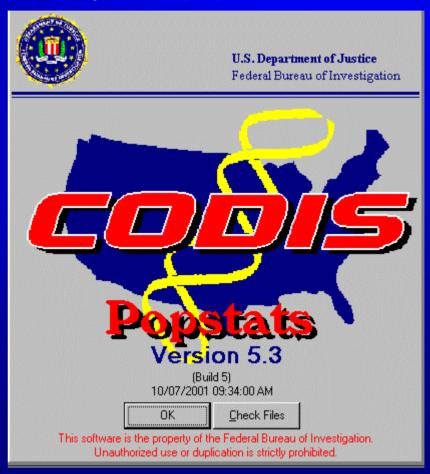
Popstats Unplugged

14th International Symposium on Human Identification

John V. Planz, Ph.D. UNT Health Science Center at Fort Worth



Forensic Statistics

From the ground up...

Why so much attention to statistics?

Exclusions don't require numbers

Matches do require statistics

Problem of verbal expression of

numbers

Transfer evidence

Laboratory result

- 1. Non-match exclusion
- 2. Inconclusive- no decision
- 3. Match estimate frequency

Statistical Analysis

Focus on the question being asked...

About "Q" sample

"K" matches "Q"

Who else could match "Q"

partial profile, mixtures

Match - estimate frequency of:

Match to forensic evidence

NOT suspect DNA profile

Who is in suspect population?

So, what are we really after?

Quantitative statement that expresses the rarity of the DNA profile

Estimate genotype frequency

- 1. Frequency at each locus

 Hardy-Weinberg Equilibrium
- 2. Frequency across all loci
 - Linkage Equilibrium

Terminology

Genetic marker variant = allele

DNA profile = genotype

Database = table that provides frequency of alleles in a population

Population = some assemblage of individuals based on some criteria for inclusion

Where Do We Get These Numbers?



POPULATION DATA and Statistics



DNA databases are needed for placing statistical weight on DNA profiles

vWA data (N=129)

	14	15	16	17	18	19	20	freq
14	9							75
15	3	0						6
16	19	1	1					46
17	23	1	14	9				72
18	6	0	3	10	4			31
19	6	1	7	3	2	2		23
20	0	0	0	3	2	0	0	5
								258

Because data are not available for every genotype possible,

We use allele frequencies instead of genotype frequencies to estimate rarity.

Estimate allele frequencies by "gene counting"

$$A_1$$
 n_1 $n_1 + n_2 + n_3 + n_4 + ... n_k = 2N$
 A_2 n_2
 A_3 n_3 $p_1 = n_1/2N$
 A_4 n_4 $p_2 = n_2/2N$
 \vdots \vdots $p_3 = n_3/2N$

Population database

Look up how often each allele occurs at the locus in a population (or populations)

AKA looking up the "allele" frequency

TECHNICAL NOTE

Bruce Budowle, ¹ Ph.D.; Tamyra R. Moretti, ¹ Ph.D.; Anne L. Baumstark, ¹ B.S.; Debra A. Defenbaugh, ¹ B.S.; and Kathleen M. Keys, ¹ B.S.

Population Data on the Thirteen CODIS Core Short Tandem Repeat Loci in African Americans, U.S. Caucasians, Hispanics, Bahamians, Jamaicans, and Trinidadians*

REFERENCE: Budowle B, Moretti TR, Baumstark AL, Defenbaugh DA, Keys KM. Population data on the thirteen CODIS core short tandem repeat loci in African Americans, U.S. Caucasians, Hispanics, Bahamians, Jamaicans, and Trinidadians. J Forensic Sci 1999;44(6):1277–1286.

markers are required, and all laboratories that contribute to the database should use the same genetic loci. Short tandem repeat (STR) loci are the most informative PCR-based genetic markers available to date for attempting to individualize biological material (2–5). The 13 STR loci CSF1PO, FGA, TH01, TPOX, vWA, D3S1358, D5S818,

CSF1PO.SWH - Notepad Edit Format Help Bin Range (alleles) Fraction Count <6̃− 0.0000 <6 6 0 6-0.0000 1 7-0.0024 0.0000 9-0.0072 10 -10 106 0.2536 10.3-10.3 0 0.0000 11 – 11 111 0.2656 9 12-164 12 0.3923 12.1-12.1 10 0.0000 11 27 13 -13 0.0646 12 14-14 0.0096 13 15-15 0.0048 14 >15- >15 0.0000 Totals 418 1.0001 Minimum allele frequency = 0.0120 Null allele frequency = 0.0120 Min Allele Frequency and Null Allele Frequency calculated as 5/2N where N = 209 single-allele patterns are entered twice in database Provided by B. Budowle, FBI Academy

FBI population Data

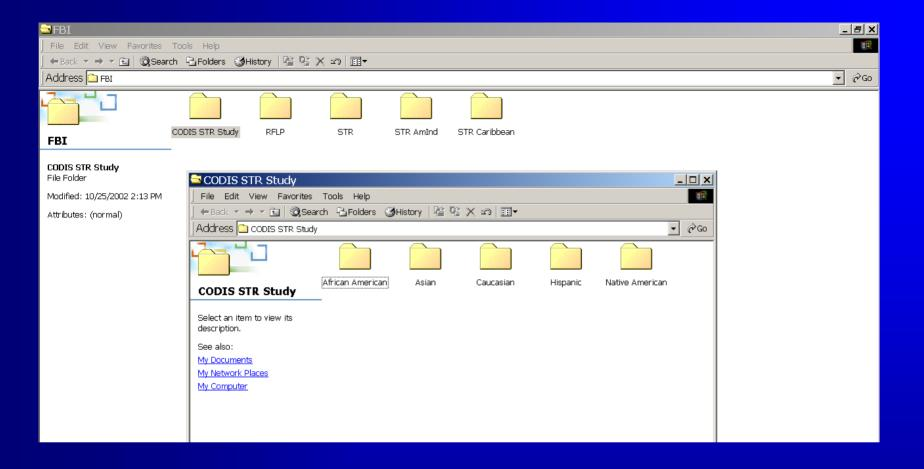
Bruce Budowle, ¹ Ph.D.; Brendan Shea, ² M.S.; Stephen Niezgoda, ² M.B.A.; and Ranajit Chakraborty, ³ Ph.D.

CODIS STR Loci Data from 41 Sample Populations*

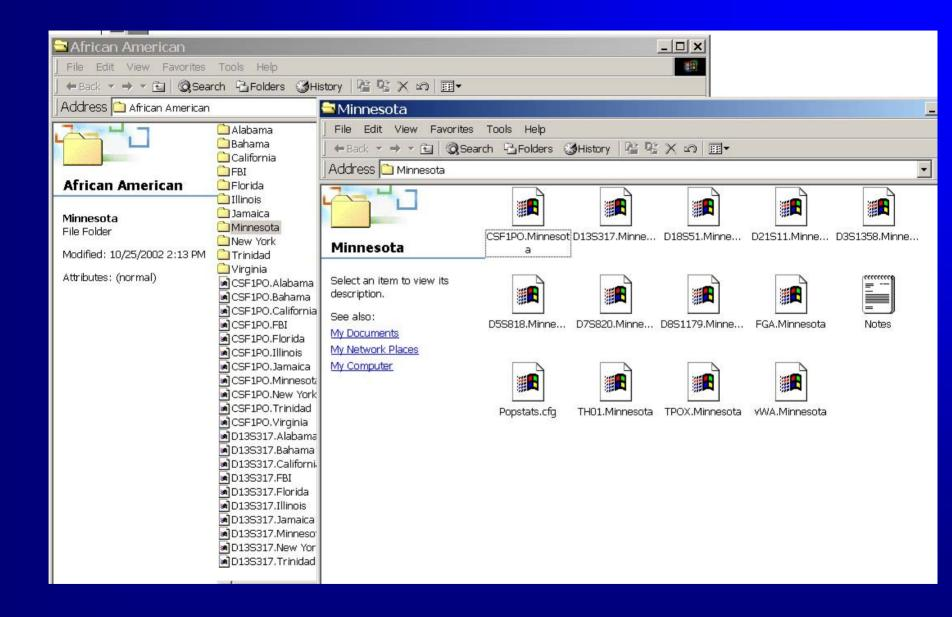
REFERENCE: Budowle B, Shea B, Niezgoda S, Chakraborty R. CODIS STR loci data from 41 sample populations. J Forensic Sci 2001:46;(3):453–489.

