

## Three Types of DNA Forensic Issues

- Single Source: DNA profile of the evidence sample providing indications of it being of a single source origin
- Mixture of DNA: Evidence sample DNA profile suggests it being a mixture of DNA from multiple (more than one) individuals
- Kinship Determination: Evidence sample DNA profile compared with that of one or more reference profiles is to be used to determine the validity of stated biological relatedness among individuals

• Interpretation of a result:

• 1. Non-match - exclusion

• 2. Inconclusive - no decision

• 3. Match - estimate frequency

### What is an Exclusion?

- Single Source: DNA profiles of the evidence and reference samples differ from each other at one or more loci; i.e., barring sample mix-up and/or false identity of samples, reference individual is not the source of DNA found in the evidence sample
- DNA Mixture: Reference DNA profile contains alleles (definitely) not observed in the evidence sample for one or more loci; i.e., reference individual is excluded as a part contributor of the mixture DNA of the evidence sample
- Kinship: Allele sharing among evidence and reference samples disagrees with the Mendelian rules of transmission of alleles with the stated relationship being tested

### What is an Inclusion?

- Single Source: DNA profiles of the evidence and reference samples are identical at each interpretable locus (also called DNA Match); i.e., reference individual may be the source of DNA in the evidence sample
- DNA Mixture: Alleles found in the reference sample are all present in the mixture; i.e., reference individual can not be excluded as a part contributor of DNA in the evidence sample
- Kinship: Allele sharing among evidence and reference samples is consistent with Mendelian rules of transmission of alleles with the stated relationship being tested; i.e., the stated biological relationship cannot be rejected

# When is the Observation at a Locus Inconclusive?

- Compromised nature of samples tested failed to definitively exclude or include reference individuals
- May occur for one or more loci, while other loci typed may lead to unequivocal definite inclusion/exclusion conclusions
- Caused often by DNA degradation (resulting in allele drop out), and/or low concentration of DNA (resulting in alleles with low peak height and/or area) for the evidence sample

#### Statistical Assessment of DNA Evidence

- Needed most frequently in the inclusionary events
- (Apparent) exclusionary cases may also be sometimes subjected to statistical assessment, particularly for kinship determination because of genetic events such as mutation, recombination, etc.
- Loci providing inconclusive results are often excluded from statistical considerations
- Even if one or more loci show inconclusive results, inclusionary observations of the other typed loci can be subjected to statistical assessment

## Exclusion vs Match

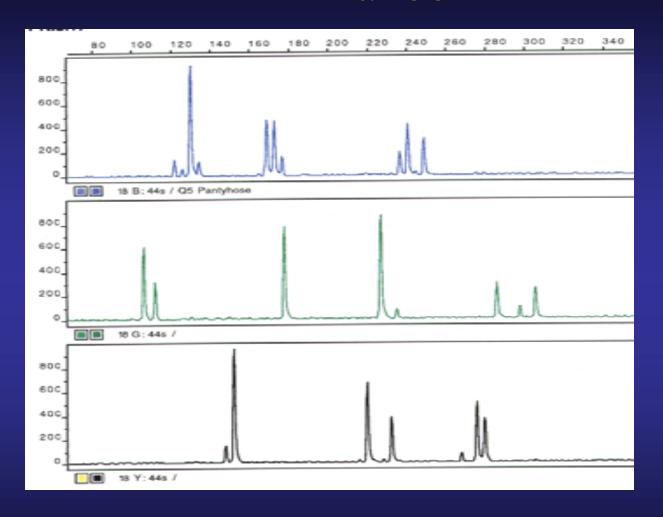
- Exclusion numbers are not needed
- Match requires a numerical estimate (weight of evidence)

# Statistical Analysis

### About "Q" sample

- "K" matches "Q"
- Who else could match "Q"
- partial profile, mixtures

# Mixtures



### What Constitutes a DNA Mixture?

• Presence of three or more alleles at several loci typed

Rationale: Typically a single individual displays one or two alleles

However, exceptions do occur – e.g., three allele profiles at the TPOX locus

### What Constitutes a DNA Mixture?

Notable imbalance in alleles at a locus

Rationale: Peak height/area difference, intensity difference, etc. caused by different contributions of amount of DNA from contributors of the sample.

# Common Mixture Interpretations

- Profiles of the tested persons can explain all alleles detected in the mixture sample
- Profiles of the tested persons do not explain the alleles detected in the mixture sample
- Mixture contains unexplained alleles beyond the ones present in the profiles of the known persons profiled
- Major and minor contributors
- Interpretable and uninterpretable

# Mixture Interpretation

Alleles considered first must meet interpretation guidelines for data analysis

- > RFU Threshold
- > Peak ratios/percent stutter
- > "Match" criteria

Calculations for mixed profiles are based on same statistical assumptions used for calculating single contributor profiles

### DAB Statistics

Alternate methods for assessing weight Rarely is there only one approach Philosophy, experience, legal system Practicality, available data, assumptions Simplistic approaches are acceptable Convey to fact finder Probability of Exclusion or LR Bayesian inferences NRC II Report

### Two common questions can be asked:

- How often would a random person be excluded as a contributor of the observed mixture? (Exclusion Probability)
- What statistical support is there for postulated hypotheses on the origin of the mixture? (Likelihood Ratio)

