

# Introduction to Real-Time PCR: Basic Principles and Chemistries

Leta Steffen, PhD

Applications Scientist

# Outline

---

- I. Real-Time PCR overview
  - Basics of Real-Time PCR
  - Understanding the data
- II. Chemistries and Instrumentation
  - Dye-based chemistries (e.g. SYBR)
  - Label-based chemistries (e.g. TaqMan)
  - Reverse Transcription qPCR
  - Instrumentation
- III. Quantification – an intro to the math
- IV. The MIQE Guidelines

# PCR Refresher

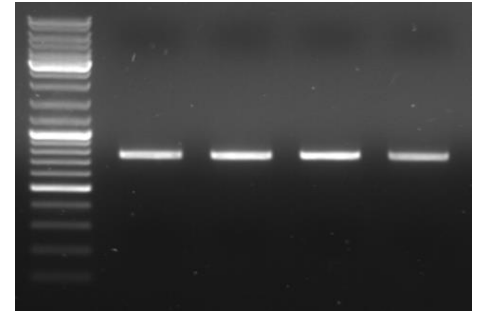
Taq  
dNTPs  
MgCl<sub>2</sub>  
Buffer  
Primers  
Template



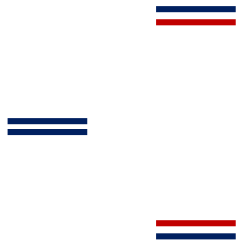
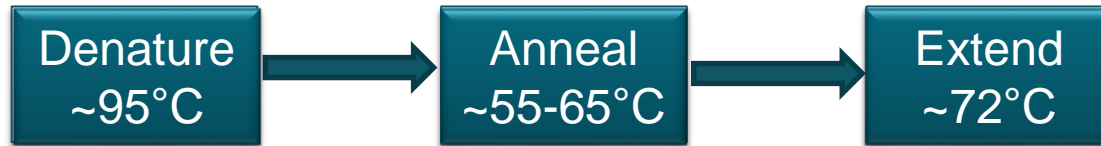
Stage	Temp	Time	Cycles
Activation	95°C	2 min	1
Dissociation	95°C	15 s	25 - 40
Annealing	60°C	15 s	
Extension	72°C	1 min	



Gel electrophoresis

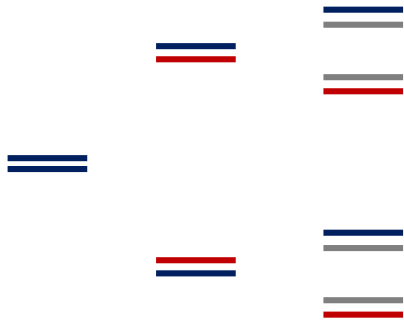


# PCR Refresher



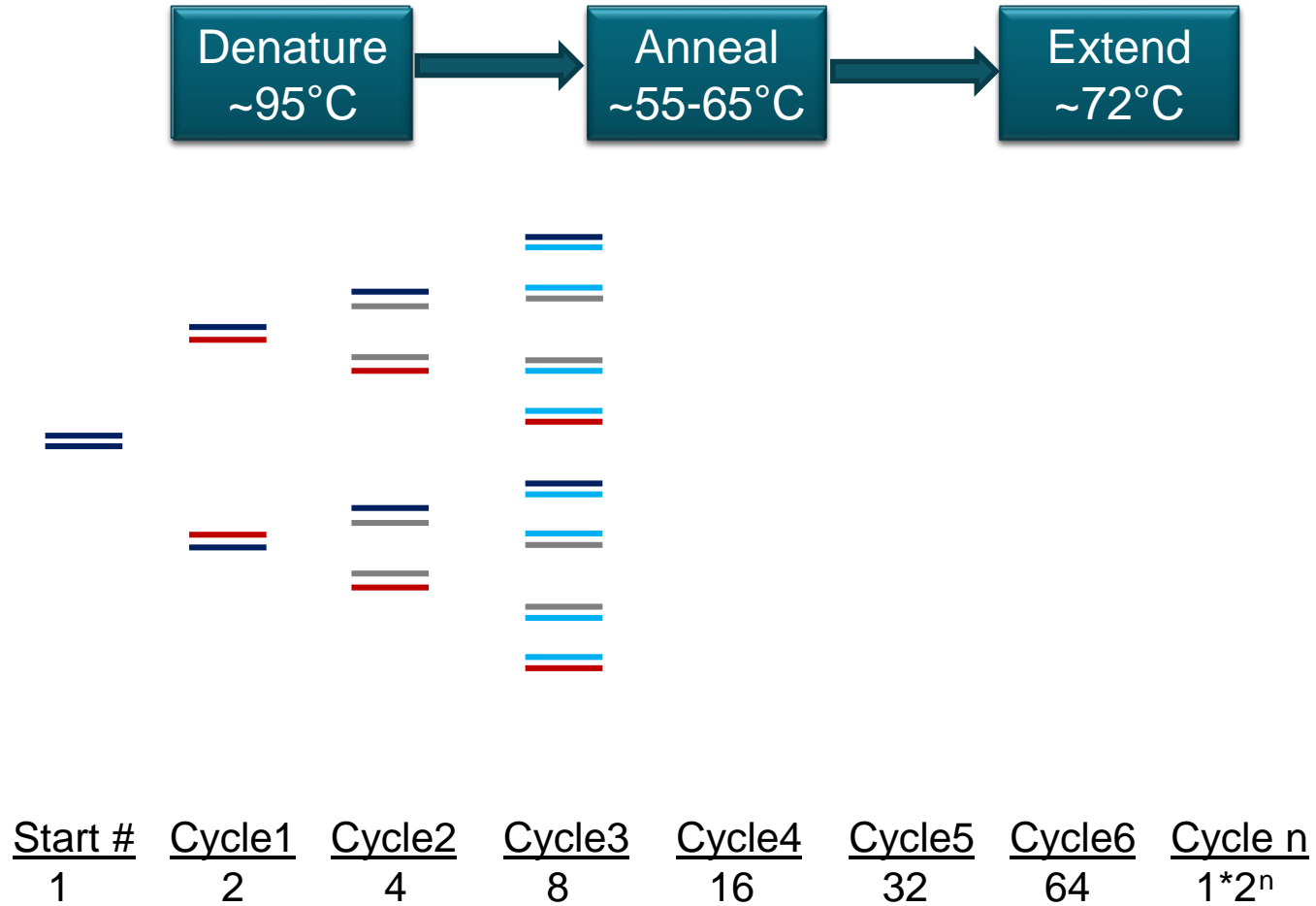
<u>Start #</u>	<u>Cycle1</u>	<u>Cycle2</u>	<u>Cycle3</u>	<u>Cycle4</u>	<u>Cycle5</u>	<u>Cycle6</u>	<u>Cycle n</u>
1	2	4	8	16	32	64	$1 \cdot 2^n$

# PCR Refresher

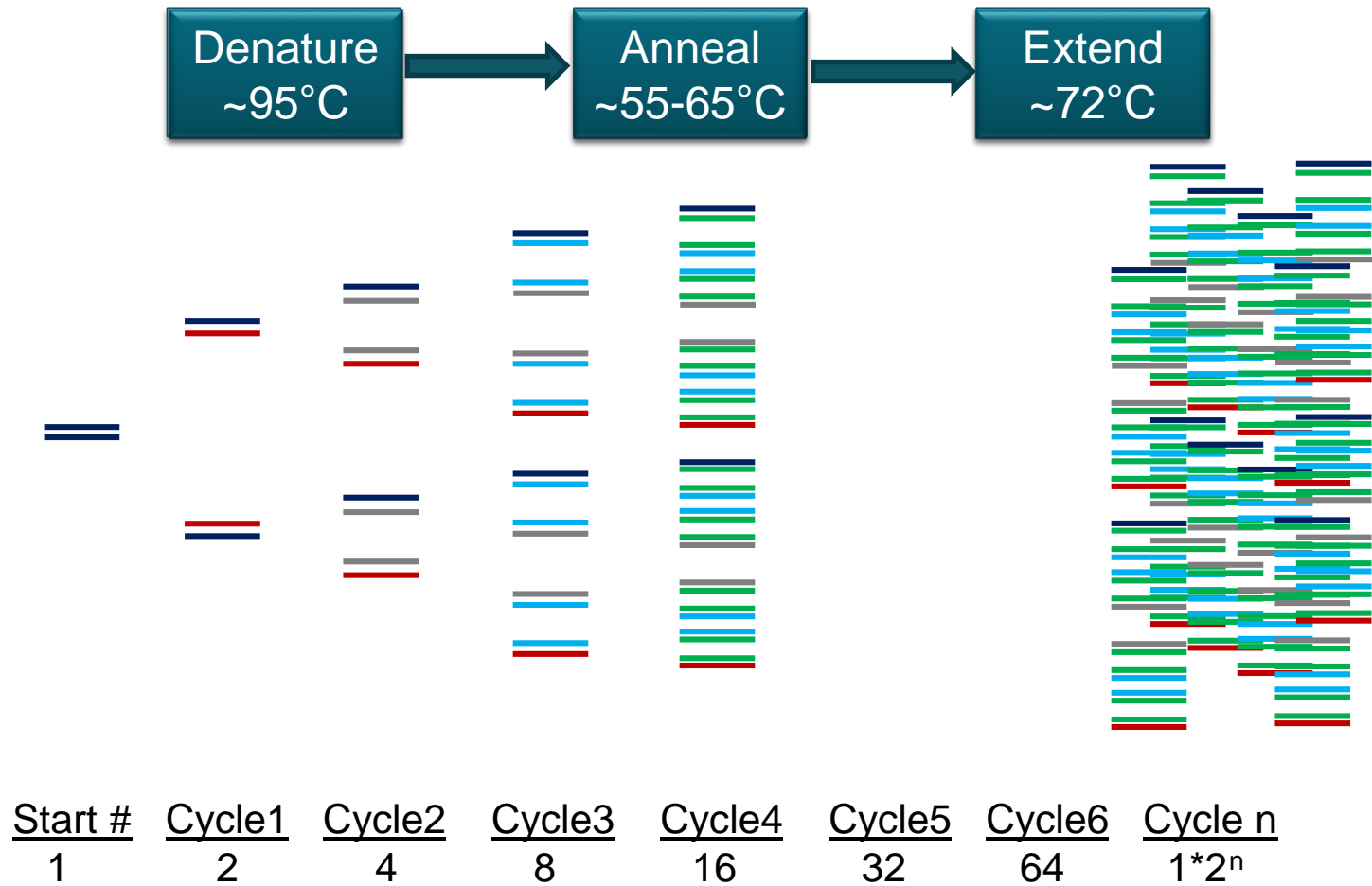


<u>Start #</u>	<u>Cycle1</u>	<u>Cycle2</u>	<u>Cycle3</u>	<u>Cycle4</u>	<u>Cycle5</u>	<u>Cycle6</u>	<u>Cycle n</u>
1	2	4	8	16	32	64	$1 \cdot 2^n$