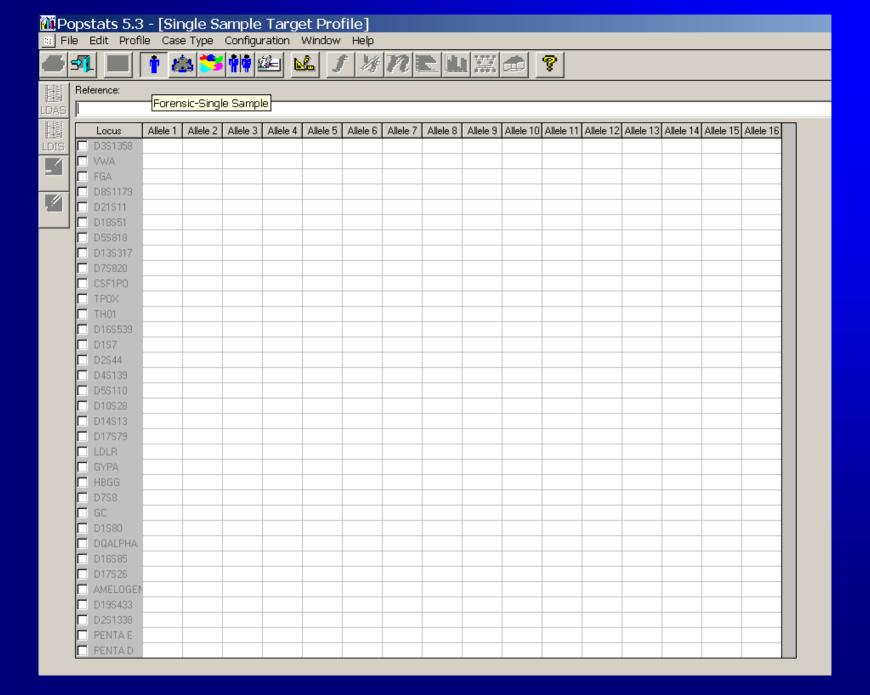
Materials and Methods

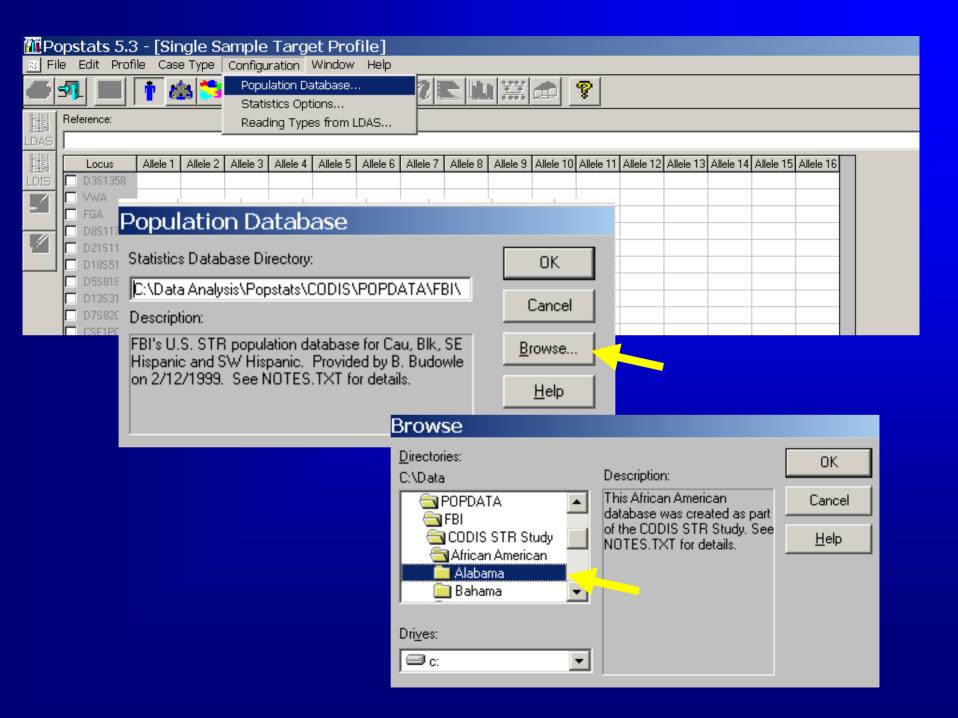
Samples



You can get other population databases to be used by Popstats if needed







OK

So now we have our population databases

How do they work for us in Popstats??

Anchor principle

Analysis of genetic makeup in individuals is based on the *Genotype* at the locus being queried

To remove "individual variation" so that we can focus on population-wide variation we must meld all the genotypes into a pool...separated as alleles

Alleles in populations The Hardy-Weinberg Theory

Basis: Allele frequencies are inherited in a Mendelian fashion and frequencies of occurrence follow a predictable pattern of probability

The Hardy-Weinberg principle states: that single-locus genotype frequencies after one generation of random mating can be represented by a binomial (with two alleles) or multinomial (with multiple alleles) function of the alleles frequencies.

This mathematical model serves as our *Null Hypothesis* when comparing populations, demes etc.

Hardy - Weinberg Equilibrium

$$A_1A_1$$
 A_1A_2 A_2A_2
 p_1^2 $2p_1p_2$ p_2^2 A_1 A_2

freq $(A_1) = p_1$ A_1 A_1A_2 A_1A_1 A_1A_2

freq $(A_2) = p_2$ A_2 A_1A_2 A_1A_2 A_2A_2

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

A Hardy-Weinberg Population

LARGE POPULATION
NO NATURAL SELECTION
NO MUTATION
NO IMMIGRATION / EMIGRATION
RANDOM MATING

Estimate genotype frequency:

1. Frequency at each locus

2. Frequency across all loci

Product Rule

Product Rule

The frequency of a multi-locus STR profile is the product of the genotype frequencies at the individual loci

 $f locus_1 \times f locus_2 \times f locus_n = f_{combined}$

				Profller F	Plus				
Item	D3S1358	vWA	FGA	D8S1179	D21S11	D18S51	D5S818	D13S317	D7S820
Q1	16,16	15,17	21,22	13,13	29,30	16,20	8,12	12,12	8,11

CoFller

Item	D3S1358	D16S539	TH01	TPOX	CSF1P0	D7S820
Q1	16,16	10,12	8,9.3	9,10	12,12	8,11

D3S1358 = 16, 16 (homozygote)

Frequency of 16 allele = ??

TABLE 1—Observed allele distributions (as %) for 13 STR loci in six population groups.

D3S1358	African American (N=210)	Bahamian (N=157)	Jamaican (N=194)	Trinidad (N=80)	Caucasian (N=203)	Hispanic (N=209)
<12	0.476	0.000	0.000	0.000	0.000	0.000
12	0.238	0.000	0.515	0.000	0.000	0.000
13	1.190	0.000	1.546	0.000	0.246	0.239
14	12.143	7.643	6.701	5.625	14.039	7.895
15	29.048	31.847	33.763	31.250	24.631	42.584
15.2	0.000	0.318	0.258	0.000	0.000	0.000
16	30.714	33.758	30.670	31.875	23.153	26.555
17	20.000	19.745	21.134	20.000	21.182	12.679
18	5.476	6.369	4.639	11.250	16.256	8.373
19	0.476	0.318	0.773	0.000	0.493	1.435
>19	0.238	0.000	0.000	0.000	0.000	0.239
Homozygosity (Obs.)	21.4%	25.5%	27.8%	16.3%	19.2%	26.3%
Homozygosity (Exp.)	23.5%	26.2%	25.8%	25.0%	20.3%	28.0%
(p)	0.482	0.838	0.513	0.070	0.691	0.595
Exact Test	0.797	0.758	0.270	0.222	0.084	0.333
PD	0.903	0.885	0.886	0.878	0.920	0.880
PE	0.543	0.499	0.508	0.511	0.589	0.492

D3S1358 = 16, 16 (homozygote)

Frequency of 16 allele = 0.3071

When same allele:

Frequency = genotype frequency (p²)
(for now!)

Genotype freq = $0.3071 \times 0.3071 = 0.0943$

				Profller F	Plus				
Item	D3S1358	vWA	FGA	D8S1179	D21S11	D18S51	D5S818	D13S317	D7S820
Q1	16,16	15,17	21,22	13,13	29,30	16,20	8,12	12,12	8,11

CoFller

Item	D3S1358	D16S539	TH01	TPOX	CSF1P0	D7S820
Q1	16,16	10,12	8,9.3	9,10	12,12	8,11

VWA = 15, 17 (heterozygote)

Frequency of 15 allele = ??

Frequency of 17 allele = ??

