

# Dry amorphization of itraconazole: twin-screw processing and crystallinity

Eric Berthier, Dr. Simon Welzmler, Thermo Fisher Scientific, Chemin de Verney 2, 1024 Ecublens, Switzerland

Dr. Margarethe Richter, Thermo Fisher Scientific, Dieselstraße 4, 76227 Karlsruhe, Germany

Corresponding author: [eric.berthier@thermofisher.com](mailto:eric.berthier@thermofisher.com)

## INTRODUCTION

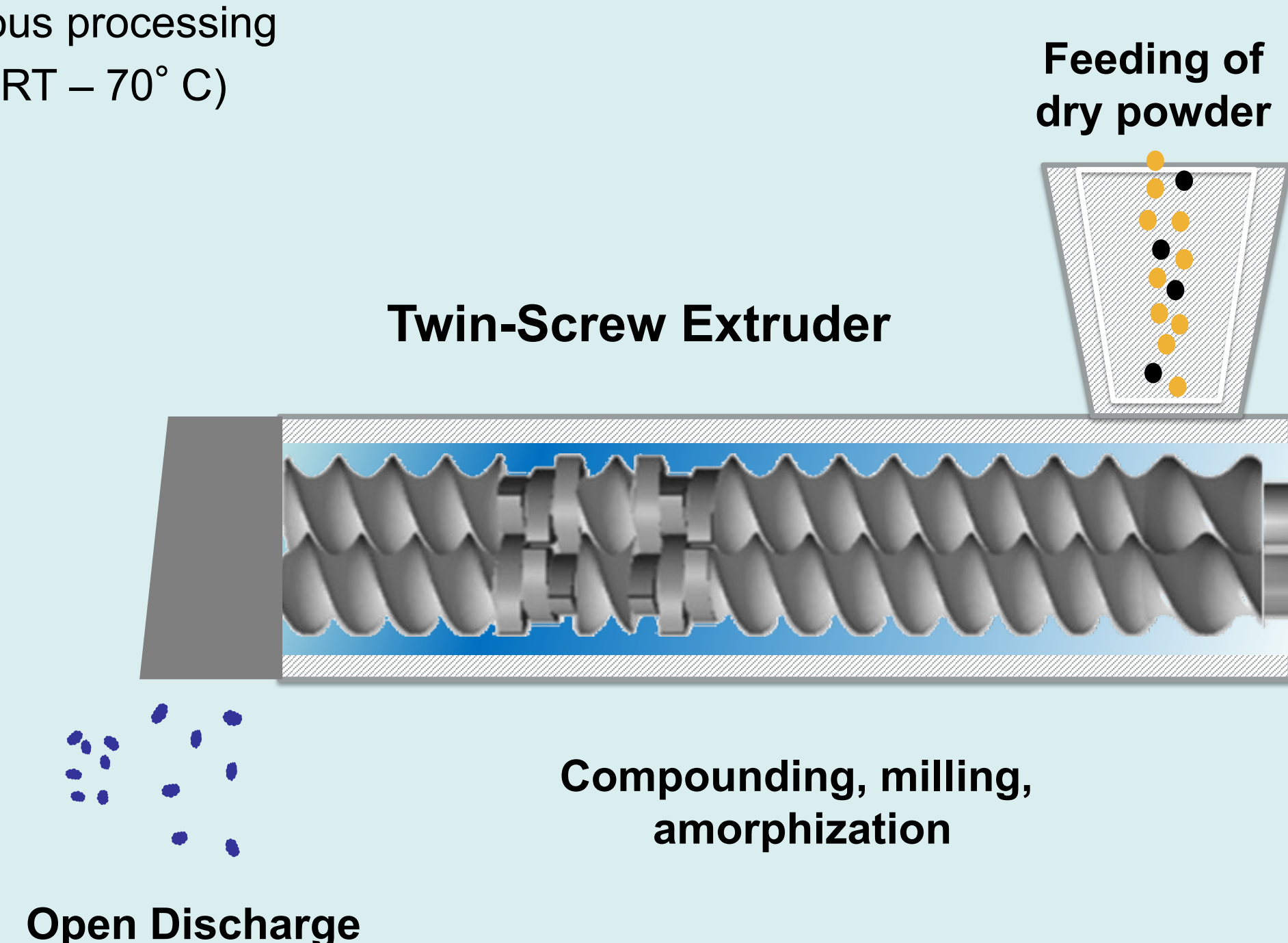
In the development of active pharmaceutical ingredients (APIs), their bioavailability is of main concern. The metastable amorphous state of APIs can reduce the dissolution time and therefore increase bioavailability, but it needs to be stabilized by a matrix material. Polymers are predominantly used to form solid dispersions with APIs and therefore stabilize the amorphous state. Solid dispersions are mostly prepared by hot melt extrusion (HME) and spray drying. HME requires temperatures above the glass transition temperature of the polymer which often can lead to decomposition or structural change of the API, whereas spray drying requires tedious handling of a solvent. An alternative route to stabilize the amorphous state of APIs is the usage of porous silica with a well-defined pore size distribution. Recent studies [1]-[3] have shown that solvent-free ball-milling of API/porous silica mixtures at moderate temperatures can improve the dissolution rate. Nonetheless, extensive milling times and batch processing is required. Twin-screw compounding technology is a mature process which can synthesize material in a constant flow with increased reproducibility and flexibility.[4][5]

