

PowerPlex® Fusion

Overview & Developmental Validation Preliminary Summary

Ann MacPhetridge, Marketing Manager

October 2012



PowerPlex® Fusion System

Webinar outline

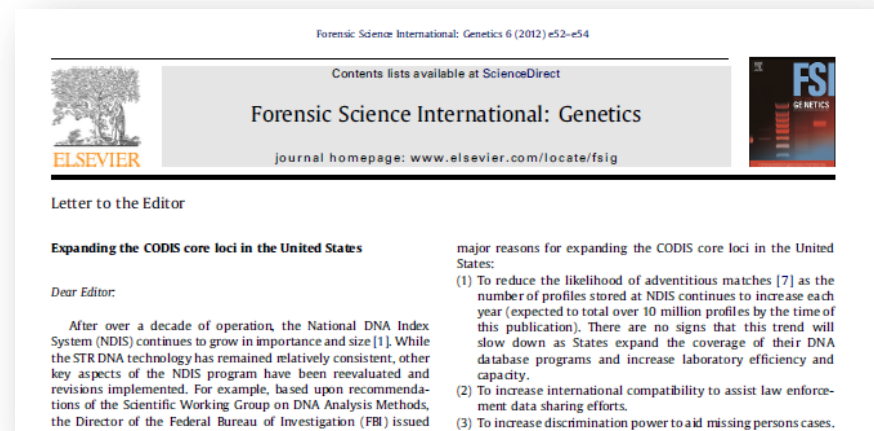
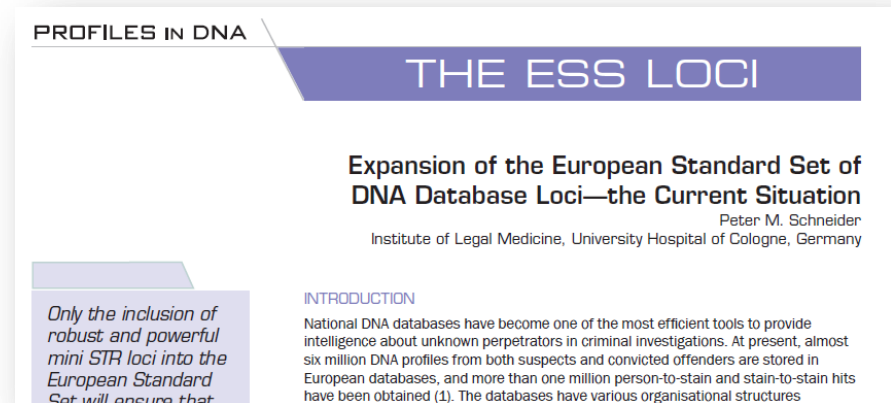


- Expansion of CODIS
- Overview of the PowerPlex® Fusion System
- PowerPlex® Fusion Developmental Validation Data – a summary of work performed to-date by Promega

Expansion of Databases... Expansion of Multiplexes



- ✓ Need for common loci is recognized
- ✓ European Standard Set has expanded
- ✓ CODIS will expand in the coming years
- ✓ Goals:
 - More data sharing
 - Fewer adventitious matches



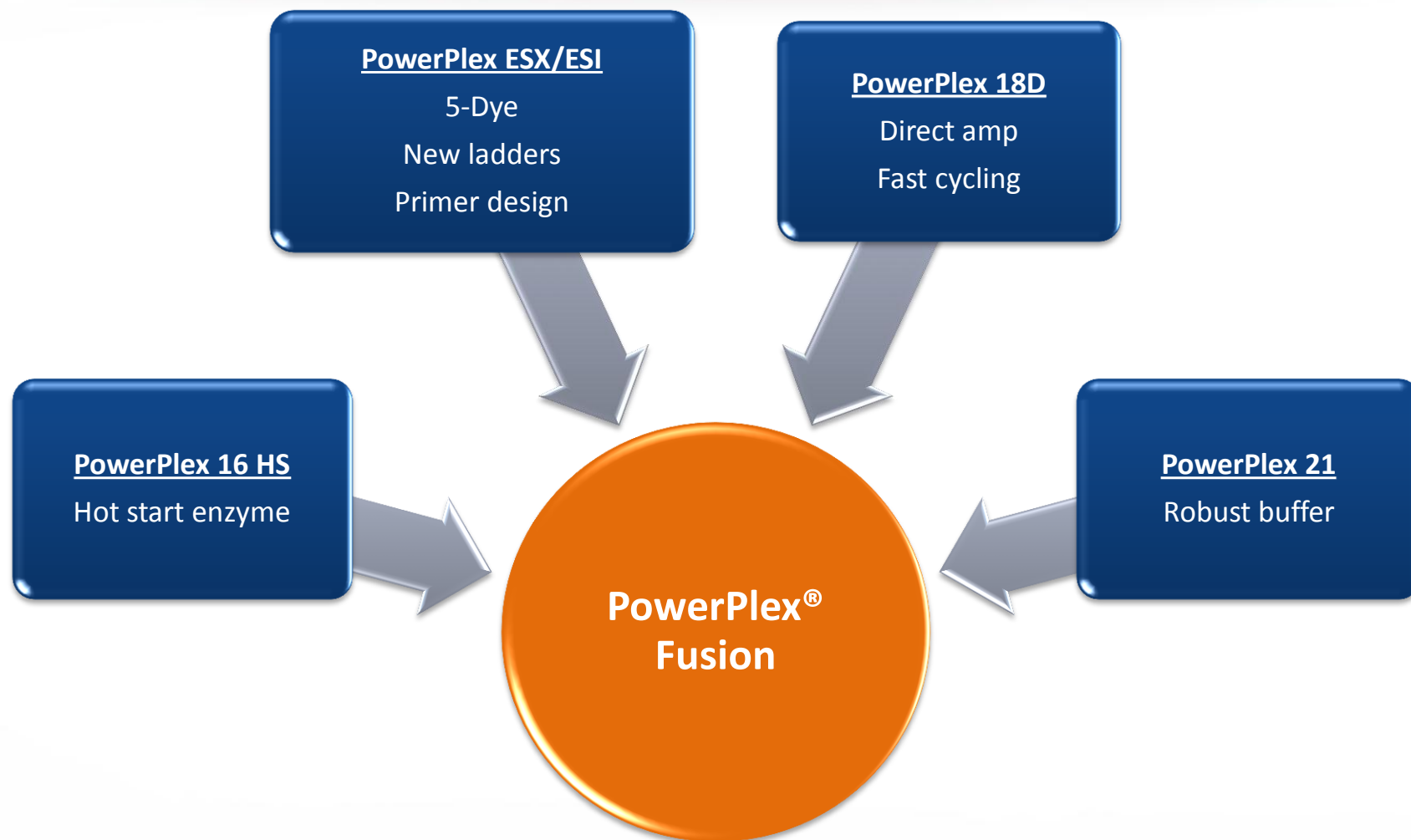
Supporting Expansion of Core Loci Product Requirements



1. Include all current CODIS and ESS loci, plus Amelogenin
 - US SWGDAM: include DYS391
2. Support current CE instruments and software
 - 5-dye chemistry proven on the 31xx and 3500 Genetic Analyzer
3. Enable fast cycling
4. Create one kit for all labs to simplify validation, purchasing, etc.
 - Provide database (i.e. direct amp) and casework protocols

Development Approach

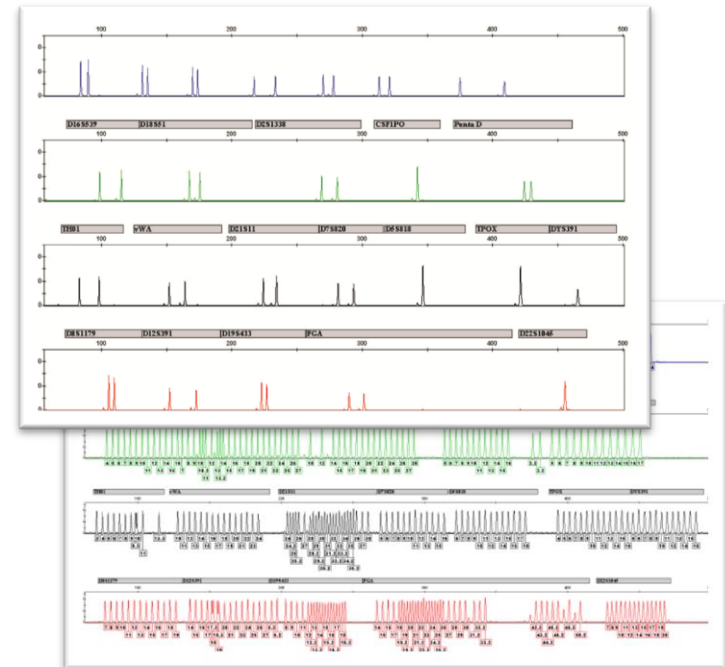
Building on Proven Technology



What is the PowerPlex® Fusion System?

