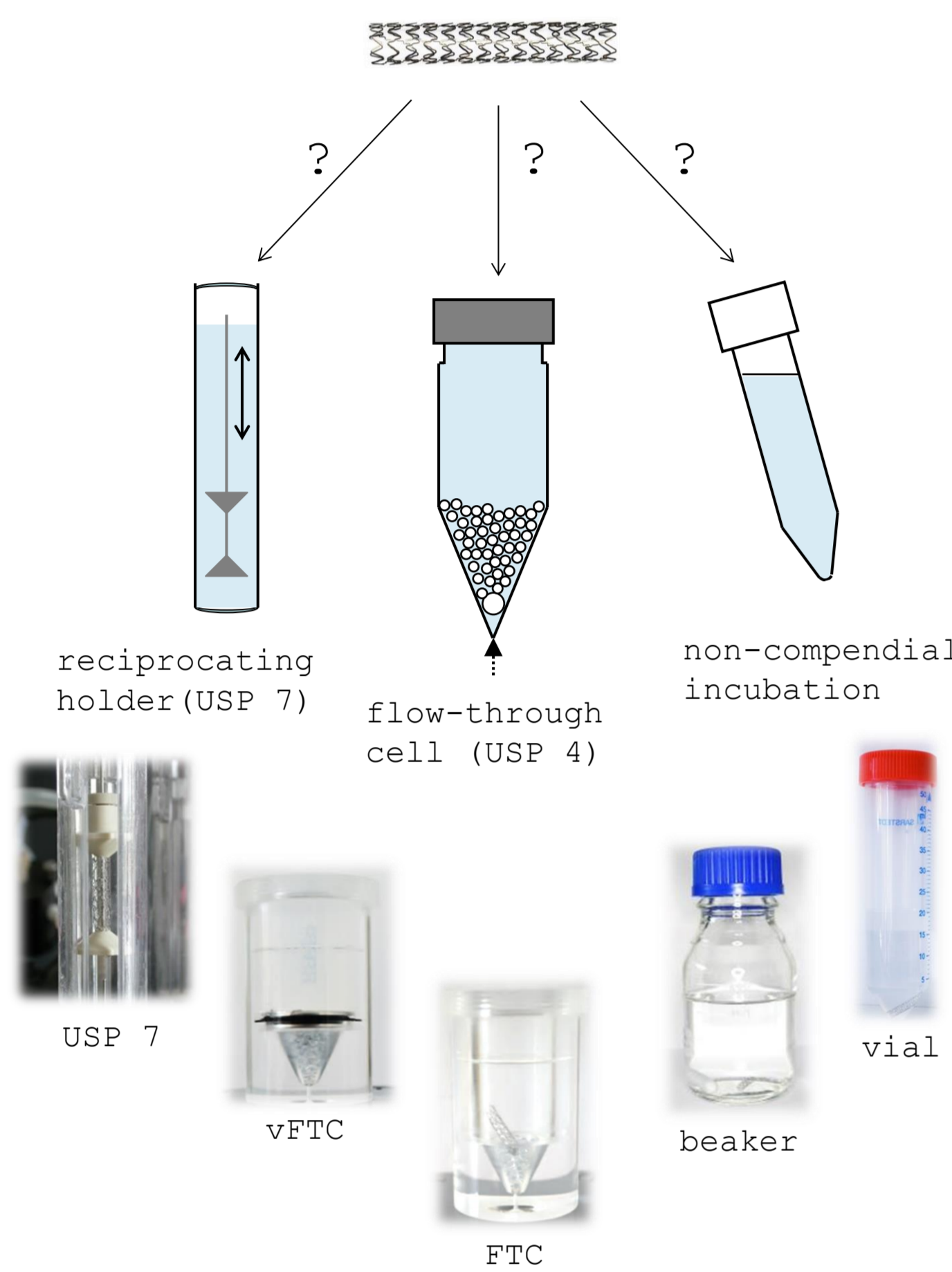




A. Seidlitz<sup>1</sup>, K. Burnashov<sup>1</sup>, T. Reske<sup>2</sup>, V. Senz<sup>2</sup>, N. Grabow<sup>2</sup>, and W. Weitschies<sup>1</sup>  
<sup>1</sup>Center of Drug Absorption and Transport, Institute of Pharmacy, University of Greifswald, Germany  
<sup>2</sup>Institute for Biomedical Engineering, University of Rostock, Germany  
 anne.seidlitz@uni-greifswald.de

## Purpose



### Drug-eluting stents

- Dosage form for local drug delivery (extended release) to stenosed portions of the vessel wall

