

# A powerful combination

Advancing powder materials  
characterization from particle  
to performance



## Agenda

| Time          | Subject                                                                                                                                                                                                |
|---------------|--------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| 8:30 - 9:00   | Registration                                                                                                                                                                                           |
| 9:00 - 9:05   | Introduction                                                                                                                                                                                           |
| 9:05 - 9:50   | <b>Session 1 - part 1.</b> <ul style="list-style-type: none"><li>• Particle size, stability, concentration, and shape</li><li>• MasterSizer, SediGraph, Zetasizer, SAS, Empyrean, Morphologi</li></ul> |
| 9:50 - 10:10  | Demo & Break                                                                                                                                                                                           |
| 10:10 - 10:55 | <b>Session 1 - part 2.</b> <ul style="list-style-type: none"><li>• Particle size, stability, concentration, and shape</li><li>• MasterSizer, SediGraph, Zetasizer, SAS, Empyrean, Morphologi</li></ul> |
| 10:55 - 11:15 | Demo & Break                                                                                                                                                                                           |
| 11:15 - 12:00 | <b>Session 2.</b> <ul style="list-style-type: none"><li>• Powder flow, density and porosity</li><li>• FT4, AccuPyc, GeoPyc</li></ul>                                                                   |
| 12:00 - 1:30  | Demo & Lunch                                                                                                                                                                                           |
| 1:30 - 2:15   | <b>Session 3.</b> <ul style="list-style-type: none"><li>• Porosimetry, surface area, and structure</li><li>• AutoPore, AccuPore, TriStar, ASAP, 3Flex, Gemini, Empyrean, Aeris</li></ul>               |
| 2:15 - 2:35   | Remote Demo                                                                                                                                                                                            |
| 2:35 - 3:20   | <b>Session 4.</b> <ul style="list-style-type: none"><li>• Activity</li><li>• AutoChem, ASAP 2020 Chemi, 3Flex Chemi, BTA, FR and ICCS</li></ul>                                                        |
| 3:20 - 3:30   | Break                                                                                                                                                                                                  |
| 3:30 - 4:15   | <b>Session 5.</b> <ul style="list-style-type: none"><li>• Elemental composition</li><li>• Epsilon, Revontium, Zetium, SciAps</li></ul>                                                                 |
| 4:15 - 4:30   | Demos                                                                                                                                                                                                  |
| 4:30          | Wrap up                                                                                                                                                                                                |