

Application and modeling of a lab scale Twin Screw Extruder for the continuous mixing of Li-Ion battery slurries

Juan Meza¹, Annika Völp², Hermann Nirschl¹

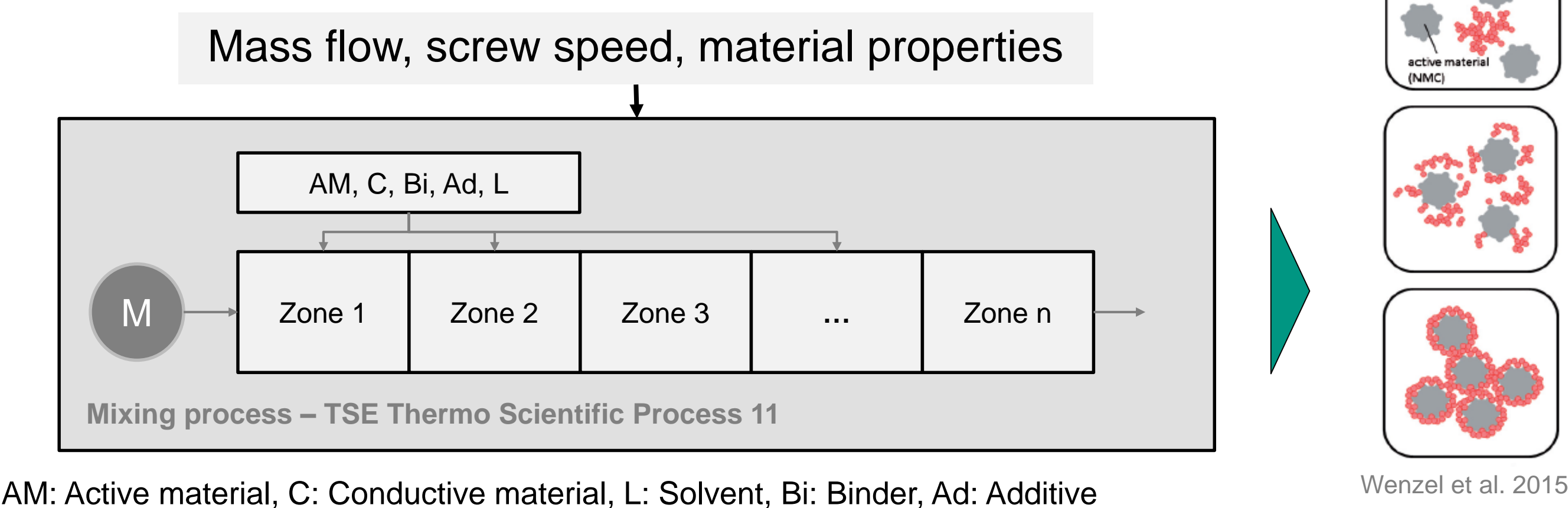
¹ Karlsruhe Institute of Technology (KIT), Institute of Mechanical Process Engineering and Mechanics (MVM), Karlsruhe, Germany

