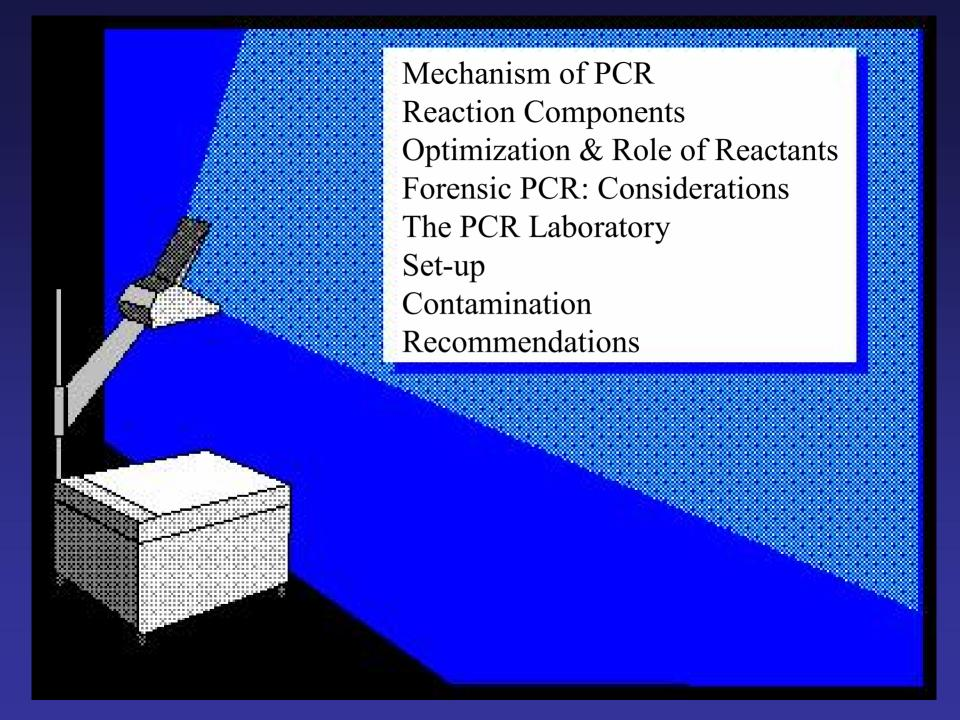
### Polymerase Chain Reaction

Workshop
Phoenix, AZ
October 4, 2004









Medicine
HIV detection
Disease diagnosis
Prenatal & carrier detection



#### Development of PCR

- 1969 Brock & Freeze: Thermus aquaticus
- 1976 Characteristics of Taq polymerase
- 1980 Temperature optimum for DNA synthesis
- 1985 Mullis: PCR using thermostable polymerase
- 1985 Cetus: automated thermocycling

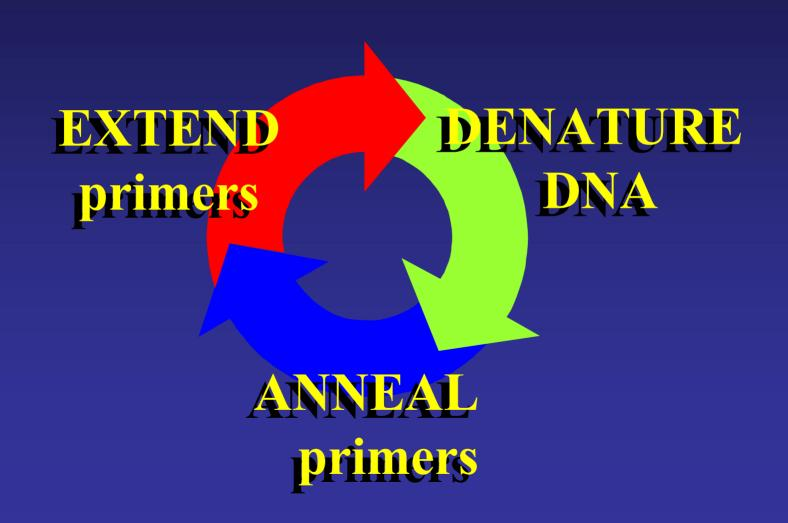
#### PCR Defined...

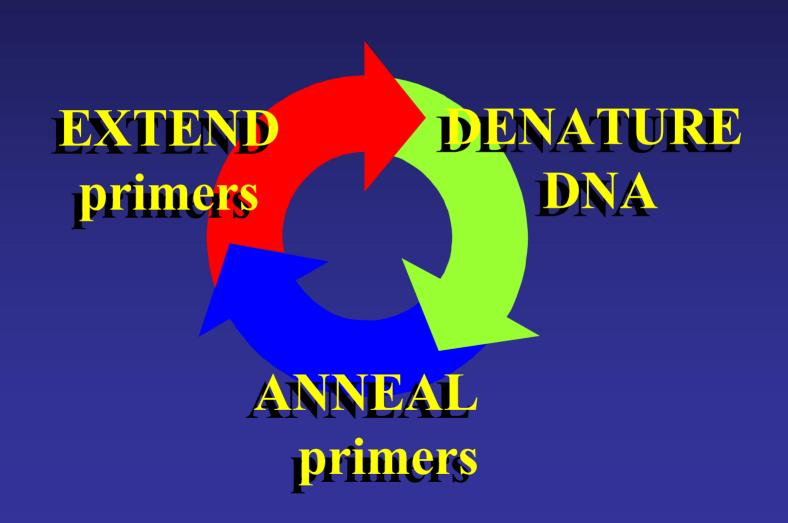
- An in vitro process
- Produces millions/billions of copies of target DNA fragments (subanalytical to analytical)
- Cyclical enzymatic reaction
- Replicates DNA products of previous cycles are used as templates in subsequent cycles

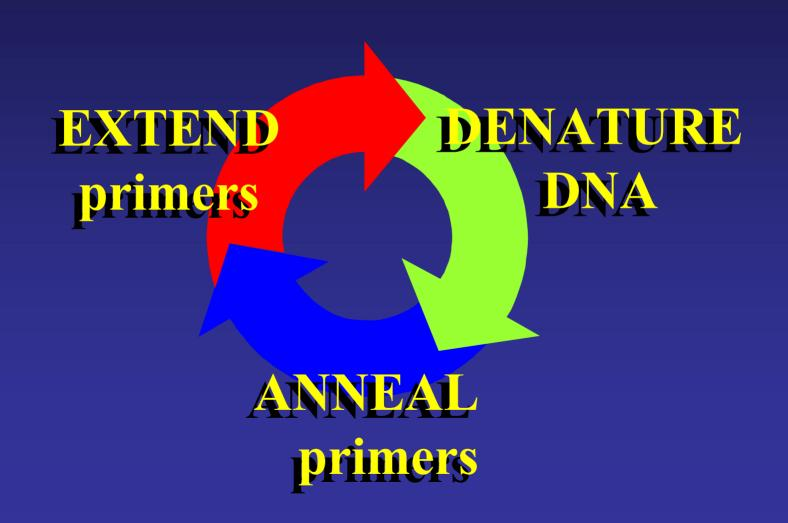
# 2<sup>30</sup> Copies PCR!

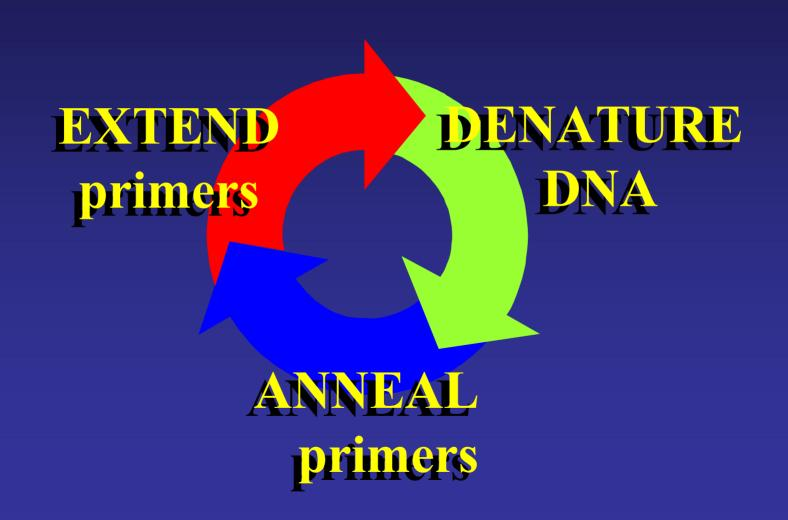
#### POWER OF PCR

- Subanalytical levels of DNA amplified to analytical levels
- Exponential increase in the amount of product:  $2^n$  where n = the number of cycles
- Theoretically:
  - -n = 20 yields a million-fold increase
  - -n = 30 yields a billion-fold increase

