

# How NOT to mess up your qPCR

Promega GmbH



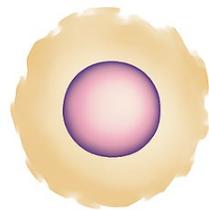
# Stimulus-Triggered Acquisition of Pluripotency (STAP)

ARTICLE

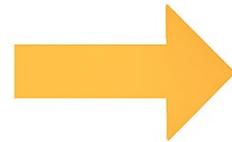
doi:10.1038/nature12968

## Stimulus-triggered fate conversion of somatic cells into pluripotency

Haruko Obokata<sup>1,2,3</sup>, Teruhiko Wakayama<sup>3†</sup>, Yoshiki Sasai<sup>4</sup>, Koji Kojima<sup>1</sup>, Martin P. Vacanti<sup>1,5</sup>, Hitoshi Niwa<sup>6</sup>, Masayuki Yamato<sup>7</sup> & Charles A. Vacanti<sup>1</sup>



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 <sup>a,\*</sup> Jan M. Ruijter,<sup>b</sup> Maurice J.B. van den Hoff,<sup>b</sup> Mikael Kubista,<sup>c</sup> Michael W. Pfaffl,<sup>d</sup>  
Gregory L. Shipley,<sup>e</sup> Nham Tran,<sup>f</sup> Stefan Rödiger,<sup>g</sup> Andreas Untergasser,<sup>h</sup> Reinhold Mueller,<sup>i</sup> Tania Nolan,<sup>j</sup>  
Mojca Milavec,<sup>k</sup> Malcolm J. Burns,<sup>l</sup> Jim F. Huggett <sup>l</sup> Jo Vandesompele <sup>m</sup> and Carl T. Wittwer<sup>n</sup>

---

# What is Real-Time PCR?

## End-point PCR

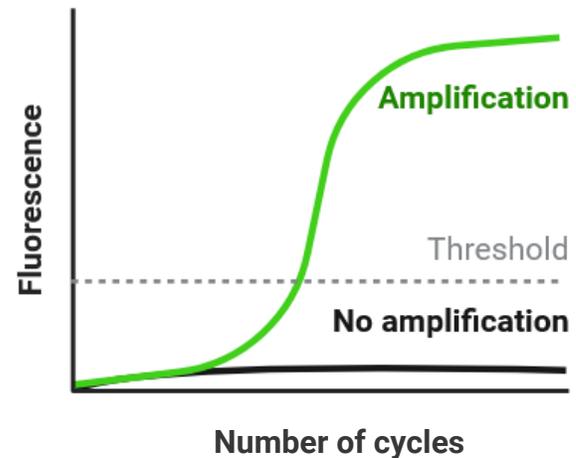
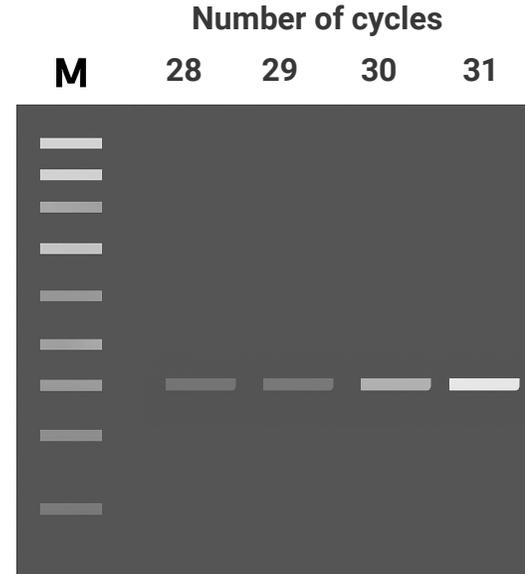
Product formation measured