# PCR Refresher

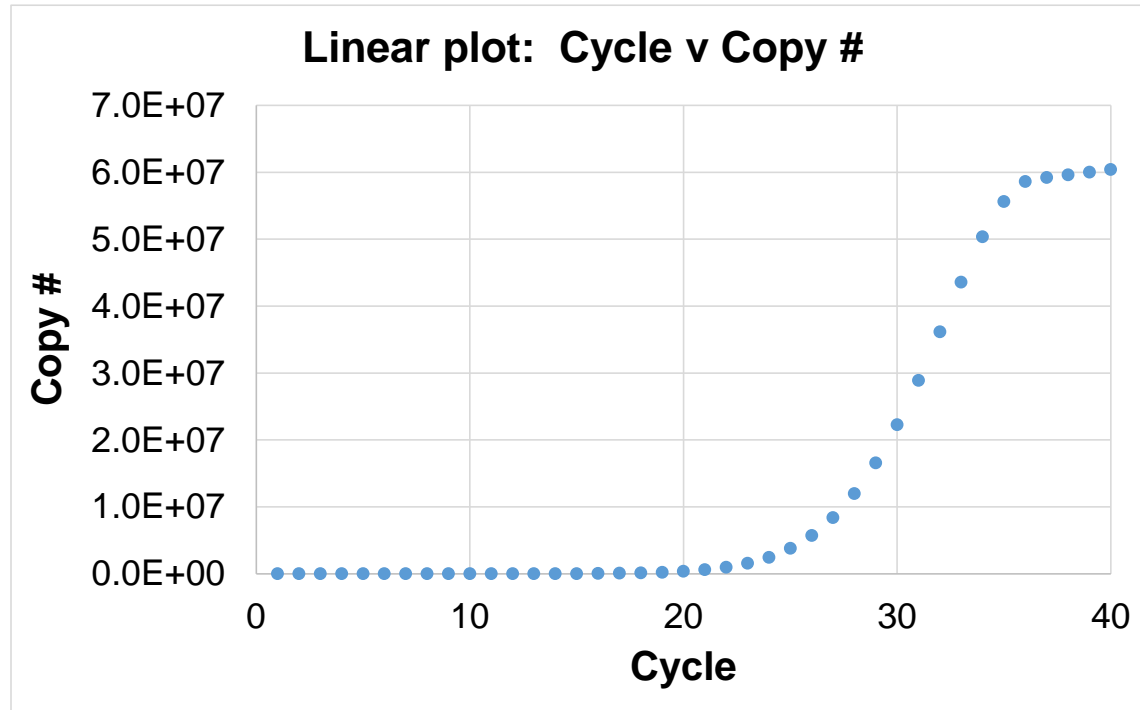


# PCR Refresher



# Traditional PCR

## *Dynamics of amplification*

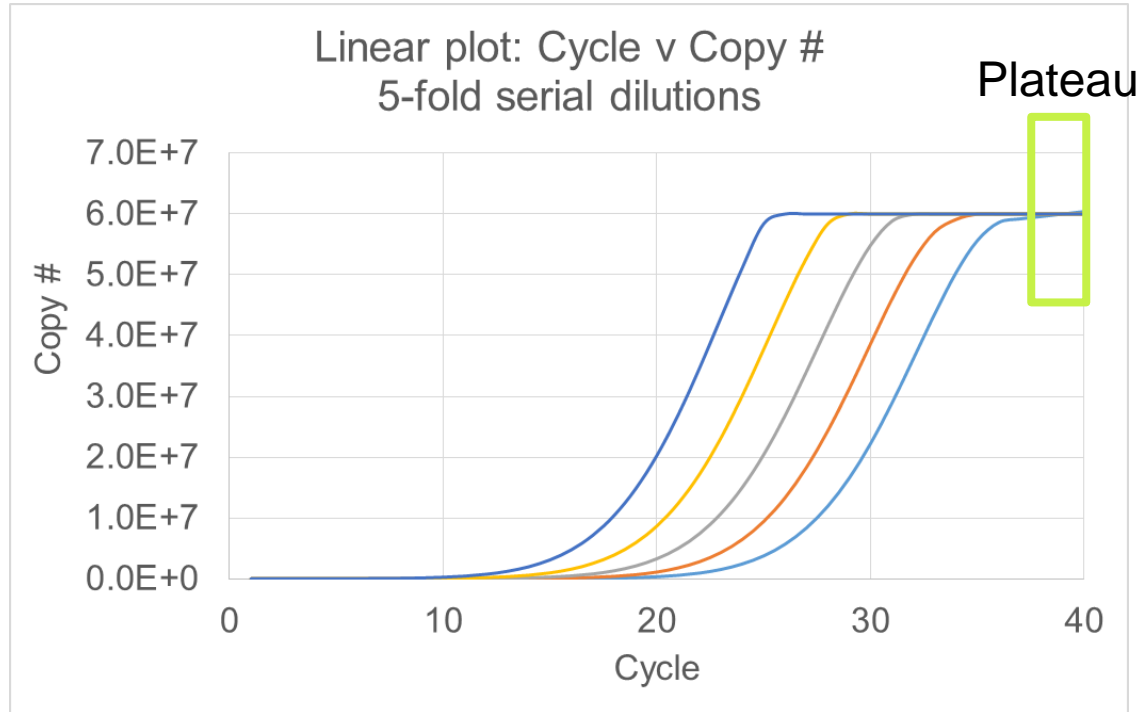


<u>Start #</u>	<u>Cycle1</u>	<u>Cycle2</u>	<u>Cycle3</u>	<u>Cycle4</u>	<u>Cycle5</u>	<u>Cycle6</u>	<u>Cycle n</u>
1	2	4	8	16	32	64	$1 \cdot 2^n$



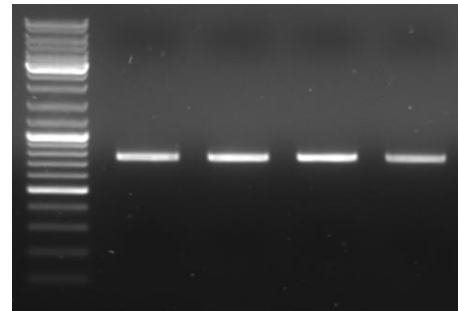
# Endpoint PCR

*Analyze reactions at plateau*



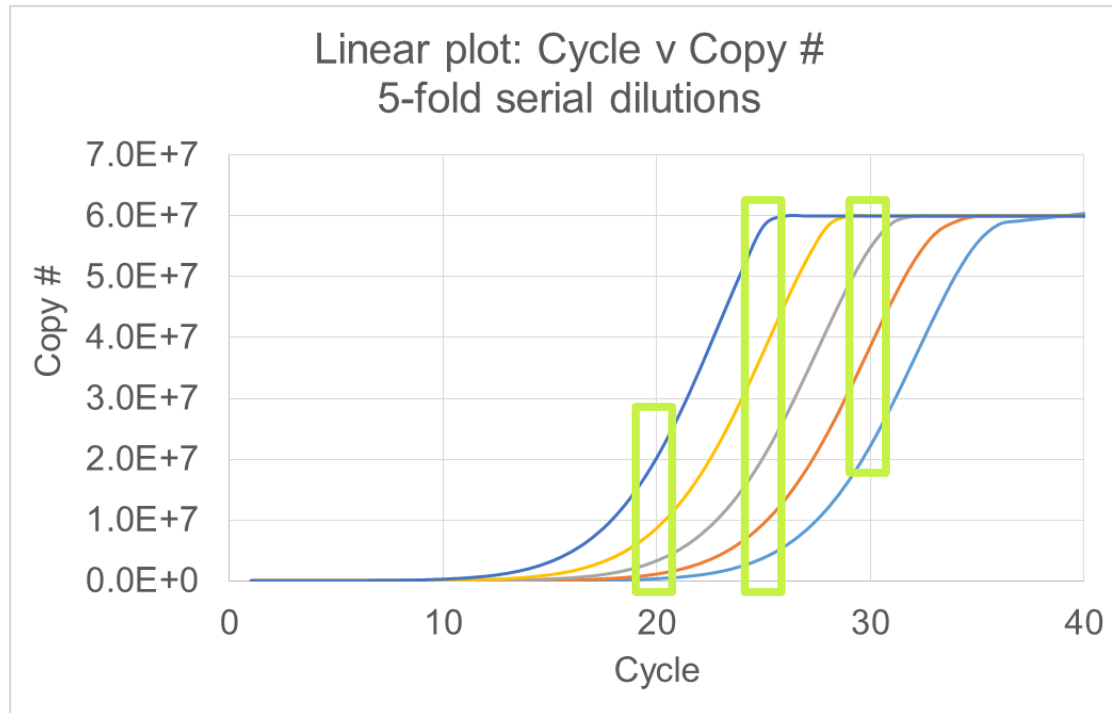
## Endpoint PCR

- Assayed at plateau
- Not quantitative
- Size & number
- Limited by gel sensitivity



# Semi-Quantitative PCR

*Analyze reactions at an intermediate cycle*



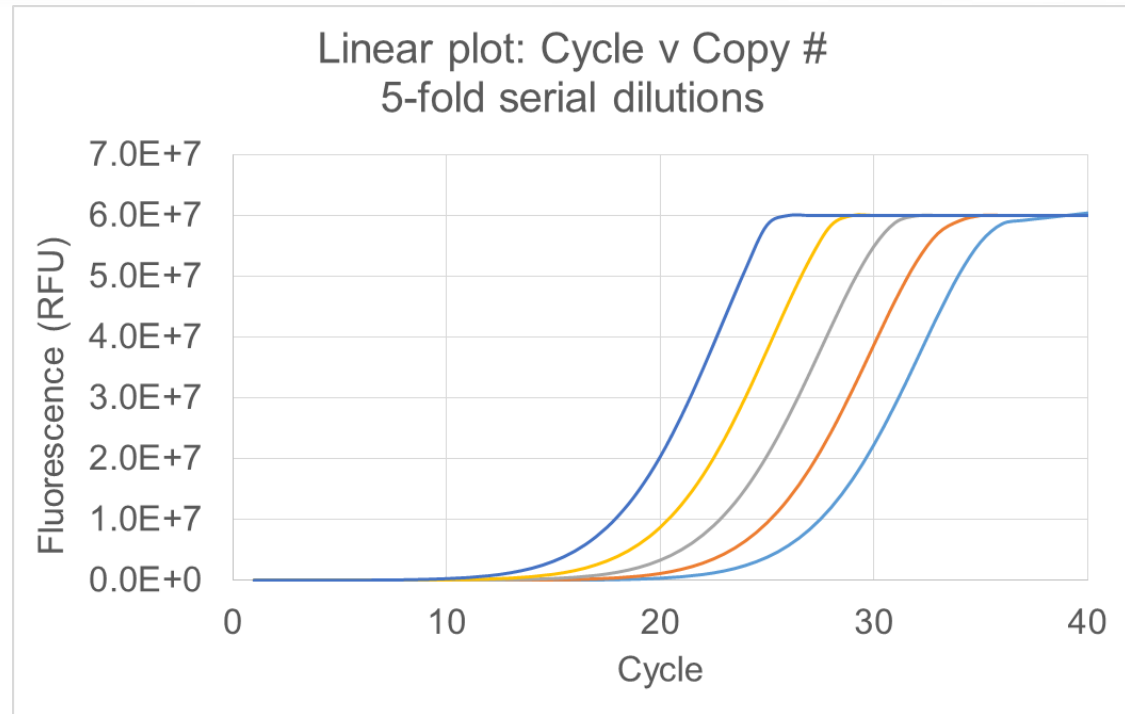
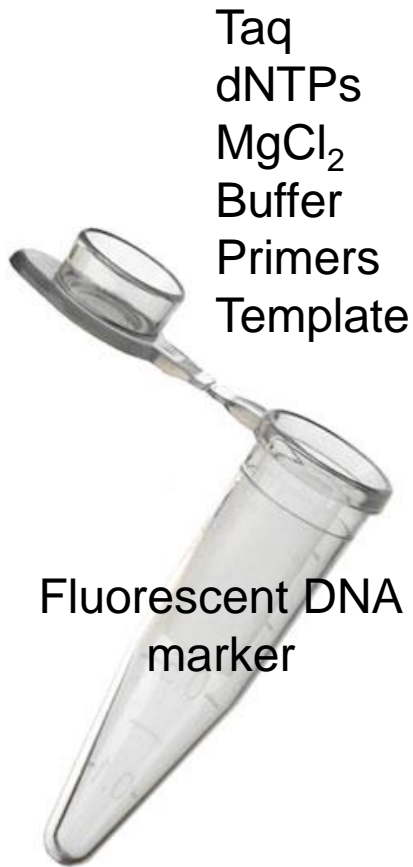
## Semi-quantitative PCR

