

GeneMarker® HID

**Expert System Human Identification Software with Linked
Post-Identification Database Search, Kinship, and DNA
Mixture Applications**

Promega Webinar
July 24, 2013



SoftGenetics LLC

- Founded: 2001

- Mission:

“Provide Biologists with highly accurate, sensitive & user-friendly analysis tools.”

- Founders:

CS Liu, PhD. VP Development.
VP Software Development, Spectrumedix
17 Issued Patents

John Fosnacht, VP Sales & Marketing VP Sales
& Marketing, Spectrumedix ; 2 years
President, AccuStandard, 4 years General
Manager EMD Reagents USA, 17 years

Product Portfolio:

GeneMarker®HID STR Analysis

- Human Identity Expert System
- Mixture Applications
- Familial and Exact Match Database Search
- Parentage and Kinship Testing

Mutation Surveyor® for Sanger Sequence

- mtDNA Analysis and SNP detection
- Detect Heteroplasmy
- Patented Core Technology

NextGENe® analysis of NGS Sequencing

- Patent Pending Core Technology
Condensation of “short reads”
- Analysis of mtDNA, STRs and mixtures

[Croat Med J.](#) 2011 Jun;52(3):299-313.

Second generation sequencing allows for mtDNA mixture deconvolution and high resolution detection of heteroplasmy.

[Holland MM](#), [McQuillan MR](#), [O'Hanlon KA](#).

GeneMarker®/MTP Genotyping

AFLP®/TRFLP
Microsatellite, Trisomy
MLPA®, MS-MLPA
LOH/ MSI
SNaPshot®

ChimerMarker™

STR Analysis for Chimerism Testing
Single Donor Chimerism
Double Donor Chimerism
Long-Term Monitoring
MCC Analysis

GeneticistAssistant™

NGS Web-based visualization tool

JelMarker®

Reading and converting gel images

GeneMarker® HID (Human Identity)

- **Concordant, Validated Expert System**
 - Alternative to Genotyper®, GeneMapper® IDX
 - Compatible with major chemistries, CE and Rapid Systems output; Windows® XP, Vista, 7 and 8
- **Single Source casework and databasing**
 - Streamlined Workflow with Audit Trail
 - CODIS compatible reporting
- **Reference Sample and Familial Search**
 - Mass Disaster, Crime Scene Response --- No Data Transfer Needed
 - Positive Matches Ranked by Likelihood Ratio
- **Kinship and Paternity Testing**
 - Identity by Descent (IBD) Calculations with pre-loaded or custom allele frequency tables
 - Automated Pedigree drawing and AABB trio/single parent PI calculations
- **Mixture Analysis**
 - Detects Mixtures and calculates Probability of Inclusion and Exclusion (PI, PE) for all mixtures
 - Performs Statistics two contributor including LR and searches database for deduced contributor (with or without reference sample)

Run Wizard

Template Selection

Run Wizard

Data Process - HID Analysis

Run Wizard

Additional Settings - HID Analysis

Set additional options related to the different analysis types

Allelic Ladder:

Positive Control Template:

Allele Evaluation

Peak Score:

Reject < Check

☐ Mixture Evaluation

Valid Mixture Peak Percentage:

Min Mixture Marker Number:

Access Rights of User Types

Access Rights:

- ☒ Edit Alleles
- ☒ Insert Alleles
- ☒ Delete Alleles
- ☒ Undelete Alleles
- ☒ Confirm Alleles
- ☒ Unconfirm Alleles
- ☒ Comment Alleles
- ☒ Recall Alleles
- ☒ Re-Run
- ☐ **Set Run Option**
- ☒ Change Template
- ☒ Edit Calibration
- ☒ Recover Data
- ☒ Enable/Disable Samples
- ☒ Comment Samples
- ☒ Add Samples to Project
- ☒ Comment Project
- ☒ Save Project
- ☐ Edit Panel
- ☐ Edit Size
- ☒ Export Trace
- ☐ Update Software
- ☐ Edit Mixture Analysis Parameters

User Type:

Reviewer
Analyst
 Lab Manager
 Administrator

Users:

Set Default

Ok

Cancel

Review Allele Calls In Main Analysis Screen or All Color Browser

Any edits or comments are recorded as part of the audit trail

GeneMarker HID - Analyst_A_w_patDiagInclusionExclusionExamplePaternityTests.SGF

File View Project Applications Tools Help

Analyst_A_w_patDiagInclusionExclusionExamplePaternityTests.SGF

Raw Data

Allele Call

- LD: LADDERD09.hid
- LD: LADDERD09.hid
- LD: LADDERE09.hid
- LD: LADDERE09.hid
- Fusion_TrioB_AF_500pg_1E02.h
- Fusion_TrioB_C_500pg_1C02.hic
- Fusion_TrioB_M_500pg_1A02.hic
- Fusion_TrioE_AF_500pg_2B05.h
- Fusion_TrioE_C_500pg_2D05.hic
- Fusion_TrioE_M_500pg_2F05.hic

Fusion_TrioE_M_500pg_2F05.hid (Ref Ladder: LADDERE09.hid)

TH01 vWA D21S11 D7S820 D5S818 TPOX DYS391

9.3 16 28 32 10 11 12 8

No.	Size	Height	Marker	Allele	Allele Comments	Quality Reasons
1	97.8	13717	TH01	9.3		
2	152.0	9327	vWA	16		
3	220.2	3882	D21S11	28		
4	236.5	4488	D21S11	32		
5	289.3	4575	D7S820	10		
6	293.3	4056	D7S820	11		
7	342.6	4525	D5S818	11		
8	346.6	3888	D5S818	12		
9	410.0	9242	TPOX	8		
10	449.2	89	DYS391	6	[<Deleted>]extreme pull up	IHO, LS

Report Table Actions:

- Insert an Allele
- Delete an Allele
- Delete Alleles
- Edit an Allele
- Confirm an Allele
- Confirm Alleles
- Sort Alleles

Peak Quality Reasons:

- LS = Low Score
- OL = Off Ladder
- OB = Out of Bin
- BC = Bin Conflict
- SR = Saturated (Repaired)
- SP = Saturated (Pull- up)
- PL = Beyond Ploidy
- LO = Low Intensity
- HI = High Intensity
- IMB = Heterozygote Imbalance
- IHO = In Homozygote Inconclusive
- IHE = In Heterozygote Inconclusive

Project Panel .XML Recognizes Variant Alleles and Shifts Bins to Fit Allelic Ladders Marker Specific Settings to Customize to the Lab Standard Operating Procedure

The screenshot displays the 'Panel Editor' software interface. On the left, a 'Project Panel' tree shows a list of markers. The main window shows a gel electrophoresis image with peaks labeled with marker names and sizes. Below the gel image is a table of marker data.