after reaction is complete

## Real-Time PCR

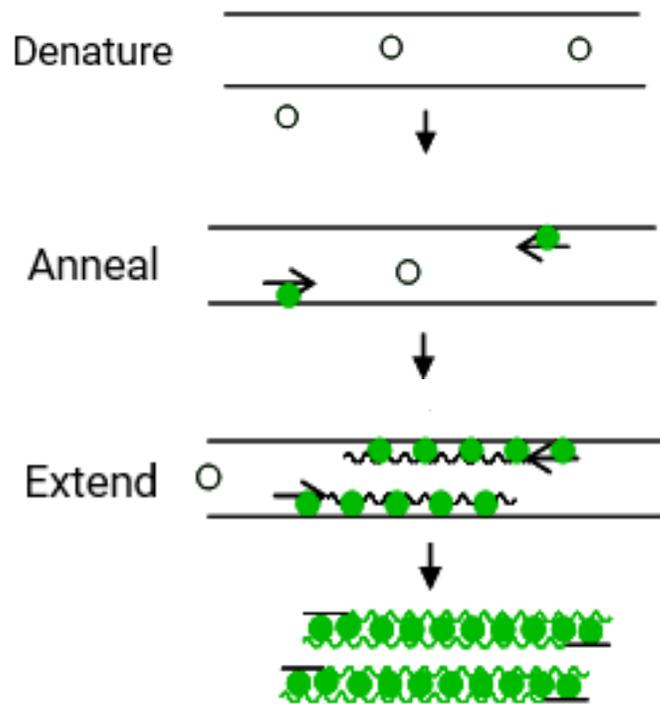
Product formation measured

at each cycle, during the reaction

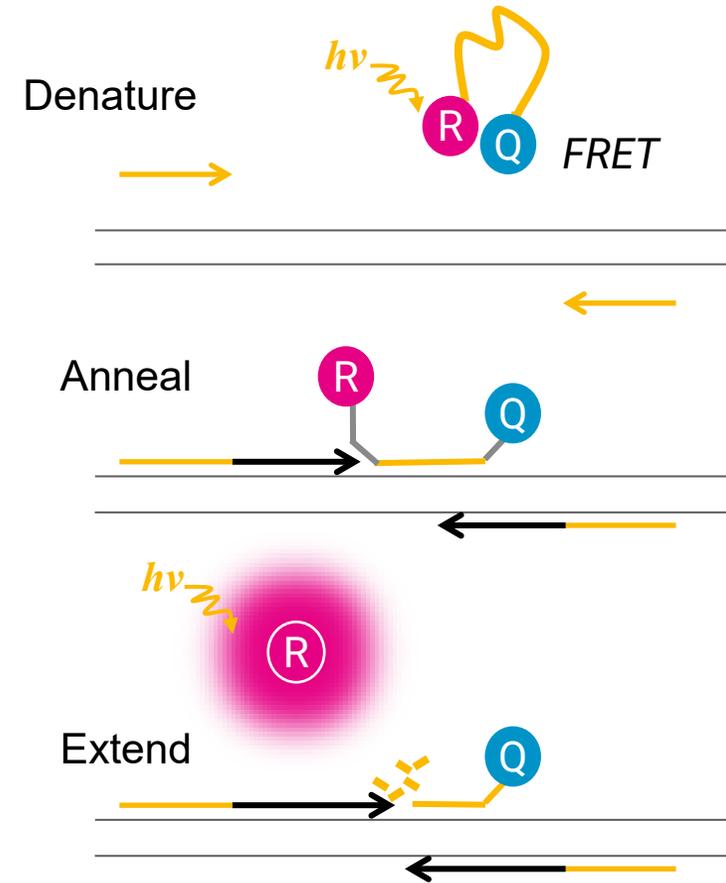


# Real-time PCR Chemistries: **Dye-based** vs. **Probe-based** qPCR

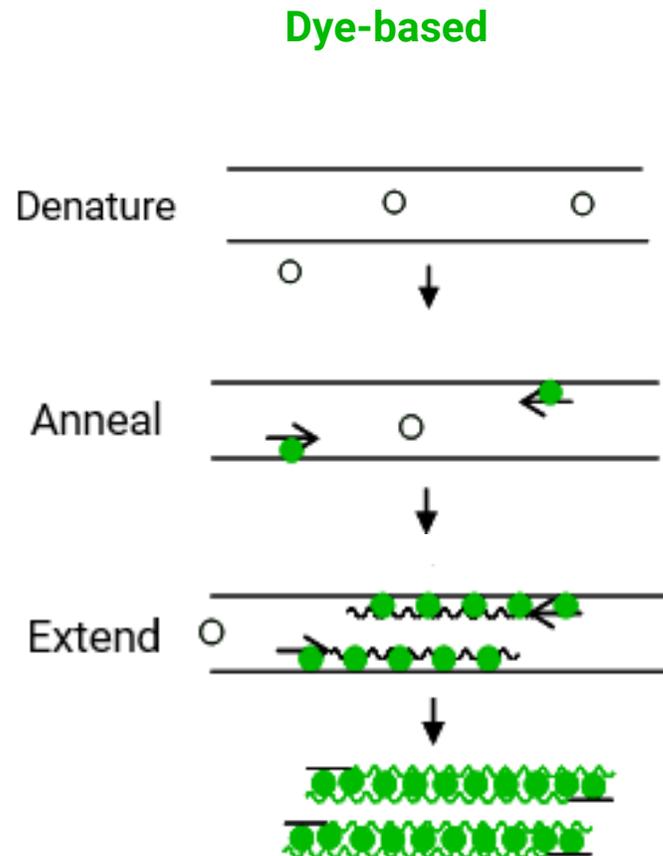
## Dye-based



## 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

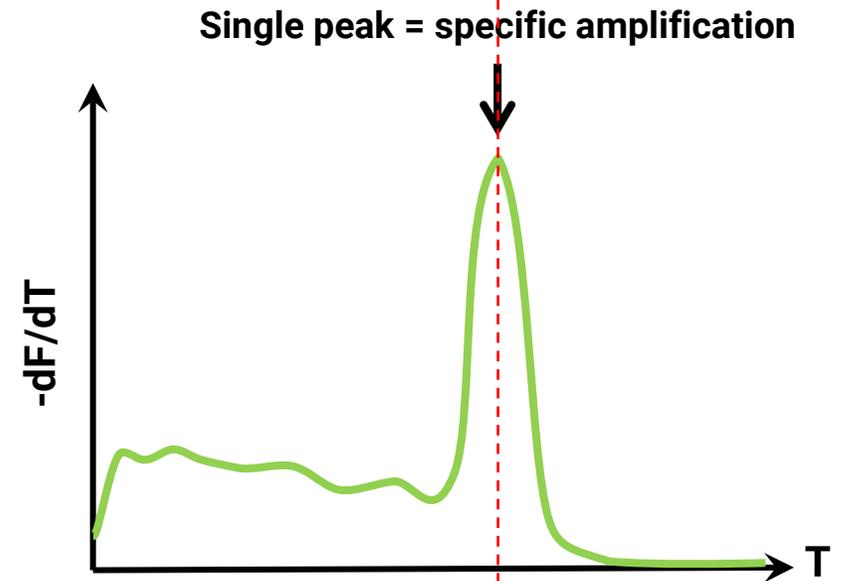
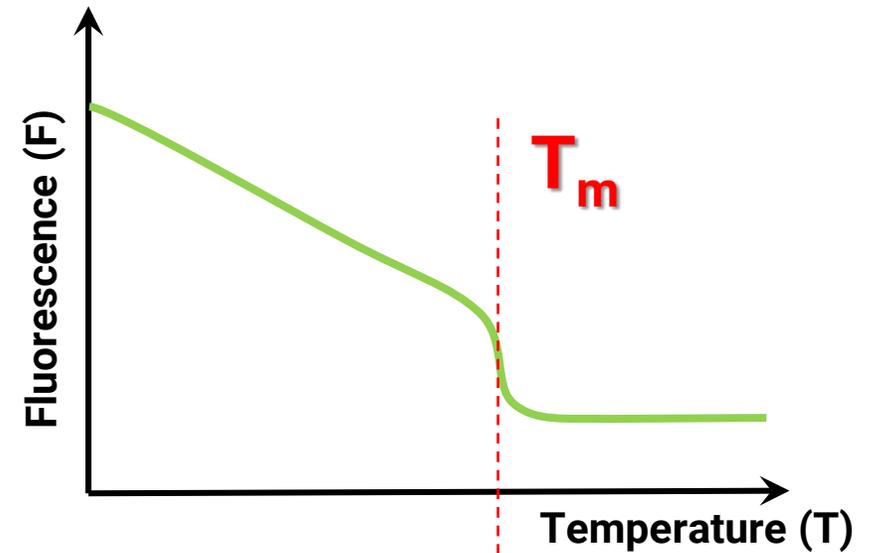
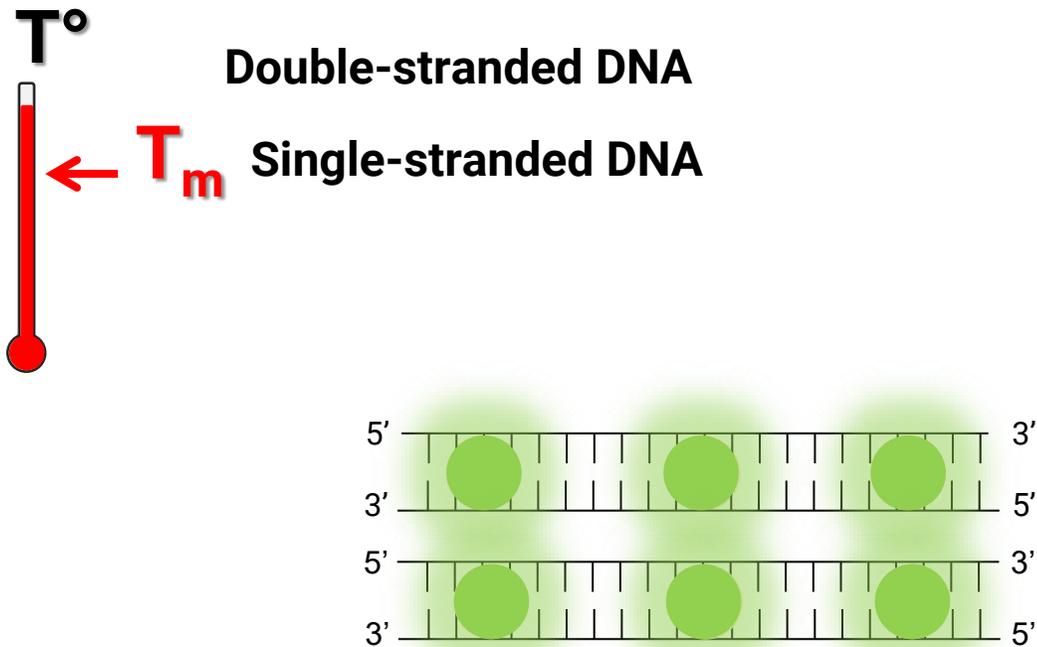
- Free Dye → low fluorescence
- Bound Dye → high fluorescence

