



Challenging biological samples: stabilization and simultaneous purification of DNA and RNA

Stefanie Schröer

1 October 10, 2012

■ Requirements for analysis of precious samples

- Precious and inhomogeneous samples

- Different analytes from the same sample

■ Stabilization of RNA and DNA

■ Simultaneous purification of different analytes

- Simultaneous analysis of DNA, RNA, and protein

- DNA, RNA, and miRNA from challenging starting material

- Analysis of DNA and RNA from FFPE samples

■ Downstream applications

■ Summary



Precious and nonhomogeneous samples

- Clinical samples that are irreplaceable
- Samples of limited quantity
- Tumor samples that often nonhomogeneous
- Highly variable starting material (e.g., different tissue types)
- Samples from biobanks



- Requirements for analysis of precious samples
 - Reliable comparison of mRNA/miRNA expression profiles
 - Reliable comparison of methylation status of samples
 - Different analytes from the same sample

- Stabilization of RNA and DNA
 - Stabilization is required in biomedical research
 - QIAGEN offers *RNA/later* and Allprotect products

- ▶ Need for reliable stabilization and parallel purification of nucleic acids from even the most challenging samples

- Requirements for analysis of precious samples

- Precious and inhomogeneous samples

- Different analytes from the same sample

- **Stabilization of RNA and DNA**

- Simultaneous purification of different analytes

- Simultaneous analysis of DNA, RNA, and protein

- gDNA, RNA, and miRNA from challenging starting material

- Analysis of DNA and RNA from FFPE samples

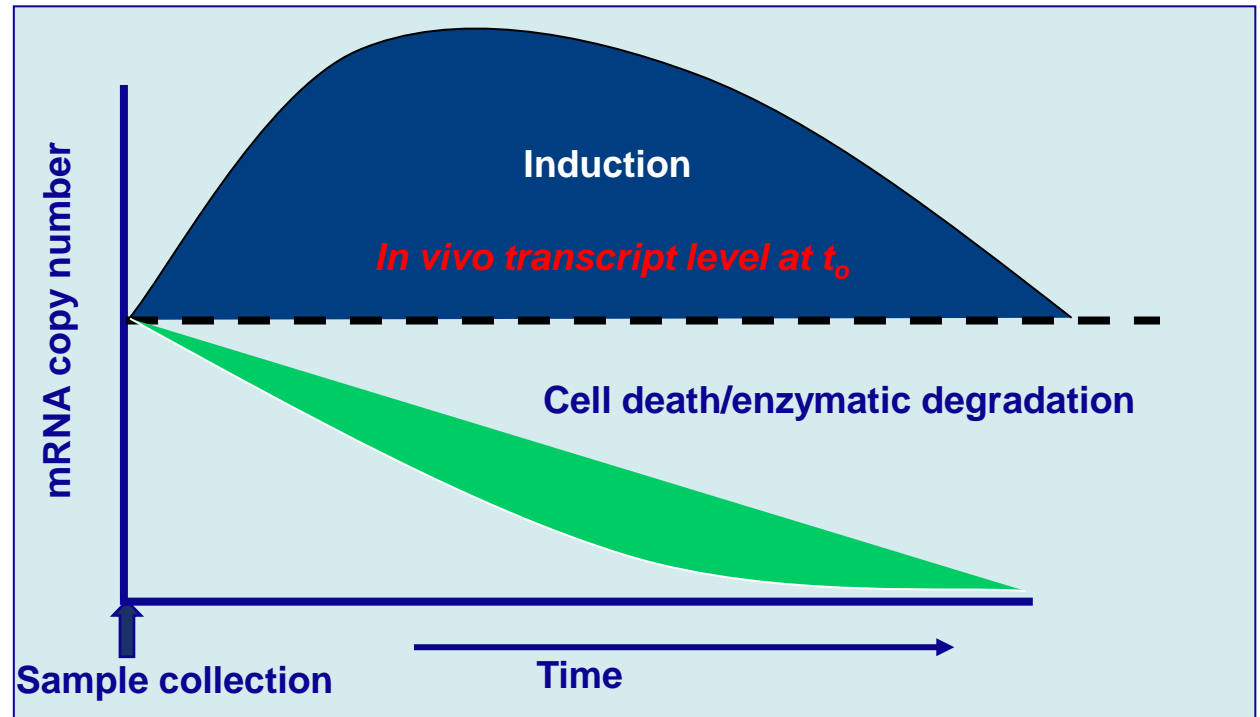
- Downstream applications

- Summary

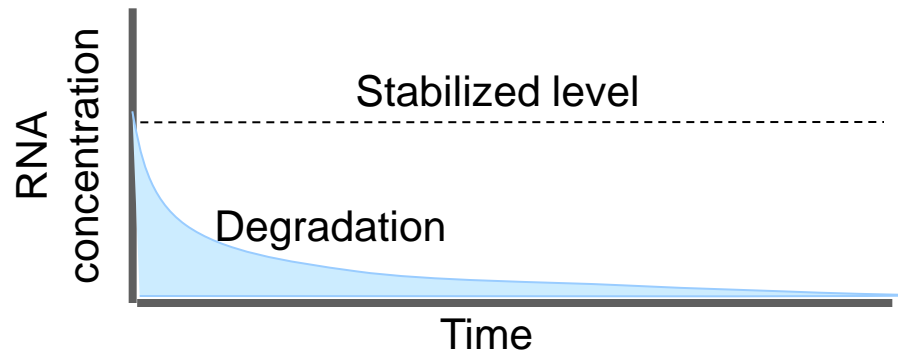
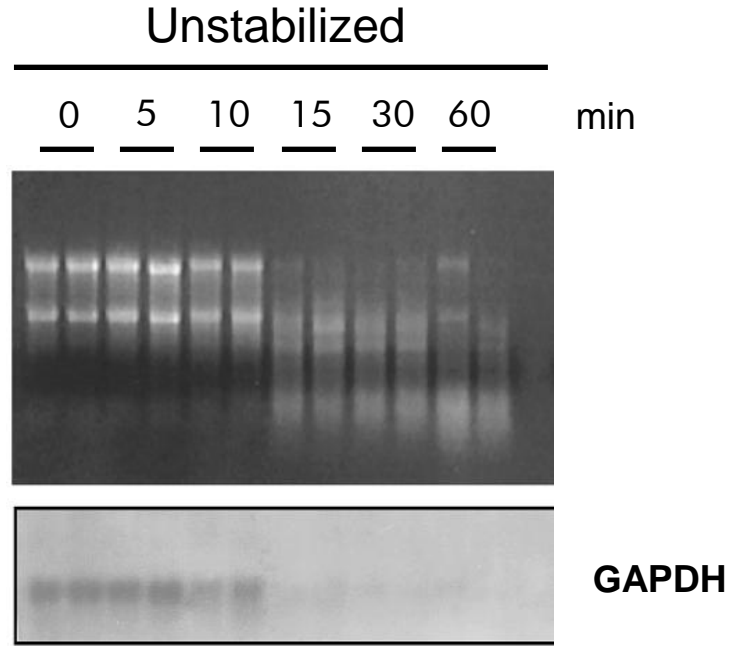
Stabilization of RNA and DNA

Stabilization:

- Protection of RNA from degradation
- Prevention of induction or downregulation of mRNA
- Prevention of degradation of genomic DNA




Degradation of RNA without stabilization



Guidelines for handling RNA

- Work in an RNase-free environment
- Always wear latex or vinyl gloves
- Only use RNase-free water and buffers
- Only use new tips
- Use sterile, disposable plasticware

 Prevent introduction of RNases into the RNA sample during or after isolation procedure

Stabilization reagents:

- Depends on analyte (DNA, RNA, protein)
- Depends on sample type

Products:

