

How NOT to mess up your qPCR

Promega GmbH



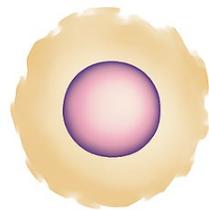
Stimulus-Triggered Acquisition of Pluripotency (STAP)

ARTICLE

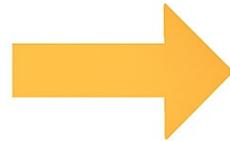
doi:10.1038/nature12968

Stimulus-triggered fate conversion of somatic cells into pluripotency

Haruko Obokata^{1,2,3}, Teruhiko Wakayama^{3†}, Yoshiki Sasai⁴, Koji Kojima¹, Martin P. Vacanti^{1,5}, Hitoshi Niwa⁶, Masayuki Yamato⁷ & Charles A. Vacanti¹



Lymphocytes



Mild acid bath



Pluripotent cells

„STAP Cell Scandal“

News | Published: 18 March 2014

Stem-cell method faces

The STAP case scandal: Researcher Haruko Obokata
resigns after failing to

ta

News | Published: 01 April 2014

Stem-cell scientist found guilty of misconduct

STAP Cell Scandal: Japan's Biggest Science Fraud

The Final Word on STAP

Researchers fail to replicate STAP study; computational analysis reveals genomic inconsistency

Lessons Learned: Why Standards Are Crucial

MIQE 2.0 Guidelines

Standardized set of recommendations for

- Designing
- Optimizing
- Validating
- Analysing
- Reporting

of qPCR results

Clinical Chemistry 00:0
1–18 (2025)

Special Report

MIQE 2.0: Revision of the Minimum Information for Publication of Quantitative Real-Time PCR Experiments Guidelines

Stephen A. Bustin ^{a,*} Jan M. Ruijter,^b Maurice J.B. van den Hoff,^b Mikael Kubista,^c Michael W. Pfaffl,^d
Gregory L. Shipley,^e Nham Tran,^f Stefan Rödiger,^g Andreas Untergasser,^h Reinhold Mueller,ⁱ Tania Nolan,^j
Mojca Milavec,^k Malcolm J. Burns,^l Jim F. Huggett ^l Jo Vandesompele ^m and Carl T. Wittwerⁿ

What is Real-Time PCR?

End-point PCR

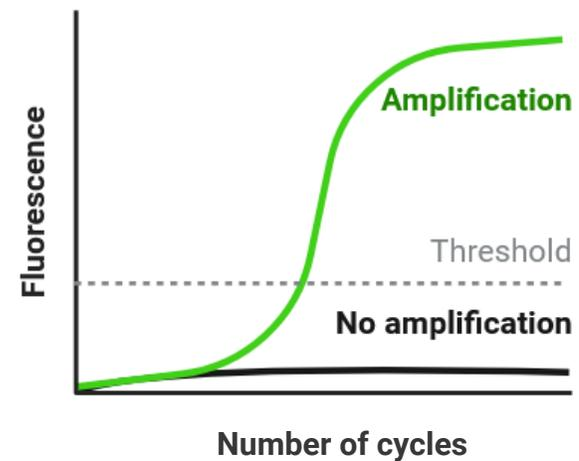
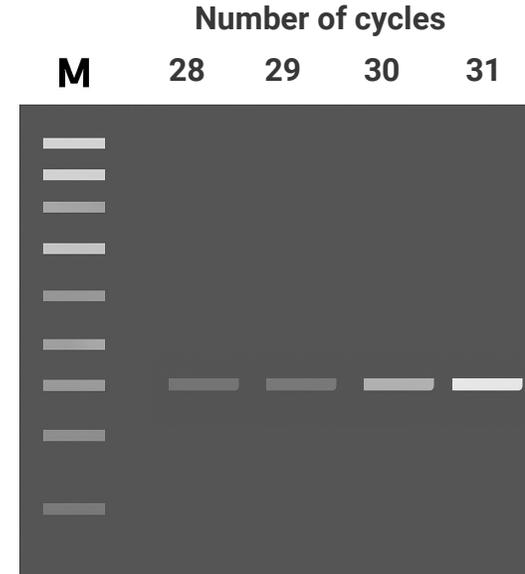
Product formation measured