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

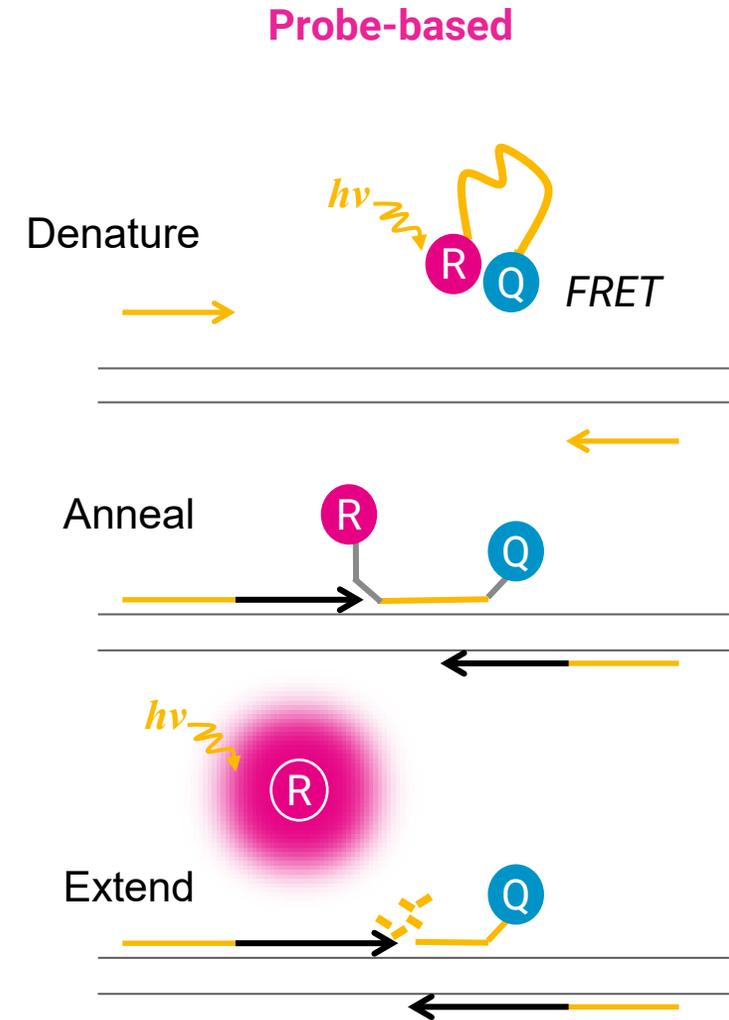
**TaqMan<sup>®</sup> is the most familiar type:**

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

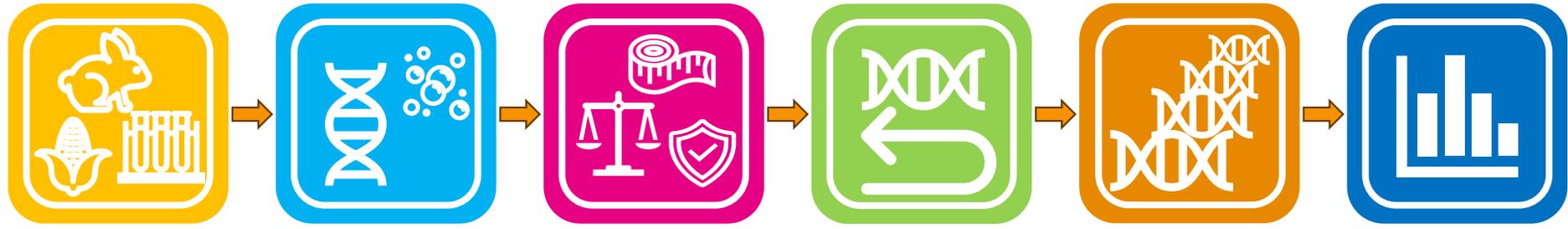
**Free probe → FRET occurs**

**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

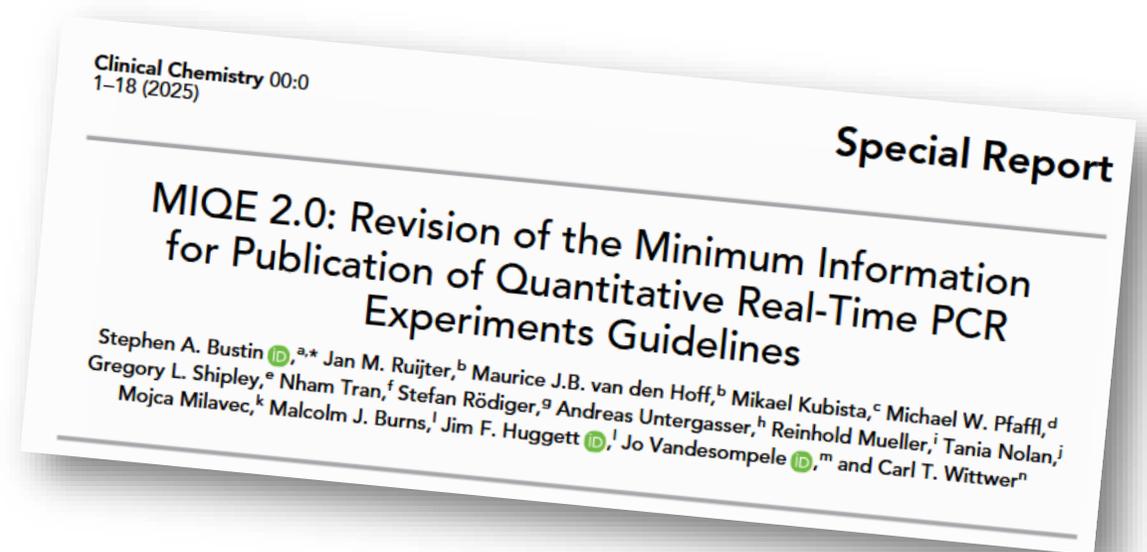


# Ask Yourself 5 Questions Before You Publish Your qPCR Results

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

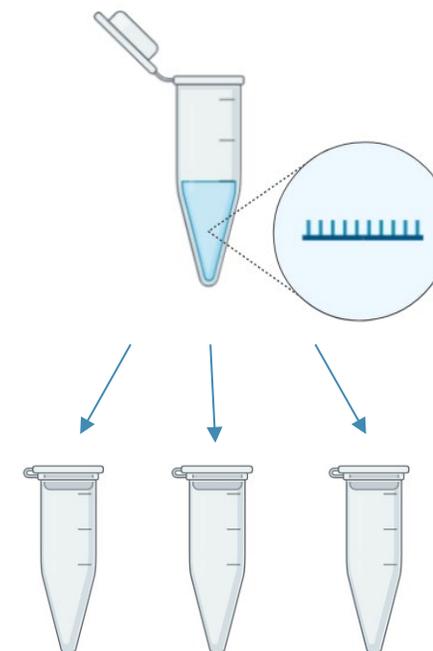
# 1. Can I Trust This Sample As A Valid Starting Point For My Quantification?



