

William Kang, Thermo Fisher Scientific, Carlsbad, CA, 92008, USA

© 2026 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. CyTek™ Aurora and SpectroFlo™ are trademarks of CyTek Biosciences. Brilliant™ is a trademark of Becton Dickinson and Company and is under license. Cy™ is a trademark of GE Healthcare. This information is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others.