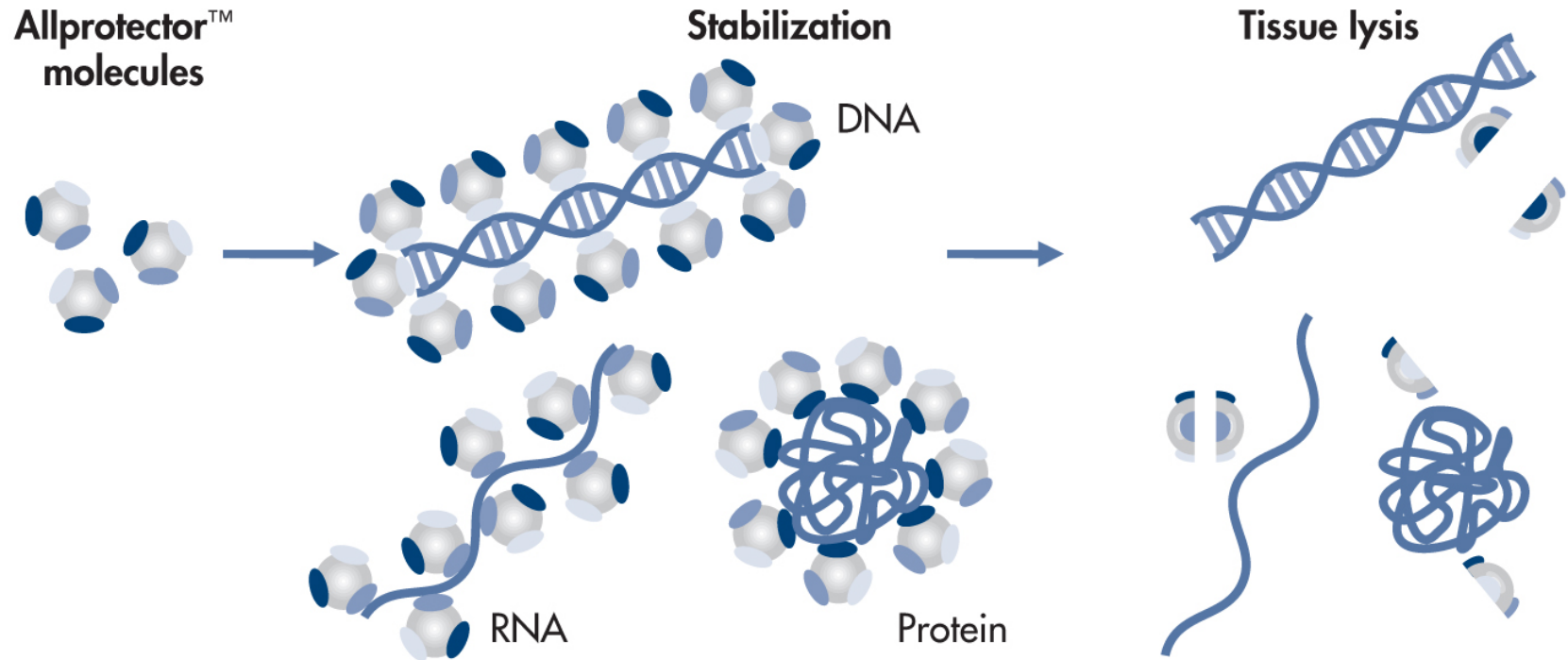
- Tissue: Allprotect Tissue Reagent (DNA/RNA/protein),
RNA_{later} RNA Stabilization Reagent (RNA)
- Cells: RNAprotect Cell Reagent
- Bacteria: RNAprotect Bacteria Reagent
- Saliva: RNeasy Protect Saliva Mini Kit
- Animal blood: RNeasy Protect Animal Blood System

Allprotect Tissue Reagent ensures simultaneous stabilization of DNA, RNA, miRNA and proteins in tissue samples

- Immediate, convenient stabilization of harvested tissue
- Streamlined workflow with no need for liquid nitrogen or phenol
- No toxicity
- No degradation, induction, or modification
- Storage up to 1 day at 37°C, 7 days at 25°C, and 1 year at 2–8°C
- Long-term storage of samples at –20°C or –80°C
- Conveniently used in combination with AllPrep Kits



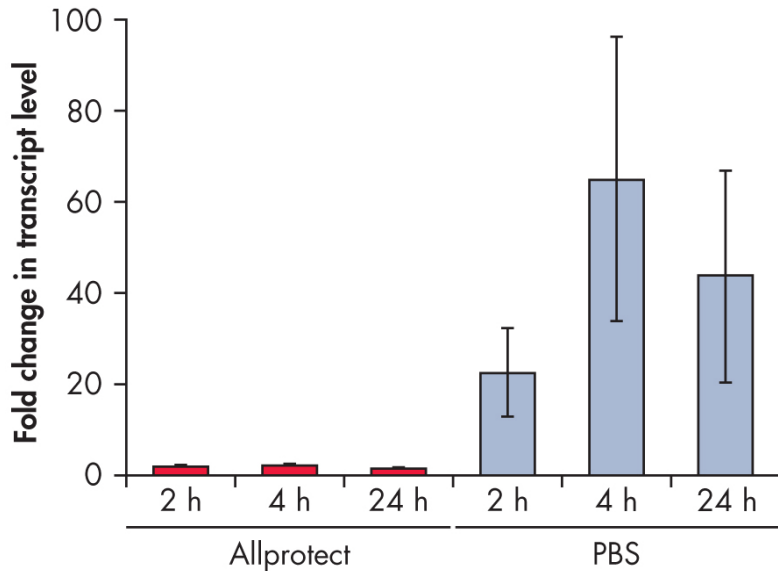
Principle: Allprotector molecules



- Allprotector molecules with contact points to envelop and protect DNA, RNA, and protein
- Upon lysis, Allprotector molecules diluted and release of DNA, RNA, and protein

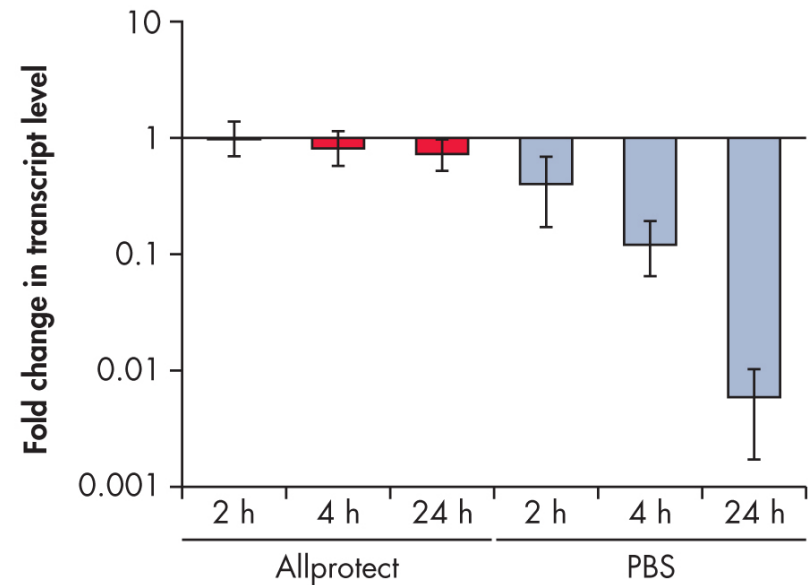
$\Delta\Delta C_T$ analysis of transcriptional induction

Transcript levels of c-fos relative to N₂



➤ Prevention of induction

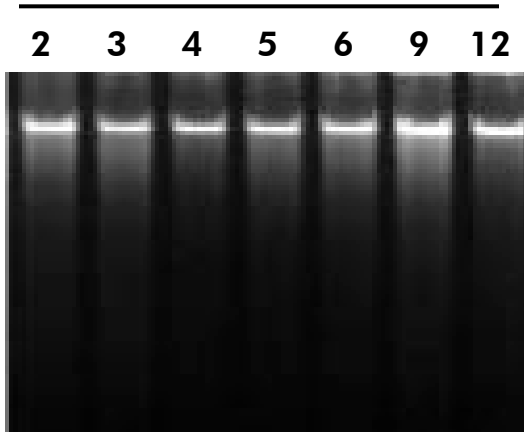
Transcript levels of Madh7 relative to N₂



➤ Prevention of degradation

Genomic DNA

Storage (months)



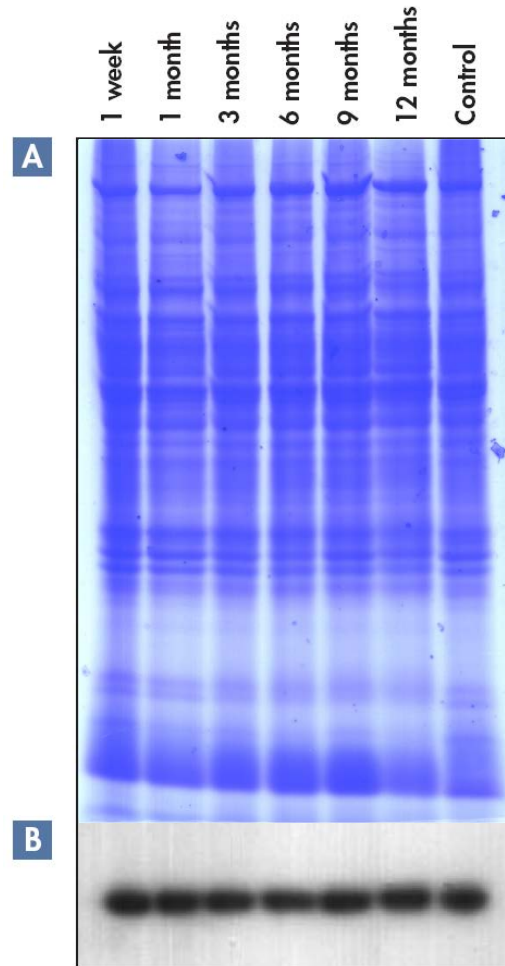
- Rat intestine tissue (10 mg) analyzed
- EZ1 RNA Tissue Mini Kit on the BioRobot EZ1
- Agarose gel