## How do I store my sample correctly?



- Extracted RNA:
  - 20°C → days,
  - 80°C → months,
  - 180°C (liquid N<sub>2</sub>) → 5 years
- Store DNA/RNA in small aliquots
- Avoid multiple freeze-thaw cycles
- Check nucleic acid concentration after each thawing



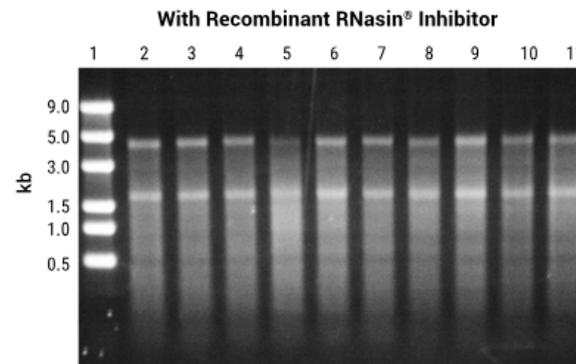
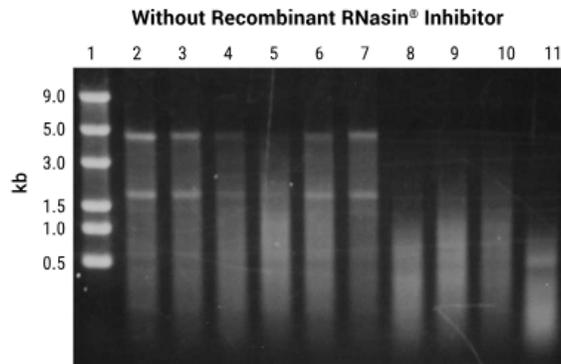


# 1. Can I trust this sample as a valid starting point for my quantification?

Is my sample intact or degraded?

## Electrophoresis

- Size
- Integrity
- Quantity



## Dye Staining (QuantiFluor)

- Concentration (very sensitive)

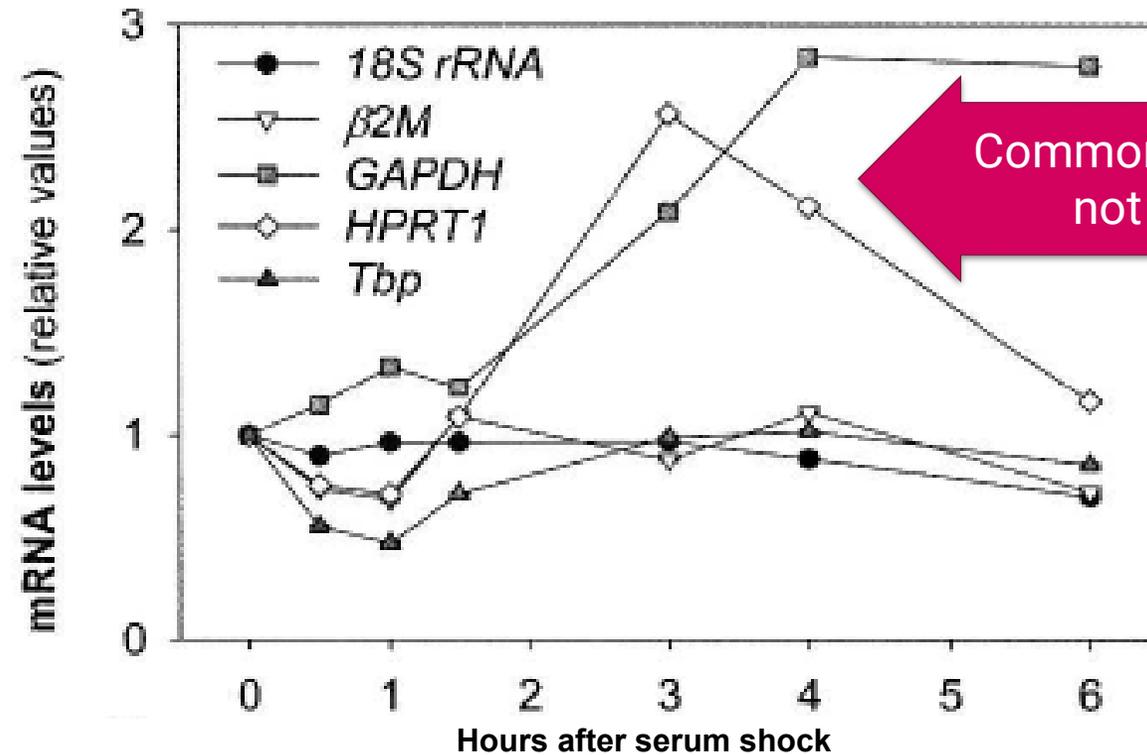


MIQE suggests **combination of electrophoresis & dye staining** to assess **size, quantity, concentration and integrity** (although this does not reveal contamination)