- Assayed before plateau
- Not linear
- Limited dynamic range



# Real-time PCR

*Measure product at every cycle!*

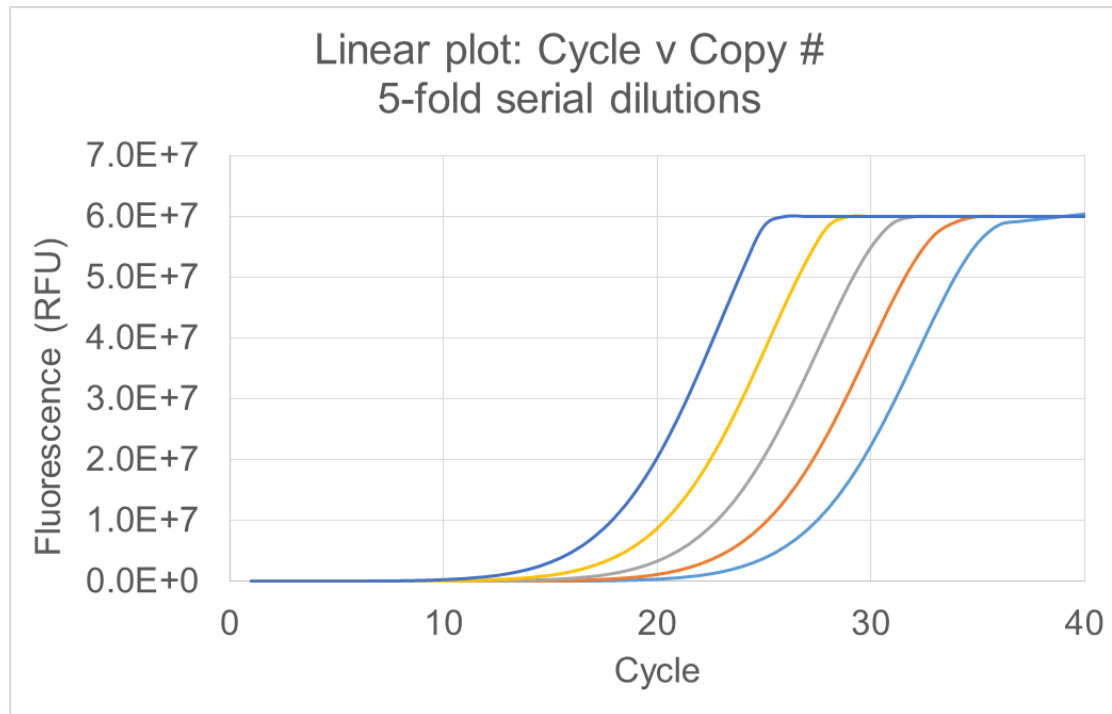


## Real-time PCR, quantitative PCR, qPCR

- Assayed at every cycle
- Requires specialized instrument
- No additional sample handling
- Broad dynamic range ( $10^6$ - $10^8$ )
- Quantitative
- High-throughput capable
- Multiplex capable
- Typically used for <250bp

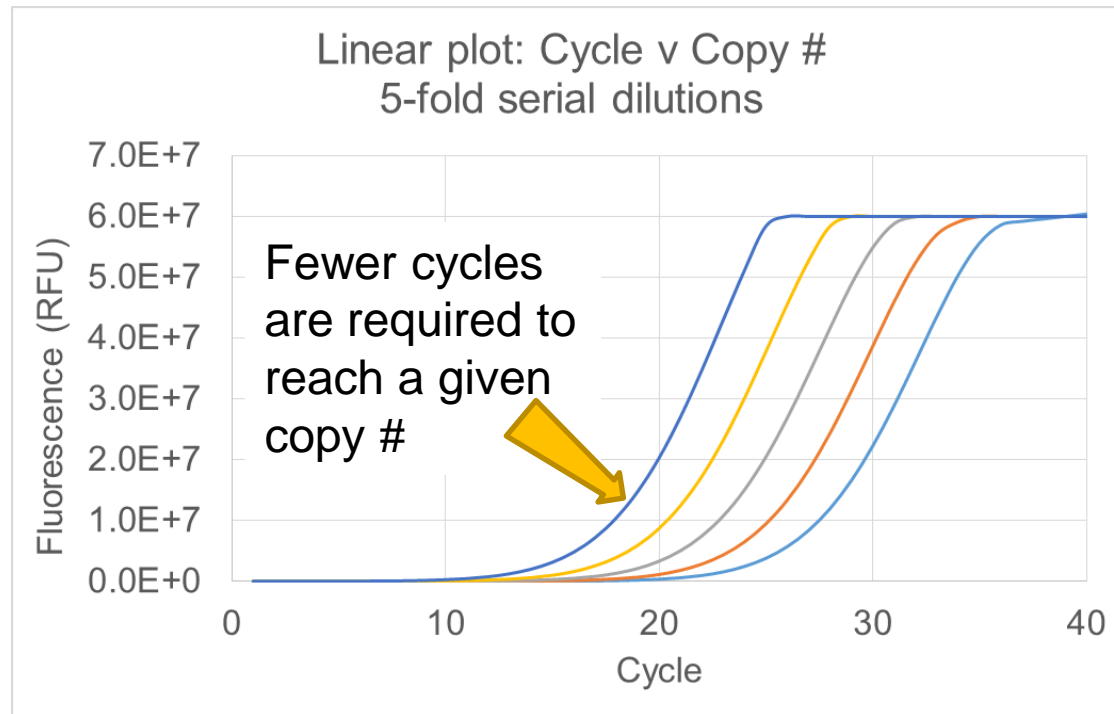
# Think About the Data

QUESTION 1: Which sample has a higher target concentration? Explain.



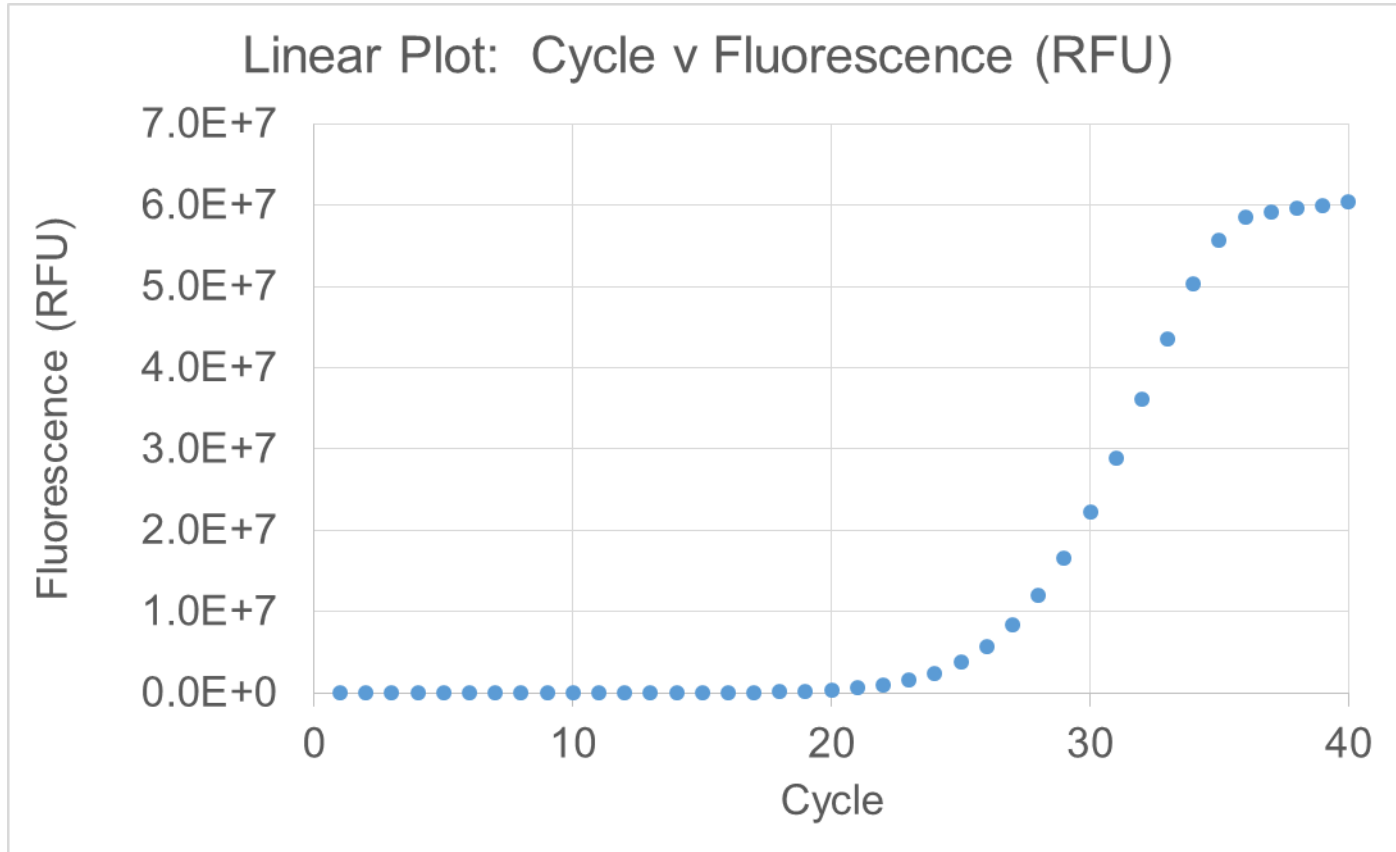
# Think About the Data

QUESTION 1: Which sample has a higher target concentration? Explain.