# Frequentist Approach of Statistical Assessment for DNA Mixture

- When the evidence mixture DNA profile fails to exclude a reference sample as a part contributor, and more commonly a set of reference samples together explains all alleles seen in the mixture, one or more of the following questions are answered:
- How often a random person would be excluded as a part contributor of the mixture sample? also called Exclusion Probability, the complement of which is the inclusion probability, giving the expected chance of Coincidental Inclusion
- (Note: This answer is based on the data on the evidence sample alone, without any consideration of the profiles of the reference samples)
- With a stipulation on the number of contributors, how often a random person's DNA, mixed with that of one or more of the reference persons, would provide a mixture profile as seen in the evidence sample, given that the reference persons are also part contributors of the DNA mixture
- (Note: This answer considers data on the profiles of evidence sample as well as those of the reference samples stipulated to be part contributors)

# Hardy - Weinberg Equilibrium

$$\begin{array}{c|cccc} A_1A_1 & A_1A_2 & A_2A_2 \\ \hline p_1^2 & 2p_1p_2 & p_2^2 \end{array}$$

$$freq(A_1) = p_1$$

$$freq(A_2) = p_2$$

	$A_1$	$A_2$
$A_1$	$p_1^2$ $A_1A_1$	$ \begin{array}{c} p_1p_2\\A_1A_2 \end{array} $
$A_2$	$ \begin{array}{c} p_1p_2\\A_1A_2 \end{array} $	$p_2^2$ $A_2A_2$

$$(p_1 + p_2)^2 = p_1^2 + 2p_1p_2 + p_2^2$$

D3S1358 = 16, 16 (homozygote)

Frequency of 16 allele = 0.3071

When same allele:

Genotype Frequency =  $p^2$  (for now!)

Genotype freq =  $0.3071 \times 0.3071 = 0.0943$ 

VWA = 15, 17 (heterozygote)

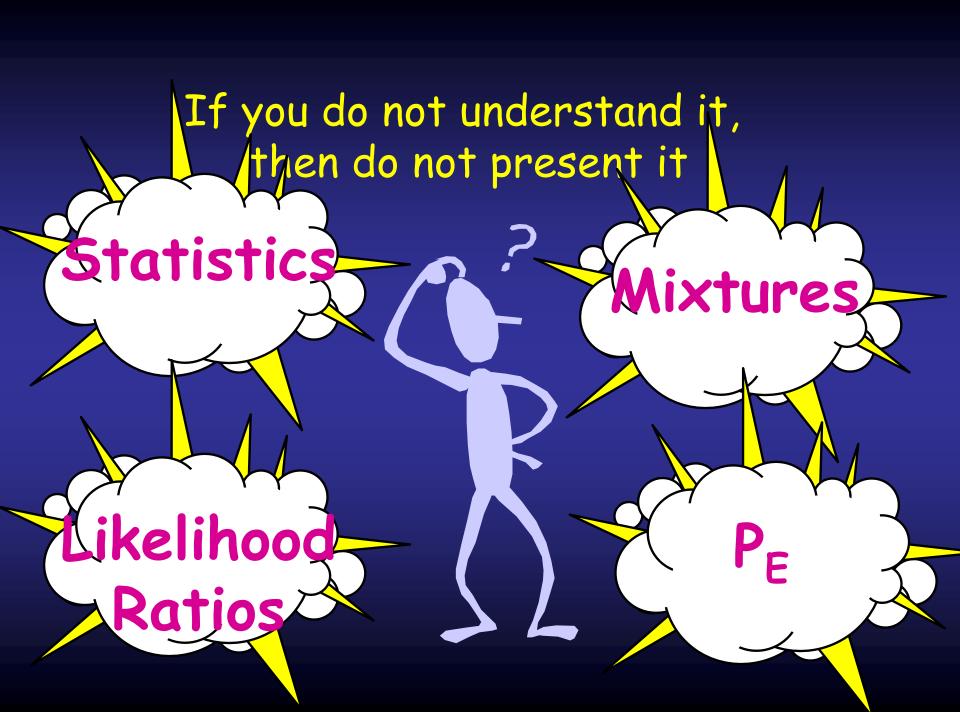
Frequency of 15 allele = 0.2361

Frequency of 17 allele = 0.1833

When heterozygous:

Frequency = 2 X allele 1 freq X allele 2 freq (2pq)

Genotype freq =  $2 \times 0.2361 \times 0.18331 = 0.0866$ 



# Probability of Exclusion

Not as powerful as LR

# **Exclusion Probability**

How often a random person would be excluded as a contributor of the observed DNA mixture?

# **Exclusion Probability**

 Power of the DNA testing panel for excluding potential noncontributing individuals from the profile

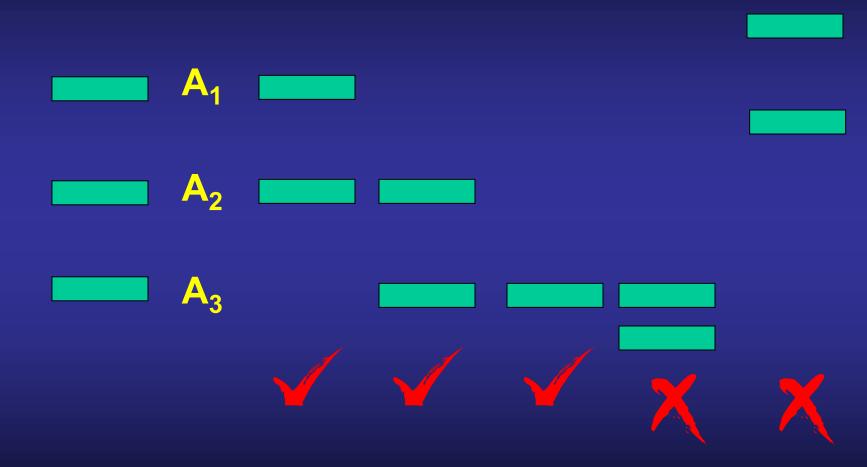
• Based on the alleles in the mixture...not the profiles of potential contributors