after reaction is complete

Real-Time PCR

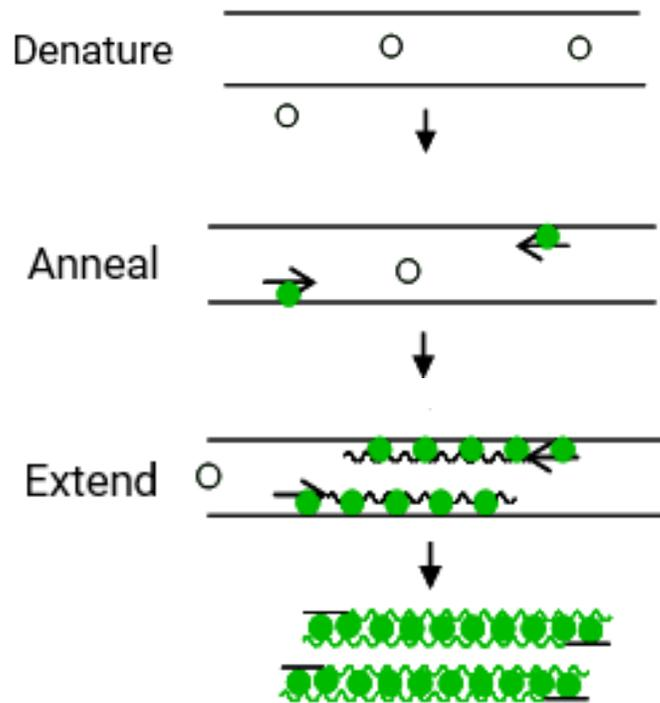
Product formation measured

at each cycle, during the reaction

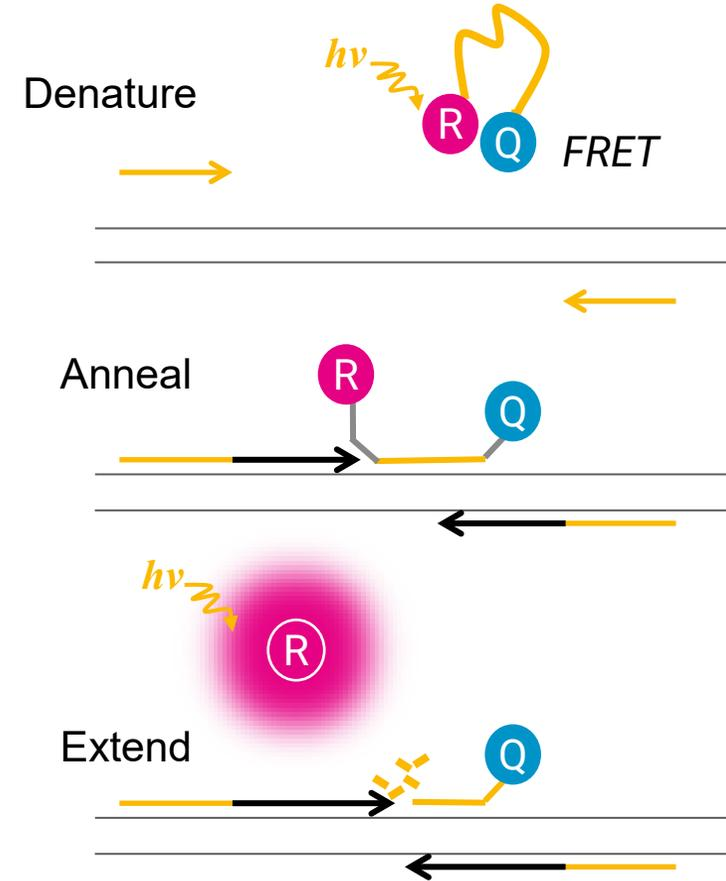


Real-time PCR Chemistries: **Dye-based** vs. **Probe-based** qPCR

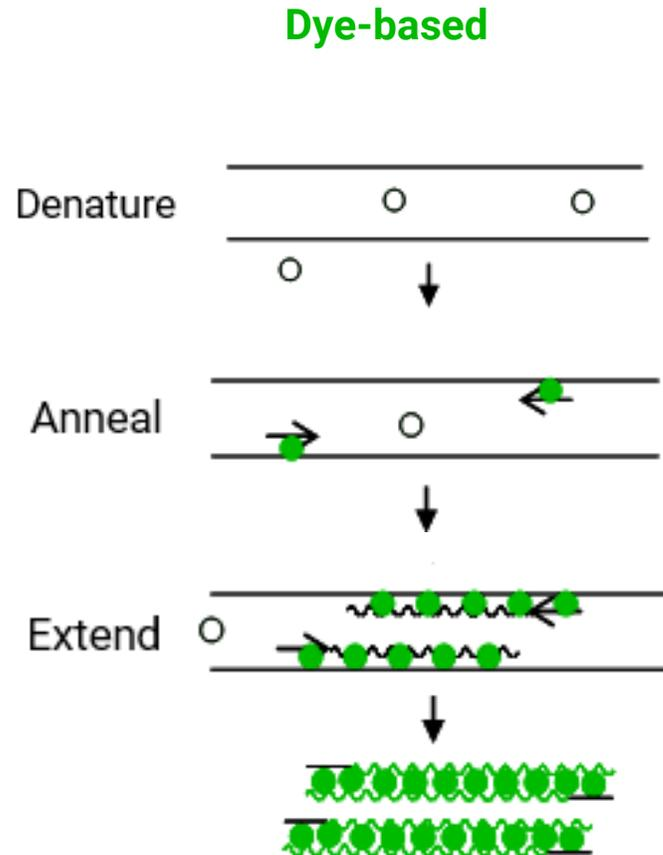
Dye-based



Probe-based



Real-time PCR Chemistries: **Dye-based**



- dsDNA-binding dye is included in PCR master mix
- Standard primers used
- Dye associates with PCR product

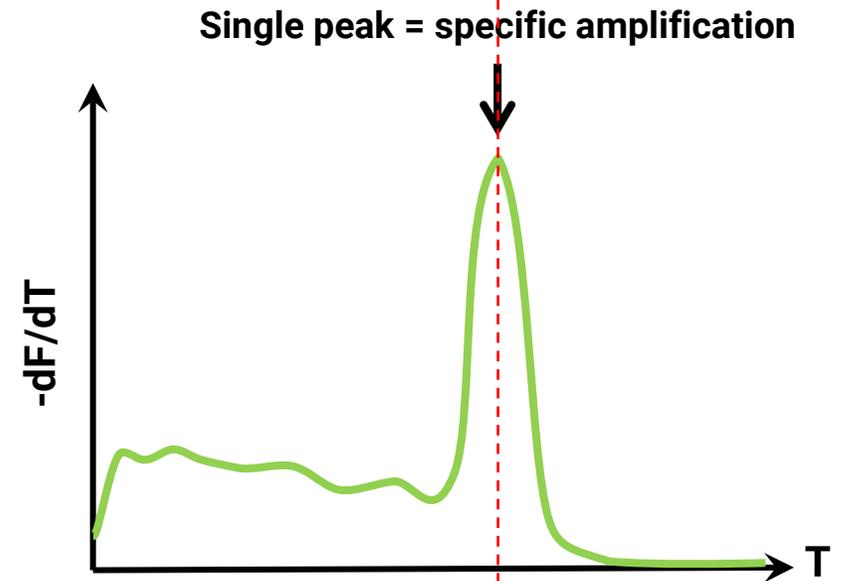
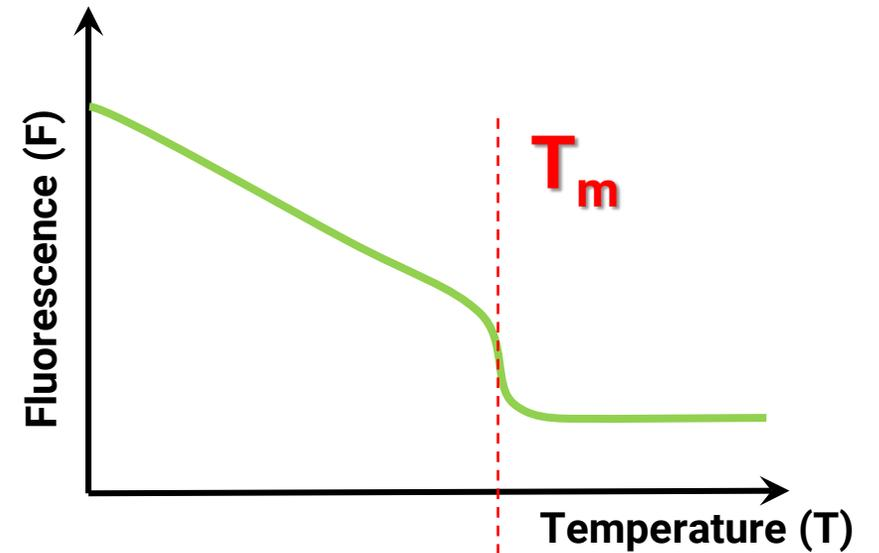
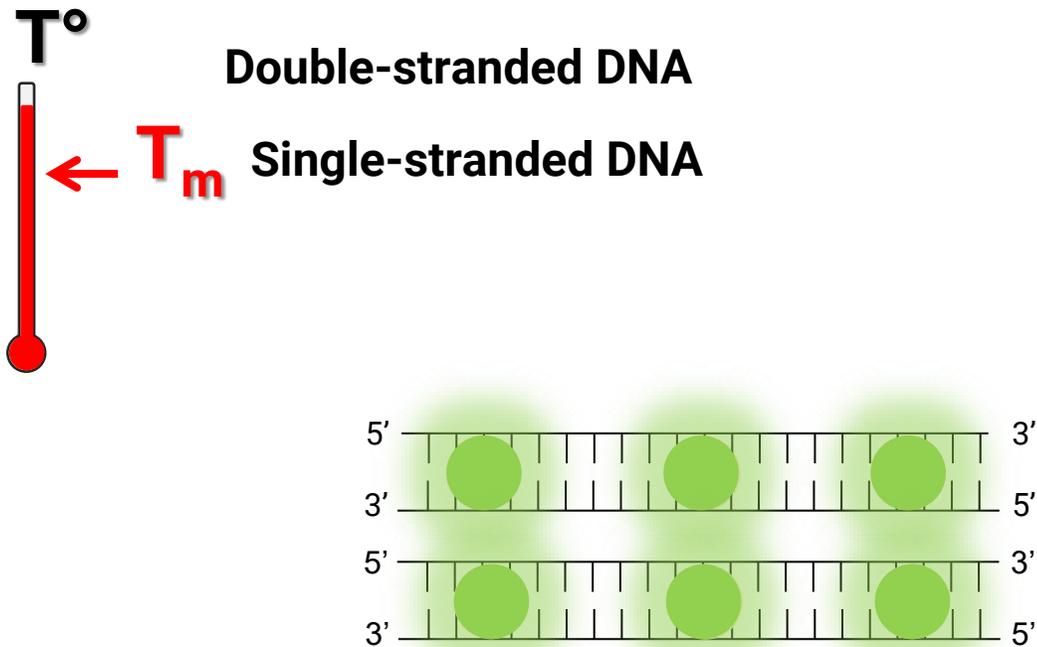
- Free Dye → low fluorescence
- Bound Dye → high fluorescence

Fluorescence is proportional to the amount of PCR product

As more PCR product is produced, more dye is bound

Melting Curve

- Produced in a second, linked thermal profile performed after amplification
- Product is heated slowly, signal is continually measured