### In vitro release testing

- No stent-specific method established
- Implementation of flow-conditions and stent embedding in a gel compartment have been shown to impact on in vitro release behavior [1]
- Drug loss may also occur prior to expansion during passage towards site of application

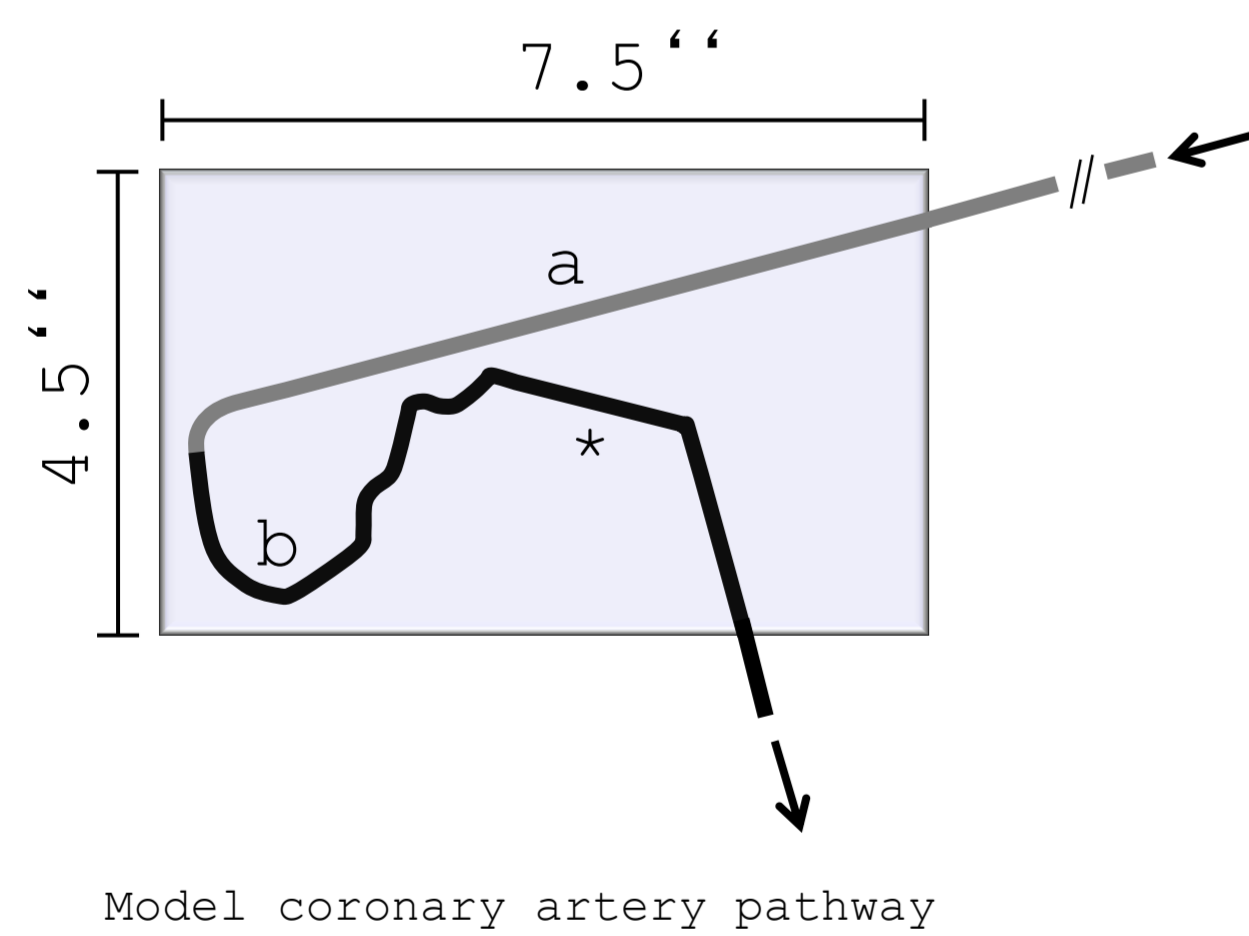
### Concept

- Estimation of potential drug loss during the passage to the site of implantation using an in vitro model
- In vitro drug release testing from a commercially available drug-eluting stent using different test methods and evaluation whether the used method influences the obtained test results

### Commercial Coroflex® ISAR stent

- Coated with Sirolimus (SIR, also known as Rapamycin)
- Release controlling agent Probucol (bioresorbable)
- Abluminal coating location

