

## Circulating Nucleic Acids in Cancer

26.10.2012

Marco Polidori

# Agenda



- Background on fcNA
  - What is fcNA
- Principle
  - Where does fcNA come from in cancer
- Known Alterations
  - That may be associated with cancer
- Advantages of analyzing fcNA
  - What makes fcNA so interesting as diagnostic marker
- fcNA associated biomarkers
  - What examples exist already
  - What may be preferred detection methods
- Challenges of finding Biomarkers
  - What should be considered when looking for new biomarkers
  - Questions about the baseline
- Extraction methods
  - What we can learn from pre-natal diagnostics
  - Our experience on extraction methods

## Background

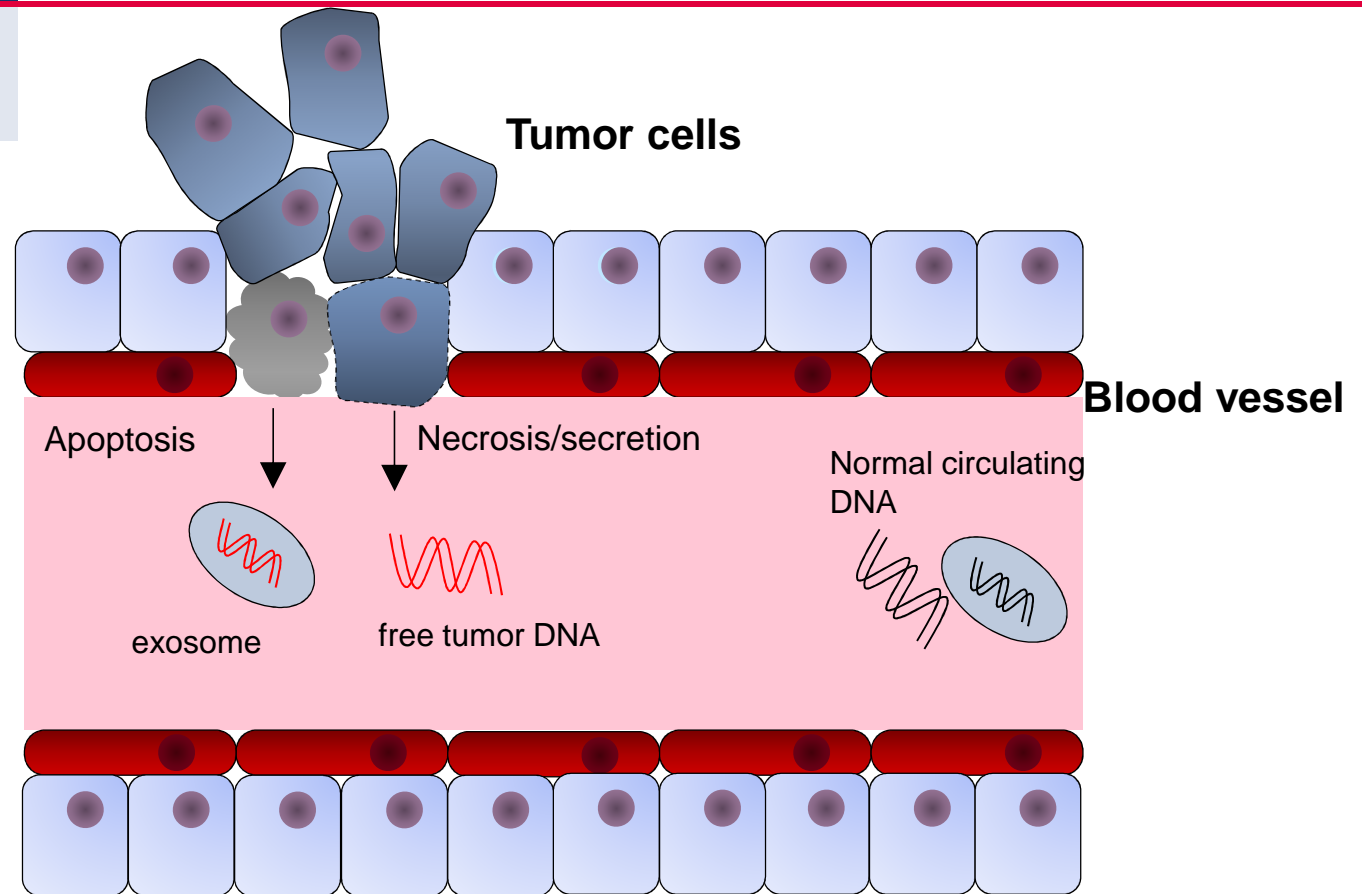
What are free circulating nucleic acids (fcNA) ?