### Suspect Evidence

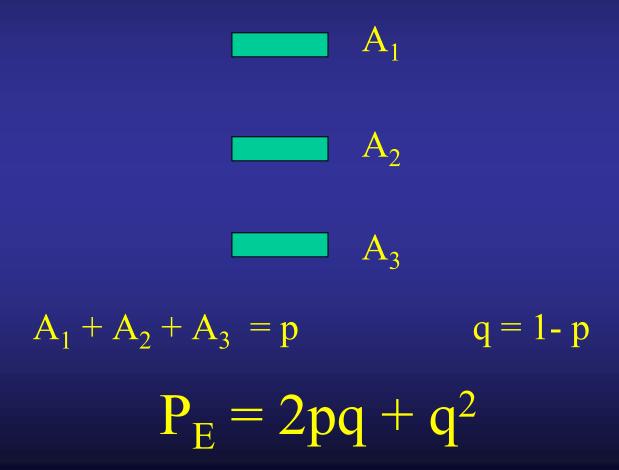


Three Allele Scenario

### Evidence



#### Evidence



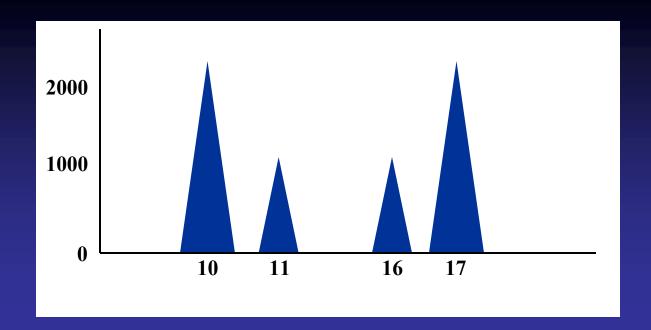
$$p_1 + p_2 + p_3 = p_c$$

$$p_1$$
 $p_2$ 
 $p_3$ 
 $(p_1 + p_2 + p_3)^2 = (p_c)^2$ 

$$p_1^2 + 2p_1p_2 + 2p_1p_3 + p_2^2 + 2p_2p_3 + p_3^2 = p_c^2$$

# Combined Probability of Exclusion

$$P_{Ec} = 1 - (1 - P_{E1}) (1 - P_{E2}) (1 - P_{En})$$



What is approach for P<sub>E</sub>?

Even though we can interpret profiles as a 10,17 and 11,16

### Likelihood Ratio

- compares the probabilities of a given observation under two different hypotheses
- a relatively subtle concept and LR computations should be explained with care
- a useful concept and used widely in statistics

### Likelihood Ratio

With two (mutually exclusive) hypotheses, say  $H_1$  and  $H_2$ , the likelihood ratio (LR) is the ratio of probabilities of observing the same data under  $H_1$  and  $H_2$ , giving

$$LR = Prob. (Data \mid H_1) / Prob. (Data \mid H_2).$$

#### Meaning of LR:

LR < 1: Data less well supported by  $H_I$ , compared with  $H_2$ 

LR = 1: Data equally well supported by  $H_1$  and  $H_2$ 

LR > 1: Data better supported by  $H_1$ , compared with  $H_2$ 

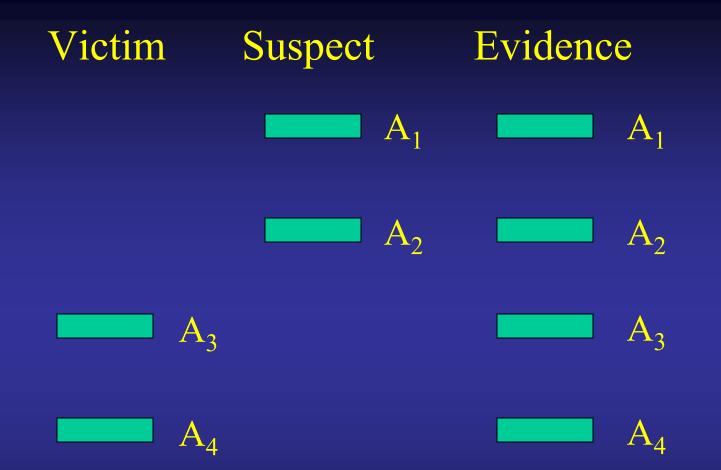
# Assumptions for LR

Independence No appreciable substructure All contributors of the same race All unrelated No allele dropout No intensity differences Defined hypotheses

# Mixtures Likelihood Ratio

- 1.  $H_p$  prosecution hypothesis
- 2.  $H_d$  defense hypothesis
- 3. Mutually exclusive
- 4. Probabilities ----  $LR = H_p/H_d$