## Methods



Characteristics of in vitro test methods

method	agitation condition	media volume	media exchange
USP 7	40 dpm	10 mL	yes
USP 7	5 dpm	10 mL	yes
vial	shaking incubator	20 mL	yes
beaker	magnetic stir bar	150 mL	no
FTC	perfusion 35 mL/min	150 mL	no
vFTC	perfusion 35 mL/min	150 mL	no

### Stent examination

- Environmental scanning electron microscopy (ESEM)
- Drug elution with MeOH and SIR content determination via HPLC
- Microscopic determination of coating thickness distribution via spectral reflectometry

### In vitro estimation of potential drug losses

- Model coronary artery pathway adapted from ASTM F-2394-07
- Includes a guiding catheter (a) and a tube (b) perfused with dissolution media, flow-rate 35 mL/min
- Rapid advancement of balloon-mounted stent through perfused system
- Resting at the marked position (\*) until completion of perfusion time (5 min including advancement)

### In vitro release testing

- Use of different test apparatuses: incubation in shaken vial, incubation in stirred beaker, flow-through cell (FTC), reciprocating holder (USP 7), vessel-simulating flow-through cell (vFTC) including gelled acceptor compartment and central perfusion through stent lumen
- Dissolution media 0.9 % saline solution containing 0.05 % polyoxyethylene (23) laurylether (Brij 35) and 0.0003 % 3,5-di-tert-4-butylhydroxytoluene (BHT)
- Sink conditions in all setups

## Results

### Stent examination

- 147 ± 12 µg drug load on the stent and 154 ± 7 µg on the balloon surface
- Mean coating thickness of abluminal side 6.1 ± 3.6 µm and 1.3 ± 1.5 µm on luminal side

### Drug loss during simulated passage to site of application

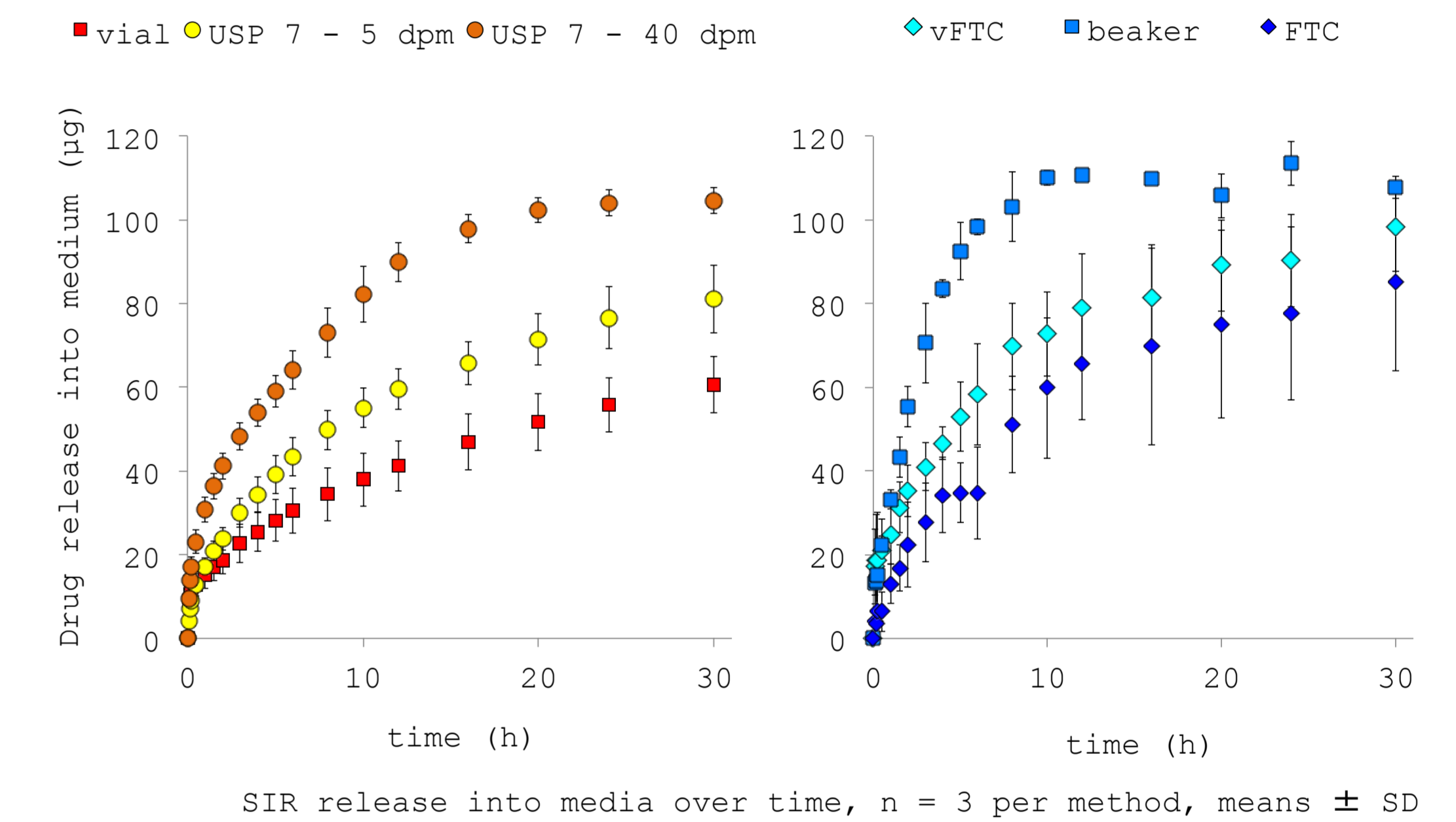
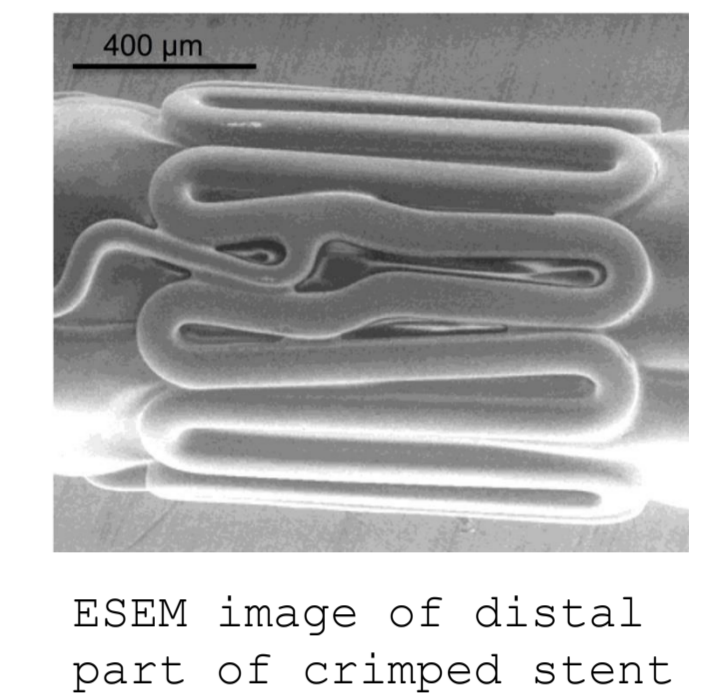
- Mean drug loss of 9.4 ± 7.9 % during simulated passage incl. 5 min perfusion

