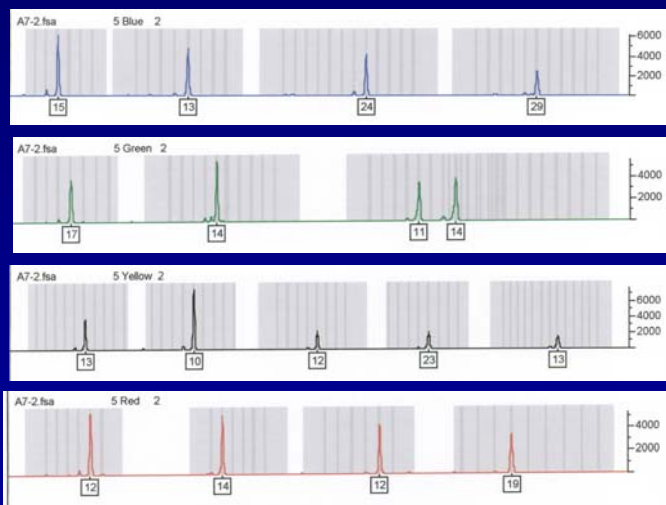


Demonstrating the Utility of the AmpFISTR Yfiler PCR Amplification Kit While Conducting Internal Validation Studies



Future Trends in
Forensic DNA Technology
Phoenix, AZ
March 22-23, 2006

Joseph Varlaro
Spencer Laboratories, Inc.

Presentation Overview

- Evaluation of the utility of Y-STRs (Yfiler)
- Development of validation strategy
- Highlights of internal validation studies
- Validation summary
- Casework example
- “Future trends”

Evaluation of Y-STRs

- **Major Forensic Advantage**
 - Useful with mixtures of DNA containing a small quantity of male DNA in a larger background of female DNA (i.e. oral assaults, fingernail clippings from female victim)
 - These types of mixtures often yield complex, ambiguous or inconclusive results with autosomal STRs
- **Major Forensic Limitation**
 - Sharing of Y-STR haplotype among paternal male relatives; moderate discriminatory power among unrelated men

Development of Validation Strategy

Establish analysis parameters to generate optimal data

Perform standard validation studies to confirm performance (Reproducibility, Sensitivity, Mixture)

Focus on analyzing mock casework samples to demonstrate utility of Y-STRs (Yfiler)

Sample Preparation for Simulated Oral Assault Evidence

- **Collected swabbings of neck and breasts of 2 females 30-60 minutes after each area was licked by male**
- **Performed DNA extraction according to standard organic extraction procedure**
- **Expecting high female: low male DNA mixture**

Analysis Strategy

Amplify 1 ng input DNA
in **Identifiler**

Evaluate Amelogenin
for the presence of
male DNA in mixture

Male DNA in Mixture
Amplify 0.75 ng male DNA
(if possible) in **Yfiler**

No male DNA (Female only)
Amplify *range* of input DNA
in **Yfiler**

Identifiler Results for Simulated Oral Assault Evidence

- **Single source male profile detected; no presence of mixture**
- **Yfiler analysis of samples would not demonstrate the utility of Y-STRs; samples were not amplified for Yfiler!**

Sample Preparation for Simulated Sexual Assault Evidence

- Collected oral swab from 2 females
- Dipped each oral swab in very dilute semen
- Performed DNA extraction (w/o differential) to simulate a sperm cell fraction with a high degree of epithelial cell carryover