In this study, itraconazole is used as a model drug which is stabilized by a mesoporous silica matrix. Processing is carried out by a laboratory sized pharmaceutical twin screw extruder with varying screw configurations and process parameters. To analyze the quality of the processed material, powder X-ray diffraction (XRD) is used to quantify the amorphous to crystalline ratio by means of standard-less combined whole pattern refinements.

## Dry Amorphization - Synthesis

Dry amorphization using continuous twin-screw technology at low temperatures

- Consistent continuous processing
- Low temperatures (RT – 70° C)
- No solvent needed



## Analytical Instrumentation

### Twin-Screw Extruder

- Thermo Scientific™ Pharma 11 Twin-Screw Extruder
- Residence time in extruder
- Torque of extruder
- Specific mechanical energy



### Differential Scanning Calorimetry (DSC)

- NETZSCH DSC 204F1 Phoenix (240-12-0023-L)
- Heating/cooling rate: 10 K/min
- 10 mg sample size



### Powder X-RAY Diffractometer (XRD)

- Thermo Scientific™ ARL EQUINOX 100
- Cu Ka Radiation @ 36W Power
- Unique CPS detector technology
- 5 min measurement in reflection



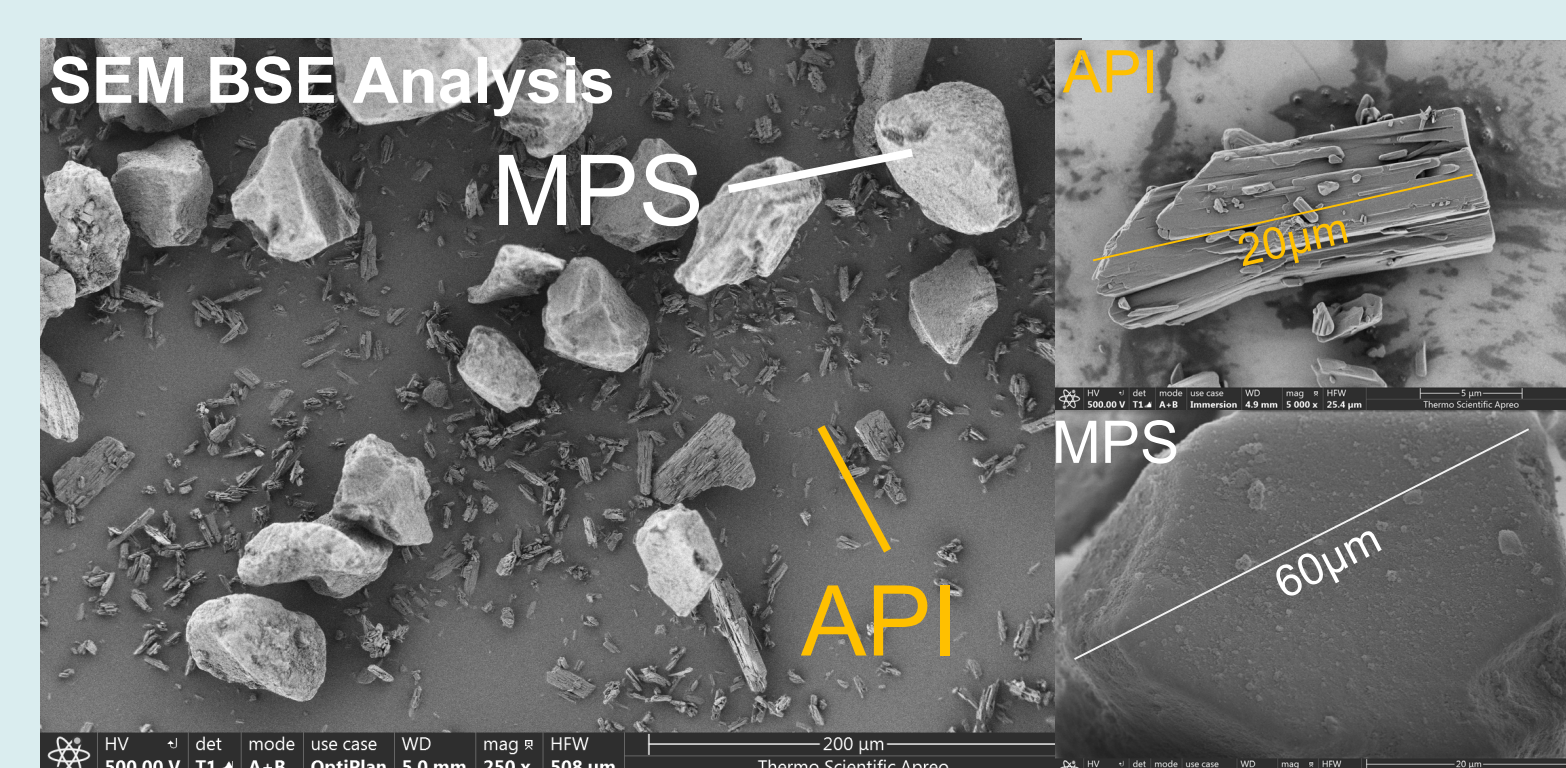
### Scanning Electron Microscope (SEM)