**Cell-free nucleic acids present in body fluids like serum/plasma, urine, CSF etc.**

### Properties

- Cell free DNA, mRNA and miRNA
- Mostly present as shorter fragments (less than 1000 bp or nt)
  - Circulating DNA: predominantly in the range of 140-100 bp
- Can also be several kb long
- Low concentration in plasma and serum (1-100 ng/mL)
- Can be protein-bound (nucleosomes)
- Can be “vesicular” (exosomes)
- Half life of minutes to several days (EDTA plasma)

# Principle of cancer associated fcDNA



Release of DNA is thought to occur through

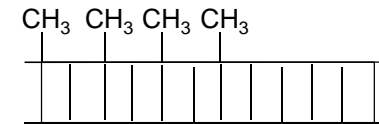
- Apoptosis
- Necrosis
- Active secretion

# Alterations of fcNA

Various alterations of fcNA that may indicate disease MH3

Main alterations known

- Methylation (epigenetic alterations)



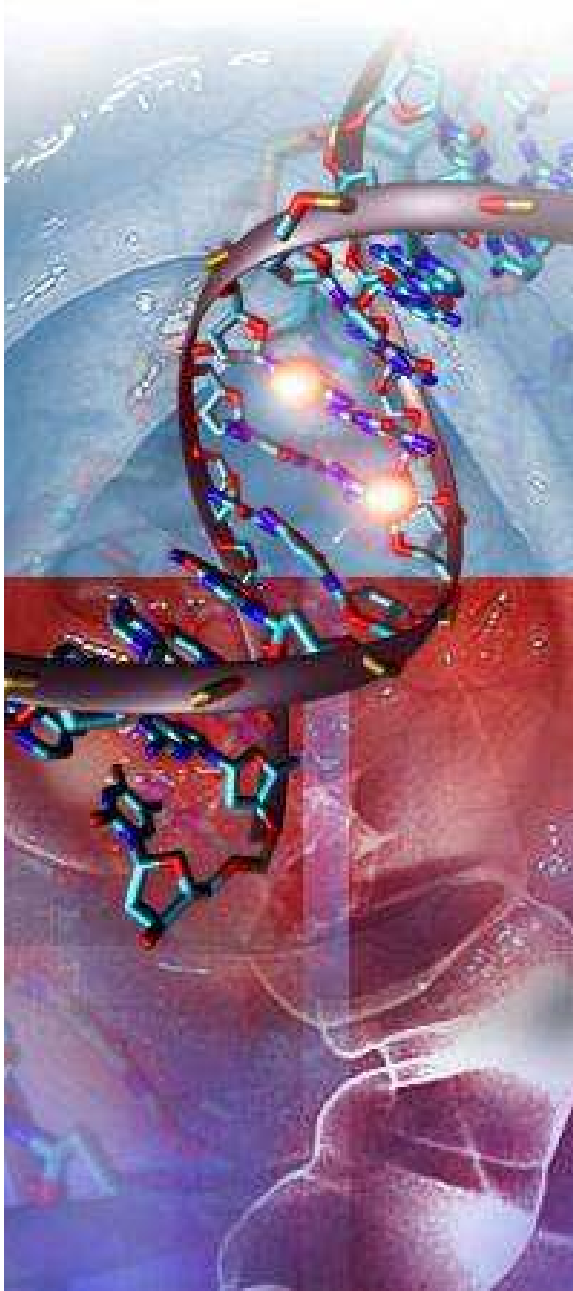
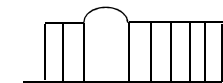
- Mutation (i.e. oncogenes like *KRAS*)



- Microsatellite alterations



- miRNA expression levels



## Folie 5

---

### MH3

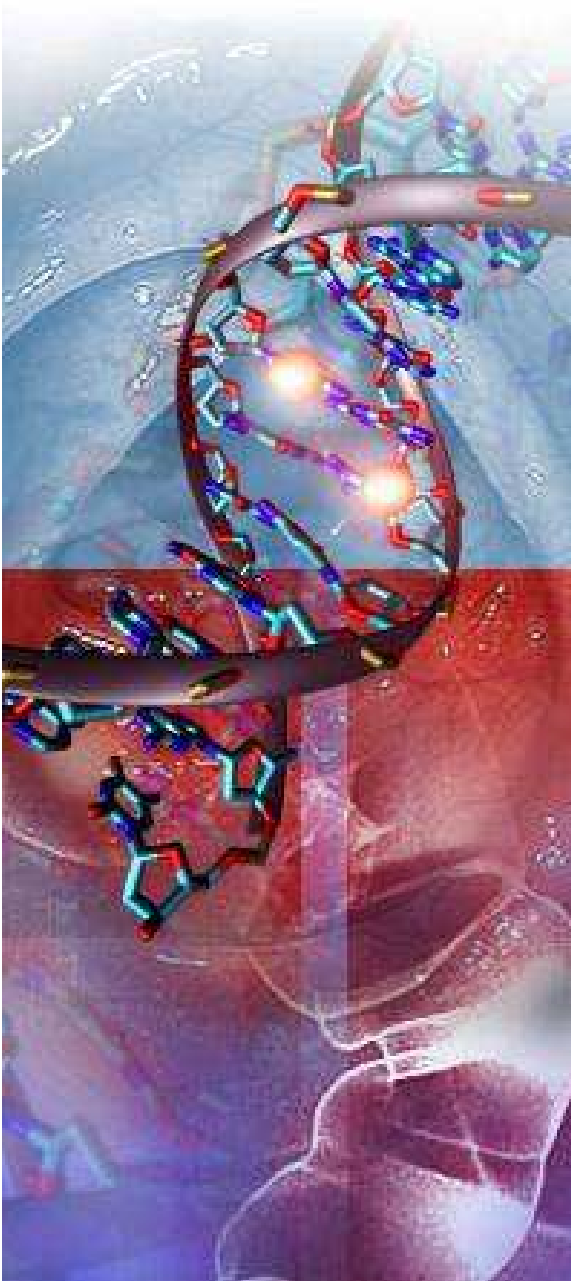
also:

-- copy number variations

-- chromosomal aneuploidies detectable in fetal ccfDNA in maternal plasma

Martin Horlitz; 23.10.2012

## Advantages of fcNA as analyte



- Minimally invasive
- Blood is a readily available sample material
- DNA from distant metastatic sites is also being captured
- Can also be used recurrently during treatment monitoring

## fcNA associated biomarkers

Many biomarkers have been found to be associated with cancer

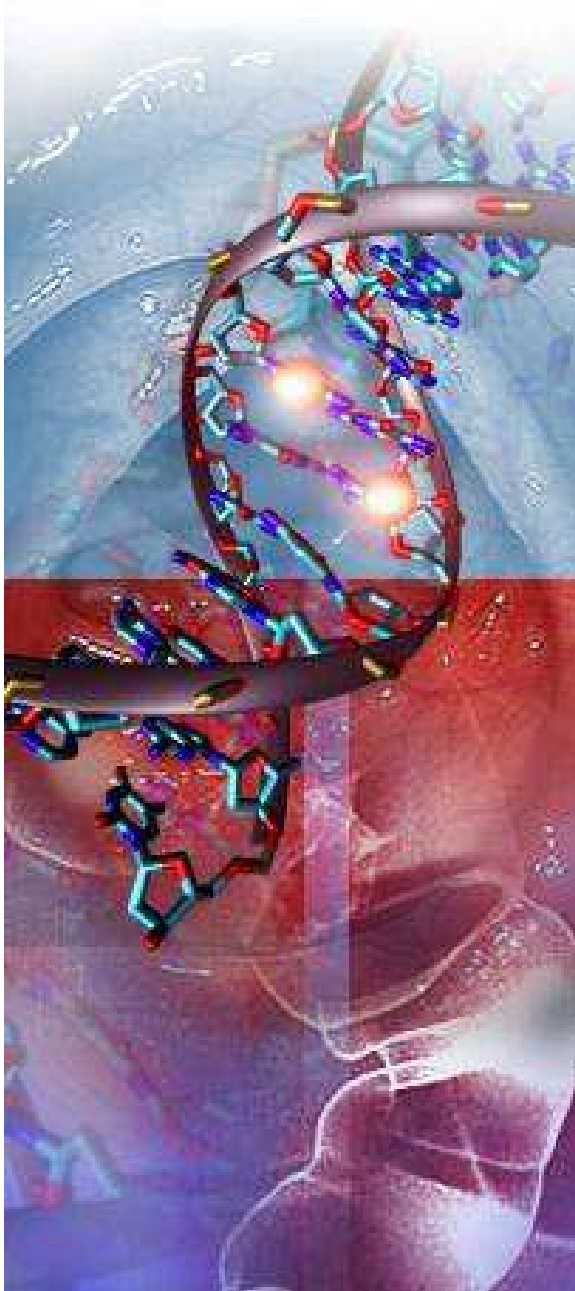
Examples include

- Breast (e.g. PIK3CA)
- Prostate (e.g. miR 21/141, GSTP1 methylation)
- Lung (e.g. **SHOX2** methylation, EGFR)
- Ovarian (e.g. TP53)
- Colorectal (e.g. **SEPT9** methylation, KRAS)
- Pancreatic (methylation pattern of several markers including SOCS1, CCND2 and others)
- Melanoma (e.g. BRAF)

And several others

Don't forget about

- CNVs
- Chromosomal rearrangements



- “The needle in the haystack”
  - What markers are already known from biopsies for a certain cancer type?
  - Can it also be detected in fcNA
- fcNA concentrations are usually very low
  - assays have to be extremely sensitive
  - Have to be detectable against high background of WT DNA
- fcNA levels increase in advanced stages of cancer
  - Early diagnostics can be challenging using fcNA since concentrations are too low in these stages
- High diversity of cancer genomics may require novel approaches to detect
  - e.g. mutations with low allelic frequency
- Variable sample processing methods are being used
  - Uniform blood collection & extraction method needed

More studies needed with large cohorts

## Do we know enough about the general biology of fcNA?

More studies are required to understand the natural occurrence of fcNA:

- Which factors influence formation of cNA in healthy humans?
- General health, implications of other common diseases (e.g. infections)
- Circadian rhythms
- Age
- Way of life (nutrition, sports, etc.)

Shall help finding better guidelines

- When is it best to draw blood for diagnostic testing?
- Take other implications into account (when to not use fcNA)

Reduce background as much as possible

## Extraction methods



Uniform extraction method needed

Blood stabilization agents?

serum or plasma?

How much serum/plasma?

Size cut-offs?

## What we can learn from pre-natal DX

Looking beyond the horizon of cancer could be beneficial. Pre-natal Diagnostics is already well established.

K<sub>3</sub>EDTA tubes (Barrett et al.)

Use Plasma (if applicable)

As much sample as possible (>1 ml)

- Whole Genome Sequencing shown to work from maternal plasma (Kitzman et al.; Fan et al.)

Also applicable for cancer research!

- Massively parallel sequencing (targeted sequencing using plasma from cancer patients (Forsheew et al.; Morgan et al.)
- WGS approach on plasma from cancer patients (Chan et al.)

