

# Maintaining Scientific Integrity Through Cell Line Authentication

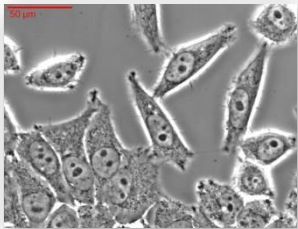
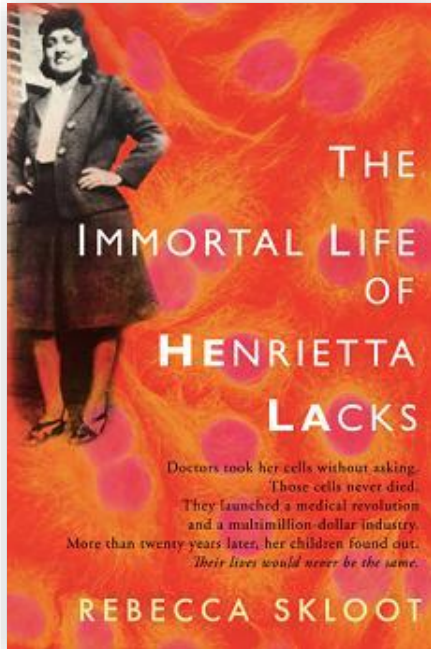
Gabriela Saldanha  
Genomics Business Unit  
Life Sciences Research

# Presentation Outline

---

- Background on cell line authentication (CLA)
- Basics of STRs and CLA protocols

# The legend of HeLa



HeLa cells



- HeLa- first immortal cancer line to be established in 1951 by George Gey
  - Cervical cancer derived from patient Henrietta Lacks
  - Widely used in research
- According to Rebecca Skloot, “More than 60,000 scientific articles had been published about research done on HeLa, and that number was increasing steadily at a rate of more than 300 papers each month.”*
- Rapid growing; can contaminate and overtake other cell types

# Stanley Gartler describes HeLa contamination of cell lines

1968: Gartler publishes first evidence of HeLa contamination based on G6PD isoenzymes

NAME	DESCRIPTION	ORIGIN	G6PD VARIANT
HeLa	Cervical adenocarcinoma, human	African	Type A (fast)
KB	Oral epidermoid carcinoma, human	Caucasian	Type A (fast)
HEp-2	Larynx epidermoid carcinoma, human	Caucasian	Type A (fast)
Chang liver	Liver, human	Caucasian	Type A (fast)
Int-407	Embryonic intestine, human	Caucasian	Type A (fast)



Nature Reviews, 2010

**Conclusion: 90% (18/20)  
human cell lines are 'HeLa'**

# Poor culture condition leads to contamination

- Poor tissue culture environment
- No disposable, plastic culture dishes
- No commercial media
- Cells grown on bench-top
- Bunsen burners and steam used for sterilization
- Technicians wore surgical masks, coats, gloves, booties, hair covers



# How does false cell line identity occur?

- Cross-contamination during original cell line development
- Cross-contamination in the lab from which the cell line was acquired
- Cross-contamination in your lab
- Mislabeling

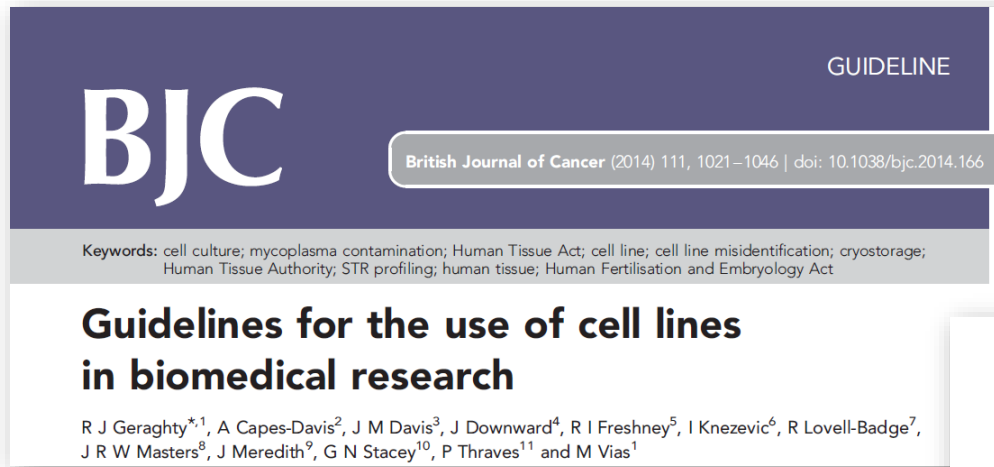
1/3

OF RESEARCHES

obtain their cell lines  
from another lab



# Good cell culture practice



# Misidentification of cell lines persists...

Year	Title of Article	Reference
2004	<u>LCC15-MB cells are MDA-MB-435: a review of misidentified breast and prostate cell lines</u>	Clinical & Experimental Metastasis 21 (6): 535, 2004
2007	<u>MDA-MB-435: the questionable use of a melanoma cell line as a model for human breast cancer is ongoing</u>	Cancer Biology & Therapy 6(9): 2007
2007	<u>A case study in misidentification of cancer cell lines: MCF-7/AdrR cells (re-designated NCI/ADR-RES) are derived from OVCAR-8 human ovarian carcinoma cells</u>	Journal: Cancer Letters, 245 (1-2): 350, 2007
2008	<u>Persistent use of “false” cell lines</u>	Journal: International Journal of Cancer 122 (1): 1, 2008
2010	<u>Verification and unmasking of widely used human esophageal adenocarcinoma cell lines</u>	NCI J Natl Cancer Inst 102(4): 271, 2010
2011	<u>MCF-7/ADR cells (re-designated NCI/ADR-RES) are not derived from MCF-7 breast cancer cells: a loss for breast cancer multidrug-resistant research</u>	Journal: Medical Oncology 28 (1): 135, 2011
2013	<u>Misidentification of putative medullary thyroid cancer cell lines RO-H85-1 and RO-D81-1</u>	J Clin Endocrinol Metab, 98(3):954, 2013
2013	<u>Beware imposters: MA-1, a novel MALT lymphoma cell line, is misidentified and corresponds to Pfeiffer, a diffuse large B-cell lymphoma cell line</u>	Genes, Chromosomes and Cancer (10): 986,2013



# Impact of misidentified cell lines on applied research

## Experimental results based on contaminated cell lines ...

- Clinical trial recruiting EAC patients
- 100 scientific publications
- At least 3 NIH cancer research grants
- 11 US patents

**CONTAMINATED**  
CELL LINES ARE EXPENSIVE

\$	\$\$\$
TO PROTECT	TO REPEAT ANALYSIS

# Consequences of using misidentified cell lines

- Loss of cell lines
- Loss of time and money
- Misinformation in the public domain
- Discordant or irreproducible results
- Tarnished reputation



© DAN PAGE COLLECTION/THEISPOIT

*“If we’re not using what we think we are using, we’re not testing our hypotheses. We’re just gumming up the literature. I’m not sure what we’re doing, but that’s not science.”*

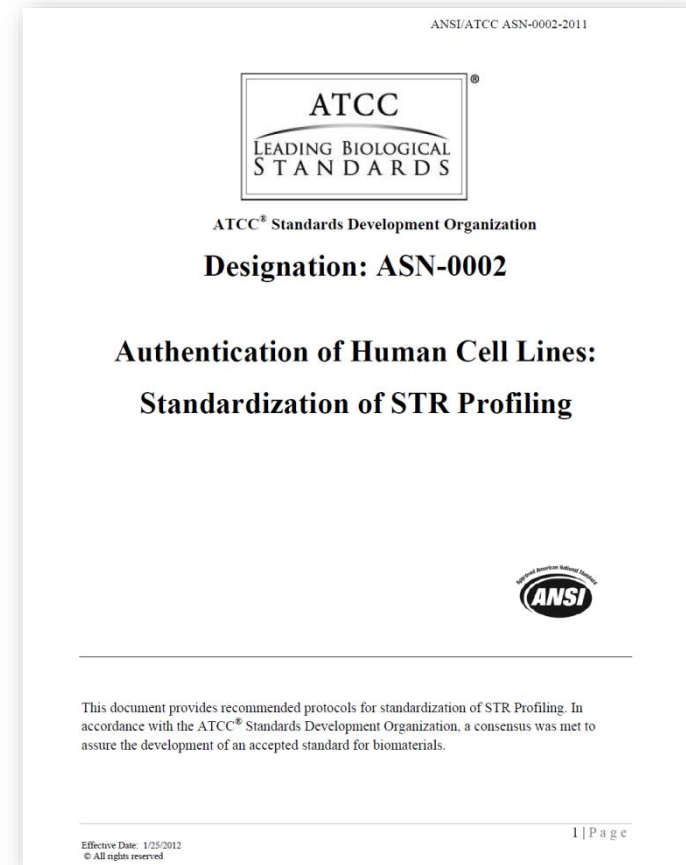
Jeffrey Boatright, Emory University, [The Big Clean-Up. The Scientist Magazine. September 2015](#)

# New ANSI/ATCC Standards Development Organization Guidelines

ASN-0002 consensus standard describing the use of **short tandem repeat (STR) analysis** for cell line authentication

With STR profiling authentication may include:

- (i) verifying the cells are of human origin
- (ii) evaluating the consistency of profiles between isolates/passages
- (iii) comparing STR profile to a database
- (iv) detecting contaminating human DNA



# Journals taking action

**nature** International weekly journal of science

Home | News & Comment | Research | Careers & Jobs | Current Issue | Archive | Audio & Video | RSS

Archive > Volume 520 > Issue 7547 > Editorial > Article

NATURE | EDITORIAL

## Announcement: Time to tackle cells' mistaken identity

15 April 2015

PDF Rights & Permissions

The differences between a cow and a monkey are clear. It is easy to tell a moth from a mosquito. So why are there still scientific studies that mix them up? The answer is simple: hundreds of cell lines stored and used by modern laboratories have been wrongly identified. Some pig cells are labelled as coming from a chicken; cell lines advertised as human have been shown to contain material from hamsters, rats, mice and monkeys.