Analysis Strategy

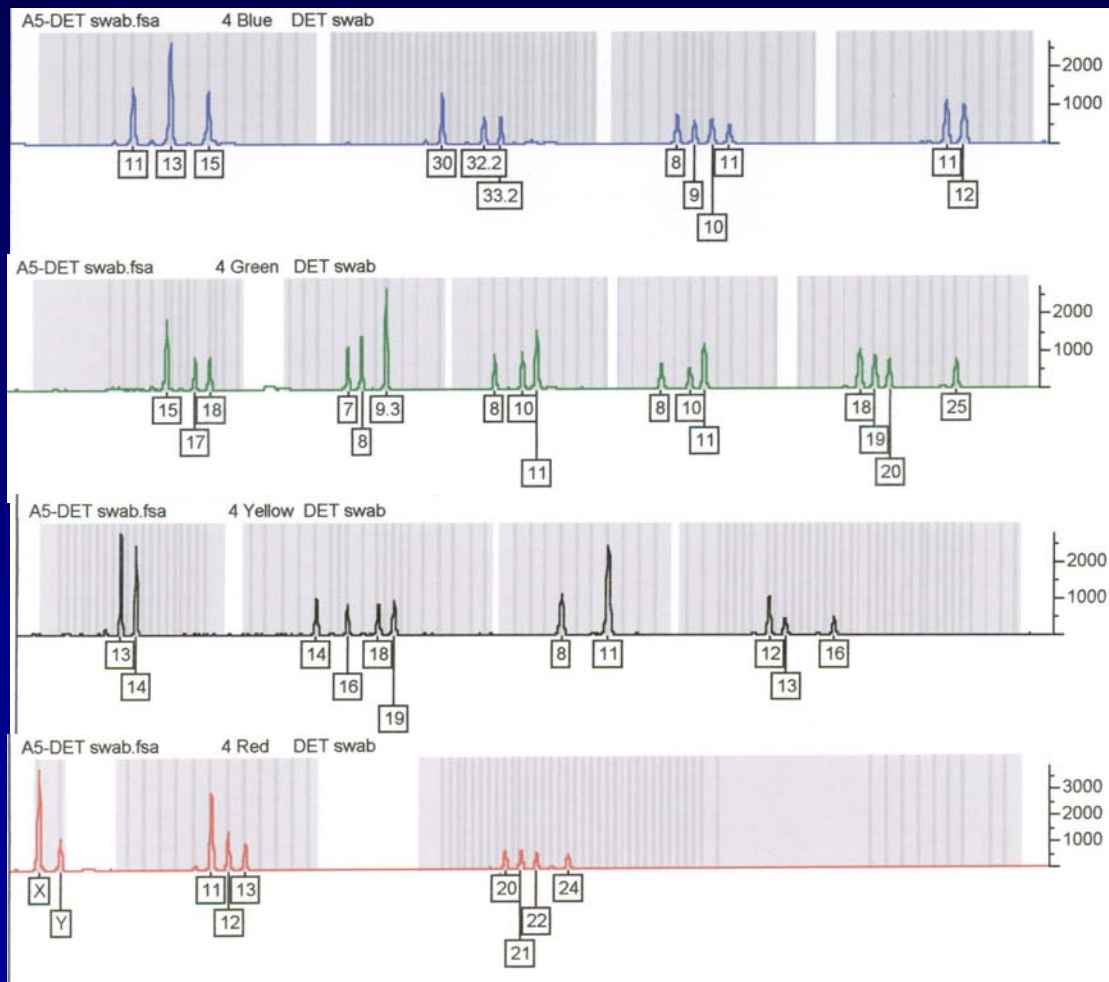
Amplify 1 ng input DNA
in **Identifiler**