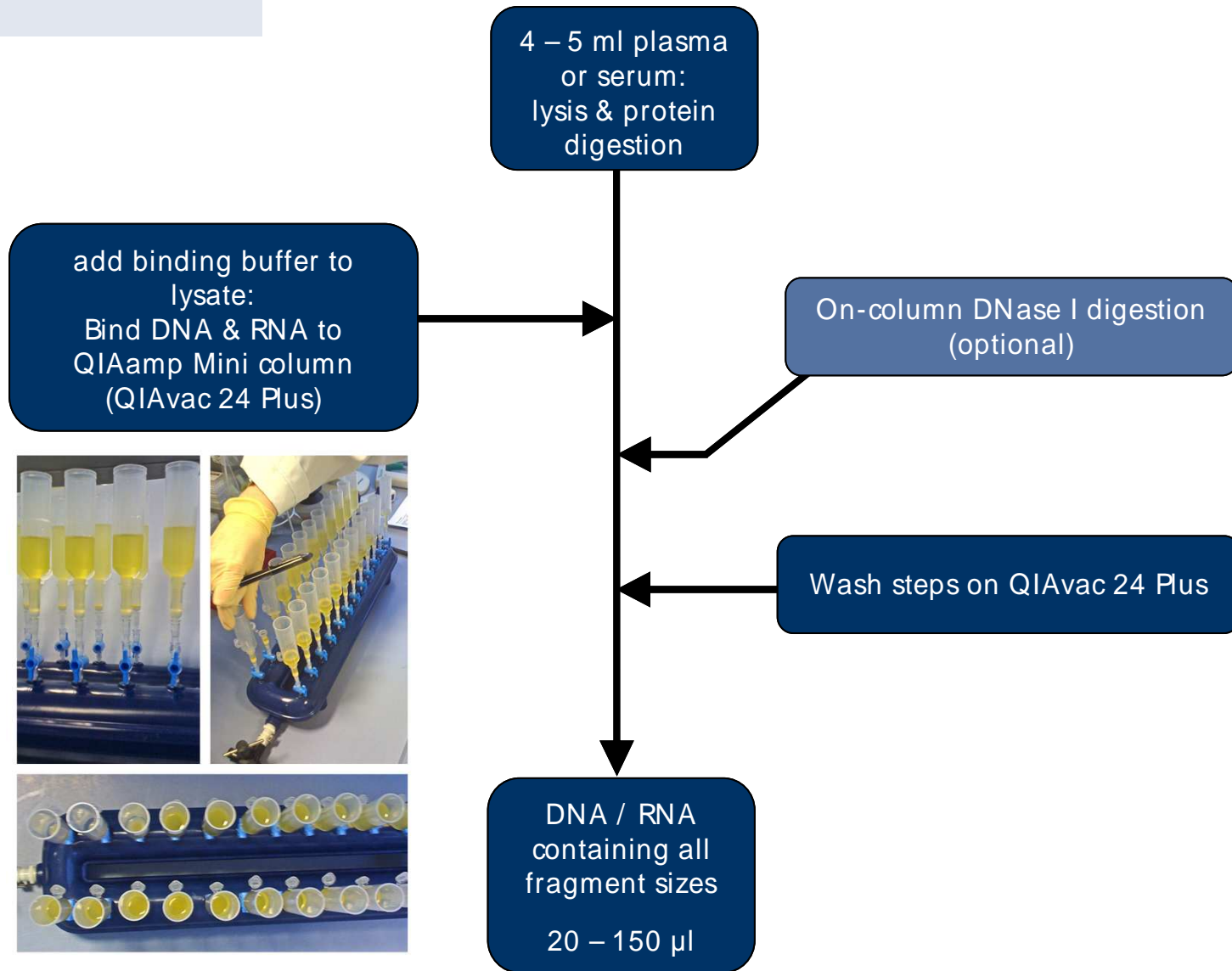


## QIAamp circulating NA Kit

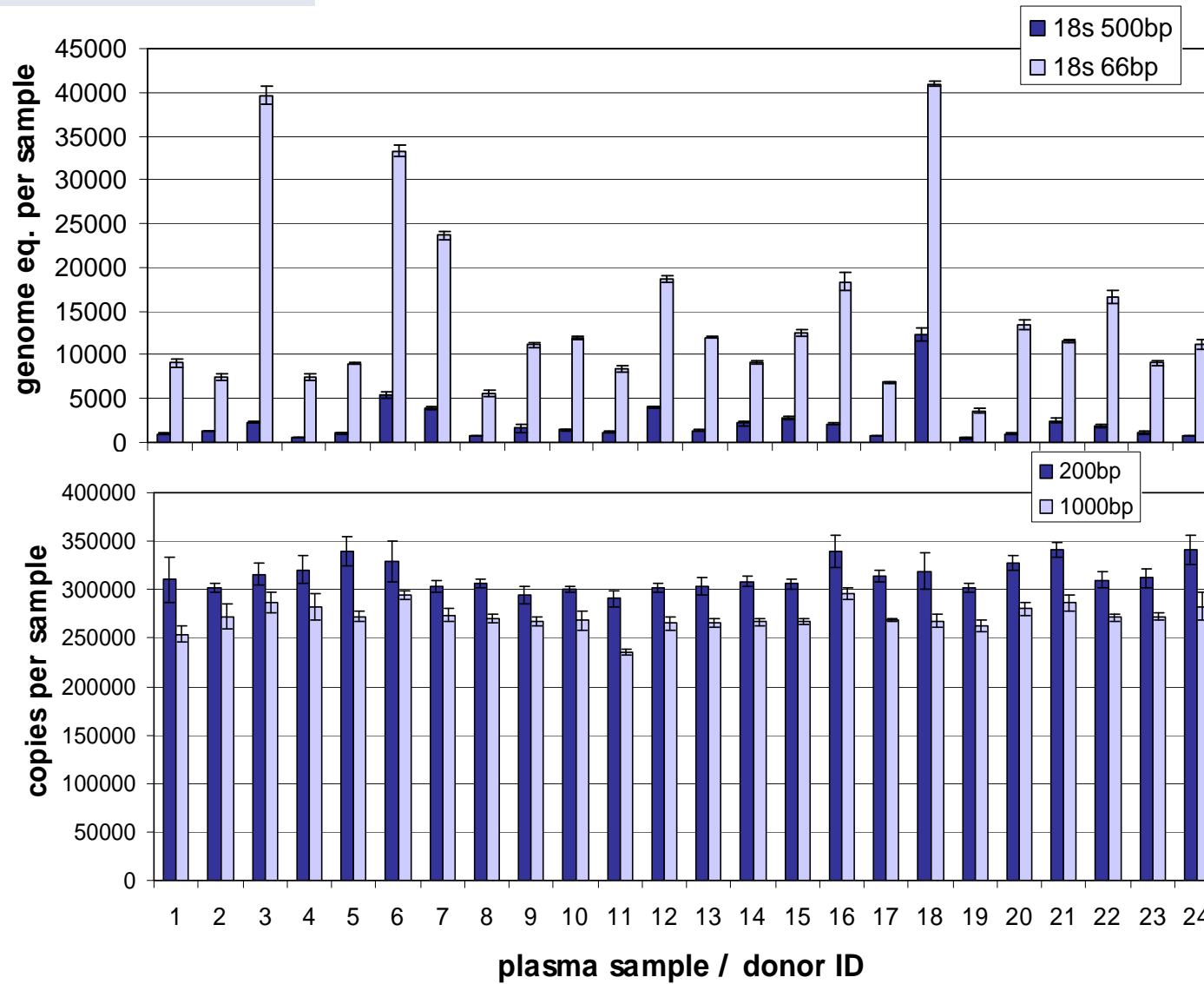


- High input volumes possible (up to 5 ml sample)
- No size bias: small and large fragments are recovered