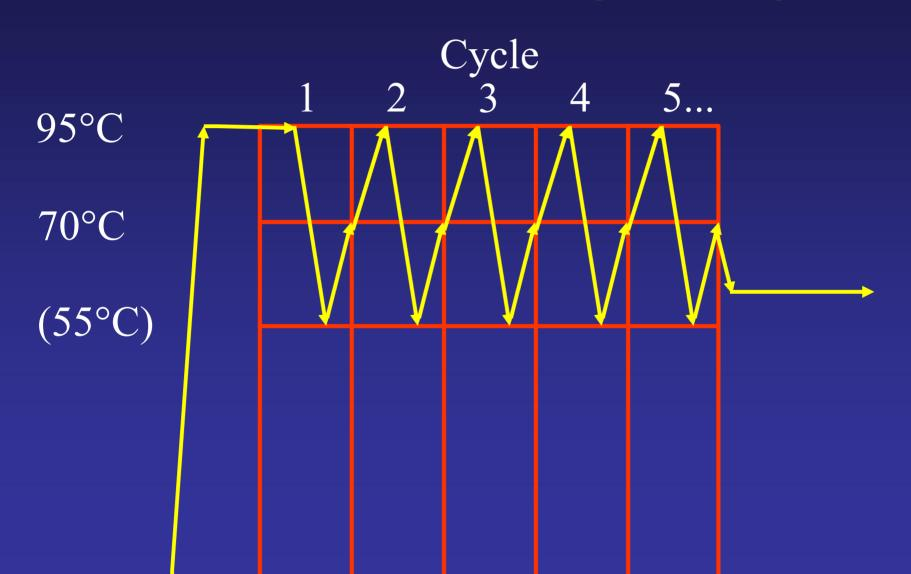








#### THREE-TEMPERATURE PCR



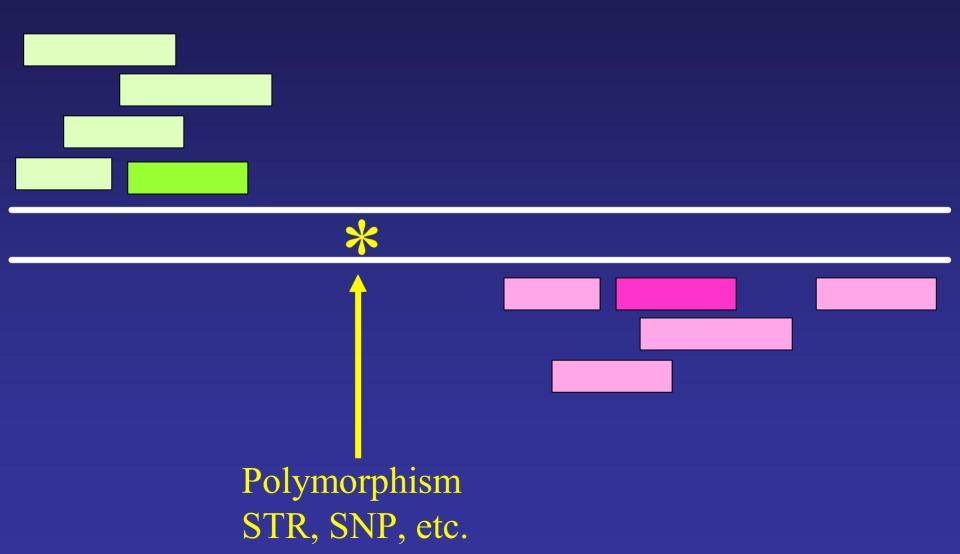
#### **Primers**

- Short oligonucleotides
- Linear
- Single Stranded
- Sequence complementary to target

5'-ATACC-3'

3'-GATGG-5'

#### Primer Selection



#### Primer Annealing

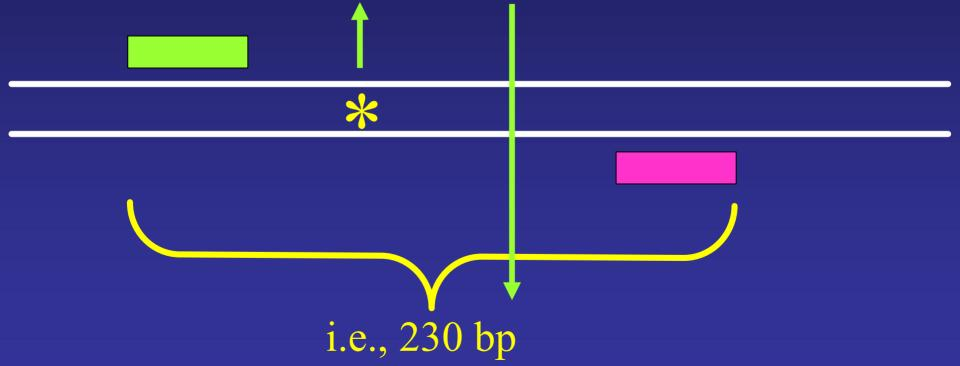
5' CTAAGT 3' 3' GATTCA 5'



5' **AATCTT** 3' 3' **TTAGAA** 5'

#### Defined Target Region

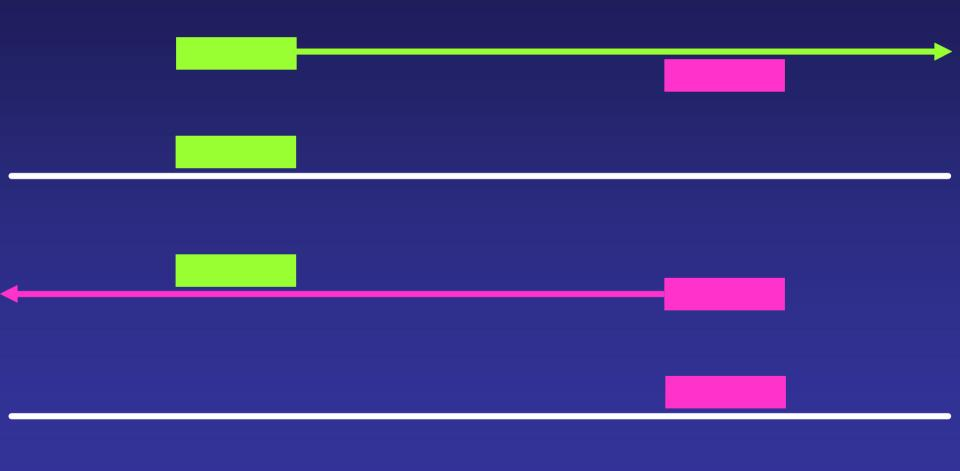
nucleotide difference, insertion, deletion, VNTR