VWA	African American (N=180)	Bahamian (N=162)	Jamaican (N=244)	Trinidad (N=85)	Caucasian (N=196)	Hispanic (N=203)
11 13 14 15 16 17 18 19 20	0.278 0.556 6.667 23.611 26.944 18.333 13.611 7.222 2.778	0.926 2.778 6.173 15.123 26.235 20.679 18.210 7.099 2.778	0.410 0.820 7.377 22.746 29.098 18.238 13.115 5.328 2.254	0.588 0.588 8.824 14.118 29.412 26.471 13.529 4.706 1.765	0.000 0.510 10.204 11.224 20.153 26.276 22.194 8.418 1.020	0.246 0.000 6.158 7.635 35.961 22.167 19.458 7.143 1.232
Homozygosity (Obs.)	0.000	0.000	0.615	0.000	0.000	0.000
Homozygosity (Obs.) Homozygosity (Exp.) (p)	18.9% 0.014	17.5% 17.6% 0.928	19.4% 0.557	20.0% 20.991	18.7% 0.179	22.9% 0.564
Exact Test	0.328	0.790	0.655	0.229	0.063	0.928
PD PE	0.926 0.624	0.942 0.648	0.933 0.617	0.917 0.602	0.932 0.625	0.914 0.563

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

Overall profile frequency =

Frequency D3S1358 X Frequency vWA

 $0.0943 \times 0.0866 = 0.00817$

This is basically what Popstats does for us in it's simplest task

Steps – Single Sample Target Profile

- enter alleles of target profile
- look up allele frequencies at all loci for all populations
- determine if homozygous or heterozygous at each locus
- calculate genotype frequency at each locus
- calculate profile frequency with product rule

But this doesn't address all of the issues!

What if...

We encounter alleles not represented in the population database...

...or alleles that are extremely rare in the database???

Minimum allele frequency

The first NRC report proposed a minimum allele frequency based on NO empirical data and without any statistical basis!

10 % or 0.1

What...you are surprised??

Ceiling Principle

Minimum allele frequency

Weir, B.S. 1992. minfreq = 1 - $\alpha^{1/2N}$

Budowle, B., K. Monson, R. Chakraborty, 1996. minfreq = $1 - [1 - (1 - \alpha)^{1/c}]^{1/2N}$

NRC II, 1996. minfreq = 5/2N

Minimum allele frequency

This method requires a minimum of 5 copies of an allele before the allele frequency can be used for calculation of genotype frequency

Total number of alleles at locus

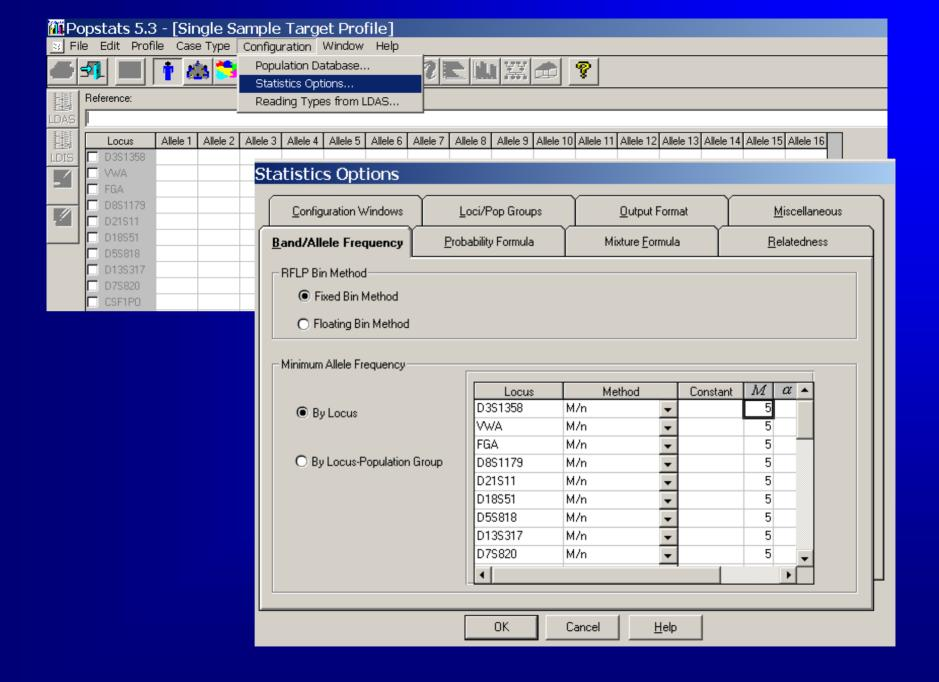
For the 13 allele at vWA: Actual Freq = 2 / 392 = 0.0051Minimal Freq = 5 / 392 = 0.0128

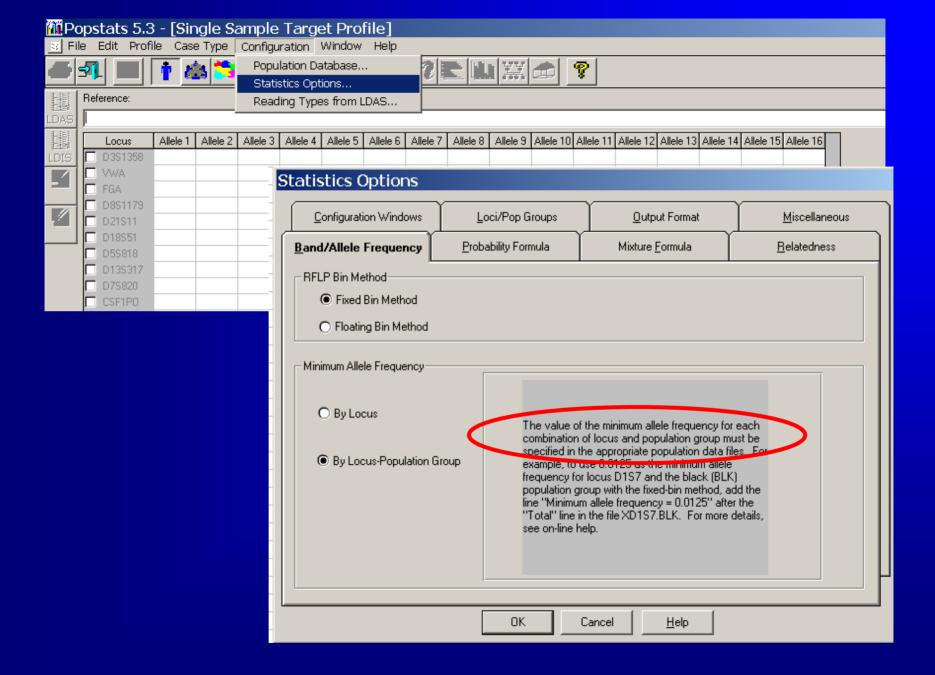
Conservatism & also addresses some substructure effects

This estimate is strictly driven by database size:

<u>N</u>	min allele freq					
100	2.50 % (0.025)					
150	1.67 % (0.0167)					
200	1.25 % (0.0125)					
250	1.00 % (0.01)					
300	0.83 % (0.0083)					

Where N is the number of individuals in database





CSF1PO.SWH - Notepad

File	Edit	Format F	łelp		
File	Edit Bin 1 2 3 4 5 6 7 8 9		(alleles) <6 6 7 8 9 10 10.3 11	Count 0 0 1 0 3 106 0 111 164 0	Fraction 0.0000 0.0000 0.0024 0.0000 0.0072 0.2536 0.0000 0.2656 0.3923 0.0000
	10 11 12 13 14	12.1- 13- 14- 15- >15-	13 14 15 >15	27 4 2 0	0.0000 0.0646 0.0096 0.0048 0.0000
			imum allele	418 e frequency = equency = 0.	

Min Allele Frequency and Null Allele Frequency calculated as 5/2N where N = 209

Note: single-allele patterns are entered twice in database

Provided by B. Budowle, FBI Academy

So the only other real thing left to consider regarding the NRC concerns is population subdivision.

Population Structure

Racial, ethnic subgroups

Excess of homozygotes

What is "theta" θ

Why modify just homozygous calculation?

NRC Formula 4.1 vs 4.4 vs 4.10

Population Subdivision

We've always surmised...

Racial / ethnic group composed of distinct sub-groups within the sample population

Only a concern if sub-groups differ substantially at allele frequencies at the loci

Human Genetic Variation

between populations within racial groups ...

between racial groups

within populations within racial groups

- Barbujani, Magagni, Minch, Cavalli-Sforza. 1997. An apportionment of human DNA diversity. *PNAS* 94:4516-4519.