*Quality is more important than yield!*

## 2. Did I Normalize the Right Way?

### Reference genes

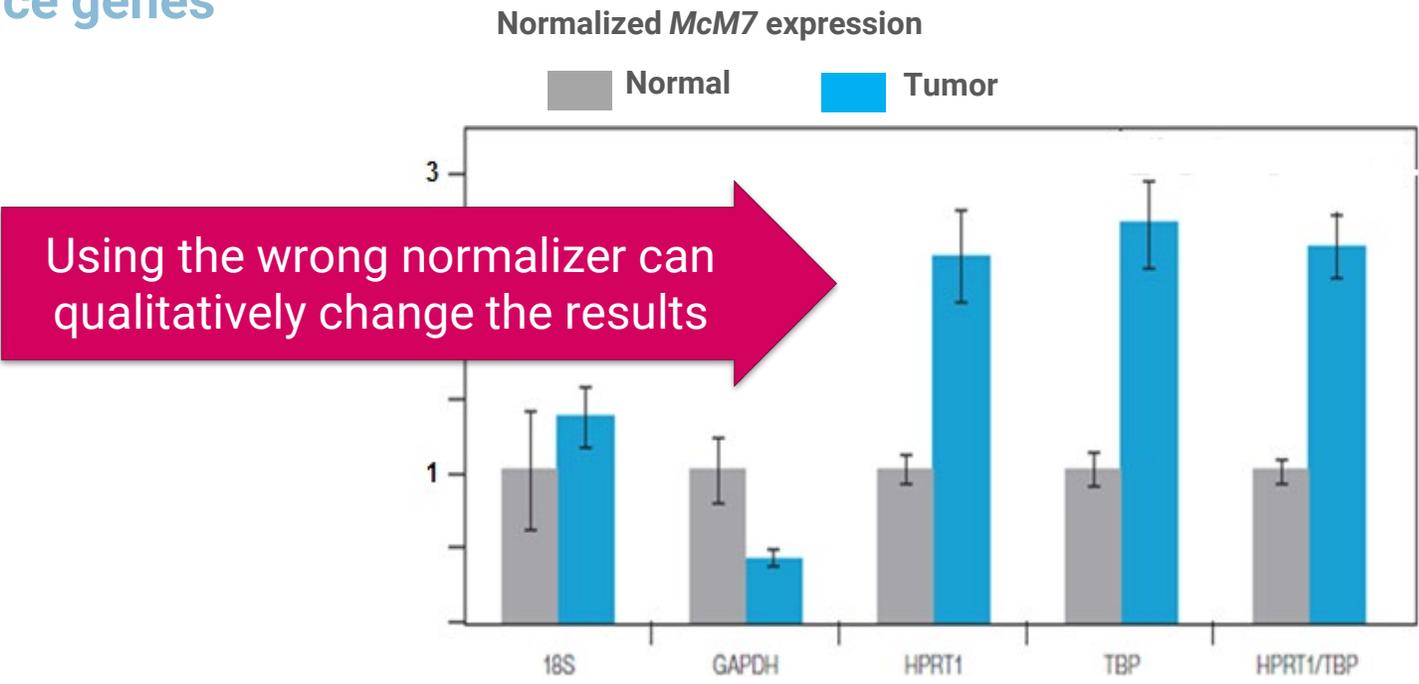


Commonly used normalizers are not always constitutive

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

## 2. Did I Normalize the Right Way?

### Reference genes

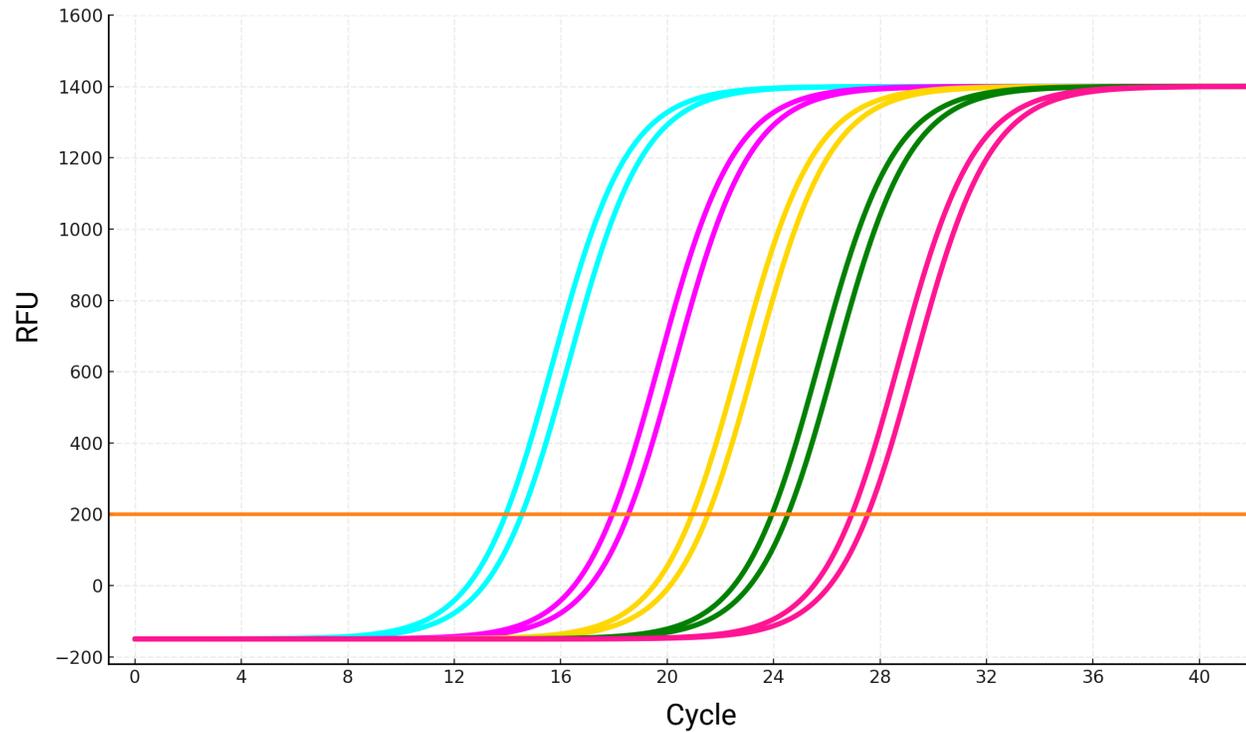


Using the wrong normalizer can qualitatively change the results

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?