# Cycle 1, Step 2 ANNEALING

### Cycle 1, Step 3 EXTENSION

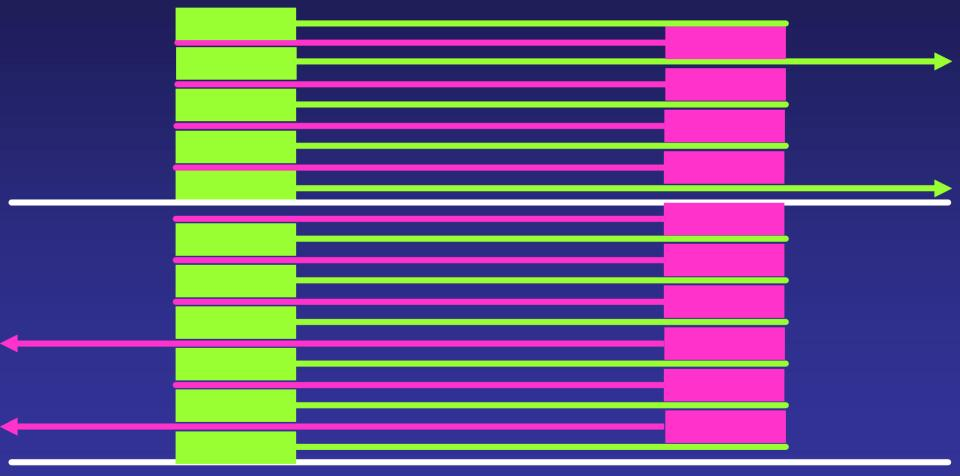
### Cycle 2, Step 2 ANNEALING



### Cycle 2, Step 3 EXTENSION

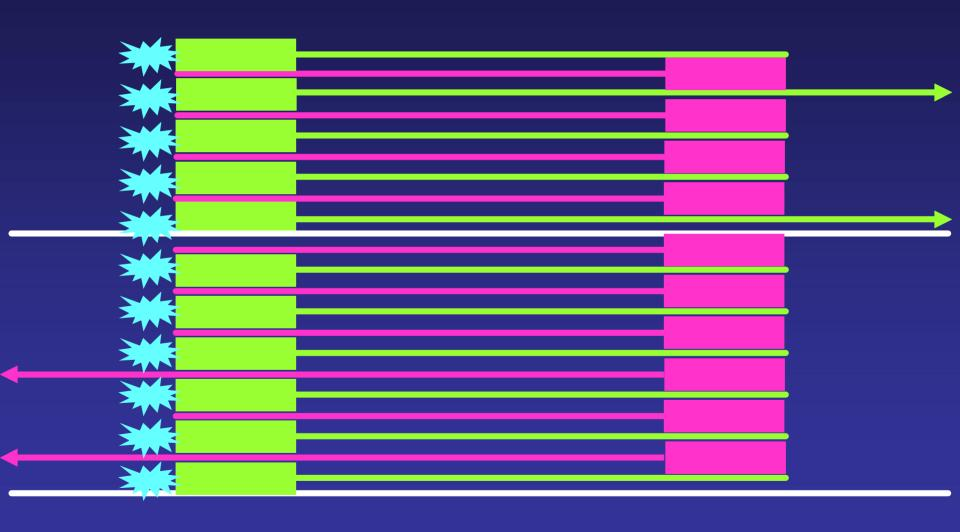


# Accumulation of Target Fragment



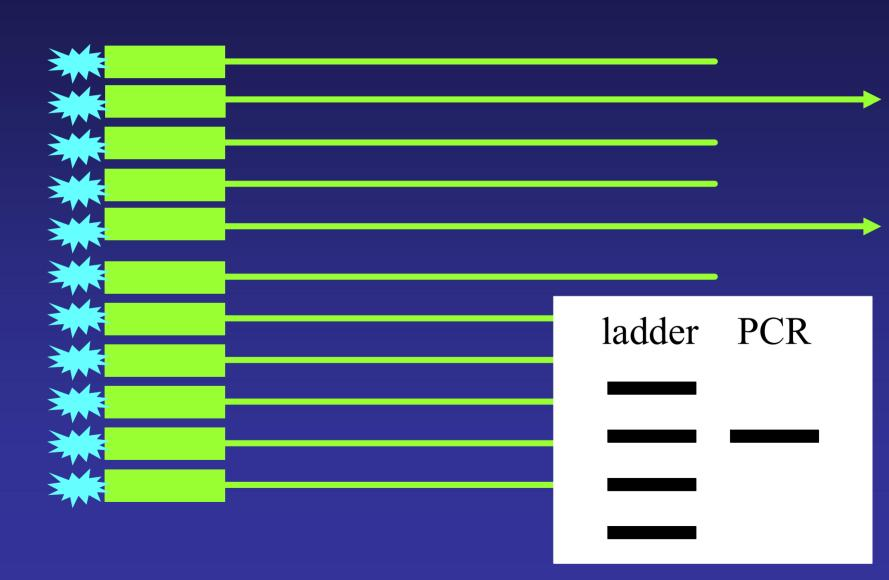


#### Fluorescent Detection





#### Fluorescent Detection



### EXAMPLE: 3-Temperature PCR

• Pre-Cycling Denaturation:

- Taq 95°C 1 min

AmpliTaq Gold 95°C 11 min

• Cycle 28 times:

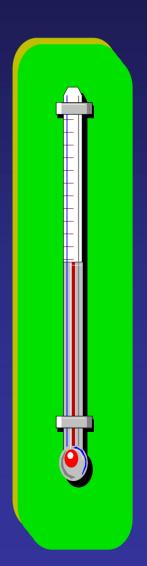
– Denaturation 95°C 1 min

- Annealing 60°C 1 min

- Extension 72°C 1 min

• Final Extension

- Extension 60°C 45 min



#### Final Extension



Complete truncated strands



Template-independent nucleotide addition

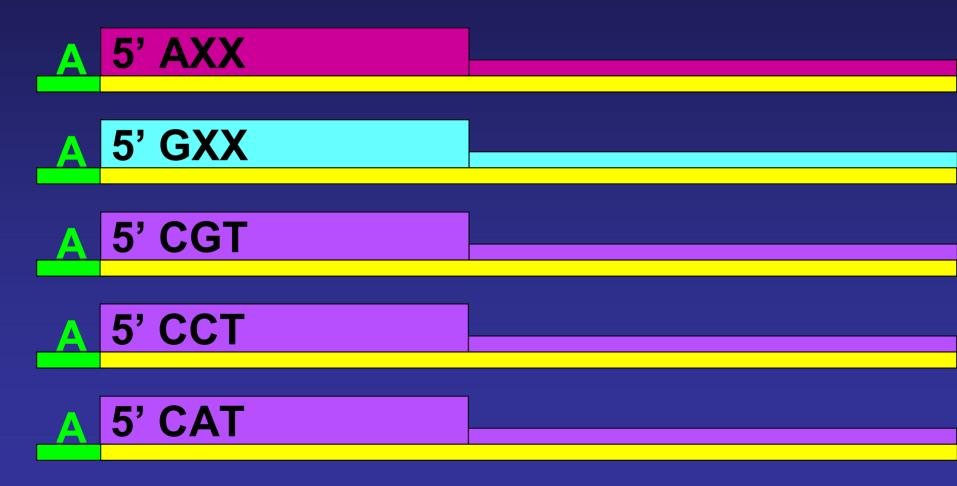
### Non-Templated Nucletide Addition ("Plus A")

- Many DNA polymerases add an extra nucleotide to the blunt end of a DNA product *without* the assistance of a template
- Usually A is added
- The resulting band/peak is 1-bp larger in size than that predicted from the DNA sequence and primer locations

### Non-Templated Nucletide Addition ("Plus A")

- Occurs during each PCR cycle and during final extension
- A is not added as well to some blunt ends
  - The sequence at the end of the DNA has an influence

# 5' end of primer affects adenylation



# 5' end of primer affects adenylation

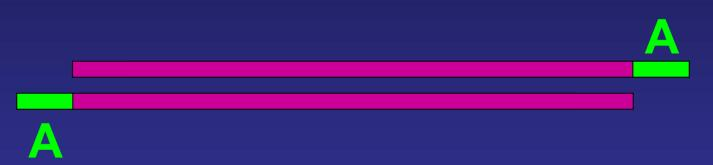
**5' TXX** 

5' CTT

Adenylation reduced

## Non-Templated Nucletide Addition ("Plus A")

Double-stranded DNA / PCR Product



## Non-Templated Nucletide Addition ("Plus A")

Temperature effects whether the ends remain together



## Non-Templated Nucletide Addition ("Plus A")

Final Extension time can been increased (e.g., 30 min to 60 min)

A



#### Taq DNA Polymerase

- Thermostable DNA Polymerase
- Isolated from *Thermus aquaticus* YT1
- 94 kDa protein
- $T_{1/2} = 40$  minutes at  $95^{\circ}C^{***}$
- 5'-3'exonuclease activity
- no 3'-5' exonuclease activity
- processivity = 50-60 nucleotides
- extension rate = 4-26kb/ min at 70-80°C

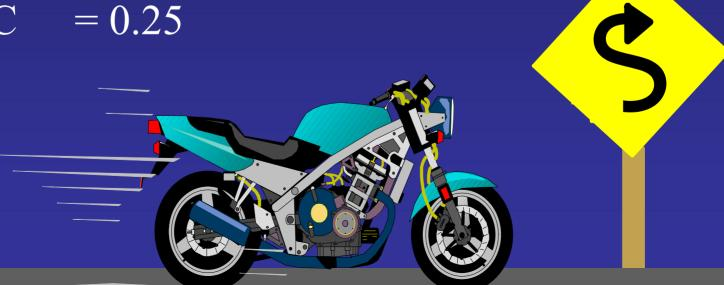
#### Taq Polymerase Activity

#### Extension Rate at:

- 70 °C = 60 nt/sec
- 55 °C = 24
- $36 \, ^{\circ}\text{C} = 1.5$
- $\overline{22 \, ^{\circ}\text{C}} = 0.25$