Interpretation of mixture depends on circumstances of the case



Four Allele Scenario

#### Four Alleles

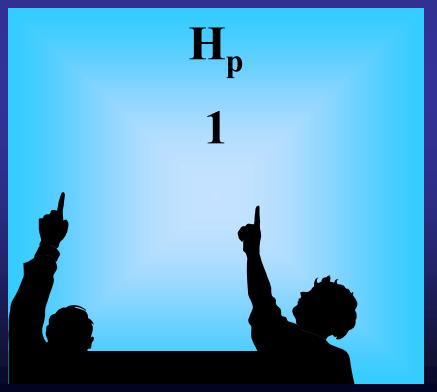
Two Match the victim - A<sub>3</sub> A<sub>4</sub>

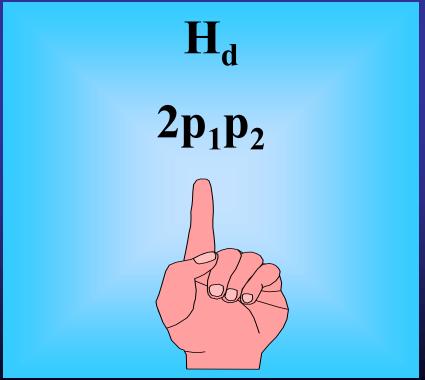
Two Match the suspect - A<sub>1</sub> A<sub>2</sub>

# Mutually Exclusive Hypotheses



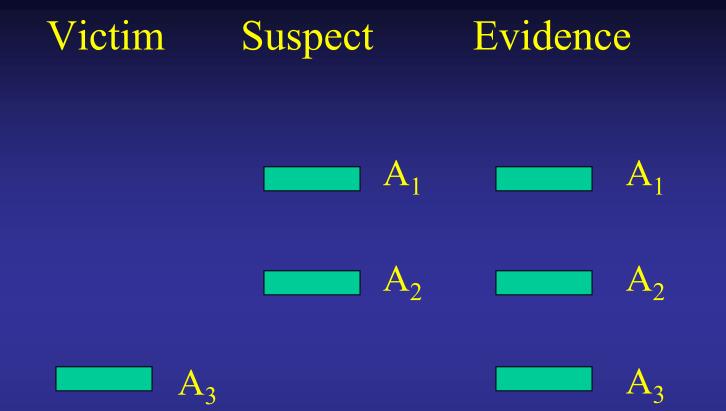
## Explain the evidence by the hypotheses





$$LR = H_p/H_d$$

 $1/2p_1p_2$ 



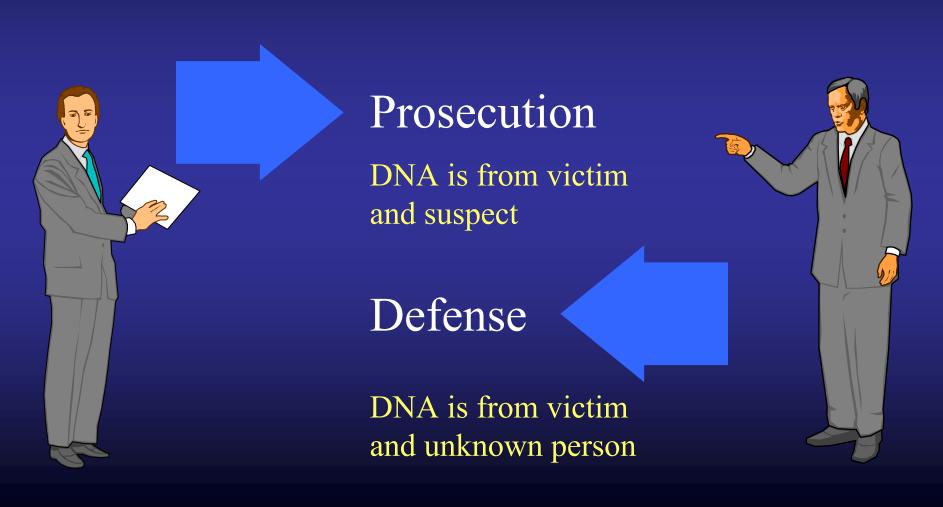
Three Allele Scenario

#### Three Alleles

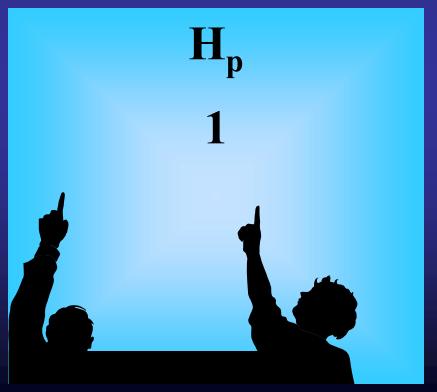
Victim is homozygote - A<sub>3</sub>A<sub>3</sub>

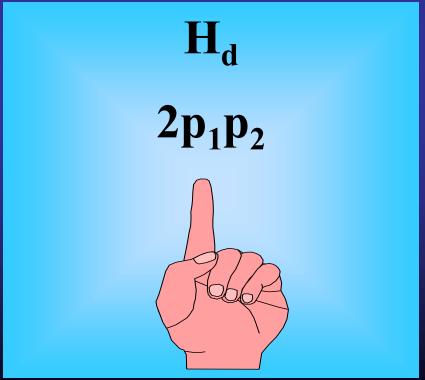
Two Match the suspect - A<sub>1</sub>A<sub>2</sub>