Problems created by population subdivision

Genotype frequencies calculated from population average allele

frequencies COUIC lead to:

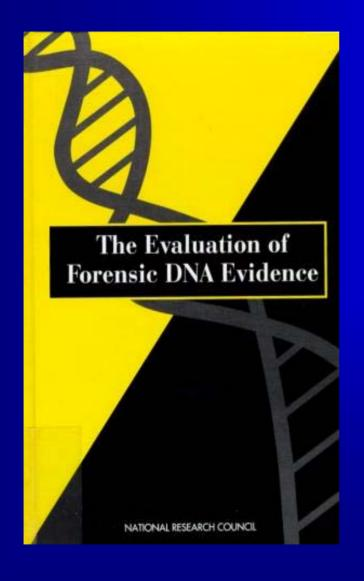
Wrong estimates!

Employ a Theta (θ) Correction

 θ is used as a measure of the effects of population subdivision (inbreeding)

How many Great, Have?

National Research Council Report II



National Academy of Sciences

Data support the recommendation that F_{ST} of 0.01 is conservative

Issued in May 1996

National Research Council Report II

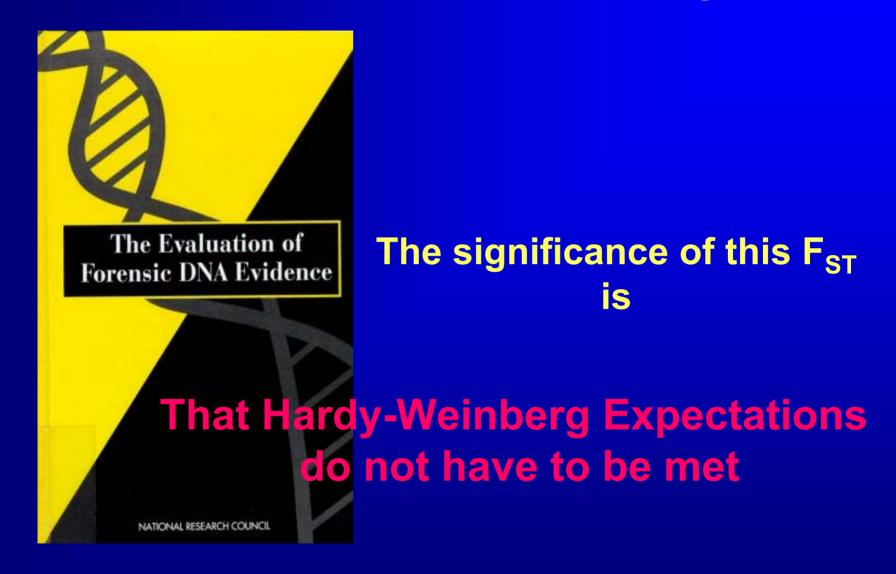


TABLE 6— F_{ST} values for the thirteen CODIS core STR loci.

Locus	African American	Caucasian	Hispanic	Asian	Native American
CSF1PO D3S1358 D5S818 D7S820 D8S1179 D13S317 D16S539 D18S51 D21S11 FGA TH01 TPOX	-0.0009 -0.0005 0.0010 0.0000 -0.0001 0.0029 -0.0013 0.0012 0.0005 0.0004 0.0015 0.0021	-0.0007 -0.0009 -0.0001 -0.0005 0.0000 -0.0008 -0.0005 0.0001 0.0008 -0.0004 -0.0012 -0.0015	-0.0003 0.0014 0.0010 0.0010 0.0005 0.0047 0.0067 0.0011 0.0013 0.0008 0.0041 0.0024	-0.0012 0.0035 0.0028 0.0039 0.0025 0.0071 0.0017 0.0046 0.0056 0.0029 0.0058 0.0100	0.0244 0.0764 0.0656 0.0201 0.0125 0.0157 0.0132 0.0268 0.0371 0.0168 0.0356 0.0164
vWA F _{ST} over all loci	0.0011	-0.0011 -0.0005	0.0029	0.0027	0.0172

Modifying the pro Intermediate to the

$$P^2 + P(1-P)\theta$$

 $\theta = 0.01$

F_{ST} that you would Use correction factor f find in populations with 1st and 2nd cousin matings

0.03 for Native populations

use $2p_ip_i$ for heterozygotes (ie: no correction)

Really, this is more than ten fold more conservative

Modifying the product rule

Formula 4.1 - HW

Formula 4.4 - Simple subdivision

Formula 4.10 - assumption of population

Conditional vs Unconditional Probability

$$p^2 + p(1 - p)\theta$$

$$\frac{[2\theta + (1-\theta)p_{i}][3\theta + (1-\theta)p_{i}]}{(1+\theta)(1+2\theta)}$$

This last formula addresses a conditional probability of the suspect genotype, given that of the perpetrator, $P(A_iA_i \mid A_iA_i)$, considering the person contributing the evidence and the suspect are from the same subgroup.

Although we **CAN** correct the heterozygote genotype estimate...it is not generally necessary.

HWE:

2pq

NRC II, 4.4a:

 $2pq(1-\theta)$

$$\frac{2[\theta + (1 - \theta)p_{i}][\theta + (1 - \theta)p_{j}]}{(1 + \theta)(1 + 2\theta)}$$