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

Real-time PCR Chemistries: **Probe-based** qPCR

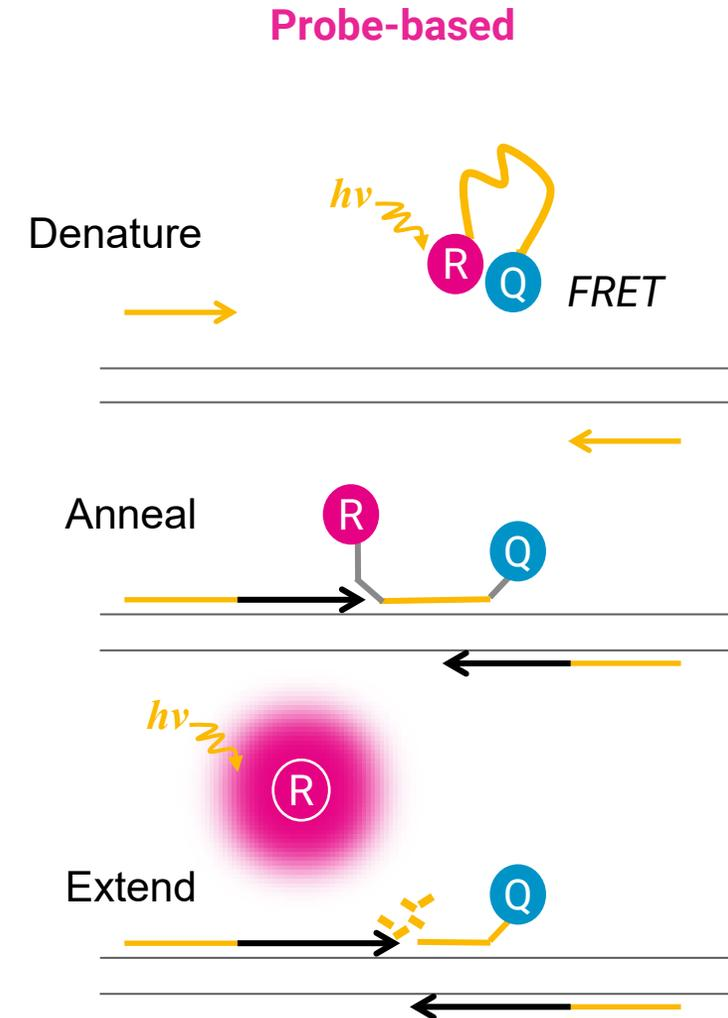
TaqMan[®] is the most familiar type:

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

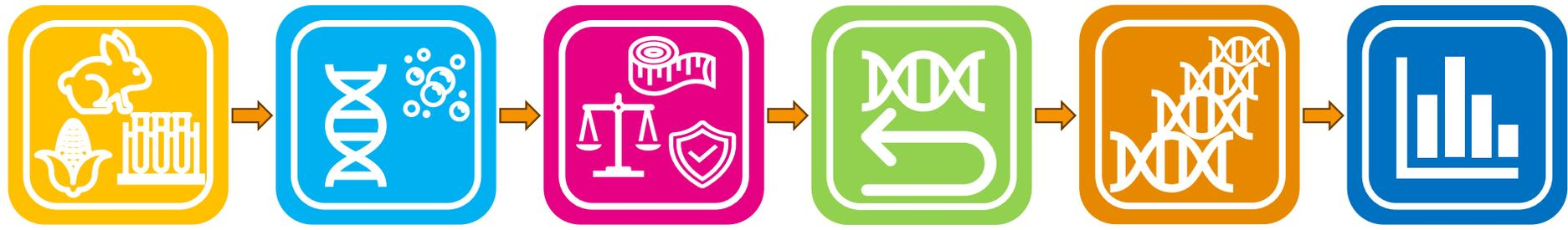
Free probe → FRET occurs

Degraded probe → reporter unquenched

Fluorescence is proportional to the amount of proper PCR product



Experimental Workflows



RT-qPCR

Sample collection & Processing

RNA Extraction

RNA QC & Protection

Reverse Transcription

qPCR

Data Analysis

qPCR

Sample collection & Processing

DNA Extraction

DNA QC

qPCR

Data Analysis

What general precautions should be taken when working with nucleic acids?



- Change **gloves** regularly
- **RNase-** and **DNase-free environment**

- Avoid **cross-contamination**:
 - Use sterile, certified RNase- and DNase-free **pipette tips**
 - **Separate work areas** pre- & post-amplification
 - Uracil N-glycosylase (**UNG**), **dUTP** instead of dTTP
- Avoid gDNA contamination:
 - Extended incubation with **DNase I**
 - **gDNA shearing** with syringe

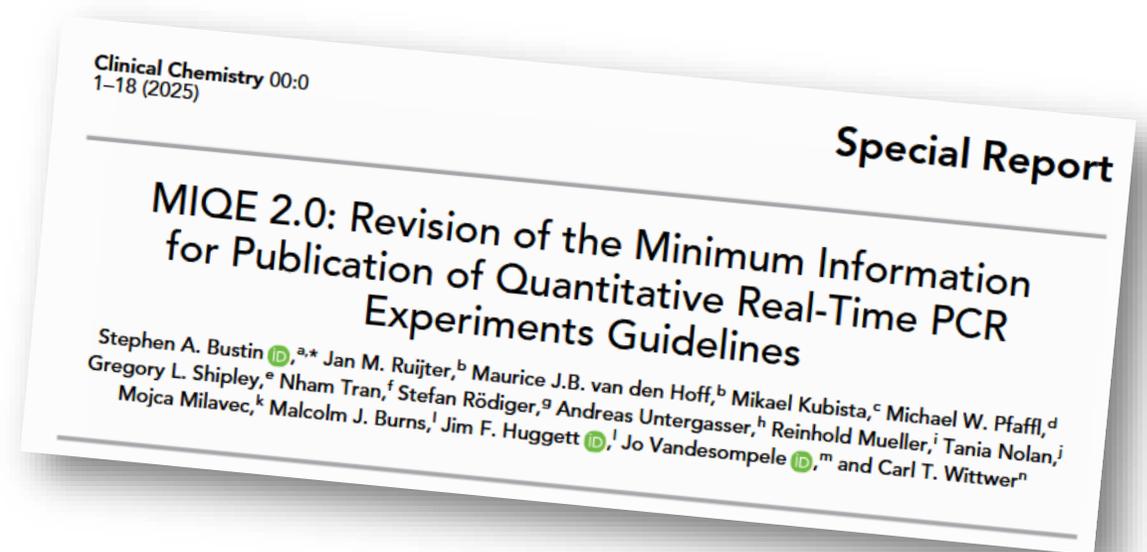


Ask Yourself 5 Questions Before You Publish Your qPCR Results

1. Can I Trust This Sample As A Valid Starting Point For My Quantification?
2. Did I Normalize The Right Way?
3. What Is My Result And What Does It Mean? How Do I Interpret Cq Values?
4. What Controls Have My Back?
5. Can Someone Else Re-Run My Analysis?

Keep in mind: **MIQE isn't paperwork, it's guardrails.**

It ensures that you work cleanly, remain confident, and don't lose valuable time.





1. Can I Trust This Sample As A Valid Starting Point For My Quantification?

How Do I Prepare My Sample Correctly?

Maxwell® RSC Instruments

→ Research Use Only



Efficient

Up to 16 or 48 samples simultaneously



Isolation of high-quality DNA, RNA or TNA

Flexible

Wide range of samples types
From Aspergillus to Zebrafish

Maxwell® CSC Instruments

→ CE-IVD certified



Time-saving

On average < 45 minutes per run

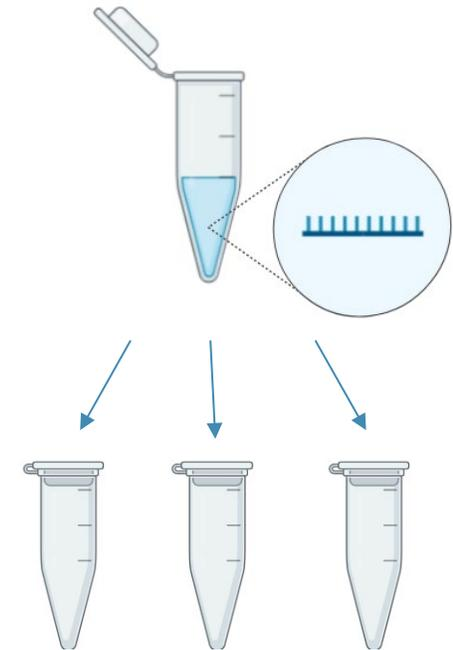
1. Can I Trust This Sample As A Valid Starting Point For My Quantification?