- Optimized serum/plasma protocol
- Compatible with urine
- miRNA protocols available

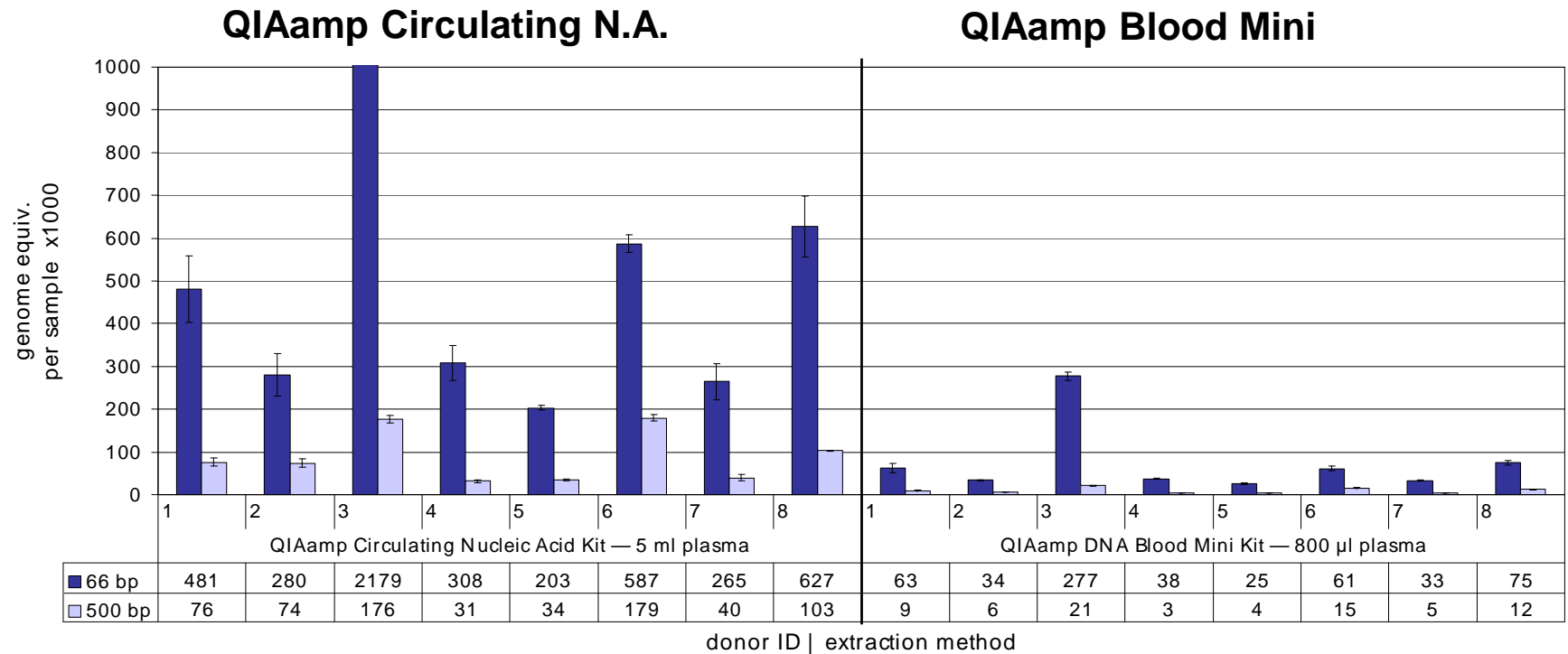
# QIAamp circulating Nucleic Acid Kit workflow