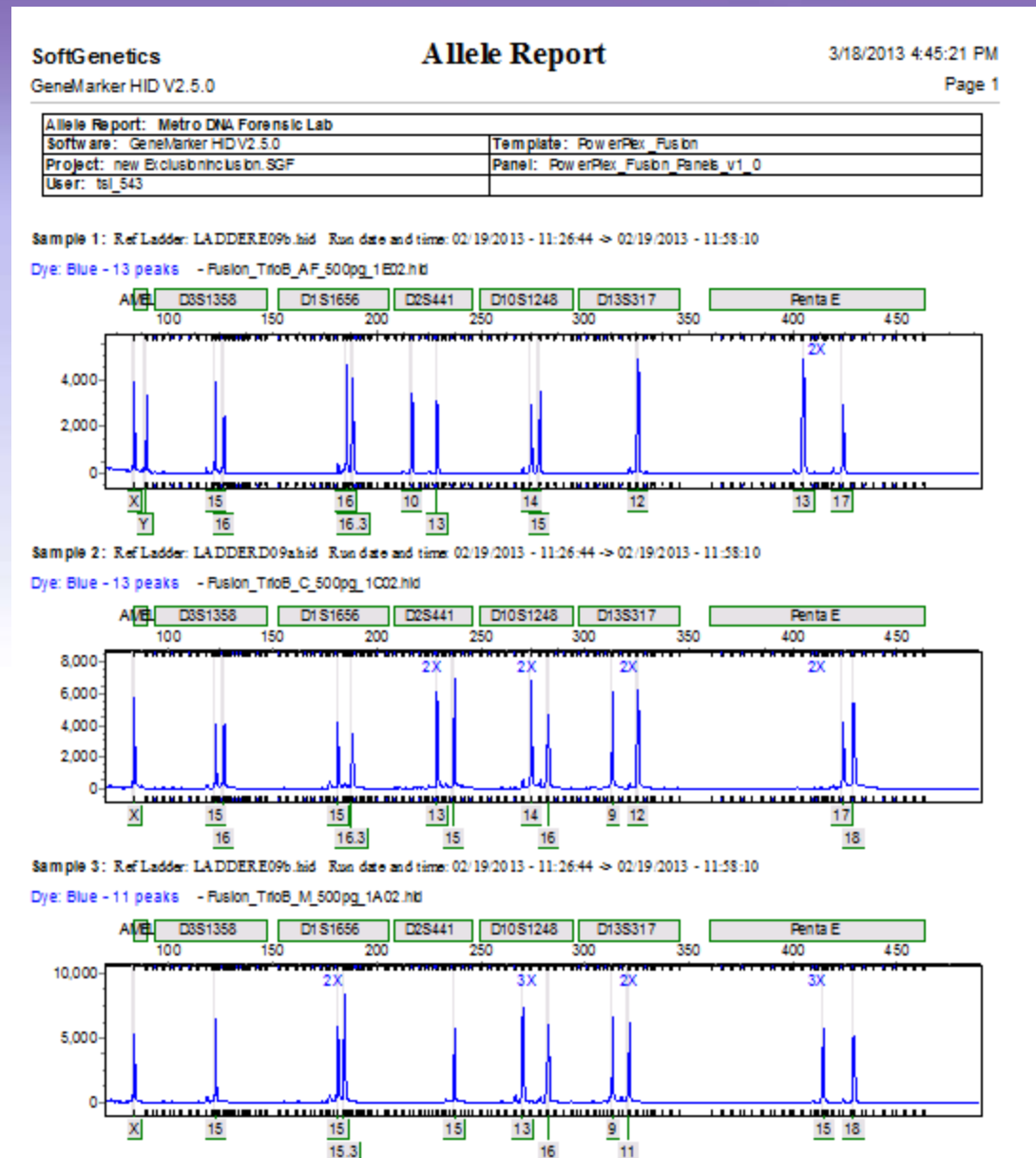
No.	Dye	Marker	Size	Left Range	Right Range
90	Blue	D13S317	332.6	0.5	0.5
91	Blue	D13S317	333.6	0.4	0.5
92	Blue	D13S317	337.7	0.5	0.5
93	Blue	D13S317	341.7	0.5	0.5
94	Blue	D13S317	345.8	0.5	0.5
95	Blue	Penta E	361.3	0.5	0.5
96	Blue	Penta E	366.2	0.5	0.5
97	Blue	Penta E	371.0	0.5	0.5
98	Blue	Penta E	375.8	0.5	0.5
99	Blue	Penta E	380.7	0.5	0.5
100	Blue	Penta E	385.6	0.5	0.5
101	Blue	Penta E	390.5	0.5	0.5
102	Blue	Penta E	395.4	0.5	0.5

The 'Edit Marker' dialog box is open, showing the following settings for marker D3S1358:

- Marker Name: D3S1358
- Nucleotide Repeats (x): 4
- Boundary: 93.5 To 148.2
- Min Homozygote Intensity: 50
- < = Inconclusive < =: 200
- Min Heterozygote Intensity: 50
- < = Inconclusive < =: 100
- Max Heterozygote Imbalance: 60 %
- Min Heterozygote Imbalance: 20 %
- ☐ Apply Homo/Hetero Settings to All Markers
- Stutter Filter:
 - N - x: 12 %
 - N - 2x: 1 %
 - N + x: 1 %
- ☐ Apply Stutter Settings to All Markers

The background shows a gel electrophoresis image with peaks labeled with marker names and sizes. The 'Penta E' marker is highlighted in the gel image.