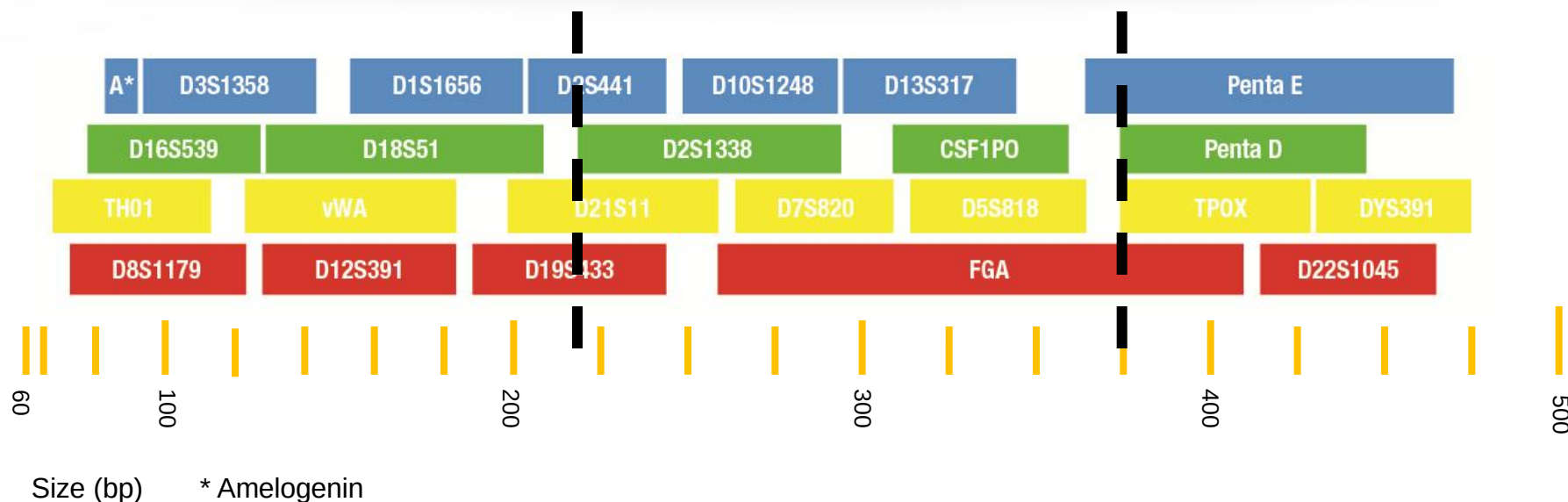


- Analyzes 24 loci in a single reaction
 - CODIS + ESS + Pentas + D2S1338 + D19 +DYS391 + Amelogenin
 - Highest discrimination
- Uses rapid PCR technology (~85 min)
- Offers casework *and* database protocols ***in one kit***
 - One kit to validate, purchase & QC
- Designed to work with 31xx CE
 - No 6-dye upgrade required



PowerPlex® Fusion System

Locus placement



Most essential loci located below ~400bp; nine under 220bp

- Excluding “extended” FGA alleles, the most informative required loci are <375bp
- All required “new CODIS loci” (“Panel A”)
- 2 of 3 “recommended” loci, TPOX and D22 (“Panel B”) -- no SE33
- Pentas

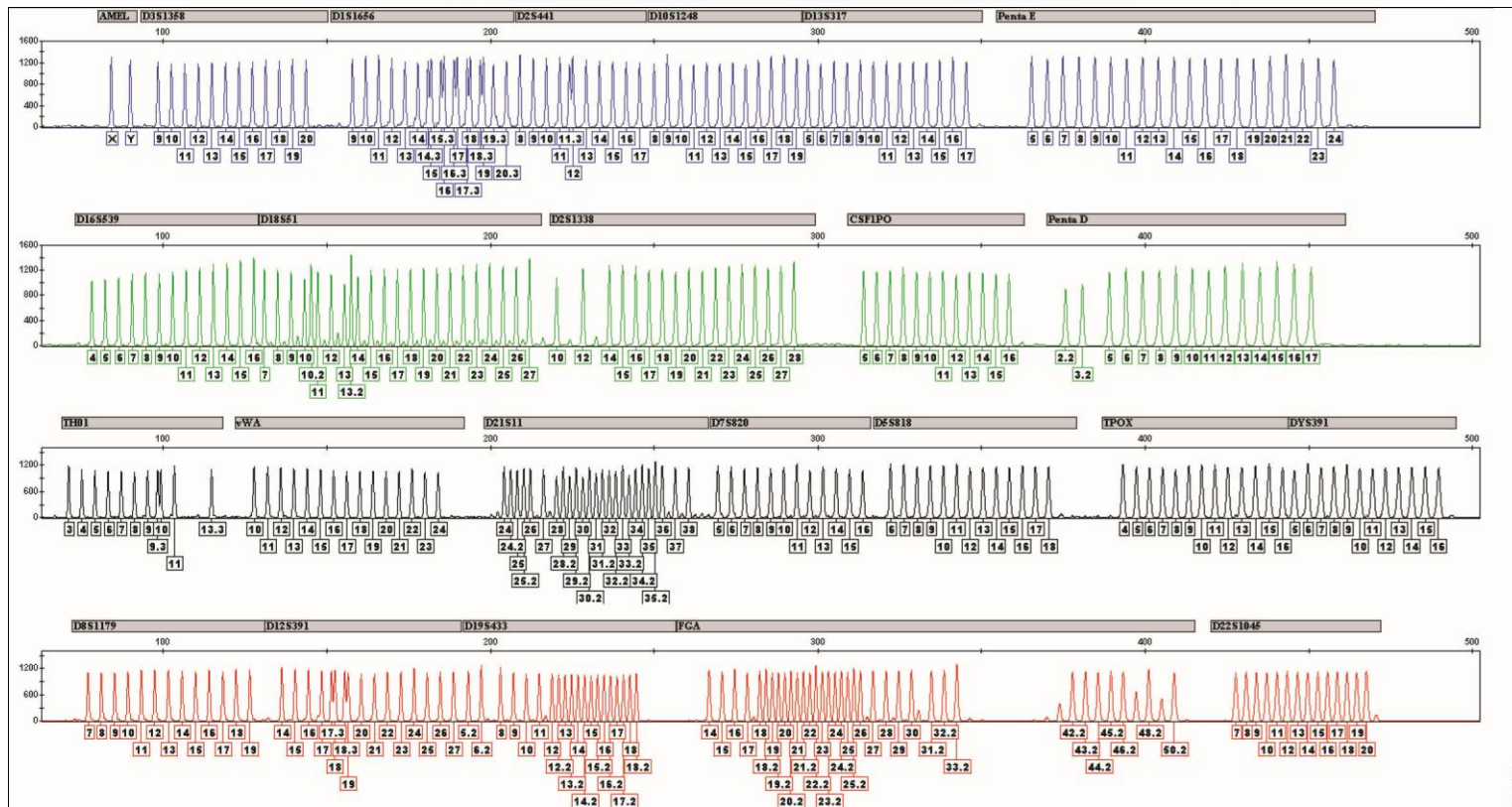
Locus Selection

Inclusion of Pentas



- Straightforward inclusion in multiplexes because of “size range” (allele range)
- Commercially-available for more than a decade
- Easy mixture interpretation due to low stutter
- High discrimination added by two highly polymorphic loci
- Used routinely in Asia, Europe, North and South America
 - >1M samples in US NDIS contain Pentas

Enhanced allelic ladder



- 36 new alleles compared to PowerPlex® 16
- New virtual bins to increase genotyping efficiency
- Fewer inconclusive calls

Configuration and protocol



Product Overview—Configuration



Kit sizes

- 200 reactions (DC2402)
- 800 reactions (DC2408)



Components

- PRE-AMP: 5X Master Mix (w/ hot start *Taq*), 5X Primer Mix, 2800M Control DNA, Amplification Grade Water
- POST-AMP: Size standard (ILS 500), Allelic Ladder

Related Products

- 5-dye Matrix for AB 31xx & 3500 (DG4700)
- GeneMapper ID & ID-X panel/bin/stutter file downloads (free)

Supported Instruments



Thermal Cycler:

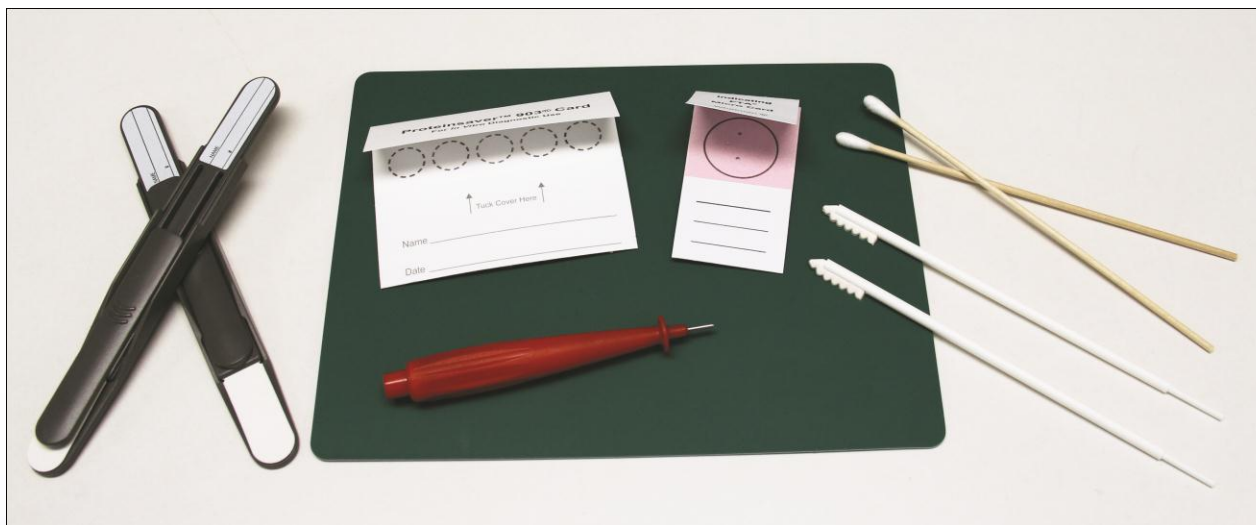
- GeneAmp 9700

Capillary Electrophoresis:

- 31xx Genetic Analyzer
- 3500 Genetic Analyzer

Usable Sample Types

- Extracted DNA
- Direct amplification from solid support materials
 - FTA[®] and Indicating FTA[®] cards
 - Swabs – Cotton, Omni Swab [treated with SwabSolution[™]]
 - Bode Buccal DNA Collector[™] [treated with PunchSolution[™]]
 - S&S 903 Specimen Collection paper [treated with PunchSolution[™]]



PowerPlex® Fusion Protocol



PowerPlex® Fusion System

Extracted DNA
OR punch
OR swab extract

Reaction Assembly

- 5X Master Mix
- 5X Primer Mix
- 15µl volume available

Amplification: <1.5 hours

1 cycle	96°C for 1 min
30	94°C for 10 seconds
cycles:	59°C for 1:00min
	72°C for 30 seconds
	60°C for 10 minutes
	4°C soak

Detection and Analysis

- 5 Dye Chemistry (G5)
- 3130/3130xl or 3500/3500xl

Protocols

Extracted DNA, FTA[®] punches, NonFTA punches, Swabs



Extracted DNA

Extract DNA
from sample

Add up to 15µl
of sample into
reaction

FTA punches

Punch FTA[®]
into plate

Add reaction
mix

NonFTA Punches

Punch card into
plate, treat with
PunchSolution[™]

Add reaction
mix to dried
punch

Swabs

SwabSolution[™]
incubation of
swab

Add 2µl of
extract into
reaction mix

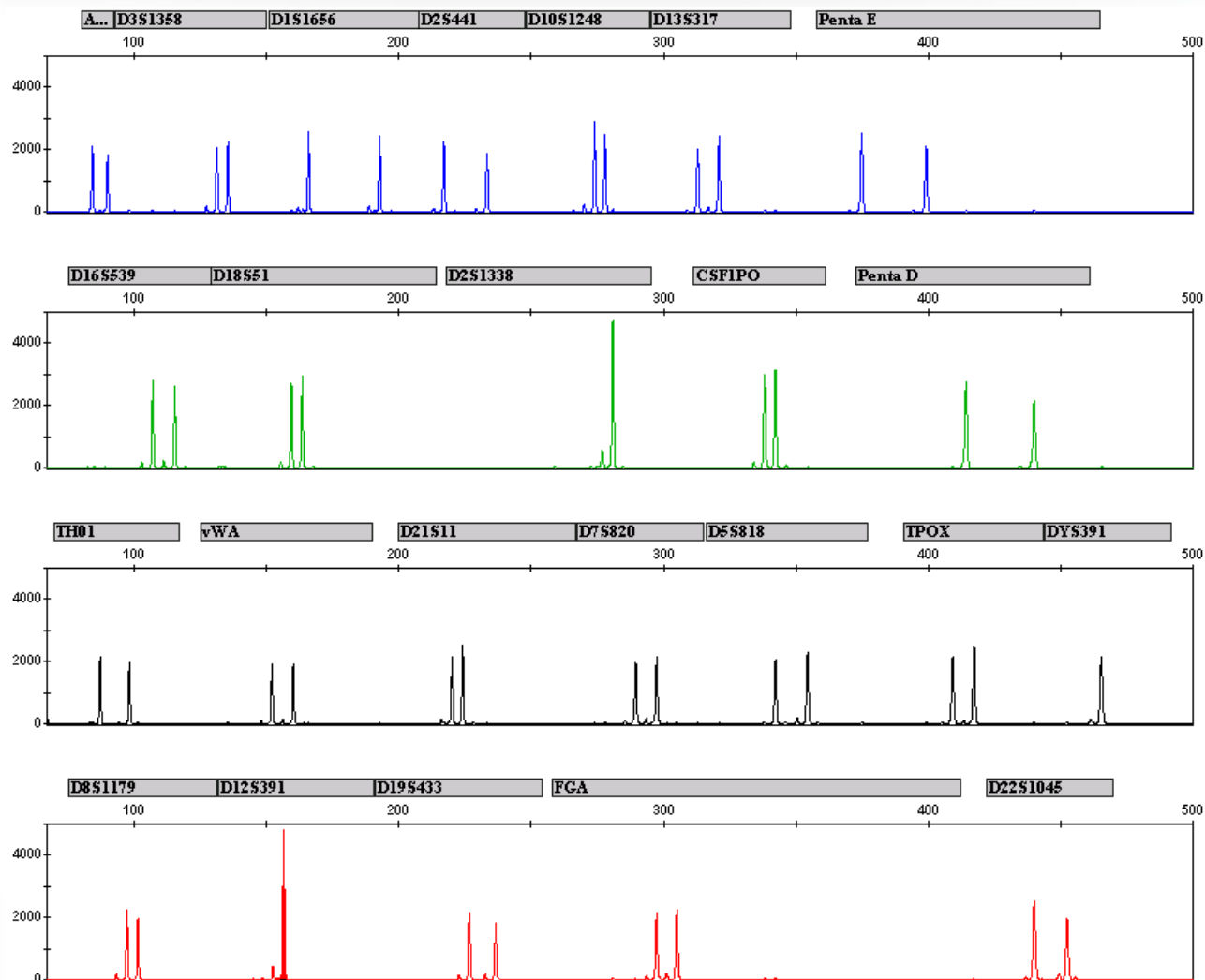
Cycling time ~85 minutes for all applications

Extracted DNA

0.5ng

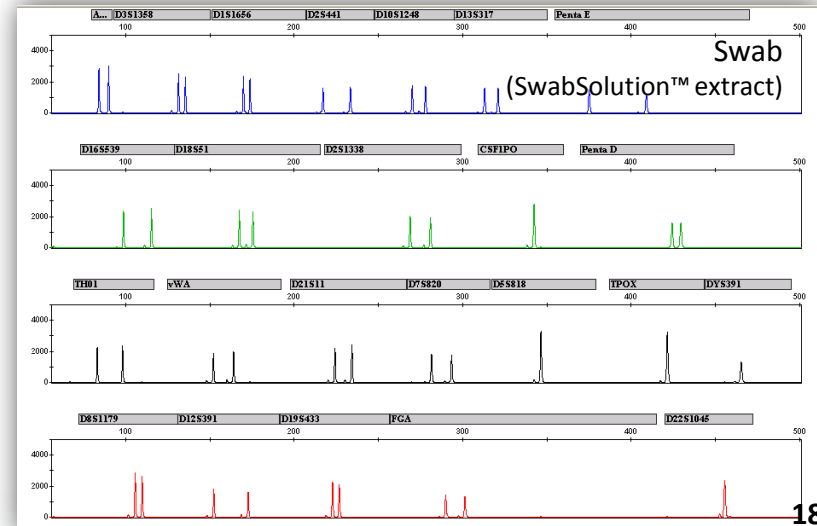
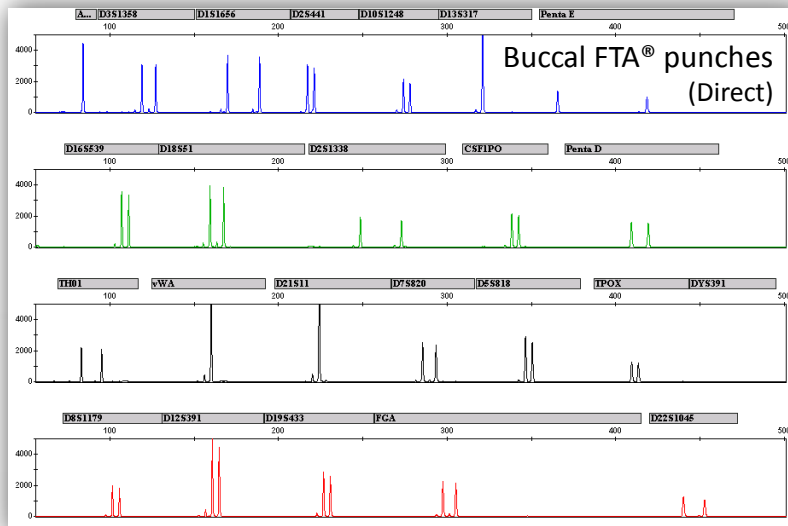
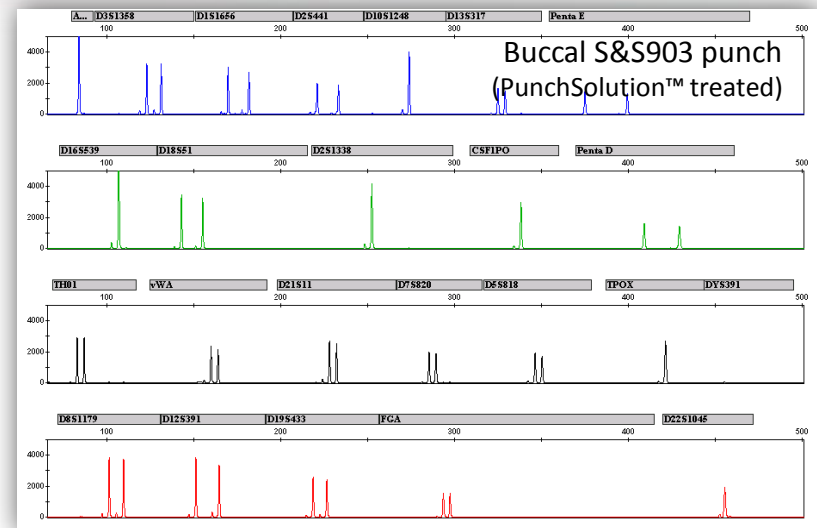
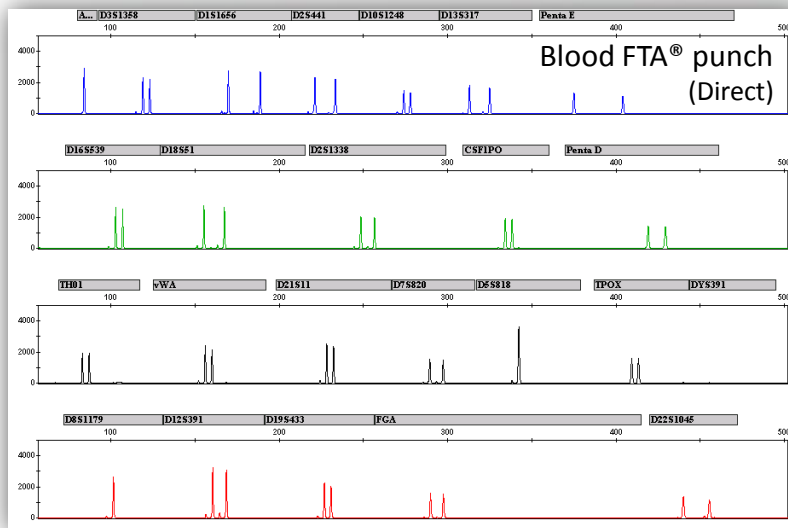


