

Maximize Your Reverse Transcription-qPCR (RT-qPCR) Assays

Carl A. Strayer, PhD February 2012



*If you see this icon in the upper left hand corner of a slide, **DOUBLE CLICK IT** to view the speaker notes.*

Outline



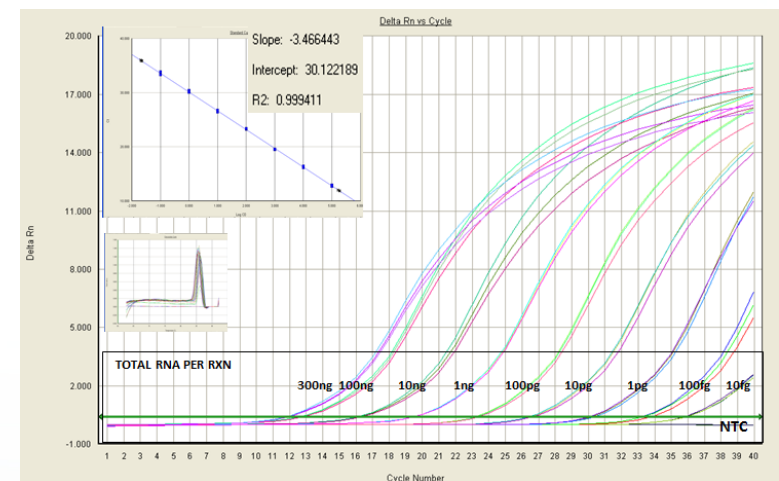
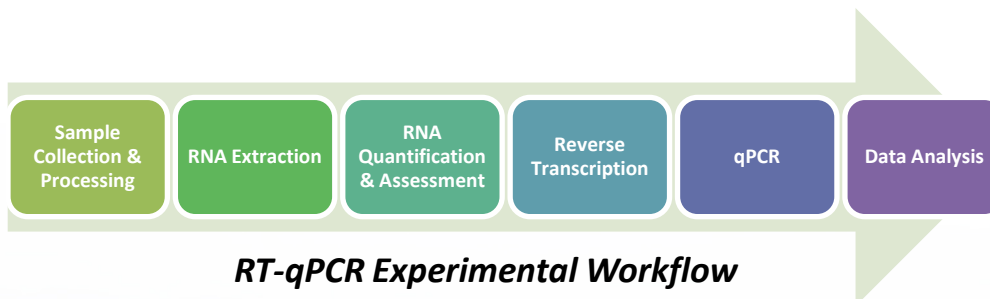
I. Fundamentals of Real-Time PCR

- What is Real-Time PCR?
- Real-Time PCR Metrics & Analysis

II. RT-qPCR Assay Design

- Primer Design
- Assay Validation

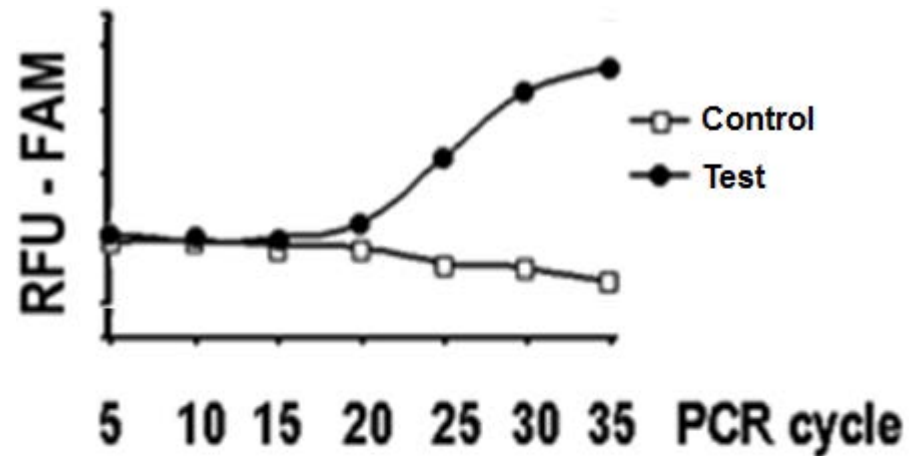
III. Tips & Troubleshooting



What is Real-Time PCR?

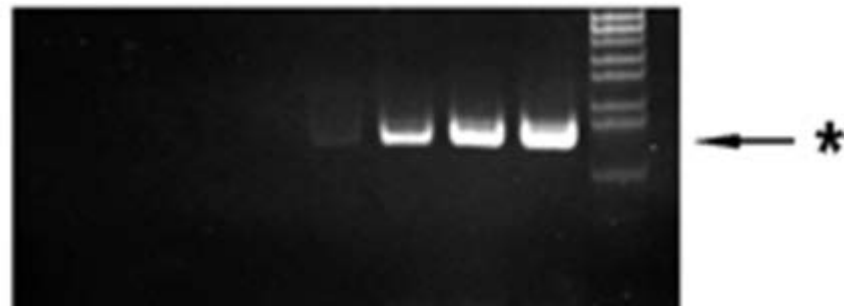
Real-Time PCR

Product formation
measured
at each cycle,
during the reaction



Endpoint PCR

Product formation
measured
after reaction is
complete



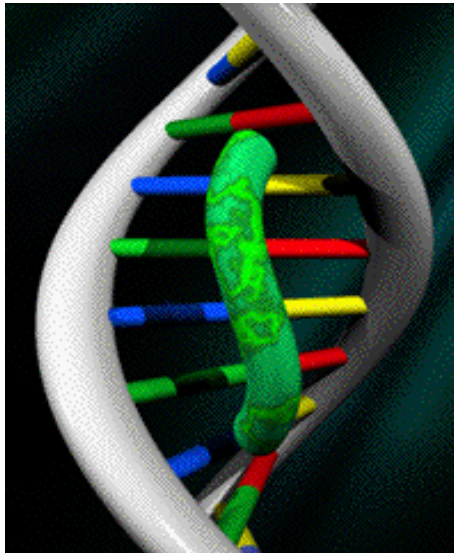
Adapted from Sherrill et al., JACS 2004

Real-Time PCR Chemistries

A fluorescent **Reporter** is used to detect product formation

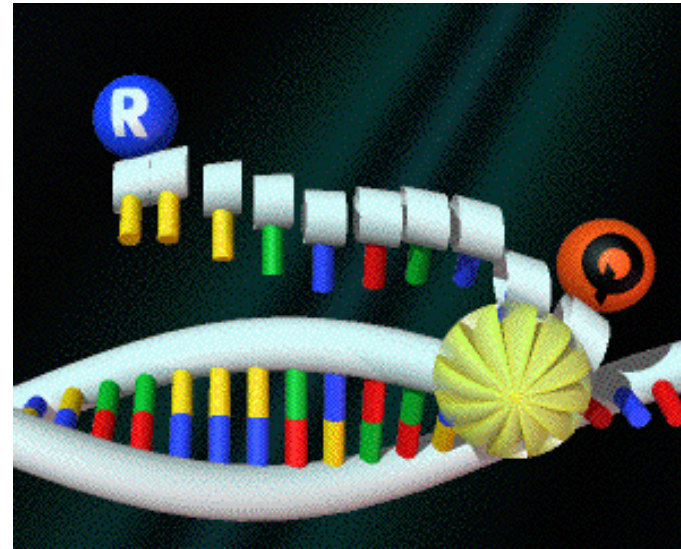
- part of the reaction mix
- **two general types...**

dsDNA binding dye



or

Labeled primer or probe



Real Time PCR Instruments



Thermal cycler + fluorescence detection module

Many manufacturers, many models, for example...



ABI 7500
Real Time System



Bio-Rad MyiQ2



Stratagene Mx3005P®

Hardware differences - determine reporter compatibility, multiplexing capability, cost

Excitation source

Lamp, laser, LEDs

Detection method

CCD Camera or PMT

Fluorescence Filters

Type & number

Real-Time PCR Chemistries – Dye-based

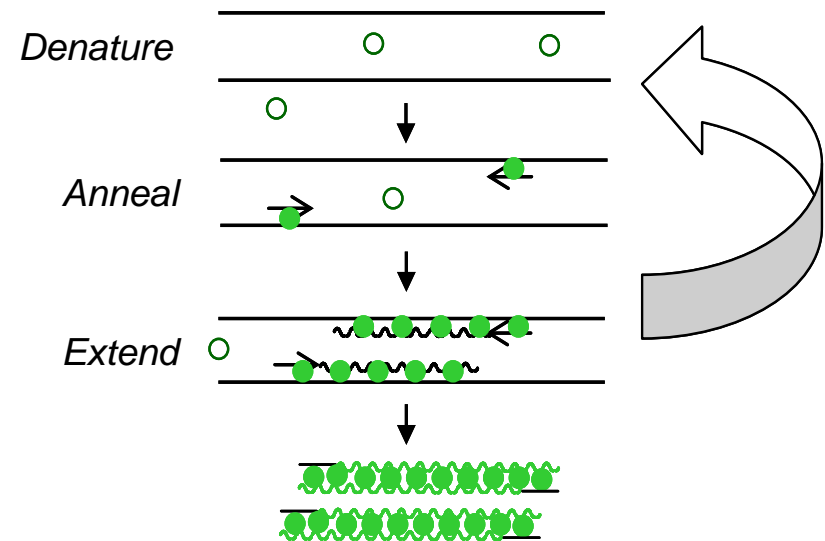


- *dsDNA-binding dye is included in PCR mastermix*
- *Standard primers used*
- *Dye associates with PCR product*

- *Free Dye -> low fluorescence*
- *Bound Dye -> high fluorescence*

As more PCR product is produced,
more dye is bound

*Fluorescence is proportional to the
amount of product*



*SYBR® Green is familiar...
...improved dyes now available...*

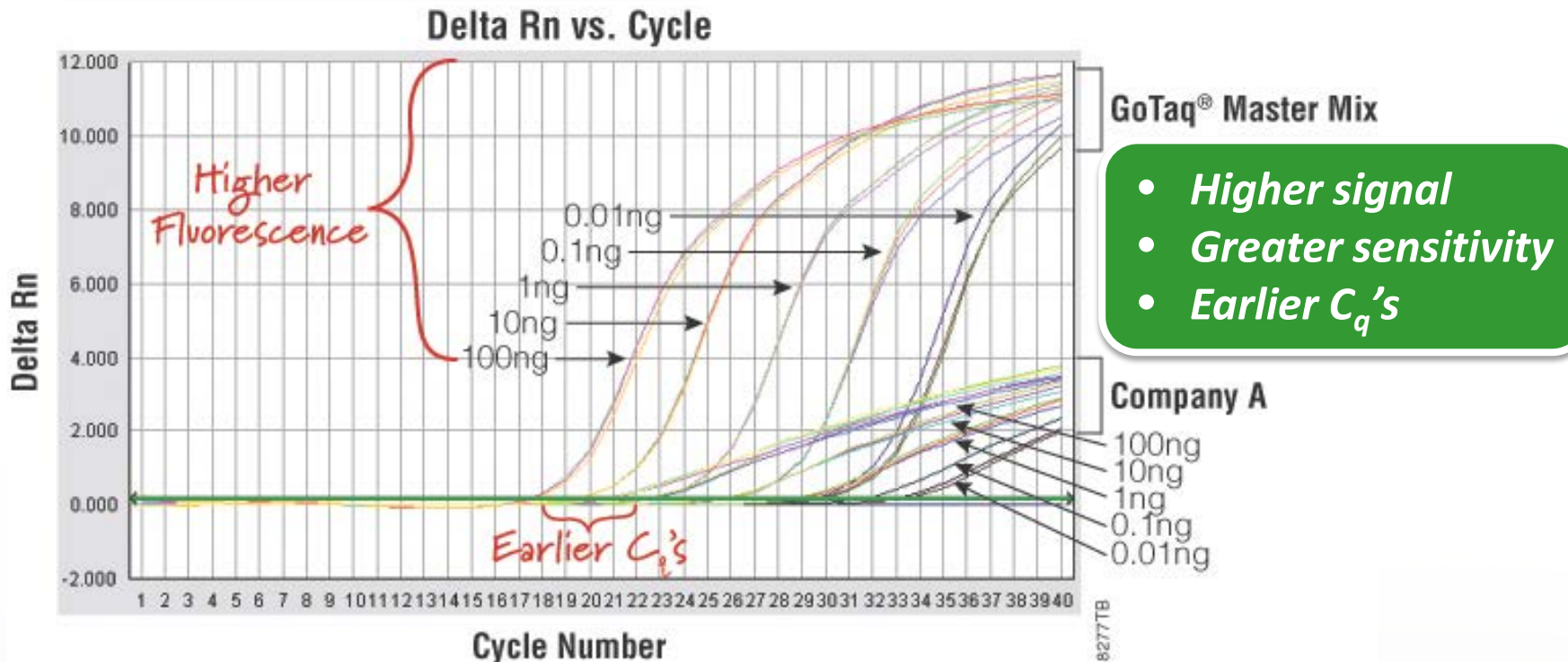
New generation of Dye-based Real-Time systems offer performance advantages



GoTaq[®] qPCR Master Mix

- BRYT Green[®] Dye
- Hot-start Taq
- Optimized reaction buffer

- **BRYT Green[®]** is a new dsDNA binding dye developed by Promega
- Spectra nearly identical to SYBR[®] Green I, detected using same filters