### In vitro drug release testing

- Very different results in dependency of applied test method (range of 61-108 µg released into media within 30 h)
- Fastest release upon incubation in stirred beaker
- Slowest release upon incubation in shaken vial in spite of higher volume and same sampling times as USP 7
- Distinct differences between USP 7 at 5 dpm and 40 dpm

SIR loss during simulated passage to site of application

stent #	SIR loss 1 <sup>st</sup> min (µg)	SIR loss total (µg)	SIR loss total (% of detec. amount)
1	10.4	15.7	5.6
2	28.3	47.8	18.5
3	9.0	11.9	4.1



## Conclusion

### Conclusions

- Microscopic examination and drug load distribution indicate coating of the balloon-mounted stent
- Potential drug loss during passage to site of application variable with higher loss during the advancement (abrasion forces) opposed to the resting time (dissolution)
- In vitro drug release dependent on used test method and most likely also on media (not examined here)
- Difference in release in the USP 7 in dependency of the dip rate in combination with the highest release in the stirred beaker with harsh stirring conditions may indicate a strong influence of hydrodynamics in the test system on the release profile of the tested stent system

### References

[1] Neubert et al., J Control Release 2008; 130 (1): 2-8.

### Acknowledgement

The authors thank Agilent Technologies for generously providing the USP 7 apparatus. Financial support by the European Regional Development Fund (ERDF), the European Social Fund (ESF) and the Federal Ministry of Education and Research (BMBF) within REMEDIS is gratefully acknowledged. Anne Seidlitz thanks Galenus Privatstiftung for a personal grant.

In vitro release profiles of drug-eluting stents are highly dependent on the employed in vitro test method



REMEDIS  
 IMPLANTATTECHNOLOGIE DER ZUKUNFT

SPITZENFORSCHUNG INNOVATION  
 IN DEN NEUEN LÄNDERN

Galenus Privatstiftung