Which is worse: that such crude mix-ups exist, or that, every day, researchers use cell lines that somebody, somewhere has already found to be mislabelled, misidentified or contaminated? To solve the first problem is a huge challenge. To address the second is a more manageable task, and one that researchers, journals, universities and funders must take seriously.

*Nature* and the Nature research journals are strengthening their policies to improve the situation. From next month, we will ask authors to check that they are not working on cells known to have been misidentified or cross-contaminated, and will ask them to provide more details about the source and testing of their cell lines.

NATURE METHODS | METHAGORA

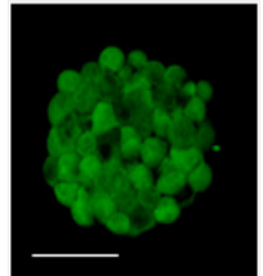
## A retraction resulting from cell line contamination

09 Sep 2013 | 2:09 AM | Posted by Daniel Evanko | Category: General Interest, Journal Policy, Nature Methods papers

After nine years in print, *Nature Methods* today published its first retraction; one that could have been prevented by cell line authentication. What does this mean for journal-mandated cell line testing?

In a *Nature Methods* paper published in 2010, Ivan Radovanovic and colleagues described a method to isolate cancer-initiating cells in human glioma without the need for molecular markers. Based on morphology and on a green autofluorescence, the authors reported they could use FACS to sort cancer-initiating cells from gliomasphere cultures (which had been derived from primary tumors). They also detected autofluorescence in cells from fresh glioma specimens, but at a much lower level.

Cells from the autofluorescent fraction could self renew clonogenically in vitro and were tumorigenic when transplanted into mouse brains, the authors reported, and in both cases performed better than non-autofluorescent cells from the rest of the culture or tissue. The origin of this autofluorescent signal was not understood at the time. The authors speculated it may have been related to the unique metabolism of the cancer-initiating cells.



Two-photon fluorescence image of live primary gliomasphere from retracted manuscript.

# Media outlets are paying attention

Up to 36% of all cell lines are contaminated...

[View this email in your browser](#)

## What is in your cell lines?

### The word is out!

National Public Radio is [reporting today](#) on the prevalence of contaminated and mis-identified cell lines being used in research. The Genomics core is hosting a [seminar](#) about this today. Find out how to verify your cell lines, join us TODAY at 1pm.

[Listen to the story](#)

npr news arts & life music topics programs listen

shots HEALTH NEWS FROM NPR

your health treatments & tests health inc. policy-ish public health

treatments

## Scientists Often Skip A Simple Test That Could Verify Their Work

DECEMBER 10, 2014 4:54 AM ET

*"There's a simple test that scientists could use to make sure the cells they're studying in the lab are what they think they are. But most of the time, academic scientists don't bother."*

WISCONSIN PUBLIC RADIO NPR NEWS & ...

Catch up with the NPR Hourly Newscast

shots HEALTH NEWS FROM NPR

your health treatments & tests health inc. policy-ish public health

TREATMENTS

## Mistaken Identities Plague Lab Work With Human Cells

Updated January 4, 2015 - 4:23 PM ET  
Published December 9, 2014 - 5:00 AM ET

# The culture of cell culture practices and authentication—Results from a 2015 survey

Concerns continue to grow regarding irreproducible basic biological and preclinical research.

Accurate documentation of cell line tissue of origin (i.e., identity), sex, and species are critical to ensure the credibility, reproducibility, and translation of data and results from cell culture based experiments.

The financial implications of misidentified or contaminated cell lines can be profound; as much as \$700 million dollars per year in research that could be at risk.





## ...Results continued

**18-36%**

OF CELL LINES

are misidentified

**52%**

“never” perform authentication or quality control

**74%**

Never conduct STR DNA profiling

**61%**

Cost

**53%**

Time

**35%**

Delay in research

**24%**

Reported “I don’t see the necessity; I am careful”

**22%**

laboratory manager or principal investigator is “unaware of or ignores the issue”

# Cell line authentication to improve reproducibility in cancer research



NIH revised guidelines for funding – [Enhancing Reproducibility through Rigor and Transparency \(effective Jan 25, 2016\)](#)

- Authentication of key biological and/or chemical resources

*“NIH plans to enhance reproducibility in multiple ways, by promoting greater scientific rigor and transparency in funding applications and publications, encouraging robust peer review, providing adequate training, removing perverse incentives, and emphasizing overlooked areas such as the consideration of both sexes in research **and the use of authenticated cell lines.**”*

# Cell line authentication to improve reproducibility in cancer research



Prostate Cancer Foundation  
*Cell Line Authentication Initiative*

Perform genetic and pathogenic tests  
for cell lines for future funding



Duke University  
*Free Human Cell Line Authentication Assays*  
**Free of charge**, a human cell line  
authentication service to Duke University  
School of Medicine investigators.

## CELL LINE AUTHENTICATION POLICY

### PURPOSE

The contamination or misidentification of cell cultures can seriously compromise research performed with such cell lines. The purpose of this policy is to ensure that information generated using cultured cells at The University of MD Anderson Cancer Center (MD Anderson) is obtained from samples that have been authenticated by DNA validation techniques, to rule out misidentification and inter- or intra-species contamination. **This will prevent the necessity for retraction of papers in which cell lines are found not to be of the reported lineage. It has been indicated by NIH that grant applications that fail to use acceptable experimental practices “would not fare well in the review process.”**

### POLICY STATEMENT

Validation by DNA analysis will be performed on cell lines used to produce scientific information at MD Anderson. This validation process **will enhance the quality of research at MD Anderson and maintain the scientific integrity and reputation of the institution.**

### SCOPE

This policy **applies to all workforce members utilizing cell lines for research.**

# Champions who are dedicated to advancing the #Authenticate goals



# Which journals ask for cell line authentication?\*

- AACR journals:
  - Cancer Discovery
  - Cancer Research
  - Clinical Cancer Research
  - Cancer Epidemiology, Biomarkers & Prevention
  - Molecular Cancer Research
  - Molecular Cancer Therapeutics
  - Cancer Prevention Research
- Carcinogenesis
- Cell Biochemistry and Biophysics
- Cell Biology International
- Endocrine Society journals:
  - Endocrinology
  - Endocrine Reviews
  - Journal of Clinical Endocrinology & Metabolism
  - Molecular Endocrinology
  - Hormones and Cancer
- International Journal of Cancer
- In Vitro Cellular & Developmental Biology - Animal



- Journal of Molecular Biology
- Journal of the National Cancer Institute
- Molecular Vision
- Nature Publishing Group:
  - Nature
  - Nature Reviews Molecular Cell Biology
  - Nature Genetics
  - Nature Reviews Immunology
  - Nature Reviews Cancer
  - Nature Reviews Neuroscience
  - Nature Biotechnology
  - Nature Methods
- Neuro-Oncology
- Placenta
- PLoS ONE
- Society for Endocrinology journals:
  - Journal of Endocrinology
  - Journal of Molecular Endocrinology
  - Endocrine-Related Cancer

\*Requirements vary with each journal



# How researchers can prevent the use of misidentified cell lines?

**ONLY 1%**

OF AUTHORS

authenticate their  
cell lines

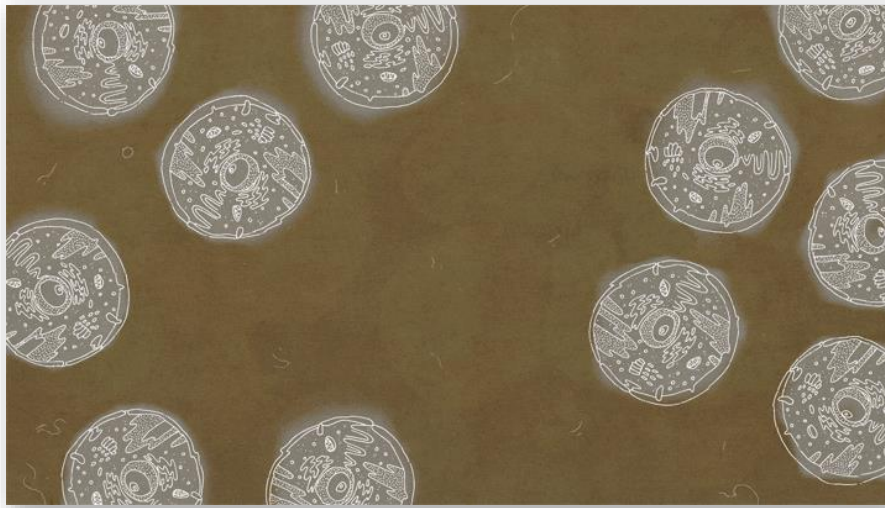
- Become aware of the magnitude of the problem
- Authenticate all cell lines you are working with!
- Eliminate laboratory practices that result in misidentification
- Notify recipients of misidentified cell lines
- Publish retractions/explanations of “false” publications