² Thermo Fisher Scientific, Karlsruhe, Germany

Motivation

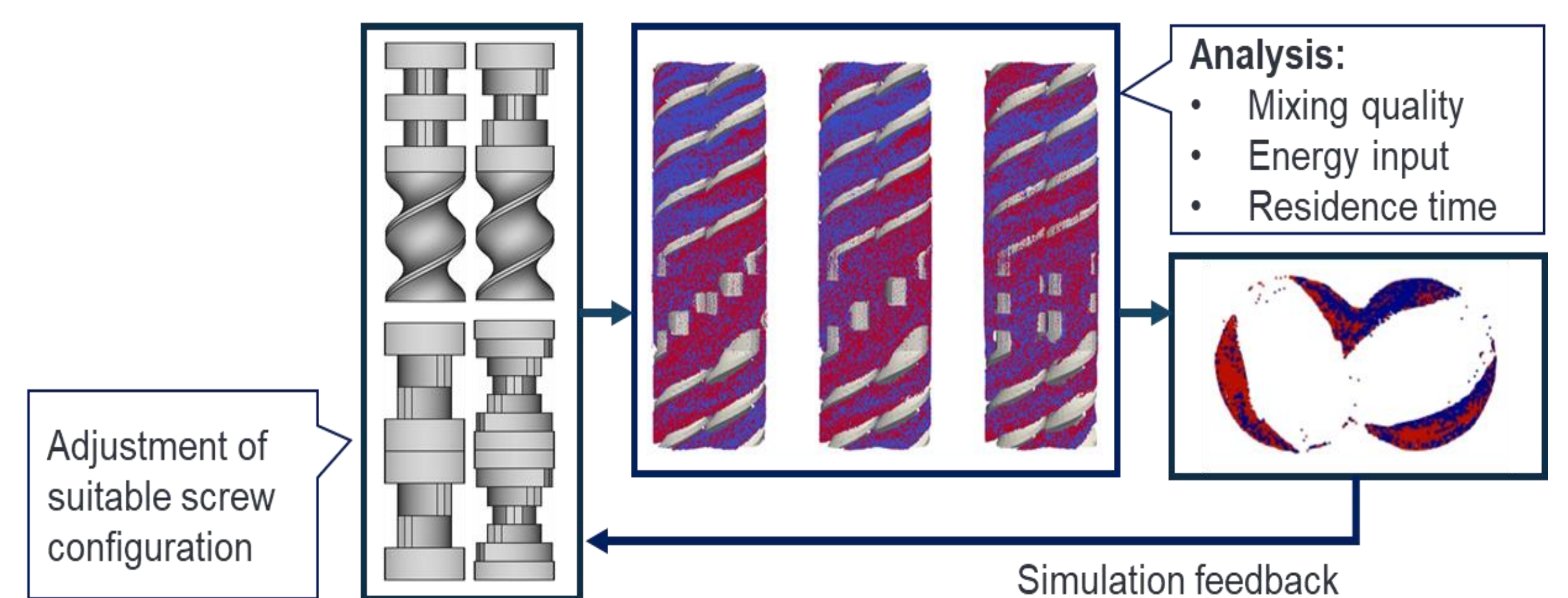
- Towards **continuous** and **agile** production of battery slurry:
 - **Versatile** dosing of slurry components during mixing
 - **Variable** slurry production for **testing** purposes or **scale-up**
 - **Reduction** of waste material after ramp up
- **Adaptive control** of production rate, temperature, energy input and viscosity

Extrusion process



Digital Twin Approach

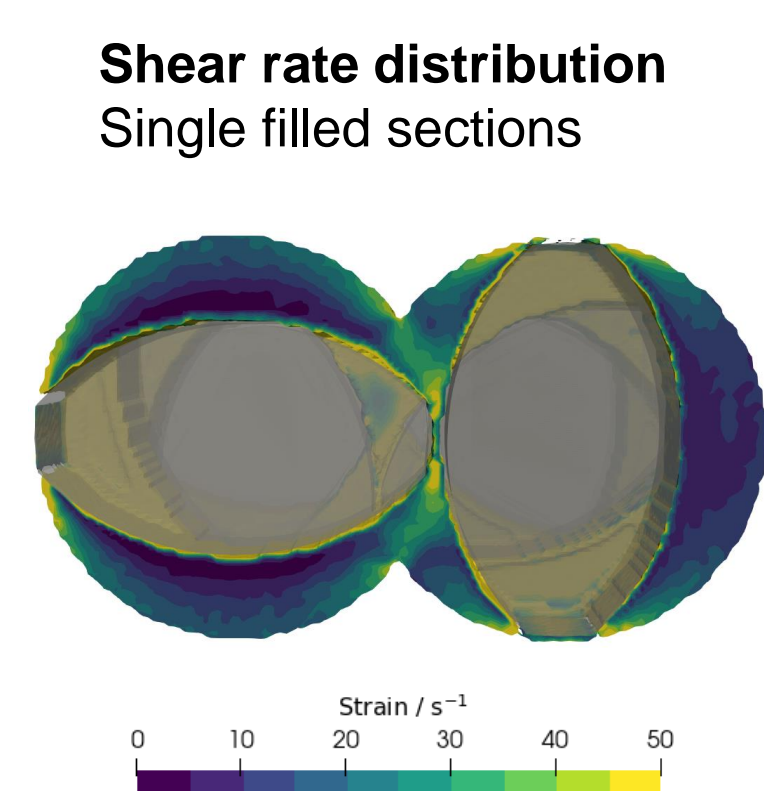
- **Simulation** based investigation of material **strain** corresponding to various process parameter variations prior to experiments
- Evaluation of feeding, kneading and conveying **performance** through screw **segmentation**
- **Iterative approach** for adaptation of process parameters and screw design to quality requirements