Real-Time PCR Chemistries – Label-based

- Primer or probe is synthesized with reporter
- Product formation alters fluorescence of the reporter

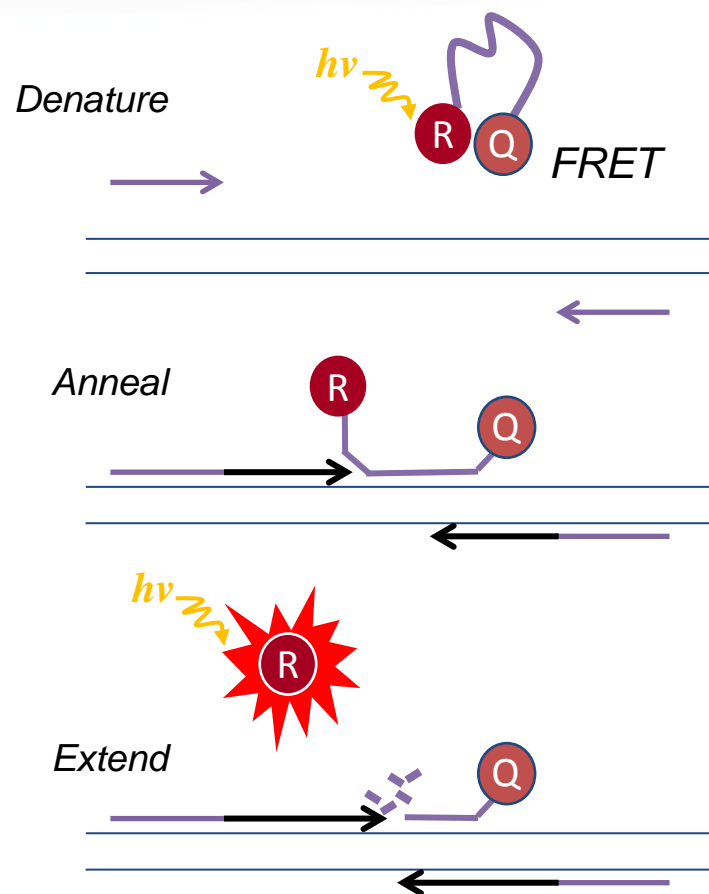
TaqMan[®] is the most familiar type:

- 2 PCR primers + 1 probe
- Probe labeled with reporter & quencher
- Primers & probe anneal to target
- During extension, 5' nuclease activity of Taq degrades probe

Free probe -> FRET occurs

Degraded probe -> reporter un-quenched

Fluorescence is proportional to the amount of product



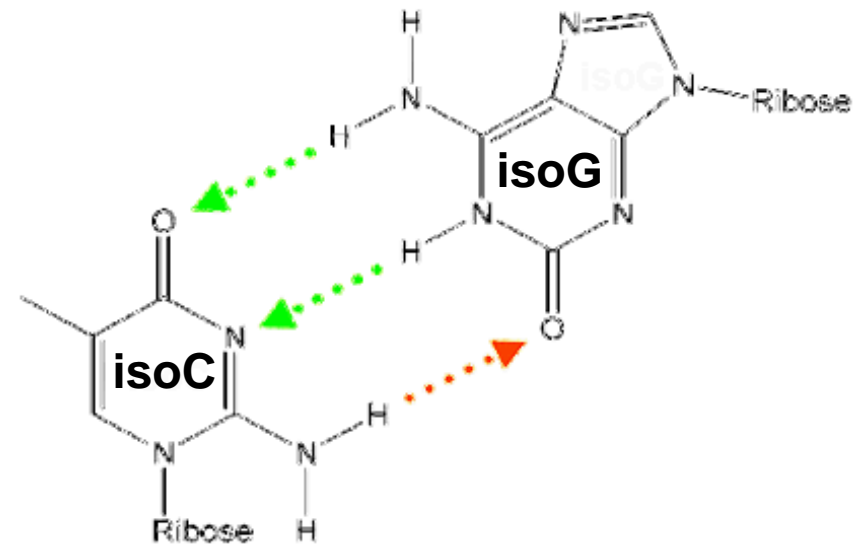
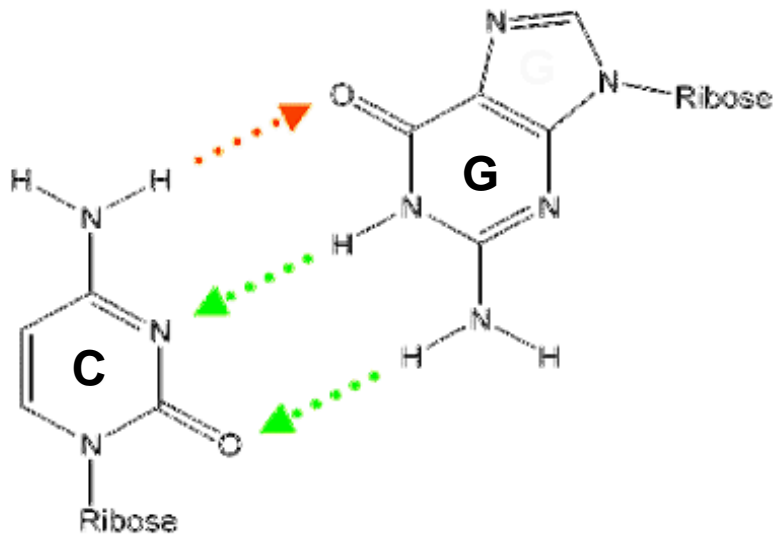
As with dye-based chemistry, other options available...

Plexor[®] Technology - Novel Label-based Chemistry



Iso-C & Iso-G dNTPs

- Pair only with each other - not with A·C·G·T·U
- Recognized by DNA Polymerase



Johnson, S.C., et al. (2004) *Nucleic Acids Res.* **32**, 1937-41.

Plexor[®] Technology - Novel Label-based Chemistry

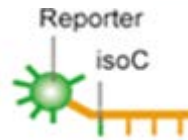


2 Primer Method

- 1 standard primer
- 1 primer with iso-C base & reporter at 5' end

Amplification Master Mix Contains

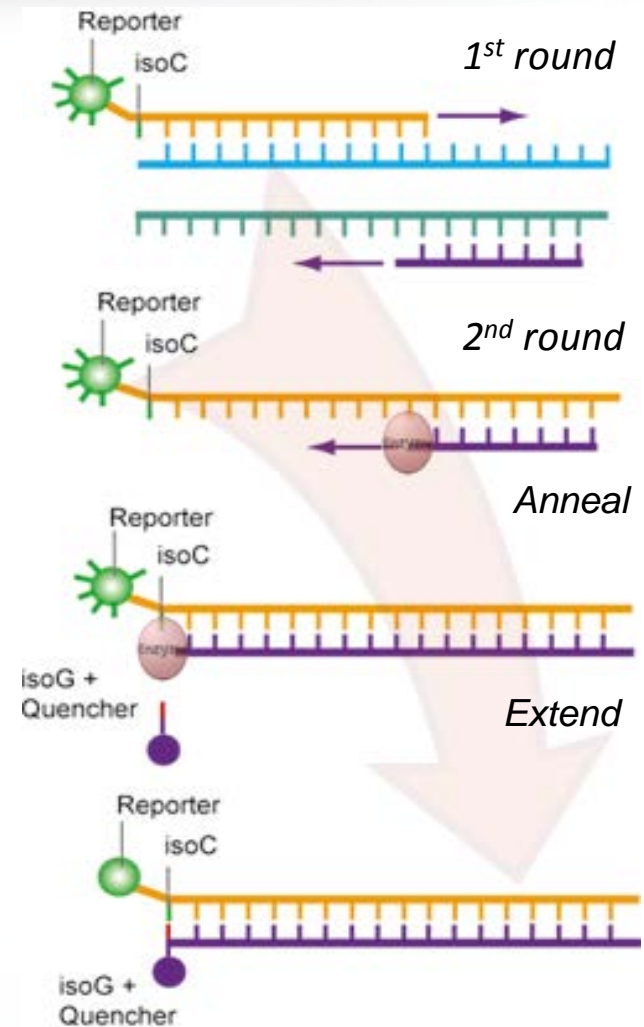
- Standard dNTPs
- Quencher-labeled iso-dGTP



Free primer-reporter & quencher -> no FRET

Incorporation of primer & quencher -> FRET occurs

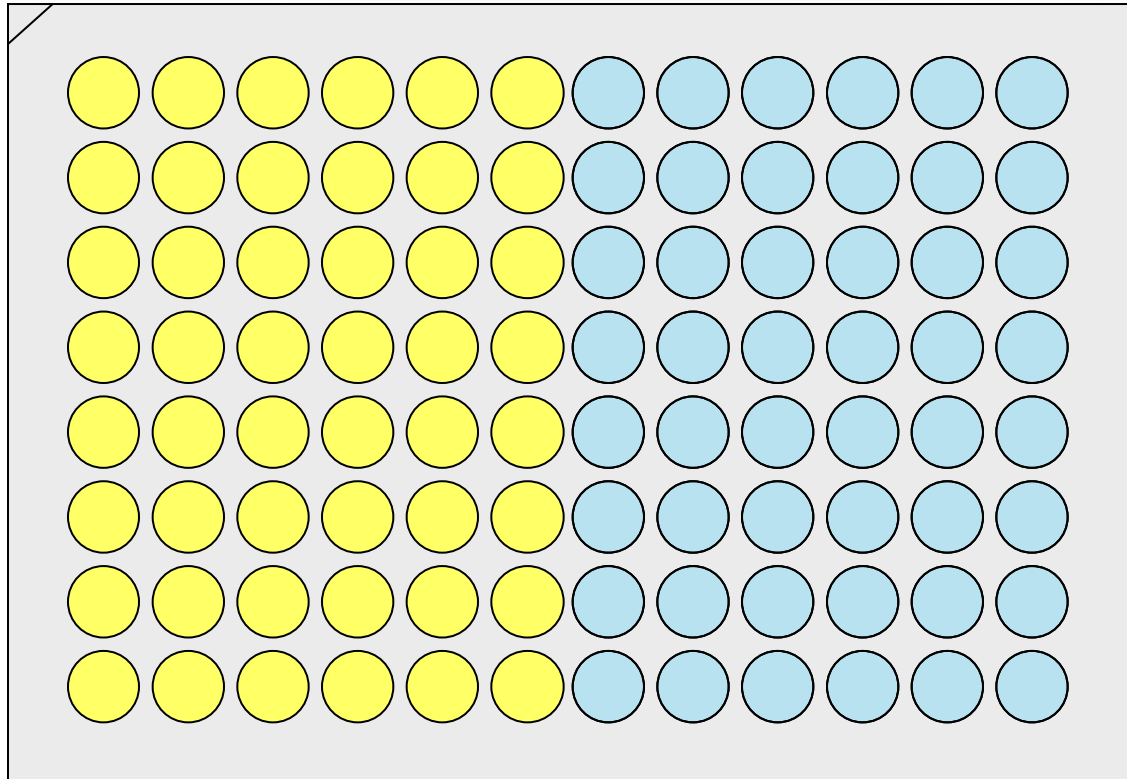
Fluorescence is inversely proportional to the amount of product



Label-based methods allow multiplexing

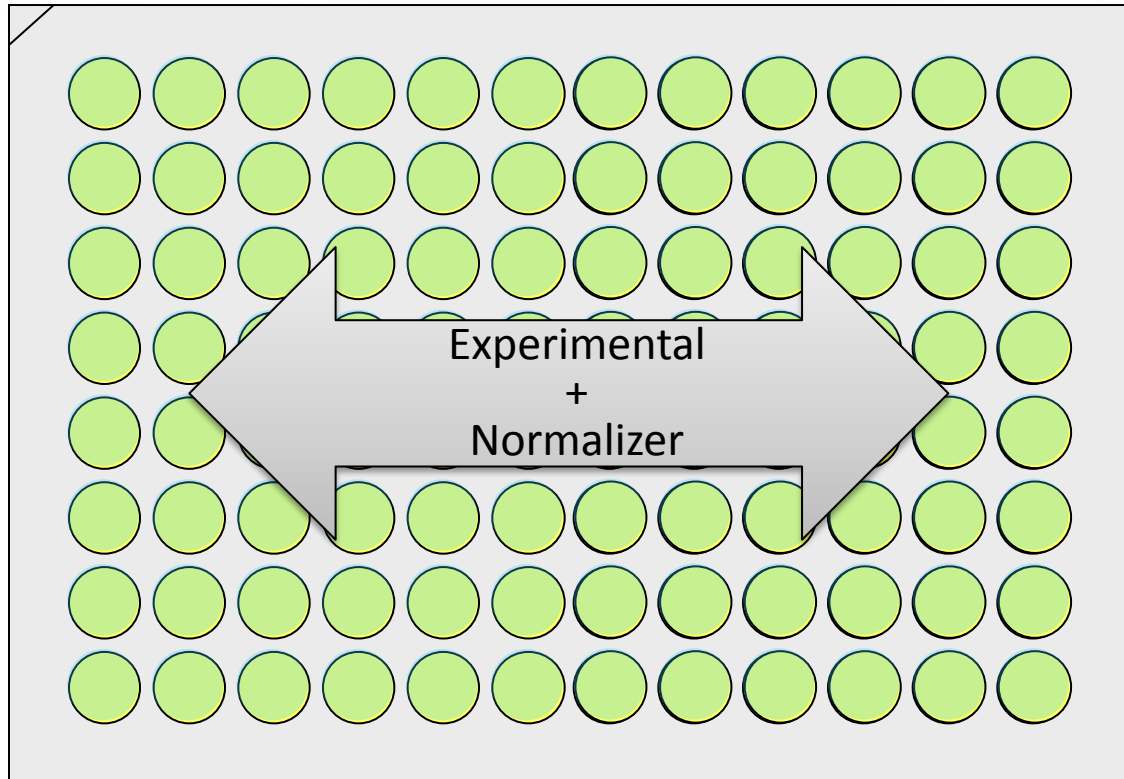


Experimental
target



Normalizer

Label-based methods allow multiplexing



- ***Better data*** - Targets & Normalizer in same reaction
- ***Economical*** - More samples analyzed per well/plate
- ***Conserves sample***