- Thermo Scientific™ Apreo S SEM

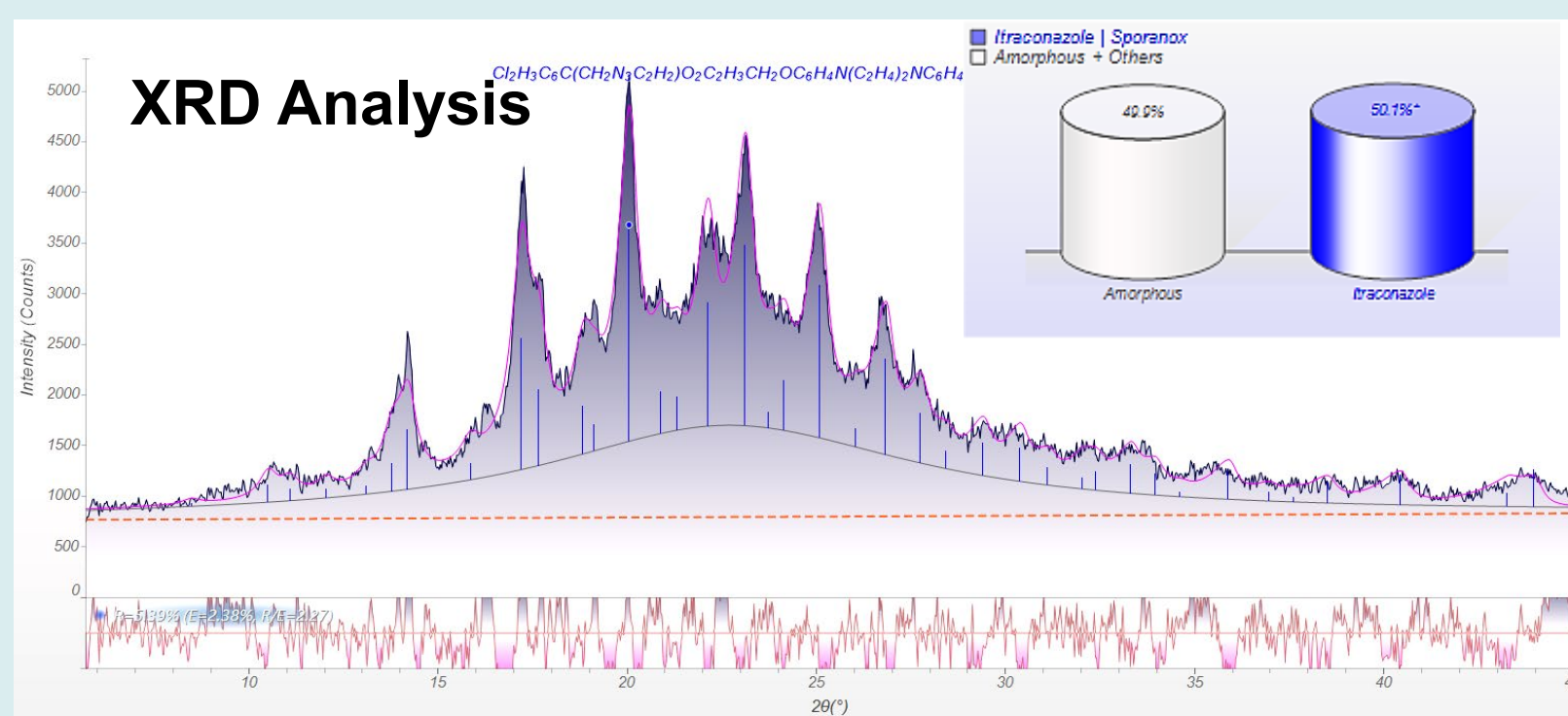