RNA

Storage (months)	RIN value		
	Liver	Intestine	Brain
1	9.4	9.3	8.5
2	9.2	9.3	8.6
4	9.4	9.3	8.4
6	9.2	9.5	8.6
9	9.4	9.7	8.2
12	9.4	9.3	8.4

- Rat tissue (10 mg) analyzed
- RNeasy, RNeasy Lipid Tissue
- Agilent BioAnalyzer

▶ Stabilization of nucleic acids for up to 12 months without freezing

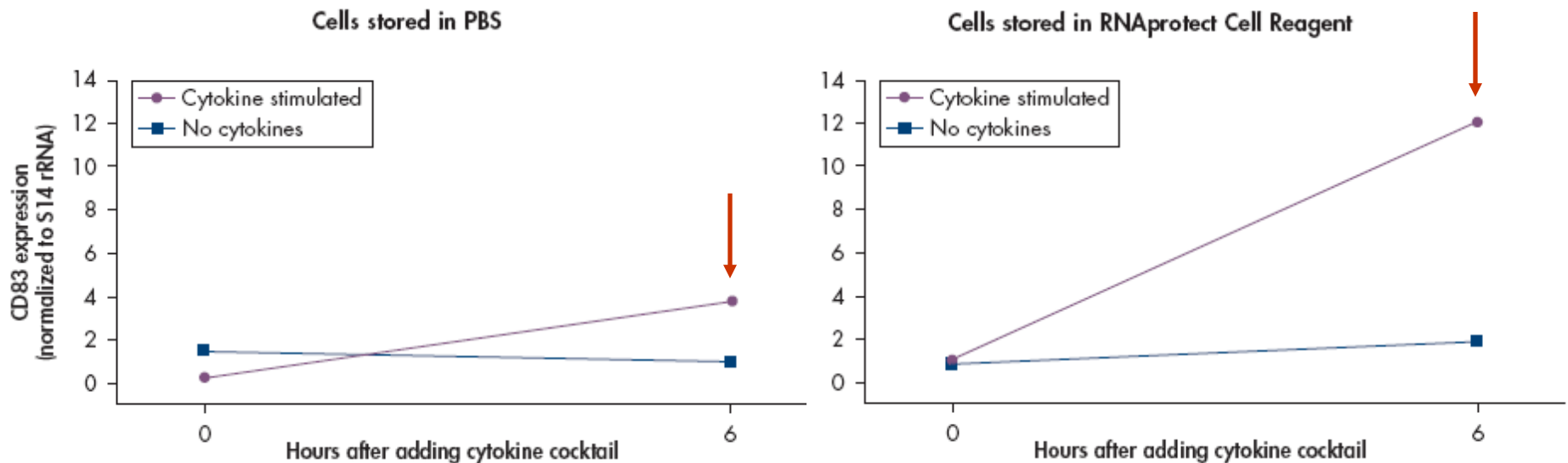


- Rat liver
- Total protein purified with Qproteome Mammalian Protein Prep Kit
- ERK2 analysis by SDS PAGE and western blotting
- Control: Liquid nitrogen

▶ Stabilization of proteins for up to 12 months without freezing

RNAprotect Cell Reagent:

- Dendritic cells stimulated with cytokines
- Stored in PBS or RNAprotect immediately at -20°C for 6 hours
- Quantification of CD83 expression
- Normalization to S14 rRNA



Data kindly provided by Dr. Alexander Prechtel, University Hospital Erlangen, Germany.

- Requirements for analysis of precious samples

- Precious and inhomogeneous samples
- Different analytes from the same sample

- Stabilization of RNA and DNA

- **Simultaneous purification of different analytes**

- Simultaneous analysis of DNA, RNA, and protein**
- DNA, RNA, and miRNA from challenging starting material**
- Analysis of DNA and RNA from FFPE samples**

- Downstream applications


- Summary

Multiple analytes from the same sample

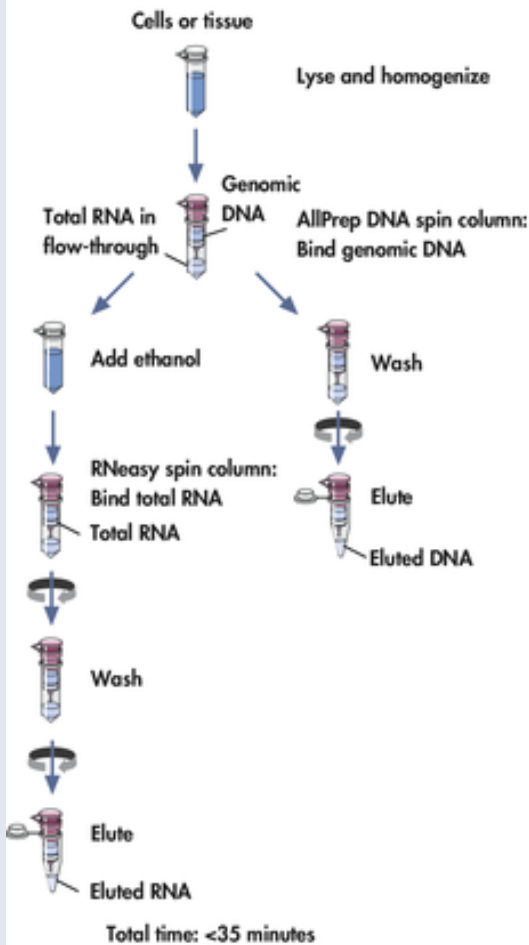
All information derived from the same sample:

- Precious, nonhomogeneous samples
 - All analytes from exactly the same part of the sample
- Complete picture needed

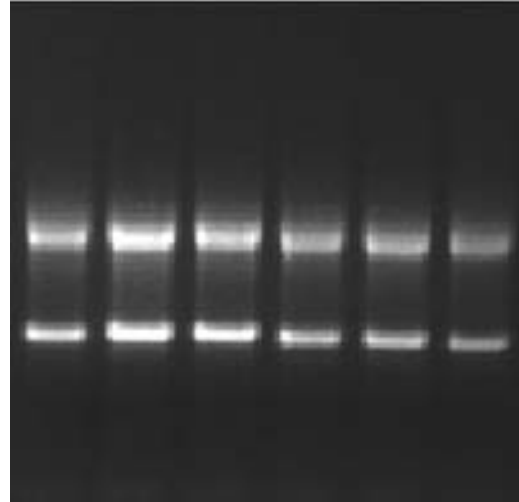


- AllPrep DNA/RNA Mini Kit
- AllPrep DNA/RNA Micro Kit
- AllPrep DNA/RNA 96 Kit
- AllPrep DNA/RNA FFPE Kit
- AllPrep DNA/RNA/Protein Kit
- AllPrep RNA/Protein Kit
- AllPrep gDNA/RNA/miRNA Universal Kit 

AllPrep DNA/RNA Procedure

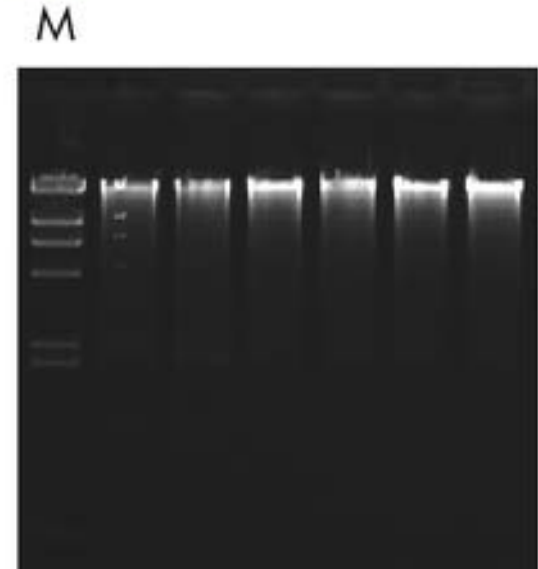


A RNA



- Rat liver (10 mg) analyzed
- 6 independent samples
- Analyzed on 1.2% formaldehyde agarose gel