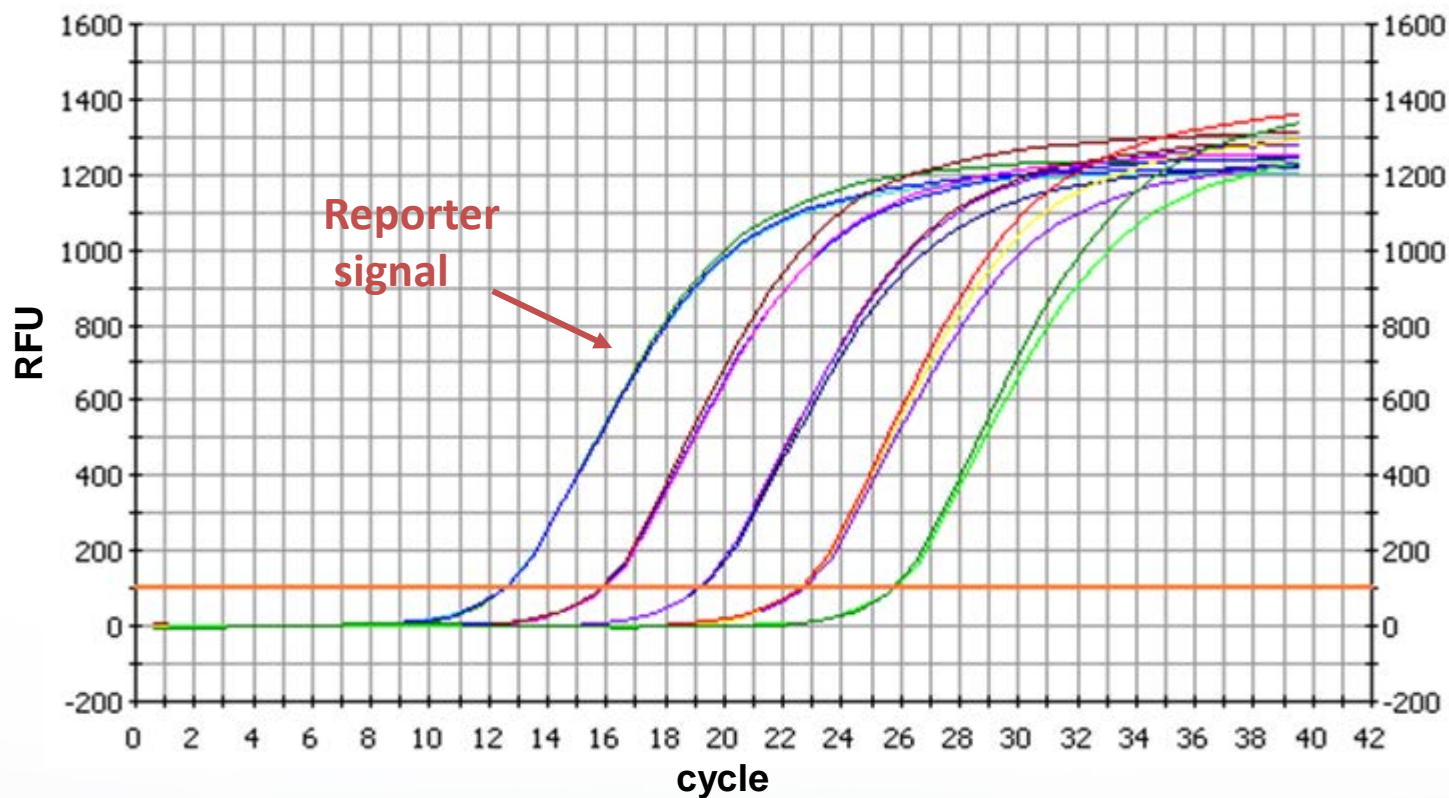
Real-Time PCR Metrics & Analysis



Primary output is the Amplification Curve

Amplification Curve – shows accumulation of product as PCR progresses

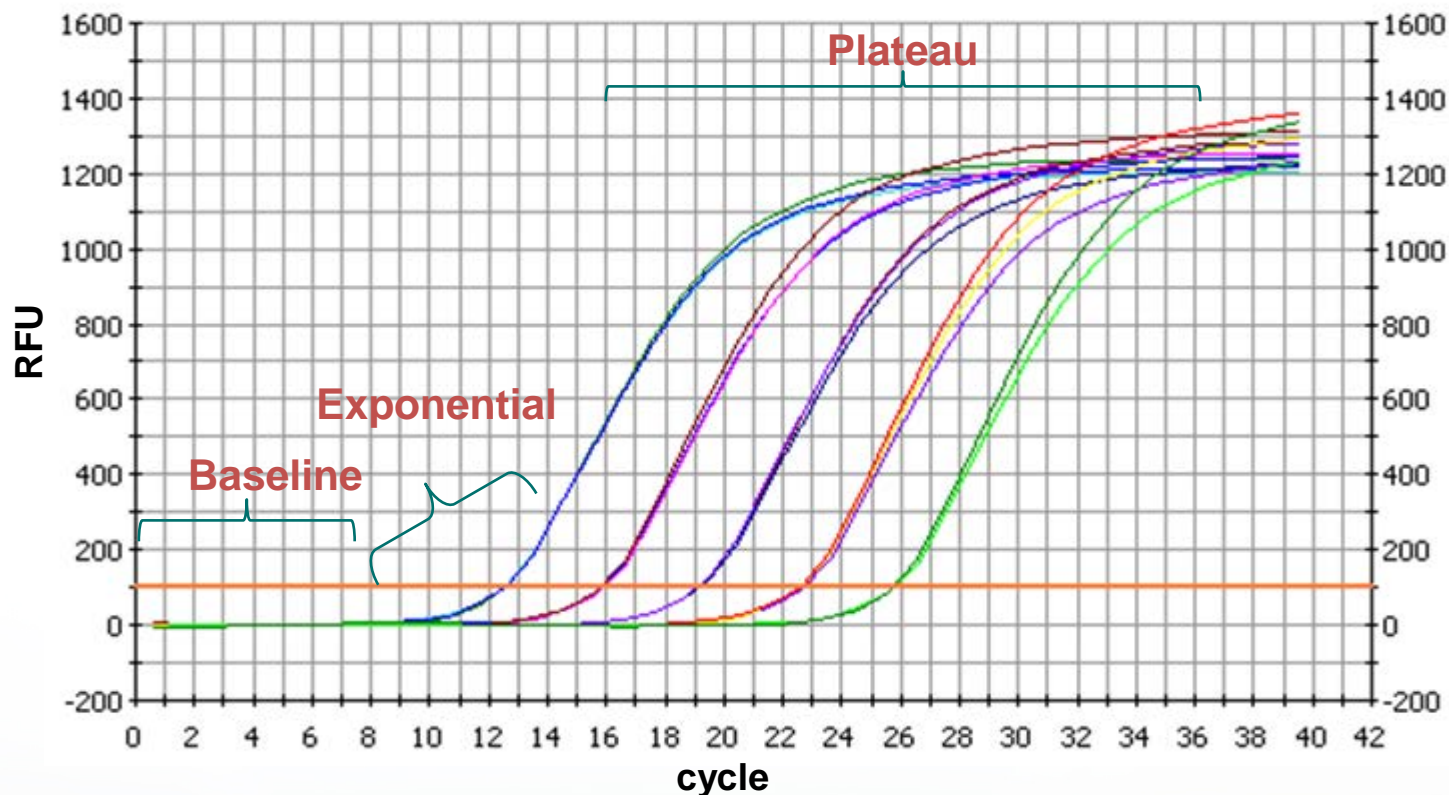
- **reporter** – fluorescent dye or label used to monitor PCR product formation
- **R** – raw fluorescence of reporter (RFU = relative fluorescence unit)



Primary output is the Amplification Curve

Amplification Curve – shows accumulation of product as PCR progresses

- **baseline** – initial reporter fluorescence, before significant product formation occurs
- **exponential phase** – stage of reaction when product is doubling with each cycle
- **plateau phase** – stage of reaction when rate of product formation is diminishing