Reporting Options -- Print/Save Allele Reports



Print Report

Print Type
☒ Normal ☐ Chart Overlay

Samples
☒ All Samples ☐ Selected Samples

Contents
☒ Electropherogram
☐ Peak Table
☐ Follow trace chart
☒ Start after all charts finished
☐ Start on separate page
☐ Forensics Table

Dyes
☒ Dye1 ☒ Dye5
☒ Dye2 ☐ Dye6
☒ Dye3
☒ Dye4
☐ Mix Dyes
☐ Hide Bins

Advance Options
☒ Print Project Comments -- on Each Page ☐ Word Wrap ☐
☒ Print Report Header -- on Each Page ☐
☐ Label Dyes & Peak Numbers ☒ Print Markers
☒ Implement Y Axis Settings ☒ Print Alleles
☐ New Page for Each Sample ☐ Abide By Panel
☒ Grouped by Dye ☒ Auto Scale Markers
☐ Mark Deleted/Edited Peaks ☐ Print Edited Peak Only
☐ Label Peak Ratio ☐ Print Samples With Grouping
☒ Label Peak Height Ratio ☐ Label Peak Area Ratio
 Chart Height (mm): 55

Report Table Format Options:

Report	Bin														
	Sample	Marker	Allele#1	Size#1	Height#1	Ht_Ratio#1	Area#1	Ar_Ratio#1	Allele#2	Size#2	Height#2	Ht_Ratio#2	Area#2	Ar_Ratio#2	
1	Fusion_TrioB_AF	AMEL	X	84.0	3994	1.00	22185	1.00	Y	89.8	3312	0.83	18839	0.85	
		D3S1358	15	122.8	3927	1.00	22348	1.00	16	126.8	2491	0.63	14175	0.63	
		D1S1656	16	185.7	4676	1.00	27201	1.00	16.3	188.7	4071	0.87	23740	0.87	
		D2S441	10	217.1	3436	1.00	21333	1.00	13	229.3	3155	0.92	19729	0.92	
		D10S1248	14	274.5	3018	0.85	20548	0.83	15	278.4	3557	1.00	24626	1.00	
		D13S317	12	325.5	4909	1.00	37273	1.00							
		Penta E	13	405.0	2436	1.00	20903	1.00	17	424.1	1491	0.61	13416	0.64	
		D16S539	9	98.5	4809	1.00	25623	1.00	11	106.8	4531	0.94	25144	0.98	
		D18S51	15	163.0	4897	1.00	27272	1.00	18	175.1	4282	0.87	24077	0.88	
		D2S1338	21	264.8	4960	1.00	33252	1.00	22	268.8	4492	0.91	30771	0.93	
		CSF1PO	10	334.8	4486	1.00	34780	1.00	11	338.8	3721	0.83	29062	0.84	
		Penta D	12	423.9	1794	0.83	16802	0.80	14	434.1	2162	1.00	20893	1.00	
		TH01	7	86.9	4685	1.00	24641	1.00	8	90.9	3735	0.80	19870	0.81	
		vWA	16	152.1	3828	0.92	22237	0.92	18	160.1	4147	1.00	24049	1.00	
		D21S11	30	228.4	3573	1.00	22960	1.00	31	232.4	3362	0.94	21250	0.93	
		D7S820	10	289.3	3206	0.83	22520	0.82	12	297.3	3875	1.00	27541	1.00	
		D5S818	8	330.4	3819	1.00	29269	1.00	13	350.6	2650	0.69	20699	0.71	
		TPOX	8	410.0	7293	1.00	63871	1.00							
		DYS391	10	464.9	3244	1.00	Allele Report Settings								
		D8S1179	13	101.2	5983	1.00	<div>Report Style</div> <div><input type="radio"/> Allele List</div> <div><input type="radio"/> Forensics</div> <div><input type="radio"/> Bin Table</div> <div><input checked="" type="radio"/> Peak Table</div> <div><input type="radio"/> Allele Count</div> <div><input type="radio"/> Sample Name <input checked="" type="radio"/> File Name</div> <div>Orientation</div> <div><input checked="" type="radio"/> Horizontal <input type="radio"/> Vertical</div> <div>Options</div> <div><input type="checkbox"/> Size Range (bps)</div> <div>From <input type="text" value="0"/> to <input type="text" value="500"/></div> <div><input checked="" type="checkbox"/> Abide By Panel</div> <div><input checked="" type="checkbox"/> Grouped by Markers</div> <div><input type="button" value="Columns..."/></div> <div><input checked="" type="checkbox"/> Show ^{xxx} when no allele call</div> <div><input type="checkbox"/> Show Only Uncertain Alleles</div> <div><input checked="" type="checkbox"/> Show Rejected Low Score Alleles</div>								
		D12S391	17	147.8	6568	1.00									
		D19S433	13	223.2	4527	1.00									
		FGA	21	294.2	4808	0.96									
		D22S1045	11	440.0	3479	1.00									
2	Fusion_TrioB_C_S	AMEL	X	84.0	5757	1.00									
		D3S1358	15	123.1	4164	1.00									
		D1S1656	15	181.7	4197	1.00									
		D2S441	13	229.3	3095	0.90									
		D10S1248	14	274.5	3441	1.00									
		D13S317	9	313.3	3098	0.99									
		Penta E	17	424.1	2097	0.77									

Allele Report Settings

Report Style:

- ☐ Allele List
- ☐ Forensics
- ☐ Bin Table
- ☒ Peak Table
- ☐ Allele Count

☐ Sample Name ☒ File Name

Orientation:

- ☒ Horizontal
- ☐ Vertical

Options:

- ☐ Size Range (bps)

From to
- ☒ Abide By Panel
- ☒ Grouped by Markers
-
- ☒ Show when no allele call
- ☐ Show Only Uncertain Alleles
- ☒ Show Rejected Low Score Alleles
- ☒ Hide Extra Sample Names

Set Peak Table Columns

All Columns:

- Dye
- Marker
- Difference
- Quality
- Score
- Start
- End
- Allele Comments
- Sample Comments
- Quality Reasons

Selected Columns:

- Allele
- Size
- Height
- Ht_Ratio
- Area
- Ar_Ratio

Expert System Summary

Accurate, Rapid Analysis of Forensic STR Data

Quality Flagging, Size Calibration Check, Automated Control Concordance

Compatible with all major file types

from CE and Rapid Instruments and Human Identity PCR Kits and Custom Chemistries

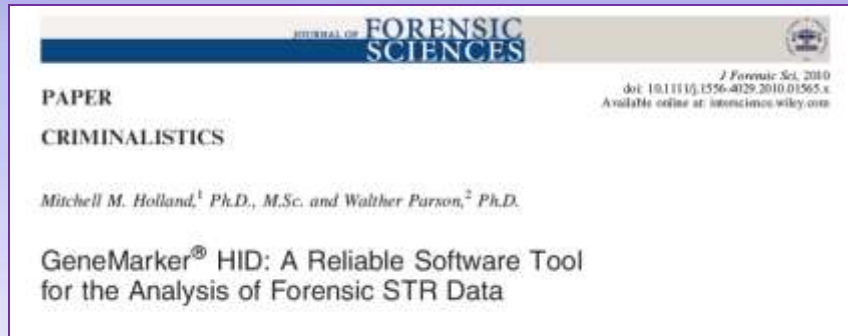
User Management, Access Rights and Audit Trail