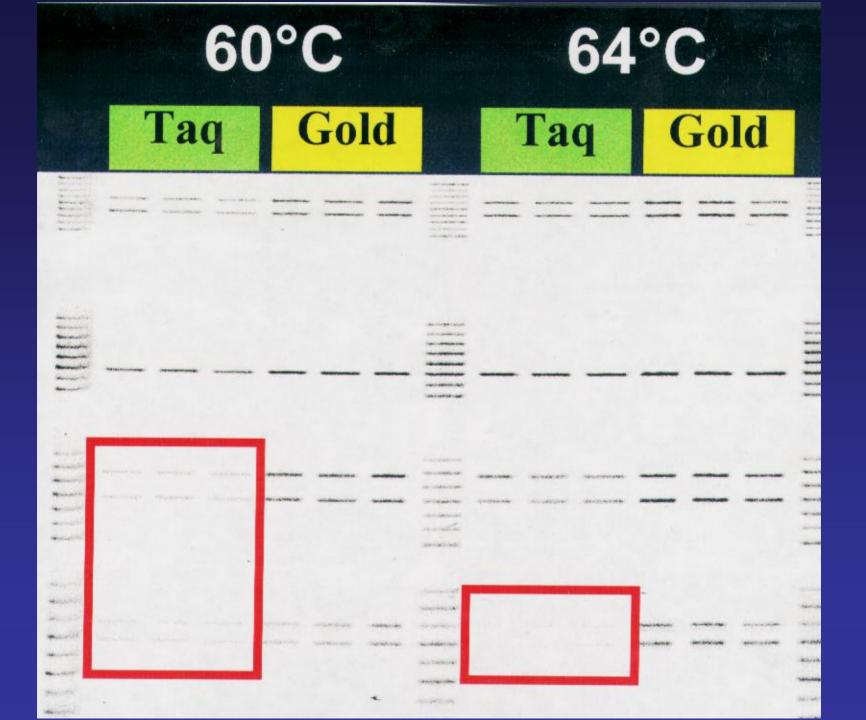
#### **Processivity:**

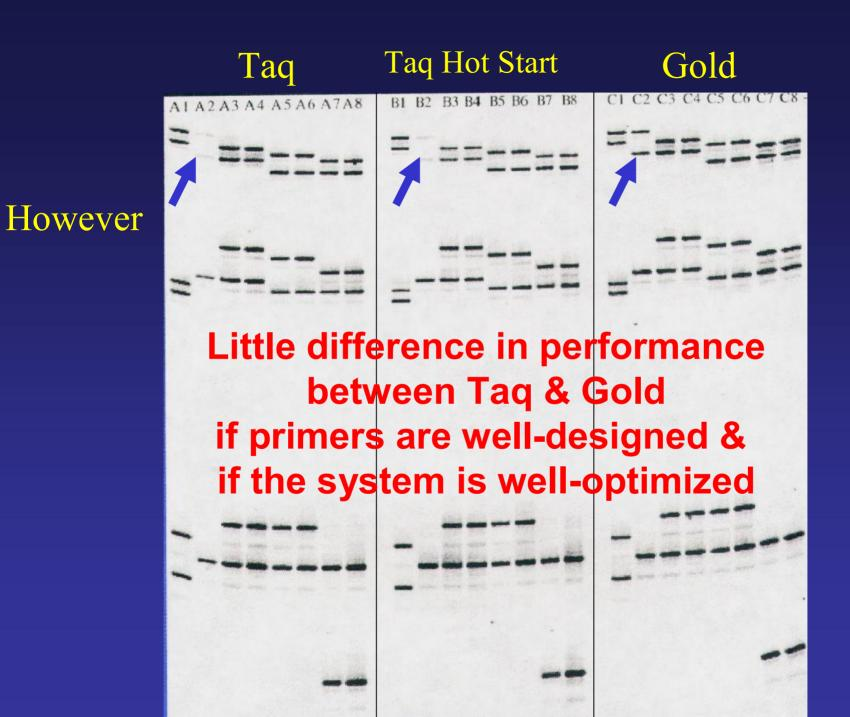
•  $70^{\circ}C = 50-60 \text{ nt}$ 

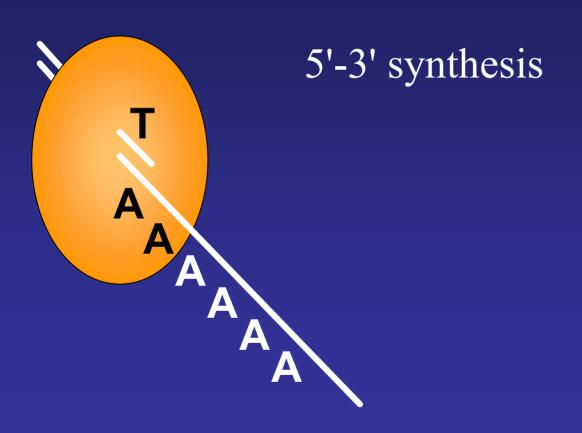


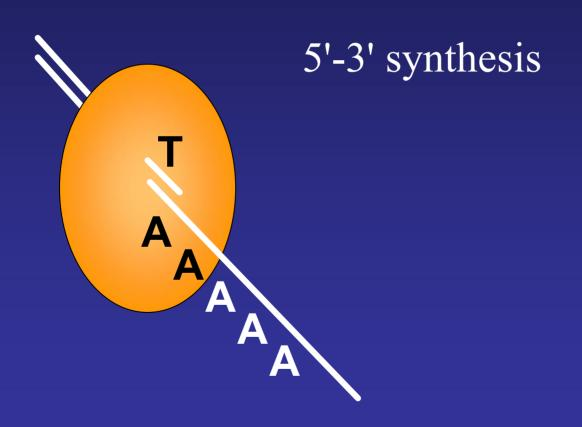
# AmpliTaq Gold DNA Polymerase

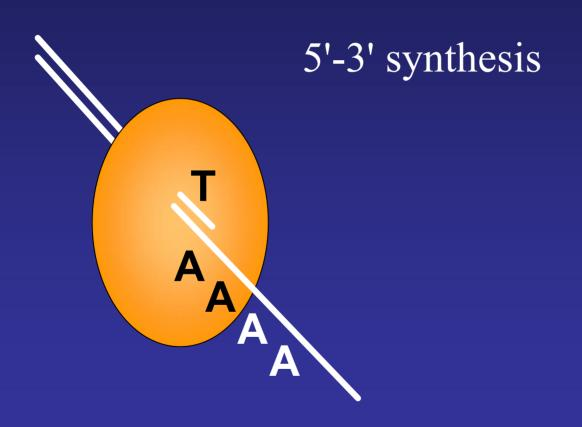
- Inactive until heated prior to thermocycling
- Simulates "Hot Start" PCR
- Increases yield & specificity of PCR
  - reduces non-specific products
  - reduces primer dimers
- Facilitates multiplexing
  - reduces primer design requirements
- Automation

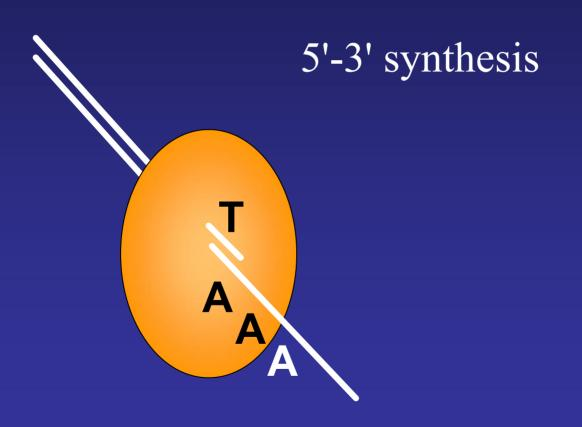


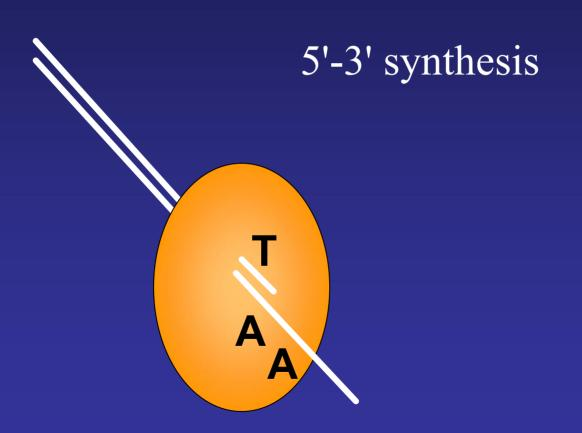




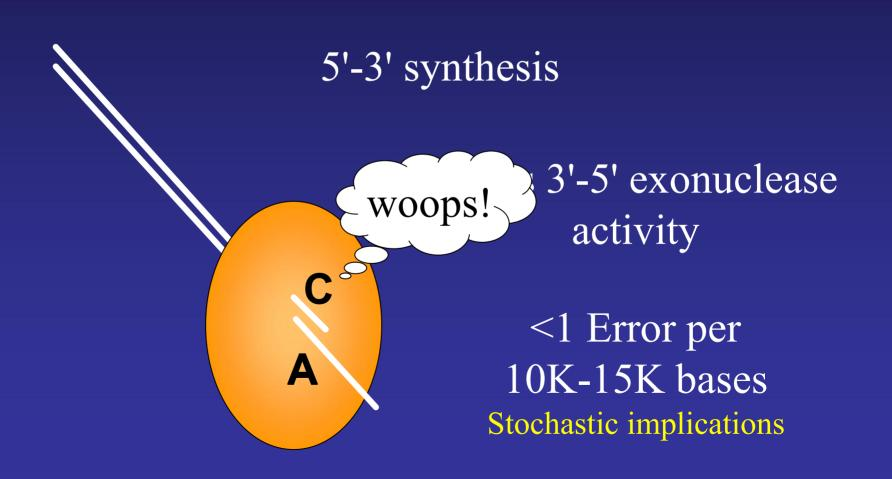






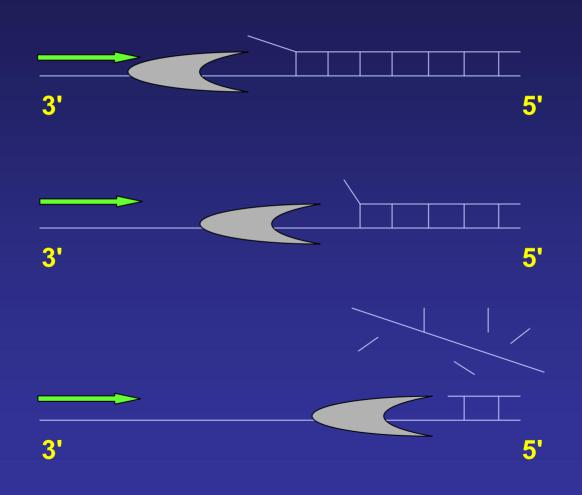


#### No Proofreading Ability



#### No 3' - 5' exonuclease activity

#### 5' --- 3' Exonuclease Activity



#### Template

- Sensitive to a single copy of template
- Various types
  - DNA: genomic, cloned, bacterial, viral, mitochondrial
  - RNA / cDNA
- Various sources
  - Blood, semen, saliva, hair, bone, teeth, etc.
  - Stained/unstained slides, parafin-embedded tissue
- Highly purified or crude lysate