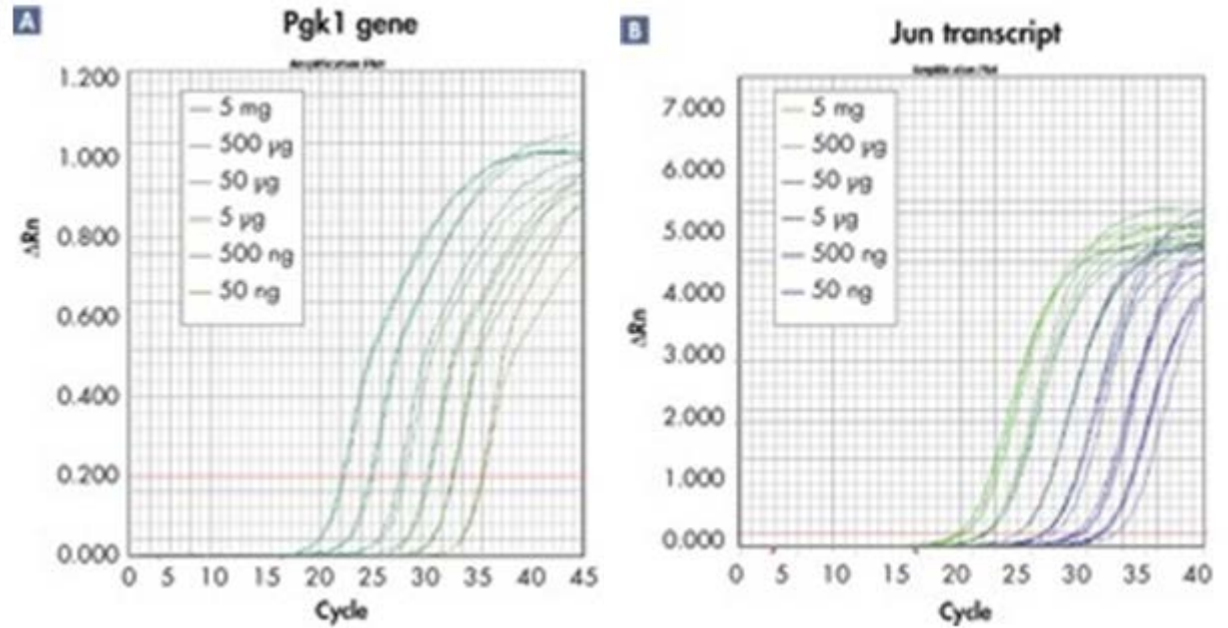
B Genomic DNA



- Rat liver (10 mg) analyzed
- 6 independent samples
- Analyzed on 0.8% agarose gel
- M: Size marker

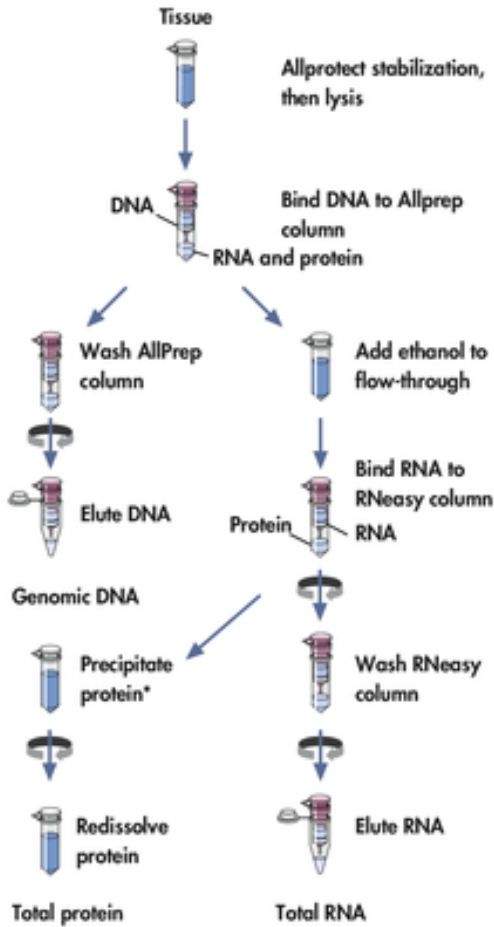
▶ High-quality of RNA and gDNA with high reliability

Serial dilutions of rat kidney homogenate



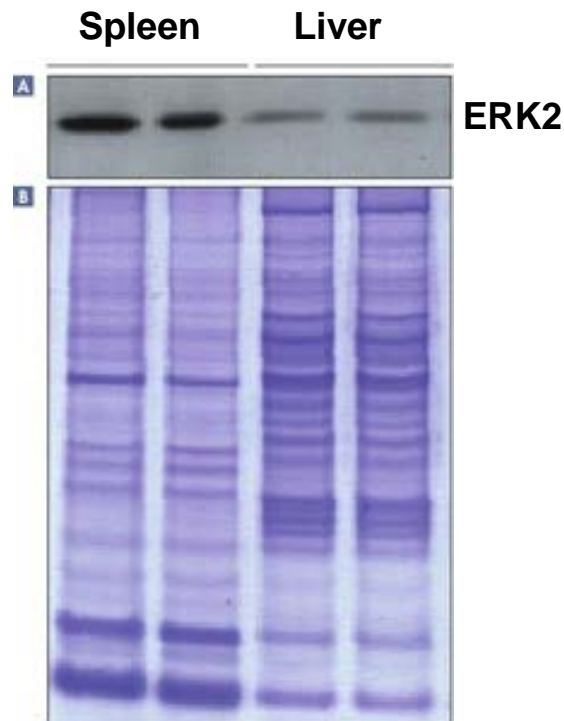
► Wide dynamic range also for small sample sizes

Allprotect/AllPrep Tissue Procedure



* Using a novel protein precipitation solution.

- Technology is based on the proven Allprep DNA/RNA technology
- Additional aqueous protein precipitation solution
- miRNA isolation possible



► Reliable western blotting and SDS-PAGE results

Kit optimized for various difficult-to-lyse starting materials

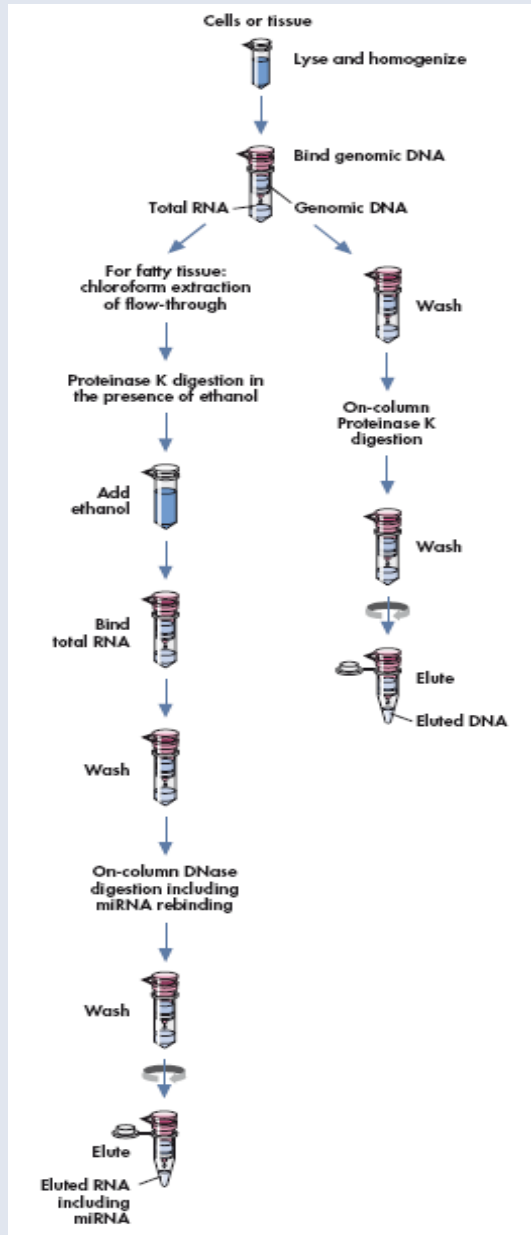
- Suitable for lipid-rich tissues (e.g., brain, adipose tissue)
- Suitable for fiber-rich tissues (e.g., muscle, heart)
- Suitable for isolated blood cells
- Lipid- and fiber- rich tissues are often very scientifically relevant
- Isolation of nucleic acids from these samples is highly difficult
- Difficult to obtain sufficient yields and quality with existing kits



