



thermo
scientific

Everything we have ever done has been focused on one thing:

**empowering the people who dedicate
their lives to microbiology.**

Those who diagnose against the clock, around the clock, research against the odds, and uncover the incredible. All of these people deserve uncompromised quality and impeccable service.

And they need a partner with the drive to accelerate innovation, enhance productivity and help shape the future. This is what we promise, and what we strive to deliver.

From the gold-standard in culture media that laboratories have relied on for decades, to the game-changing workflows of tomorrow. Or the latest testing technologies, to supply chain optimization.

And what's driving us?

**We want to help laboratories work smarter.
To enable microbiologists to focus on what
truly matters – timely, accurate diagnosis.**

When that happens, outcomes improve and we enable our customers to make the world healthier, cleaner and safer. In vital areas of microbiology, we are the catalyst for empowering your expertise.

The world leader in serving science

ThermoFisher
SCIENTIFIC

your expertise, **empowered™**

Everything we have ever done has been focused on one thing:

empowering the people who dedicate their lives to microbiology.

Those who diagnose against the clock, around the clock, research against the odds, and uncover the incredible. All of these people deserve uncompromised quality and impeccable service.

And they need a partner with the drive to accelerate innovation, enhance productivity and help shape the future. This is what we promise, and what we strive to deliver.

From the gold-standard in culture media that laboratories have relied on for decades, to the game-changing workflows of tomorrow. Or the latest testing technologies, to supply chain optimization.

And what's driving us?

We want to help laboratories work smarter. To enable microbiologists to focus on what truly matters – timely, accurate diagnosis.

When that happens, outcomes improve and we enable our customers to make the world healthier, cleaner and safer. In vital areas of microbiology, we are the catalyst for empowering your expertise.