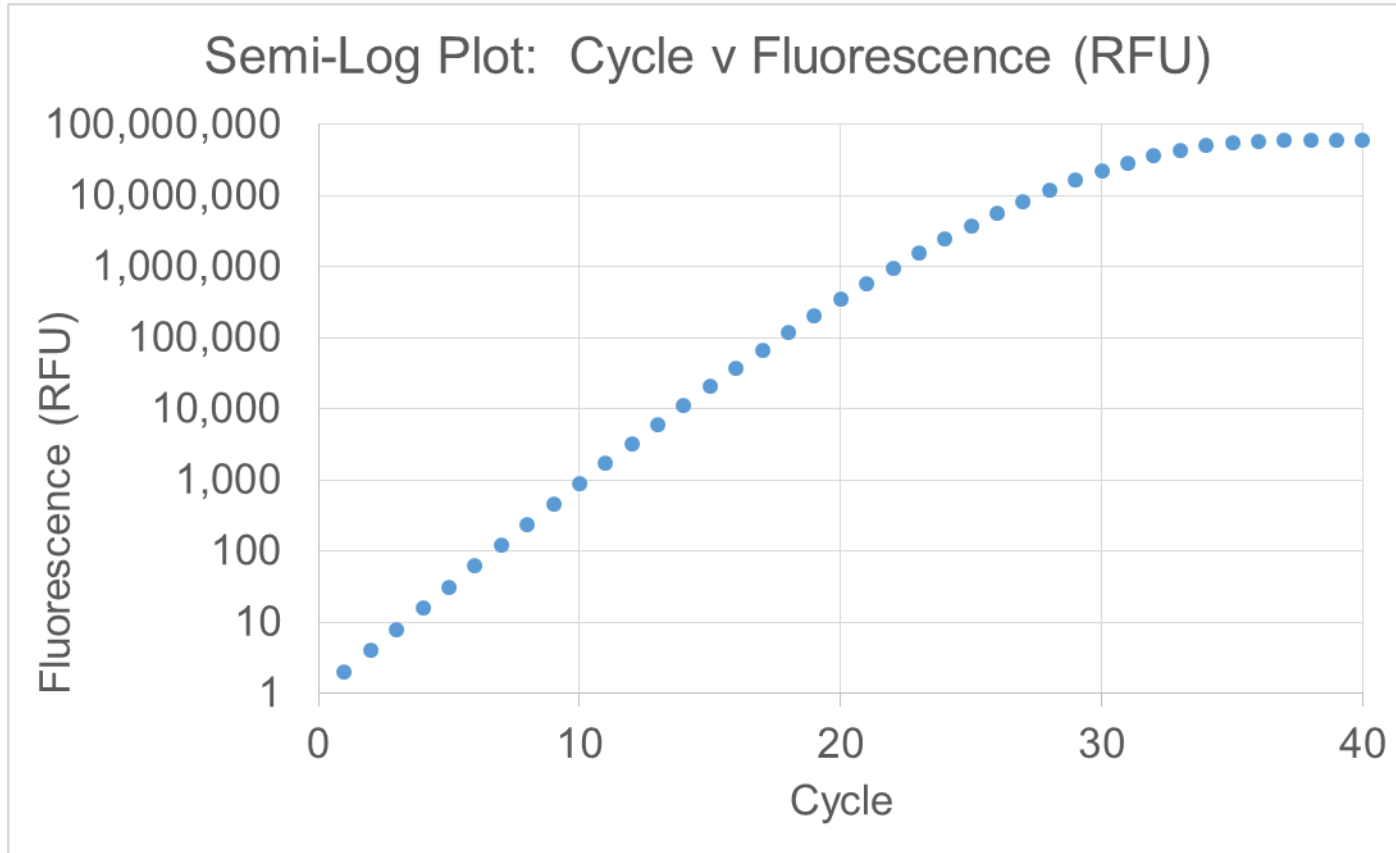


***Amplification is inversely proportional to starting concentration!***

# Viewing the Data



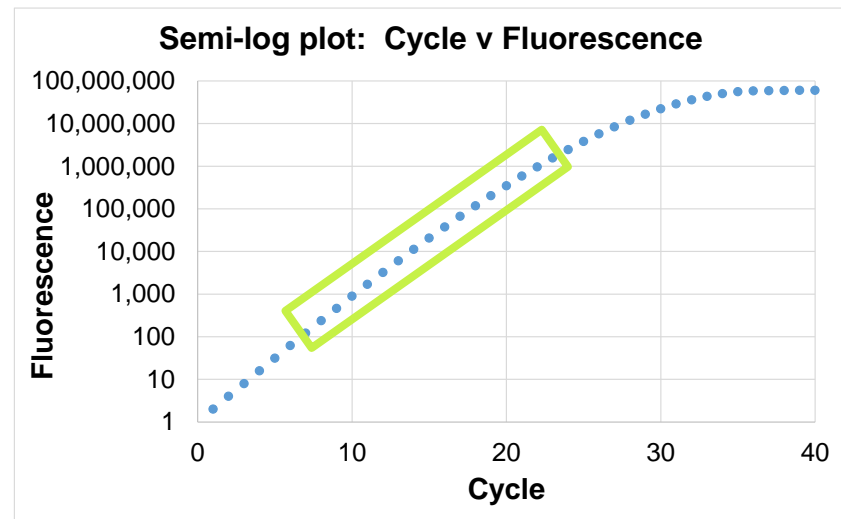
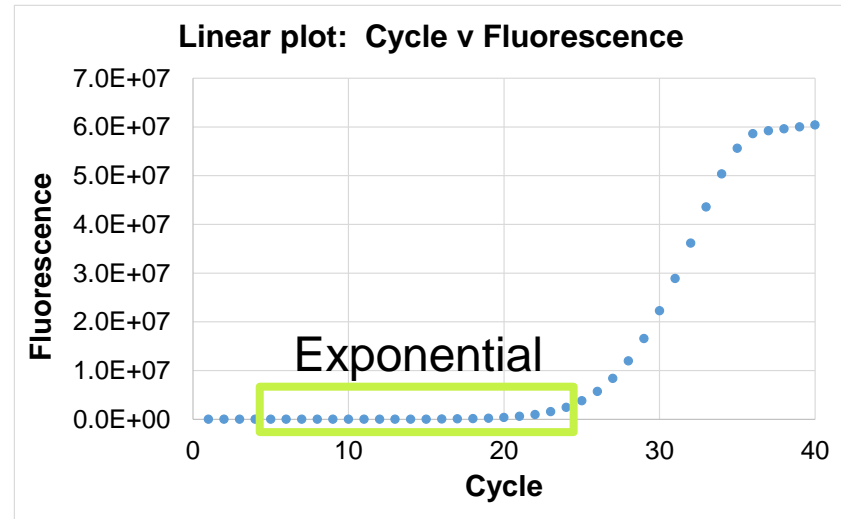
# Viewing the Data



# Viewing the Data

Quantification is most robust in the exponential phase of amplification

The exponential phase is best viewed in a semi-log plot





# From Raw Data to Quantification

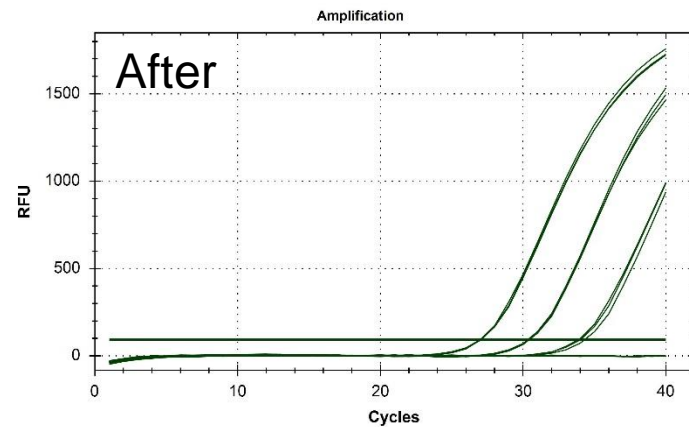
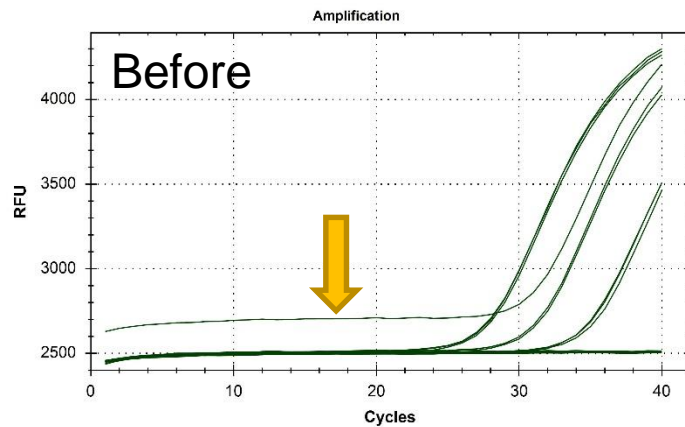
## *Software processing*

Step 0: Normalize to passive reference ( $R_n$ , normalized RFU)

- Only required on some platforms
- Use a second fluorescent dye, not influenced by DNA
- Divide reporter dye fluorescence by passive reference

Step 1: Subtract background fluorescence (RFU,  $\Delta R$ ,  $dR$ ,  $dR_n$ )

- Appropriate background cycles determined automatically
- Average background fluorescence for each reaction
- Subtract avg background fluorescence at each cycle
- De-trend data based on background fluorescence slope



# From Raw Data to Quantification

## *Software processing*

---

Step 0: Normalize to passive reference ( $R_n$ , normalized RFU)

- Only required on some platforms
- Use a second fluorescent dye, not influenced by DNA
- Divide reporter dye fluorescence by passive reference

Step 1: Subtract background fluorescence (RFU,  $\Delta R$ ,  $dR$ ,  $dR_n$ )

- Average background fluorescence for each reaction
- Subtract avg background fluorescence at each cycle
- De-trend data based on background fluorescence slope
- Appropriate background cycles determined automatically

Step 2: Fit a curve to the data

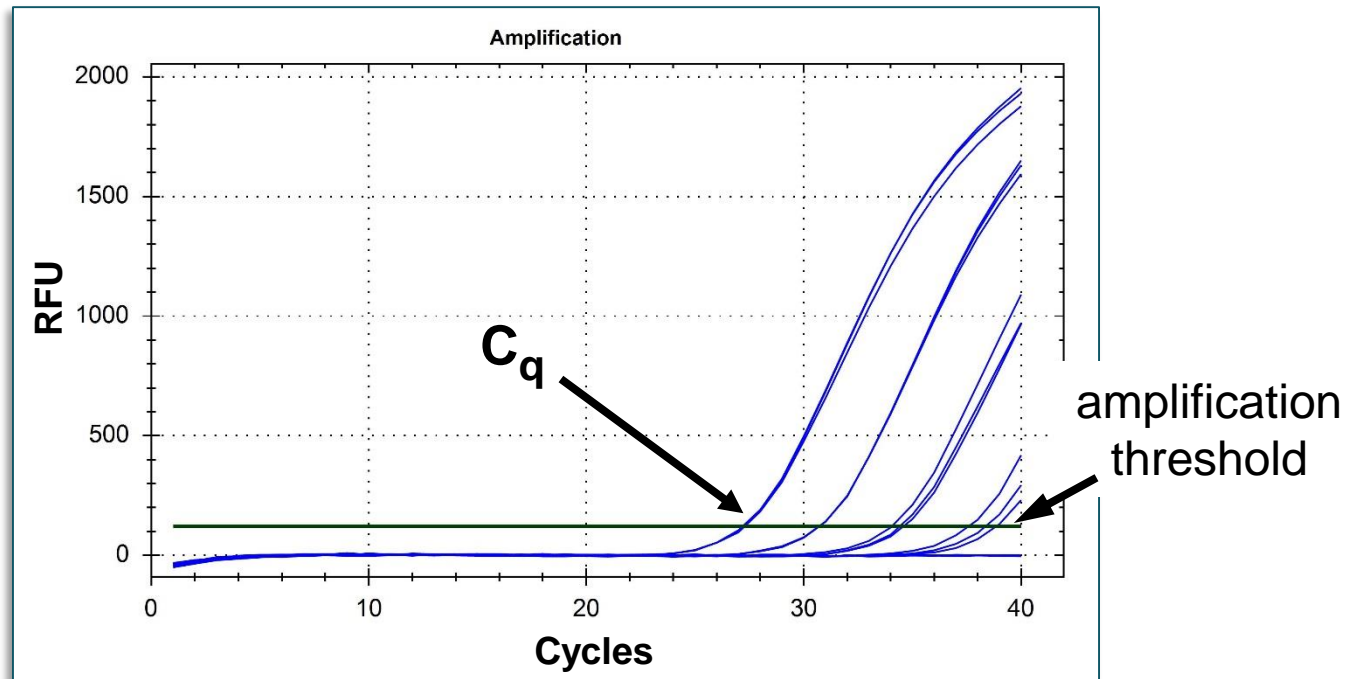
Step 3: Determine a value ( $C_q$ ) for each sample

# From Raw Data to Quantification

## *Quantification Cycle, $C_q$*

### Threshold method

- Cycle at which fluorescence crosses threshold
- Threshold set at  $\sim 10$  StDev above background
- Threshold should be drawn in exponential phase
- Threshold variable between assays; must analyze separately!

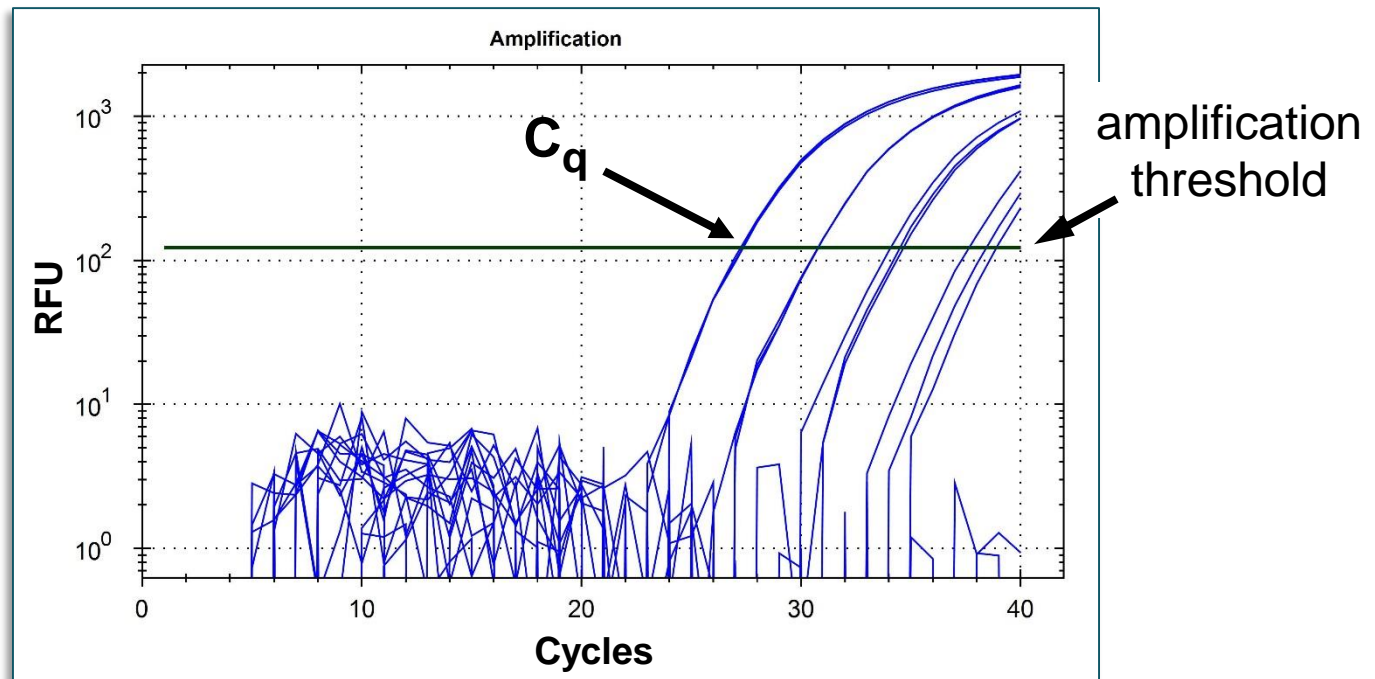


# From Raw Data to Quantification

## *Quantification Cycle, $C_q$*

### Threshold method

- Cycle at which fluorescence crosses threshold
- Threshold set at  $\sim 10$  StDev above background
- Threshold should be drawn in exponential phase
- Threshold variable between assays; must analyze separately!