New Allprep DNA/RNA/miRNA Universal Kit

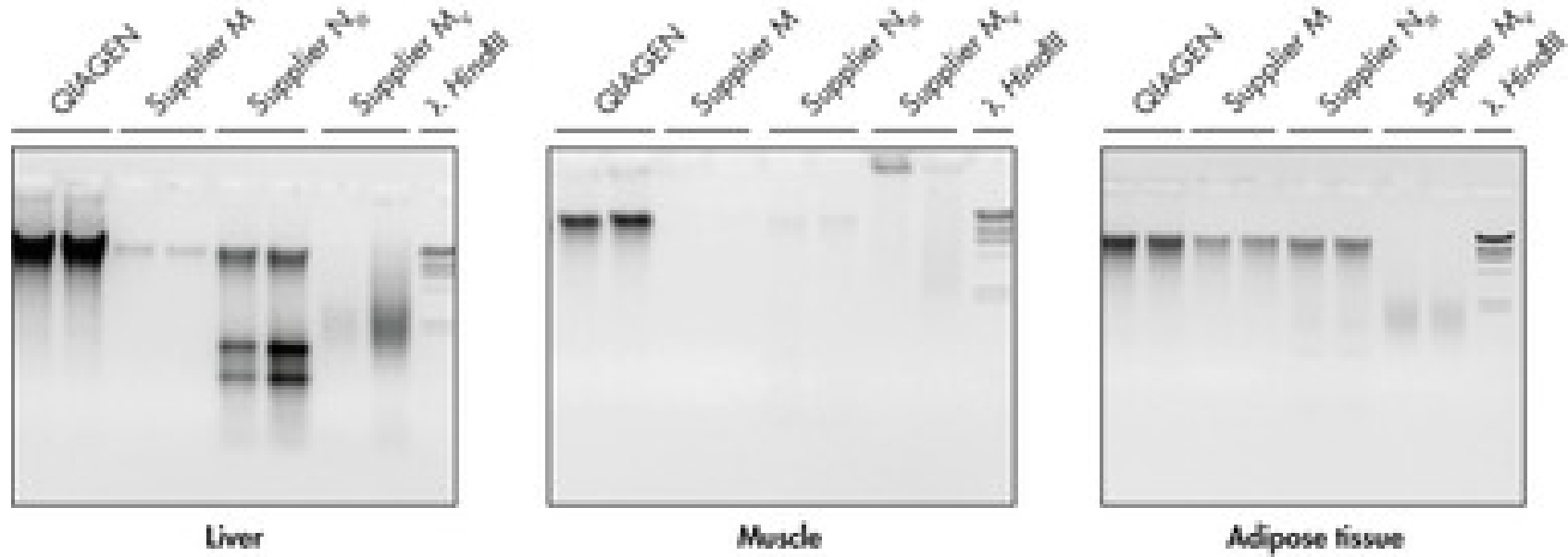
Protocol and technology

- Based on well-established AllPrep DNA/RNA technology
- Optimized proteolytic digestion steps for RNA and DNA
- No phenol required
- Co-purification of miRNA



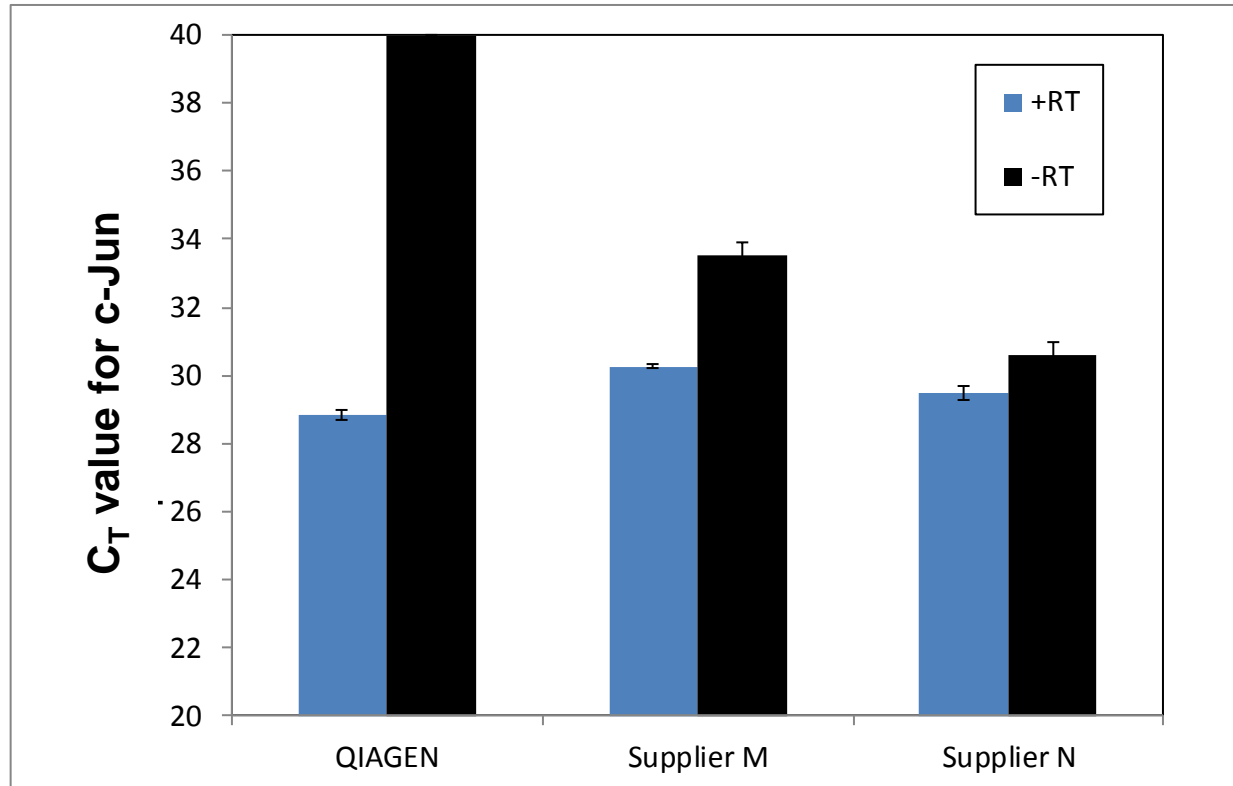
▶ High-quality results from precious and challenging samples

Purification of high yields of DNA from difficult-to-lyse tissues



▶ Intact genomic DNA from most challenging tissue types

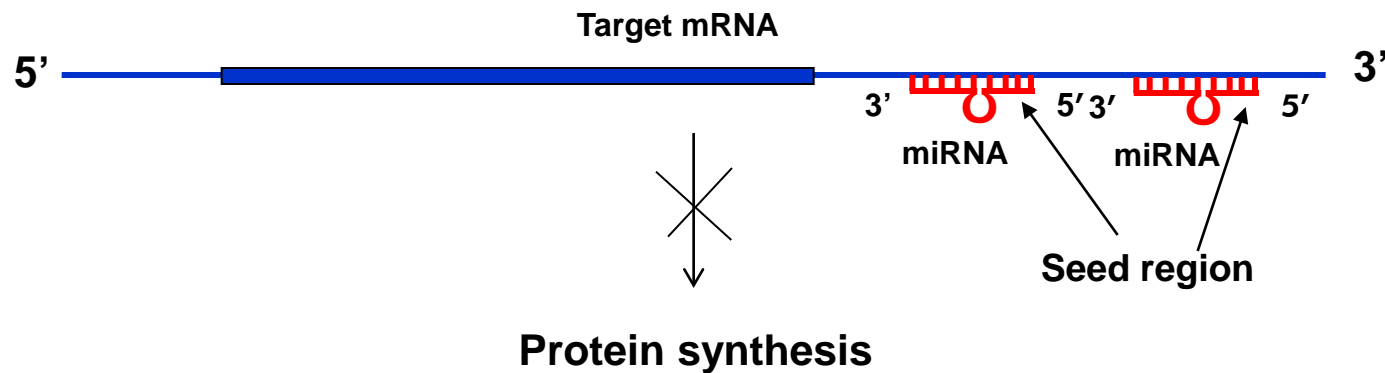
Rat muscle tissue: c-Jun expression



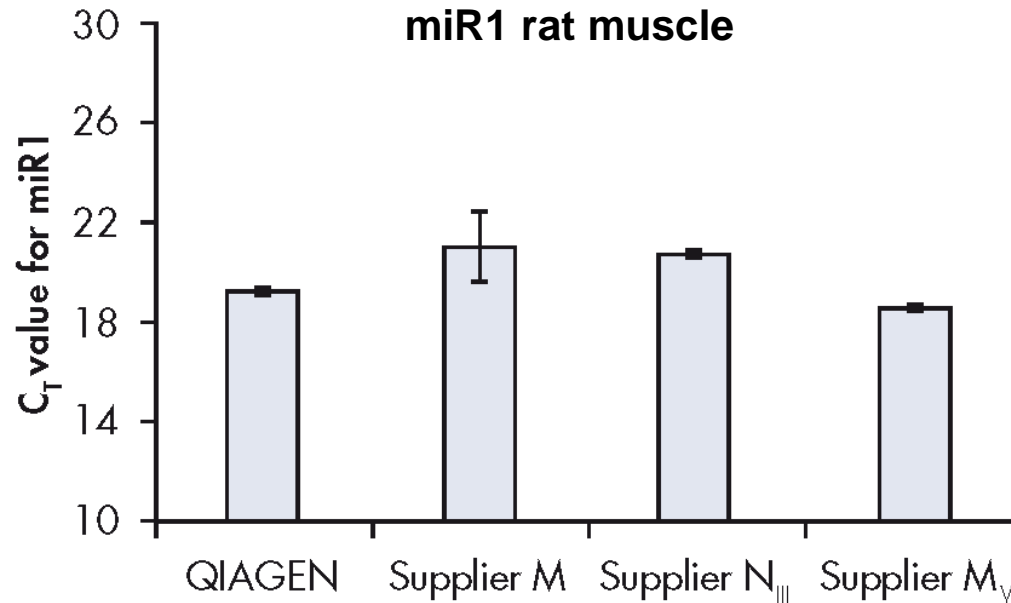
Complete removal of genomic DNA from RNA fraction

microRNAs — micromanagers of gene expression

- Naturally occurring, endogenous small RNA
- Mature miRNA approximately 22 nt long
- Regulate at least 1/3 of protein encoding genes
- miRNA binding sites typically at 3'-UTR of target transcripts
- Imperfectly base pair to target resulting in translational repression
- > 1700 miRNAs in human (miRBase V17.0)