# Primers

- Sequence
  - random base distribution
  - no poly-purine or pyrimidine stretches
  - $-\sim50\%$  G+C content
- Length
  - typically 18-28 bases
- Balance T<sub>m</sub> of all primers in a reaction
  - $-T_{\rm m} = 2 \times (A+T) + 4 \times (G+C) {\rm simplistic}$

# Primers

 Avoid 3' complementarity among primers, which can result in 'primer dimer' formation

• Primer dimer = template-independent duplex PCR product comprised of primers

# Primers

• Inter-primer 3' complementarity (two different primers)

5'-TTTTTTTCCCCC GGGGGTTTTTTT-5'

• Intra-primer 3' complementarity (one primer complementary to itself)

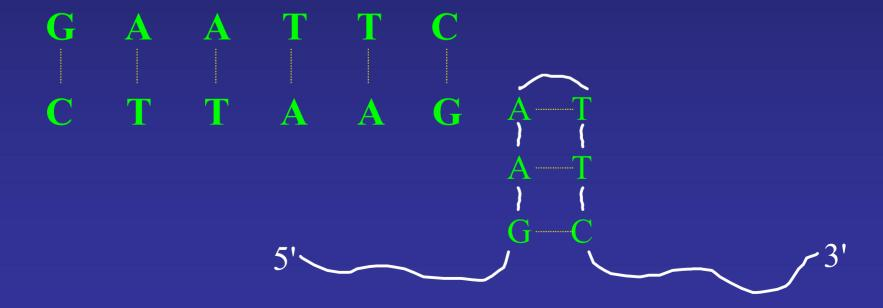
5'-TTTTTTCCCGGG GGCCCTTTTTTT-5'



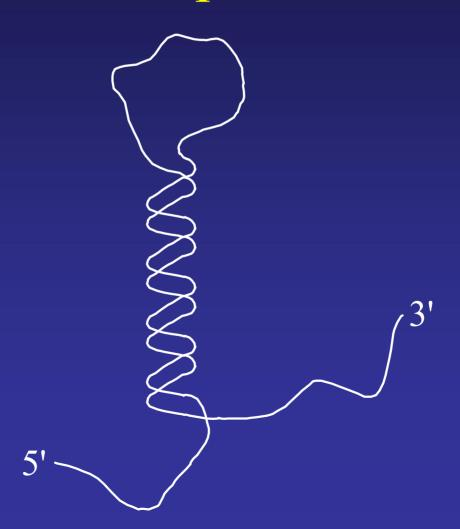
- Avoid runs of >3 G's or C's at 3' end
  - may misprime in G+C-rich regions
- Empirically determine ability to adenylate
- 5' additions to primers
  - restriction site, M13 site, promoter sequence
- biotin, fluorophore, radioisotope



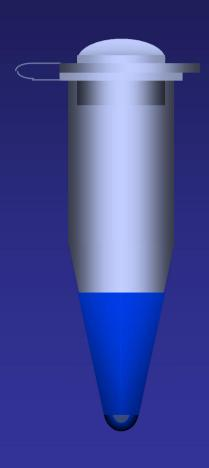
• Avoid palindromic sequences and internal complementarity



# Hairpin due to Internal Complementarity



#### Typical Reaction (25-100 ml)



- 50 mM
- 10 mM
- 1.25-1.50 mM
- 100 mg/ml
- 0.25 mM
- 200 mM
- 1-5 units

KC1

Tris-HCl

MgCl<sub>2</sub>

Gelatin

Each primer

Each dNTP

Taq Polymerase

**Yield** 

**Specificy** 

## OPTIMIZATION

**Fidelity** 

#### DNA CONCENTRATION



Depends of method of detection

Too Much DNA:



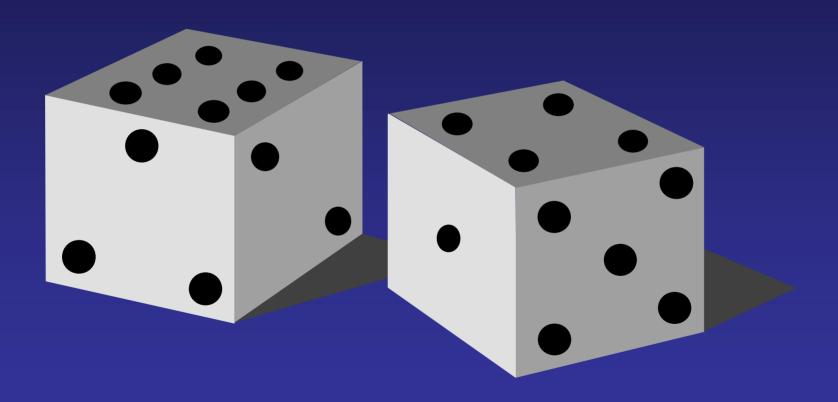
Non-specific products
Lower yield
(WHY?)

Too Little DNA:



Lower yield
Stochiastic effects

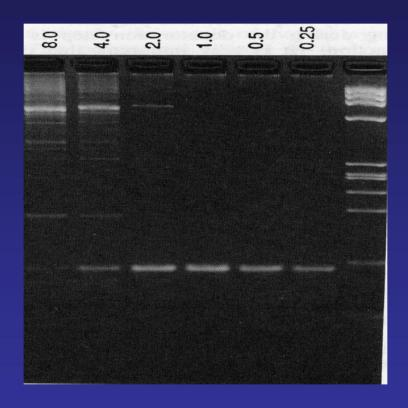
#### **Stochastic Effects**



#### TAQ CONCENTRATION

- Additional Taq may be recommended to overcome the effects of Taq polymerase inhibitors that may co-extract with DNA from forensic biospecimens
- However, too much Taq may affect specificity
- Too little Taq may result in lower yields

## VARYING TAQ POLYMERASE CONCENTRATION



5 units per reaction is typical for many forensic kits

#### **Divalent Cations**

Magnesium is the cation of choice for Taq polymerase

Cofactor in the catalytic addition of deoxynucleoside monophosphates to the 3' end of the growing DNA chain

So why 1.5 mM magnesium?

 $0.8 \text{ mM Mg}^{2+}$ 

is bound by dNTPs (200 uM x 4 = 0.8 mM)



 $0.7 \text{ mM Mg}^{2+}$ 

is needed by Taq

0.7 mM + 0.8 mM 1.5 mM



#### MgCl<sub>2</sub> Concentration



Typically
1.25 - 1.50
mM

Too High:



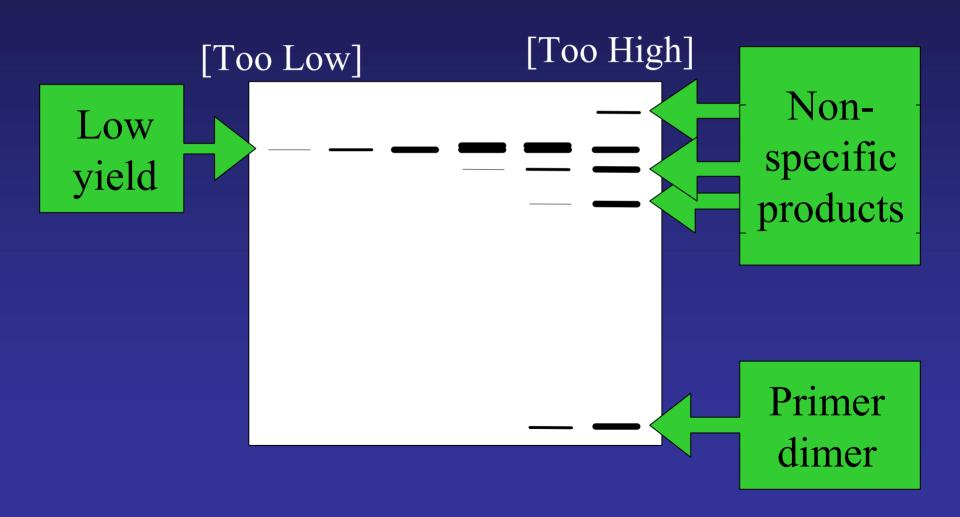
Non-specific products
Inhibits Taq
(WHY?)