Evaluate Amelogenin
for the presence of
male DNA in mixture

Male DNA in Mixture
Amplify 0.75 ng male DNA
(if possible) in **Yfiler**

No male DNA (Female only)
Amplify *range* of input DNA
in **Yfiler**

Identifiler Analysis of Swab 1



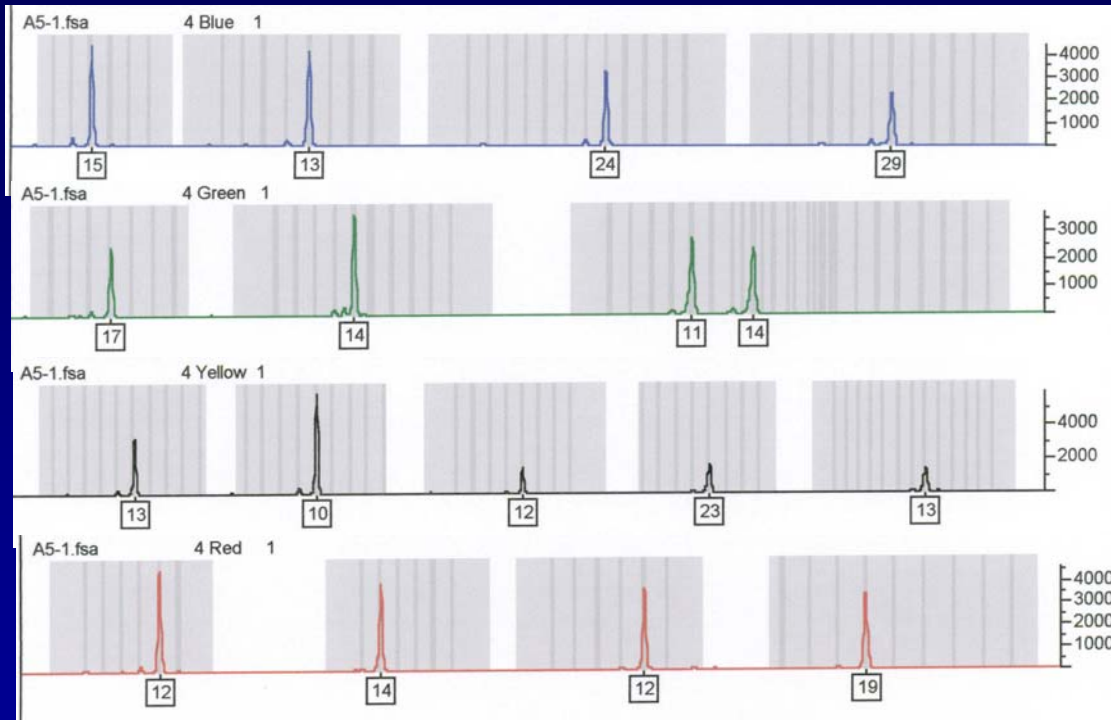
Identifiler Results for Swab 1

- **Mixture of male and female detected**
- **All alleles from male contributor were detected**

Yfiler Strategy for Swab 1

- Based on evaluation of Amelogenin, male contributor was present at approx. 50%
- TOTAL DNA concentration determined by QuantiBlot was multiplied by 50% to arrive at the approx. concentration of *male* DNA in the extract
- 0.75 ng male input DNA was targeted and amplified in Yfiler