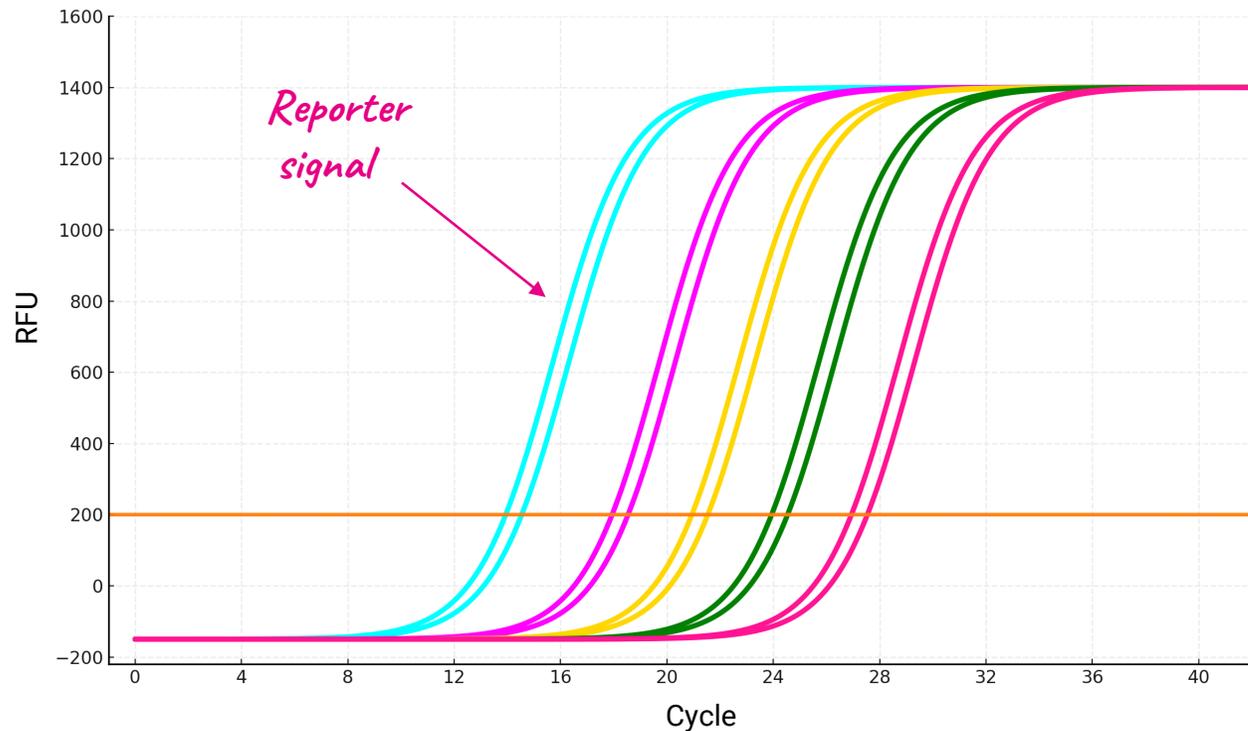


### 3. What Is My Result And What Does It Mean? How Do I Interpret Cq Values?



**Amplification Curve** – shows accumulation of product as PCR progresses

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

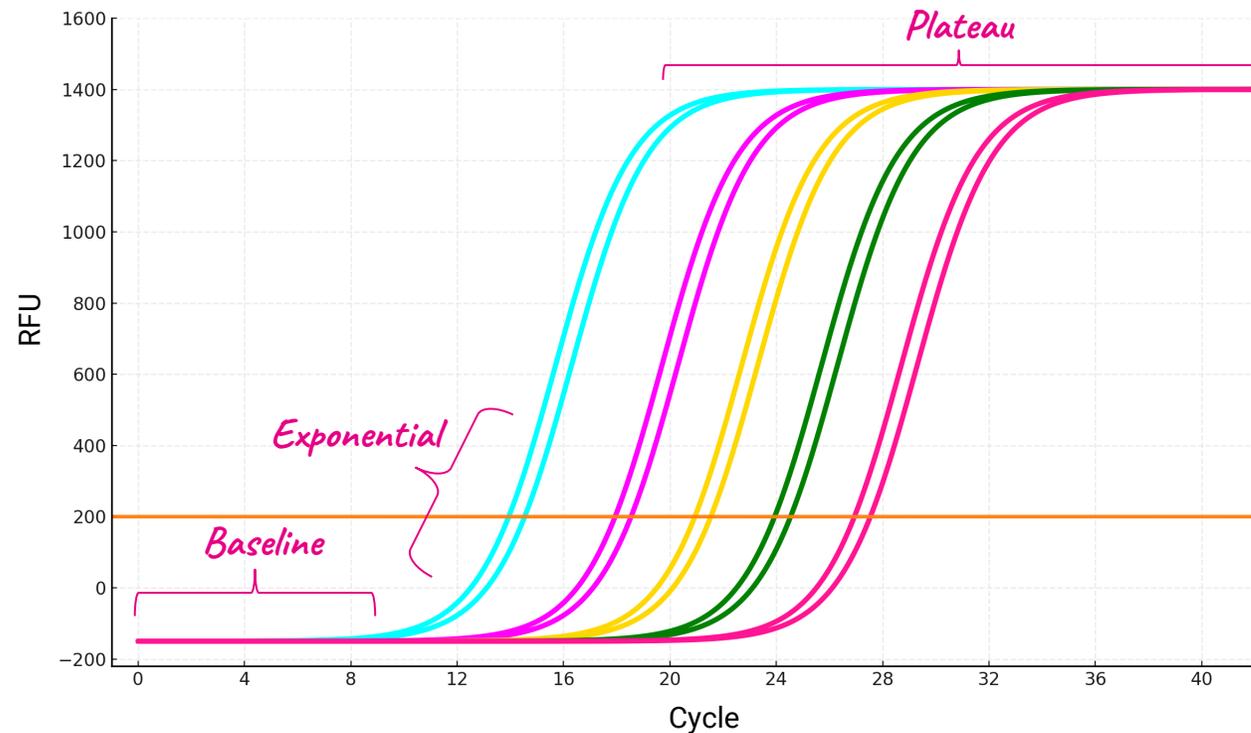




# Primary output is the amplification curve

**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



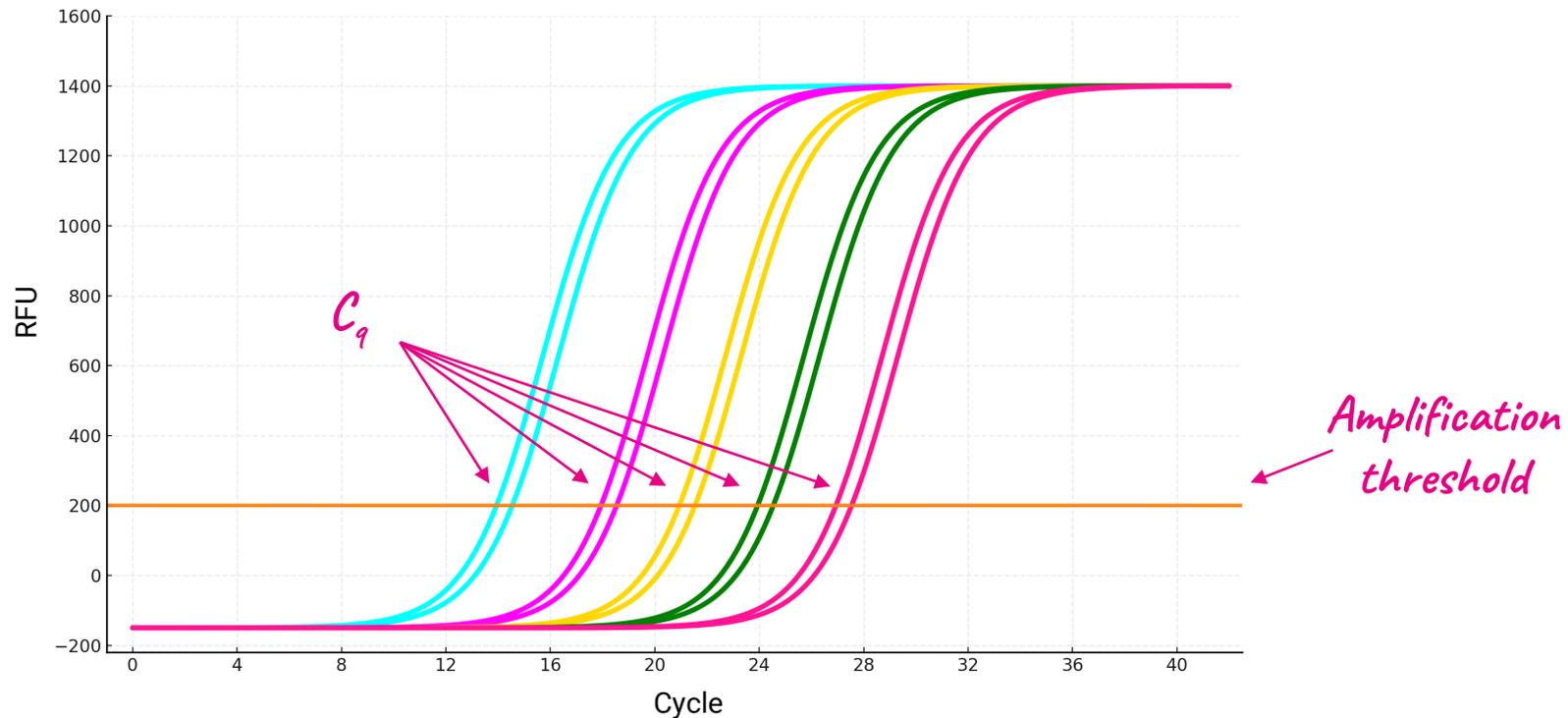


# What is a $C_q$ value?

$C_q$  = quantification cycle

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

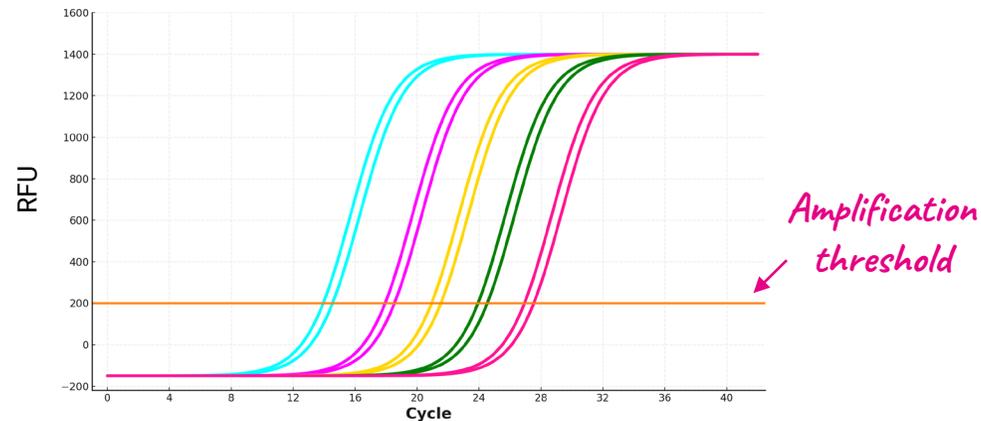
$C_q$  value is inversely proportional to amount of starting template