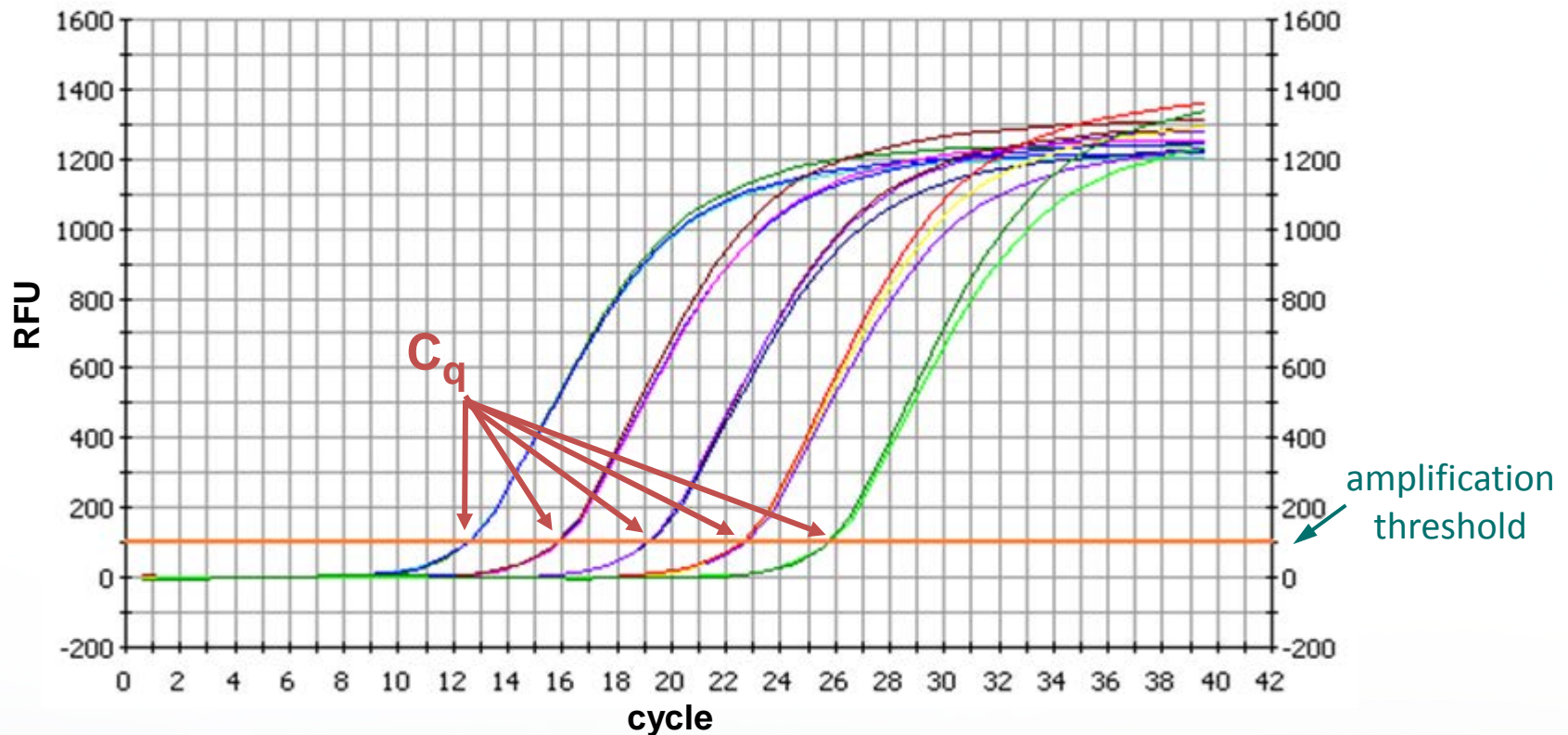


Analysis of the amplification curves gives C_q value



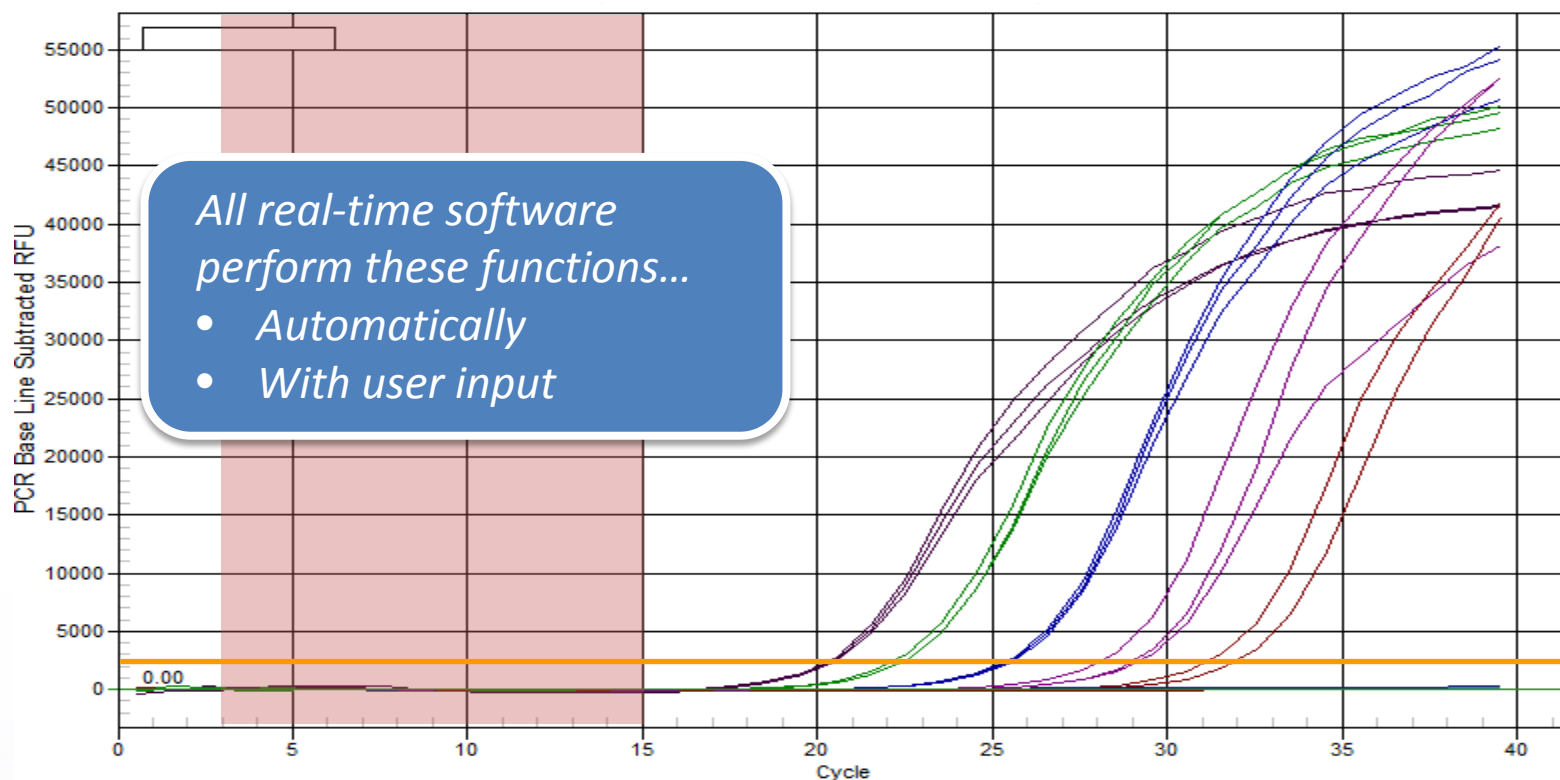
C_q = **quantification cycle** – Cycle number at which amplification curve crosses amplification threshold (aka C_t) – this is the “take-away” metric...

C_q value is inversely proportional to amount of starting template



Steps in the analysis of amplification curves

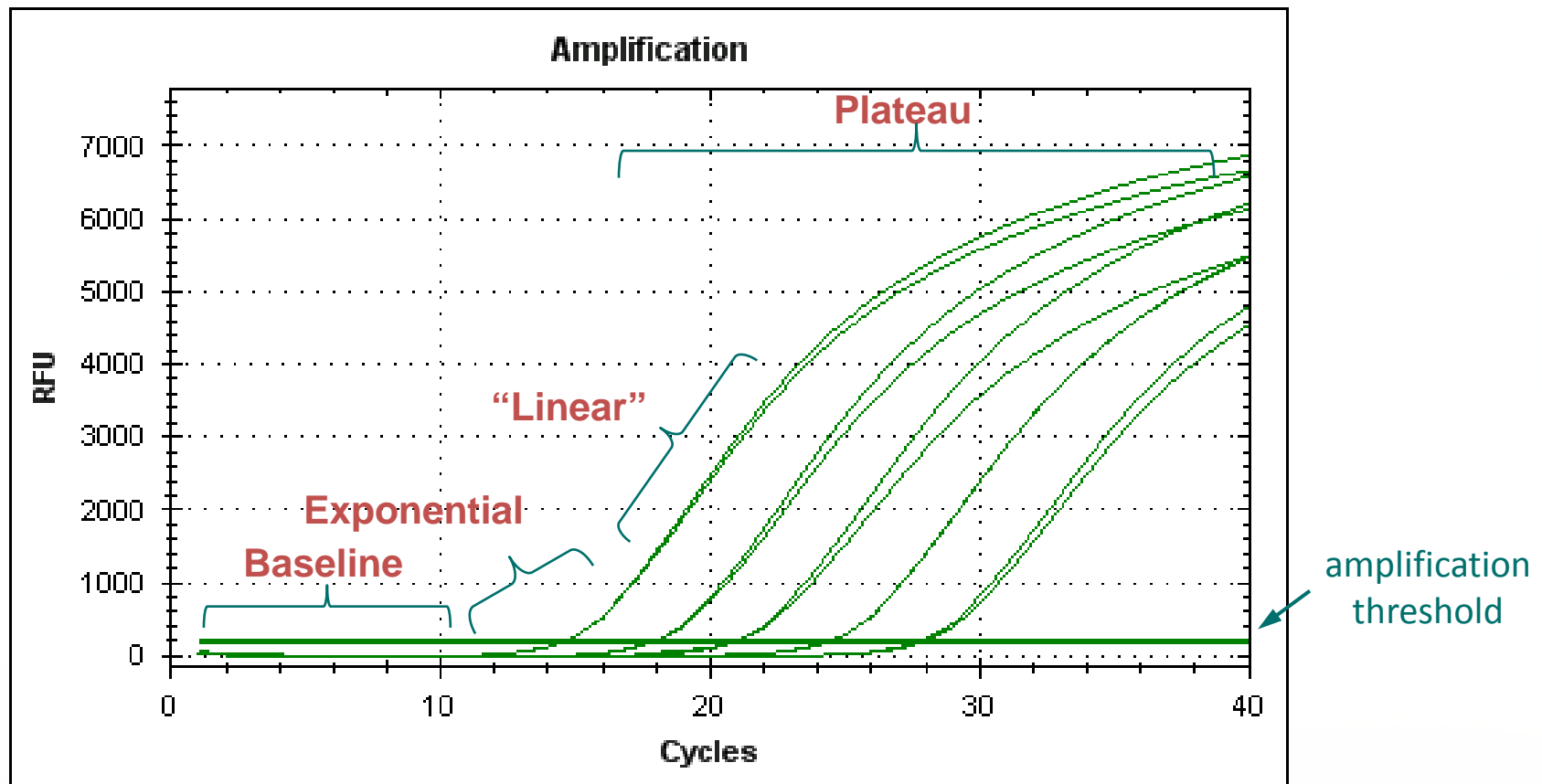
- Passive reference normalization is applied (R_n)
- Baseline regions are defined for each amplification curve
- Curves are baseline-corrected (subtracted & de-trended) (ΔR_n)
- Threshold is set (function of noise in baseline regions for all samples)



Graphs of amplification curves can be re-scaled to reveal more detail



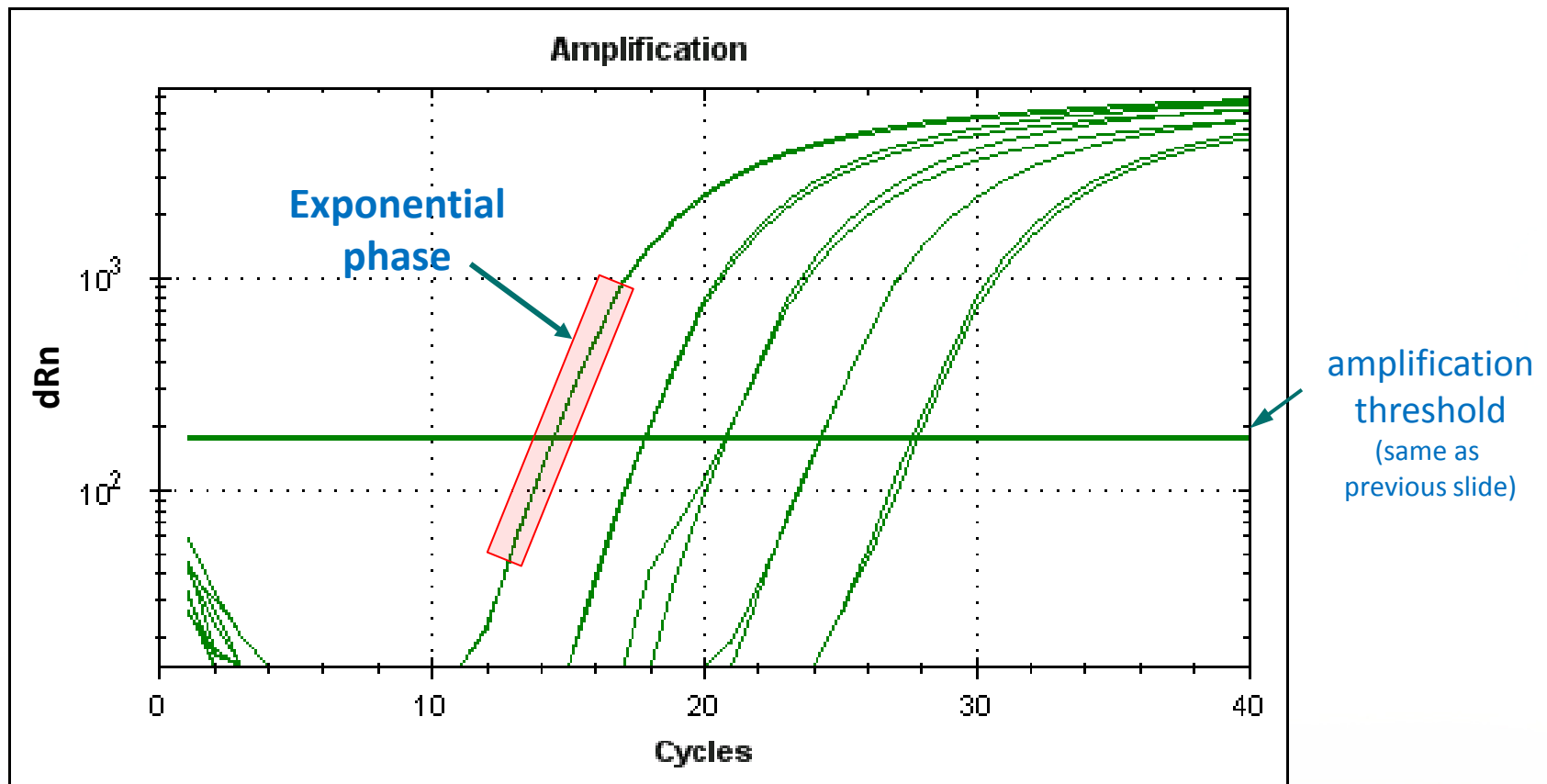
Standard plot of amplification curves allows you to see the baseline phase & plateau phase, but can't really tell anything about the exponential phase!



Graphs of amplification curves can be re-scaled to reveal more detail

Semi-log plot of amplification curves emphasizes exponential phase...