## RESULTS MATTER

- Confidence in the published data
- Meet the requirements of peer reviewed journals and grant submission
- Aid in troubleshooting undetermined or unexpected results

# Summary

- ✓ Cell authentication is more than just identity
- ✓ It consists of a number of orthogonal endpoints (identity, morphology, phenotype, ploidy, purity)
- ✓ Look for increasing demand for authentication of cells used in research and biological activity assays



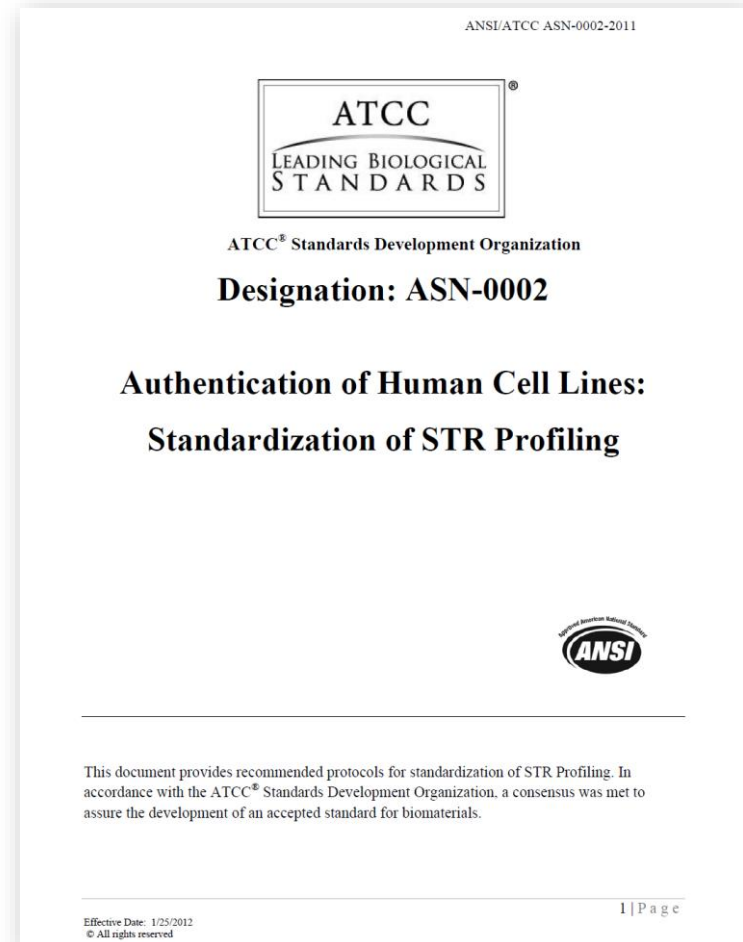
*Credit: Chris Nickels for NPR*

# STR Profiling: Authentication Standard for Human Cell Lines

# New ANSI/ATCC Standards Development Organization Guidelines

- ASN-0002 consensus standard describing the use of **short tandem repeat (STR) analysis** for cell line authentication
  - Also recommend assessing cell behavior, such as doubling time and morphology
- Plans to create a database containing STR genotypes of validated cell lines
  - Hosted by the National Center for Biotechnology Information (NCBI)
- Public database available on ATCC website

ANSI: American National Standards Institute



# ANSI authentication points

---

With STR profiling authentication may include:

- (i) verifying the cells are of human origin
- (ii) evaluating the consistency of profiles between isolates/passages
- (iii) comparing STR profile to a database
- (iv) detecting contaminating human DNA\*

\*Note that commercial STR kits are designed to detect human DNA, so other species' contaminating DNA will not be detected.

# STR analysis workflow

## Multiplex PCR through data analysis

DNA  
Sample

Multiplex PCR

- Amplification of select STR loci
- Simultaneous fluorescent labeling

Capillary Electrophoresis (CE)

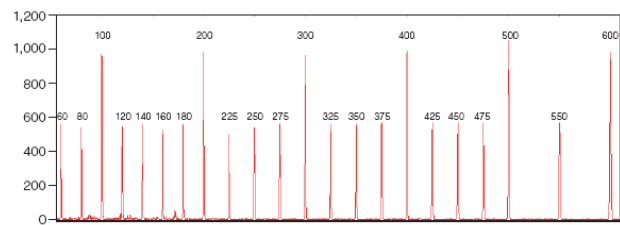
- Add Internal Lane Standard
- CE to size separate
- Fluorescent detection
- Also run Allelic Ladder in parallel

Data Analysis

- Calculate sizes based on ILS
- Compare fragment sizes to allelic ladder to determine STR alleles present in sample
- Compare to databases

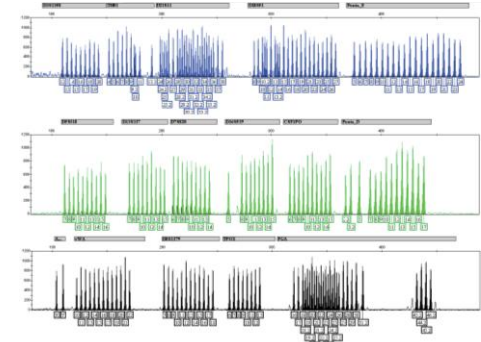


*Internal Lane Standard*



*Labeled size standards in a different color than STR fragments*

*Allelic Ladder*



*Labeled fragments of all possible alleles for each STR locus*



# STR (Short Tandem Repeat) genotyping

## STR Examples

Trinucleotide: ---CTTCTTCTTCTTCTT---

Tetranucleotide: ---AATGAATGAATG---

Pentanucleotide: -----AAAGAAAAGAAAAGAAAAGA-----

**STR = Short Tandem Repeats:** Multiple copies of a short (2-6bp), identical DNA sequence arranged in direct succession within a chromosome.