3130xl



3kv 5s

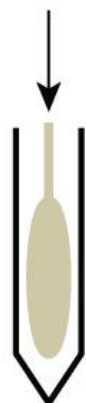
Example “Direct Amp” Data



Buccal Swab Workflow

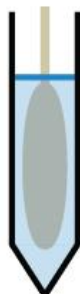


Add buccal swab
to tube or plate*.



Add 1ml of
SwabSolution™
Reagent.

Incubate 30 minutes
at 70°C**.



Transfer 2µl of
swab extract.

Add to 23µl of PCR
amplification mix‡.

Perform thermal
cycling.

Simple, rapid process for all swabs

* For plate format, use 2.2mL deep well plate (V6781)

** Use new heat transfer block (A2661)

‡ For reactions other than PowerPlex® 18D and PowerPlex® 21, add AmpSolution™ Reagent

NonFTA[®] punch workflow

S&S903, Bode Buccal DNA Collector[™] Device



Punch card
into well.



Add 10µl of
PunchSolution[™]
Reagent.



Incubate 30 minutes
at 70°C (to dryness).

Add PCR
amplification mix.

Perform thermal
cycling.

10583MB

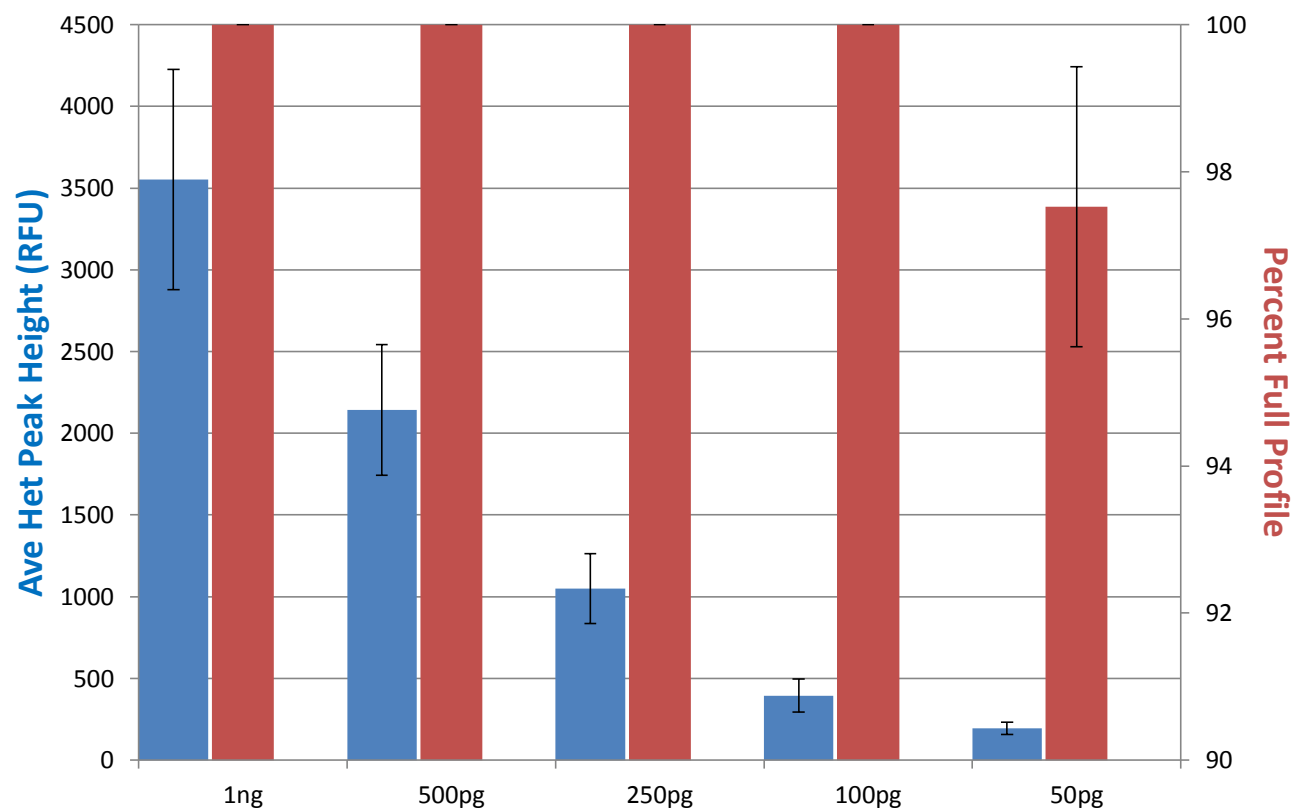
Rapid, same-plate processing of non-FTA[®] punches