Too Low:



Lower yield

#### PRIMER CONCENTRATION



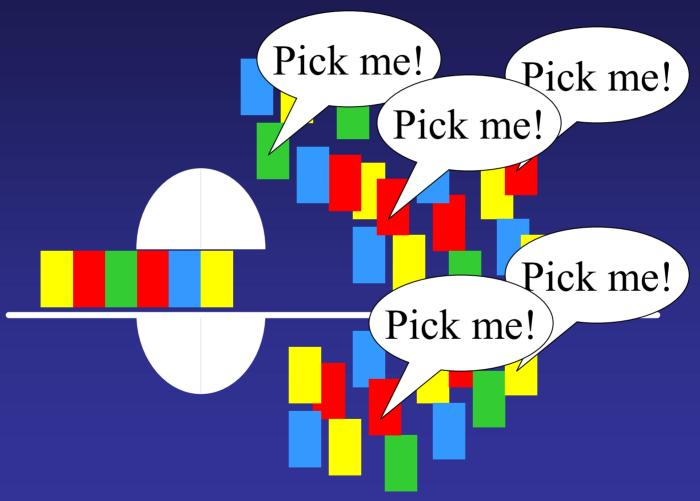
## Deoxynucleoside Triphosphates (dNTPs)

The four dNTPs are present in the PCR

Concentrations range from 20 to 200 µM each

A concentration of 200  $\mu$ M of each dNTP is theoretically sufficient to synthesize 25 $\mu$ g

#### dNTP CONCENTRATION



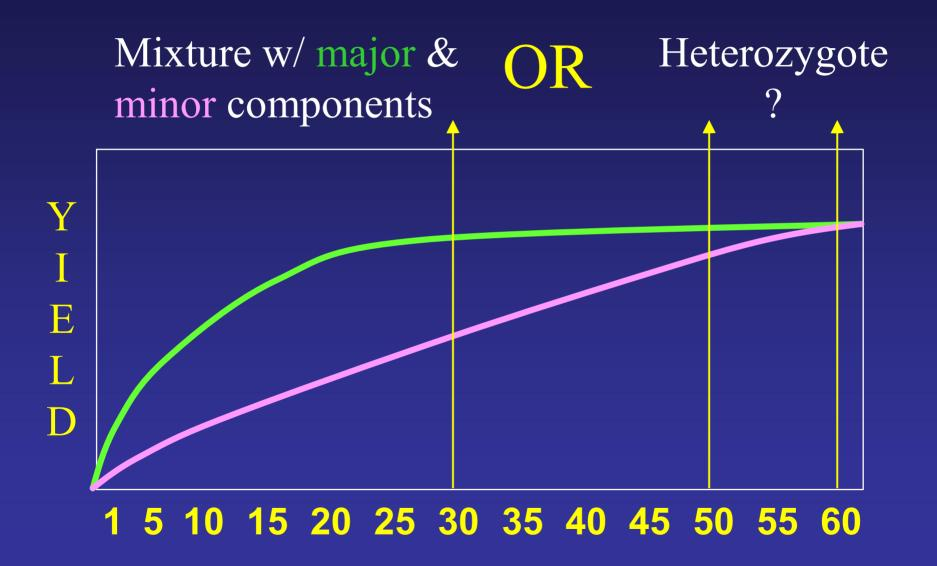
#### CYCLE NUMBER

Non-specific products and primer dimers may be detected if too many cycles are used

Depends on target DNA quantity

Copy Number	Cycle Number
300,000	25-30
15,000	30-35
1,000	35-40
50	40-45

#### CYCLE NUMBER & PLATEAU



#### DENATURATION

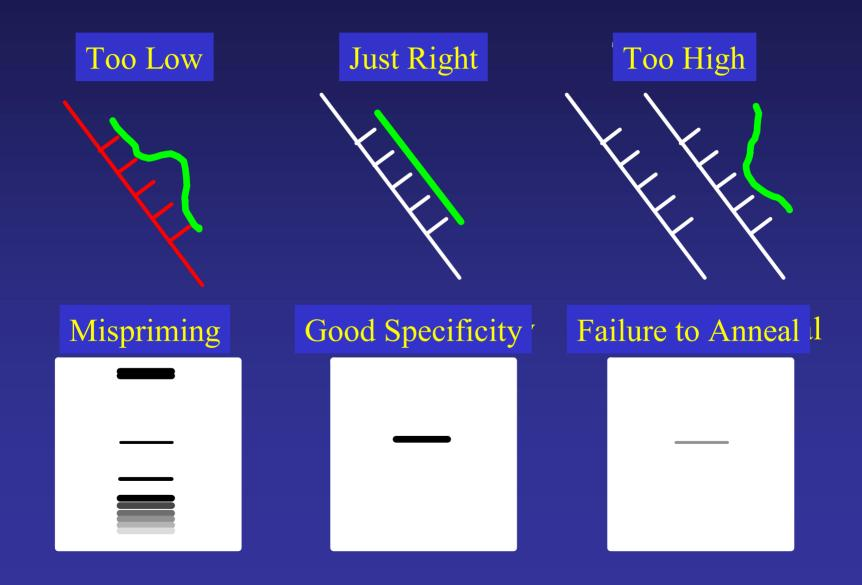
If temperature is too low:



GC-rich regions may fail to denature. Primers may not bind.

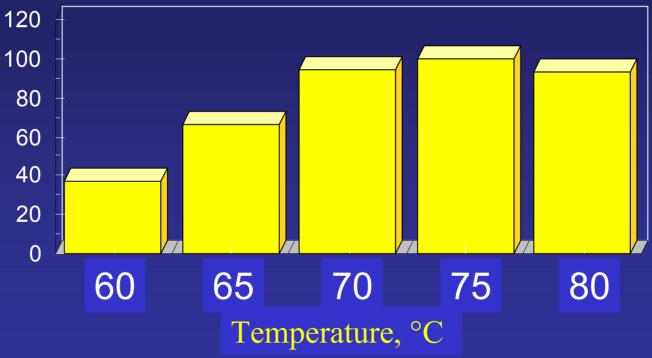


#### ANNEALING TEMPERATURE



#### **EXTENSION**





70-72°C recommended

# FORENSIC PCR Template Integrity

- Forensic samples may be:
  - old, degraded
  - exposed to environmental / chemical insults
  - contain PCR inhibitors

• However, typically in forensic PCR, small regions of DNA that can be amplified from degraded samples are targeted.

## FORENSIC PCR PCR Inhibitors



- Act by binding DNA target or inhibiting Taq.
- DNA sources & substrates that may contain inhibitors:
  - Blood... Heme

Heparin (anticoagulant)

- Hair... Melanin (pigment)
- Fabrics... Dyes (i.e., indigo)
- Soil... Humic acid (organic breakdown product)
- Urine... Metabolic substances, drugs

### Strategies for overcoming the effects of inhibitors

- Dilute out inhibitor by increasing the total PCR volume (but less DNA)
- Chelex extract DNA (or other method)
- Use microconcentrator
- Increase concentration of Taq
- Add BSA to PCR



#### Bovine Serum Albumin

• Does not affect PCR when no inhibitor

 Helpful with samples containing Taq inhibitors

 Degree of affect depends on manufacturer and fraction of BSA

#### FORENSIC PCR

Use at a concentration of 160 ug / ml

Sigma Fraction V (#A3350 or #A2153)

- = 16 ug for 100 ul reactions
- = 8 ug for 50 ul reactions

## FORENSIC PCR Multiplexing

• Amplification of more than one locus in a single PCR tube.

#### • Decreases:

- number of manipulations
- chance of sample mix-up
- chance of contamination
- DNA and reagent consumption

### FORENSIC PCR Contamination

- Sensitivity of PCR: single DNA copy
- FROM THE ENVIRONMENT:
  - nature of the crime
  - handling in the field or lab
- BETWEEN SAMPLES:
  - during preparation
- PCR-PRODUCT CARRY-OVER:
  - a serious concern because amplicons are usually in h concentration and are ideal substrates.