How do I store my sample correctly?



- Extracted RNA:
 - 20°C → days,
 - 80°C → months,
 - 180°C (liquid N₂) → 5 years
- Store DNA/RNA in small aliquots
- Avoid multiple freeze-thaw cycles
- Check nucleic acid concentration after each thawing



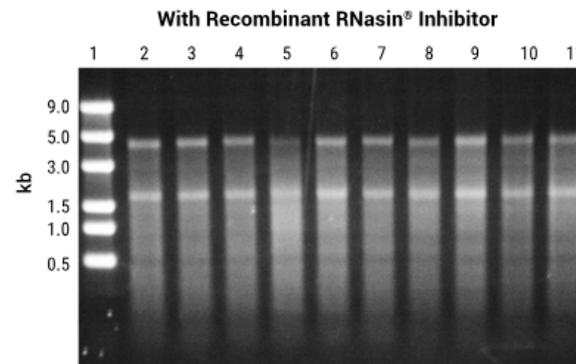
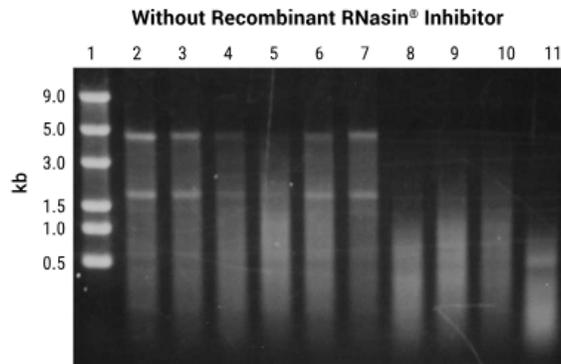


1. Can I trust this sample as a valid starting point for my quantification?

Is my sample intact or degraded?

Electrophoresis

- Size
- Integrity
- Quantity



Dye Staining (QuantiFluor)

- Concentration (very sensitive)

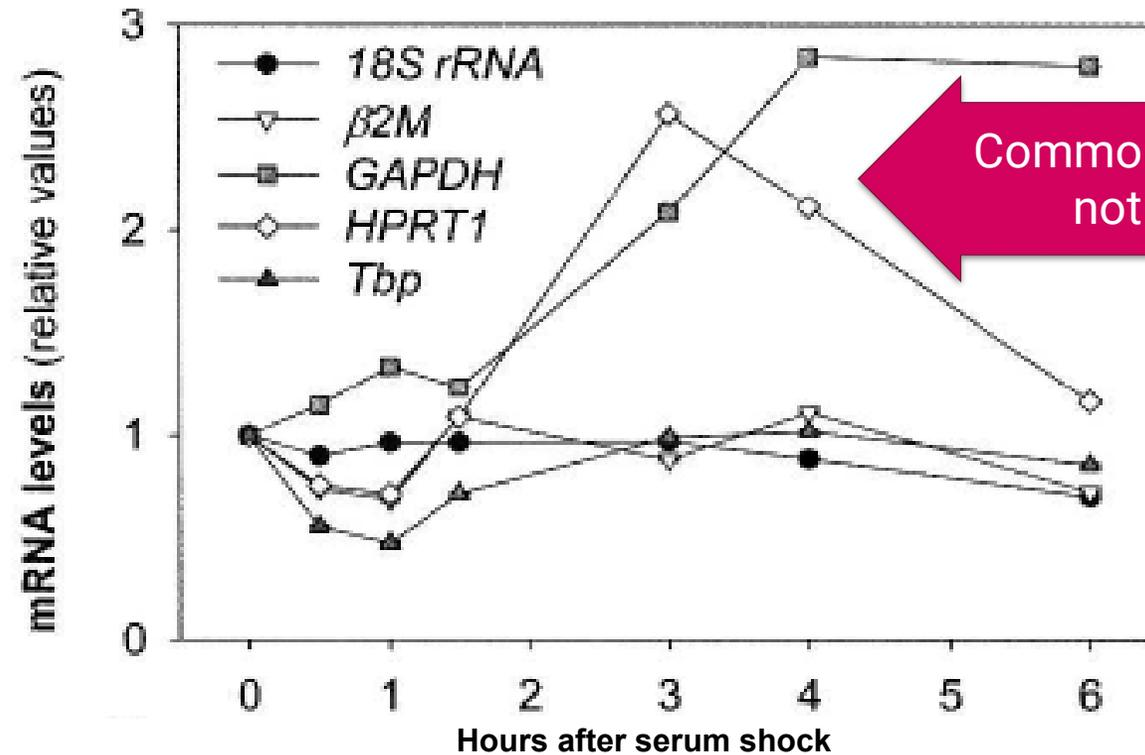


MIQE suggests **combination of electrophoresis & dye staining** to assess **size, quantity, concentration and integrity** (although this does not reveal contamination)

Quality is more important than yield!

2. Did I Normalize the Right Way?

Reference genes

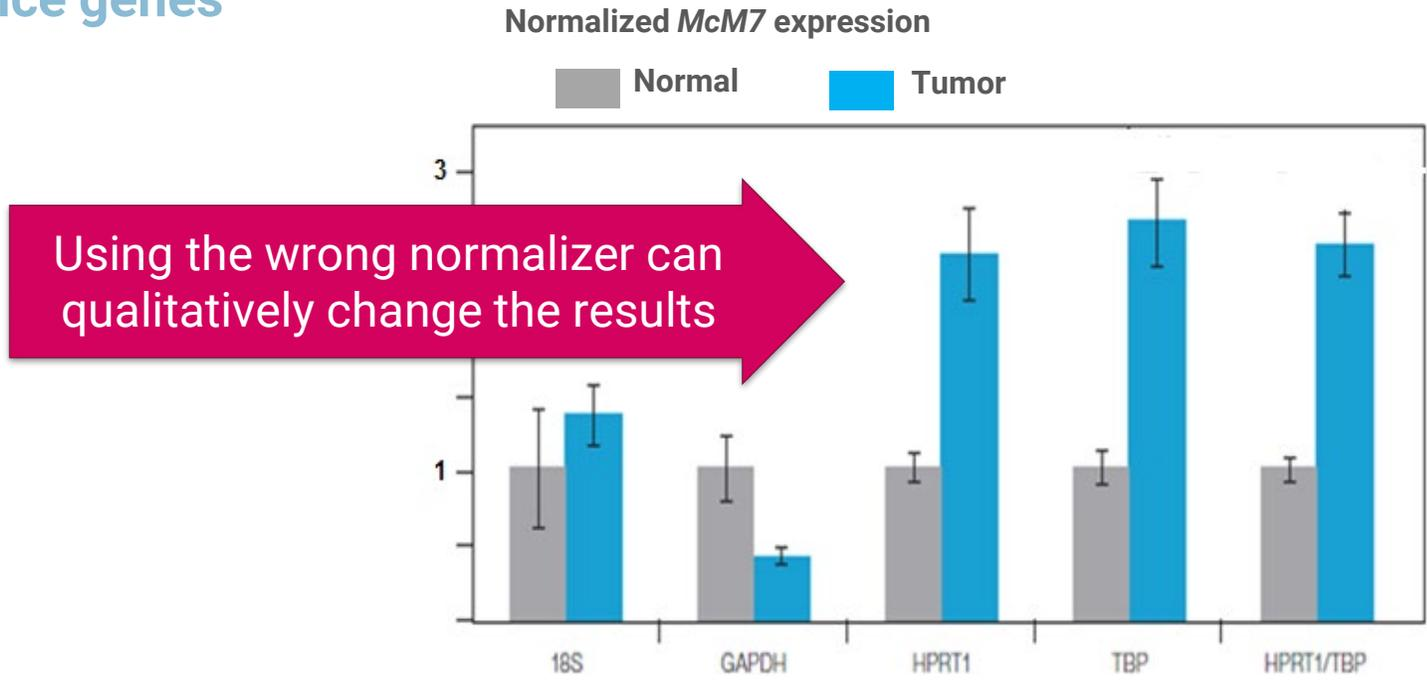


Commonly used normalizers are not always constitutive

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

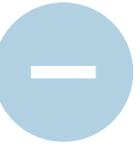
2. Did I Normalize the Right Way?