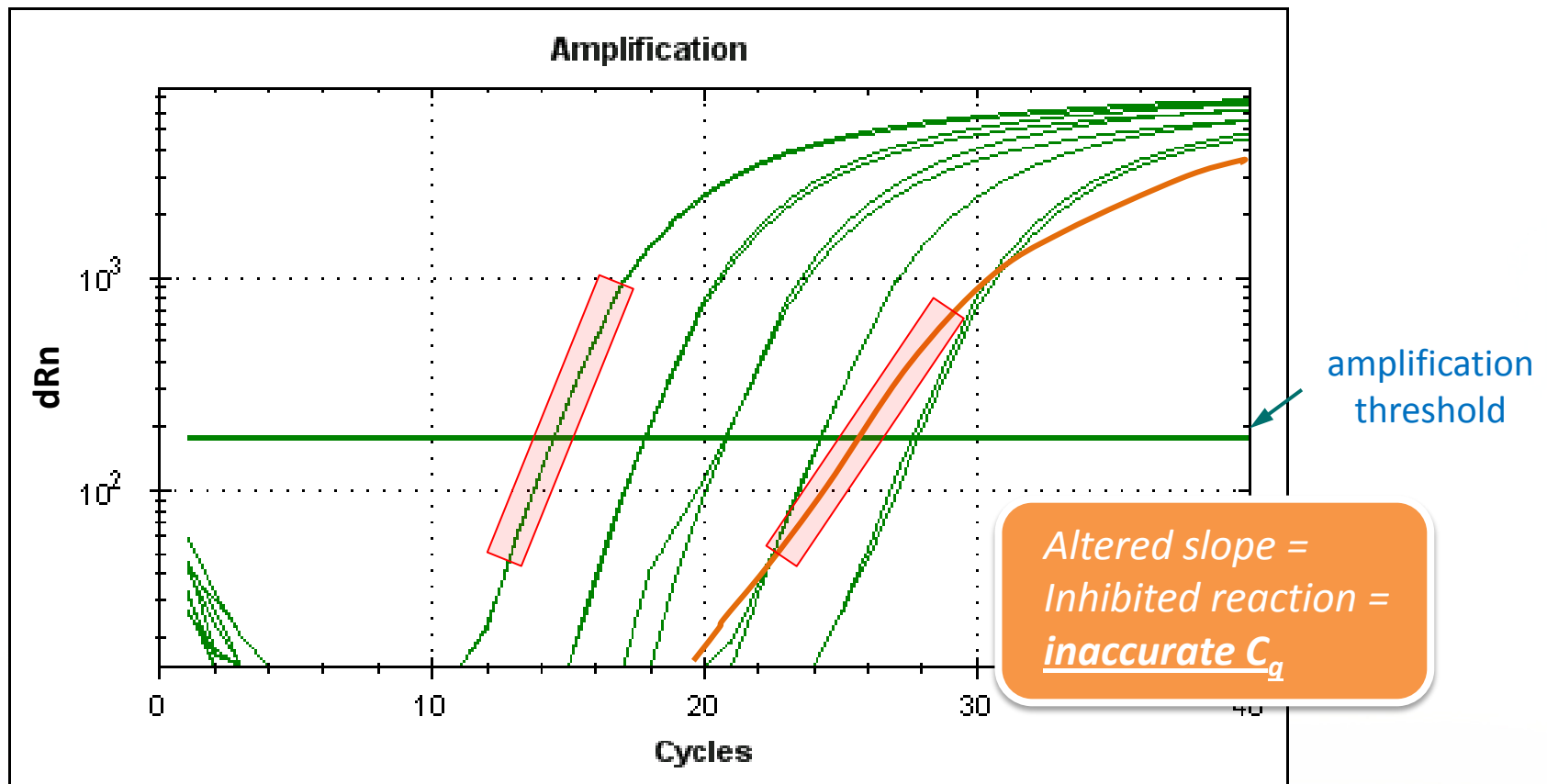
- Allows better visualization of amplification threshold crossing (C_q)



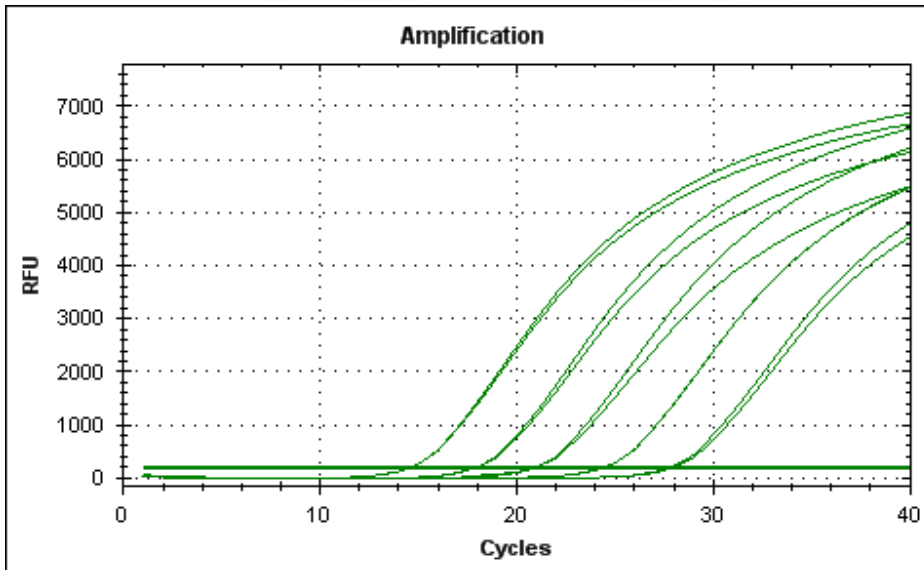
Graphs of amplification curves can be re-scaled to reveal more detail

Semi-log plot of amplification curves emphasizes exponential phase...

- Provides information about efficiency of individual reactions

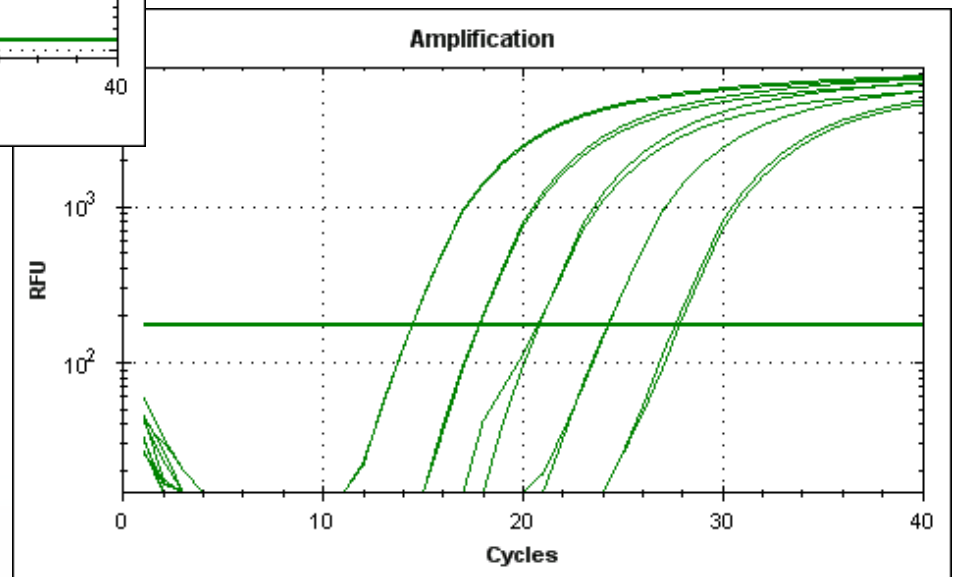


Experienced users utilize both views of amp curves



*Check baseline settings
in standard plot...*

*...check threshold setting
& amplification curve
quality in semi-log plot*



Converting C_q values to quantity - Two approaches



Now you have C_q values – how do you use them?

Absolute (Standard Curve) Quantitation

- Use C_q values to determine amount in unknown samples based on standard curve
- Normalize the amount of target relative to...
 - internal reference (another target that is always at the same level, e.g. GAPDH, beta-actin, 18S rRNA, amplified in same or parallel reaction)

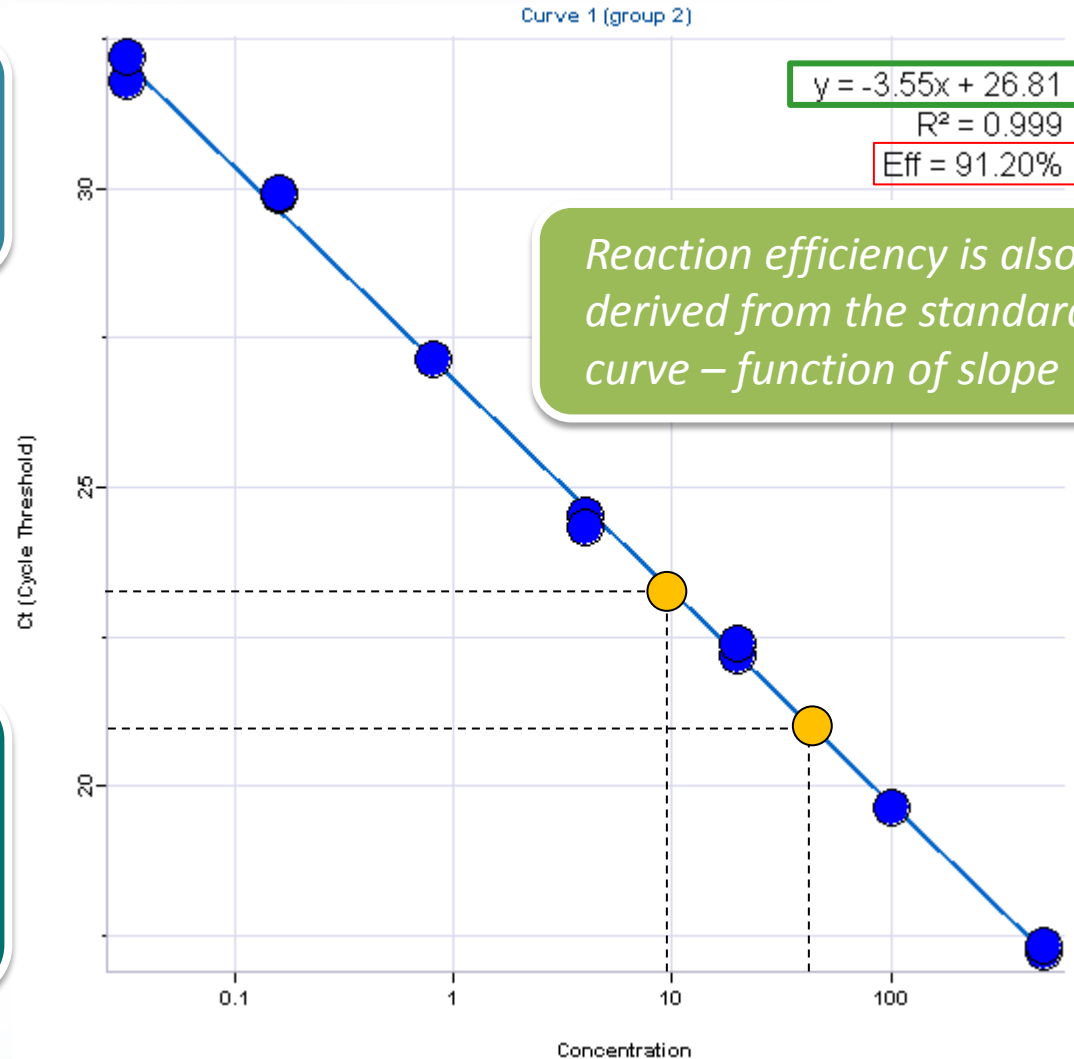
Relative Quantitation

- Compare the C_q values of target in test sample versus control samples - ΔC_t
- Can also normalize the amount of target in each sample relative to internal reference (e.g., GAPDH, beta-actin, 18S rRNA) - $\Delta\Delta C_t$

Standard Curve Quantitation

Standard curve is made
by plotting
 C_q vs (log) concentration

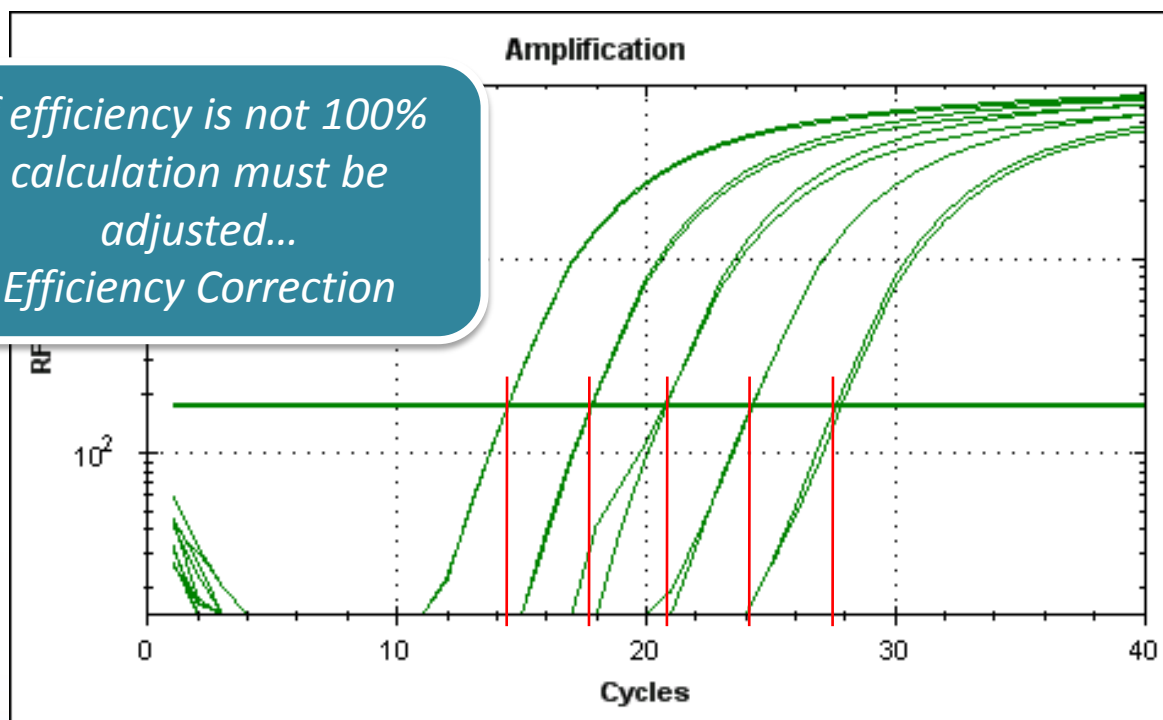
Concentrations of
unknowns
are derived from C_q
based on standard curve