Purification of miRNA from difficult-to-lyse tissues



▶ High yields of miRNA also from difficult-to-lyse tissues



Formaldehyde-fixed paraffin-embedded samples

- Formaldehyde fixation resulting in cross linking of:
 - Protein– protein
 - Protein–nucleic acid
 - Nucleic acid–nucleic acid
- Paraffin embedding resulting in:
 - Nucleic acid fragmentation
 - Deparaffinization required prior to nucleic acid preparation
- Quality of nucleic acids from FFPE samples compromised by:
 - Remaining crosslinks
 - Chemical modifications of analytes
 - Fragmentation

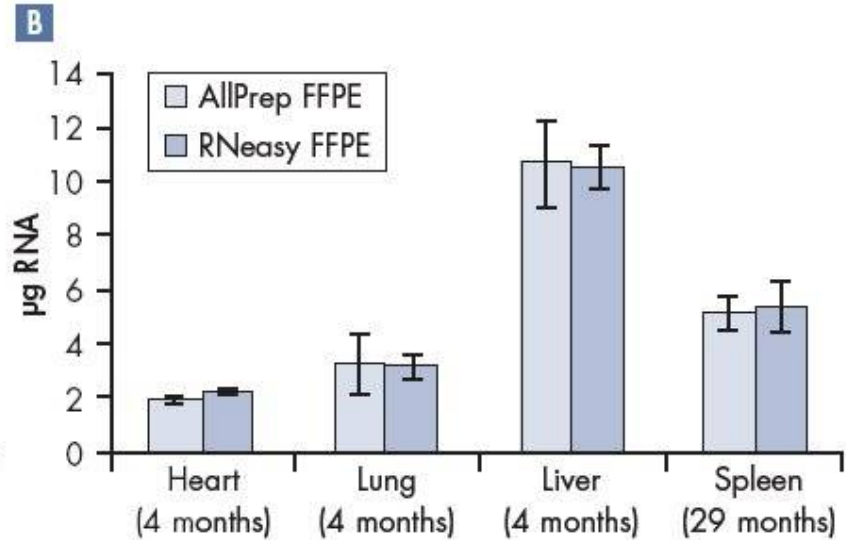
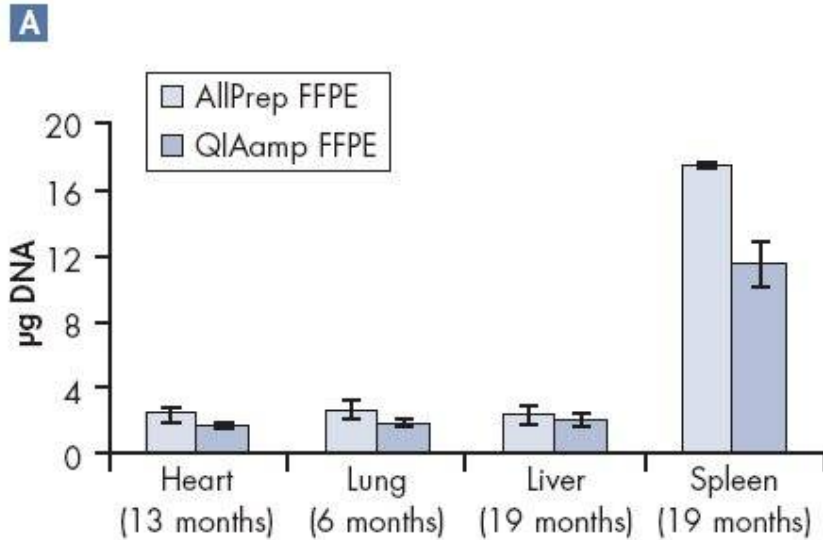
Consequences

- For effective lysis, harsh conditions required
- Poor yield (especially due to overfixation)
- Interference with enzymatic assays