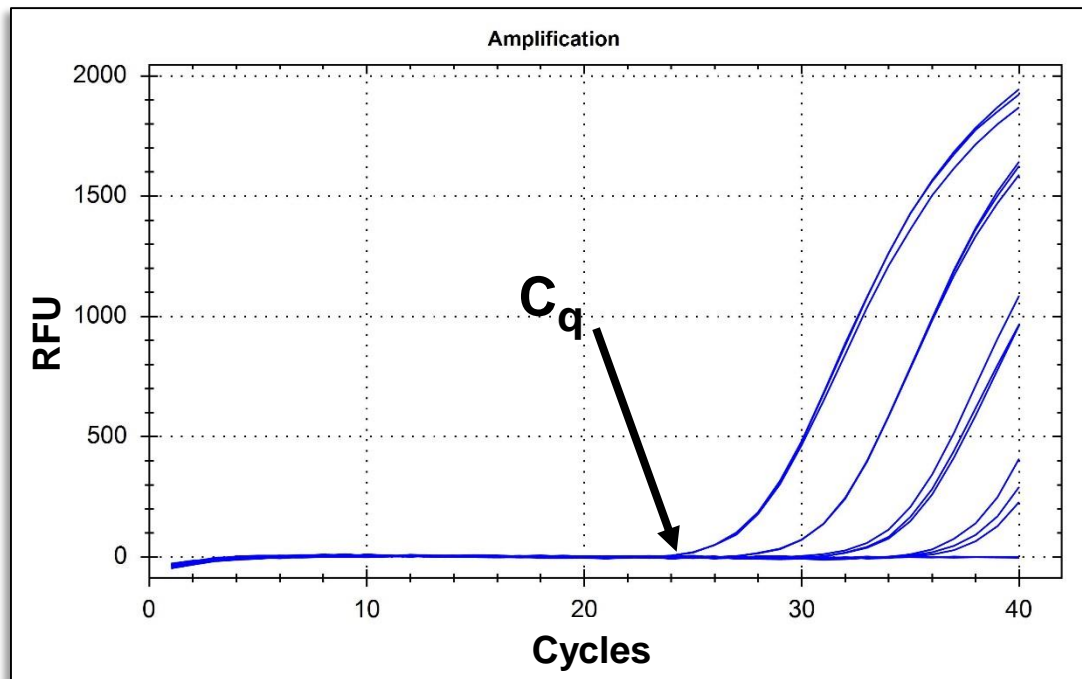


# From Raw Data to Quantification

## *Quantification Cycle, $C_q$*

### Regression method, or 2<sup>nd</sup> derivative maximum

- Point of maximum fluorescence increase
- Independent of other wells or assays on plate
- More consistent from plate to plate





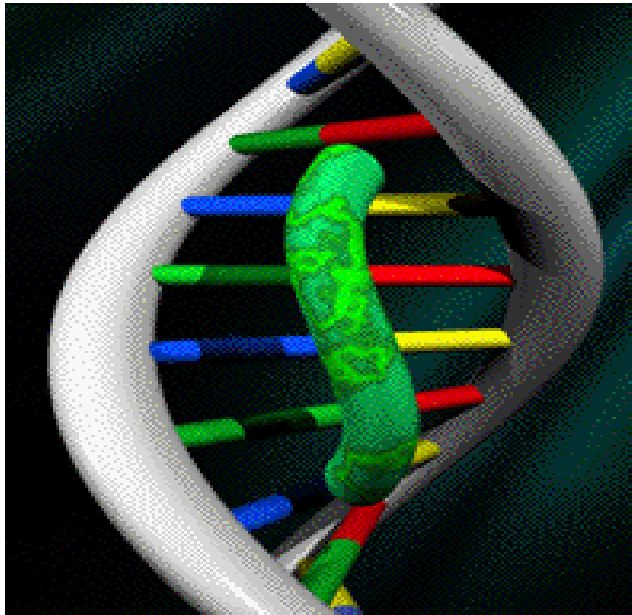
# Real-Time PCR Chemistries & Instrumentation

Promega Corporation

©2013 Promega Corporation.

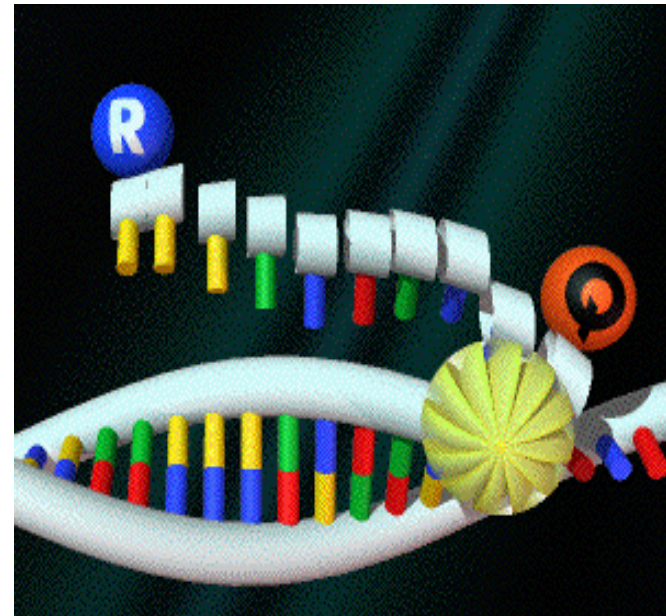
# Leading Real-Time PCR Chemistries

## Dye-based



Binds double stranded DNA

## Probe-based

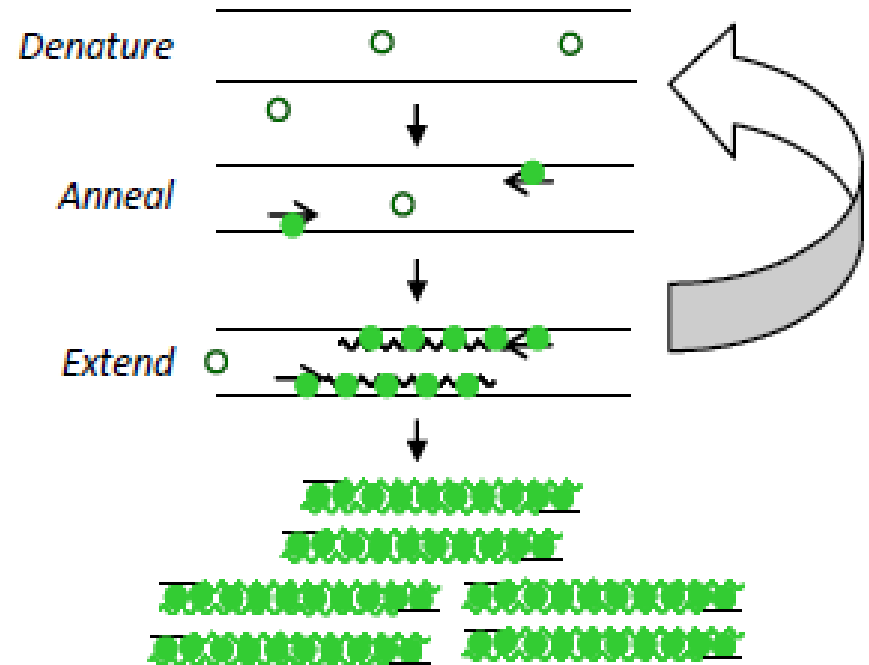


Labelled probe

# Dye-based qPCR

Uses a dsDNA-binding dye

- Low fluorescence of unbound dye
- Dye binds dsDNA stoichiometrically
- Measures all dsDNA in reaction
- Existing PCR assays can easily be adapted\*
- Less expensive
- Not multiplexed
- Allows melt analysis for QC or genotyping



## Examples:

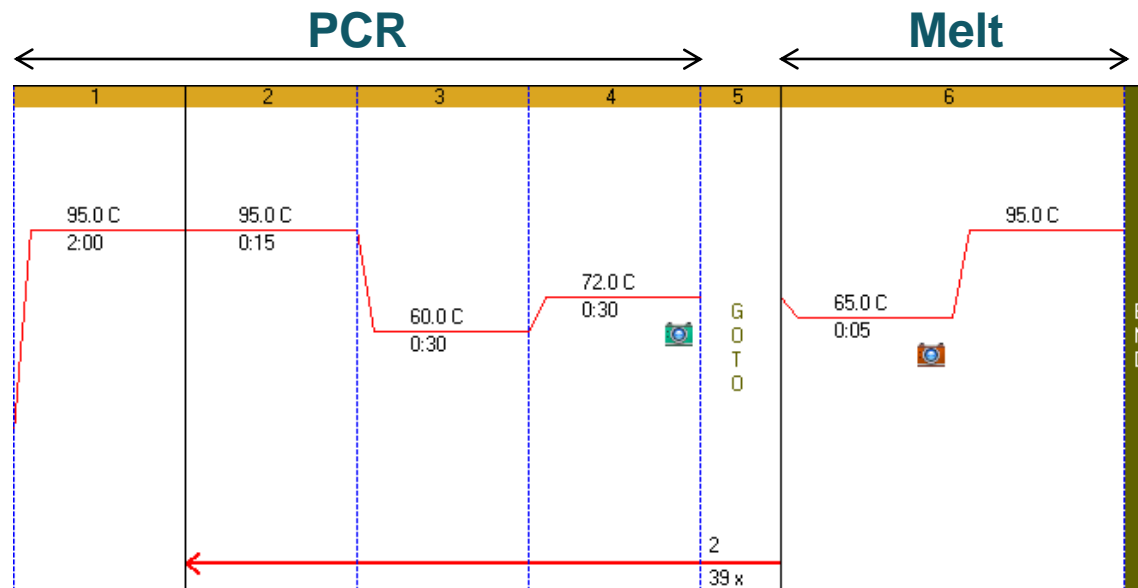
SYBR Green, BRYT Green, LC Green, Eva Green



# Melt Analysis: An internal QC

## Melt or Dissociation Curve

- After amplification, product is heated slowly, signal is continually measured
- As dsDNA amplicon denatures, dsDNA dye is displaced, signal decreases
- Melt peak  $\approx$   $T_m$  of product
- Impacted by # of amplicons, size & base composition



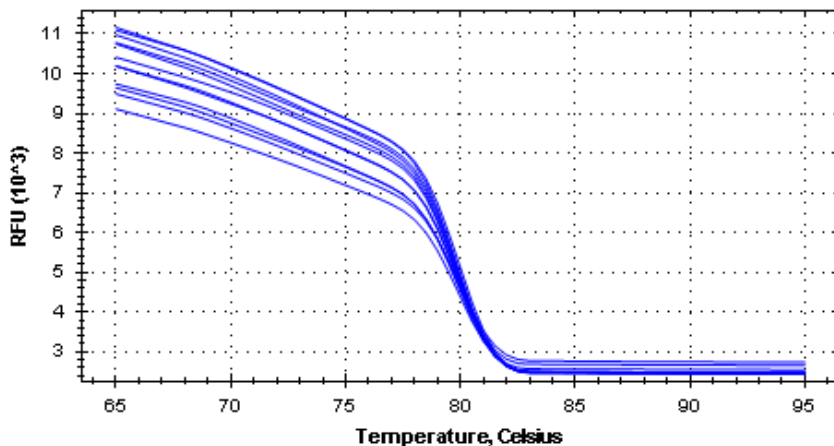
# Melt Analysis: An internal QC

## Melt or Dissociation Curve

- After amplification, product is heated slowly, signal is continually measured
- As dsDNA amplicon denatures, dsDNA dye is displaced, signal decreases
- Melt peak  $\approx$   $T_m$  of product
- Impacted by # of amplicons, size & base composition