$$P(A_iA_j | A_iA_j)$$

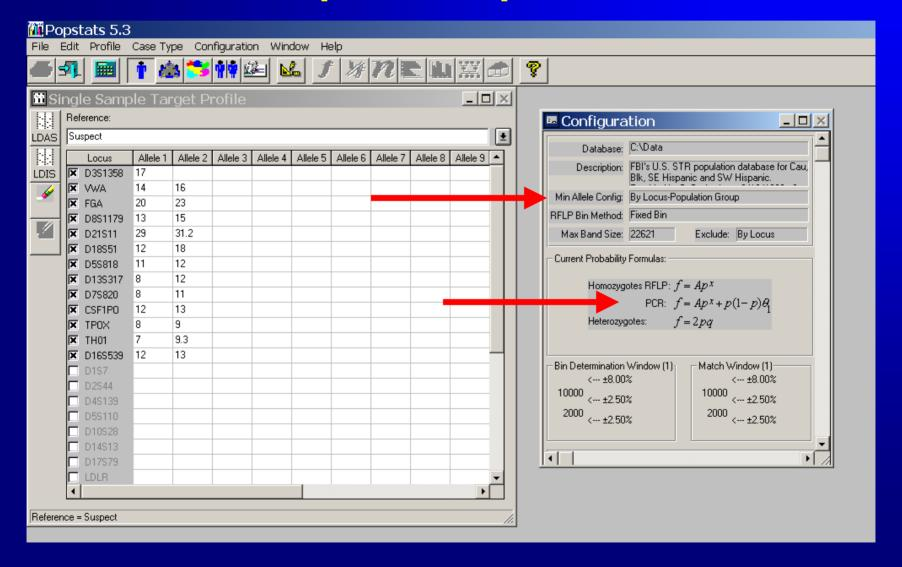
Theta Values Commonly Employed

0.01 for Cau, AA, SEH, and SWH

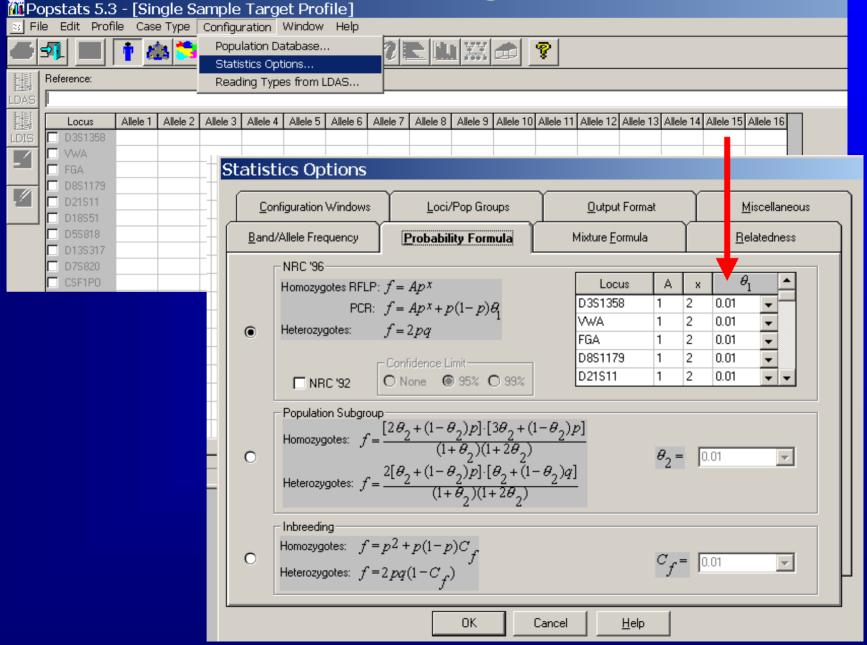
0.03 for Native American groups

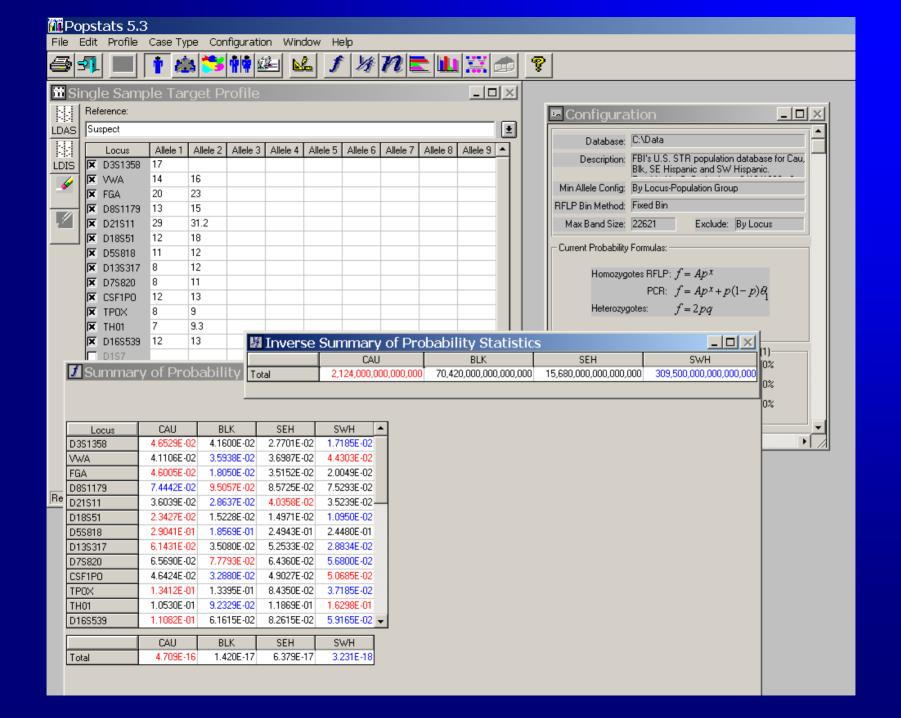
Conservative Values

So how do we deal with these concepts in PopStats??