Reference genes

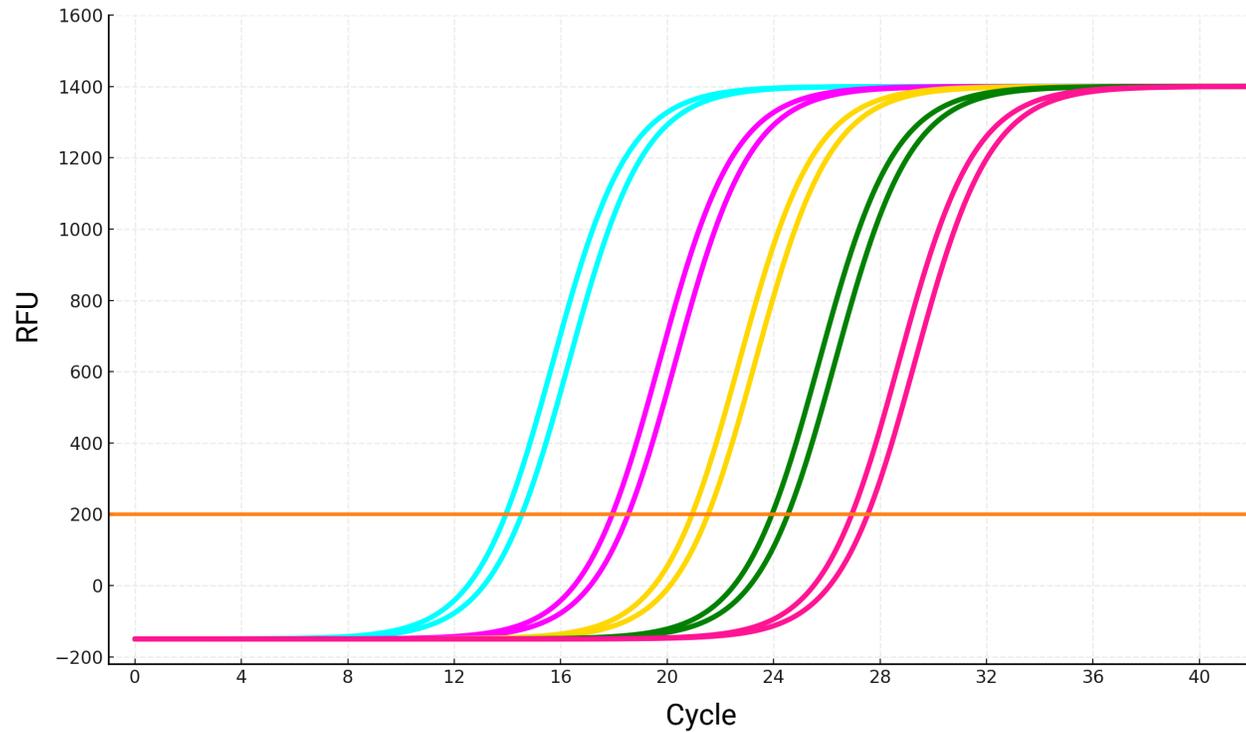


Adapted from Taylor, S. (2011) Bio-Rad tech note 6245

- MIQE suggests at least two normalizers
- Software to evaluate:
geNORM (<https://genorm.cmgg.be/>)
- Expression of normalizers should be in the range of your GOI



3. What Is My Result And What Does It Mean? How Do I Interpret Cq Values?

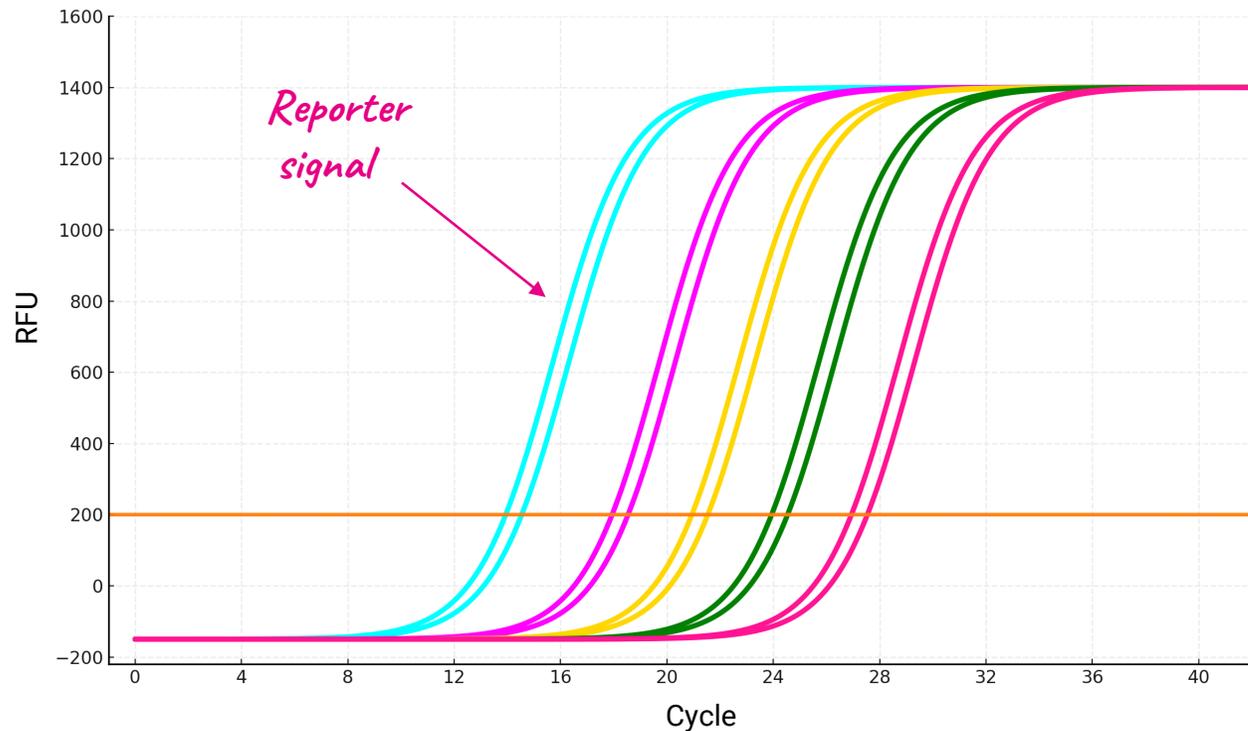


3. What Is My Result And What Does It Mean? How Do I Interpret Cq Values?



Amplification Curve – shows accumulation of product as PCR progresses

- **Reporter** – fluorescent dye or label used to monitor PCR product formation
- **RFU** – relative fluorescence unit