Performance



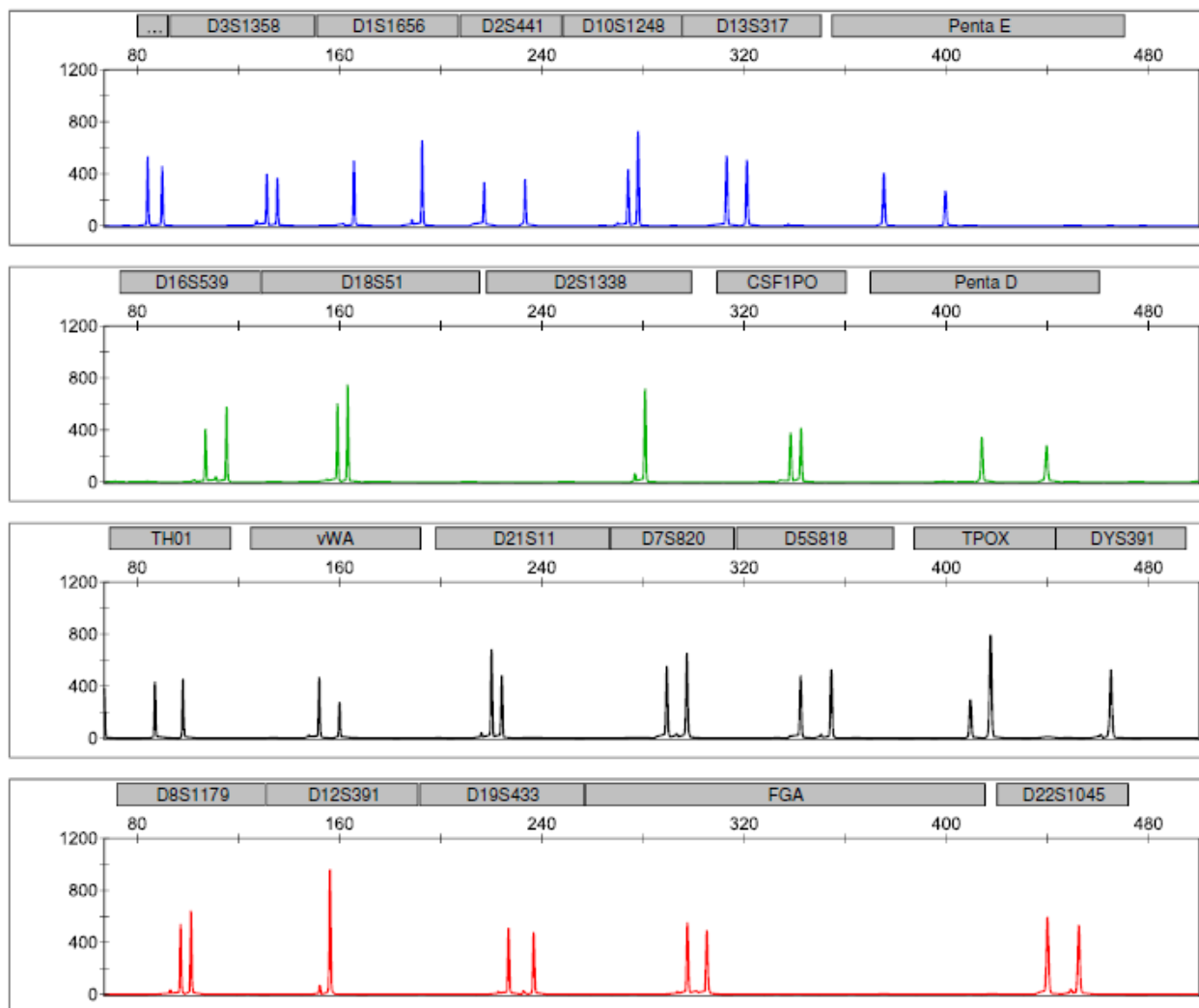
Sensitivity

Titration of Purified DNAs

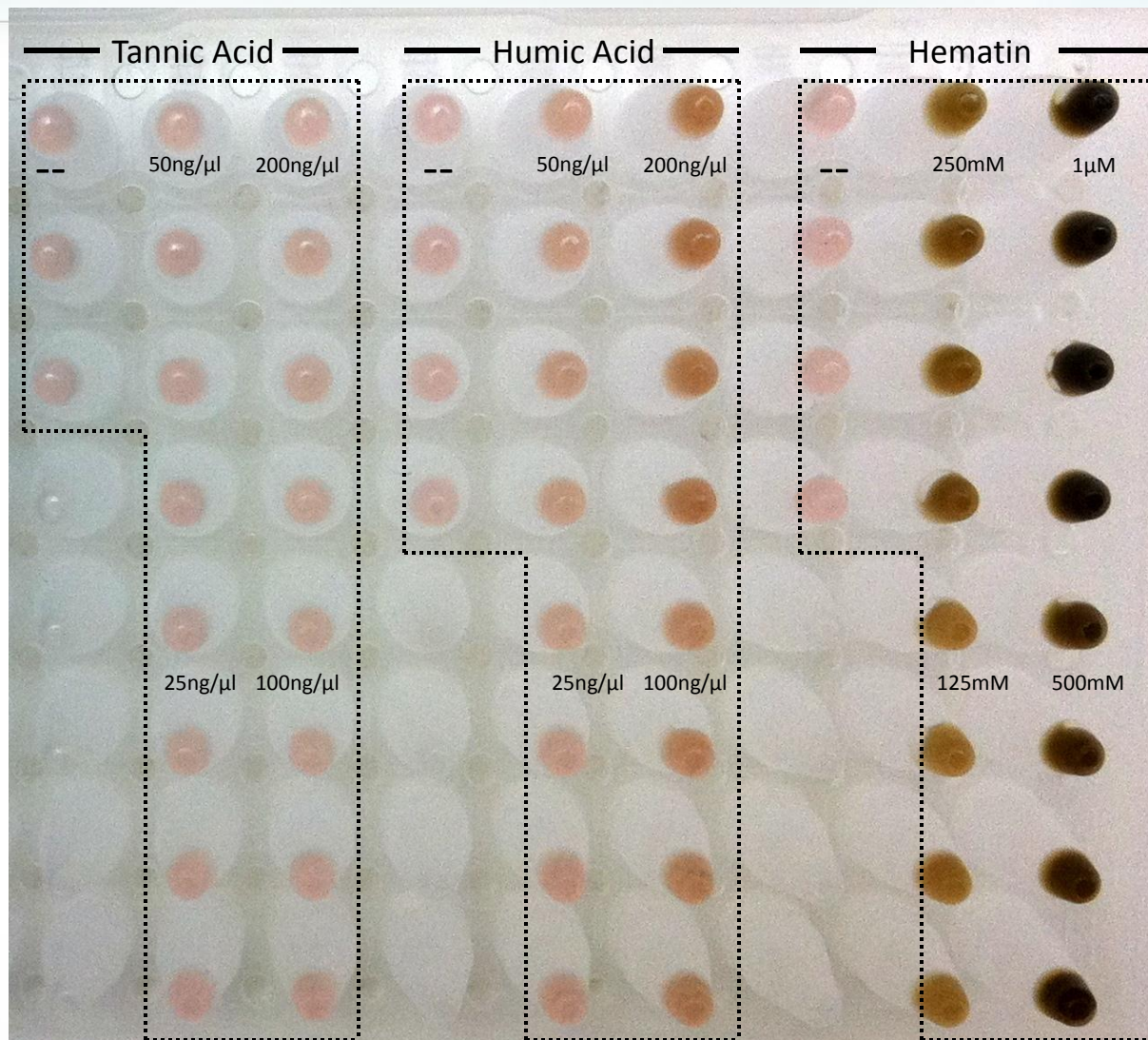


Sensitivity

100pg Amplification

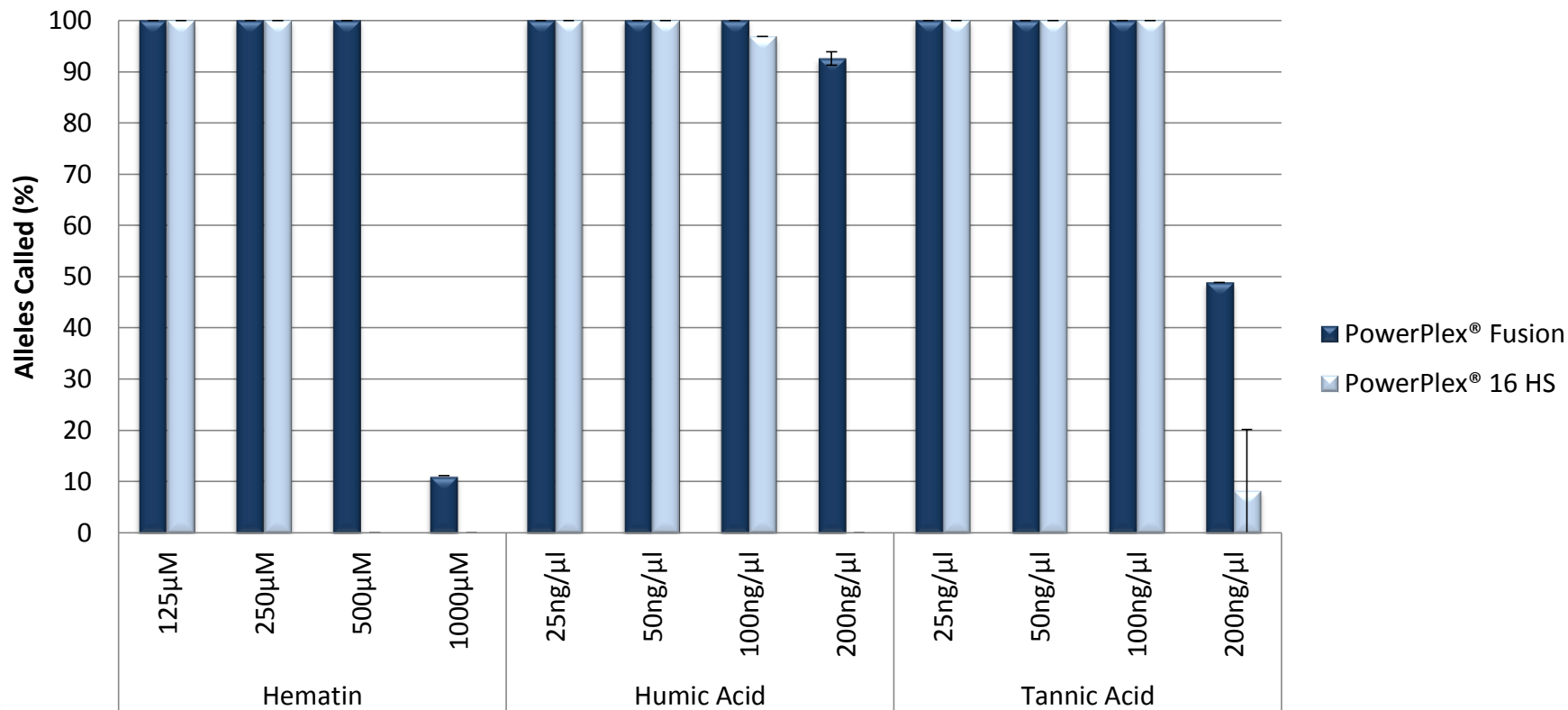


PowerPlex® Fusion Inhibitor Titration



Inhibitor Tolerance

Titration of Common Forensic PCR Inhibitors



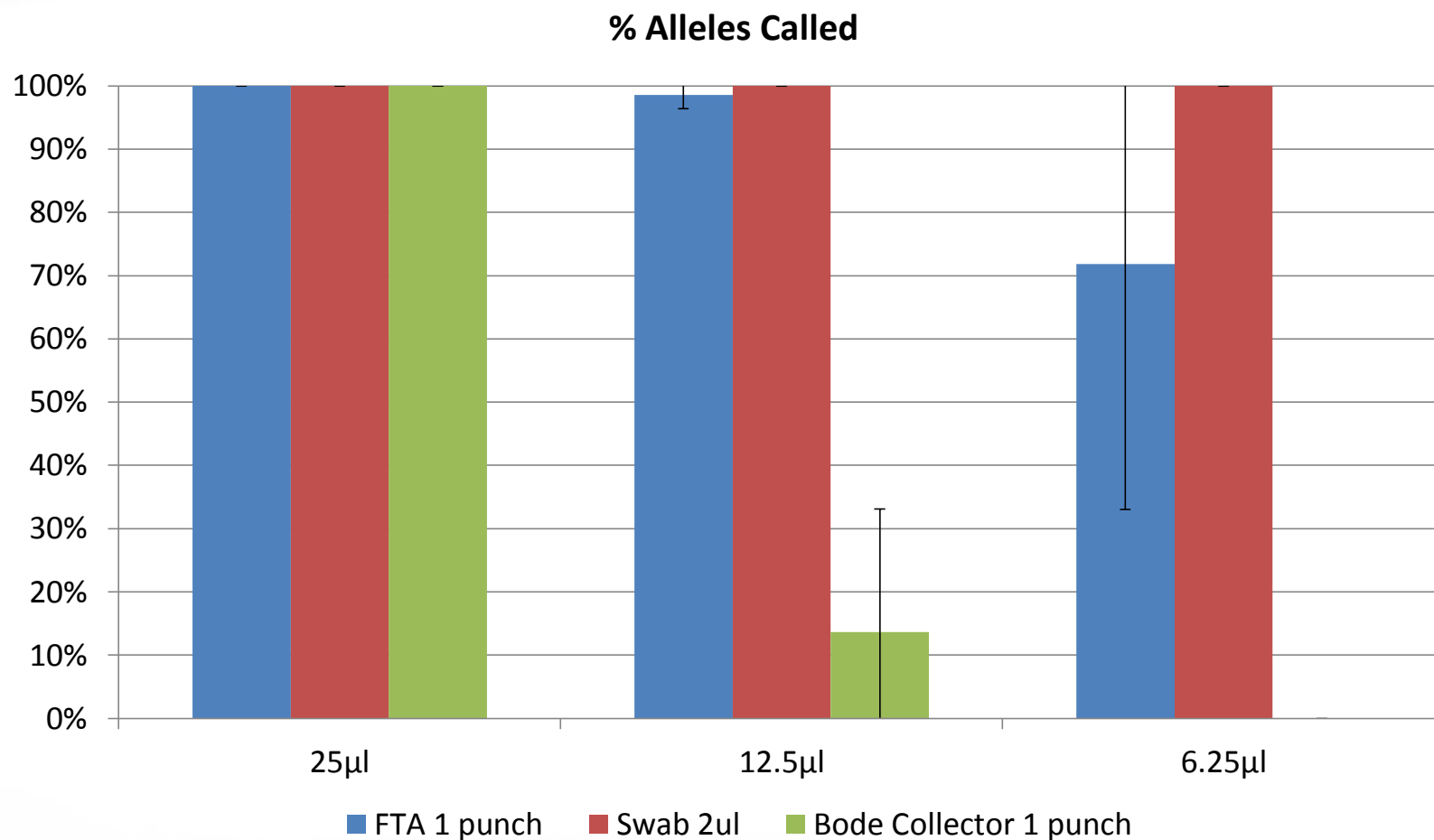
500pg of DNA (n=3) was amplified with each kit in the presence of the listed inhibitors.