## STR Variation in an Individual

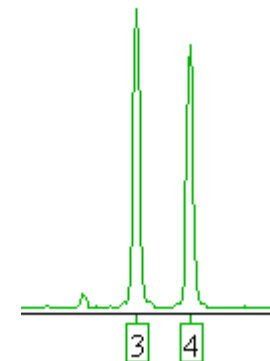
STR Locus



Homologous  
Chromosomes

**Allele 4** 1 2 3 4  
CTTCTTCTTCTT

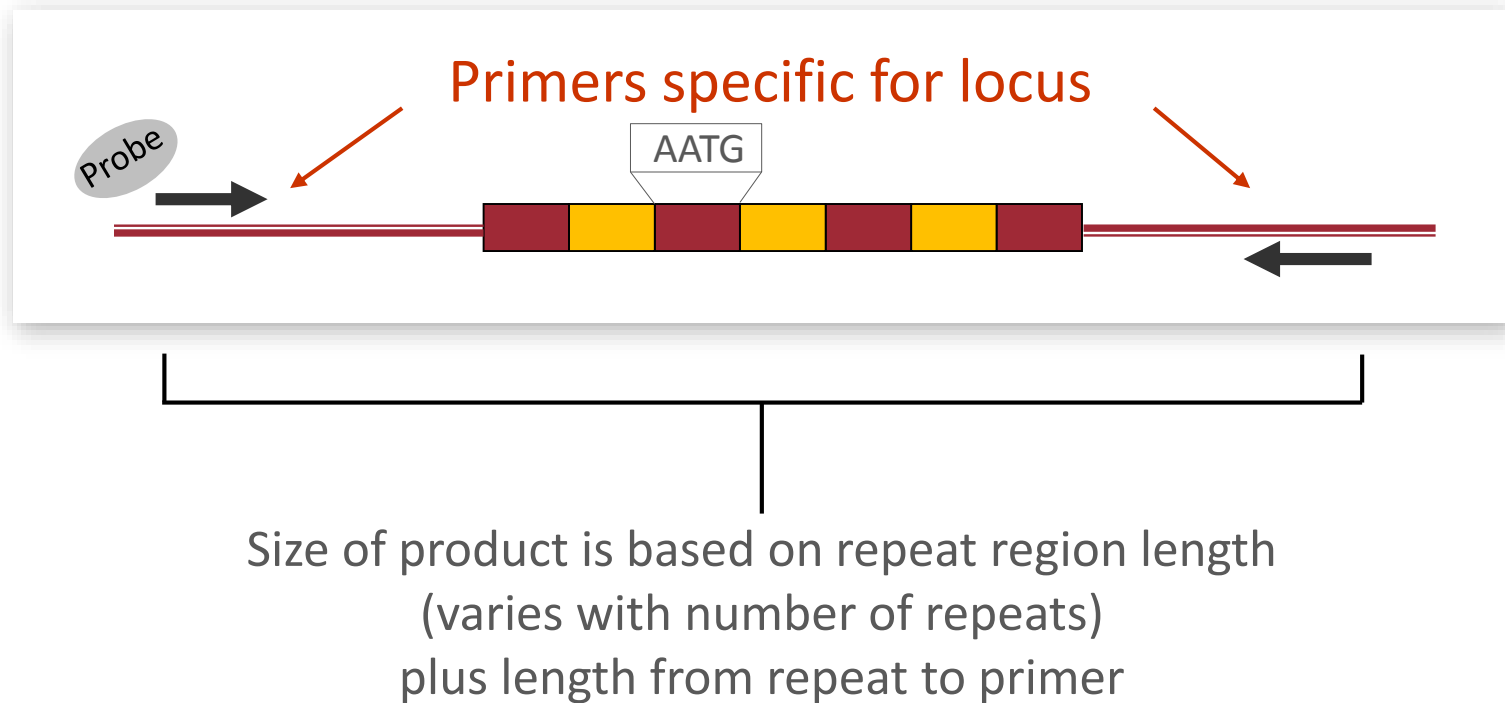
**Allele 3** 1 2 3  
CTTCTTCTT



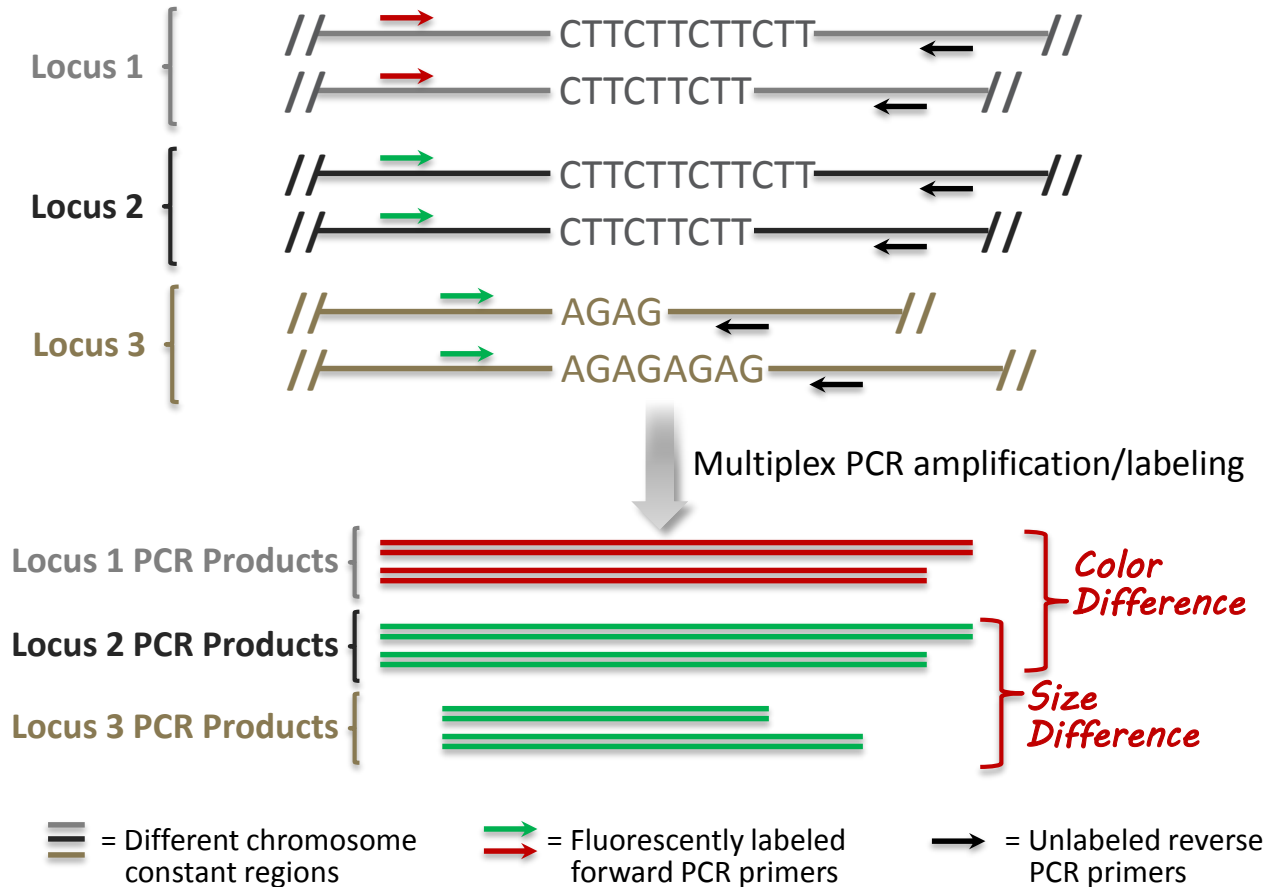
# Amplifying STRs

DNA is extracted from source material

Extracted DNA is the template for PCR



# Multiplex STR analysis combines size and color differences to analyze multiple loci



# Separating and detecting amplified STRs

## Capillary Electrophoresis

Separates amplification products based on size using Applied Biosystems' Genetic Analyzers



3130



3500

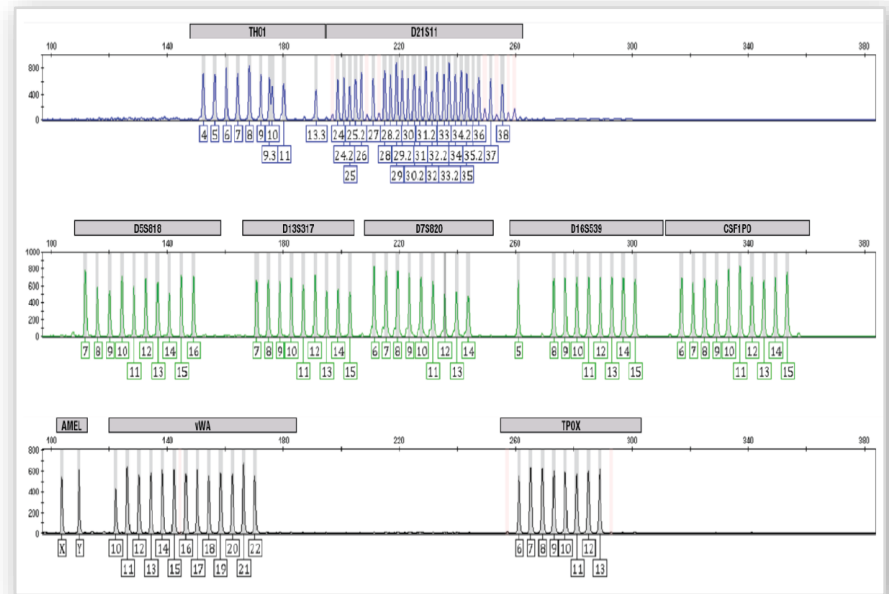
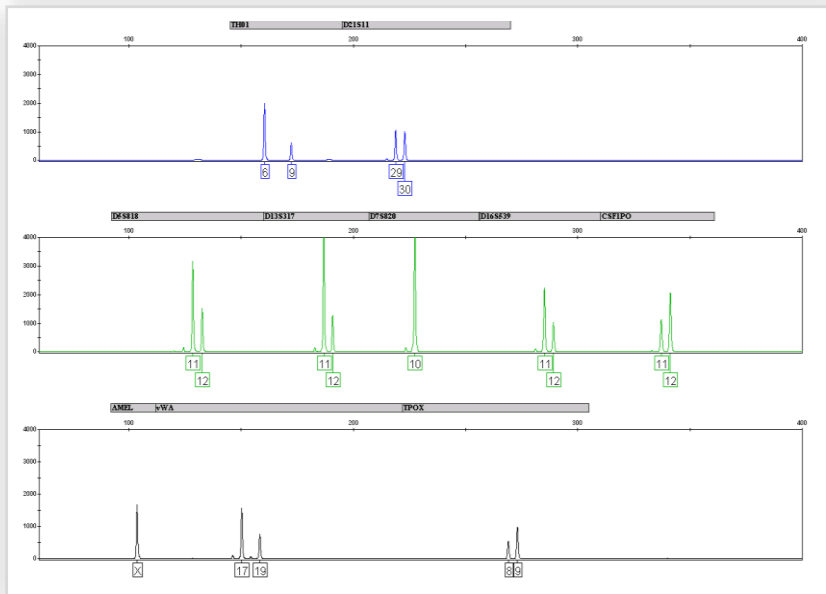