Linked Navigation – Time Savings and Automated Edit History

Report Flexibility – Allele, Peak Table and CODIS formats

Selected Presentations and Journal Articles

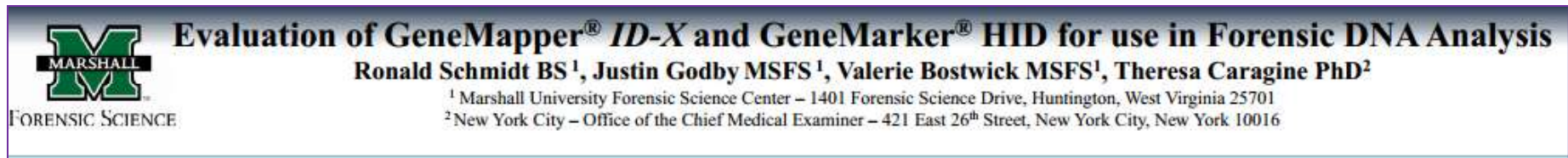
Expert System, Concordance with GeneMapper®
estimated analysis time savings 25%



Autosomal and Y STR analysis of LCN data



“..GeneMarker HID to be a more efficient software system....results in a gain of information and a significant reduction in the number of edits needed per sample..making it the preferred software package...NYC OCME”



GeneMarker HID better able to
resolve alleles with Single-Base-Pair

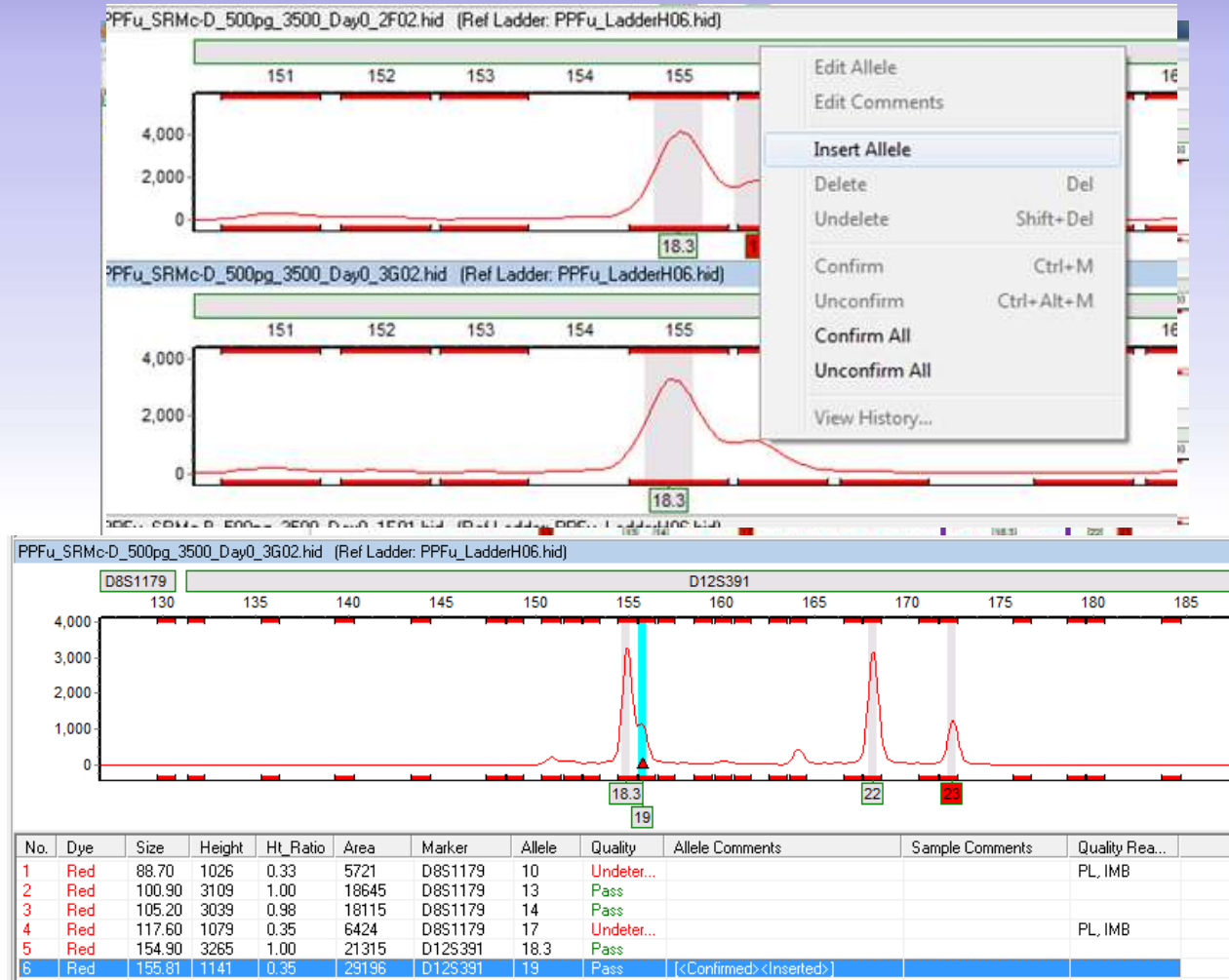


Example of Single Base Pair Resolution

Samples amplified with PowerPlex® Fusion and separated on a Applied Biosystems® 3500 CE

Minor component allele 19 is automatically called in 2 out of 3 samples using default analysis settings

Manually insert allele 19 using a right mouse click to select insert allele



Searchable Database - Relationship Testing

Match Reference Samples or Family Members

Likelihood Ratios calculated using Identity by Descent

- **Missing Persons, Mass Disasters**
- **Immigration / Human Trafficking**
- **Detect Bioterrorism – with the appropriate chemistry development**
- **Paternity/Kinship -- Three Generations (IBD)**
- **Pedigree Diagrams – Visualize Allele Conflicts**
PI Calculations AABB Recommendations

Missing Person Example

Relationship Testing Settings

Samples
☒ All Samples ☐ Selected Samples ...