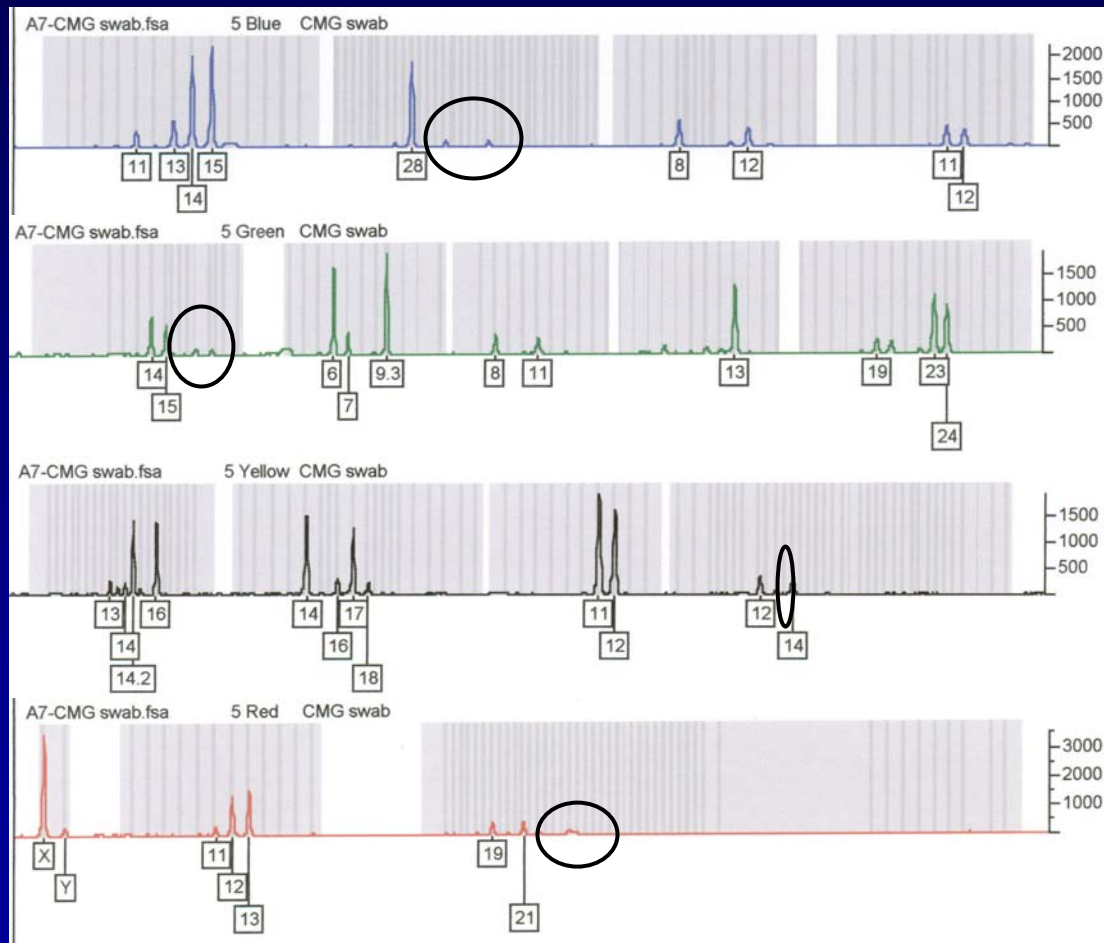
Yfiler Analysis of Swab 1



Yfiler Results for Swab 1

- **Single source Y haplotype detected**
- **Consistent with known male contributor (data from reference not shown)**
- **“Cleaned-up” the ID data; simplified data interpretation**