## Final Product analysis (1:1 ratio)

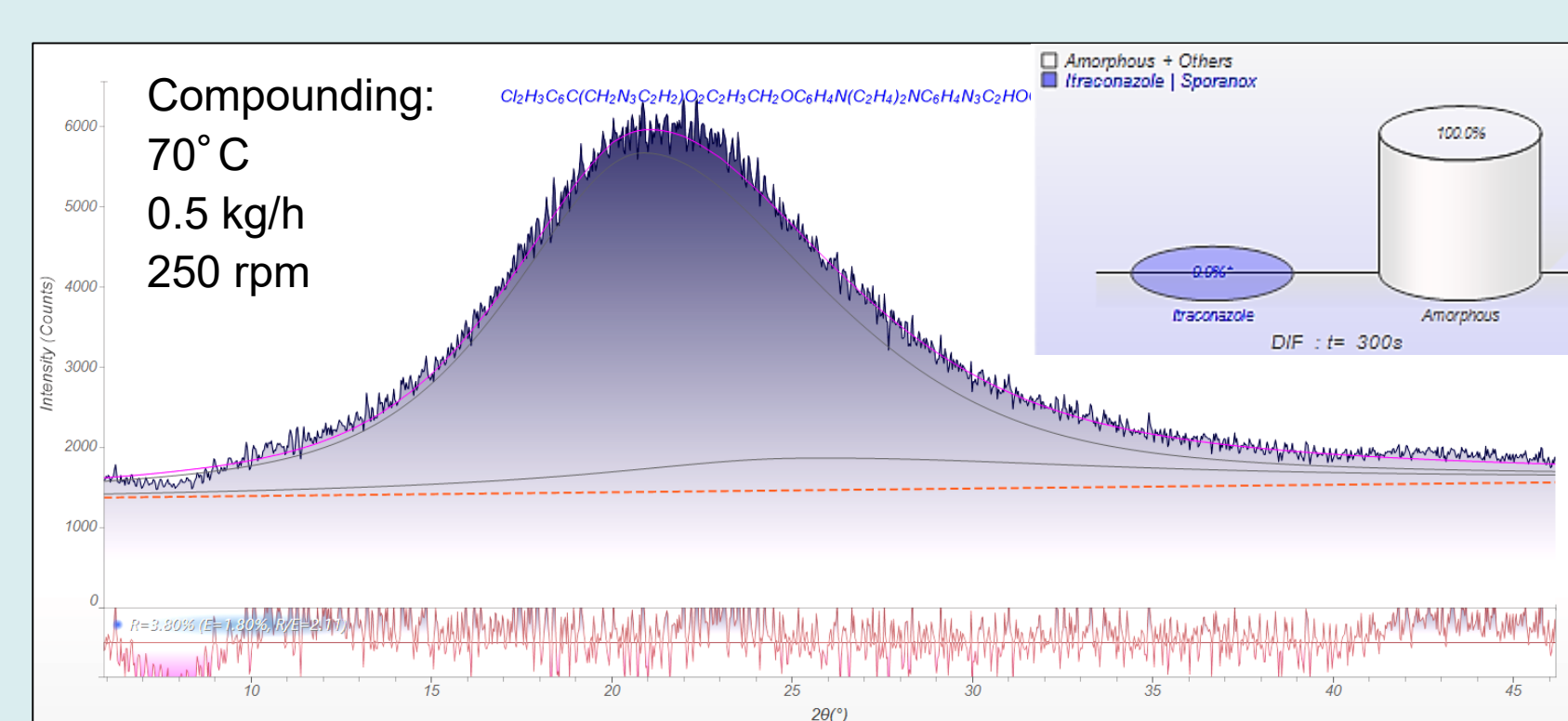