Default Allele Frequency
 Panel Name: PowerPlex Fusion
 Population: US African-American

☒ **Mutation**
 Mutation Rate: AABB 2003
 Max Mutation Markers: 2
☒ **STR Step Difference**
 One Extra Step STR Mutation Ratio: 0.100000
☐ Paternal and Maternal Mutation Difference

Prior Probability: 0.0500

Search Report Display
☒ Gender Determined
 Minimum LR Value: 10.000 Maximum File Number: 10

☒ Display Genotypes ☒ Display Locus Identifier
☐ Save Parameters when Save Report Normal

Search Scope

Relationships

- ☒ Sample-Individual
- ☒ Father/Son
- ☒ Mother/Daughter
- ☒ Full-Sibs
- ☒ Half-Sibs

Matched Ratio

Matched Ratio with Full-Sibs

Matched Ratio without Full-Sibs

Ok Cancel

Samples
Charts
Report
Calculation Details

	File Name	ID	Name	X/Y	Matched Alleles	Matched Markers	PI/KI
US African-American							
1	Same-Individual Fusion_TrioB_M_50...	1017		XX	46/46	23/23	2.59E+36
	Father/Son						
1	Mother/Daughter Fusion_TrioB_C_500...	1016		XX	26/46	23/23	2.57E+11
	Full-Sibs						
	Half-Sibs						
US Asian-American							
1	Same-Individual Fusion_TrioB_M_50...	1017		XX	46/46	23/23	3.17E+34
	Father/Son						
1	Mother/Daughter Fusion_TrioB_C_500...	1016		XX	26/46	23/23	1.20E+09
	Full-Sibs						
	Half-Sibs						
US Caucasian							
1	Same-Individual Fusion_TrioB_M_50...	1017		XX	46/46	23/23	5.73E+31
	Father/Son						
1	Mother/Daughter Fusion_TrioB_C_500...	1016		XX	26/46	23/23	1.74E+09
	Full-Sibs						
	Half-Sibs						
US Hispanic-American							
1	Same-Individual Fusion_TrioB_M_50...	1017		XX	46/46	23/23	2.26E+32
	Father/Son						
1	Mother/Daughter Fusion_TrioB_C_500...	1016		XX	26/46	23/23	4.79E+08
	Full-Sibs						
	Half-Sibs						

Likelihood Ratios of Relationship Level versus Random Person From the Population Calculated using Identity by Descent Equations and Allele Frequency Tables. Highest scoring relationship level displayed in exportable table

Kinship Analysis Across Three Generations

Kinship Analysis

Single Pair | Multi-Pair

Marker	Individual A Fusion_TrioB_C_50C		Individual B AF_500pg_1E02.hid		Parent/Child (LR)	Full-Sibs (LR)	Half-Sibs (LR)	Cousins (LR)
D3S1358	15	16	15	16	1.62584	2.38318	1.31292	1.15646
D1S1656	15	16.3	16	16.3	2.68068	1.59034	1.84034	1.42017
D2S441	13	15	10	13	5.42652	2.96326	3.21326	2.10663
D10S1248	14	16	14	15	0.91565	0.70783	0.95783	0.97891
D13S317	9	12	12		1.15207	0.82604	1.07604	1.03802
PENTA E	17	18	13	17	5.95238	3.22619	3.47619	2.23810
D16S539	9	10	9	11	1.32443	0.91222	1.16222	1.08111
D18S51	17	18	15	18	2.25225	1.37613	1.62613	1.31306
D2S1338	20	21	21	22	1.93483	1.21742	1.46742	1.23371
CSF1PO	11	14	10	11	1.22070	0.86035	1.11035	1.05518
PENTA D	12		12	14	4.31035	2.40517	2.65517	1.82759
TH01	6	7	7	8	0.63937	0.56968	0.81968	0.90984
vWA	17	18	16	18	2.17391	1.33696	1.58696	1.29348
D21S11	28	31	30	31	3.16456	1.83228	2.08228	1.54114
D7S820	11	12	10	12	2.76243	1.63122	1.88122	1.44061
D5S818	12	13	8	13	1.07759	0.78879	1.03879	1.01940
TPQX	8	10	8		1.49254	0.99627	1.24627	1.12313
D8S1179	13	14	13	15	1.20773	0.85386	1.10386	1.05193
D12S391	18.3	22	17	22	6.35647	3.42824	3.67824	2.33912
D19S433	13	14	13	14	2.24156	3.84595	1.62078	1.31039
FGA	23	24	21	23	1.40449	0.95225	1.20225	1.10112
D22S1045	11	16	11	15	1.76591	1.13296	1.38296	1.19148
Product Score:					3.76E+06	1.46E+03	1.64E+04	3.36E+02

Kinship Analysis Settings

Relationships

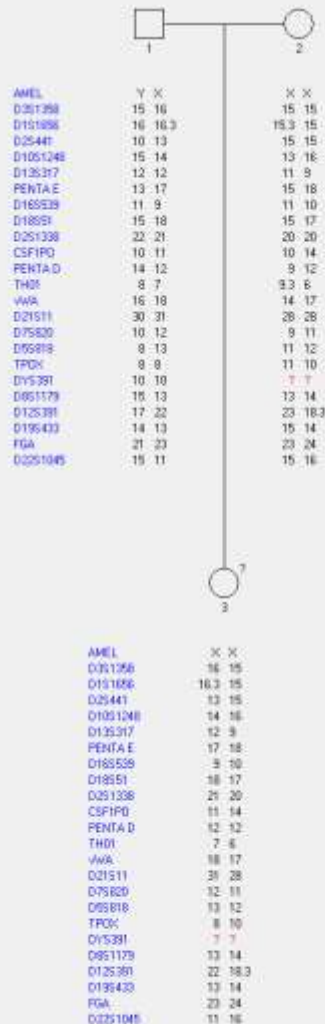
- ☒ Unrelated
- ☒ Parent/Child
- ☒ Full-Sibs
- ☒ Half-Sibs
- ☐ Uncle-Nephew
- ☒ Cousins
- ☐ Grandparent-Grandchild