Results

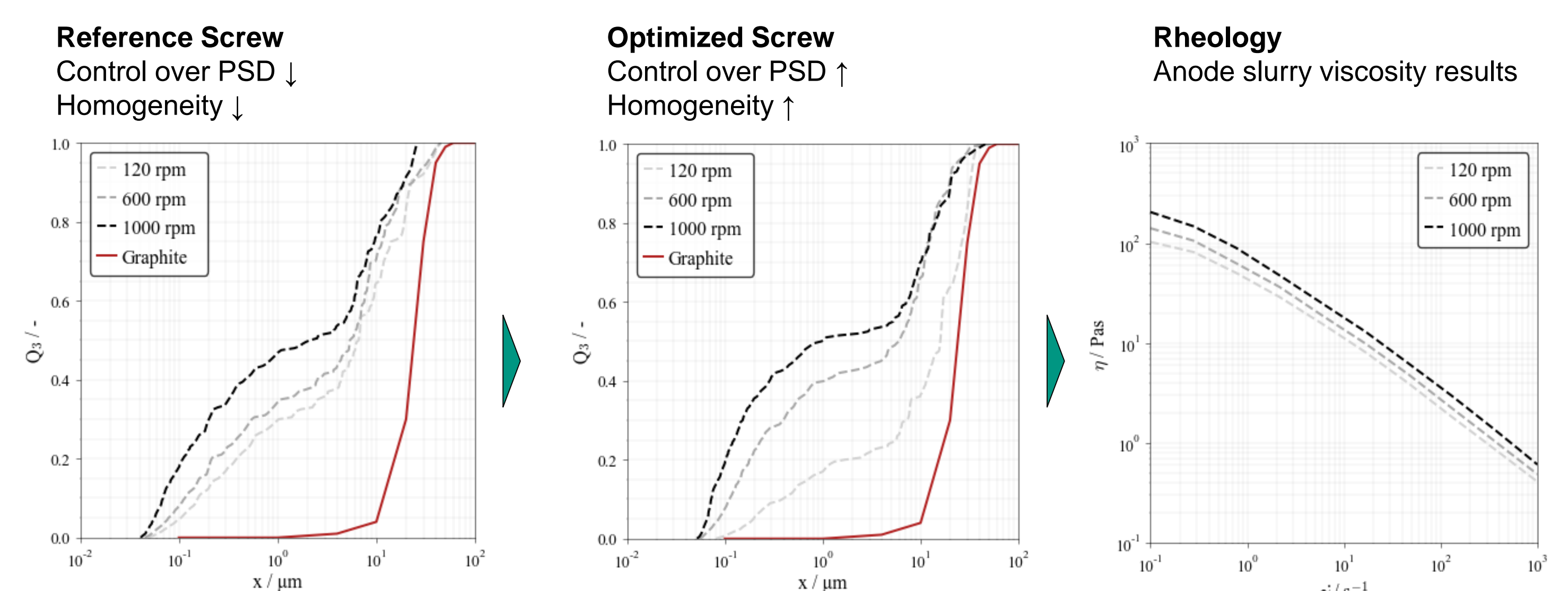
Simulation

- Analysis of the material flow behavior under experimental process parameters for efficiency evaluation
- Smoothed particles hydrodynamics (SPH) provides detailed information of slurry screw interaction
- 1D Modell enables optimization of screw configuration