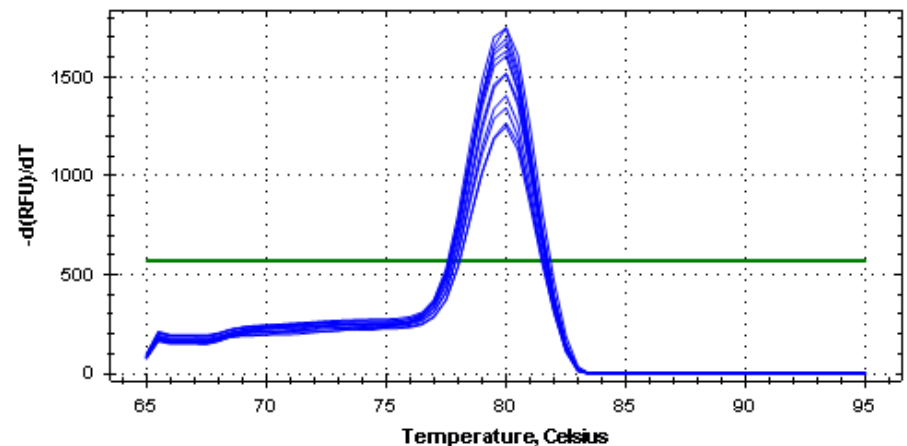
### RFU vs T(°C)

Melt Curve



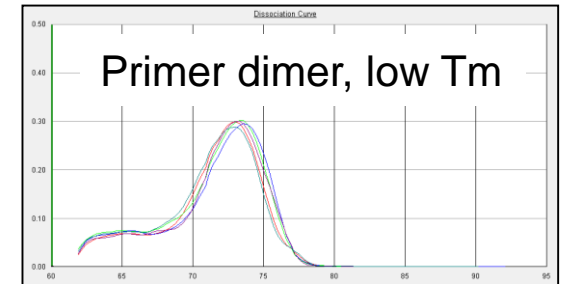
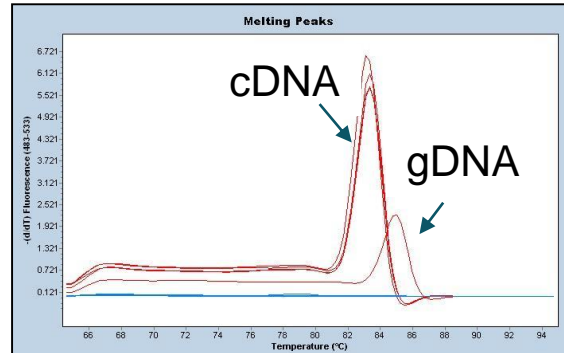
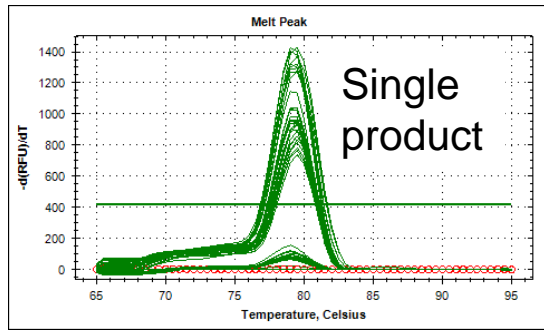
### dRFU/dT vs T(°C)

Melt Peak

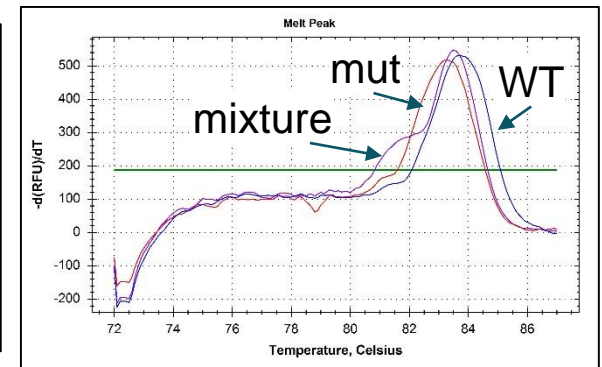
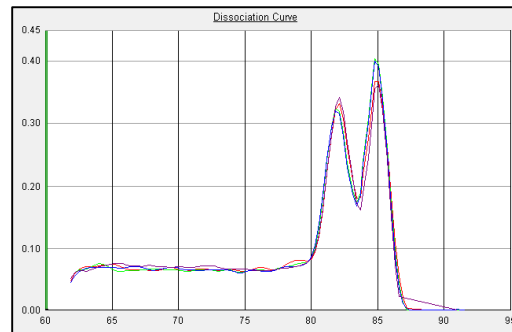
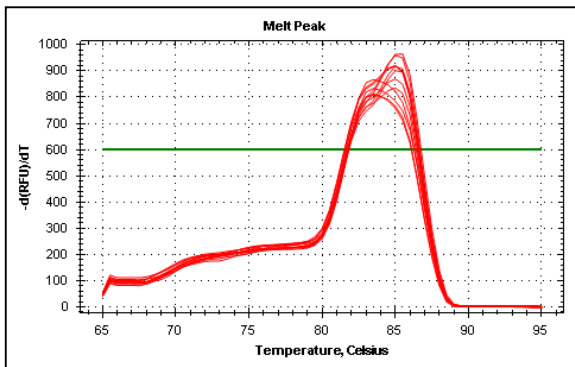


# Melt Analysis: An Internal QC

Quality control



Multiple products



High Resolution Melt (HRM)

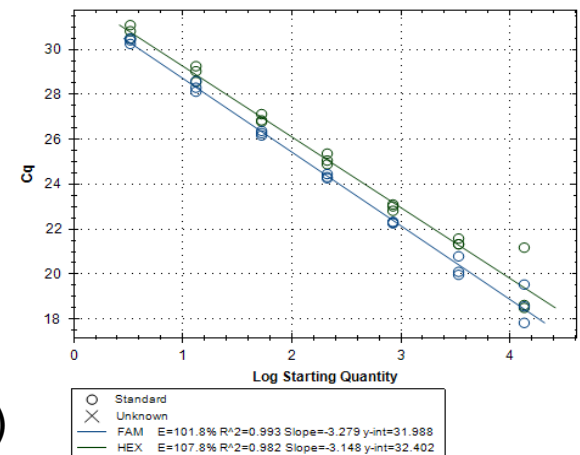
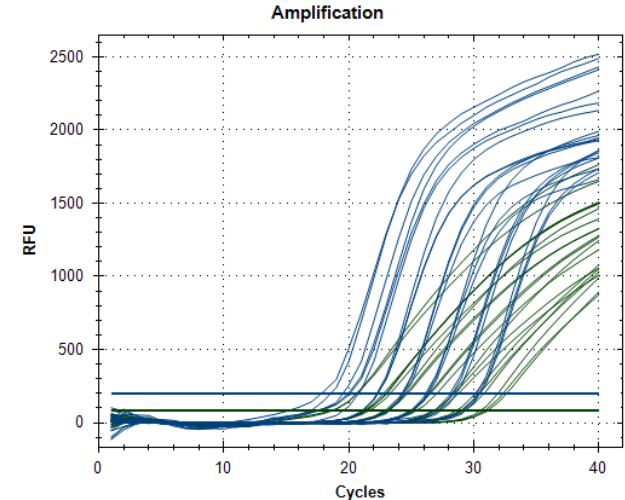
# Probe-based qPCR

Uses primers and 1-2 dye-labeled probes

- Increases specificity of target detection
- gDNA, psuedogenes still detected if amplified
- More expensive
- Allows for detection of multiple targets
- May require more optimization
- TaqMan<sup>®</sup> doesn't allow melt analysis

## Examples

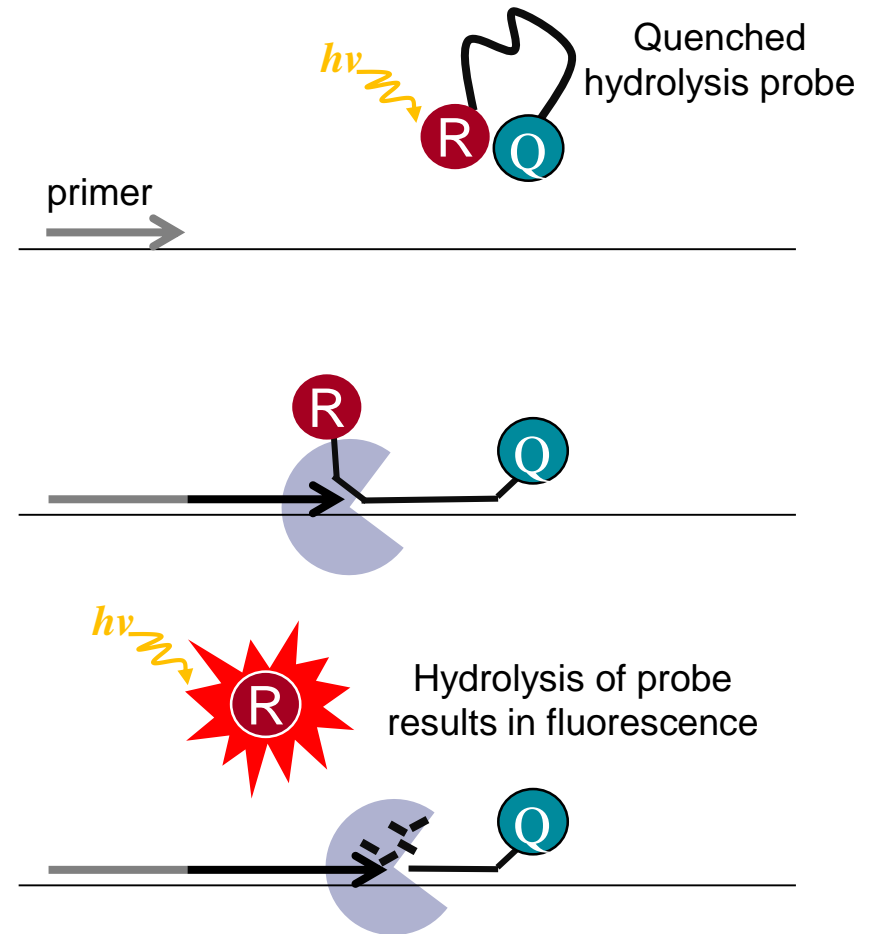
- TaqMan<sup>®</sup> (hydrolysis probe)
- Molecular Beacons<sup>®</sup> (stem-loop FRET probe)
- Dual Hybridization Probes (donor-acceptor probe)



See *Real-time PCR Detection Chemistries*, Navarro et al. (2015)

# Probe-based qPCR: *TaqMan Fluorogenic 5' Nuclease Assay*



- Probe contains reporter fluorophore and quencher
- Intact probe is quenched
- Taq degrades quenched probe during extension
- Requires 5' nuclease activity of Taq
- Irreversibly releases reporter dye from quenching



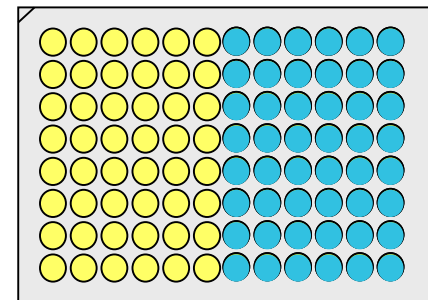
# Multiplex Analysis

*Analysis of multiple targets in the same well*

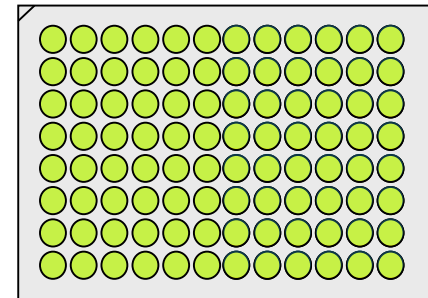
- Each target labeled with a different fluorophore
- Requires fewer wells for the same amount of data
- Requires less sample
- Better normalization
  - Same sample is assayed for all targets
- Requires more assay design & optimization
- Limited by instrumentation and dyes