# What do I do with the Cq value?

---

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



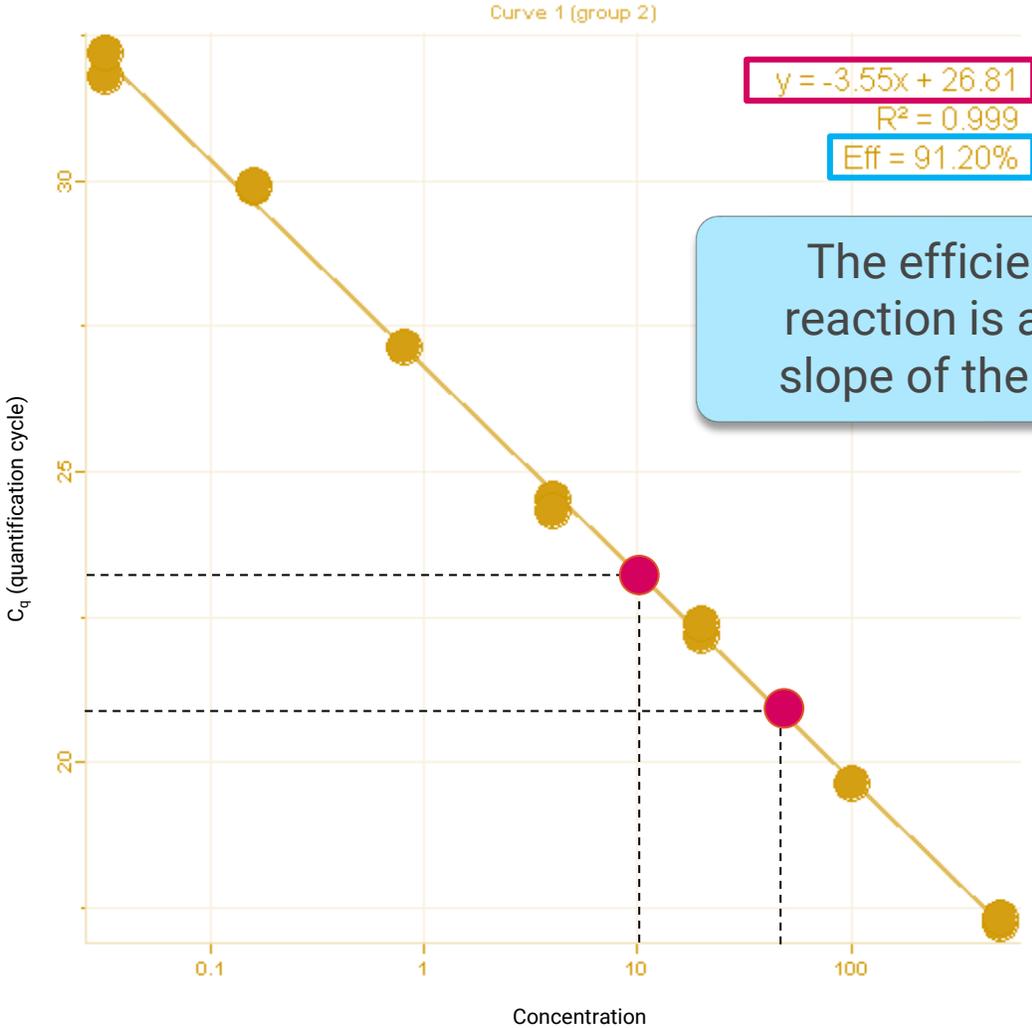
1. **Absolute quantification** = measurement of the expression of a target gene using a standard curve
2. **Relative quantification** = measurement of the expression of a target gene compared to references



# Absolute Quantifizierung über die Standardkurve

The standard curve is created by plotting C<sub>q</sub> against log (concentration).

The concentrations of unknown samples are derived from the C<sub>q</sub> 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

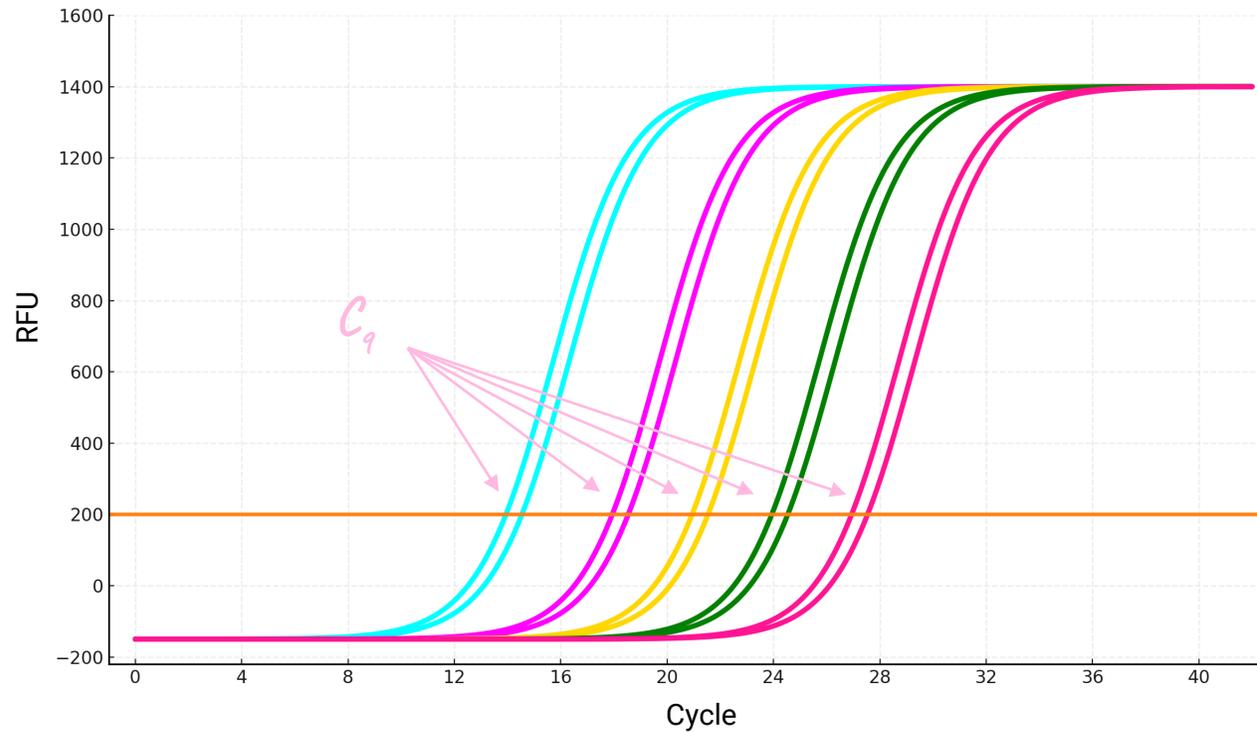
$$NRQ = \frac{E_{goi}^{\Delta C_{q,goi}}}{\sqrt[n]{\prod_i^n E_{ref_i}^{\Delta C_{q,ref_i}}}}$$