Relative Quantitation

- Based on tenet that in PCR each template is replicated at each cycle...
- Therefore, product formed after n cycles = 2^n (assuming 100% efficiency)
- *The relationship works in reverse...* if two samples have a C_q difference of 1 (threshold is reached 1 cycle apart), then they have $2^1 = \text{two-fold}$ difference in starting template concentration

*If efficiency is not 100%
calculation must be
adjusted...
Efficiency Correction*



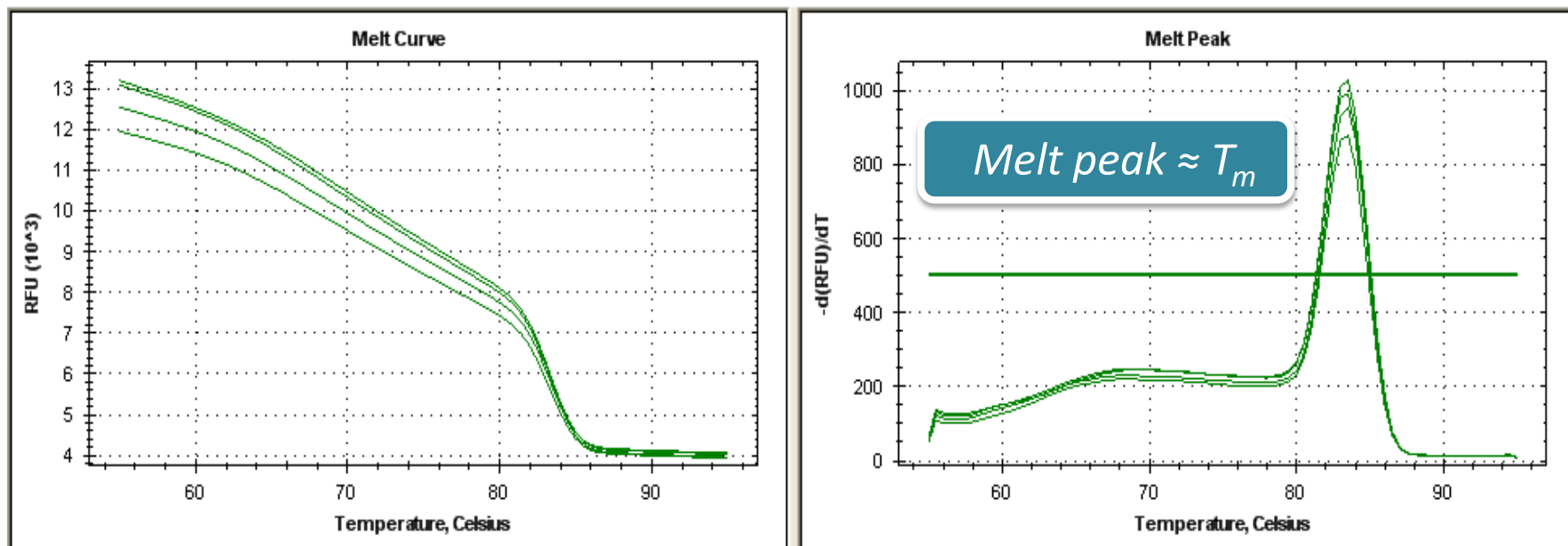
ΔC_q (n cycles)	Difference in concentration (2^n)
3.3	10
5	32
10	1024
15	~32k
20	~1M

Real Time PCR Output may include Melt Data

- Produced in a second, linked thermal profile performed after amplification
- Product is heated slowly, signal is continually measured
- As dsDNA amplicon denatures, signal changes

Provides qualitative information about PCR products – primarily, number & size

Graphed as **Melt Curve** (RFU vs T, *left panel*) or **Melt Peak** (dRFU/dT vs T, *right panel*)

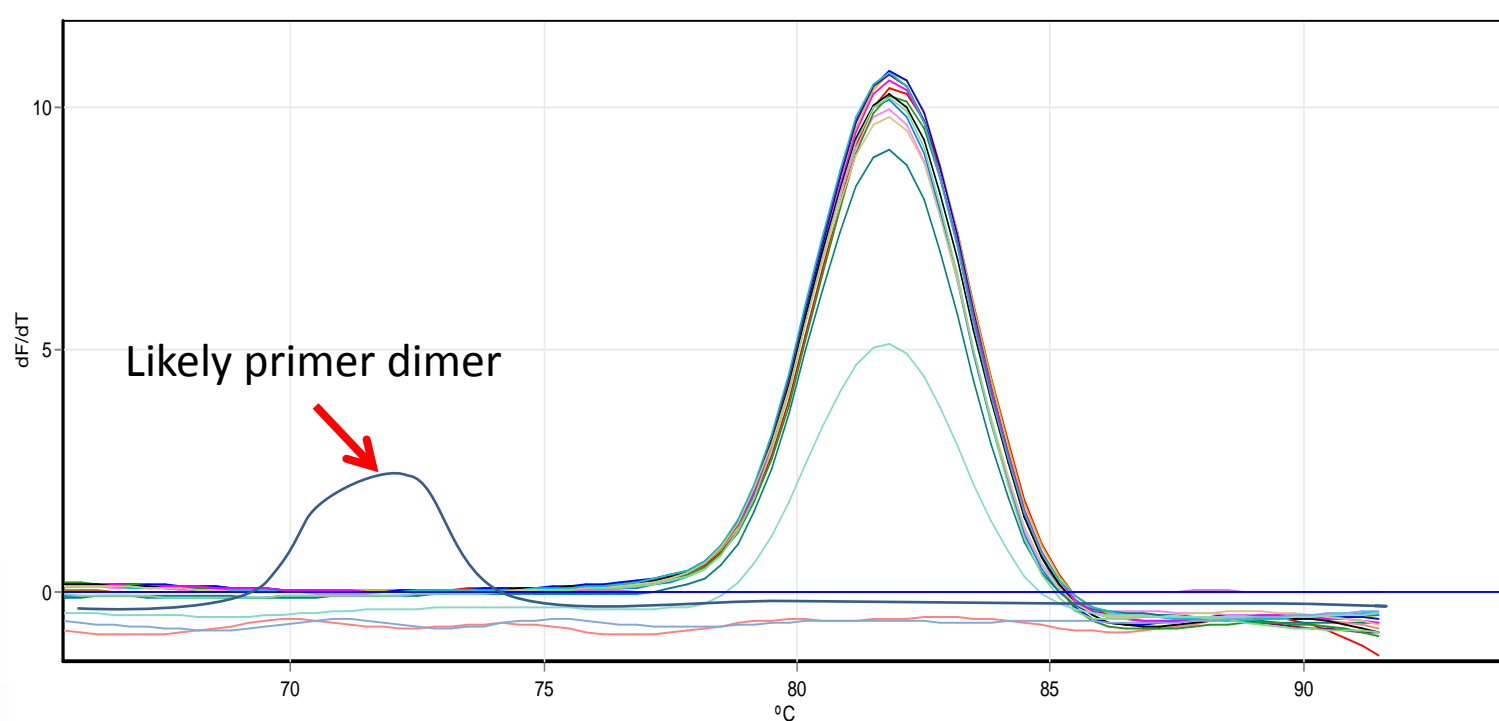


Real Time PCR Melt Data as qPCR Control



What to look for in melt data...

- Melt peak for **samples & positive controls** should be at **similar Temp (T_m)**

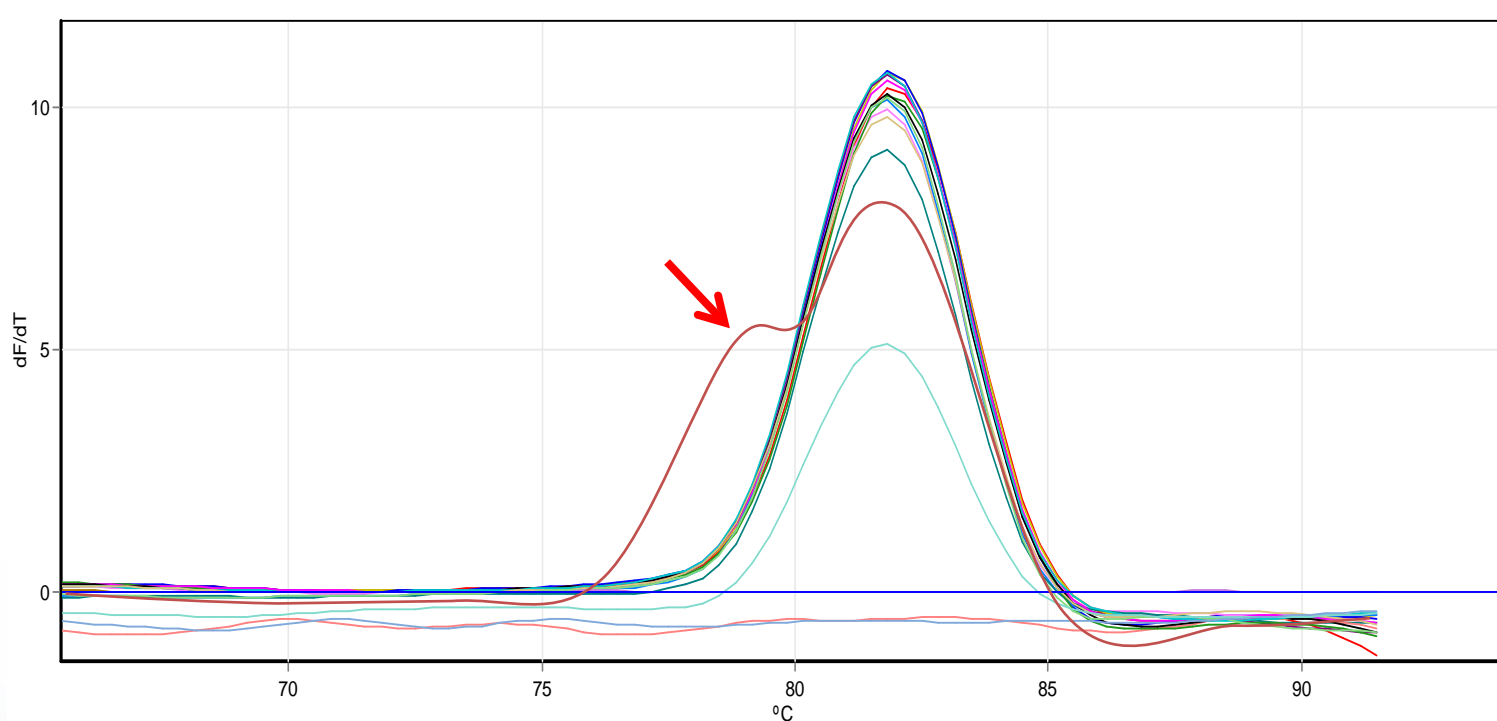


Real Time PCR Melt Data as qPCR Control



What to look for in melt data...

- Melt peak for **samples & positive controls** should be at **similar Temp (T_m)**
- **No secondary peak or shoulder** - indicates secondary (non-specific) product formation

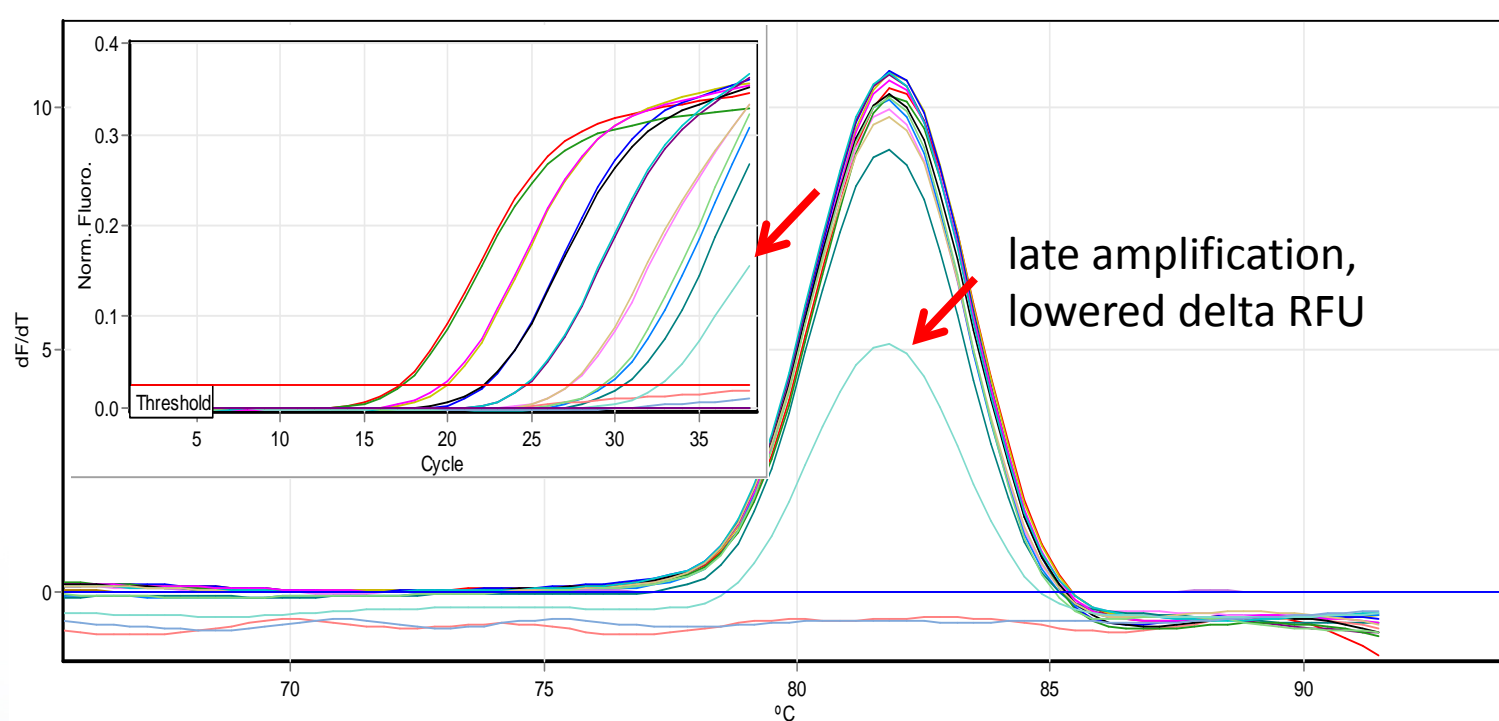


Real Time PCR Melt Data as qPCR Control



What to look for in melt data...