# Recovery rates of fcNA



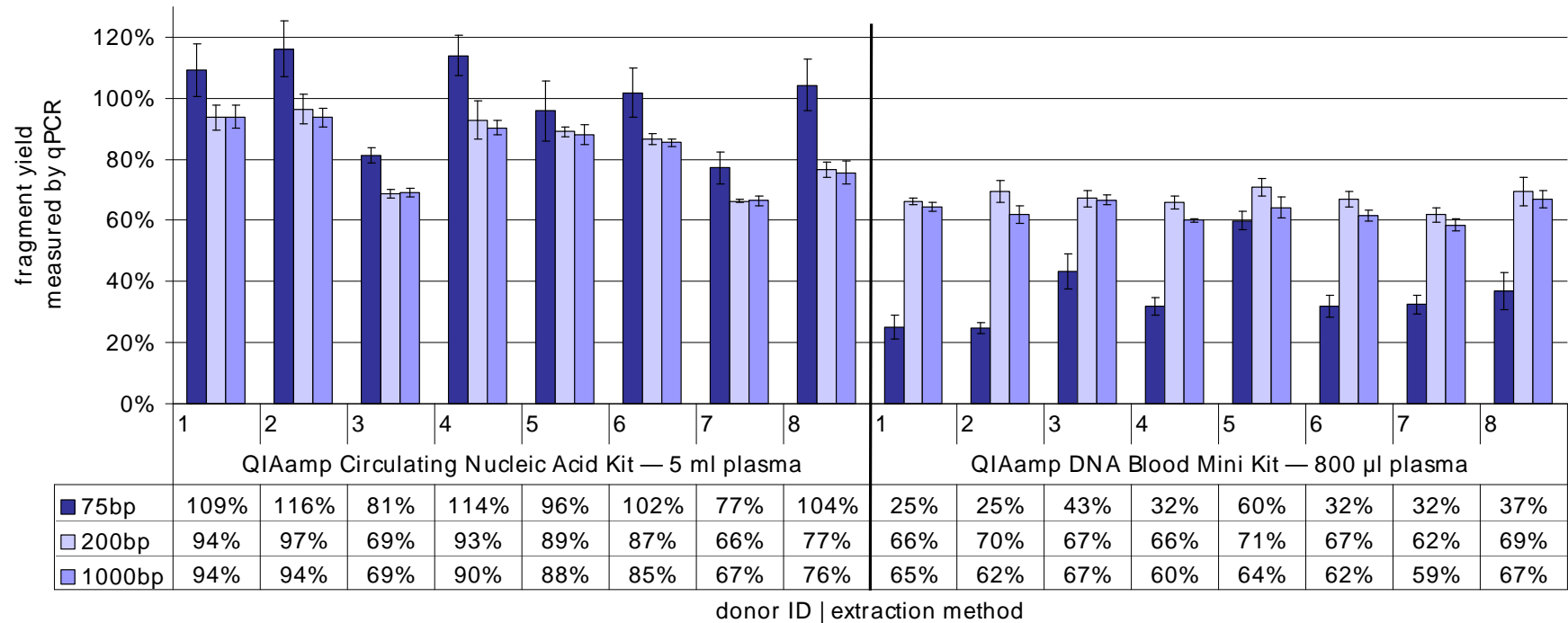
Enabling large volume preparations increases performance drastically



## Circulating DNA in EDTA plasma — individual donors

- Circulating DNA yield: 18S rDNA PCR — 66 bp and 500 bp amplicons
- Comparison to QIAamp DNA Blood Mini Kit (800 µl protocol)
  - Chemistry of QIAamp C.NA kit improves recovery of circulating DNA beyond the effect of higher sample volume