Identifiler Analysis of Swab 2



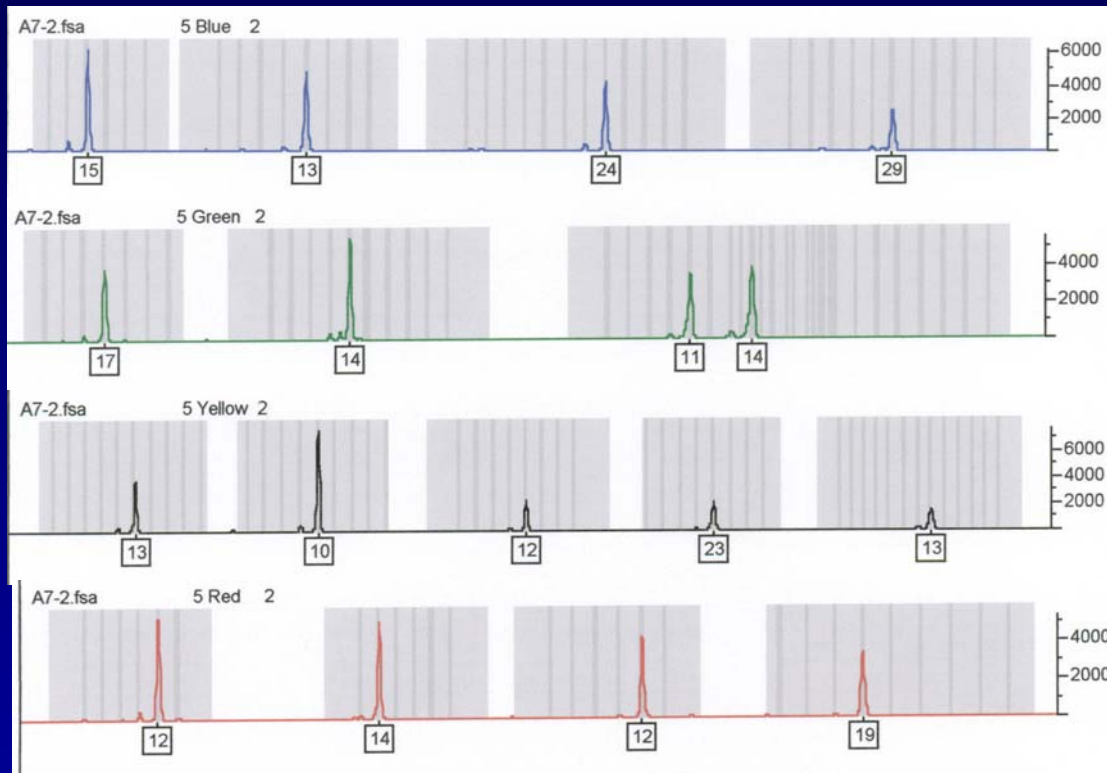
Identifiler Results for Swab 2

- **Mixture of male and female detected**
- **Numerous alleles from male contributor were not detected (partial result; subject to more complex interpretation)**

Yfiler Strategy for Swab 2

- Based on evaluation of Amelogenin, male contributor was present at approx. 20%
- TOTAL DNA concentration determined by QuantiBlot was multiplied by 20% to arrive at the approx. concentration of *male* DNA in the extract
- 0.75 ng male input DNA was targeted and amplified in Yfiler

Yfiler Analysis of Swab 2



Yfiler Results for Swab 2

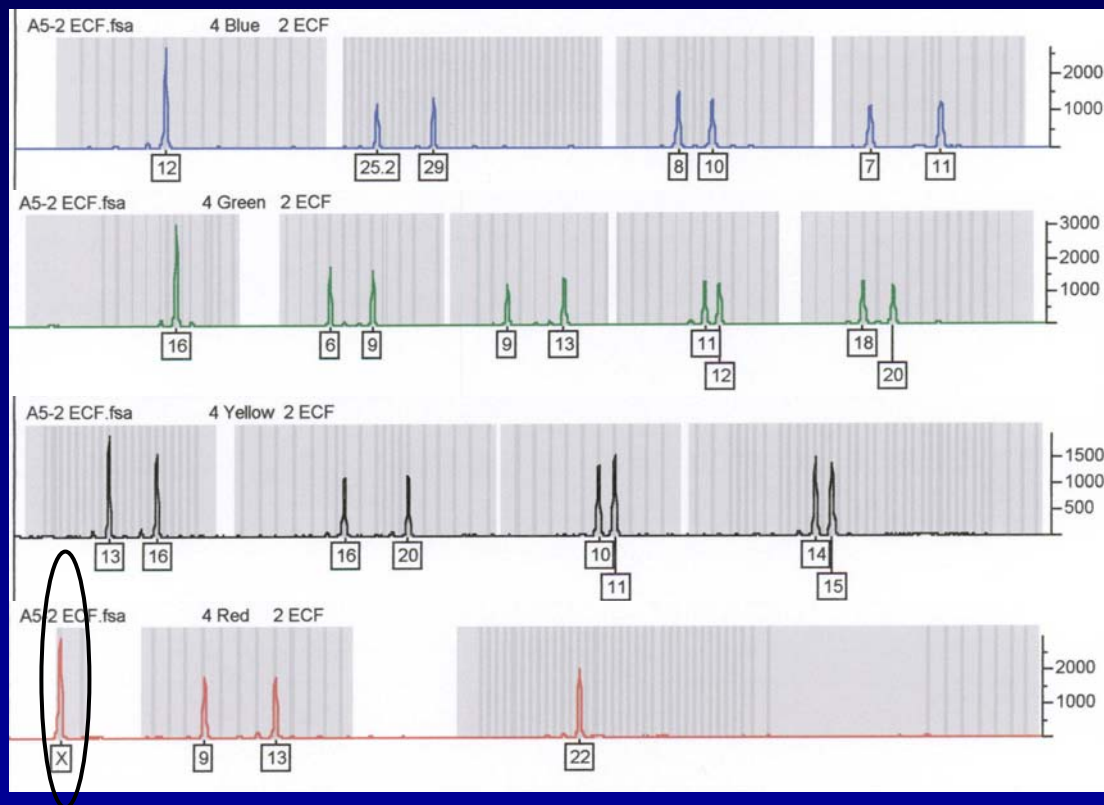
- **Single source Y haplotype detected**
- **Consistent with known male contributor (data from reference not shown)**
- **Simplified data interpretation; generated more *conclusive* result?**