# Mutually Exclusive Hypotheses



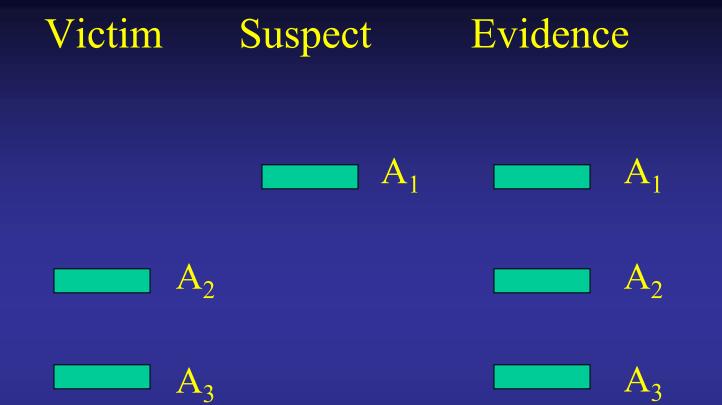
## Explain the evidence by the hypotheses





$$LR = H_p/H_d$$

 $1/2p_1p_2$ 



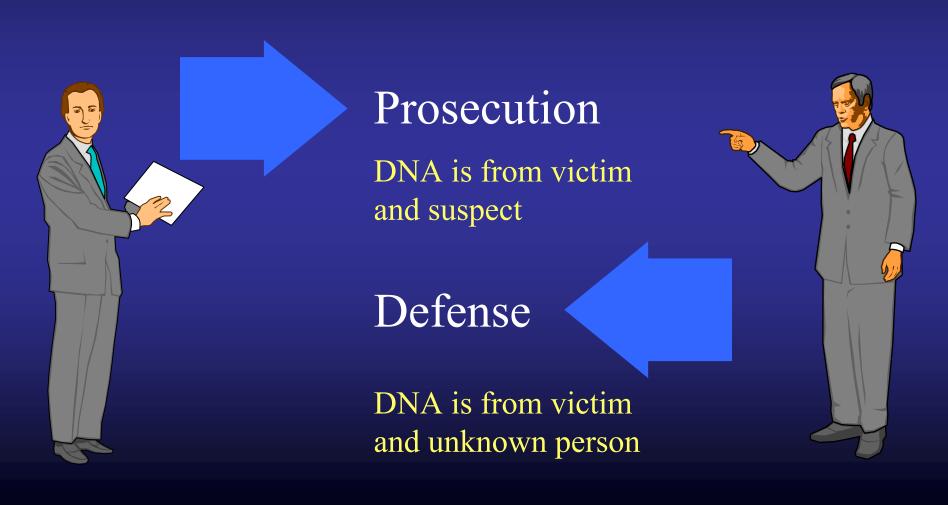
#### Three Allele Scenario

#### Three Alleles

Victim is heterozygote - A<sub>2</sub>A<sub>3</sub>

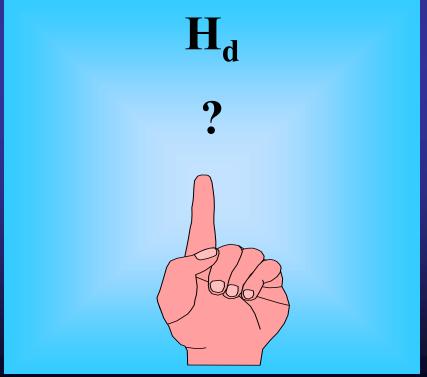
Suspect is homozygote- A<sub>1</sub>A<sub>1</sub>

# Mutually Exclusive Hypotheses



## Explain the evidence by the hypotheses





### Three possible genotypes can explain the evidence

Given that the Victim is heterozygote - A<sub>2</sub>A<sub>3</sub>

The possible genotypes to explain the evidence:

 $A_1A_1, A_1A_2, A_1A_3$ 

$$A_1A_1$$

$$A_1A_2$$

$$A_1A_3$$

$$p_1^2$$

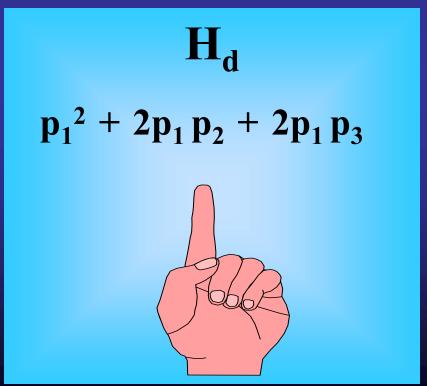
$$2p_1p_2$$

$$2p_1p_3$$

$$P_1^2 + 2p_1p_2 + 2p_1p_3$$

## Explain the evidence by the hypotheses





$$LR = H_p/H_d$$

$$1/p_1^2 + 2p_1p_2 + 2p_1p_3$$

Victim  $A_1A_1$ 

Suspect

Evidence

 $H_{\mathfrak{p}}$ 

 $H_d$ 

 $A_1A_1$ 

 $A_1A_1$ 

 $p_1^2$ 

 $A_1A_2$ 

 $A_1A_2$ 

 $A_1A_2$ 

 $p_1^2 + 2p_1 p_2 + p_2^2$ 

 $A_1A_1$ 

 $A_2A_2$ 

 $A_1A_2$ 

 $2p_1p_2 + p_2^2$ 

# HOW TO EXPRESS THE LIKELIHOOD RATIO COMPUTATIONS

- When making statements on statistical inferences remember express only the genetic data
- Not dealing with issues of "chance"
- To do so would involve Bayesian inferences which include Prior Probabilities...for which genetic data offer little assistance

Likelihood Ratio is how many times more likely it is to see this evidence under hypothesis #1 compared to hypothesis #2

$$LR = H_p/H_d$$

1.0/0.11 = 9.09

Compared with the prosecution's hypothesis  $(H_p)$ , the defense scenario  $(H_{d1})$  is

9-times less well-supported!