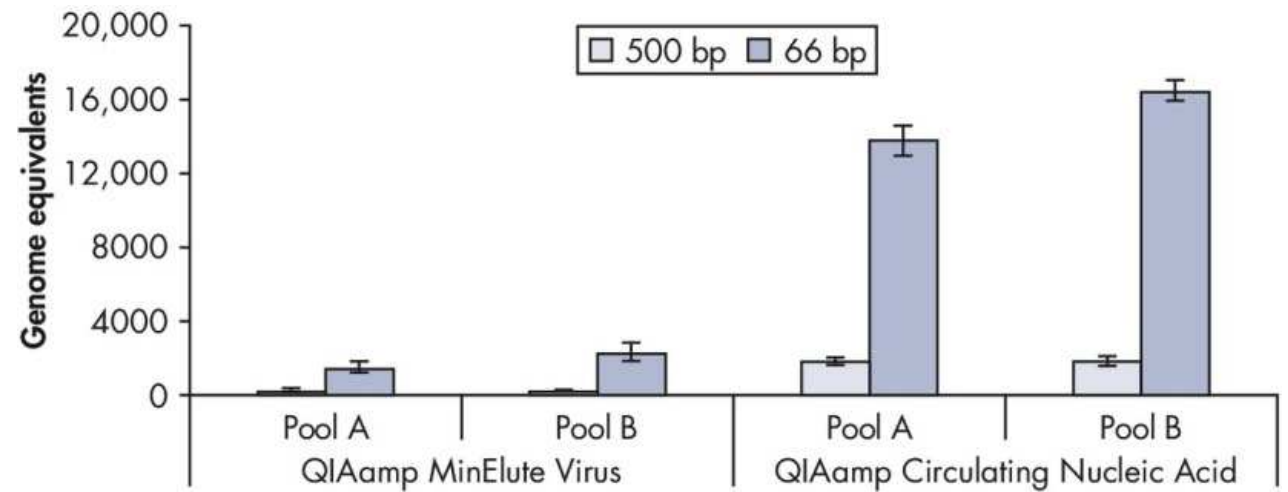
## No size bias of the QIAamp circulating NA Kit



## Recovery of spiked DNA fragments from EDTA plasma — individual donors

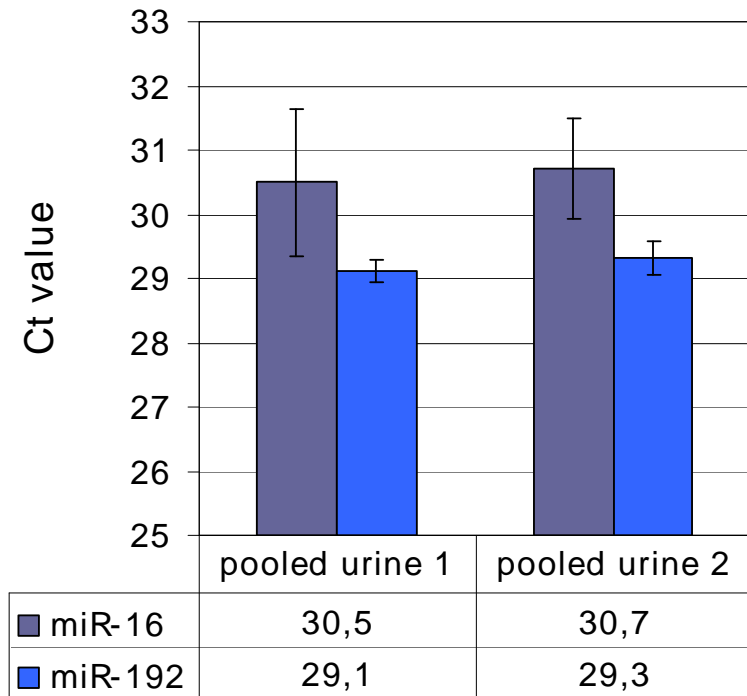
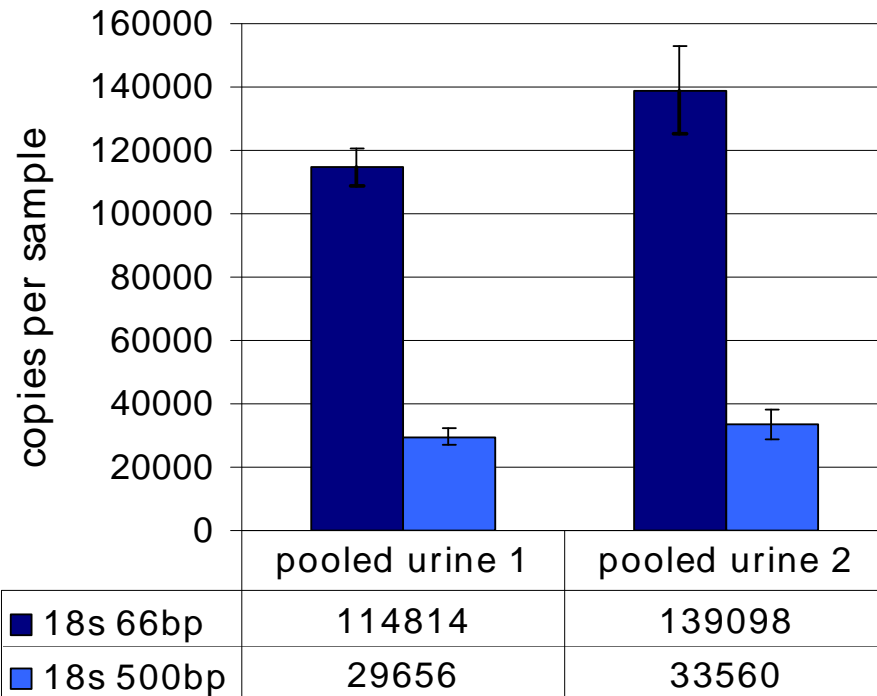
- 75 bp, 200 bp 1000 bp DNA added to plasma before extraction (200,000 copies)
- Chemistry of QIAamp C.NA kit improves recovery of DNA fragments shorter than 200 bp