*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<sub>q</sub> 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**

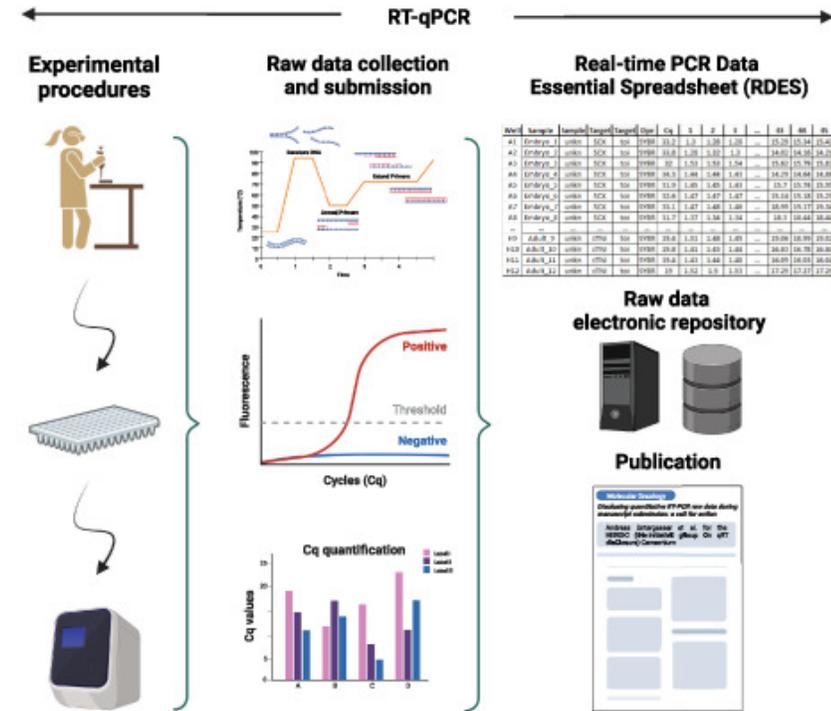
If positive:

- (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 <sup>®</sup> qPCR	GoTaq <sup>®</sup> Probe	GoTaq <sup>®</sup> Enviro	GoTaq <sup>®</sup> 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](http://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

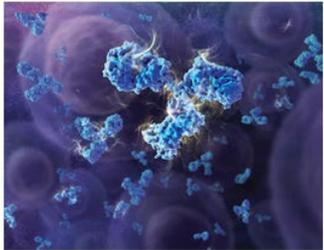


**1983**

Promega opens first international branch offices

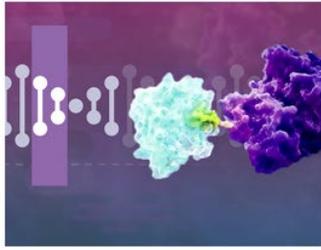
# Learning @Promega

## Popular Webinars



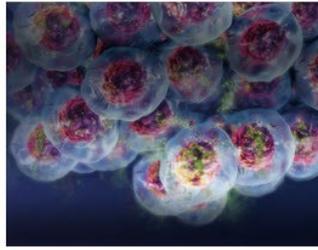
### Antibody Internalization Assay

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: Zugespitzt auf Ihre Bedürfnisse!

## Promega Connections

Thoughts, tech tips and news about science

## Nukleinsäure-Analyse

Automatisierte DNA & RNA Extraktion aus jedem Probenotyp



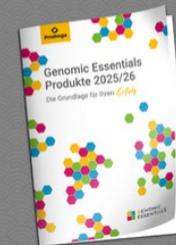
Nukleinsäure-Aufreinigung: modular und flexibel



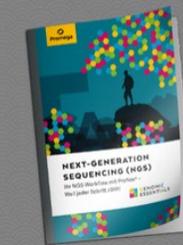
Manuelle DNA- & RNA-Aufreinigung



Genomics-Produkte



Next Generation Sequencing



Fluorometer zur Nukleinsäurequantifizierung



DNA-Sequenzierung & Fragmentanalyse



CE-Instrument für forensische STR-Analyse



Humane Identifizierung



# Technical Service

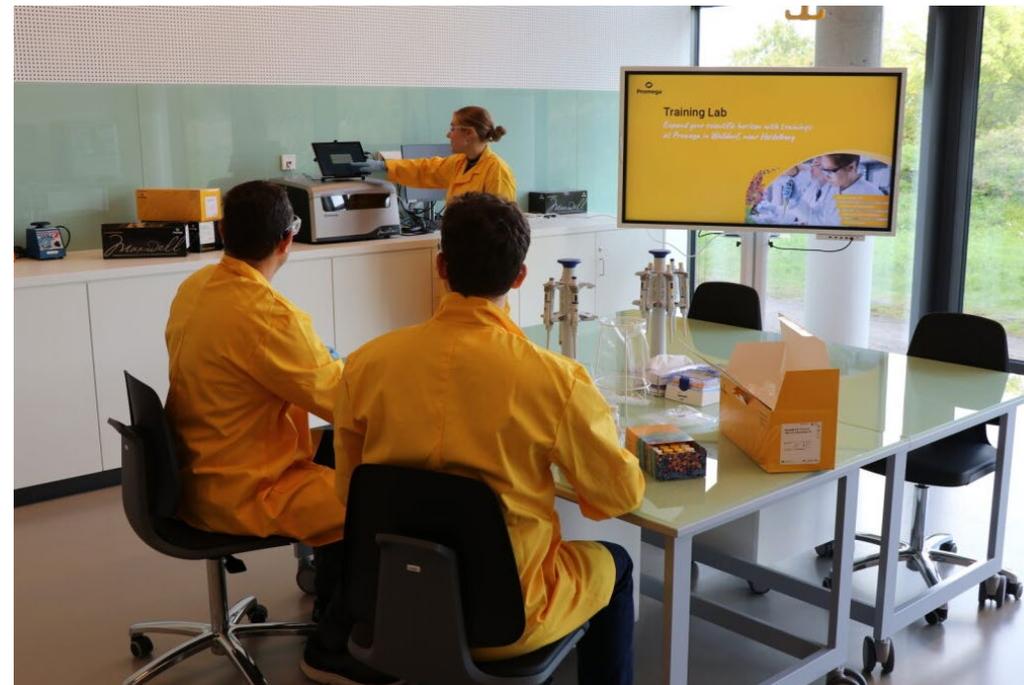
---

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](mailto: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](mailto: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/](http://www.promega.com/c/local_sales/sales_contacts/)

