

# Receiving your new media

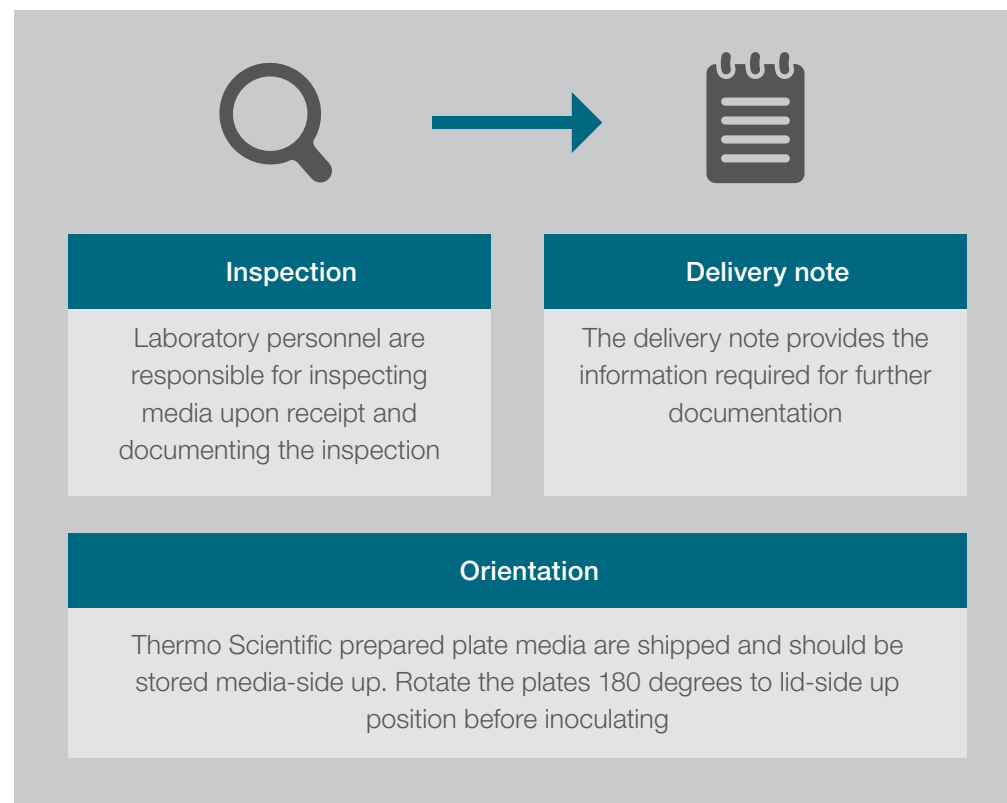
Receiving products from a new supplier can be a challenging thing if you don't know what to expect. Here are a few key reminders of what to expect and do when you receive your new Thermo Scientific™ prepared plate media.

## Orientation

Thermo Scientific prepared plate media are shipped mediaside up to protect the media from condensation and settling particulates during shipping. Continue to orientate in this way during storage and on the bench until use. 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. On each shipping box 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](https://thermofisher.com/clinical-ppm-switch)