3730

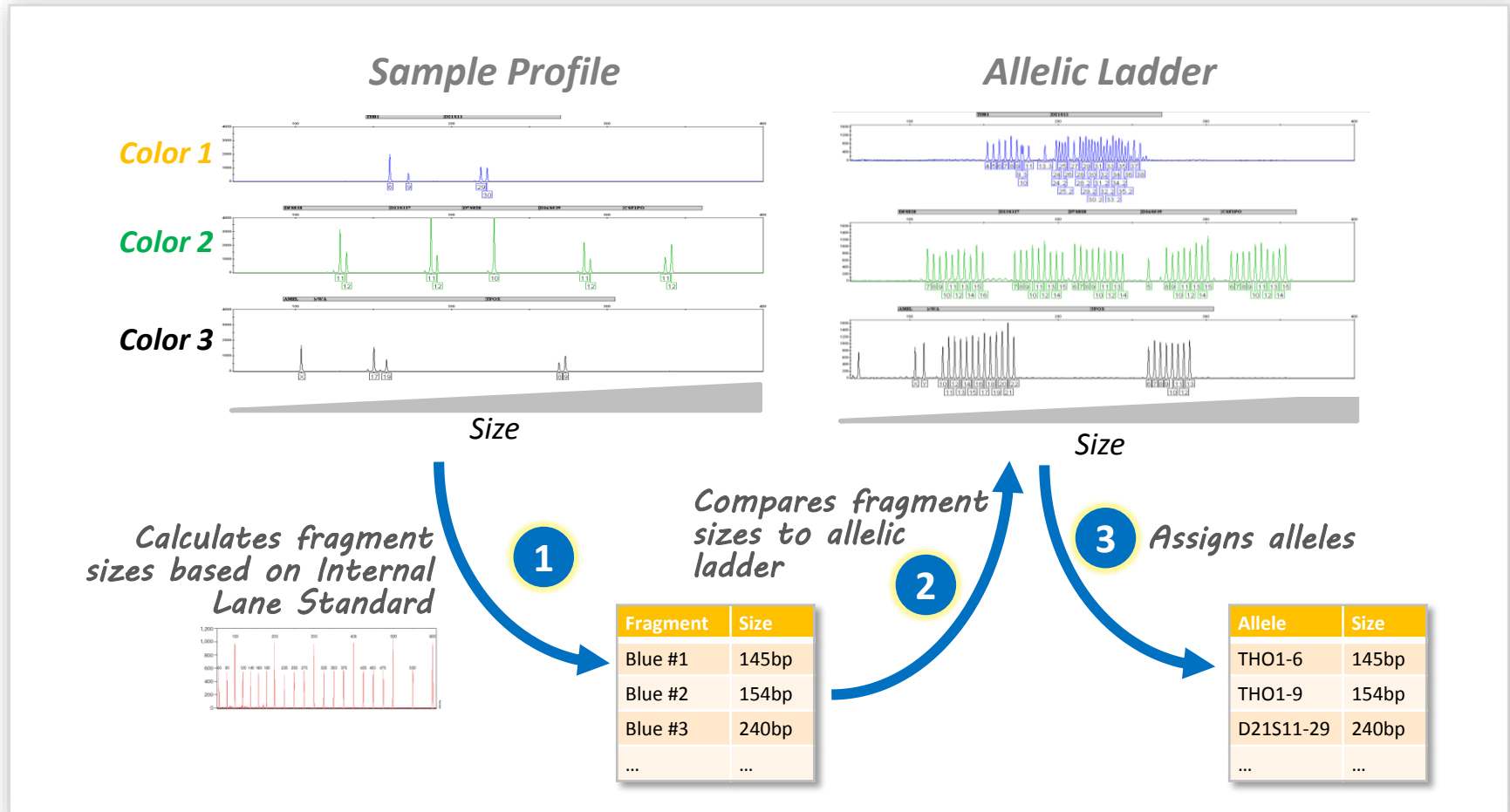
# Assigning allele types to peaks

Samples are compared to an allelic ladder to determine actual alleles

- Nomenclature is based on number of repeats
- Output is a simple text-based table



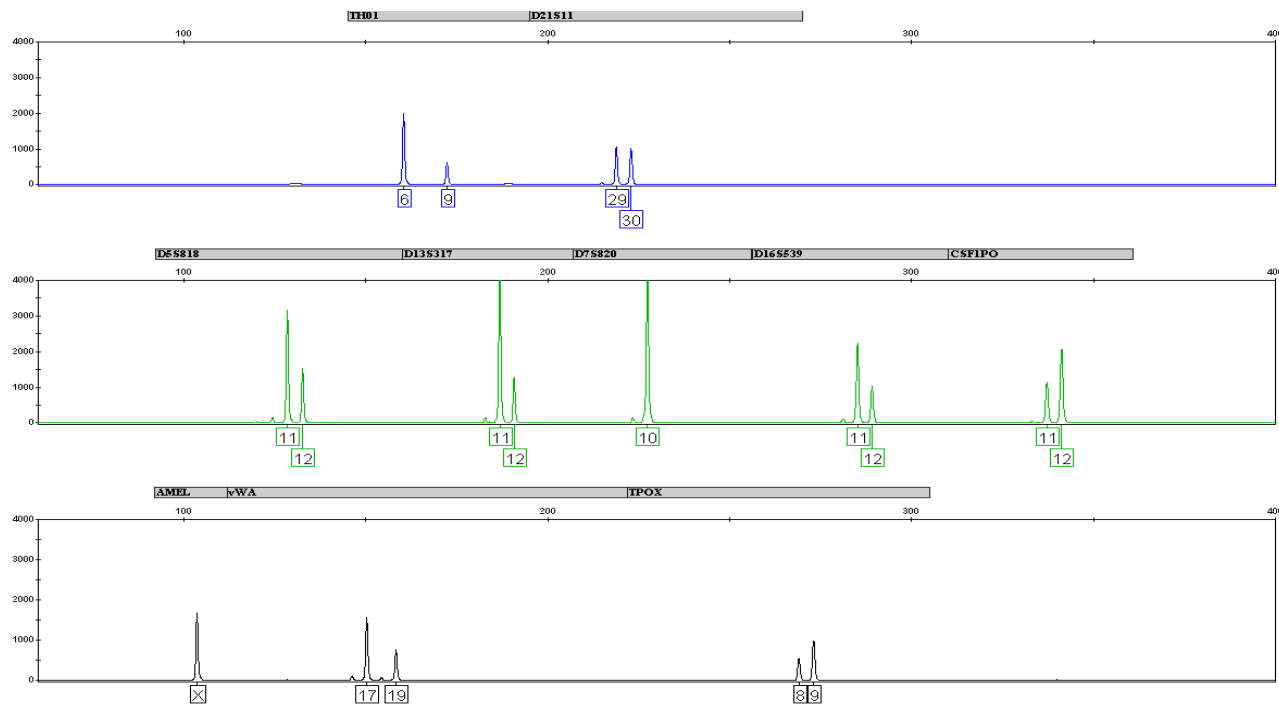
# Data analysis software identifies STR alleles





# Direct amplification profile from FTA® Paper

## HT-29 cell line with GenePrint® 10 System

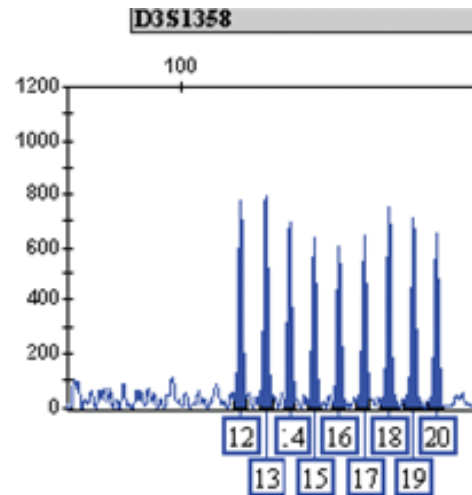


Simple text-based table

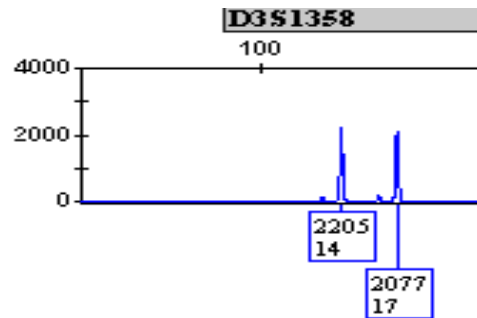
STR Locus	Alleles
TH01	6, 9
D21S11	29, 30
D5S818	11, 12
D13S317	11, 12
D7S820	10, 10
D16S539	11, 12
CSF1PO	11, 12
Amelogenin	X, X
vWA	17, 19
TPOX	8, 9

# Potential versus actual alleles at 1 locus

Alleles commonly encountered in the human population at this location

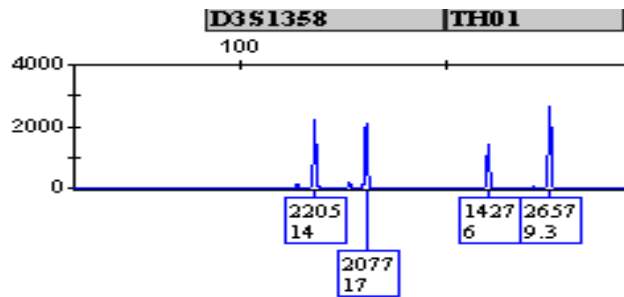
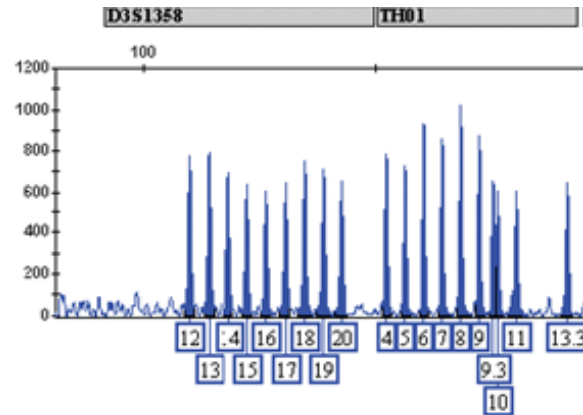