#### Alternatively

The observed mixture profile is 9-times more likely to occur under the scenario that it is a mixture of DNA from the victim and suspect, as opposed to the scenario that it originated from a mixture of DNA from the victim and an unrelated unknown person.

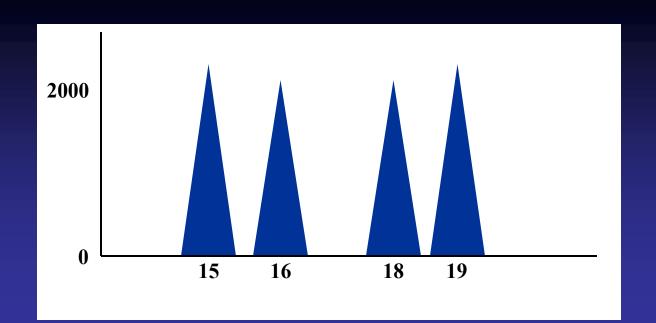
### Synthesis of Three Approaches of Statistical Assessment

- Frequency approach provides the probability of the observed DNA evidence (unconditional as well as conditional) under a given stipulated hypothesis
- Likelihood Ratio (LR) contrasts such probabilities for two mutually exclusive hypotheses
- In Bayesian approach, with the use of prior probability, LR is transformed to obtain the relative odds of one hypothesis against another given the DNA data of the evidence (and that from known persons tested)

#### Synthesis of Three Approaches (Contd.)

- The three approaches are built on one another, and hence, it is inaccurate to say one is wrong and the others are correct
- LR, without the transformation with the use of the prior probability, may be incorrectly interpreted as the answer of the Bayesian computation, but the numerator and denominator of LR can be stated with frequentist's interpretation to avoid the error of reverse conditioning
- The prior probability of the Bayesian approach generally comes from non-DNA evidence, and hence, their assumptions are untestable from DNA data

### Problems

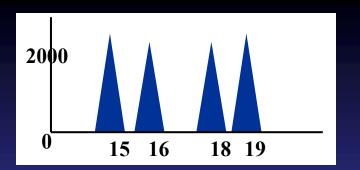


12 ... 0.010 13 ... 0.025 14 ... 0.070 15 ... 0.215 16 ... 0.230 17 ... 0.190 18 ... 0.150 19 ... 0.070 20 ... 0.030 21 ... 0.010

#### Calculate probability of exclusion for this mixture

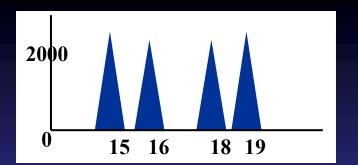
#### Calculate LR for this mixture

- a. Frame hypotheses
- b. Two contributors males
- c. Only one suspect type 16,19



Calculate probability of exclusion for this mixture

PE = 
$$1 - (0.215 + 0.230 + 0.150 + 0.070)^2$$
  
PE =  $1 - (.665)^2 = 0.558$  or  $55.8\%$ 



#### Calculate LR

### **Explain the evidence Prosecution**

Suspect – 16,19 Unknown – 15,18

 $\frac{1}{2\mathbf{p}_{15}\mathbf{p}_{18}}$ 

### What are possible genotype combinations? Defense

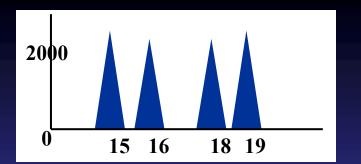
15,16 & 18,19 15,18 & 16,19

15,19 & 16,18

16,18 & 15,19

16,19 & 15,18

18,19 & 15,16

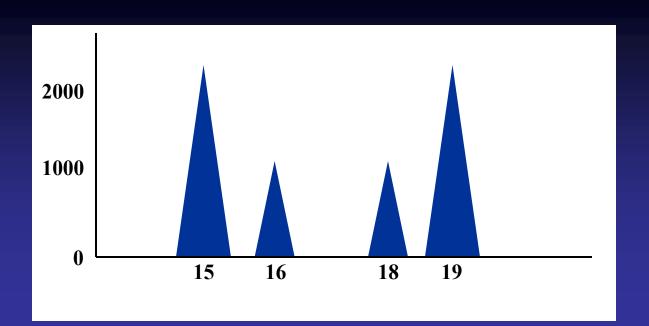


#### Calculate LR

$$LR = \frac{2p_{15}p_{18}}{24p_{15}p_{16}p_{18}p_{19}} = \frac{1}{12p_{16}p_{19}}$$