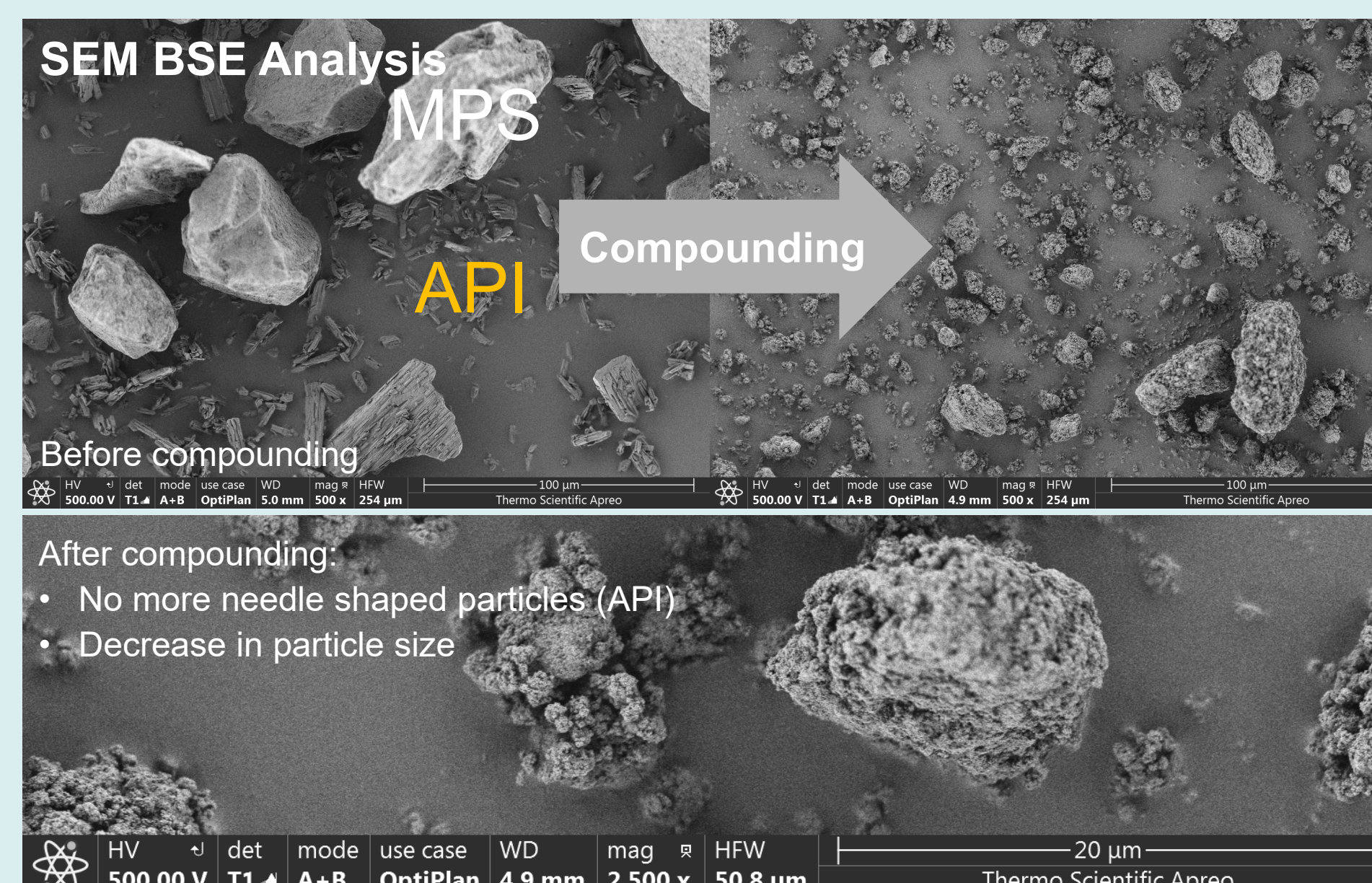
### Pre-blend analysis (1:1 ratio)