Profile generated from an individual



1 of 9 from mom  
+ 1 of 9 from dad  
1 of 18 possibilities

## ...using 2 loci



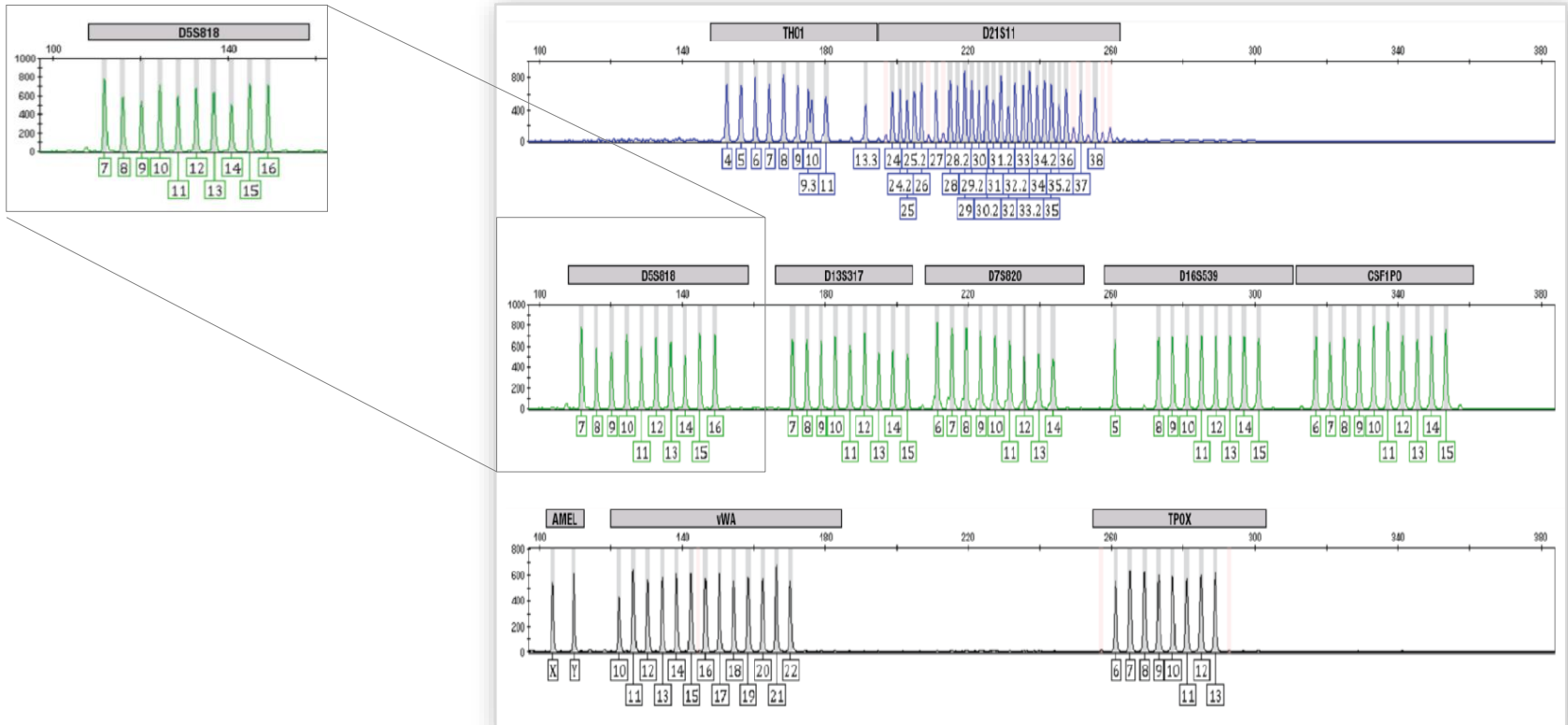
1 of 9 from mom  
+ 1 of 9 from dad  
1 of 18 possibilities  
at first location (D3S1358)

X

1 of 10 from mom  
+ 1 of 10 from dad  
1 of 20 possibilities  
at second location (TH01)

$$= 1:18 \times 1:20 = 1:360$$

...using 10 loci



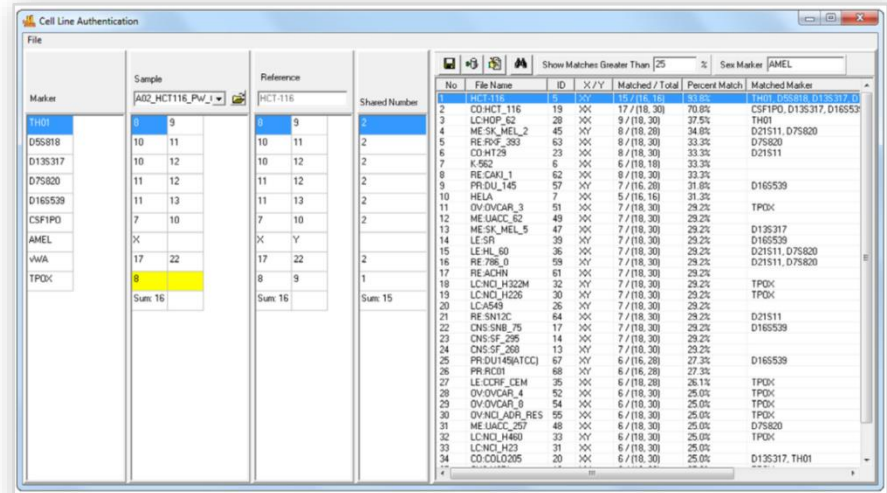
Matching probability: 1 in 18 x 1 in 20 x 1 in 50...

**1 in  $3 \times 10^9$  = 1 in 3,000,000,000 (*GenePrint*® 10)**

# Data analysis software options

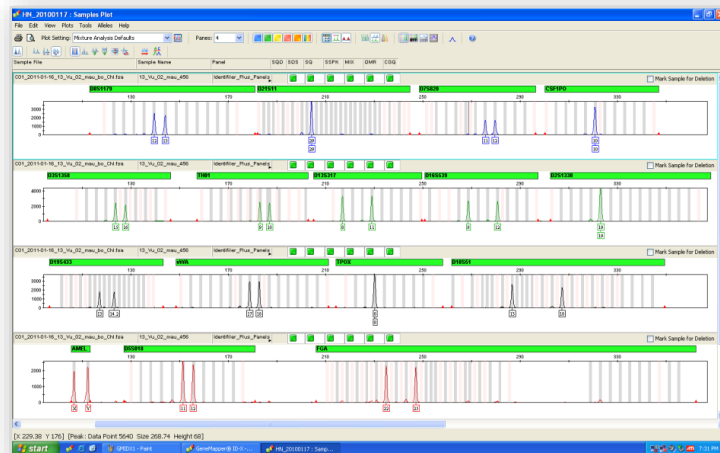
**GeneMarker® HID**  
Human Identity Software

By SoftGenetics



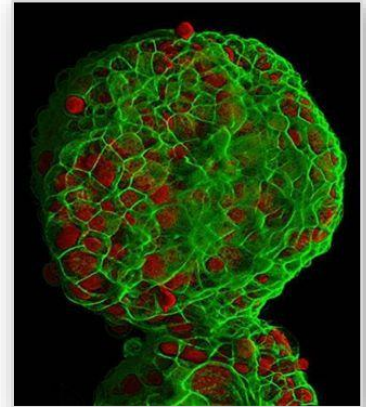
**GeneMapper® Software**  
**GeneMapper® ID-X Software**  
By Life Technologies

There are several versions. Some instrument files are not compatible with some versions. Additional files may be needed to perform full genotyping analysis depending upon the version used.



