

Find out more at thermofisher.com/nalgenelabware

thermoscientific

© 2018, 2025 Thermo Fisher Scientific, Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. Alconox is a registered trademark of Alconox, Inc. Ficol-Hypaque and Percoll are registered trademarks of GE Healthcare Bio-Sciences AB. Nucleomix is a registered trademark of QIAGEN Laboratories, Inc. Tiltrol is a registered trademark of The Dow Chemical Company. Tween is a registered trademark of ICI Americas, Inc. **STU010-25-1321**