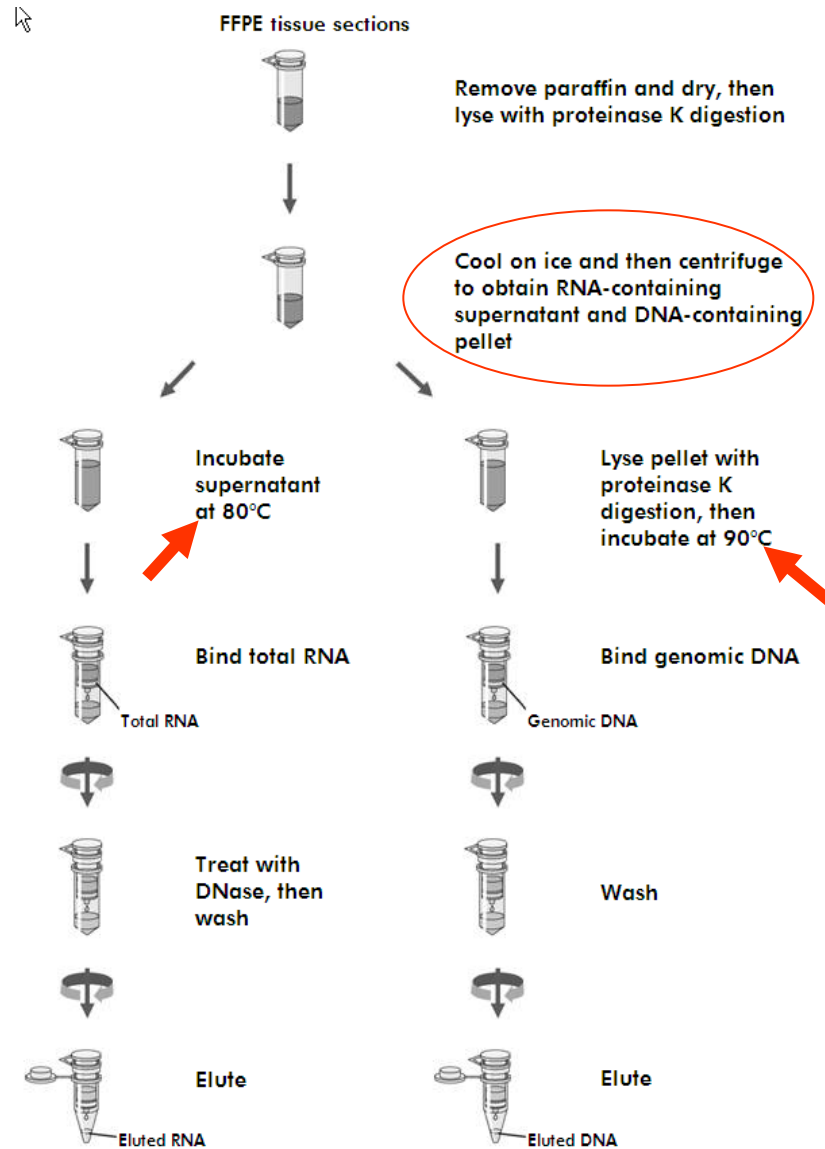


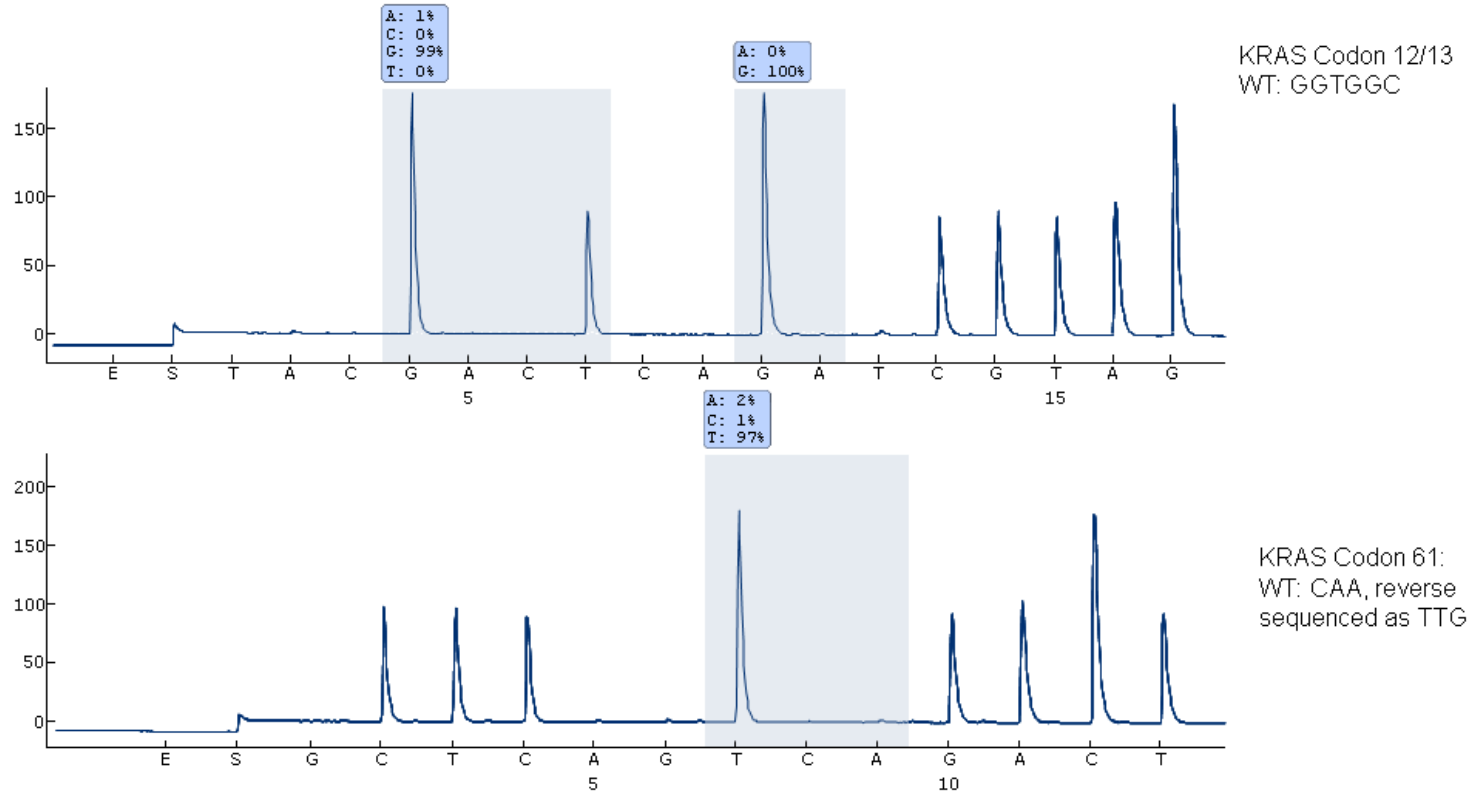
AllPrep DNA/RNA FFPE Kit:

- Simultaneous purification of DNA and RNA from FFPE samples
- Differential solubilization methodology
- Efficient removal of crosslinks
- Yield/quality comparable to that obtained with specialized kits
- Includes efficient genomic DNA removal
- Protocol including miRNA available



AllPrep DNA/RNA FFPE Kit - protocol





- DNA purified from breast cancer FFPE tissue using AllPrep DNA/RNA FFPE Kit
- Pyrosequencing to identify mutation in KRAS gene
 (*therascreen* KRAS Pyro Kit)

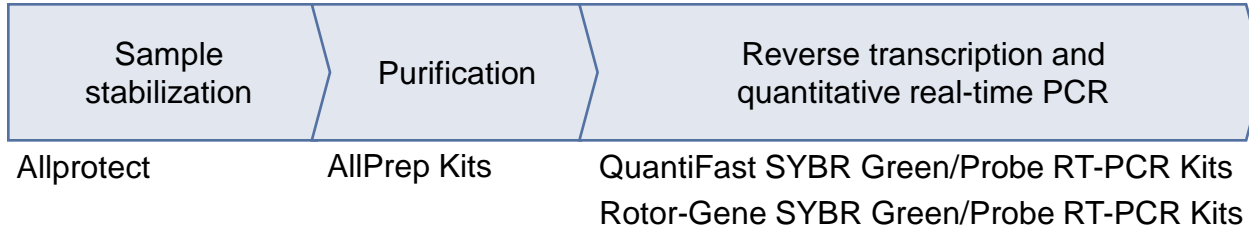
- Requirements for analysis of precious samples
 - Precious and inhomogeneous samples
 - Different analytes from the same sample
- Stabilization of RNA and DNA
- Simultaneous purification of different analytes
 - Simultaneous analysis of DNA, RNA, and protein
 - gDNA, RNA, and miRNA from challenging starting material
 - Analysis of DNA and RNA from FFPE samples

■ **Downstream applications**

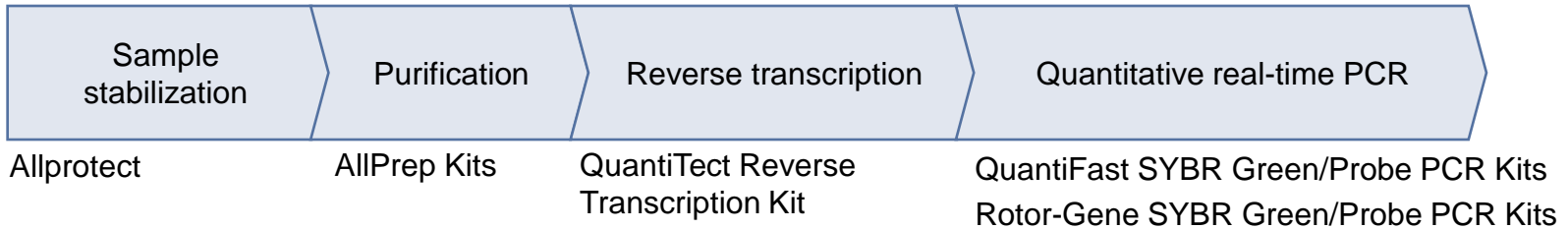
- Summary

One-step or two-step RT-PCR

One-step RT-PCR



Two-step RT-PCR



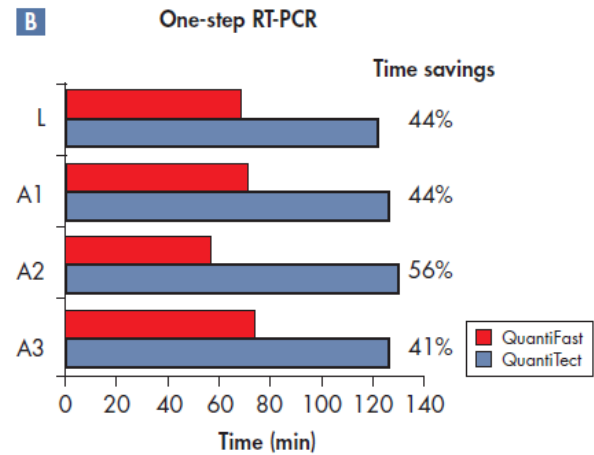
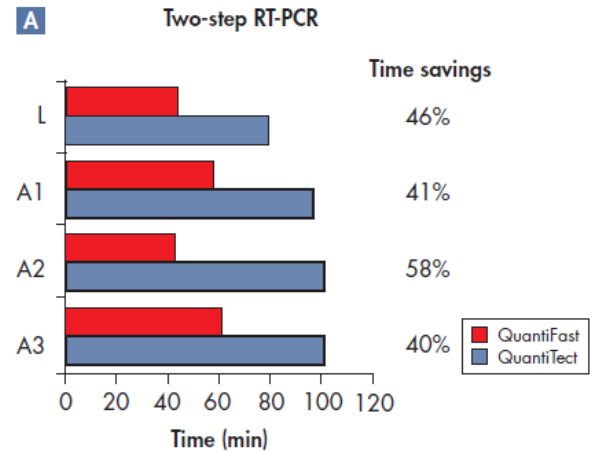
▶ Various kits for probe-based and SYBR Green detection and one-step and two-step RT-PCR

Unique composition for success at the first attempt

Enzymes

- HotStarTaq *Plus* DNA Polymerase
 - Reactivation within 5 minutes
 - Unmatched specificity and sensitivity

- QuantiFast RT Mix (for one-step kits)
 - Optimized combination of Sensiscript and Omniscript
 - High sensitivity