Report Content

- ☒ Likelihood Ratio
- ☐ Probability
- ☐ Both

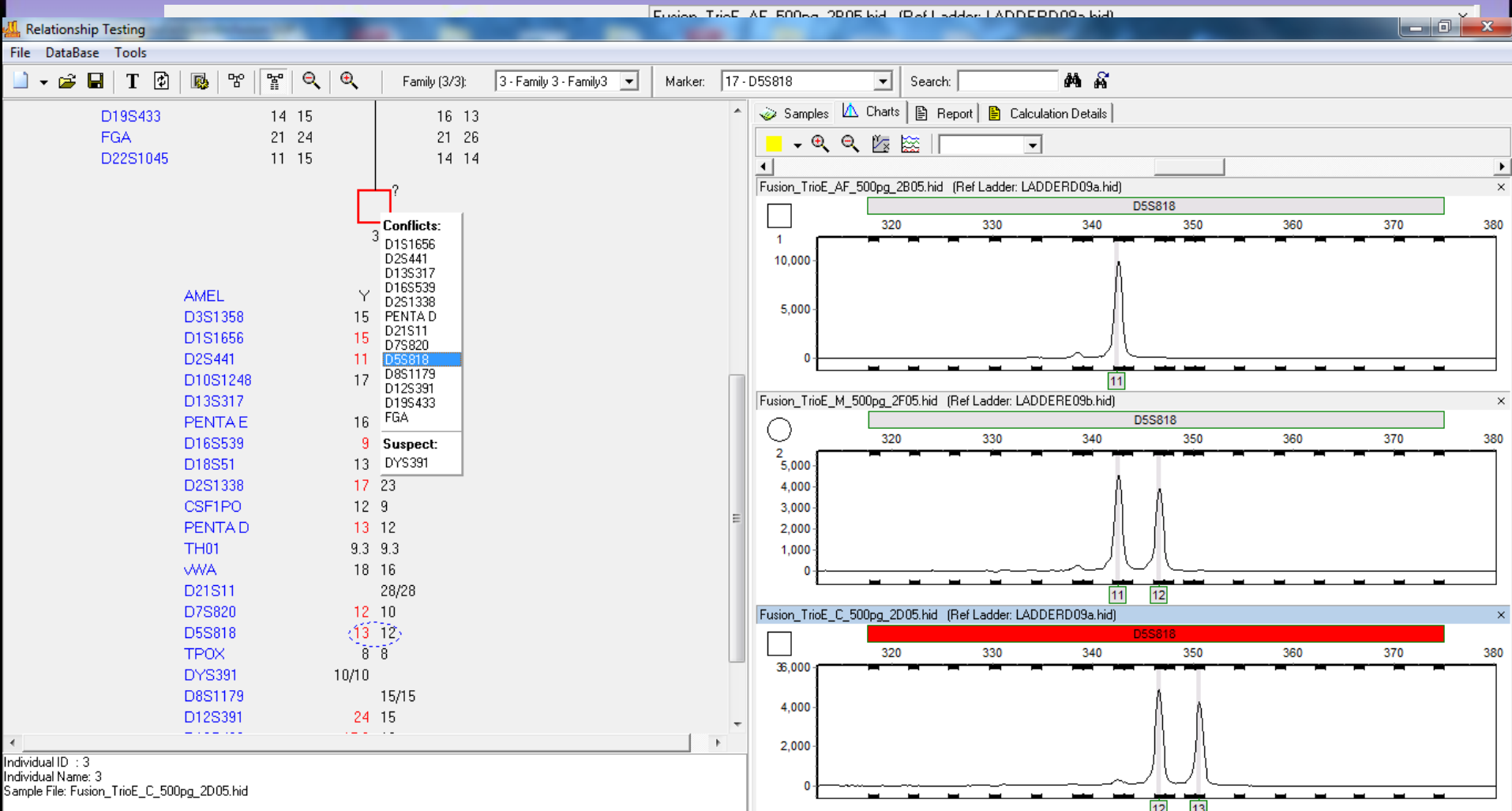
Ok Cancel

Automated Pedigree Drawing – Inclusion Example



Samples Charts Report Calculation Details							
	PI	PE	1	2	3	Numerator	Denominator
Total	3.59E+11	1.0000000					
D3S1358	1.57E+00	1.00E+00	15:16	15	15:16	9.47E-02	6.02E-02
D1S1656	5.36E+00	1.00E+00	16:16.3	15:15.3	15:16.3	5.45E-03	1.02E-03
D2S441	1.09E+01	1.00E+00	10:13	15	13:15	4.14E-03	3.82E-04
D10S1248	1.83E+00	1.00E+00	14:15	13:16	14:16	2.62E-02	1.43E-02
D13S317	2.30E+00	1.00E+00	12	9:11	9:12	9.42E-02	4.09E-02
PENTA E	1.19E+01	1.00E+00	13:17	15:18	17:18	2.56E-03	2.15E-04
D16S539	2.65E+00	1.00E+00	9:11	10:11	9:10	2.97E-02	1.12E-02
D18S51	4.50E+00	1.00E+00	15:18	15:17	17:18	8.38E-03	1.86E-03
D2S1338	3.87E+00	1.00E+00	21:22	20	20:21	1.83E-02	4.73E-03
CSF1PO	2.44E+00	1.00E+00	10:11	10:14	11:14	2.78E-02	1.14E-02
PENTA D	4.31E+00	1.00E+00	12:14	9:12	12	1.33E-03	3.09E-04
TH01	1.28E+00	1.00E+00	7:8	6:9.3	6:7	4.09E-02	3.20E-02
vWA	4.35E+00	1.00E+00	16:18	14:17	17:18	1.56E-02	3.58E-03
D21S11	6.33E+00	1.00E+00	30:31	28	28:31	1.38E-02	2.18E-03
D7S820	5.52E+00	1.00E+00	10:12	9:11	11:12	1.47E-02	2.65E-03
D5S818	2.16E+00	1.00E+00	8:13	11:12	12:13	6.61E-03	3.07E-03
TPDX	2.99E+00	1.00E+00	8	10:11	8:10	5.61E-02	1.88E-02
D8S1179	9.56E-01	1.00E+00	13:15	13:14	13:14	2.17E-02	2.27E-02
D12S391	1.27E+01	1.00E+00	17:22	18.3:23	18.3:22	2.87E-03	2.26E-04
D19S433	1.97E+00	1.00E+00	13:14	14:15	13:14	2.53E-02	1.28E-02
FGA	1.52E+00	1.00E+00	21:23	23:24	23:24	9.70E-03	6.38E-03
D22S1045	3.53E+00	1.00E+00	11:15	15:16	11:16	1.81E-02	5.11E-03

Automated Pedigree Drawing – Exclusion Example



Searchable Database - Relationship Testing Summary

Applications Include:

- Missing Persons, Mass Disaster, Paternity, Immigration, Human Trafficking**
- Linked to Main Analysis – No Data Transfer**
- Search Database for Exact Matches and Familial Search**
- Automates Repetitive Calculations LR using Identity by Descent**
- Kinship Analysis Across Three Generations**
- Immediate Visualization of Allele Conflicts in Pedigree Drawings**
- Paternity Index Calculations using Standards for Relationship Testing, AABB**

