



## PARTICLE COUNTING AND PARTICLE CHARACTERIZATION

From techniques spanning the Coulter principle, light scattering and laser diffraction to polarized intensity differential scattering (PIDS), we offer a variety of training options and formats designed to deliver the real-world knowledge and technical know-how you need to optimize your equipment investment.

Visit [beckman.com/training](http://beckman.com/training) to view all training options and register for courses.

For training inquiries or customized training requests, email us at [beckmancoulterlearningcenter@beckman.com](mailto:beckmancoulterlearningcenter@beckman.com)



© 2018 Beckman Coulter Life Sciences. All rights reserved. Beckman Coulter, the stylized logo, Biomek, Agencourt and AMPure are trademarks of Beckman Coulter, Inc. and are registered with the USPTO. All other trademarks are the property of their respective owners.

For worldwide office locations and contact detail, visit Contact Us at [beckman.com](http://beckman.com)

MULTI-3365FLY02.18