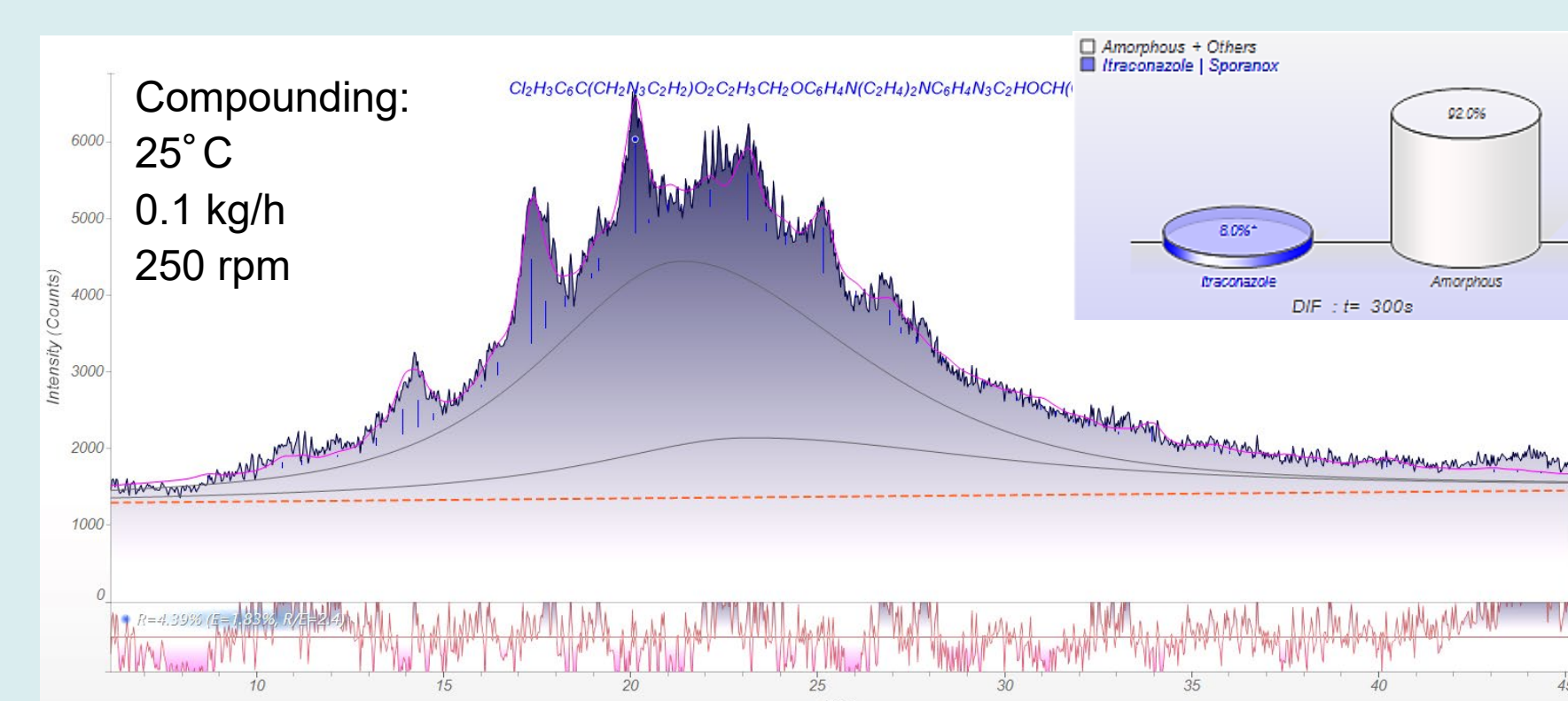
API (itraconazole): needle shaped particles  
MPS (meso-porous silica): lump like particles