Mixture Analysis Application – Assists Forensic Experts

- No Data Transfer – Linked to Results of Genotyping Analysis
- ISFG and SWGDAM Guidelines and Recommendations
- Customize Analysis Parameters to Lab Standards
- Automates Repetitive Calculations: PI, PE, CPI, CPE, RMNE, LR

Run Wizard

Additional Settings - HID Analysis
Set additional options related to the different analysis type

Allelic Ladder:

Positive Control Template:

Allele Evaluation

Peak Score:

Reject < Check < Pass

☒ Mixture Evaluation

Valid Mixture Peak Percentage: %

Min Mixture Marker Number:

☒ Auto Select Best Ladder

☒ Auto Panel Adjustment

<< Back Ok Cancel



Good_edits_male_female_mixture.SGF
 Raw Data
 Allele Call
 ✓ MX: 0.6ng_1to1_MtoF_B06.fsa
 ? MX: 0.6ng_1to3_MtoF_E06.fsa
 ? MX: 0.6ng_1to9_MtoF_B07.fsa
 ? MX: 0.6ng_3to1_MtoF_H05.fsa
 ? MX: 0.6ng_9to1_MtoF_F05.fsa
 ✓ 0.6ng_Female_F07.fsa
 ✓ 0.6ng_Male_G04.fsa

Mixture Analysis

- One Contributor
- Two Contributors
 - 0.6ng_1to1_MtoF_B06
 - 0.6ng_1to3_MtoF_E06
 - 0.6ng_Female_F07
 - 0.6ng_Male_G04
 - 0.6ng_1to9_MtoF_B07
 - 0.6ng_3to1_MtoF_H05
 - 0.6ng_9to1_MtoF_F05
- Three or more Contributors

No.	Marker	Major	Minor	Major Mx	Residual	Major Hit
1 -- 1	D3S1358	16,17	14,15	0.69	0.0034	0.97
2 -- 1	TH01	6,9,3	6,9,3		0.0002	0.96
2 -- 2	TH01	6,9,3	6,Q	0.94	0.0251	0.98
2 -- 3	TH01	6,9,3	9,3,Q	0.94	0.0309	0.90
2 -- 4	TH01	6,9,3	6,6	0.98	0.0334	
3 -- 1	D21S11	31,32,2	27,28	0.69	0.0012	0.94
4 -- 1	D18S51	12,12	13,14	0.74	0.0008	0.85
5 -- 1	Penta_E	10,21	7,16	0.71	0.0069	0.72
6 -- 1	D5S818	11,12	11,11	0.70	0.0002	
6 -- 2	D5S818	11,11	12,12	0.65	0.0100	
6 -- 3	D5S818	11,12	11,Q	0.90	0.0270	0.60
7 -- 1	D13S317	12,13	10,10	0.75	0.0025	0.89
7 -- 2	D13S317	10,12	13,13	0.65	0.0198	0.62
7 -- 3	D13S317	12,13	10,12	0.59	0.0219	0.66
7 -- 4	D13S317	10,13	12,12	0.60	0.0267	0.70
7 -- 5	D13S317	12,13	10,13	0.62	0.0339	0.55
8 -- 1	D7S820	10,12	8,11	0.62	0.0113	0.88
9 -- 1	D16S539	11,14	9,13	0.68	0.0037	0.84
10 -- 1	CSF1PO	10,11	11,13	0.72	0.0010	0.91
11 -- 1	Penta_D	9,9	13,13	0.74	0.0007	
11 -- 2	Penta_D	9,13	9,9	0.52	0.0192	
11 -- 3	Penta_D	9,9	9,13	0.48	0.0298	
12 -- 1	AMEL	X,X	X,Y	0.66	0.0020	
12 -- 2	AMEL	X,X	Y,Y	0.83	0.0236	
13 -- 1	vWA	18,18	14,14	0.78	0.0080	
13 -- 2	vWA	18,18	14,18	0.57	0.0117	
14 -- 1	D8S1179	12,14	10,12	0.75	0.0034	0.86
14 -- 2	D8S1179	12,14	10,14	0.77	0.0198	0.68
14 -- 3	D8S1179	12,14	10,10	0.87	0.0336	0.87
15 -- 1	TPOX	8,8	8,10	0.66	0.0019	
15 -- 2	TPOX	8,8	10,10	0.83	0.0239	
16 -- 1	FGA	20,21	20,24	0.77	0.0025	0.93
16 -- 2	FGA	20,21	21,24	0.80	0.0252	0.67



Contributor 1: 0.6ng_Female_F07 (Major) ☐ Contested

Contributor 2: 0.6ng_Male_G04 (Minor) ☒ Contested

Mixture	PI	PE	Contributor 1	Contributor 2	LR
14,15,16,17	0.84456	0.15544	16,17	14,15	18.814
6,9,3	0.04368	0.95632	6,9,3	6,9,3	4.910
27,28,31,32,2	0.17556	0.82444	31,32,2	27,28	56.922
12,13,14	0.03497	0.96503	12	13,14	126.582
7,10,16,21	0.03423	0.96577	10,21	7,16	122.730
11,12	0.36240	0.63760	11,12	11	6.384
10,12,13	0.37946	0.62054	12,13	10	41.254
8,10,11,12	0.72932	0.27068	10,12	8,11	12.566
9,11,13,14	0.44756	0.55244	11,14	9,13	20.222
10,11,13	0.29702	0.70298	10,11	11,13	40.980
9,13	0.07453	0.92547	9	13	23.737
X,Y	--	--	X	X,Y	--
14,18	0.03276	0.96724	18	14,18	51.188
10,12,14	0.20340	0.79660	12,14	10,12	47.456
8,10	0.16322	0.83678	8	8,10	19.611
20,21,24	0.10240	0.89760	20,21	20,24	19.593
Cumulative:	8.58E-13	1-8.58E-13			4.96E+21
PowerPlex 16: US Asian-American					
14,15,16,17	0.79924	0.20076	16,17	14,15	46.296
6,9,3	0.07508	0.92492	6,9,3	6,9,3	3.931
27,28,31,32,2	0.08703	0.91297	31,32,2	27,28	347.222
12,13,14	0.13396	0.86604	12	13,14	26.966
7,10,16,21	0.03725	0.96275	10,21	7,16	163.827
11,12	0.32376	0.67624	11,12	11	4.806
10,12,13	0.38316	0.61684	12,13	10	8.334
8,10,11,12	0.79388	0.20612	10,12	8,11	10.437
9,11,13,14	0.40960	0.59040	11,14	9,13	18.222
10,11,13	0.27248	0.72752	10,11	11,13	100.827
9,13	0.21437	0.78563	9	13	11.505
X,Y	--	--	X	X,Y	--
14,18	0.10240	0.89760	18	14,18	16.235
10,12,14	0.16810	0.83190	12,14	10,12	9.561
8,10	0.06656	0.93344	8	8,10	15.081
20,21,24	0.08880	0.91120	20,21	20,24	22.522
Cumulative:	6.25E-12	1-6.25E-12			1.30E+20
PowerPlex 16: US Caucasian					
14,15,16,17	0.68558	0.31442	16,17	14,15	14.162