Target 1   
Target 2 

Single target



Multiple targets

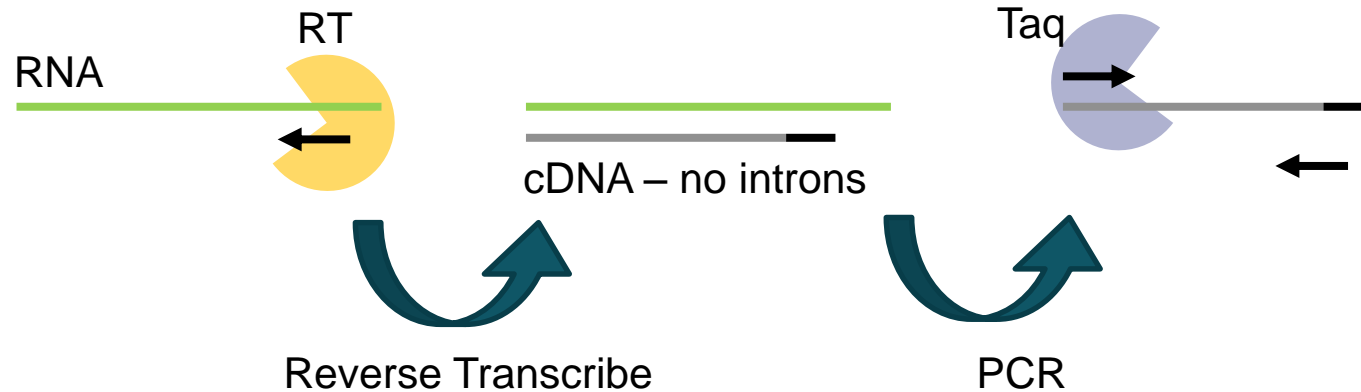


# qPCR Chemistries

## *Dye-based vs. Probe-based*

	Dye-based qPCR	Probe-based qPCR
Amplicon labeling	dsDNA-binding dye	Fluorescently labeled probes
Cost	Lower cost	Higher cost
Flexibility	All optimized assays	Single assay
Instrumentation	All qPCR instruments	Must match probes to filters
Specificity	Measures all dsDNA	Measures amplicon with probe sequence
Multiplexing	No	Yes – different dyes/filters
Melt analysis <i>QC and genotyping</i>	Yes	No (TaqMan)
Throughput	High	Highest (multiplexed)
Sample required	Low	Lowest (multiplexed)
Requires validation	Yes	Yes

# Reverse Transcription qPCR (RT-qPCR)



## Applications

- Gene expression
- Biomarker discovery
- RNA Sequencing
- RNA viruses
- cDNA cloning

## Reverse Transcriptase

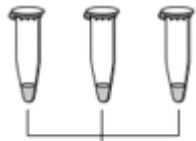
- RNA-directed DNA polymerase
- Requires priming,  $Mg^{2+}$  or  $Mn^{2+}$  cofactor
- RNase H activity in wildtype
- Inhibits Taq polymerase
- AMV, MMLV



# Reverse Transcription

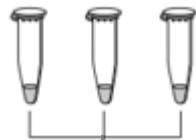
## Two-step RT-qPCR—Making a pool of cDNA

### Two Step RT-qPCR



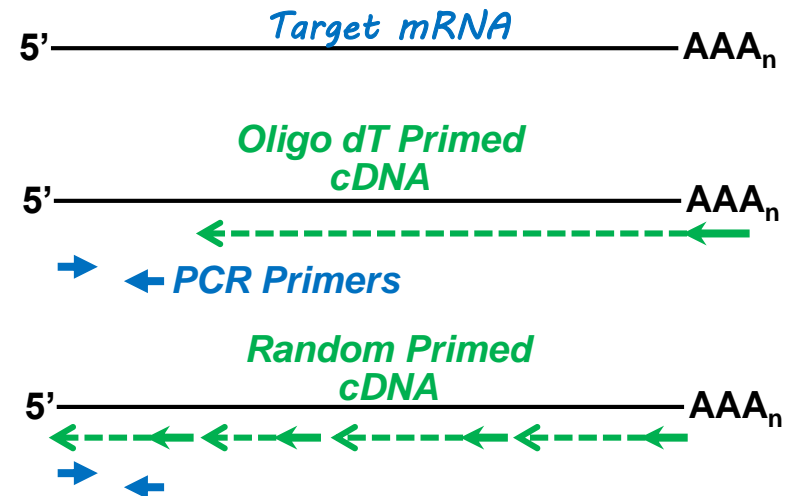
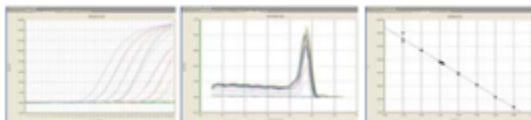
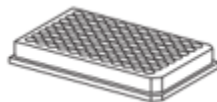
#### Reverse Transcription

Heat denature RNA w/ primers  
Add RT, Buffer, dNTPs & RNasin  
Anneal & extend  
Heat inactivate RT



#### PCR or qPCR

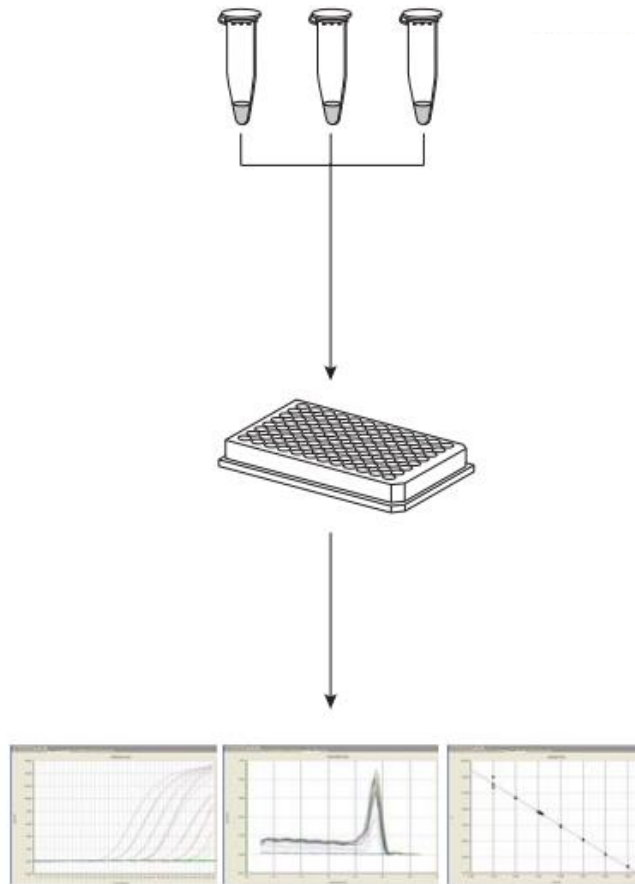
Gene-specific primers



# Reverse Transcription

## *One-step RT-qPCR—Amplification of a single target*

### *One Step RT-qPCR*



### RT & qPCR

Set up as for qPCR

Add RT and RNasin

Use gene-specific primers

### Cycling considerations

Perform RT first

Inactivate RT/ Activate Taq

Standard qPCR cycling

### Benefits

Uses less sample

Replicates over both steps

Quant & QC for FFPE RNA

# Real-Time Instruments

*Examples:*



Bio-Rad  
CFX96 Touch



ABI  
7500 Fast



Roche  
LightCycler 480



Stratagene  
Mx3005p

## Hardware Differences

- Excitation source (lamp, LED, LASER)
- Detection method (CCD camera, PMT)
- Filters (Filter wheel, # of filters)
- Specialized (Gradient block, # of wells, interchangeable blocks, fast cycling)

## Software Differences

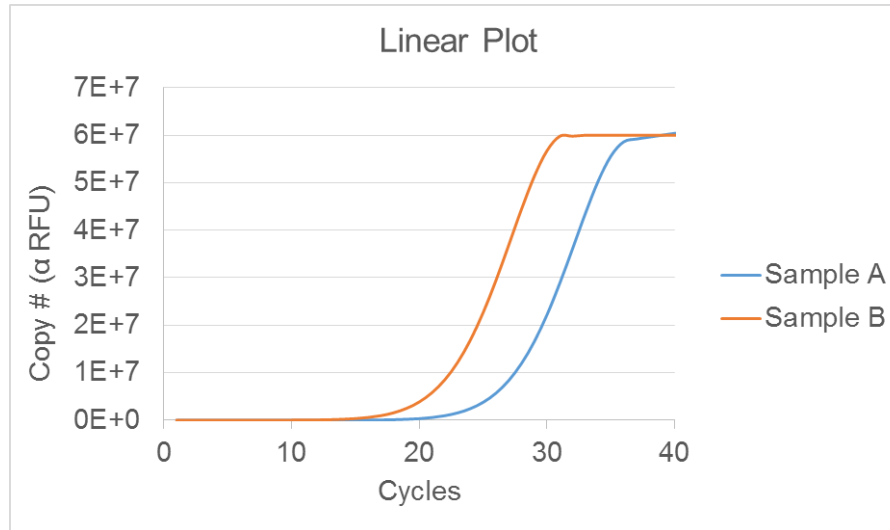
- Analysis methods
- Analysis flexibility
- Ease of use
- Specialized (HRM, bar-coding, etc.)
- Traceability, *in vitro* Diagnostic use



# Introduction to Quantification

# Comparing Samples

*Assuming identical input amounts*



<u>Sample</u>	<u>Copy#</u>	<u>Cycle1</u>	<u>Cycle2</u>	<u>Cycle3</u>	<u>Cycle4</u>	<u>Cycle5</u>	<u>Cycle6</u>	<u>Cycle n</u>
A	1	2	4	8	16	32	64	$1 \cdot 2^n$
B	10	20	40	80	160	320	640	$10 \cdot 2^n$

Quantity at a given cycle “n” is expressed as  $a * 2^n$   
 where  $a$  is starting quantity

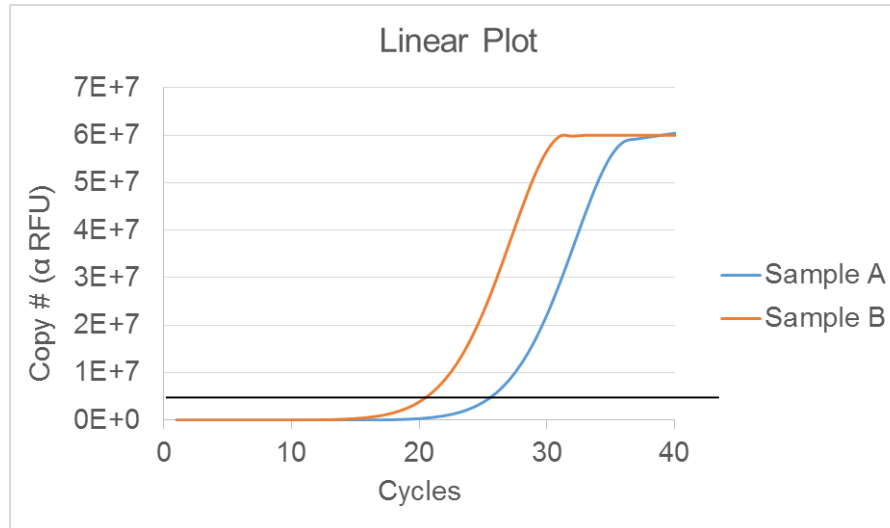
# Comparing Samples

*Assuming identical input amounts*

$\Delta Cq$  method

$$\frac{b}{a} = 2^{\Delta Cq}$$

Not recommended!