# Strengths and limitations

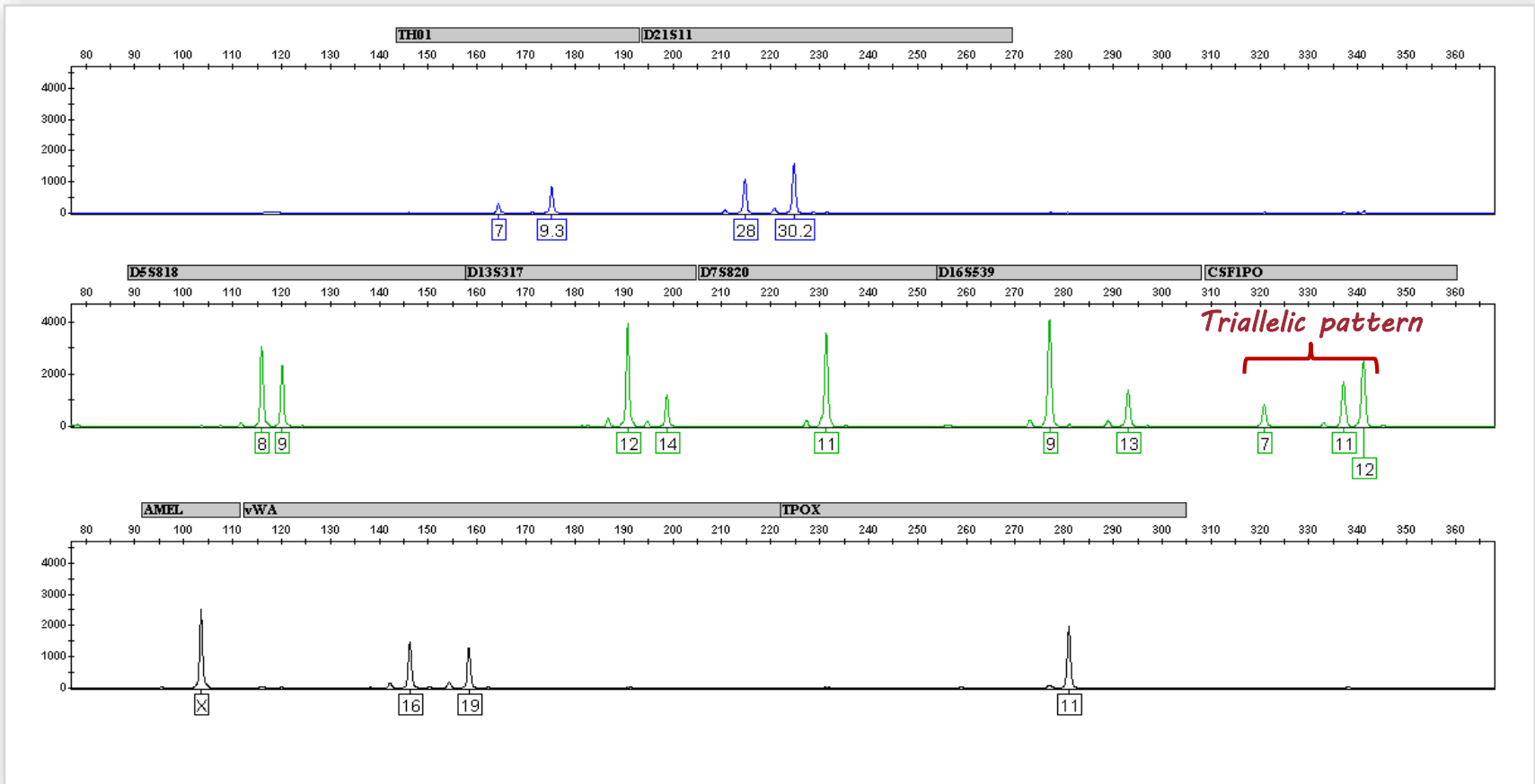
- STR systems can determine...
  - Relatedness of a cell line to a reference standard
  - Cell line cross-contamination
  - Gender (sex)
- STR systems cannot distinguish...
  - Different cell lines created from the same individual (donor)
  - Cell lines created from identical twins
- Promega STR Systems are designed to be human-specific
  - Cannot be used to genotype non-human species



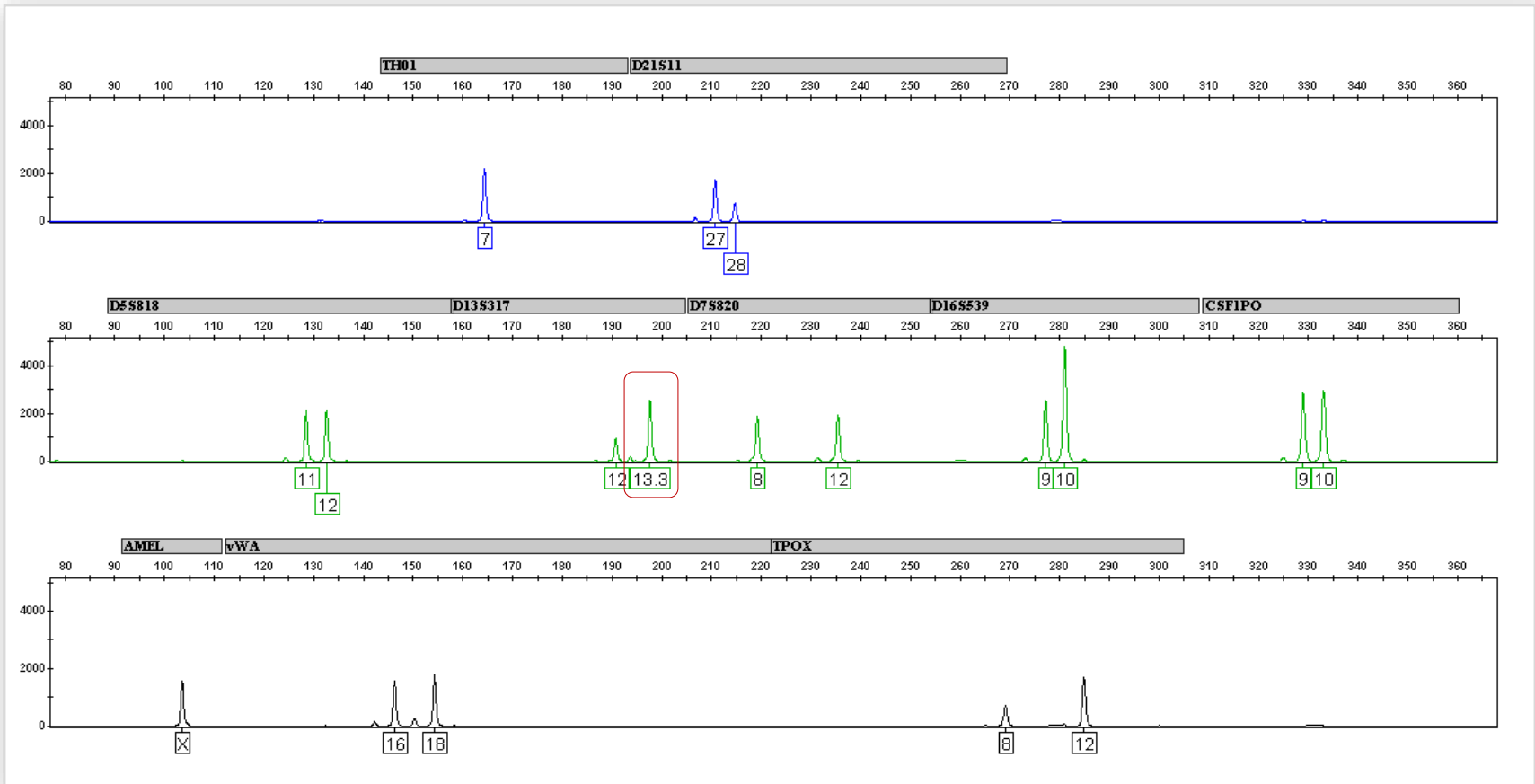
3D cell culture of breast cancer cells (MCF-7 cell line)



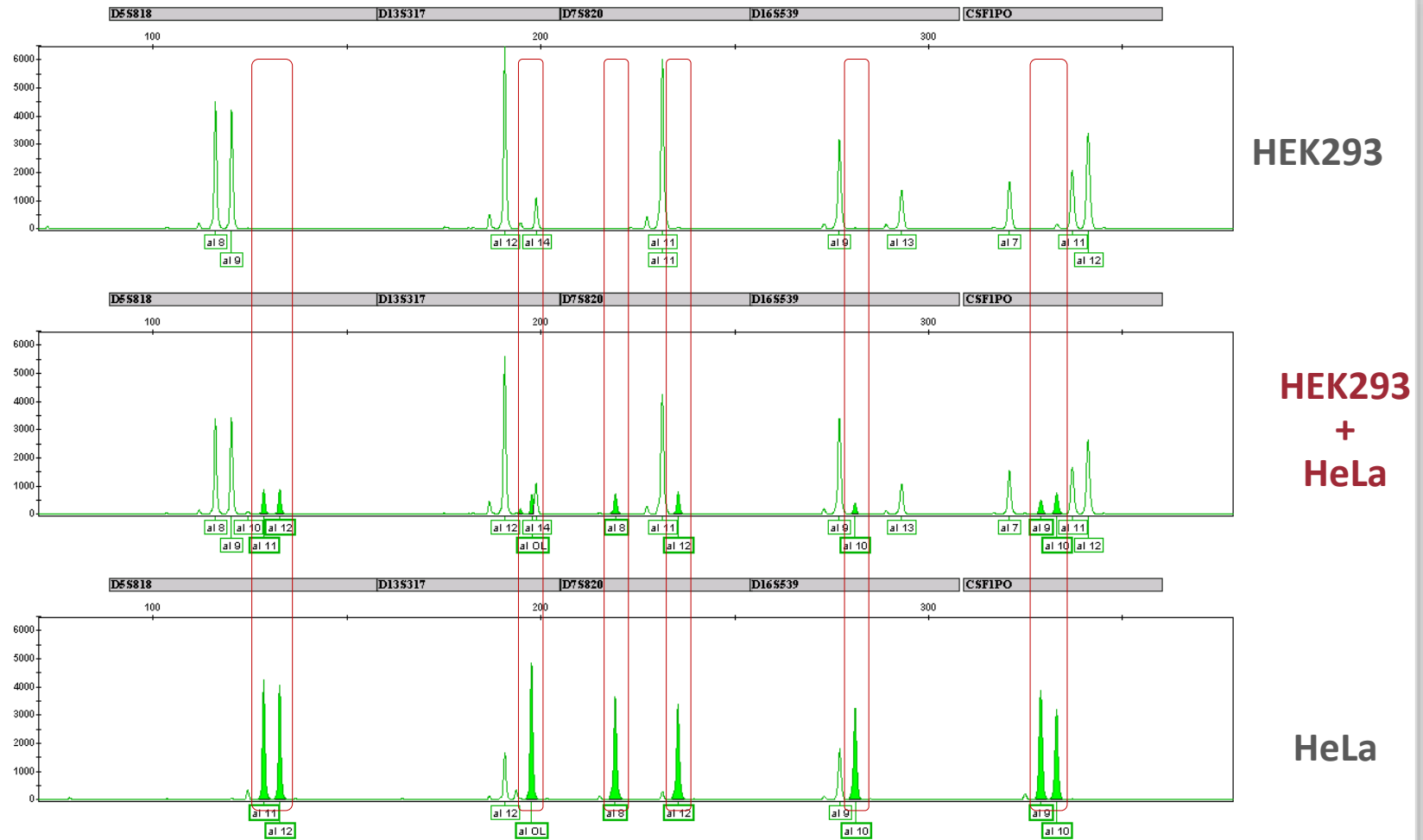
# HEK293 STR profile demonstrates triallelic pattern often found in immortalized cell lines