Comment:

Contesting minor contributor (Male sample) LR displayed for results using each of the population allele frequencies

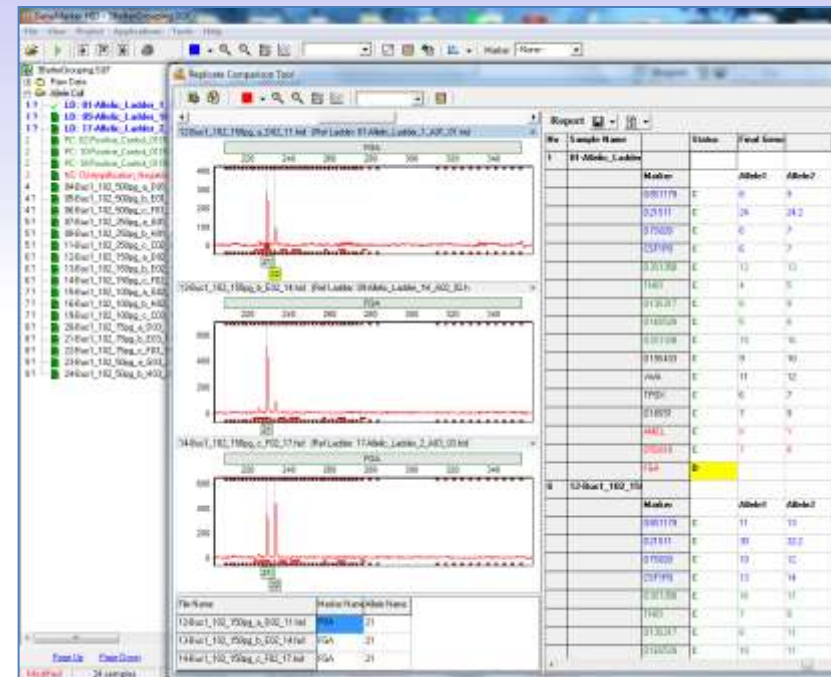
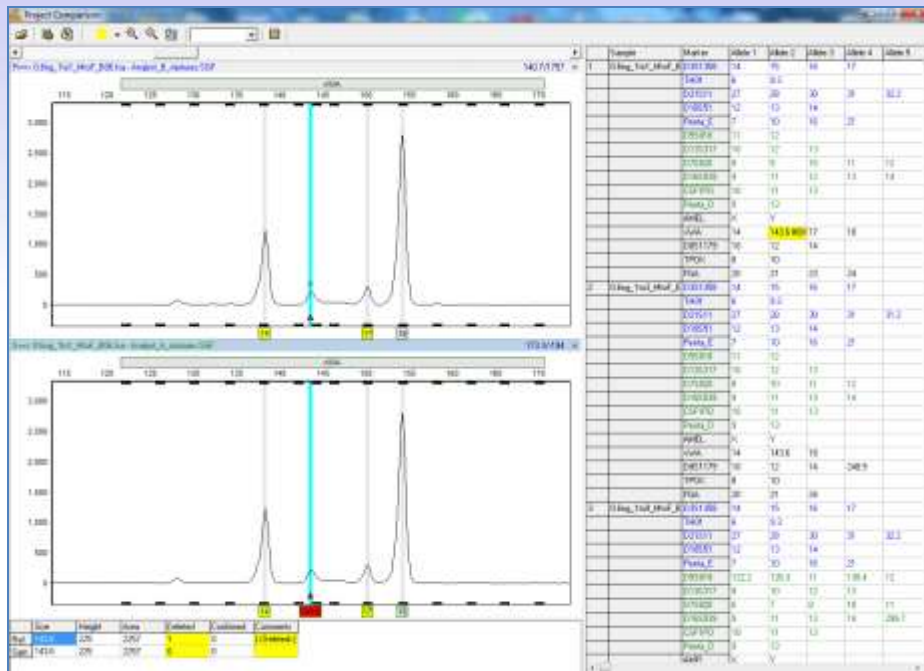
Mixture Analysis Assistant Summary

- **Developed using recommendations of the DNA Commission of the International Society of Forensic Genetics, compatible with SWGDAM Guidelines**
- **Linked to Main Analysis – No Data Transfer Needed**
- **Automates Repetitive Calculations**
- **Calculates PI, PE, LR, RMNE**
- **Database search for contributor; with or without a reference sample**

Quality Control and Validation Tools:

Project Comparison

Replicate Comparison



Highlights any Differences – Table is Linked to Electropherograms

IN CONCLUSION, GeneMarker HID:

Expert System

- Documented Time Savings
- Decrease Backlogs
- Concordant / Validated

Searchable Database / Relationship Testing

- Linked to Main Analysis – No Manual Data Transfer
- Missing Persons, Mass Disaster Applications
- Locate Same Individual or Family Members (IBD)
- Kinship Across Three Generations (IBD)
- Paternity/Immigration Pedigrees and Paternity Index (AABB relationship testing guidelines appendix 8, trios and single parent cases)

Mixture Analysis

- Linked to Main Analysis – No Manual Data Transfer
- Customize Analysis Parameters
- Automate Repetitive Calculations, PE, PI, RMNE, LR
- Search Database – Calculate LR

Thank you for attending this presentation on GeneMarker HID!

Acknowledgments: we are grateful to the following people for sharing their thoughts and experience to assist us during development

Dr. Mitchell Holland, Penn State University Forensics Program

Drs. Thomas Reid and Michael Baird, DNA Diagnostics Center

Drs. John Butler, Kristen O'Connor, Pete Valone and Michael Coble, NIST

Drs. Rhonda Roby, John Planz, Arthur Eisenburg, and Nicole Philips, UNT CHI

Drs. Andy Hopwood and Bob McLaren, Promega, for providing data