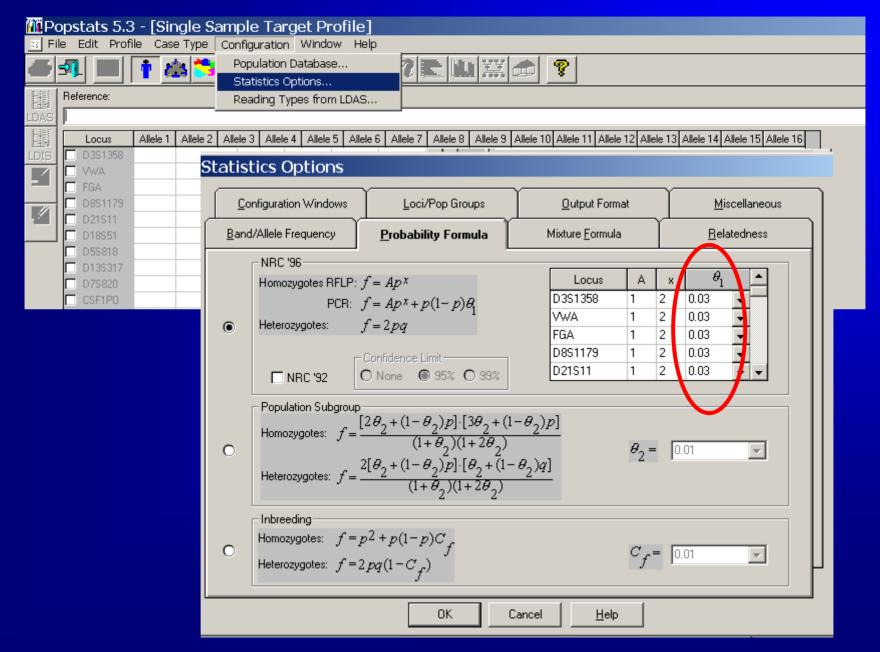


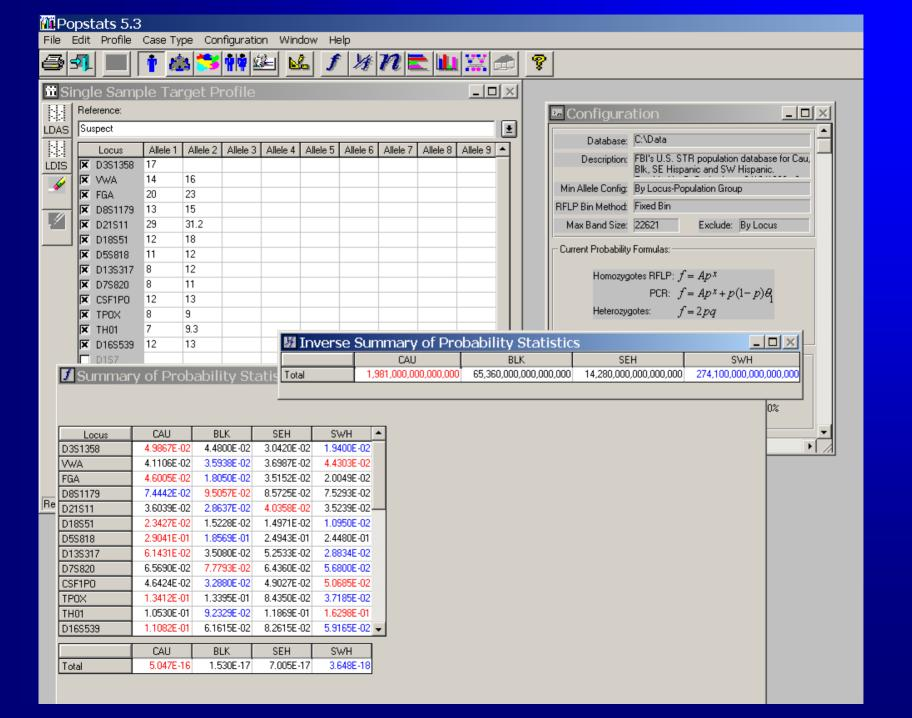
This is the default using 0.01 θ





If we want to use 0.03θ





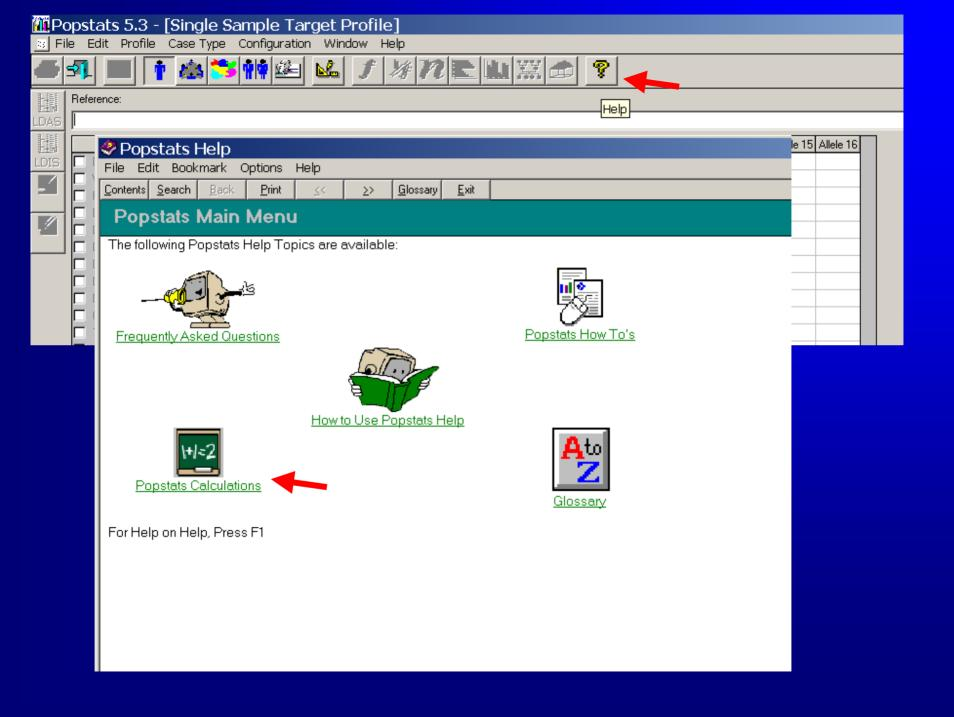
$\theta = 0.01$

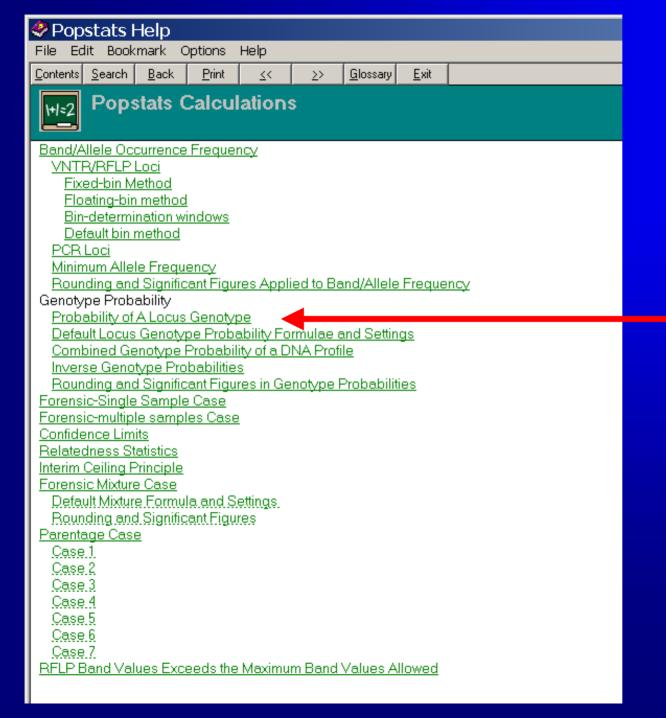
牙	Inverse Summary of Probability Statistics ■□×							
			CAU		BLK	SEH	SWH	
Tota	al	2,1	124,000,000,0	00,000	70,420,000,000,000,000	15,680,000,000,000,000	309,500,000,000,000,000	

$\theta = 0.03$

Inverse Summary of Probability Statistics ■ □ >						
	CAU	BLK	SEH	SWH		
Total	1,981,000,000,000,000	65,360,000,000,000,000	14,280,000,000,000,000	274,100,000,000,000,000		

WHERE CAN WE FIND INFORMATION ON WHAT POPSTATS CAN DO?





Probability of A Locus Genotype



Where

 ρ and q are band/allele occurrence frequencies

 \mathcal{A} and x are user-configurable constants for each locus.

For VNTR/RFLP loci, the default settings are:

$$A = 2$$

 $x = 1$

For PCR loci, the default settings are:

$$A = 1$$
 (A must be positive)

$$x = 2$$
 (x must be non-negative).

2. NRC '96 random-match probability formulae with θ adjustment for homozygotes of PCR loci:

$$f = Ap^{x}$$

$$f = Ap^{x} + p(1-p)\theta_{1}$$

$$f = 2pq$$

- Where p, q, A, and x are defined as above. The parameter θ_1 is only available for PCR loci. The user can enter any value for θ_1 . The default value is $\theta_1 = 0.01$. For some small, isolated populations like the American Indians, a value of 0.03 is suggested for θ_1 by the *National Research Council. (NRC) Report 1996*.
- 3 The conditional probability formulae with population subgroup correction:

hom caygotes:
$$f = \frac{[2\theta_2 + (1-\theta_2)p] \cdot [3\theta_2 + (1-\theta_2)p]}{(1+\theta_2)(1+2\theta_2)}$$

heterozygotes:
$$f = \frac{2[\theta_2 + (1 - \theta_2)p] \cdot [\theta_2 + (1 - \theta_2)q]}{(1 + \theta_2)(1 + 2\theta_2)}$$

- These equations calculate the probability of one person having the DNA profile genotype under the condition that another person has that DNA profile genotype. θ_2 is a user-configurable parameter that represents the degree of population subdivision. Its value is configured globally for all loci and is equal to 0.01 by default.
- 4. The random-match probability formulae with inbreeding correction are

homozygotes:
$$f = p^2 + p(1-p)C_f$$

heterozygotes:
$$f = 2pa(1 - C_s)$$

Where p and q are defined as above. The inbreeding coefficient ($^{C}_{\mathbf{f}}$) is a measure of the degree of inbreeding. The value of $^{C}_{\mathbf{f}}$ is configured globally and can be 0.01, 0.03 or other decimal number. The default value for $^{C}_{\mathbf{f}}$ is 0.01.

Ok...so lets look at our other main option

NRC II, 4.10a:
$$\frac{[2\theta + (1-\theta)p_{i}][3\theta + (1-\theta)p_{i}]}{(1+\theta)(1+2\theta)}$$

NRC II, 4.10b:
$$\frac{2[\theta + (1 - \theta)p_{i}][\theta + (1 - \theta)p_{j}]}{(1 + \theta)(1 + 2\theta)}$$

Note: both homozygotes and heterozygotes are treated in this application

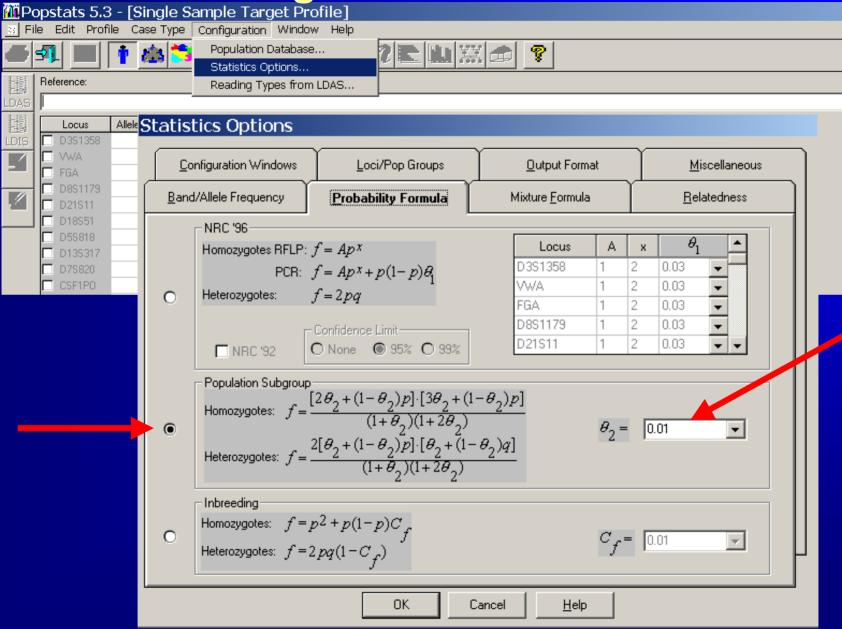
When and why should we consider this??

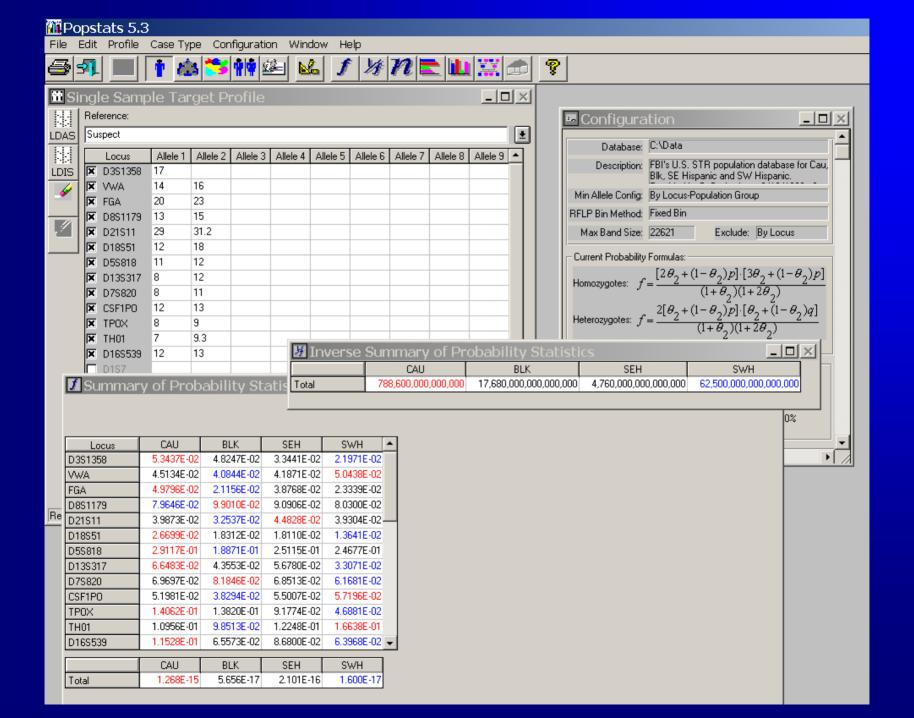
Takes into account the assumption that the person contributing the evidence and the suspect are from the same subgroup

What it gives us is a conditional probability of the suspect genotype given that of the perpetrator.