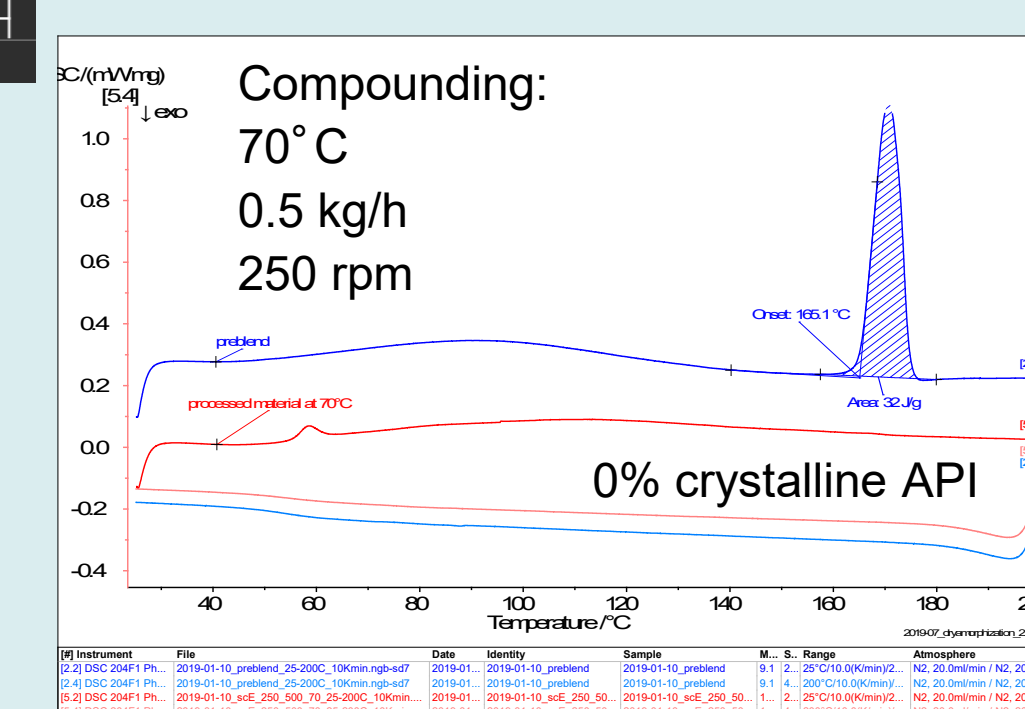
Quantitative XRD analysis of amorphous (MPS: grey) and crystalline (API: blue) components; Analysis shows 1:1 ratio and proves the principle



