

Preliminary agenda

8:30 - 9:00 Registration	Session 1:
9:00 - 9:15 Introduction	Particle size, stability, concentration, and shape
9:15 - 10:00 Session 1	Mastersizer, SediGraph, Zetasizer, NTA, SAS, Empyrean, Morphologi
10:00 - 10:15 Break	Session 2:
10:15 - 11:00 Session 2	- Powder flow, density and porosity - FT4, AccuPyc, GeoPyc
11:00 - 11:15 Break	Session 3:
11:15 - 12:00 Session 3	Porosimetry, surface area, and structure
12:00 - 1:00 Lunch	AutoPore, AccuPore, TriStar, ASAP, 3Flex, Gemini, Empyrean, Aeris
1:00- 2:00 Demos	Session 4:
2:00- 2:45 Session 4	Activity
2:45 - 3:00 Break	AutoChem, ASAP 2020 Chemi, 3Flex Chemi, BTA, FR and ICCS
3:00 - 3:45 Session 5	Session 5:
3:45 - 4:00 Wrap-up	Elemental composition
4:00 - 4:30 Tear down & finish	Epsilon, Revontium, Zetium, SciAps