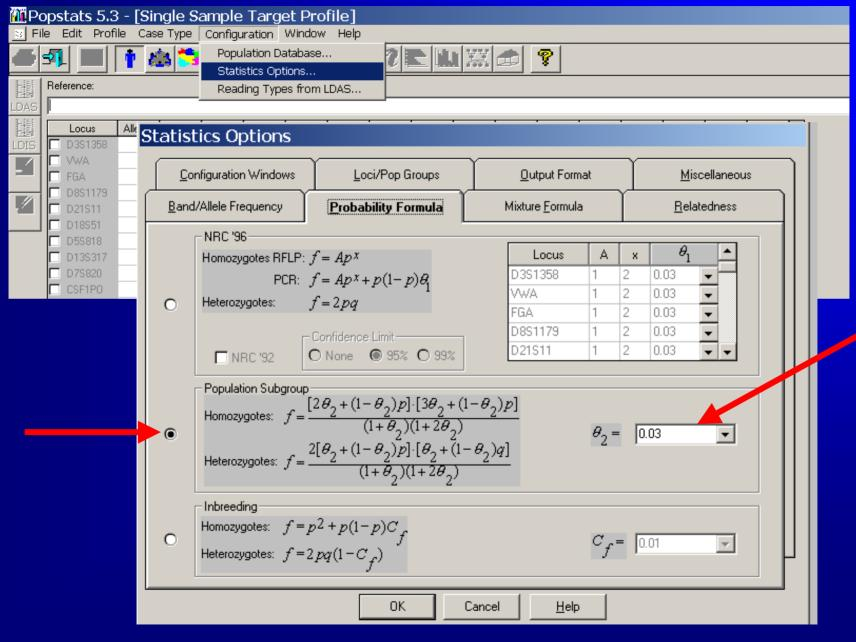
Example... use if the suspect and all possible perpetrators are from the same small isolated town i.e. religious sects, native communities

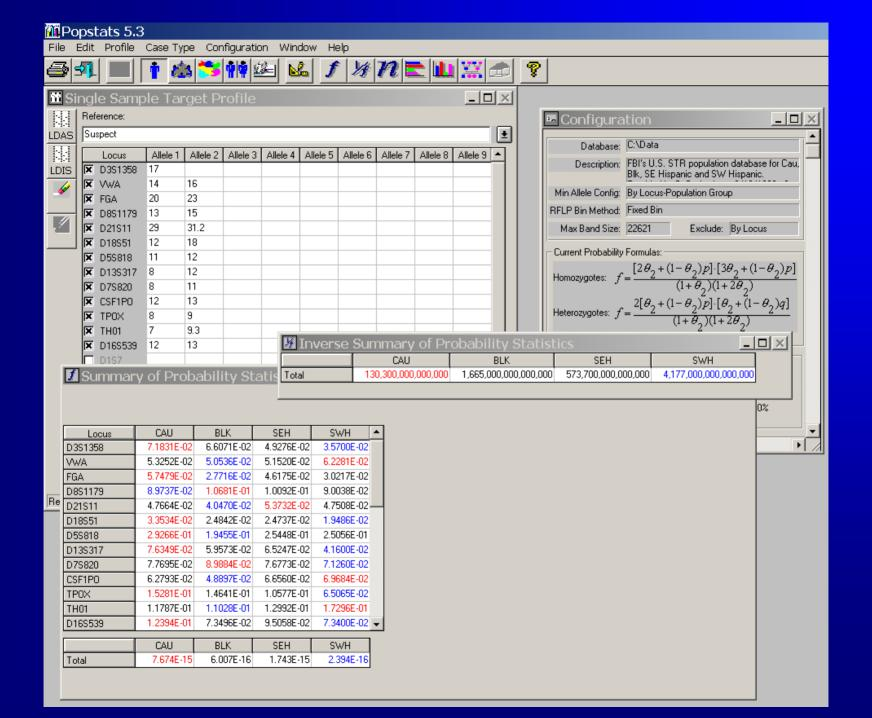
Here it is using the default of 0.01 θ





If we want to use 0.03θ





$\theta = 0.01$

Inverse		×			
	CAU	BLK	SEH	SWH	
Total	788,600,000,000,000	17,680,000,000,000,000	4,760,000,000,000,000	62,500,000,000,000,000	

$\theta = 0.03$

Inverse	Summary of Pr	×			
	CAU	BLK	SEH	SWH	
Total	130,300,000,000,000	1,665,000,000,000,000	573,700,000,000,000	4,177,000,000,000,000	

So,

PopStats has given us the numbers we desired...

What do we do with them???

Well,

We report them of course!

But we should consider what we are reporting and the information we are conveying in our "statistics"

Source attribution

Hot topic for statistical debate

With the current panel of genetic markers available to forensic testing, it is not uncommon for the reciprocal of the random match probability determined for a genetic profile to exceed the worlds population several fold.

So, how do you want to express this fact in your reports and testimony?

What do these numbers mean to you?

the prosecutor?

the defense?

the judge?

the jury?

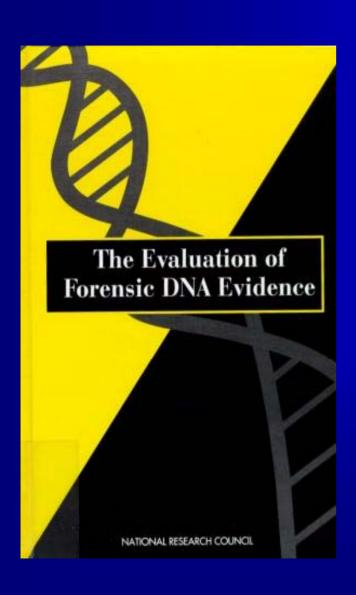
This is what really matters!!!

Big Number Names:

million	1,000,000
billion	1,000,000,000
trillion	1,000,000,000
quadrillion	1×10^{15}
quintillion	1×10^{18}
sextillion	1×10^{21}
septillion	1×10^{24}
octillion	1×10^{27}
nonillion	1×10^{30}
decillion	1×10^{33}

Even Bigger Number Names:

1×10^{36}	undecillion
1×10^{39}	duodecillion
1×10^{42}	tredecillion
1×10^{45}	quattordecillion
1×10^{48}	quindecillion
1×10^{51}	sexdecillion
1×10^{54}	septendecillion
1×10^{57}	octodecillion
1×10^{60}	novemdecillion
1×10^{63}	vigintillion



NRC II May 1996

"...that profile might be said to be unique if it is so rare that it becomes unreasonable to suppose that a second person in the population might have the same profile."

To address uniqueness we are back to the same old question... population sample size

Here the population size differs from what we discussed when calculating allele frequencies...

The relevant population is at issue here

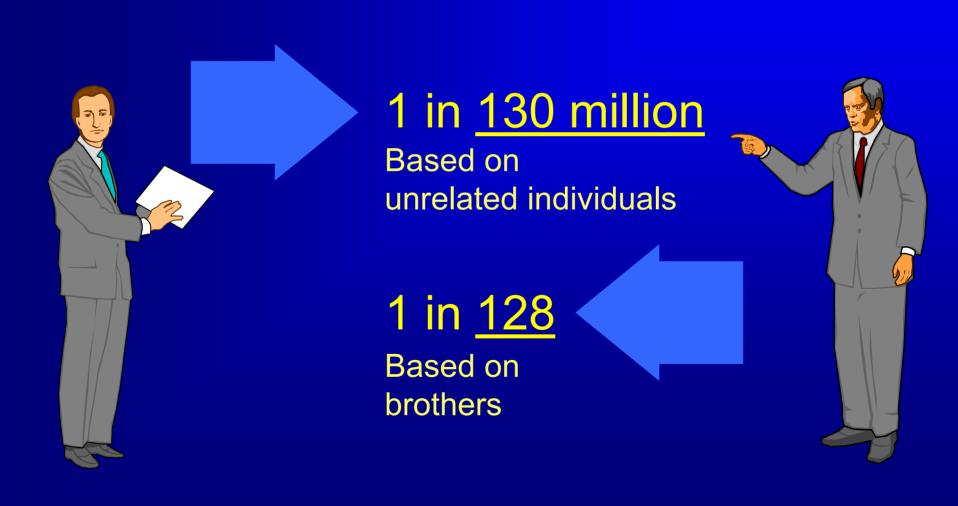
Define the Question

(or at least make sure you know what question you are answering)



Define the Question

Estimates of the Rarity of a DNA Profile:



Uniqueness / Source Attribution Webster's Definitions

only one

Unusual

Some [circumstance] that is the only one of it kind

Uniqueness / Source Attribution Webster's Definitions

Attribution evaluated within context of case

Rarely is the world's population the appropriate context

Thus, a circumstance that is the only one of its kind is appropriate context

Uniqueness

A profile that exists in one person and no other (excluding identical twins)

Context?