- Melt peak for **samples & positive controls** should be at **similar Temp (T_m)**
- **No secondary peak or shoulder** - indicates secondary (non-specific) product formation
- *Melt peaks height of samples may not be the same – that's not necessarily bad*



Real Time PCR Output – Melt Data

Some Real-Time PCR Chemistries can produce melt data...

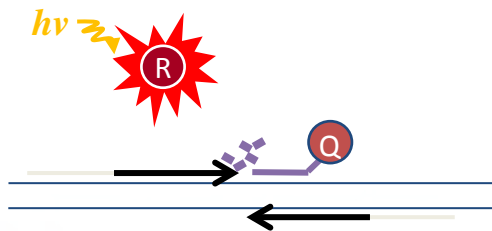


dsDNA dye & Plexor®

fluorescence signal modulation is reversible

Some Real-Time PCR Chemistries can't ...

In this case, gel analysis is crucial during validation



TaqMan® assay

because signal generation is irreversible...
melt analysis is *NOT* possible

RT-qPCR Assay Design



Which reverse transcription (RT) strategy?

- **1-step RT-qPCR**: RT & qPCR in one tube, one reaction set-up.
 - *Gene-Specific Primers (GSP) used for both RT & qPCR*
 - *Reverse primer is the RT primer*
 - *1 aliquot of RNA sample is consumed for 1 qPCR reaction*
 - *With dye-based chemistry this is necessarily a monoplex*
 - *With label-based chemistry can multiplex >1 target*
 - *May be the most sensitive approach*
- **2-step RT-qPCR**: RT reaction done separately from qPCR
 - *Oligo-(dT) &/or random primers used for RT (GSP used for qPCR)*
 - *Either primes all transcripts in an unbiased way (theoretically)*
 - *Random primers will prime all RNA; Oligo-(dT) only poly-(A) RNA (mRNA)*
 - *1 aliquot of RNA can be used for multiple qPCR reactions,*
 - *to quantify multiple targets; or for technical replicates*

RT-qPCR primer design considerations

- Target sequence is from the correct organism
- RefSeq is used (or validated mRNA)
- Paralogs, or conserved motifs in other genes
- Species or strain variation (SNPs or INDELs)
- Amplicon (or primer) should span Exon:Exon junction
- Alternate transcripts
- RT primer position
 - if Oligo-d(T) used, 3'-target may be more sensitive*
- Size of amplicon
 - 75-125 bp is typical range*

*NCBI mRNA accession pre-fixes
NM_*, XM_* = reference mRNA*

RT-qPCR Primer Design Resources



Pre-designed qPCR Primers:

Primer Bank - <http://pga.mgh.harvard.edu/primerbank/>

RTPrimerDB - <http://www.rtprimerdb.org/>

Primer design software:

Primer3 - http://www-genome.wi.mit.edu/cgi-bin/primer/primer3_www.cgi

Primer-BLAST - <http://www.ncbi.nlm.nih.gov/tools/primer-blast/>

IDT PrimerQuest - <http://www.idtdna.com/Scitools/Applications/Primerquest/>

Sequence Resources

Entrez Gene



- Gene Name & Aliases
- Reference Sequence
- Gene organization
- Evidence Viewer – cDNA alignments

Validation of RT-qPCR Primers is Essential

- *You have a new primer design – what next?*



BLAST[®] - *in silico* check to see that they are specific!

If there are matches to unintended genes, evaluate:

- *Match at 3' end?*
- *Percent identity*
- *Predicted T_m of interaction*

- *The in silico analysis looks good, & you order them – now what?*

Experimental Validation! - *Test them on a dilution series of positive control sample...*

- *Determine efficiency*
- *Determine linear dynamic range*

*This is essential ...
even if primers are from a
bank or previously published!*

MIQE is a valuable reference



Minimum Information for Publication of Quantitative Real-Time PCR Experiments

Stephen Bustin et al. (2009) Clinical Chemistry, 55:4

- Recommendations & rational for:
 - *qPCR experimental design, validation, execution, controls, analysis, & presentation.*
 - *Sample handling, nucleic acid extraction, & characterization (quantification & integrity)*
 - *Real-Time PCR terms & nomenclature.*

MIQE

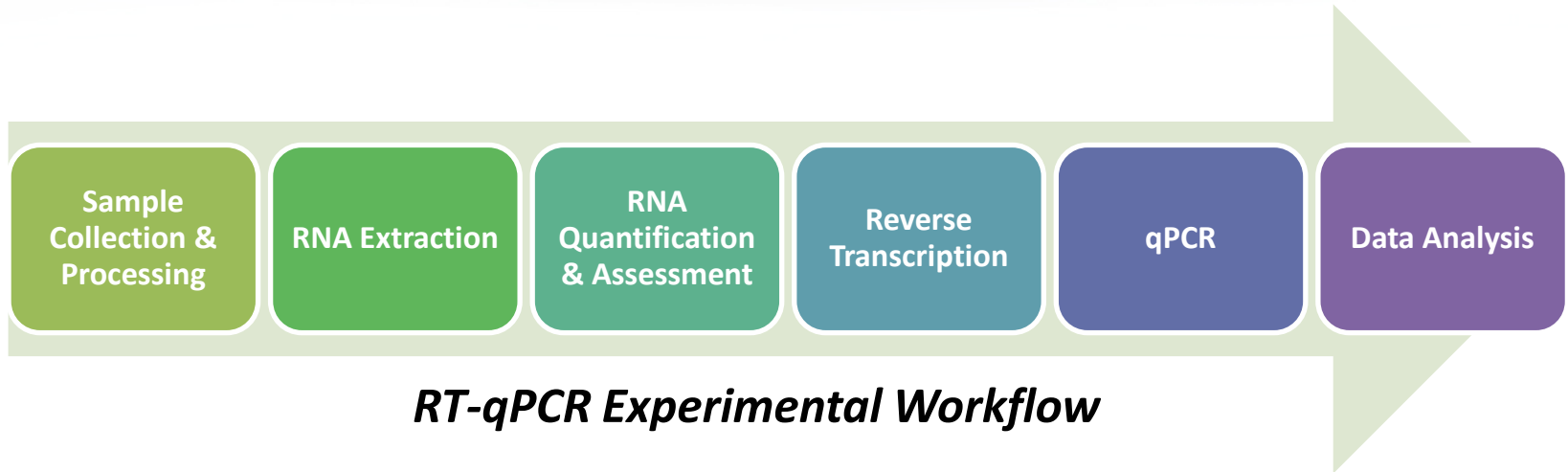
<http://www.rdml.org/miqe.php>

- *pdf of paper*
- *Checklist*

Tips & Troubleshooting



RT-qPCR Tips & Troubleshooting



- Success depends on proper tools & technique at each step
- Problems in any step of the experiment can cause experimental failure, or worse, inaccurate results

Proper handling of samples is the first step to ensure good RNA yield & quality



Degradation & loss of RNA often occurs during sample collection & processing

- Temperature abuse of samples before/during /after collection
 - *Process or store immediately*
 - *Snap freeze on liquid N or dry ice*
- Dissection takes too long
 - *If dissection is difficult, do gross dissection first (quickly)... then fine dissection in a preservative, or after preservation (e.g., RNAlater®)*
- Sample dimensions too large – takes too long to freeze & thaw
 - *Cut into smaller chunks during dissection, before further processing*
- Insufficient tissue disruption
 - *Rotor-stator is generally the best approach*
 - *Dounce homogenizer or blue pestle - may need to chop/mince first*

Consider all parameters when choosing – and using - RNA extraction method



Yield – often the primary consideration

- *Most methods give similar yield - organic extraction often perceived as best...but often at a cost*
- *Exceeding processing recommendations usually does not increase yield (& may compromise purity)*

gDNA removal – high amounts can cause quantitation error; even low amounts can cause problems in qPCR

- *gDNA contamination is often an issue with organic separation methods*
- *DNase can be added after any method*
- *Efficient removal is important, but impossible to remove 100% (& may not be necessary)*

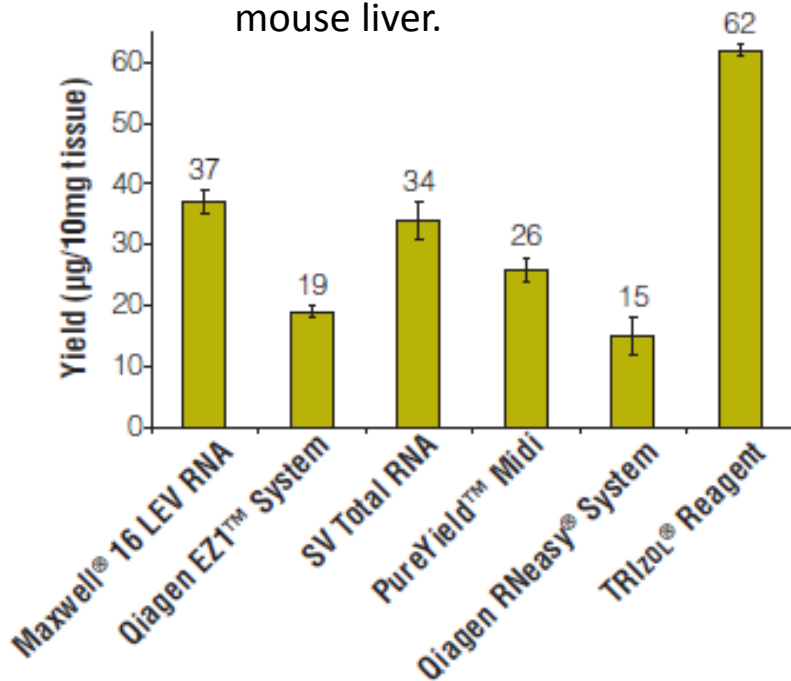
Inhibitor carryover – may lead to variation in Cq's rather than reaction failure.

- *Can be an issue with organic separation methods*
- *Exceeding processing recommendations can compromise purity*

Increased yield at the cost of purity is a poor trade

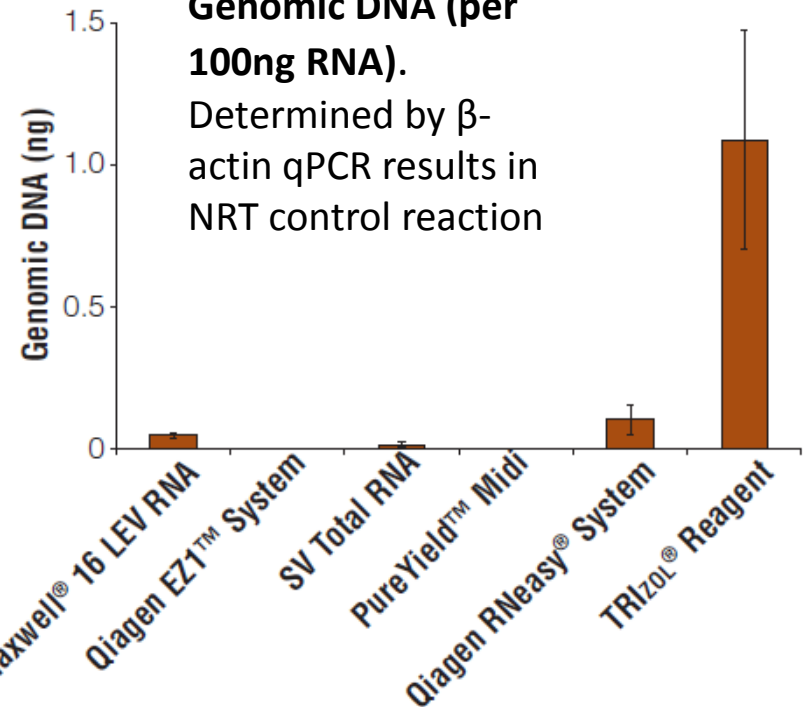


RNA yield.
mouse liver.



Genomic DNA (per 100ng RNA).

