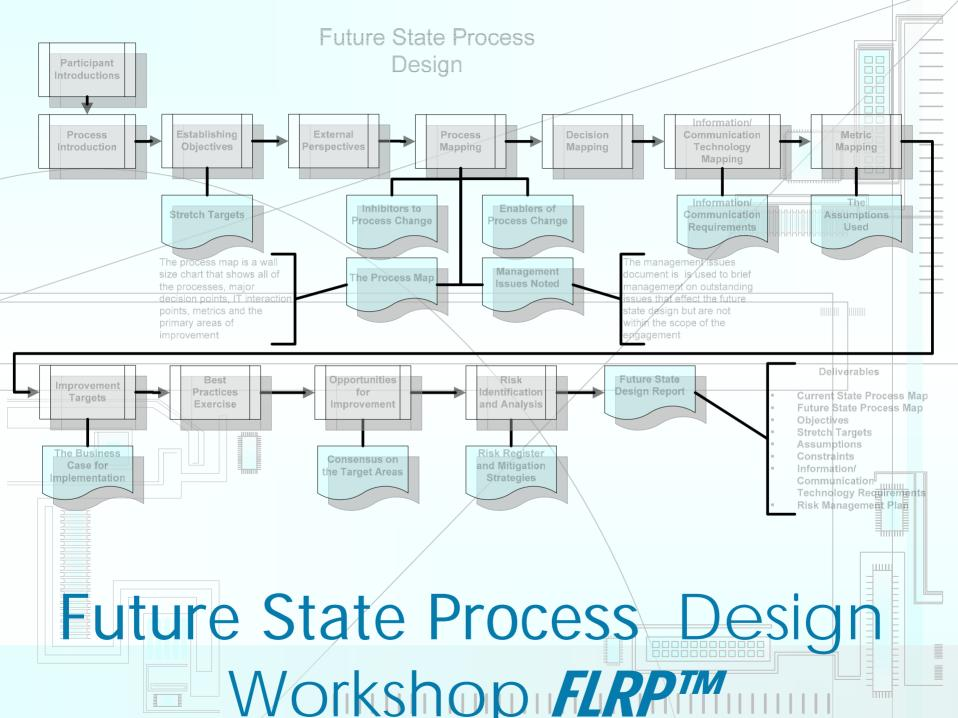


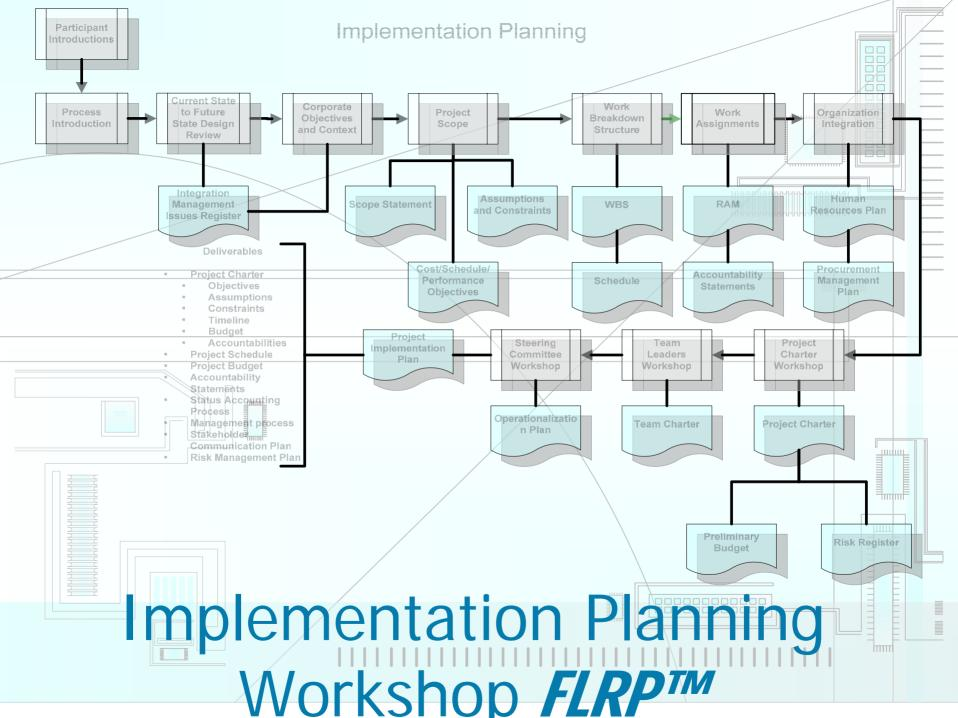
## **Four Primary Events**

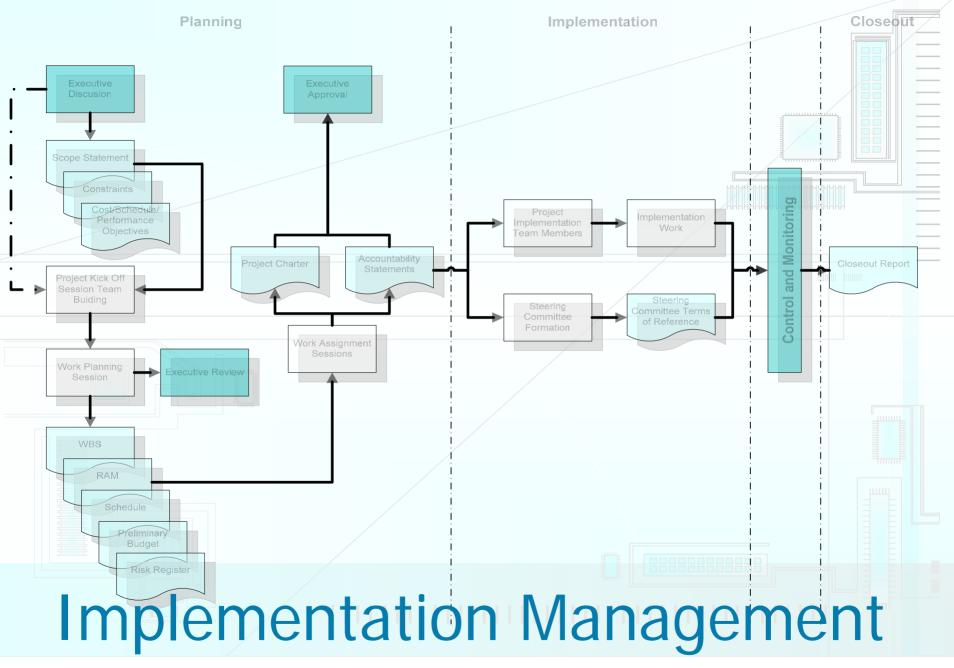
- Current State Review Workshop
- Future State Design Workshop
- Implementation Planning Workshop
- Implementation Management Process











Process FLRPTM



## **Knowledge Sources**

Determine Where You Are

Determine Where You Want to Be

Plan How Your Going to Get to Where You Want to Be

Implement Your Plan

Knowledge Workers
Process Specialists
Educational Work-Shops
OA Standards

I/C Technology Standards
Best Practices
QA Standards
Governance/Accountability

**Project Management Standards** 

Management Process Standards
Measurement Standards



## **Information Received**

- Fiscal Budget, Cash Flow, Return of Investment
- Schedule Timeline, Milestones,
   Dependencies, Tracking, Resource
   Usage
- Performance Improvement Metrics, Improvement Targets, Governance, Growth



## **Quantifiable Results**

- Rework The cost of non value added activities to correct things not done right the first time
- Throughput The number of items going through the process with no rework being performed
- Case and Sample Throughput Time Length of time it takes to get a case or a sample through the process
- Cost per Sample Processed Sample Processing cost divided by number of Samples analyzed by the process
- Investigations Aided Review the increase CODIS hits developed due to increased efficiency
- Customer Satisfaction The degree to which customers are satisfied with the performance of the process



# Acknowledgments

- We would like to thank:
  - Mike Kvasnik
  - Rick Mandy the Baintree Group



## More information

Lucy Davis Houck 304-685-6612

Lucy@QualityForensics.com

www.QualityForensics.com