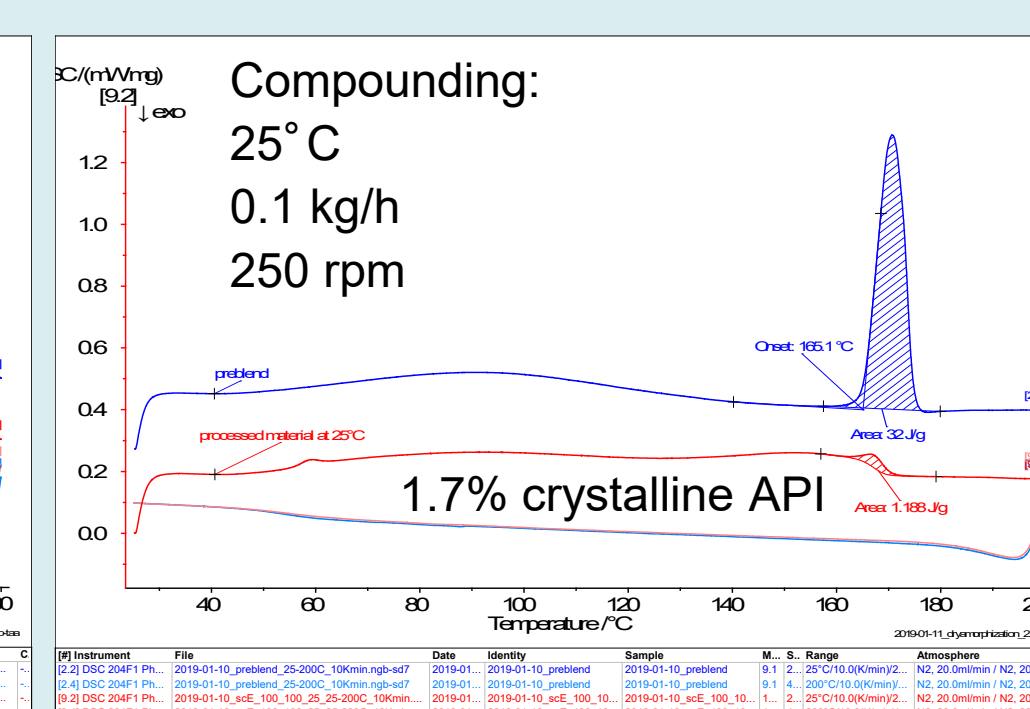
Quantitative XRD analysis of amorphous (MPS: grey) and crystalline (API: blue) components; Sample is completely amorphous



Quantitative XRD analysis of amorphous (MPS: grey) and crystalline (API: blue) components; Sample contains 8% crystalline API



DSC analysis of pre-blend (blue) and processed material (red); no melting heat in processed material indicated no remaining crystalline API phase



DSC analysis of pre-blend (blue) and processed material (red); melting heat in processed material quantifies to 1.7% remaining crystalline API phase

## CONCLUSION

- Decrease in particle size after twin-screw processing
- Full amorphization succeeded at 70° C and 175° C (according to XRD and DSC analysis)
- Up to 92% of the API in a 1:1 blend with silica reached amorphous state being processed at room temperature. (according to XRD analysis); Finding supported by DSC analysis with slightly different ratio
- Proof of concept of dry amorphization with continuous twin-screw process; No need of solvents or high temperatures
- The key factors: mean residence time and specific mechanical energy

## REFERENCES

- [1] A. Baán *et al.*, "Dry amorphisation of mangiferin, a poorly water-soluble compound, using mesoporous silica," *Eur. J. Pharm. Biopharm.*, vol. 141, no. March, pp. 172–179, 2019.
- [2] D. Bahl, J. Hudak, and R. H. Bogner, "Comparison of the ability of various pharmaceutical silicates to amorphize and enhance dissolution of indomethacin upon co-grinding," *Pharm. Dev. Technol.*, vol. 13, no. 3, pp. 255–269, 2008.
- [3] F. Monsuur *et al.*, "Solvent free amorphisation for pediatric formulations (minitables) using mesoporous silica," *Int. J. Pharm.*, vol. 511, no. 2, pp. 1135–1136, 2016.
- [4] J.C. Vervaeke and J. P. Remon, "Continuous granulation in the pharmaceutical industry," *Chem. Eng. Sci.*, vol. 60, no. 14, pp. 3949–3957, 2005.
- [5] J. Vercruysee *et al.*, "Stability and repeatability of a continuous twin screw granulation and drying system," *Eur. J. Pharm. Biopharm.*, vol. 85, pp. 1031–1038, 2013

## ACKNOWLEDGEMENTS

Joachim Quadflieg from GRACE GmbH is greatly acknowledged for scientific support of the project

**ThermoFisher**  
SCIENTIFIC