# QIAamp cNA vs. QIAamp MinElute Virus



## Circulating DNA in EDTA plasma — pooled plasma

- Circulating DNA yield: 18S rDNA PCR — 66 bp and 500 bp amplicons

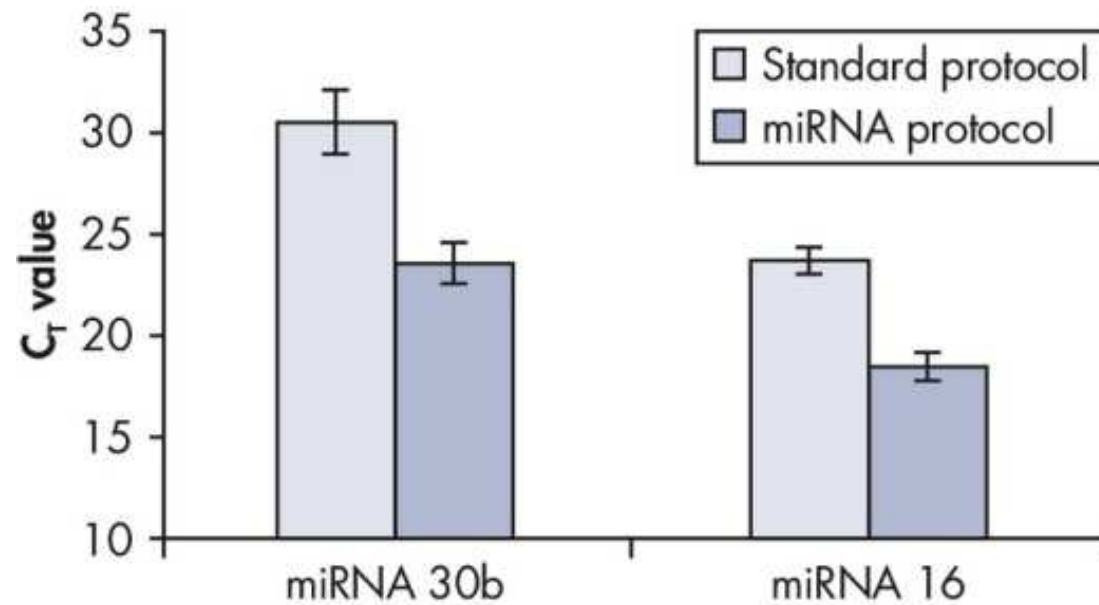


### Circulating DNA and miRNA in pooled urine samples:

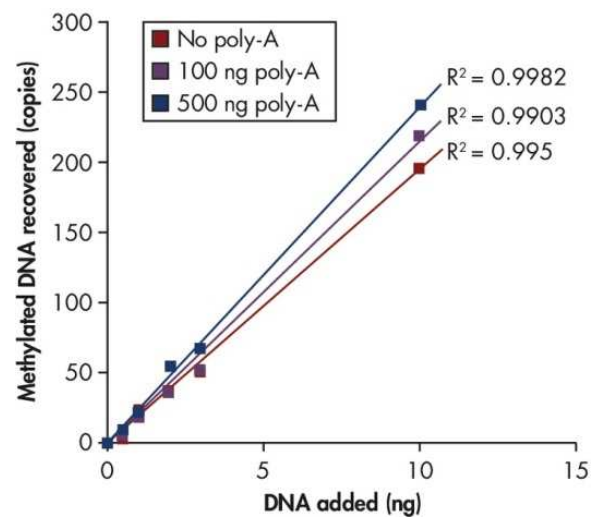
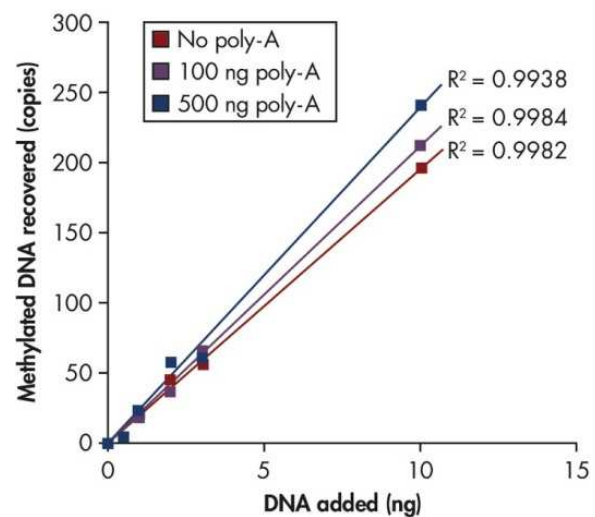
- ccfDNA is present in urine as fragments mostly <500 bp
- miRNAs can be easily detected in samples of 3 ml urine

## Purification of free circulating miRNA

Purification of miRNA can be enhanced using a specialized protocol



# Recovery of methylated DNA





## We are looking for you

---

Please reach out to us if you are interested in collaborations:

[Marco.Polidori@qiagen.com](mailto:Marco.Polidori@qiagen.com)

Asia-Pacific region

[Tina.Li@qiagen.com](mailto:Tina.Li@qiagen.com)

Europe

[Nicole.neubauer@qiagen.com](mailto:Nicole.neubauer@qiagen.com)

North America

[Ray.Ahmed@qiagen.com](mailto:Ray.Ahmed@qiagen.com)

**Thank you for your attention!**



**Marco.Polidori@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](http://www.qiagen.com) or can be requested from QIAGEN Technical Services or your local distributor.