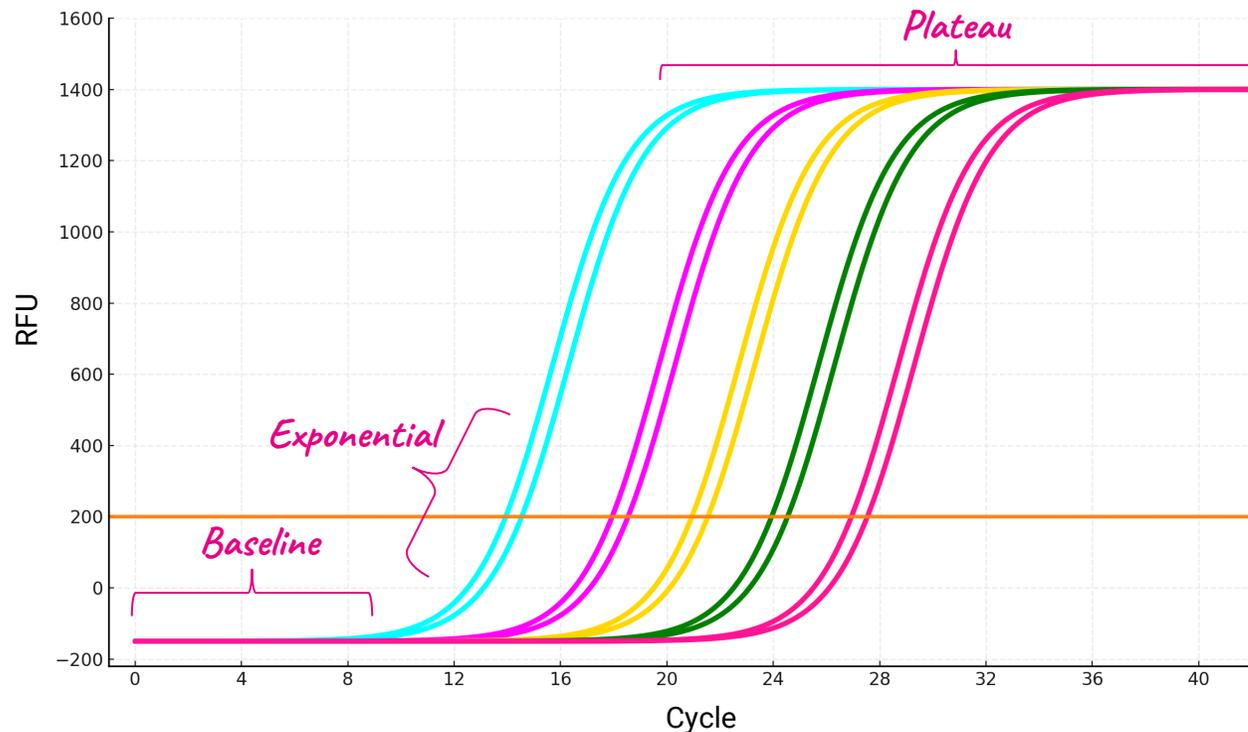




Primary output is the amplification curve

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



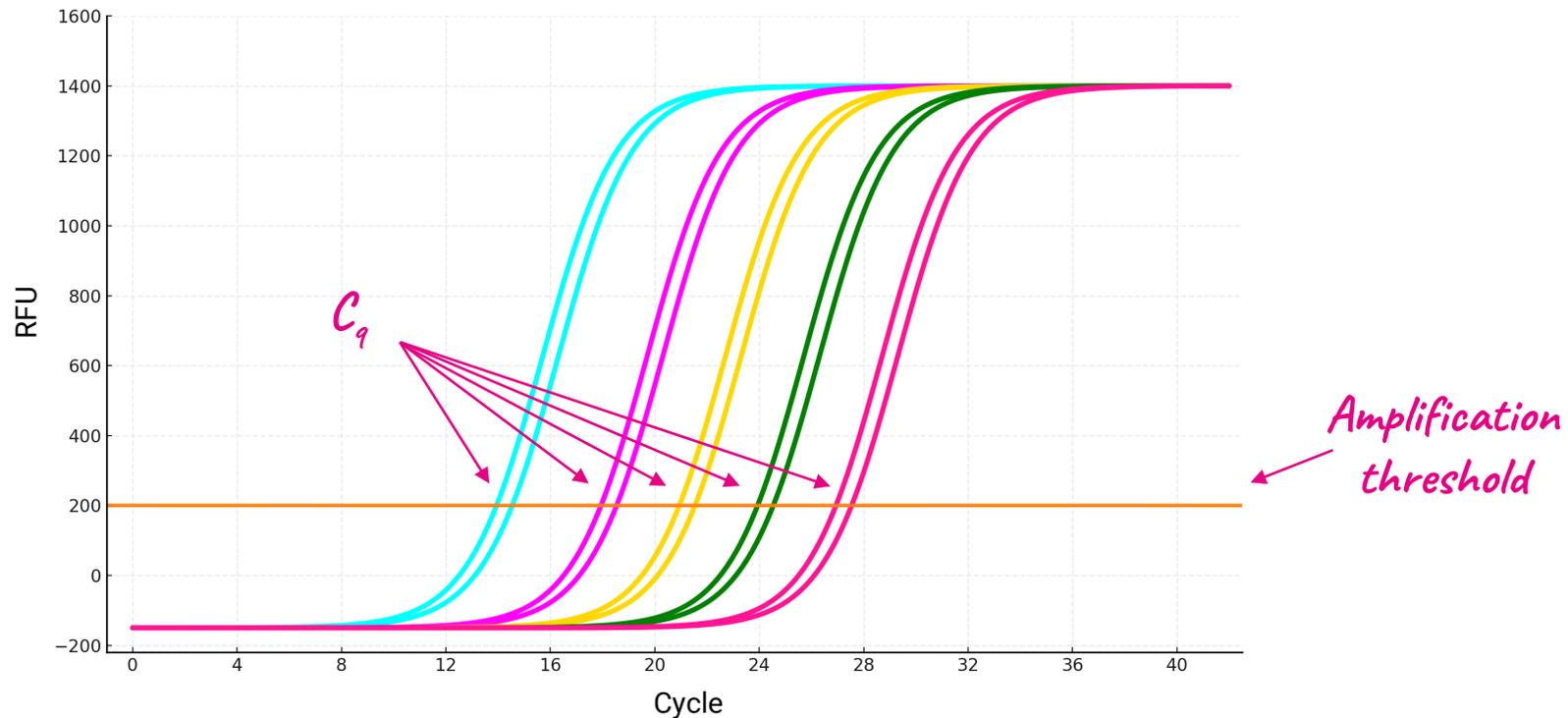


What is a C_q value?

C_q = quantification cycle

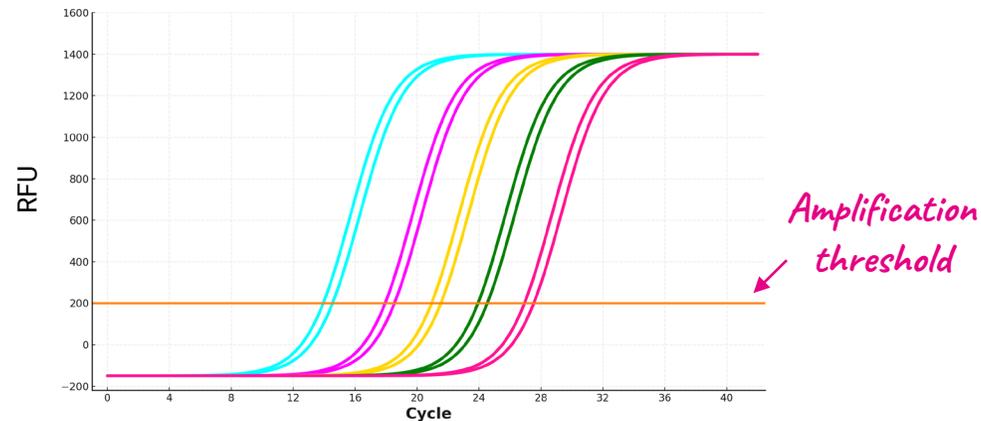
Cycle number at which amplification curve crosses amplification threshold – this is the “take-away” metric...

C_q value is inversely proportional to amount of starting template



What do I do with the Cq value?

Cq values depend on how the threshold is set and are therefore not directly comparable.