#### PCR Product Carry-Over

0.1 µl carry-over of amplicon can contain a MILLION COPIES of DNA



#### Real-Time PCR

• Monitor the accumulation of PCR product during amplification

• Detection of changes in fluorescent signal generated during cycles of PCR

\* The fewer cycles to detect fluorescence signal the greater amount of DNA in the sample

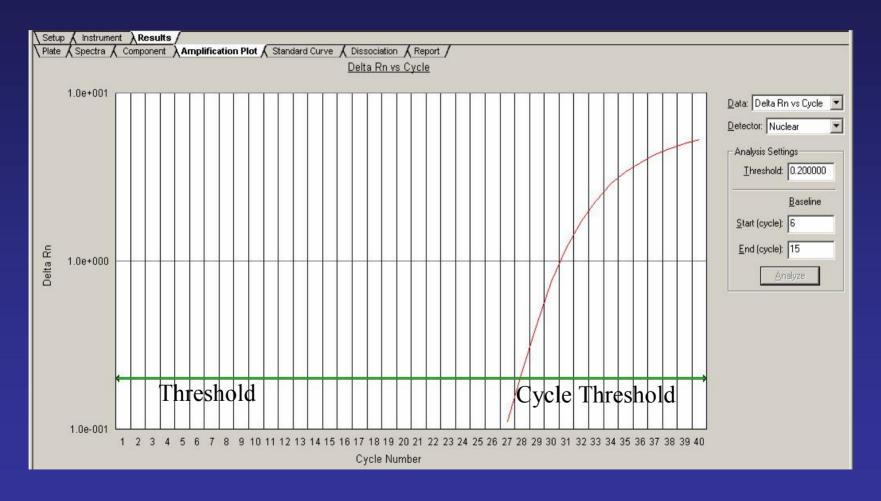
Simultaneous amplification and detection in the same well

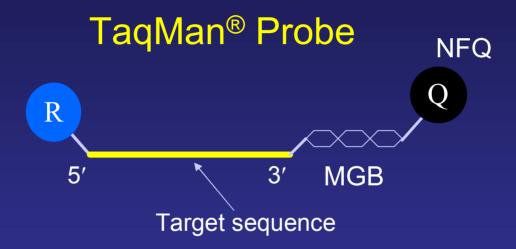
#### Real Time Definitions

Threshold – The level of detection or the point at which a reaction reaches a fluorescent signal above baseline (7000 and 7900HT SDS default settings are 0.2)

 $\underline{C_T}$  (Cycle Threshold) – The cycle at which a sample amplification curve crosses the threshold

# Example of Accumulation of PCR Product (fluorescence) per Cycle

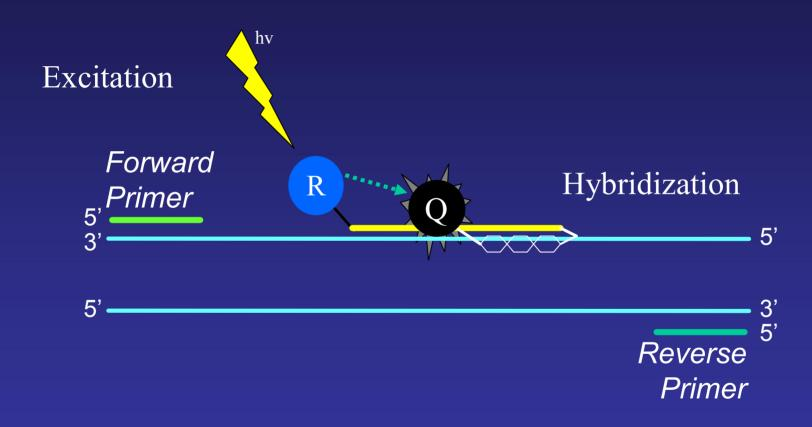


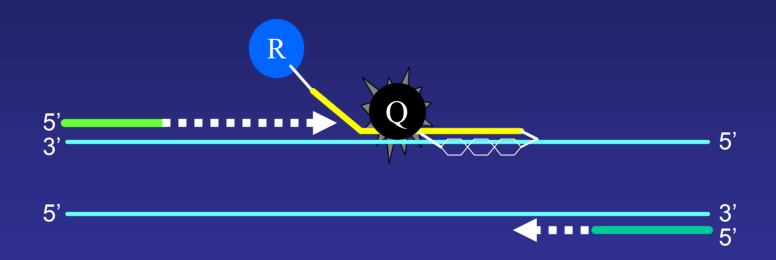


R = Reporter (FAM<sup>TM</sup> or VIC<sup>®</sup> Dyes)

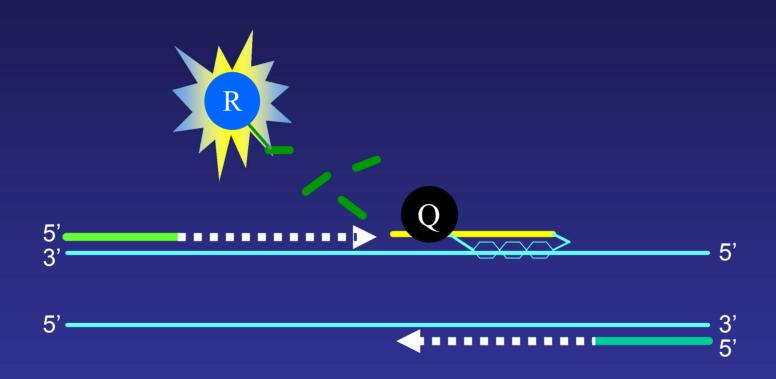
Q = Non Fluorescent Quencher (NFQ)

Acts as energy transfer acceptor that does not emit a detectable fluorescent signal

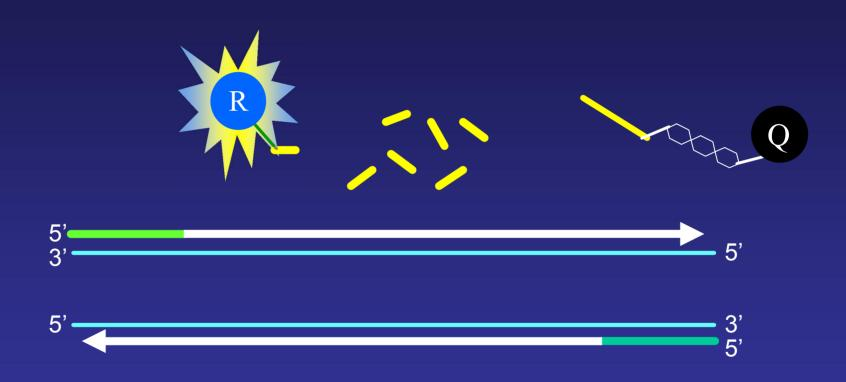




Displacement



Cleavage

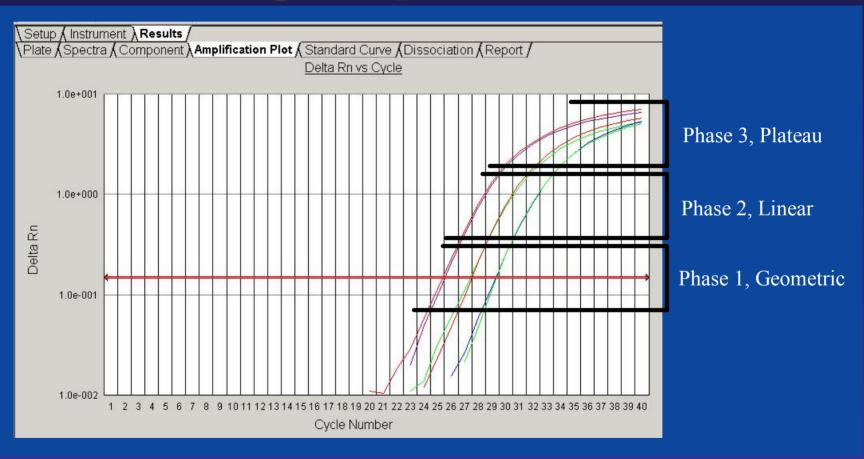


Polymerization completed

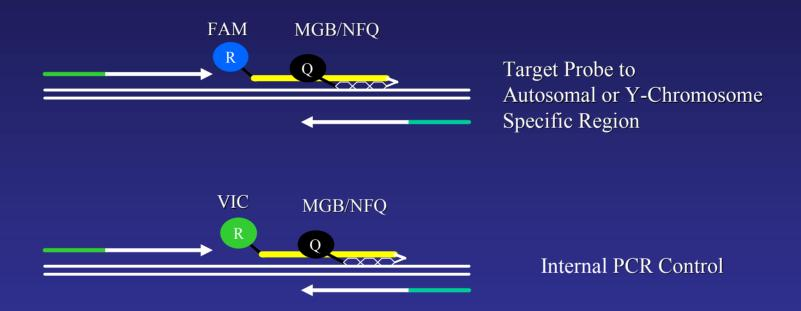
# Advantages of Using the 5'-Nuclease Assay

- 1. The probe provides specificity
- 2. Low background noise increases sensitivity
- 3. The fluorescent signal is directly proportional to the number of amplicons generated