Validation Summary

- **Yfiler is a highly effective tool with complex mixtures of male and female DNA**
- **Studies with simulated casework samples (although difficult to make) were most helpful in determining true utility of assay**
- **Yfiler is currently best used in conjunction with Identifiler**

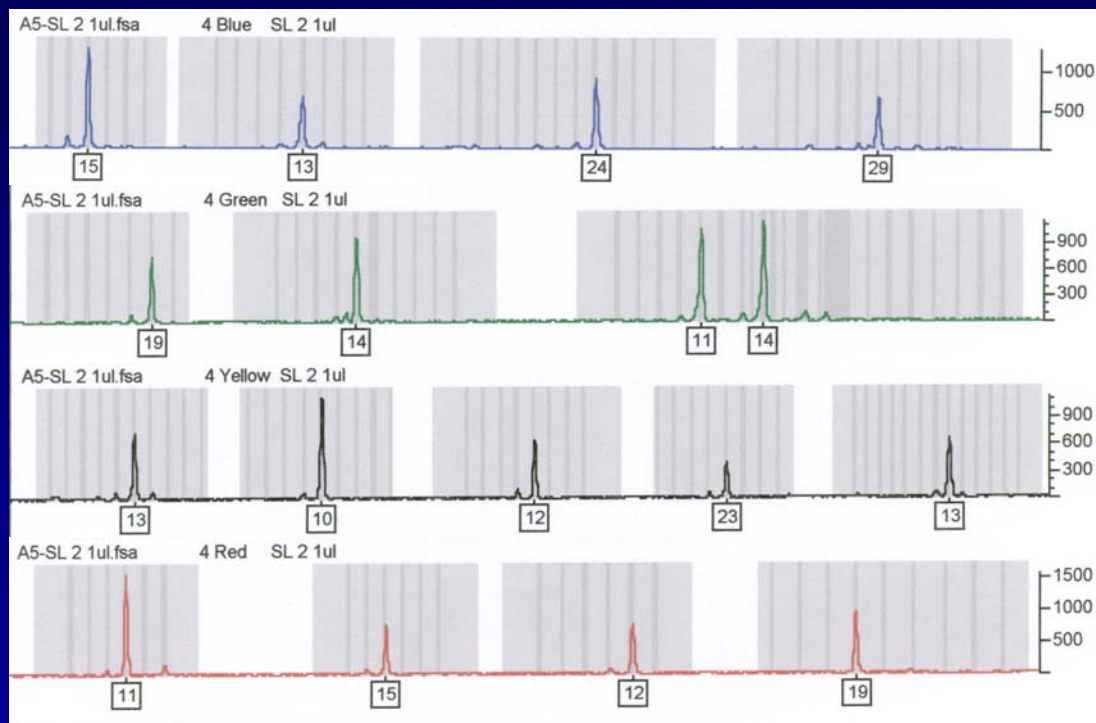
Identifiler Casework Sample

* Lip/breast swabbing from female victim of oral assault



* Since no male DNA detected, a range of input DNA will be amplified in Yfiler

Yfiler Analysis of Casework Sample



* Currently awaiting reference sample from suspect

“Future Trends”

- **Use of male-specific quantitation, along with total DNA quantitation (Quantifiler, Quantifiler Y), to determine ratio of male:female DNA**
- **Choose appropriate DNA typing system**
 - **Autosomal STRs**
 - **Y-STRs**

Acknowledgements

Spencer Laboratories, Inc.

Chantel M. Giamanco

Steven Todd

Applied Biosystems

Amie Ingold

Thank you very much!