Sample A will reach the threshold  $T = a * 2^x$  at cycle  $x$   
Sample B will reach the threshold  $T = b * 2^y$  at cycle  $y$

Since the threshold  $T$  is the same for both,

$$a * 2^x = b * 2^y$$

If we rearrange to determine fold difference in starting quantities,

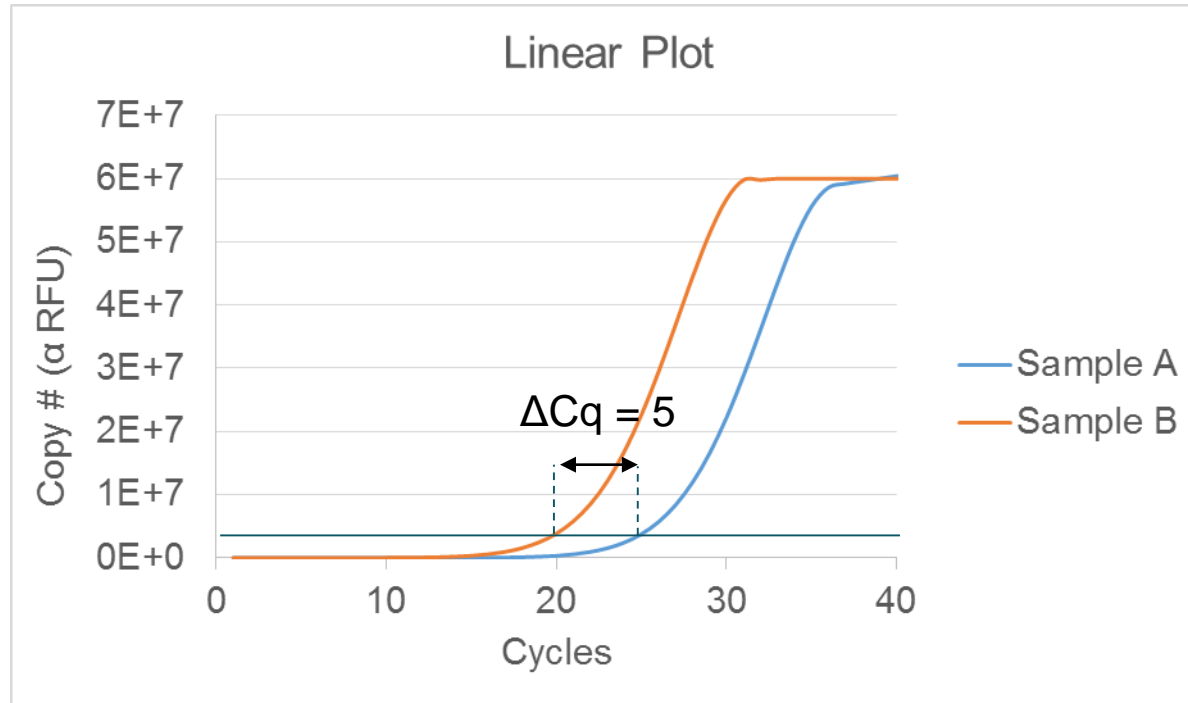
$$\frac{b}{a} = \frac{2^x}{2^y} = \frac{2*2*2*2*\dots*x \text{ times}}{2*2*2*2*\dots*y \text{ times}} = 2^{x-y} = 2^{\Delta Cq}$$

# Common Assumptions in Calculations

---

- $\Delta Cq$  method
  - **ASSUMPTION 1**: Amplification efficiency is perfect

# Common Assumptions in Calculations



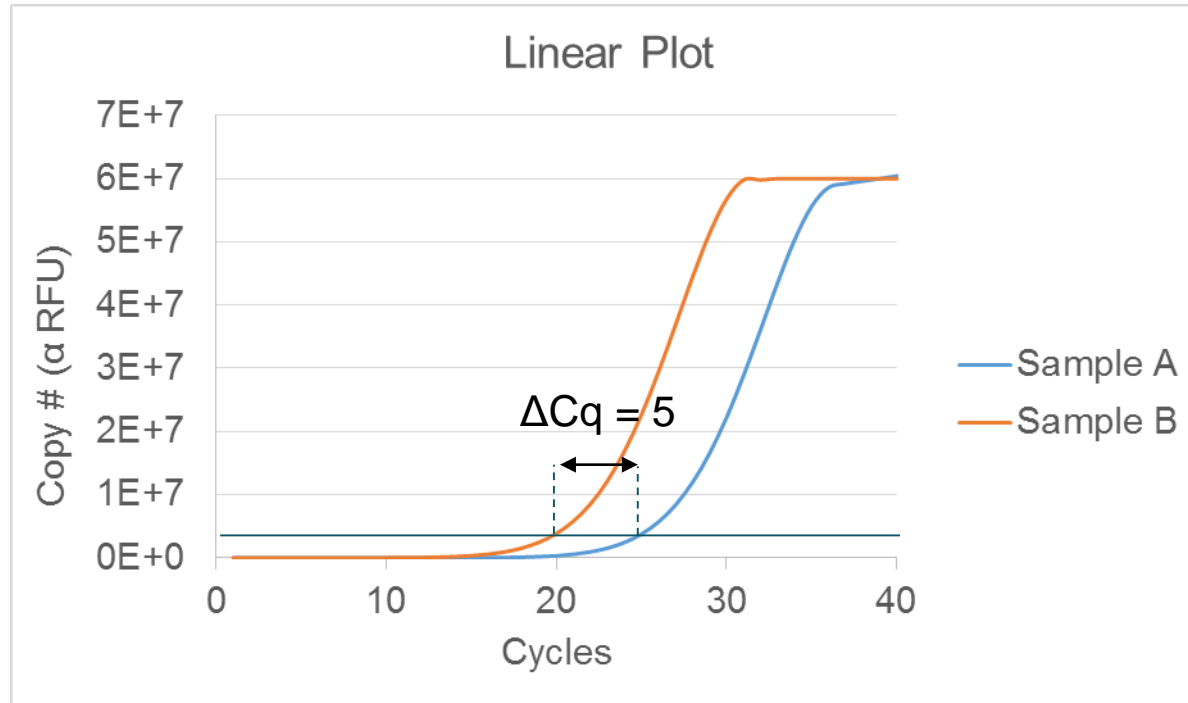
**QUESTION 2:** What is the difference in target concentration between samples A & B if the assay has 100% amplification efficiency?

$$\frac{b}{a} = 2^{\Delta Cq} = 2^5$$

**32-fold different**



# Common Assumptions in Calculations



**QUESTION 3:** What is the difference in target concentration between samples A & B if the assay has 80% amplification efficiency?

$$\frac{b}{a} = 1.8^{\Delta Cq} = 1.8^5$$

**18.9-fold different**

# Common Assumptions in Calculations

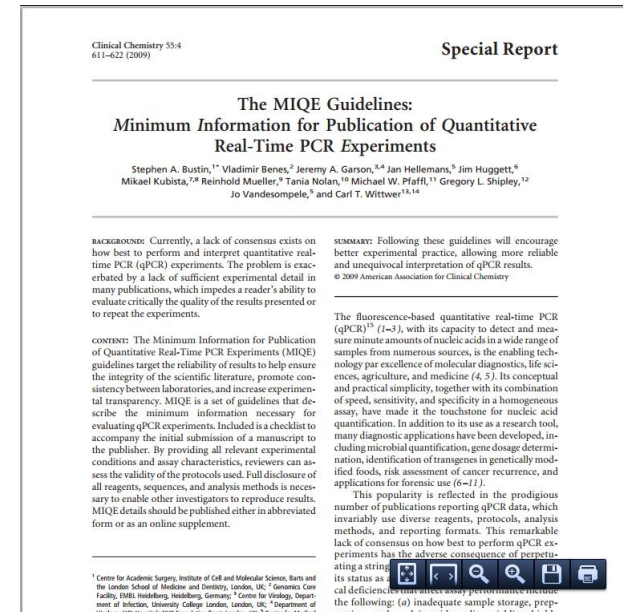
- $\Delta Cq$  method
  - **ASSUMPTION 1:** Amplification efficiency is perfect
    - Instead of  $2^{\Delta Cq}$ , measure the assay efficiency experimentally and use it!
    - Use your samples (or very similar ones) to determine efficiency
  - **ASSUMPTION 2:** Identical inputs were used from each sample
    - Quantification accuracy really matters!
    - Normalize to a reference gene that is expressed at a constant concentration
- $\Delta\Delta Cq$  method
  - Normalize  $\frac{b}{a}$  between GOI and reference gene in the same sample
  - **ASSUMPTION 3:** Reference gene expression is constant
    - Use multiple reference genes and verify expression is constant
    - See *Hellemans & Vandesompele (2014)*!

# The MIQE Guidelines—Read *BEFORE* You Plan!

## The MIQE Guidelines Minimum Information for Publication of Quantitative Real-Time PCR Experiments

Bustin, et al., *Clinical Chemistry* 55:4, 611–622 (2009)

- Goal is to improve repeatability between labs
- Outlines common vocabulary to use
- Includes a checklist for reporting & planning
- Describes minimal reporting requirements
  1. Description of sample manipulation
    - storage, extraction, quantification, integrity, etc.
  2. Description of assay reagents, protocols, and controls
    - standard curves,  $\geq 2$  reference genes, -RT controls, NTCs
  3. Assay validation—Required even for purchased assays!
    - sensitivity, specificity, and efficiency; reference gene stability
  4. Description of calculations



# Real-Time PCR Resources On-line

---

## General Real-Time PCR:

MIQE Guidelines:	Bustin <i>et al.</i> , <i>Clin Chem</i> (2009) <a href="http://miqe-press.gene-quantification.info/">http://miqe-press.gene-quantification.info/</a>
Gene Quantification:	<a href="http://www.gene-quantification.info">www.gene-quantification.info</a>
Real-time chemistries:	Navarro <i>et al.</i> , <i>Clinica Chimica Acta</i> (2015)
Reference genes:	Hellemans and Vandesompele, <i>Methods Mol Biol</i> (2014) <a href="http://www.gene-quantification.com/Bio-Rad_2008_Rethink_PCR_Conference_Hellemans_Vandesompele.pdf">http://www.gene-quantification.com/Bio-Rad_2008_Rethink_PCR_Conference_Hellemans_Vandesompele.pdf</a> geNorm and qBase+ software

## Primer design software & Pre-designed assays:

Primer3 -	<a href="http://www-genome.wi.mit.edu/cgi-bin/primer/primer3_www.cgi">http://www-genome.wi.mit.edu/cgi-bin/primer/primer3_www.cgi</a>
Primer-BLAST -	<a href="http://www.ncbi.nlm.nih.gov/tools/primer-blast/">http://www.ncbi.nlm.nih.gov/tools/primer-blast/</a>
IDT PrimerQuest -	<a href="http://biotools.idtdna.com/Primerquest/">http://biotools.idtdna.com/Primerquest/</a>
Primer Bank -	<a href="http://pga.mgh.harvard.edu/primerbank/">http://pga.mgh.harvard.edu/primerbank/</a>
RTPrimerDB -	<a href="http://medgen.ugent.be/rtpprimerdb">http://medgen.ugent.be/rtpprimerdb</a>

# Technical Services Scientists Ready to Help

## Customer & Technical Support



### Customer Support

Contact customer service to check price, availability, place an order, check delivery or check on an order status, or get a copy of your invoice. We are here to help you!

[Chat with Customer Service](#)

**Phone:** (608) 274-4330

**Toll-Free Phone:** (800) 356-9526

**Fax:** (608) 277-2516

**Toll-Free Fax:** (800) 356-1970

**Email Address:** [custserv@promega.com](mailto:custserv@promega.com)

**Hours:** 7am - 6pm CST, Monday-Friday

[Contact Us Form](#)

### Technical Support

To answer technical questions about our products, request technical seminars and for information about our new products, our team of scientists are here to support you!

[Chat with a Scientist](#)

**Phone:** (608) 274-4330

**Toll-Free Phone:** (800) 356-9526

**Fax:** (608) 277-2516

**Toll-Free Fax:** (800) 356-1970

**Email Address:** [techserv@promega.com](mailto:techserv@promega.com)

**Hours:** 7am - 6pm CST, Monday-Friday

[Technical Request Form](#)



# Questions Welcome