### Monitoring Real-Time PCR [three phases]

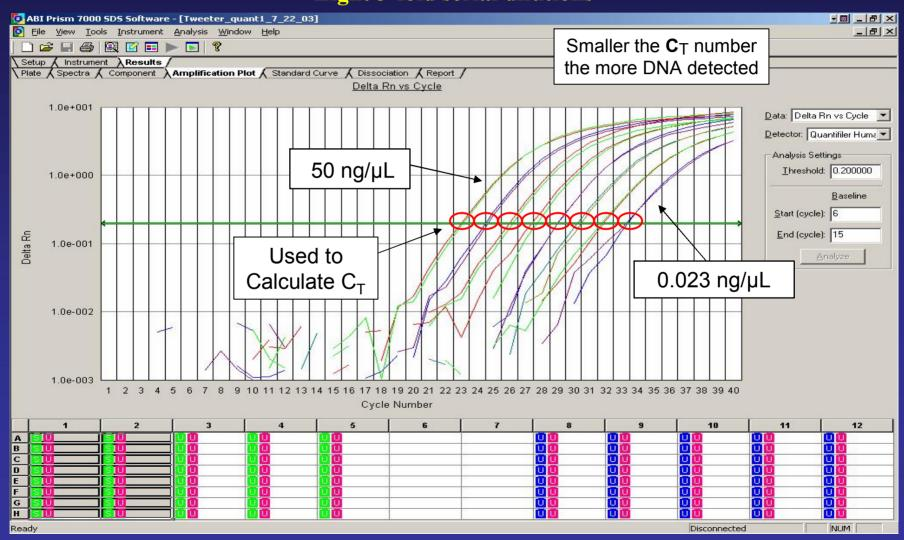


#### Quantifiler Primer Mix: Pre-Formulated as A Duplex Assay

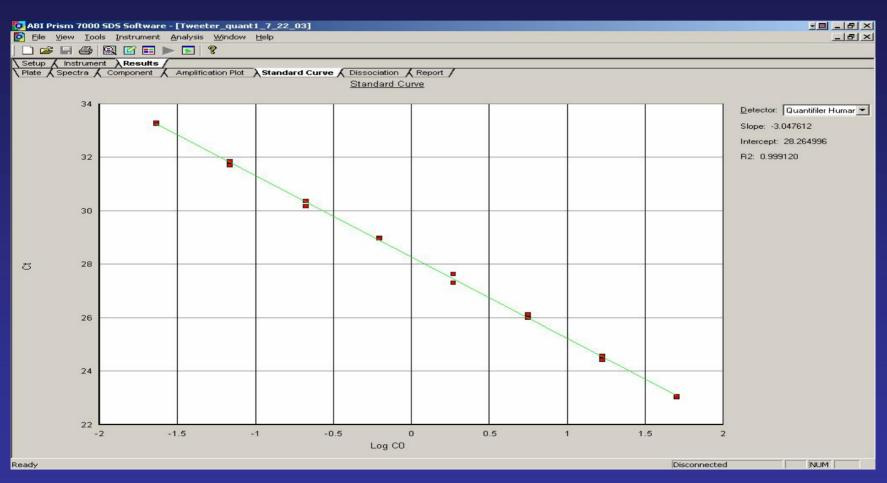


### Amplification Plots for DNA Concentration Standards

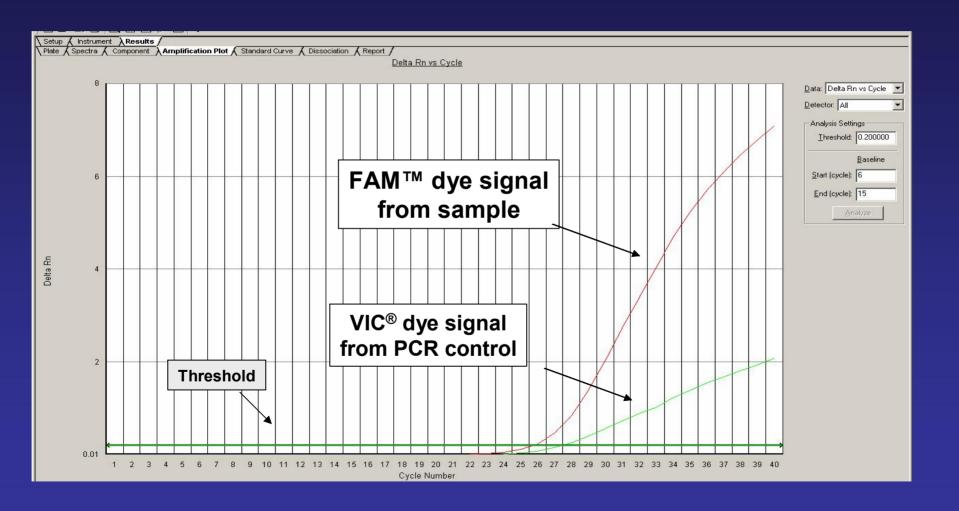
**Eight 3-fold serial dilutions** 



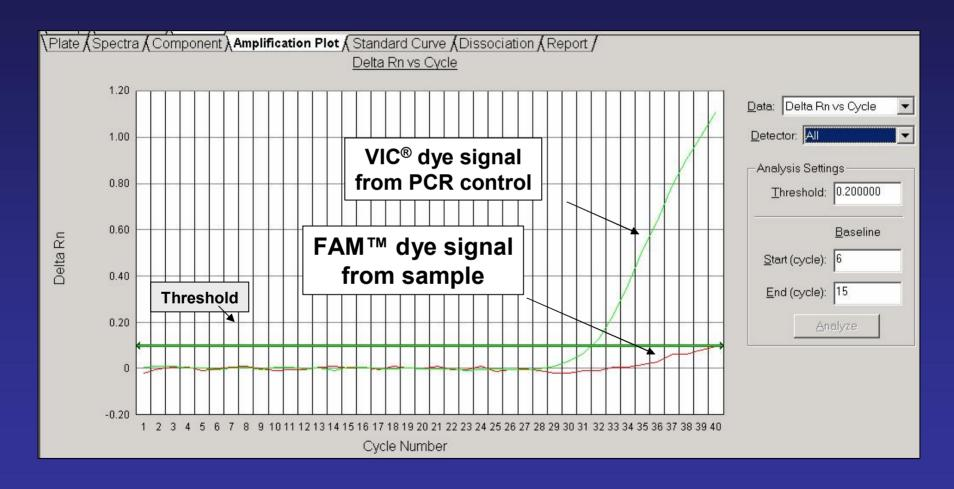
### Standard Curve from DNA Concentration Standards



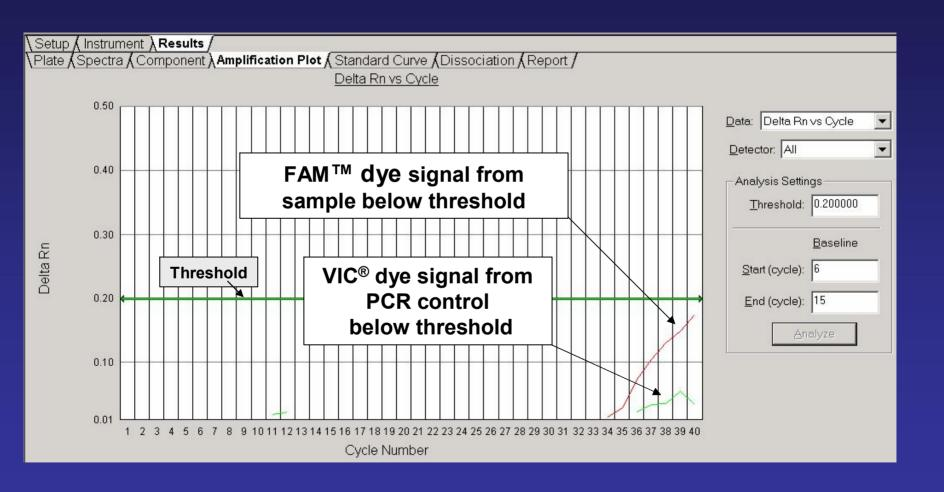
#### Example of sample positive for DNA



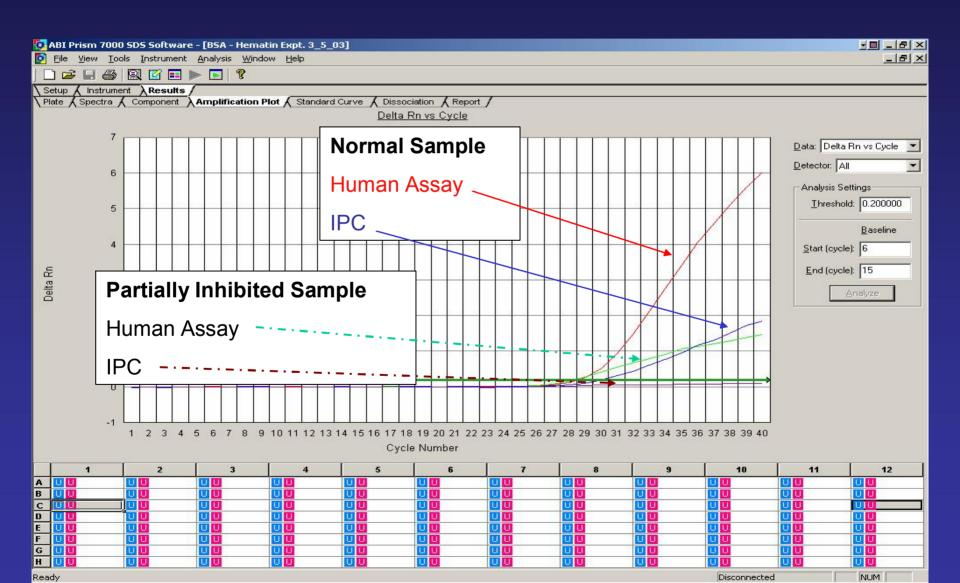
### Example of sample with no detectable DNA



#### Example of sample with inhibition



#### Example of sample with partial inhibition



### Application

- Quantitation
- SNP detection