Determined by β -actin qPCR results in NRT control reaction

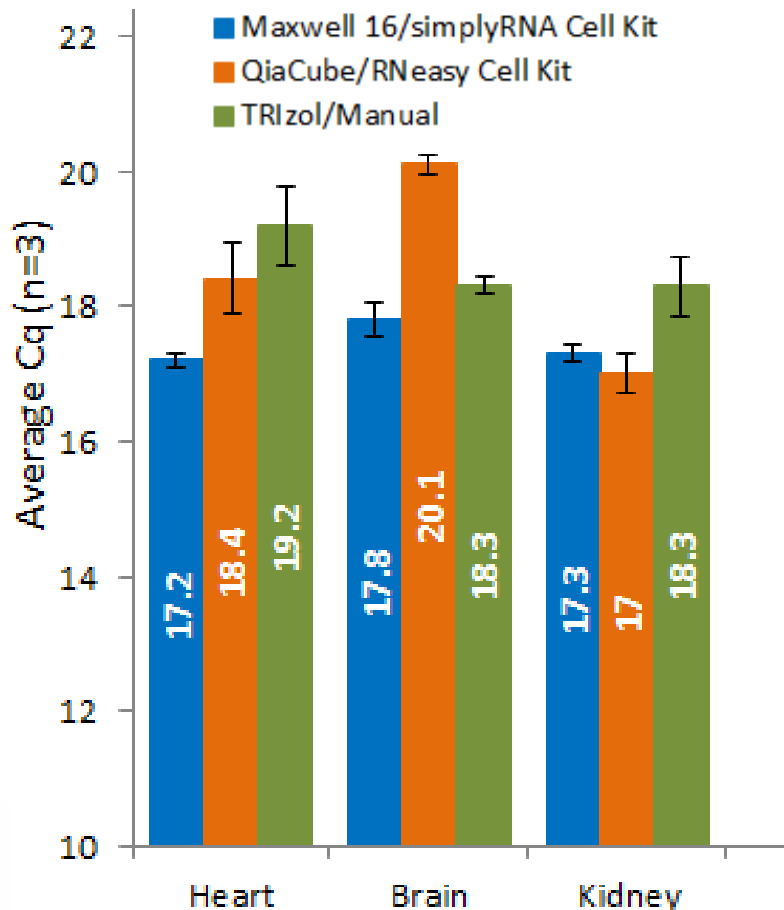


Schagat et al., 2008 Promega Notes 99 "RNA Purification Kit Comparison: Yield, Quality and Real-Time RT-PCR Performance"

75001 MA

75001 MA

Increased yield at the cost of purity is a poor trade



Lower purity may result in higher, more variable C_q s

What is advantage in 2-fold difference in yield in an assay with 10^6 -fold range?

Beware the blind pursuit of maximum yield!

There are many methods to evaluate RNA

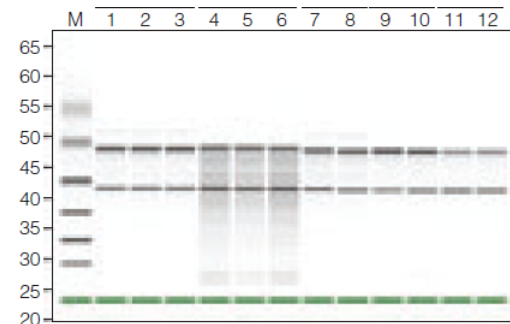


Quantification - Ideally, RNA input amount should be similar in each RT reaction

- Direct absorbance – A260; abs ratios can provide information about purity
- Fluorescent dye – greater sensitivity & dynamic range

Quality assessment –

- Gel or Bioanalyzer -
 - *Crucial to assess integrity*
- Absorbance ratios
- Spike experiment
 - *For inhibitors*
- No RT qPCR control (NRT)
 - *For gDNA contamination*
 - *Not as critical if primers span introns*



ΔC_q	Difference in concentration
3.3	10
5	32
10	1024

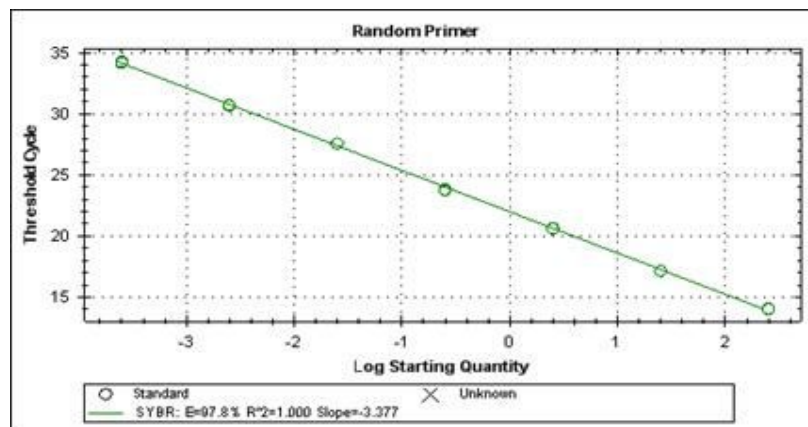
RT reaction considerations

RNA amount – current RT systems have very good range & proportional yield; however, it is still desirable to have similar amounts of total RNA within each rxn

Oligo-(dT) or random primer amount - need not necessarily be adjusted relative to mRNA input amount...

- Total Human RNA
- Diluted 1ug/ul to 1pg/ul - 1ul per rxn
- 500ng random primers per rxn
- GAPDH primers

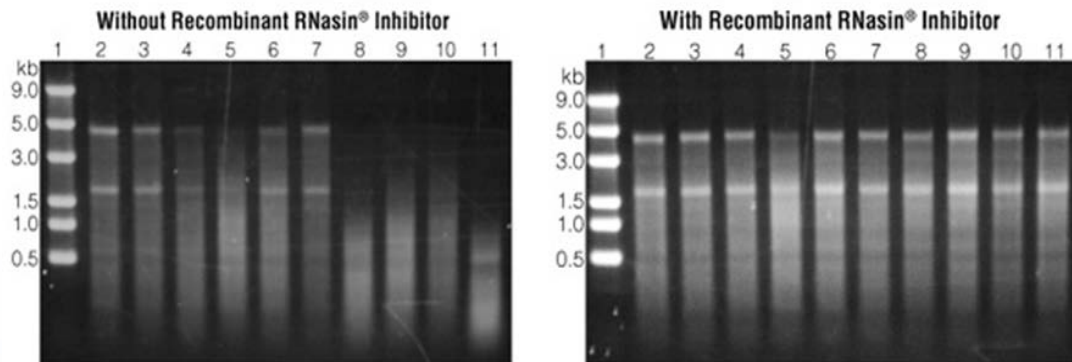
GoScript® 2-Step RT-qPCR System



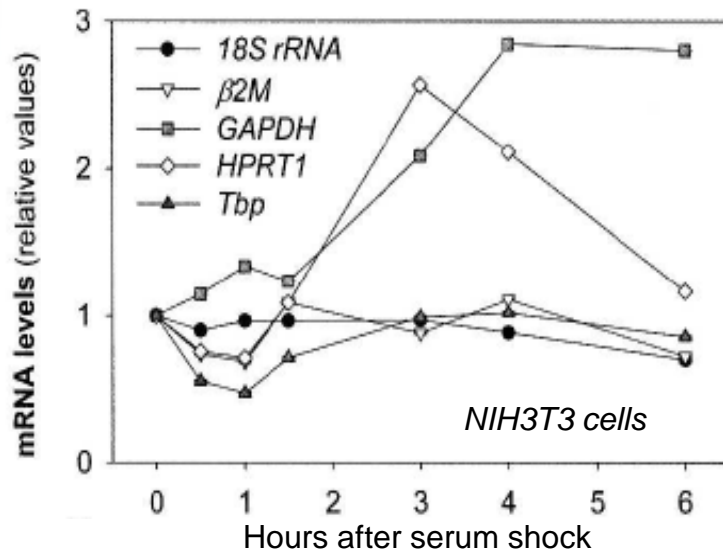
Avoid RNase contamination post-purification

- Follow best practices, e.g., gloves, barrier tips, etc
- Use **RNase inhibitor** in RT reaction

RNasin® Is included in GoScript® & GoTaq® Systems



Stability of normalization target must be verified

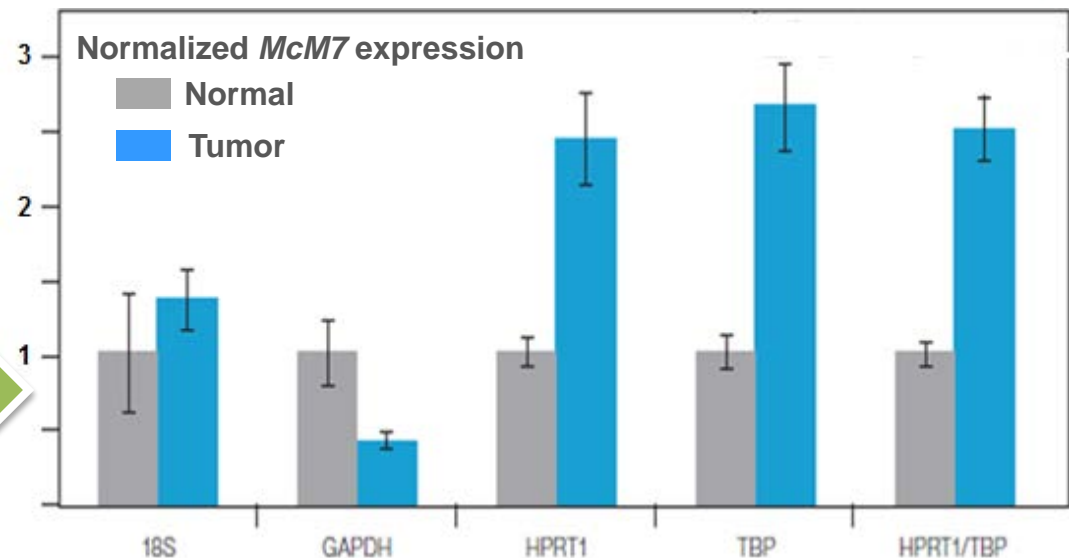


Garabino-Pico, E. et al. (2007) RNA

Commonly used normalizers are not always constitutive

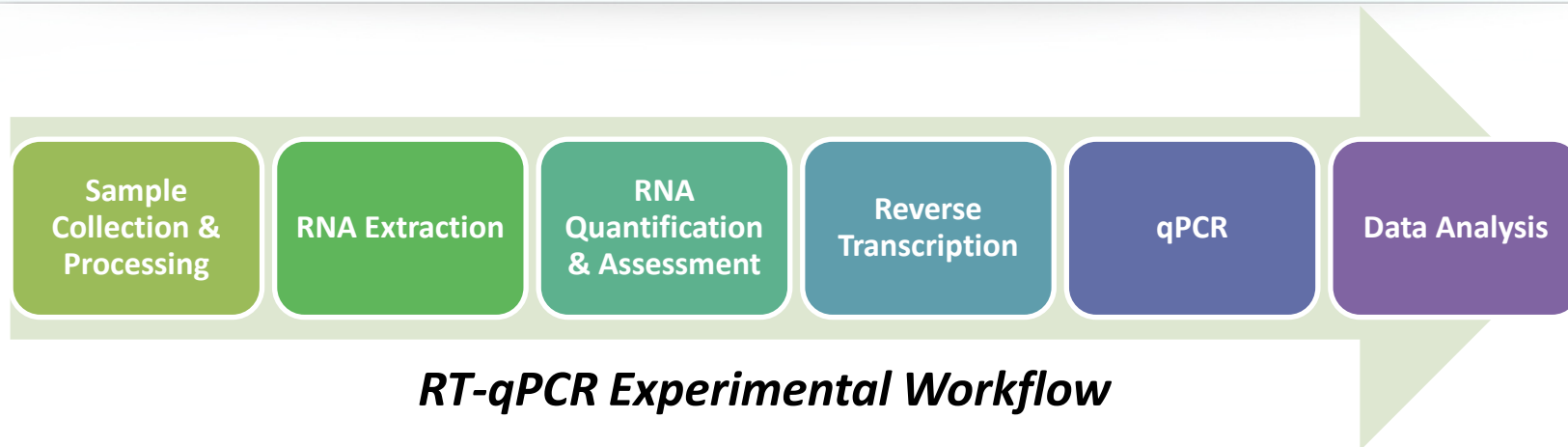
- MIQE suggests two normalizers
- Software exist to evaluate - geNORM

Using the wrong normalizer can qualitatively change the results



Adapted from Taylor, S. (2011) Bio-Rad tech note 6245

Promega Products for RT-qPCR



RNA Extraction

- SV Total RNA Isolation System
- Maxwell® 16 Instrument
- Maxwell® 16 simplyRNA Purification Kits

RNA Quantification

- QuantiFluor™ RNA System
- GloMax®-Multi+ Detection System

Reverse Transcription & qPCR

- GoScript™ Reverse Transcriptase
- GoTaq® qPCR Master Mix
- GoTaq® 1-Step & 2-Step RT-qPCR Systems
- Recombinant RNasin® Ribonuclease Inhibitor
- Plexor® qPCR & RT-qPCR Systems

Questions? Ask a Scientist!

<http://www.promega.com/support/>



Good luck with your experiments!