***PowerPlex® Fusion
Developmental Validation Data***



Developmental Validation – Preliminary data

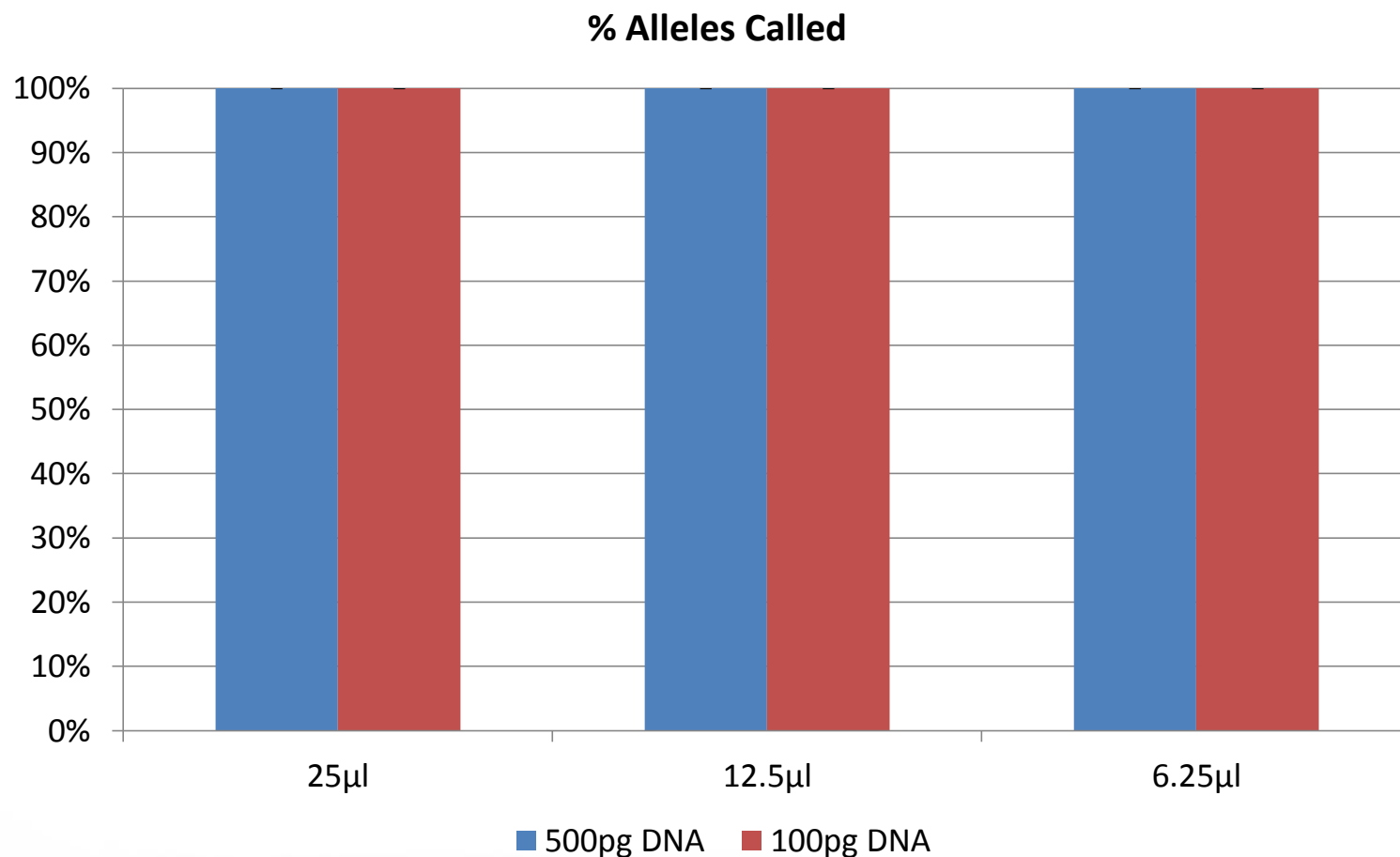
Reaction Volume – Solid Support Samples



N = 3

Developmental Validation – Preliminary data

Reaction Volume – Extracted DNA



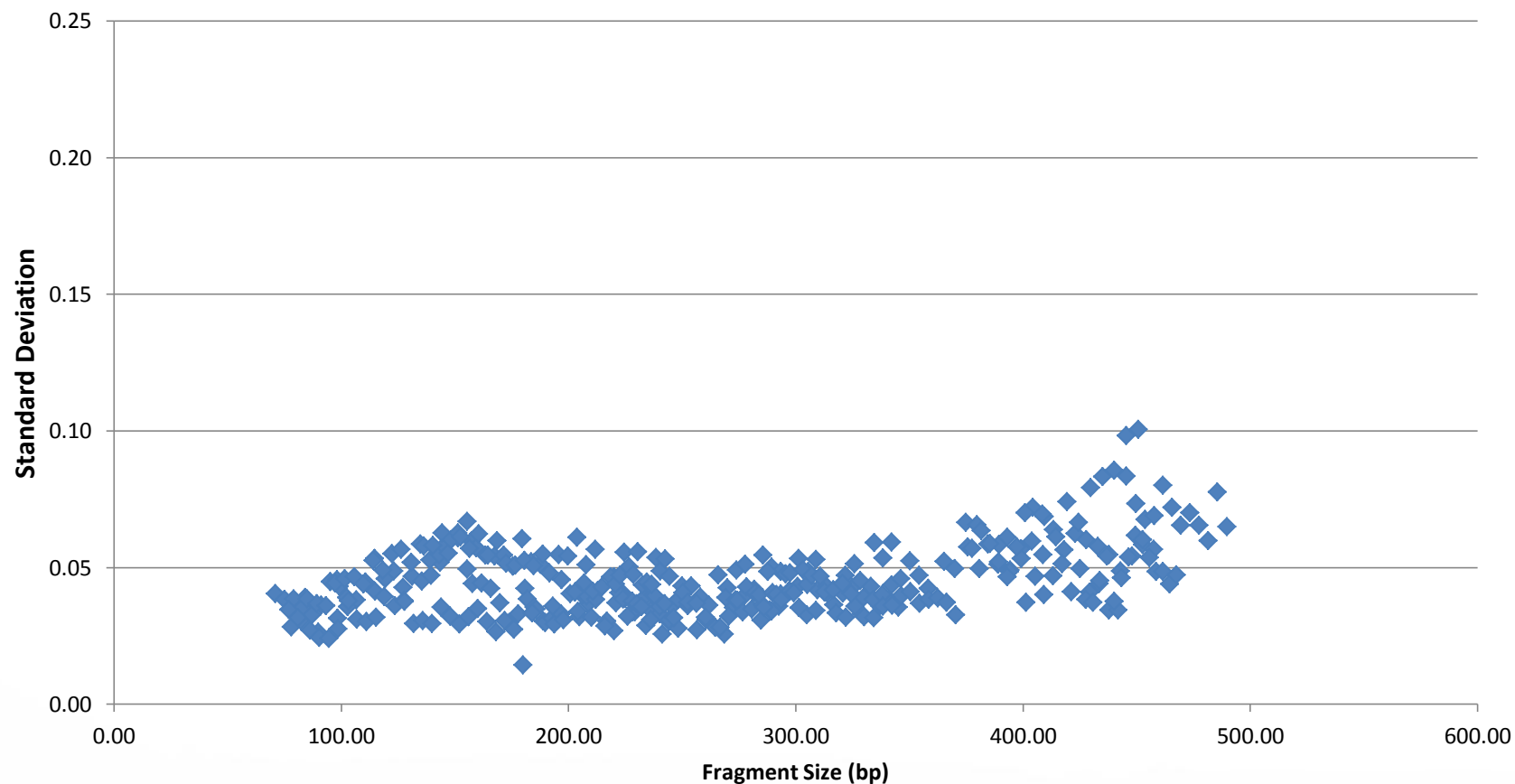
N = 3

Developmental Validation – Preliminary data

Precision Studies AB 3130xl



3130xl POP-4 Precision - 32 Ladders

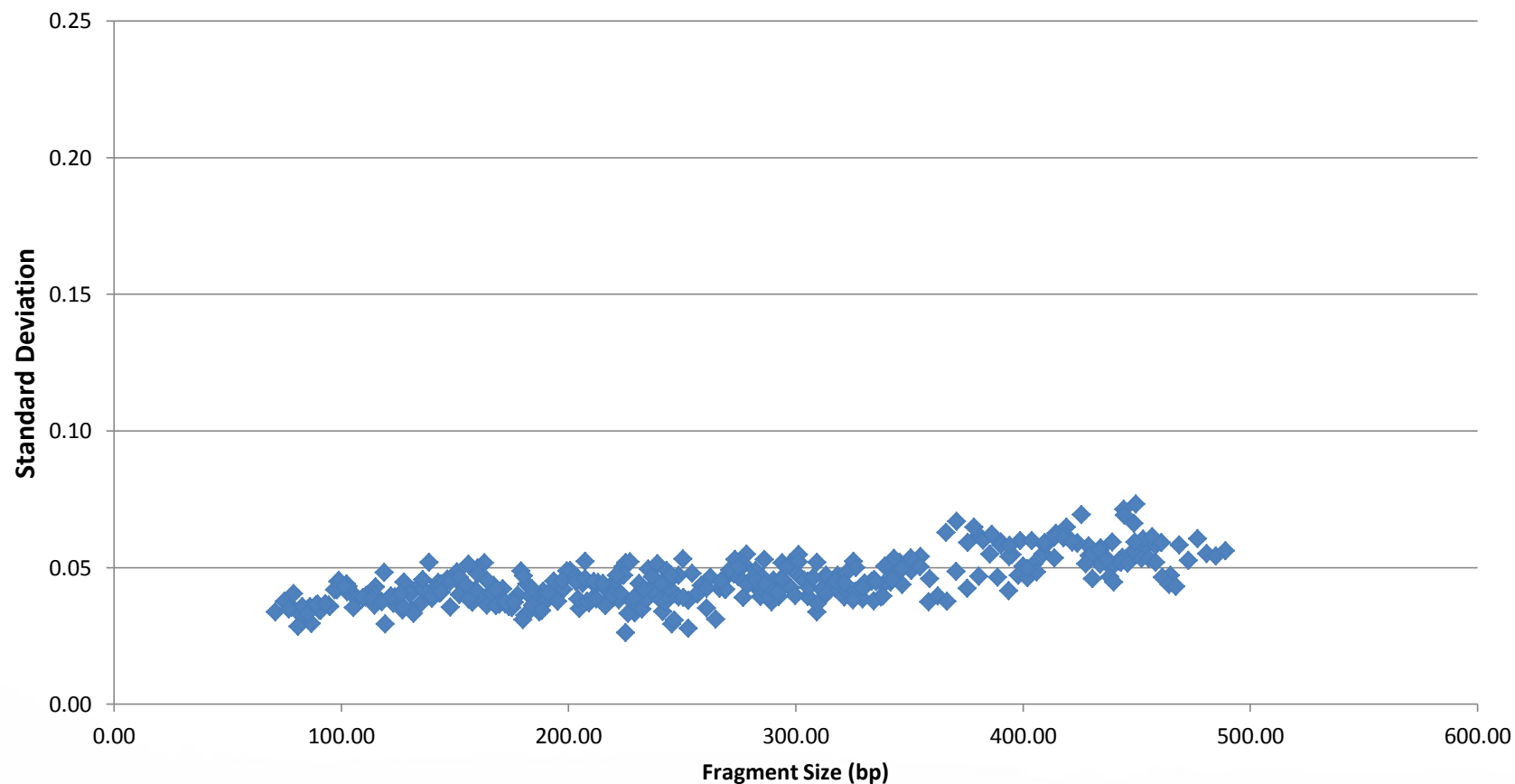


Developmental Validation – Preliminary data

Precision Studies AB 3500xl



3500xl POP-4 Precision - 48 Ladders



Developmental Validation

Population Study - NIST Concordance Testing



- PowerPlex® Fusion results compared to all other kits tested including:
Sinofiler/NGM/Identifiler/Yfiler/IDplex/ESSplex/PP16/PP21 kits with **652 unrelated individuals** (NIST U.S. population set)
- Fusion is **fully concordant with NIST SRMs 2391b&c** certified values
- No PP Fusion **null alleles**