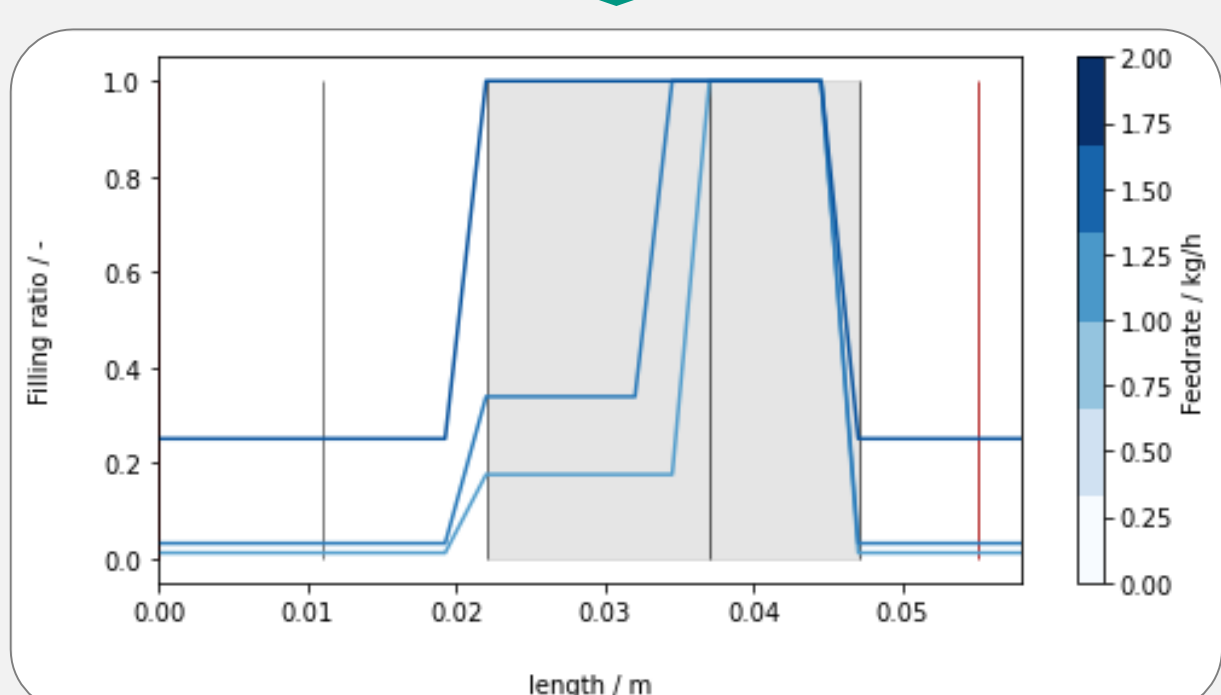
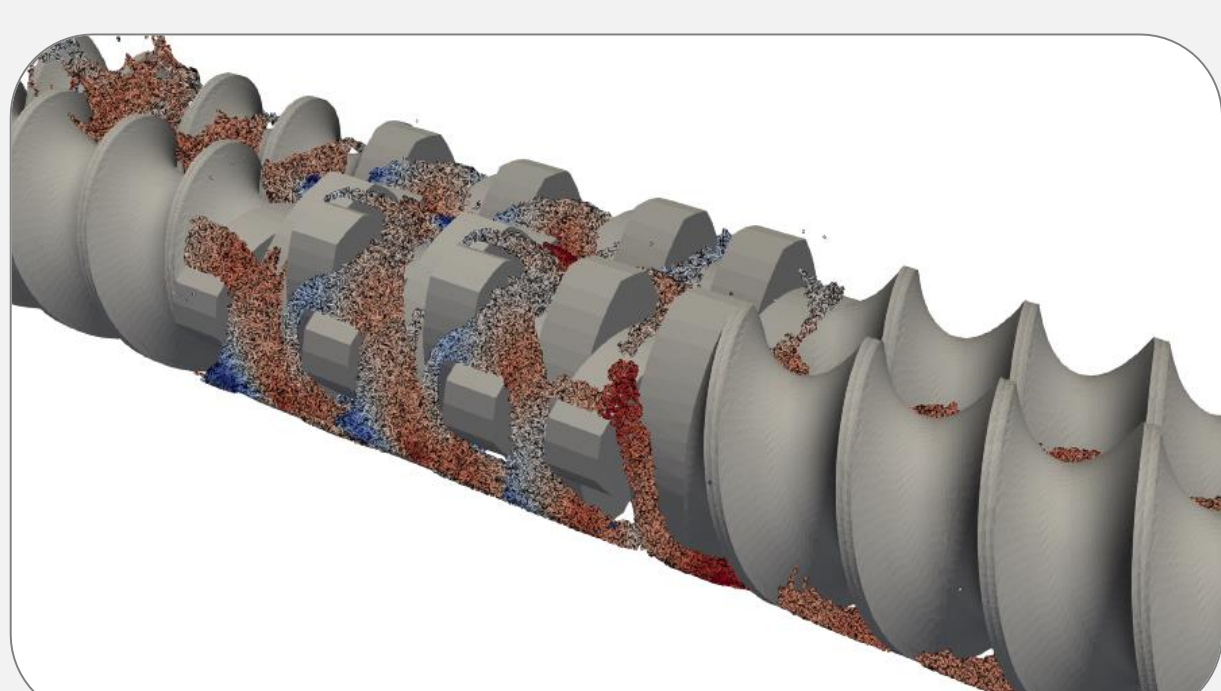