- Population of the world...maybe
- Population of the US....there is a thought!
- Population with access to a crime scene...

Uniqueness

A profile that exists in one person and no other (excluding identical twins)

Actually we are interested in source attribution, not whether the profile is unique in the world

Is it reasonable to consider the profile to be so rare that one can opine about the source of the evidence?

Let the RMP of a given <u>evidentiary</u> profile X be p_x (Calculate using NRC II Report Recommendations)

Then
$$(1-p_x)^N$$

is the probability of not observing the profile in a population of N unrelated individuals

This probability should be greater than or equal to a 1- α confidence level

$$(1-p_x)^N \ge 1-\alpha$$
 $p_x \le 1 - (1-\alpha)^{1/N}$

Source Attribution

- Specify (1- α) confidence level of 95% or 99% (uses an α of 0.05 or 0.01, respectively)
- Determine RMP threshold to assert with a specific degree of confidence that the particular evidence profile is unique with a population of N unrelated individuals

What population????

Source Attribution Values

Calculate p for major population groups

 $\theta = 0.01 \text{ or } 0.03$

Take the most common value for p Increase p by factor of 10 Determine if p \leq 1- (1- a) $^{1/N}$

What N??

The standard basis that is used here in the US is an estimate of US population of approximately 260 million people

So, taking this and if we accept an α of 0.01 (99% confidence level) with

$$p_x \leq 1 - (1 - \alpha)^{1/N}$$

A random match probability less than 3.9×10^{-11} would convey at least 99% confidence that the evidentiary profile is unique in the population

RMP thresholds for source attribution at various population sizes and confidence levels

SAMPLE		CONFIDEN	CE LEVELS	
SIZE N	0.90	0.95	0.99	0.999
2	5.1x10 ⁻²	2.5x10 ⁻²	5.0x10 ⁻³	5.0x10 ⁻⁴
3	3.5x10 ⁻²	1.7x10 ⁻²	$3.3x10^{-3}$	3.3x10 ⁻⁴
4	2.6x10 ⁻²	1.3x10 ⁻²	2.5x10 ⁻³	2.5x10 ⁻⁴
5	2.1x10 ⁻²	1.0x10 ⁻²	2.0x10 ⁻³	2.0x10 ⁻⁴
6	1.7x10 ⁻²	8.5x10 ⁻³	1.7x10 ⁻³	1.7x10 ⁻⁴
7	1.5x10 ⁻²	7.3×10^{-3}	1.4x10 ⁻³	1.4x10 ⁻⁴
8	1.3x10 ⁻²	6.4×10^{-3}	1.3x10 ⁻³	1.3x10 ⁻⁴
9	1.2x10 ⁻²	$5.7x10^{-3}$	1.1x10 ⁻³	1.1x10 ⁻⁴
10	1.1x10 ⁻²	5.1×10^{-3}	1.0x10 ⁻³	1.0x10 ⁻⁴
25	4.2x10 ⁻³	2.1x10 ⁻³	4.0x10 ⁻⁴	4.0x10 ⁻⁵
50	2.1x10 ⁻³	1.0x10 ⁻³	2.0x10 ⁻⁴	2.0x10 ⁻⁵
100	1.1x10 ⁻³	5.1x10 ⁻⁴	1.0x10 ⁻⁴	1.0x10 ⁻⁵
1x10 ³	1.1x10 ⁻⁴	5.1x10 ⁻⁵	1.0x10 ⁻⁵	1.0x10 ⁻⁶
1×10^{5}	1.1x10 ⁻⁶	$5.1x10^{-7}$	1.0x10 ⁻⁷	1.0x10 ⁻⁸
1x10 ⁶	1.1x10 ⁻⁷	5.1x10 ⁻⁸	1.0x10 ⁻⁸	1.0x10 ⁻⁹
1×10 ⁷	1.1x10 ⁻⁸	5.1x10 ⁻⁹	1.0x10 ⁻⁹	1.0x10 ⁻¹⁰
5x10 ⁷	2.1x10 ⁻⁹	1.0x10 ⁻⁹	2.0.10-10	2.0x10 ⁻¹¹
2.6x10 ⁸	4.1x10 ⁻¹⁰	2.0x10-0	3.9x10 ⁻¹¹	3.9x10 ⁻¹²
1x109	1.1x10 ⁻¹⁰	5.1x10 ⁻¹¹	1.0x10 ⁻¹¹	1.0x10 ⁻¹²
5x109	2.1x10 ⁻¹¹	1.0x10 ⁻¹¹	2.0x10 ⁻¹²	2.0x10 ⁻¹³

So with our PopStats results obtained for a θ = 0.01

		51.17	oeu	0	
Locus	CAU	BLK	SEH	SWH	
D3S1358	4.6529E-02	4.1600E-02	2.7701E-02	1.7185E-02	
VWA	4.1106E-02	3.5938E-02	3.6987E-02	4.4303E-02	
FGA	4.6005E-02	1.8050E-02	3.5152E-02	2.0049E-02	
D8S1179	7.4442E-02	9.5057E-02	8.5725E-02	7.5293E-02	
D21S11	3.6039E-02	2.8637E-02	4.0358E-02	3.5239E-02	
D18S51	2.3427E-02	1.5228E-02	1.4971E-02	1.0950E-02	
D5S818	2.9041E-01	1.8569E-01	2.4943E-01	2.4480E-01	
D13S317	6.1431E-02	3.5080E-02	5.2533E-02	2.8834E-02	
D7S820	6.5690E-02	7.7793E-02	6.4360E-02	5.6800E-02	
CSF1PO	4.6424E-02	3.2880E-02	4.9027E-02	5.0685E-02	
TPOX	1.3412E-01	1.3395E-01	8.4350E-02	3.7185E-02	
TH01	1.0530E-01	9.2329E-02	1.1869E-01	1.6298E-01	
D16S539	1.1082F-01	C.1015E 02	8.2815E 92	5 9165E-02	T
	CAU	BLK	SEH	SWH	
Total	4.709E-16	1.420E-17	6.379E-17	3.231E-18	

 3.9×10^{-11}

Profile frequency is less than 99% threshold

So with our PopStats results obtained for a θ = 0.03

Locus	CAU	BLK	SEH	SWH	_
D3S1358	4.9867E-02	4.4800E-02	3.0420E-02	1.9400E-02	
VWA	4.1106E-02	3.5938E-02	3.6987E-02	4.4303E-02	
FGA	4.6005E-02	1.8050E-02	3.5152E-02	2.0049E-02	
D8S1179	7.4442E-02	9.5057E-02	8.5725E-02	7.5293E-02	
D21S11	3.6039E-02	2.8637E-02	4.0358E-02	3.5239E-02	
D18S51	2.3427E-02	1.5228E-02	1.4971E-02	1.0950E-02	
D5S818	2.9041E-01	1.8569E-01	2.4943E-01	2.4480E-01	
D13S317	6.1431E-02	3.5080E-02	5.2533E-02	2.8834E-02	
D7S820	6.5690E-02	7.7793E-02	6.4360E-02	5.6800E-02	
CSF1P0	4.6424E-02	3.2880E-02	4.9027E-02	5.0685E-02	
TPOX	1.3412E-01	1.3395E-01	8.4350E-02	3.7185E-02	
TH01	1.0530E-01	9.2329E-02	1.1869E-01	1.6298E-01	
D16S539	1,1082F-01	0.1015E-02	0.2015E-02	F 9165F-02	Ţ
	CAU	BLK	SEH	SWH	
Total	5.047E-16	1.530E-17	7.005E-17	3.648E-18	

 3.9×10^{-11}

Profile frequency is less than 99% threshold

"To a reasonable degree of scientific certainty, _____ is the source of the DNA in specimen Q2."

"I have a high degree of confidence, is the source of the DNA in specimen Q2."

"Based on an estimate of 260 million people resident in the population of the United States, there is 99% confidence that ____ is the source of the DNA in specimen Q2."

We are not stating that ______ is the only person to possess that profile. We are stating that we would not expect to find it in a population of N individuals.

Source Attribution

- Method is simple
- Conservative because N is so large (260,000,000)
- If N = 260,000,000, then RMP threshold is 3.9×10^{-11}
- Most of the time the RMP is far less, so confidence is greater than 0.99

Source Attribution

- N can be configured to context of the case
- Two individuals to entire town, state, or whatever
- Laboratory policy to set N

Random match probability is NOT

Chance that someone else is guilty

Chance that someone else left the bloodstain

Chance of defendant not being guilty



"We are neither hunters nor gatherers. We are statisticians."

Internission