

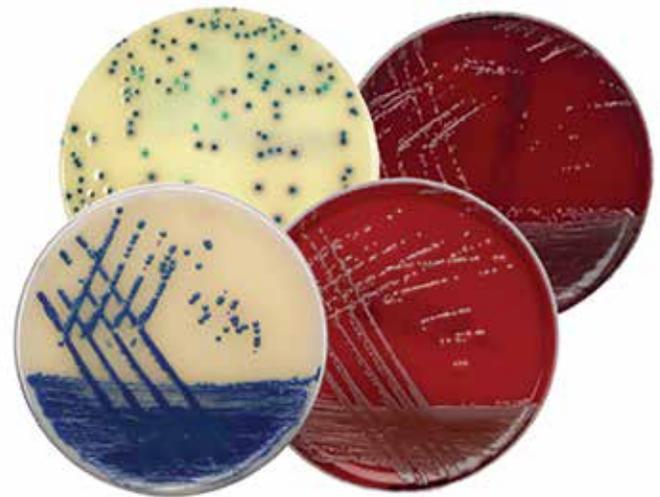
# Guide to optimizing the operation of Thermo Scientific prepared media with automated plate processing systems

Thermo Scientific™ SmartPlate media format for Thermo Scientific™ prepared plate media was recently introduced in North America. The format brings several benefits including reducing impact on the environment due to a reduction in the amount of resin used to manufacture the Petri dish and lighter weight of the finished product. The SmartPlate media format is also designed to be compatible with a wide range of automated plate processing systems.

To optimize the operation of the SmartPlate media format with your automated plate processing equipment some small adjustments will need to be made. This is due to very small variations in the dimensions of plates from different suppliers. If you are utilizing an up-front plate processing system such as the Copan WASP™ Instrument or the BD™ Kiestra™ System, we recommend that you consult your internal or contracted service teams to assist with making the necessary adjustments.

To confirm full optimization of the SmartPlate media format with your equipment, and to prevent avoidable delays to your operation, it is advisable to arrange for a service engineer to be on-hand at the point of the prepared plate media evaluation stage, ahead of a full conversion and day-to-day operation.

To find out more about converting to Thermo Scientific Prepared Media visit [thermofisher.com/clinical-ppm-switch](https://thermofisher.com/clinical-ppm-switch)



In support of this process, samples of Thermo Scientific prepared plate media can be supplied. Please speak to your local representative or get in touch using the contact information below.

© 2020 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. WASP™ is a trademark of Copan Italia SPA. BD™ and Kiestra™ are trademarks of Becton, Dickinson and Company.

**Contact Information:**

microbiology@thermofisher.com  
USA +1 800 255 6730  
International +44 (0) 1256 84114

LT2516A  
February 2020

**ThermoFisher**  
SCIENTIFIC