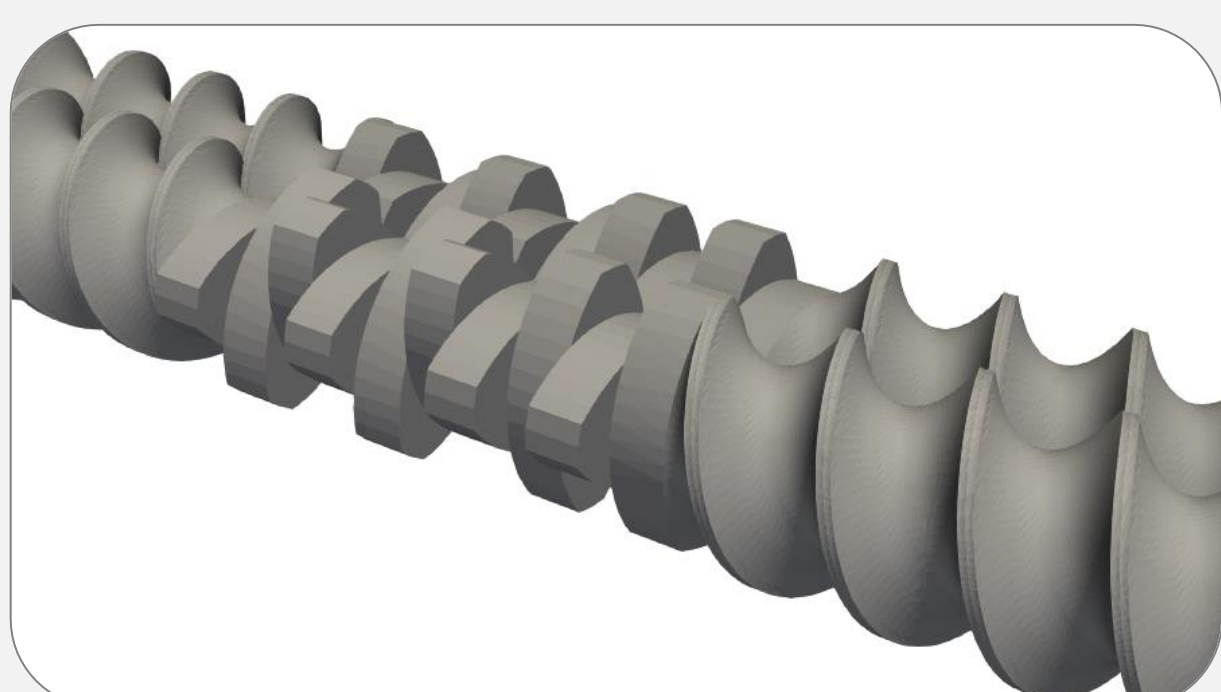