 $LR = 1/12 \times 0.230 \times 0.070 = 1/.1932 = 5.2$ 





```
12 ... 0.010

13 ... 0.025

14 ... 0.070

15 ... 0.215

16 ... 0.230

17 ... 0.190

18 ... 0.150

19 ... 0.070

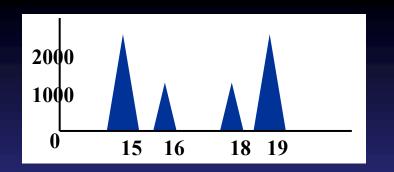
20 ... 0.030

21 ... 0.010
```

#### Calculate probability of exclusion for this mixture

#### Calculate LR for this mixture

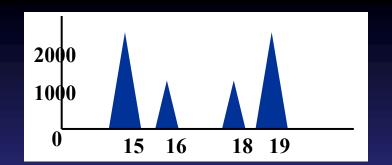
- a. Frame hypotheses
- b. Two contributors
- c. Victim type 16,18; suspect type 15,19



Calculate probability of exclusion for this mixture

PE = 
$$1 - (0.215 + 0.230 + 0.150 + 0.070)^2$$
  
PE =  $1 - (.665)^2 = 0.558$  or  $55.8\%$ 

However, these may be treated each as single source



#### Calculate LR Victim – type 16,18; suspect – type 15,19

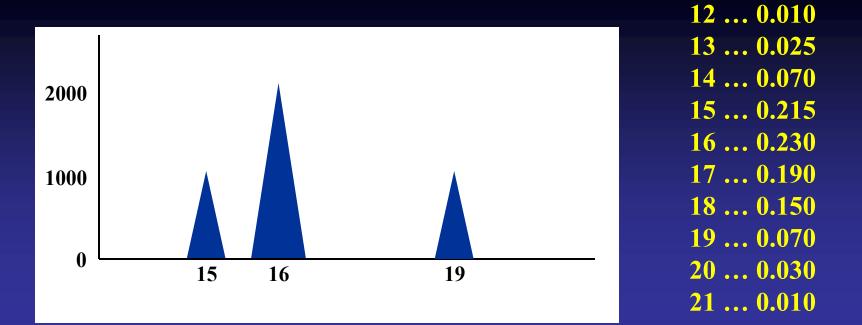
Explain the evidence

What are possible genotype combinations?

$$LR = \frac{1}{2 p_{15} p_{19}}$$

$$LR = 1/2 \times 0.215 \times 0.070 = 1/.0301 = 33.2$$

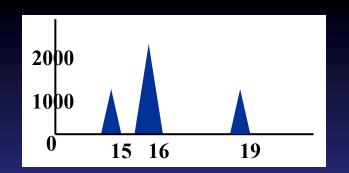




#### Calculate probability of exclusion for this mixture

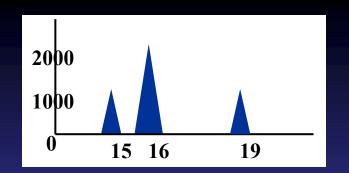
#### Calculate LR for this mixture

- a. Frame hypotheses
- b. Two contributors
- c. Victim type 15,16; suspect type 16,19



PE = 
$$1 - (0.215 + 0.230 + 0.070)^2$$

PE = 
$$1 - (.515)^2 = 0.735$$
 or  $73.5\%$ 



15 ... 0.215 16 ... 0.230 18 ... 0.150 19 ... 0.070

## Calculate LR Victim – type 15,16; suspect – type 16,19

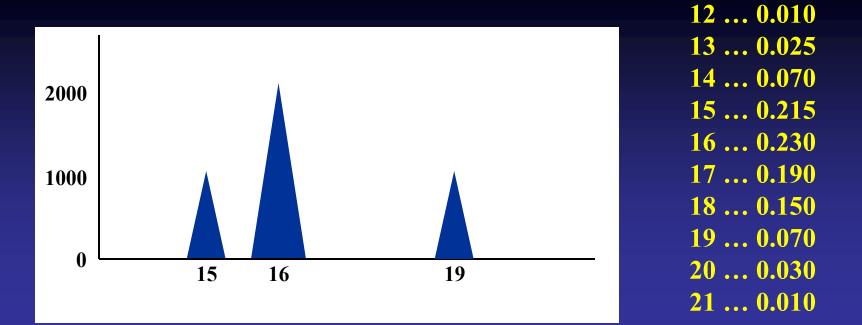
Explain the evidence

What are possible genotype combinations?

$$LR = \frac{1}{2 p_{16} p_{19}}$$

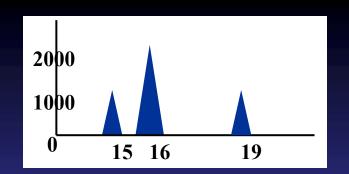
$$LR = 1/2 \times 0.230 \times 0.070 = 1/.0322 = 31.1$$





### Calculate LR for this mixture

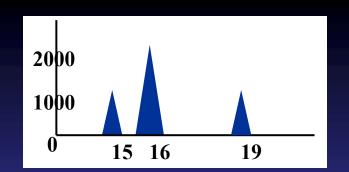
- a. Frame hypotheses
- b. Two contributors
- c. Victim type 15,16; suspect type 15,19



Victim – type 15,16; suspect – type 15,19

PE = 
$$1 - (0.215 + 0.230 + 0.070)^2$$

PE = 
$$1 - (.515)^2 = 0.735$$
 or  $73.5\%$ 



Calculate LR

15 ... 0.215 16 ... 0.230 18 ... 0.150 19 ... 0.070

Suppose defense save NA is from two unknown individuals

Victim 15,16,

Expla the evice e

Susp t – 16,19 Victio 15,16

spect-typed 65,99

What are possible genotype combinations?