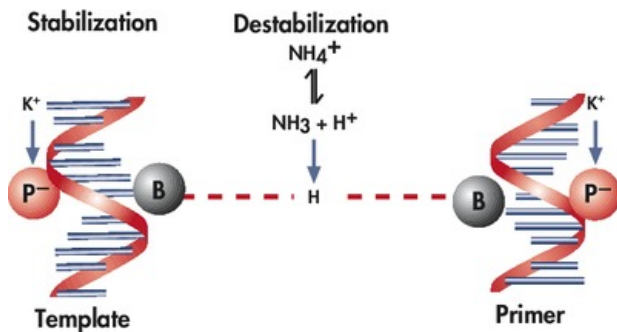


▶ Significantly reduced qPCR times

Unique composition for success at the first attempt

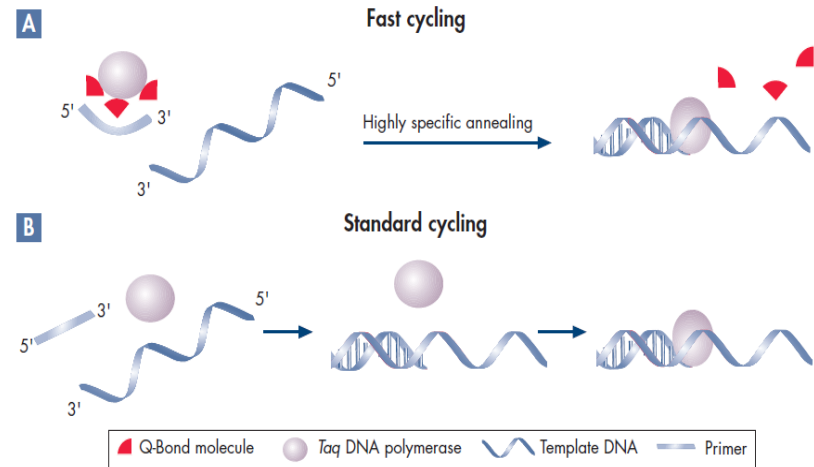
Unique qPCR reaction buffer

- Stabilizes specific bonds
- Destabilizes unspecific bonds



Q-Bond

- Increases affinity of DNA polymerase to short ssDNA
- Reduces primer annealing time to a few seconds



► Successful qPCR analysis without optimization

cDNA Synthesis from miRNA

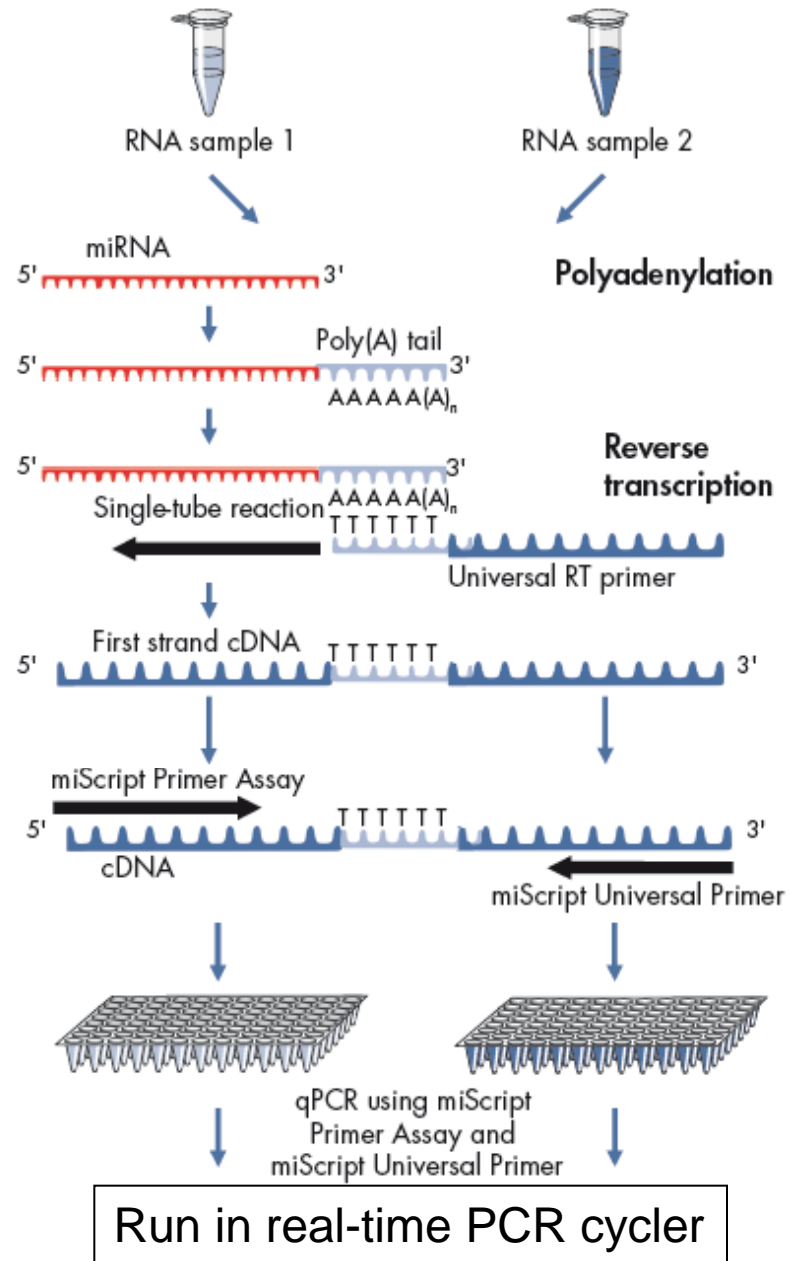
1 hour

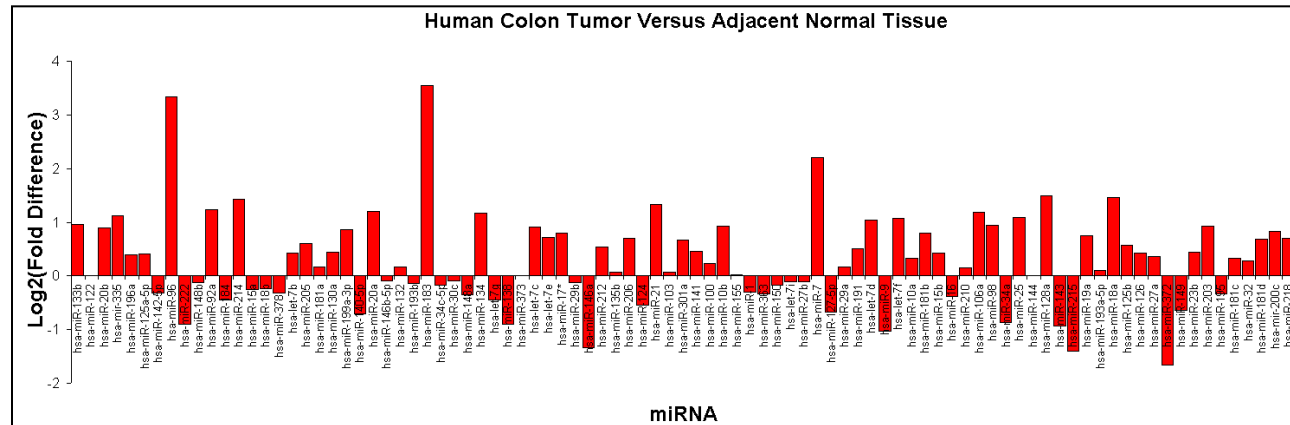
Add cDNA to master mix and load plates

2 minutes

Run 40 cycle qPCR program

2 hours





Many cancer-specific miRNA are upregulated in a colon tumor

- Requirements for analysis of precious samples
 - Precious and inhomogeneous samples
 - Different analytes from the same sample
- Stabilization of RNA and DNA
- Simultaneous purification of different analytes
 - Simultaneous analysis of DNA, RNA, and protein
 - gDNA, RNA, and miRNA from challenging starting material
 - Analysis of DNA and RNA from FFPE samples
- Downstream applications

■ Summary



- Induction and degradation RNA: a general risk
 - Careful handling of samples and stabilization needed
- Precious samples often small or nonhomogeneous
 - Co-purification of DNA, RNA, miRNA, and protein
- Lipid- and fiber-rich tissues are challenging to purify
- FFPE samples with cross-linked and fragmented nucleic acids
 - Use of specialized technologies
- Downstream applications

Thank you for your attention!



Stefanie.Schroeer@qiagen.com

For up-to-date licensing information and product-specific disclaimers, see the respective QIAGEN kit handbook or user manual. QIAGEN kit handbooks and user manuals are available at www.qiagen.com or can be requested from QIAGEN Technical Services or your local distributor.