Experimental

- Analysis of the **effect of screw speed and flowrate** on slurry rheology, particle sizes (PSD) and homogeneity



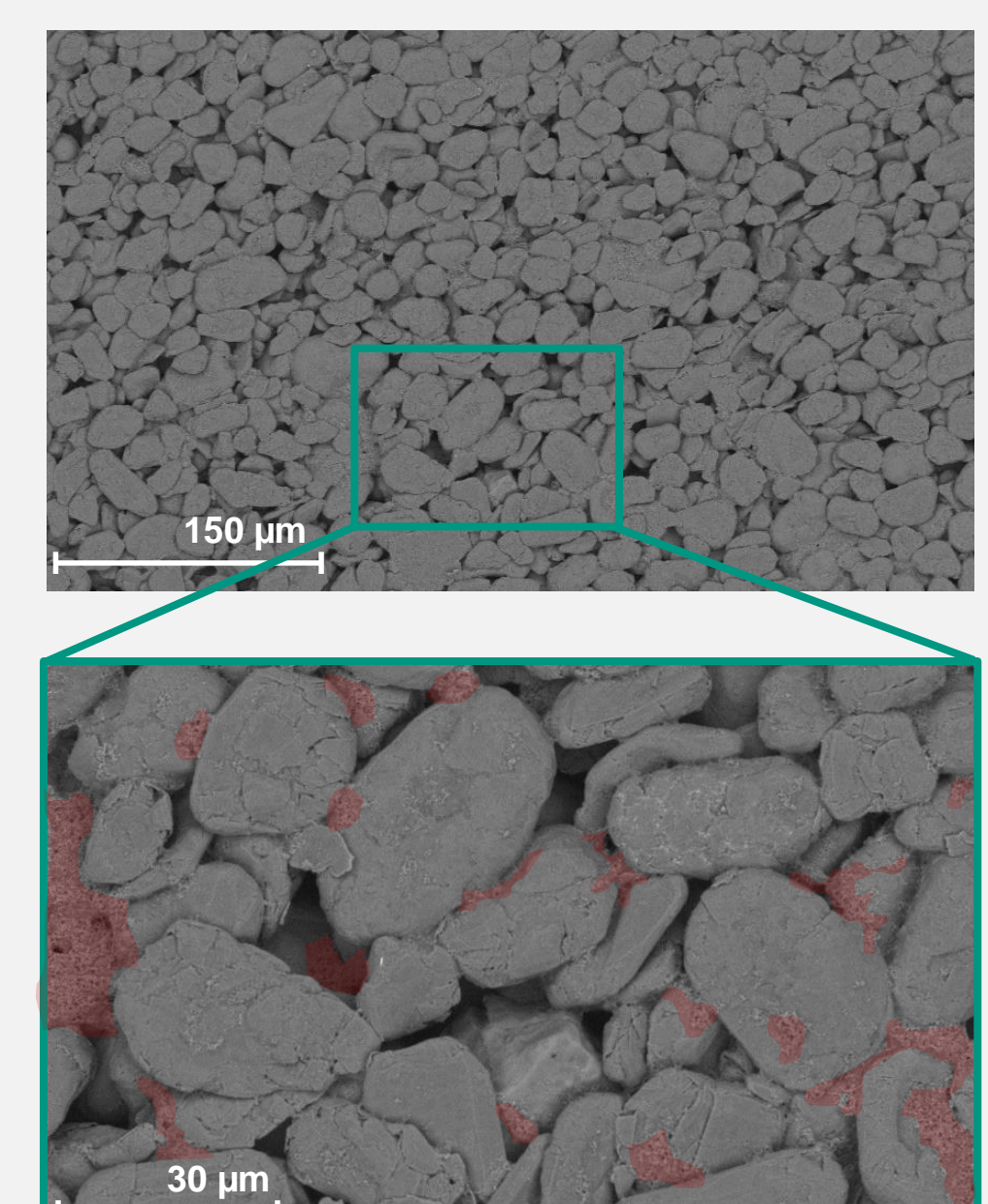
Digital Twin-Screw-Extruder (TSE)



Conclusion

- **Extrusion** process provides **continuous** slurry production
 - **Agile handling** of new recipes, slurry composition and material feasible
 - **Adaptive customization** of particle size distribution, homogeneity and stability due to variable screw design
- **Digital twin** as a key element for the **optimization** of the process
 - **Iterative modeling** supports the investigation of new screw configurations and process parameter → Waste reduction
 - Analysis of residence time, material distribution and strain feasible

Micro structures



Contact

Juan Meza, M.Sc
Karlsruhe Institute of Technology
Institute for Mechanical Process
Engineering and Mechanics
juan.meza@kit.edu

Acknowledgement

Baden-Wuerttemberg's ministry of Science, Research and Art and the Federal Ministry of Education and Research are acknowledged for funding