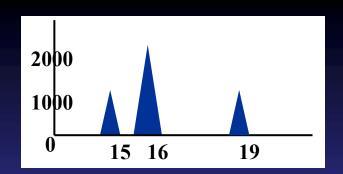
15,16<sub>u</sub> & 16,19<sub>u</sub> 15,19<sub>u</sub> & 16,16<sub>u</sub> 16,19<sub>u</sub> & 15,16<sub>u</sub> 16,16<sub>u</sub> & 15,19<sub>u</sub>

$$LR = \frac{1}{2(2p_{15}p_{16}x 2p_{16}p_{19} + 2p_{15}p_{19}x p_{16}^2)}$$

$$2 \times 2p_{15} p_{16} \times 2p_{16} p_{19}$$
  
 $2 \times 2p_{15} p_{19} \times p_{16}^{2}$ 

$$LR = 1/0.00954 = 104.8$$





Calculate LR

15 ... 0.215 16 ... 0.230 18 ... 0.150 19 ... 0.070

Suppose defense says the DNA is from two unknown individuals and one does not use quantitative data from electropherogram

Victim – type 15,16; suspect – type 16,19

Explain the evidence

**Suspect – 16,19** 

Victim – 15,16

1

What are possible genotype combinations?

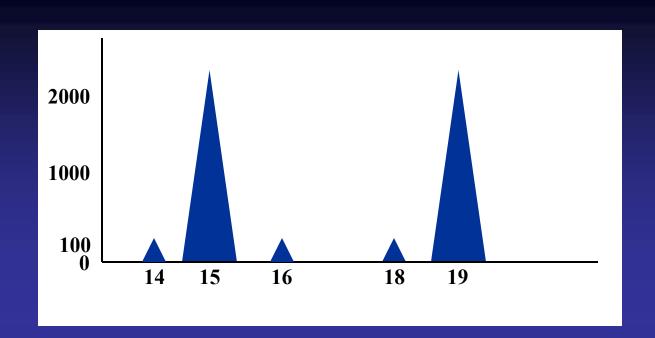
# $H_d$

<u>Unk 1</u>	Unk 2	
$A_{15}A_{16}$	$A_{15}A_{19}$	$2p_{15}p_{16} \times 2p_{15}p_{19}$
$A_{15}A_{16}$	$A_{16}A_{19}$	$2p_{15}p_{16} \times 2p_{16}p_{19}$
$A_{15}A_{16}$	$A_{19}A_{19}$	$2p_{15}p_{16} \times p_{19}^{2}$
$A_{15}A_{19}$	$A_{15}A_{16}$	$2p_{15}p_{19} \times 2p_{15}p_{16}$
$A_{15}A_{19}$	$A_{16}A_{19}$	$2p_{15}p_{19} \times 2p_{16}p_{19}$
$A_{15}A_{19}$	$A_{16}A_{16}$	$2p_{15}p_{19} \times p_{16}^{2}$
$A_{16}A_{19}$	$A_{15}A_{16}$	$2p_{16}p_{19} \times 2p_{15}p_{16}$
$A_{16}A_{19}$	$A_{15}A_{19}$	$2p_{16}p_{19} \times 2p_{15}p_{19}$
$A_{16}A_{19}$	$A_{15}A_{15}$	$2p_{16}p_{19} \times p_{15}^{2}$

$$A_{15}A_{15}$$
  $A_{16}A_{19}$   $p_{15}^2 \times 2p_{16}p_{19}$   
 $A_{16}A_{16}$   $A_{15}A_{19}$   $p_{16}^2 \times 2p_{15}p_{19}$   
 $A_{19}A_{19}$   $A_{15}A_{16}$   $p_{19}^2 \times 2p_{15}p_{16}$ 

$$12p_{15}p_{16}p_{19}(p_{15}+p_{16}+p_{19})$$





```
12 ... 0.010

13 ... 0.025

14 ... 0.070

15 ... 0.215

16 ... 0.230

17 ... 0.190

18 ... 0.150

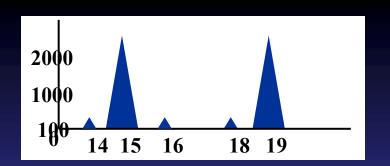
19 ... 0.070

20 ... 0.030

21 ... 0.010
```

### Calculate LR for this mixture

- a. Frame hypotheses
- b. Two contributors
- c. Victim type 15,19; suspect type 16,19



### Calculate PE

12 ... 0.010 13 ... 0.025 14 ... 0.070 15 ... 0.215 16 ... 0.230 17 ... 0.190 18 ... 0.150

Victim – type 15,19; suspect – type 16,19

19 ... 0.070

20 ... 0.030

21 ... 0.010

How to interpret?

$$PE = 1 - (p_{15} + p_{16} + p_{19})^{2}$$
or
$$PE = 1 - (p_{14} + p_{15} + p_{16} + p_{18} + p_{19})^{2}$$
or
$$PE = 1 - 2p_{16}$$
or

Treat as single source - p<sub>15</sub> p<sub>19</sub>