- No PP Fusion **discordance** with other PowerPlex kits, discordance with ABI or Qiagen kits is on their end and are previously documented

Developmental Validation

Population Study - Probability of Identity Values



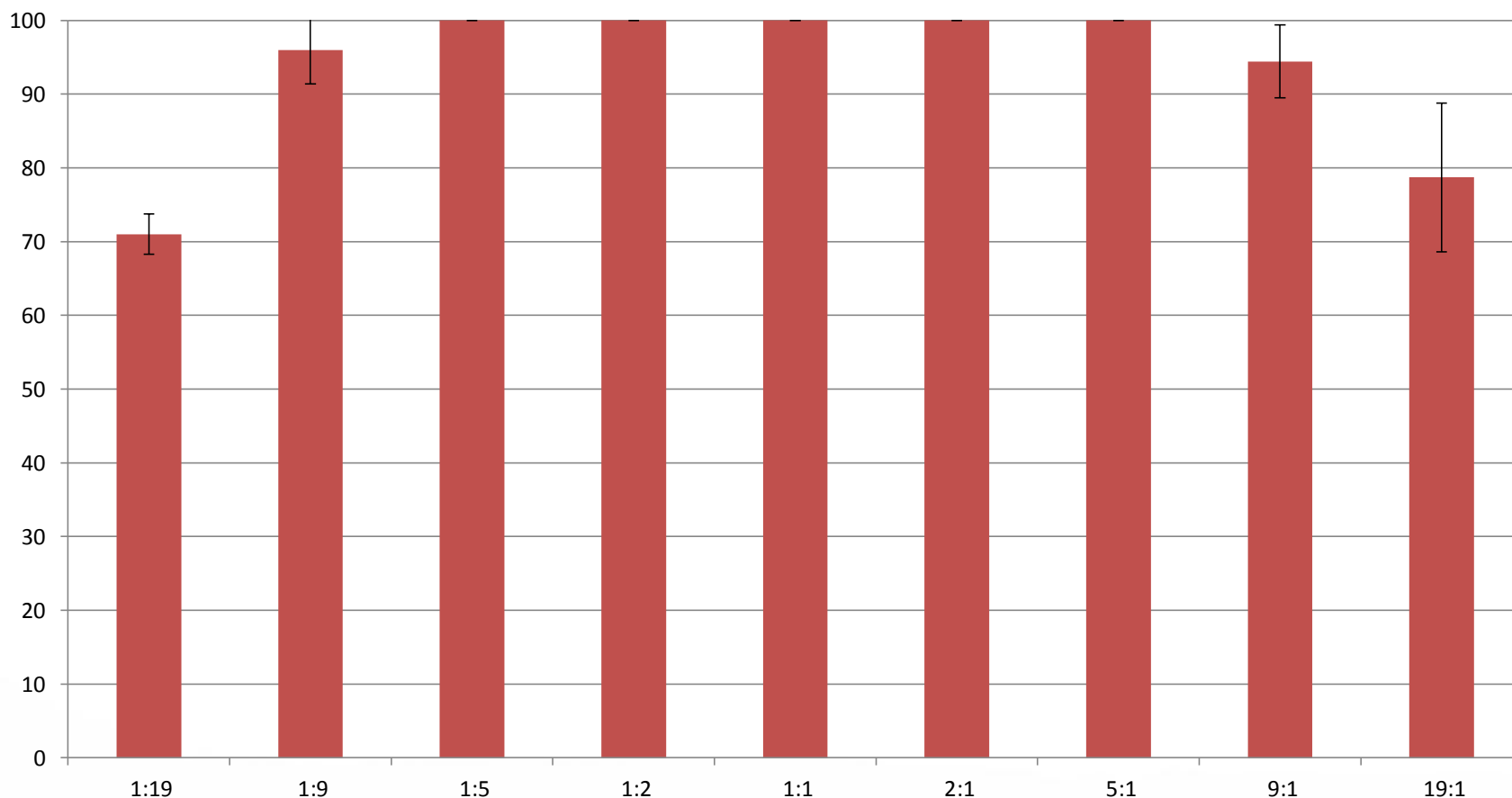
STR Typing Kits (Locus Combinations)	Total (N=1036)	African Americans (N=342)	U.S. Caucasians (N=359)	U.S. Hispanics (N=238)	U.S. Asians (N=97)
CODIS 13	5.02E-16	1.13E-15	3.01E-15	1.39E-15	1.71E-14
Identifiler	6.17E-19	1.03E-18	6.94E-18	2.76E-18	5.30E-17
PowerPlex 16	2.82E-19	6.08E-19	4.22E-18	1.29E-18	2.54E-17
PowerPlex 18D	3.47E-22	5.55E-22	9.76E-21	2.58E-21	7.87E-20
ESS 12	2.99E-16	9.14E-16	9.68E-16	2.64E-15	3.41E-14
ESI 16 / ESX 16 / NGM	2.74E-20	6.02E-20	2.21E-19	4.03E-19	9.80E-18
ESI 17 / ESX 17 / NGM SElect	1.81E-22	6.44E-22	1.74E-21	3.99E-21	1.86E-19
PowerPlex 21	6.71E-27	2.08E-26	2.51E-25	7.96E-26	5.80E-24
CODIS 19 (-DYS391)	1.95E-23	1.08E-22	1.66E-22	1.93E-22	6.01E-21
GlobalFiler (-DYS391)	1.60E-27	5.63E-27	2.95E-26	5.10E-26	2.57E-24
PowerPlex FUSION (-DYS391)	1.36E-28	2.83E-28	5.25E-27	4.81E-27	2.01E-25