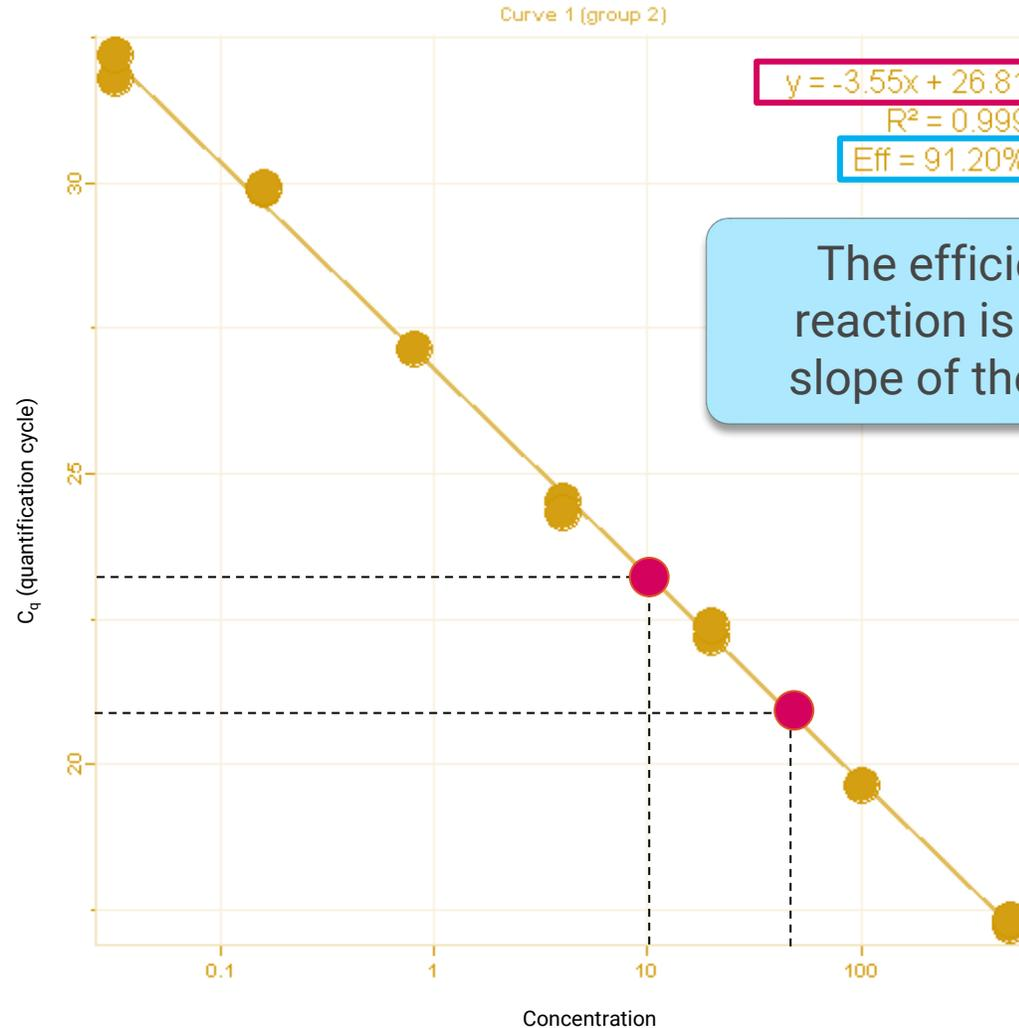
1. **Absolute quantification** = measurement of the expression of a target gene using a standard curve
2. **Relative quantification** = measurement of the expression of a target gene compared to references



Absolute Quantifizierung über die Standardkurve

The standard curve is created by plotting C_q against \log (concentration).

The concentrations of unknown samples are derived from the C_q value using the standard curve.



The efficiency of the PCR reaction is a function of the slope of the standard curve.

Versch modelle

Absolut
relativ



(Normalized) Relative Quantitation

Livak and Schmittgen (2001)

- 100% PCR efficiency
- 1 reference gene

$$NRQ = 2^{\Delta\Delta Cq}$$

$= \Delta C_{q,goi} - \Delta C_{q,ref}$

$$= \frac{2^{\Delta C_{q,goi}}}{2^{\Delta C_{q,ref}}}$$

Pfaffl (2001)

- Experimentell bestimmte PCR efficiency
- 1 reference gene

$$NRQ = \frac{E_{goi}^{\Delta C_{q,goi}}}{E_{ref}^{\Delta C_{q,ref}}}$$

reminder
 $RQ = \frac{\Delta C_q}{E}$
PCR efficiency

qBase model (2007)

- exp PCR efficiency
- Multiple reference genes

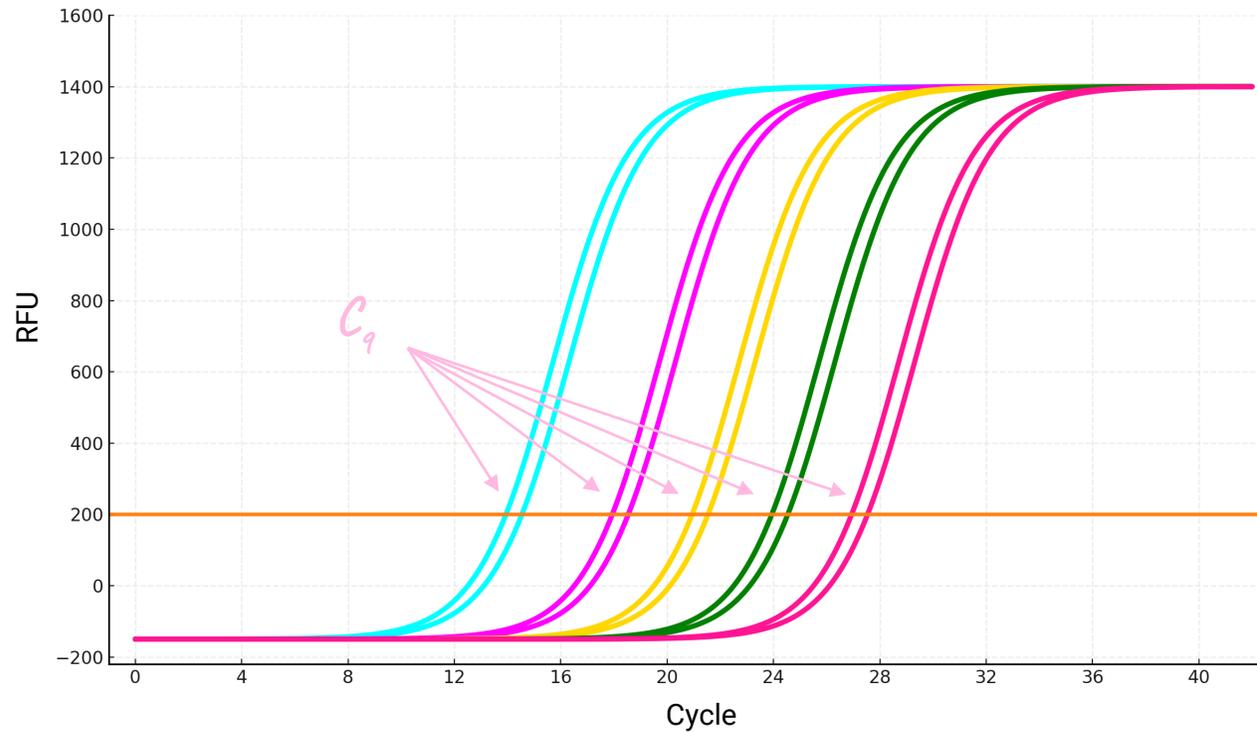
$$NRQ = \frac{E_{goi}^{\Delta C_{q,goi}}}{\sqrt[n]{\prod_i^n E_{ref_i}^{\Delta C_{q,ref_i}}}}$$

geometric mean

Don't worry!

Any real-time software can perform these functions

- Automatically
- User-defined



$$NRQ = \frac{E_{goi}^{\Delta Cq_{goi}}}{\sqrt[n]{\prod_i^n E_{refi}^{\Delta Cq_{refi}}}}$$

Amplification threshold



4. What controls have my back?

- **Positive Controls**

- **Nucleic acid target sequence**

If positive:

- helps determine LOD and LOQ,
- establishes sensitivity and linear dynamic range of the assay
- permits detection of potential inhibitors or suboptimal reaction conditions
- aids in threshold setting

- **Negative Controls**

- **Carrier nucleic acid (not target sequence)**

If positive:

- unknown sample should be considered positive only if its C_q precedes that of the negative control by at least 5 cycles

- **No RT Controls**

- **Sample without reverse transcriptase enzyme (only RT-qPCR)**

If positive:

- gDNA contamination

- **No Template Controls**

- **Water or buffer instead of nucleic acid**

If positive:

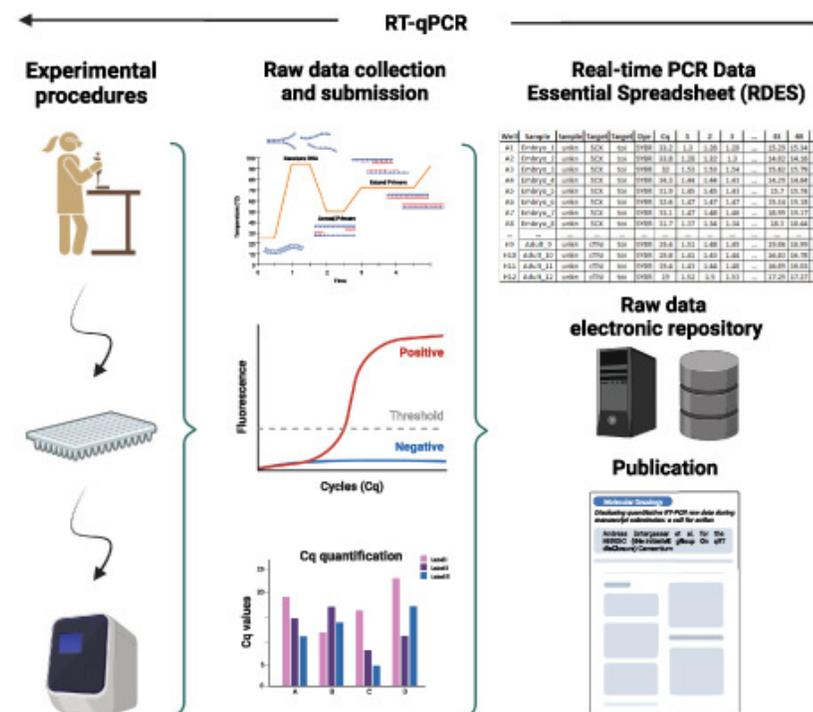
- (Cross-) contamination = results of the entire run are invalid or caution should be taken when interpreting results



5. Can someone else re-run my analysis?

Save your raw data!

- Raw data export
- Thresholding method
- Baseline settings



Adapted from Untergasser *et al.*, 2023

MIQE suggests to publish raw data in the supplementary material as RDES or RDML files



Für jede Probe der richtige Mix!

Product	GoTaq [®] qPCR	GoTaq [®] Probe	GoTaq [®] Enviro	GoTaq [®] Endure
Method of detection	Dye-based	Probe-based	Probe-based	Probe-based
Suitable for	Almost all types of samples	Almost all types of samples	Environmental samples	Samples with a lot of inhibitors
Inhibitor resistance	✓	✓	✓ ✓ ✓	✓ ✓ ✓



Probe-based GoTaq® Endure Master Mix

Inhibitors like EDTA or humic acid can distort your qPCR results!

If you have samples with many inhibitors that are difficult to remove during sample preparation, use a master mix that is resistant to inhibitors

Request a free sample!

www.promega.com/testEndure

GoTaq® Endure qPCR & RT-qPCR Master Mixes

Suitable for a variety of sample types:

- ✓ Blood
- ✓ Bacteria
- ✓ Virus
- ✓ Feces
- ✓ Soil
- ✓ Plant
- ✓ Food



Get your free sample now!

 Probe-based

 Extremely high inhibitor tolerance

 Fast Cycling

 Multiplexing

History of



1978

Bill Linton begins selling restriction enzymes at UW-Madison



1980

Promega moves operations to current location in Fitchburg, Wisconsin

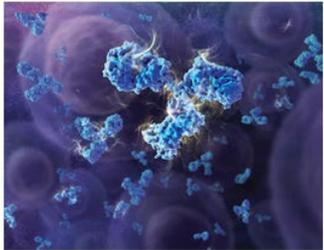


1983

Promega opens first international branch offices

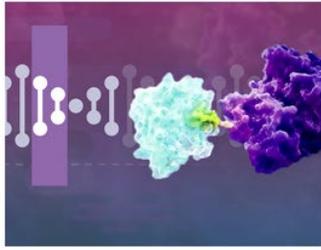
Learning @Promega

Popular Webinars



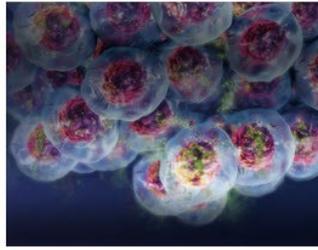
Antibody Internalization Assay

Learn how a pH sensor fluorescent dye used to identify antibodies suitable for receptor mediated internalization.



A Guide to CRISPR Mediated Gene Tagging

Learn about a simple and efficient method for CRISPR-mediated HIBIT tagging that requires no molecular cloning.



Overview of 3D Cell Culture Model Systems

Factors to consider when choosing and validating cell-based assays for use with 3D cultures.



Promega-Academy

Wissenschaftliche Seminare: Zugeschnitten auf Ihre Bedürfnisse!

Promega Connections

Thoughts, tech tips and news about science

Nukleinsäure-Analyse

Automatisierte DNA & RNA Extraktion aus jedem Probenotyp



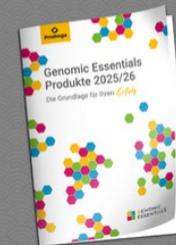
Nukleinsäure-Aufreinigung: modular und flexibel



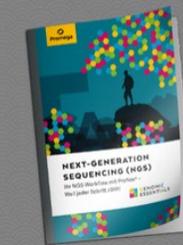
Manuelle DNA- & RNA-Aufreinigung



Genomics-Produkte



Next Generation Sequencing



Fluorometer zur Nukleinsäurequantifizierung



DNA-Sequenzierung & Fragmentanalyse



CE-Instrument für forensische STR-Analyse



Humane Identifizierung



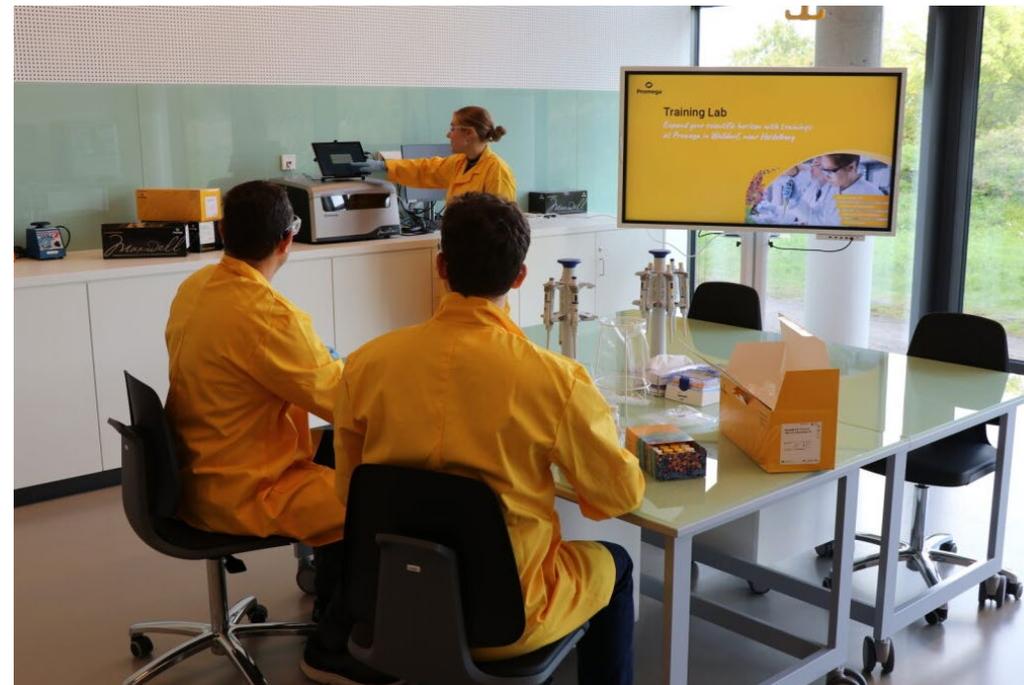
Technical Service

Visit our

- **qPCR Seminar**
- **qPCR Workshop**

Or watch our **qPCR Webinar!**

If you need help with your (PCR) reactions, feel free to contact our technical service!



Dr. Hannah Buntz

Manager Training Lab

Tel. +49 151 580 421 05

hannah.buntz@promega.com



Technical Services Team



Christof
Brückmann



Sabine
Wawro



Petra
Numrich



Thorsten
Stehlik



Jan
Adam

Contact us in case of any questions or problems

de_techserv@promega.com

+49 6227 6906 290

Request your
sample!

Thank you!

Feel free to ask questions!

 Let's connect!

Contact your local sales representative for
products & sales relevant information:
www.promega.com/c/local_sales/sales_contacts/

