

Receiving your new media

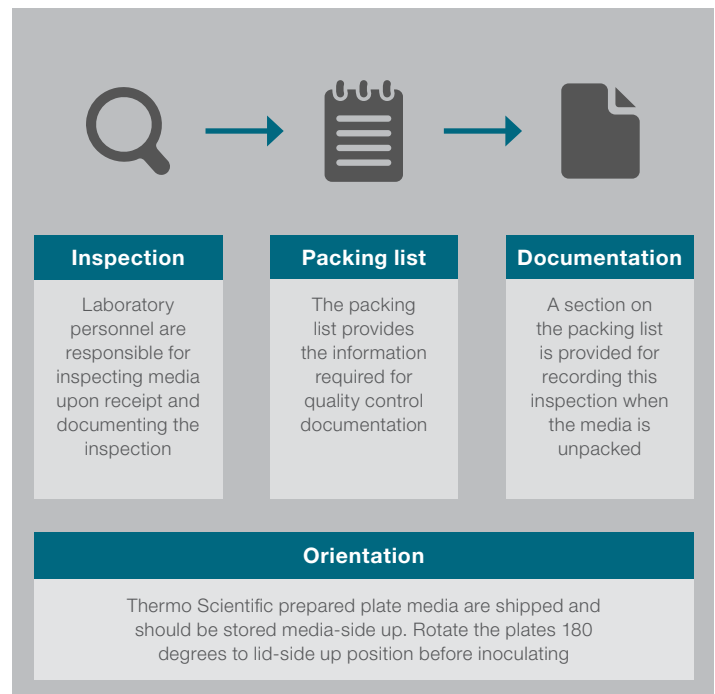
Receiving products from a new supplier can be a challenging thing if you don't know what to expect. Here are a few key reminders of what to expect and do when you receive your new Thermo Scientific™ prepared plate media.

Orientation

Thermo Scientific prepared plate media are shipped media-side up to protect the media from condensation and settling particulates during shipping. Continue to orientate prepared plate media this way during storage and on the bench until you are ready to use it. Turn the plates lid-side up before inoculating then back to media-side up for incubation.

Receiving

When you receive your Thermo Scientific prepared plate media ensure that an appropriate laboratory team member inspects the media and documents to ensure everything is present and correct. Inside each package you'll find a packing list that will describe everything that you'll need to know for your quality control documentation. Finally, on the packing list, there will be a signature line for the laboratorian to record their initial inspection.



To find out more about converting to Thermo Scientific Prepared Media visit thermofisher.com/clinical-ppm-switch

© 2020 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

Contact Information:

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

LT2516A
March 2020

ThermoFisher
SCIENTIFIC