# STR profile illustrates HeLa characteristic D13 locus 13.3 allele



# Detection of contaminating HeLa cells in HEK293 culture using GenePrint® 10 System



# Sample comparison

Test Results for Submitted Sample				ATCC Profile or Reference Profile		
Loci	Query Profile			Database (or Reference) Profile: <b>A549</b> (CCL-185)		
D3S1358	16					
TH01	8	9.3		8	9.3	
D21S11	29					
D18S51	14	17				
Penta_E	7	11				
D5S818	11			11		
D13S317	11			11		
D7S820	8	11		8	11	
D16S539	11	12		11	12	
CSF1PO	10	12		10	12	
Penta_D	9					
Amelogenin	X	Y		X	Y	
vWA	14			14		
D8S1179	13	14				
TPOX	8	11		8	11	
FGA	23					
D19S433	13					
D2S1338	24					
Number shared alleles between reference and test profile:						<b>15</b>
Total number of alleles in the database/reference profile:						<b>15</b>
Percent match between the submitted sample and the database/reference profile: =(Number shared)/(Total number in reference)*100						<b>100</b>

# Matching algorithm

1. Combine total number of alleles observed in the Test Sample and Reference (TOTAL ALLELES)
2. Count the number of alleles shared by the Test Sample and the Reference Sample (SHARED ALLELES)
3. Use the Match Algorithm to calculate a percent match result for the two samples:

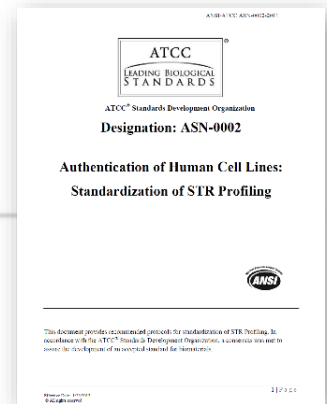
$$\text{Percent Match} = \frac{\text{SHARED ALLELES} \times 2}{\text{TOTAL ALLELES}}$$

Related samples generally yield a result in the 80-100% match range

- A percent match <80% should be investigated

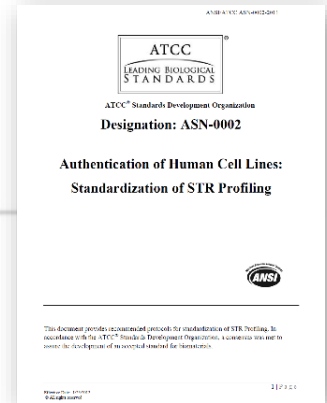
International Cell Line Authentication Committee

[http://standards.atcc.org/kwspub/home/the\\_international\\_cell\\_line\\_authentication\\_committee-iclac/](http://standards.atcc.org/kwspub/home/the_international_cell_line_authentication_committee-iclac/)



# Match criteria

**Authentic cell line**: a match at  $\geq 80\%$  of alleles across the eight (8) core STR loci.



**Match**: when two STR profiles show identical alleles. This is described as a percentage as shown above. Cell line samples matching at  $\geq 80\%$  of alleles across the eight (8) core loci are said to be related.

**Unrelated STR profiles**: when STR profiles match at  $< 55\%$  of alleles. STR profiles with alleles matching at 55-80% may be related and require further investigation.




# Cell line authentication is growing in importance

- ✓ **The number of misidentified/contaminated cells continues to grow**
  - Invalidates published and unpublished data leading to wasted time, effort and money
- ✓ **Granting agencies and journals are strongly recommending cell authentication**
  - ANS-0002 standards for cell authentication published by ANSI/ATCC
  - Check journal *Instructions to Authors* for requirements/recommendations
- ✓ **STR profiling is the gold standard for cell authentication**
  - Genotyping method that amplifies highly variable short tandem repeat loci to create a DNA fingerprint of a sample
  - Also important to observe phenotype of cells to detect possible changes
- ✓ **GenePrint® 10 System is optimized for human cell authentication**
  - Contains all loci recommended in ANS-0002 standards and available in public databases
  - Contains one additional loci for greater discriminatory power
  - Easy-to-use, complete kit compatible with multiple PCR & CE instruments

# International Cell Line Authentication Committee

**ICLAC**  
INTERNATIONAL CELL LINE AUTHENTICATION COMMITTEE

Home Resources Databases Case Studies References About ICLAC Members Partners You can Help



Cell lines are grown in laboratories worldwide to help us understand how normal cells function and how disease happens.

2014 International Cell Line Authentication Committee.

### Welcome To ICLAC

The International Cell Line Authentication Committee aims to make false or misidentified cell lines more visible, and promote awareness and authentication testing as effective ways to combat the problem.

### Resources

ICLAC resources are written to help scientists learn about misidentified cell lines and how to detect through authentication testing.

[Read more](#)

### Databases

ICLAC's list of cross-contaminated or otherwise misidentified cell lines can be found here.

[Read more](#)

### How can You Help?

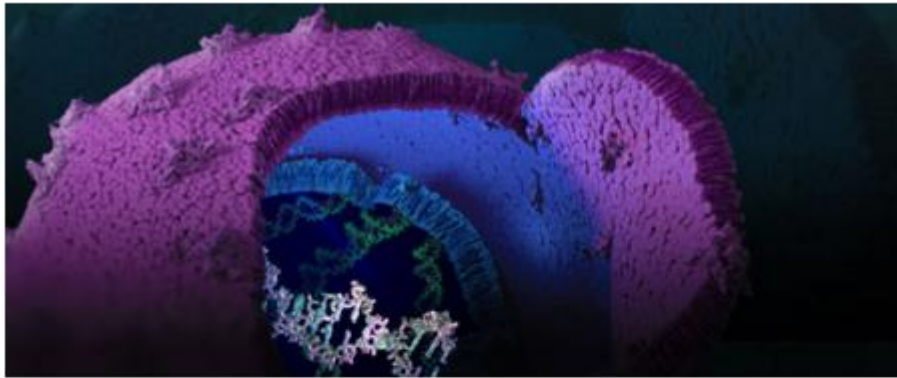
You can help to develop new resources and raise awareness of cell culture problems. Participate in the #authenticate campaign.

[Read more](#)

*ICLAC is an independent committee with support from the scientific community. Thank you to ICLAC Members and Partner Organizations for your support.*

<http://iclac.org/>

## Cell Line Authentication Testing



Cell Line Authentication Matters	How Authentication Works	Data Interpretation	Service Providers	Journals Requiring Authentication
<p>Human cultured cell lines are used in a number of biomedical research and clinical applications, including cancer research, drug discovery, genetics and biobanking. However, misidentified human and animal cell lines have continued to be used even today despite multiple and repeated warnings, articles and letters by prominent scientists in the field calling for authentication.</p> <p>The need is great for researchers to authenticate their human cell lines. The call for authentication has been made for over five decades now, yet still, many researches do not perform this vital quality assessment. Recently there have been more spotlights shining on this deficiency from National Public Radio (NPR) series last December, comments from Dr. Francis Collins of the NIH, Nature announcing changes to abstract submission and most recently, newly revised NIH grant application instructions. As a result, more institutions are coming around to offering, even requiring, human cell line authentication through their DNA core labs or service providers.</p> <p>Cell line authentication is achieved by genetic profiling using polymorphic short tandem repeat (STR) loci. STR loci consist of repetitive sequence elements 3–7 base pairs in length. These highly discriminative markers can be utilized using rapid and inexpensive multiplex-PCR based method for the identification and detection of contaminating human cells.</p>				



Share 0

Tweet 10

### EXTERNAL RESOURCES

[International Cell Line Authentication Committee \(ICLAC\)](#)

[Guide to Human Cell Line Authentication](#)

[Database of Misidentified Cell Lines](#)

[Match Criteria Worksheet](#)

[Cell Line Checklist for Manuscripts and Grant Applications](#)

[Authentication of Human Cell Lines by STR DNA Profiling Analysis](#)

[Maintaining High Standards in Cell Culture](#)

### DATABASES

[ATCC STR \(login required\)](#)

[DSMZ STR \(login required\)](#)

[NCBI BioSample \(open for sample submission\)](#)

### RELATED RESOURCES

[Short Tandem Repeat Analysis in the Research Laboratory](#)

[Doing Good Science: Authenticating Cell Line Identity](#)

### WEBINAR

[Human Cell Line Misidentification](#)

[www.promega.com/cla](http://www.promega.com/cla)

**Maintaining scientific  
integrity through cell  
line authentication**

**Thank you**