Slide kindly provided by John Butler (NIST) using data collected by Becky Hill & analysis tools developed by Dave Duewer

Developmental Validation – Preliminary data Mixture Study



Percent Minor Alleles Called

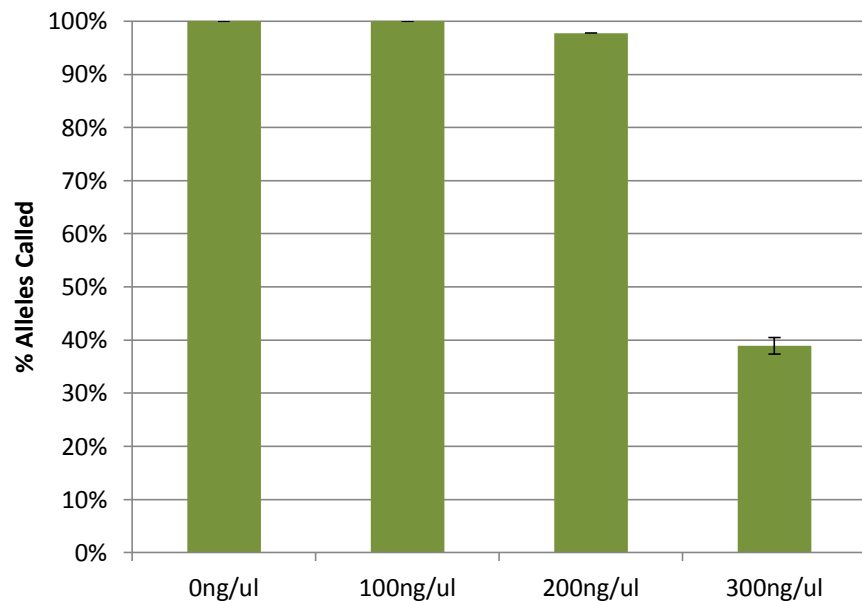


Developmental Validation – Preliminary data

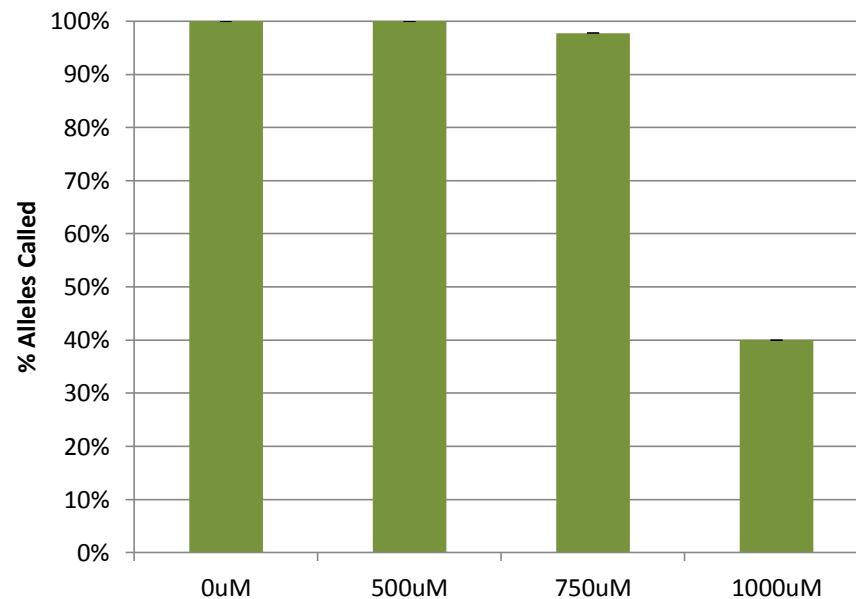
Inhibitors Study



Humic Acid



Hematin

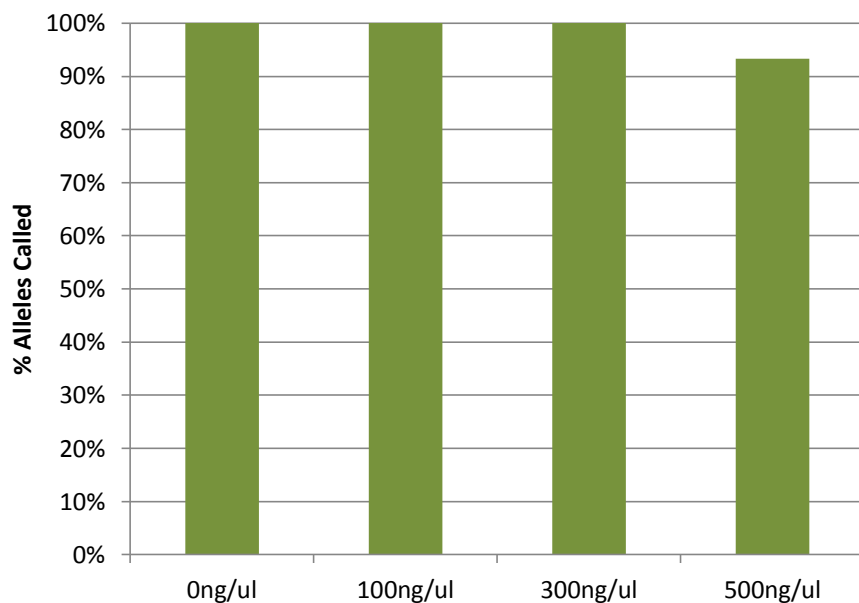


N = 2

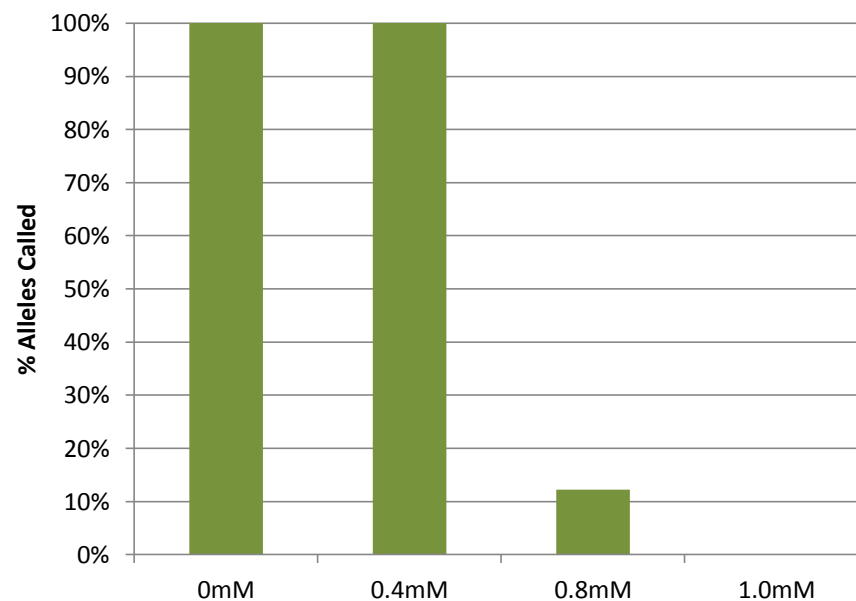
Developmental Validation – Preliminary data Inhibitors Study



Tannic Acid



EDTA



N = 2

PowerPlex® Fusion System



- Provides a rapid thermal cycling protocol
- Sensitive and robust for casework
- Allows for direct amplification for swabs and punches in databasing
- Is fully compatible with current 31xx & 3500 series Genetic Analyzers
- Includes all core CODIS and ESS loci as well as a Pentas and DYS391
 - Most discrimination available
- Will allow searching of *all* autosomal STRs historically entered in NDIS

Questions?

