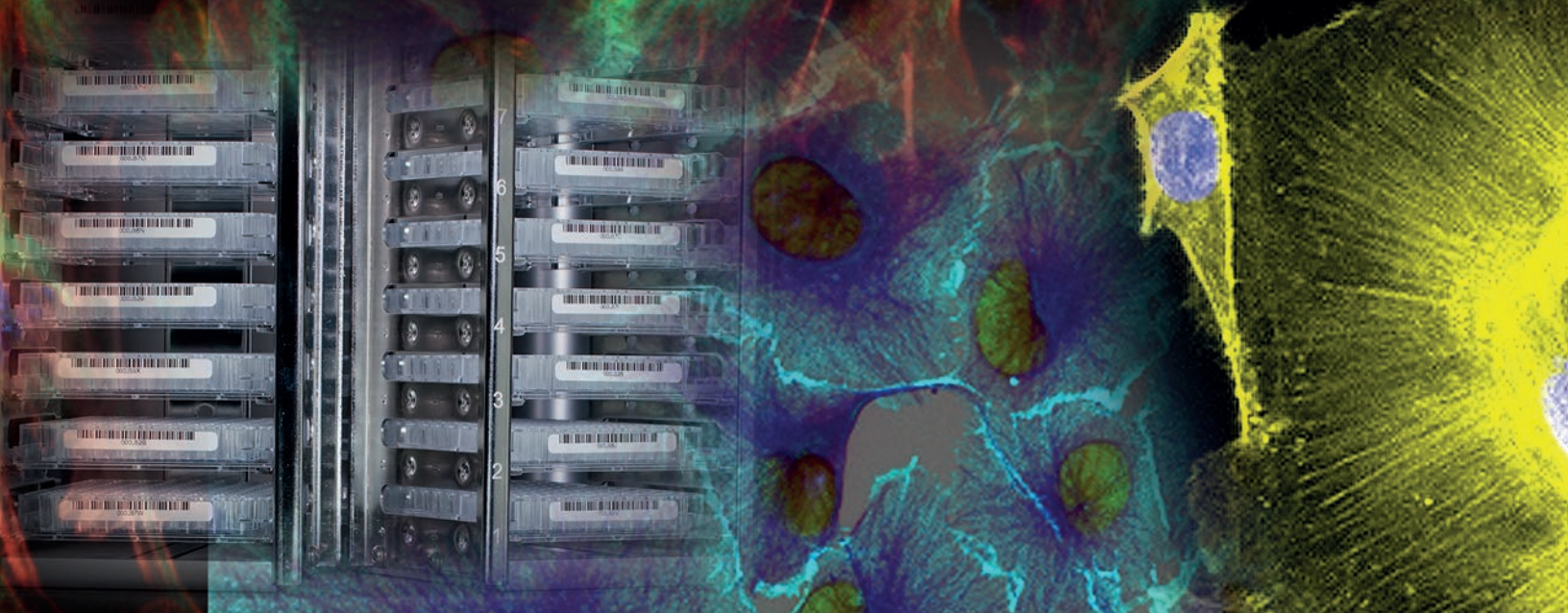


thermo scientific

Dependable sample protection
Super-fast access



Thermo Scientific™ Cytomat™ 10 Series

Automated incubators and storage systems

The next stage of advanced incubation and storage solutions, providing super-fast access, a wide temperature range and dependable sample protection

ThermoFisher
SCIENTIFIC

Dependable sample protection

High-capacity, Incubation and Storage Solutions

No compromises

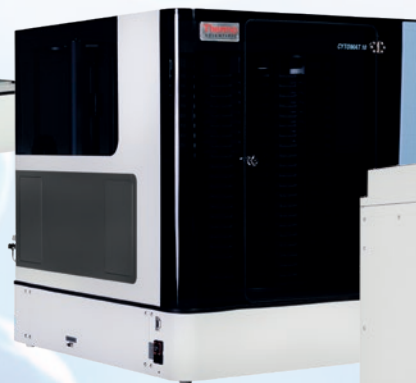
The Thermo Scientific™ Cytomat™ 10 Series provides ideal solutions for high-capacity incubation and storage, combining superior environmental conditions with super-fast access and reliable sample protection.

The range of automated incubators and ambient storage offer tightly controlled temperature regulation, as well as the flexibility to adapt to different assay and culture conditions.

- High-capacity, immediate access
- Fully automated decontamination routine
- Controlled humidity
- Wide temperature range available, from 4 to 50 °C
- Smooth and precise handling



Cytomat 10 Hotel



Cytomat 10 Hotel with
Perspex Housing

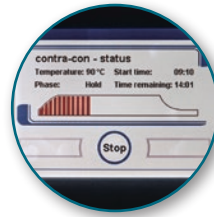


Cytomat 10 C/C450
Automated Incubator



Unrivalled Speed

The fastest incubators and hotels in their class, with a mean access time below 10 seconds. In addition, individual speed settings per plate and adjustable incremental speed settings are possible.



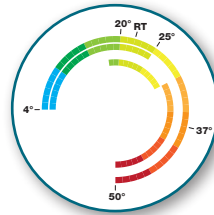
Assured Decontamination

Thermo Scientific Cytomat 10 incubators have a fully automated ContraCon thermal decontamination routine, well established in over tens of thousands of HeraCell incubators throughout the world.



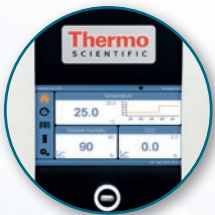
Precise Humidity Control

HydraSmart technology for controlled humidity and an external water reservoir for maximum convenience and minimal contamination risk.



Expanded Temperature Range

Incubators and storage solutions covering a range of temperatures, and including the fastest available wide temperature range unit (4 - 50 °C) in its class.



Easy Everyday Use

- LED glove-friendly display with intuitive operation
- Read-out messaging and warning alerts
- Two display sizes (5" and 7")



Flexible Solution

- Suitable for bench top and under bench
- Multiple loading/unloading location options



Cytomat 10 Hotel



Cytomat 10 Hotel with Perspex Housing



Cytomat 10 C/C450 Automated Incubator

Flexibility and versatility for all solutions

Options and accessories

Wide range of transfer stations



Standard transfer station

Supplies the microplate in portrait format. A 90° turnable version is also available for presentation in landscape format.



Swap station (180°)

180° turnable swap station to accelerate your handling process.



Linear transfer station

Available in different lengths (300, 400 and 500 mm) for extended reach. Delivers the microplate in portrait format, with further options for landscape presentation or lateral movement.



Plate lift for under-bench solution

Stepper motor driven (SMD) technology offers microplate movement to a predefined height within a 500 mm stroke.

Many more options available

General options

- Barcode reader
- Interface option for fast and easy integration
- Various stacker pitch sizes for different plate heights
- Customized stackers for different labware

Specific options

- O₂ control (1-21% or 5-90%)
- hH – high humidity up to 98% rH at 37 °C*
- CyMon – remote monitoring application, including 10" tablet (for NGD only)
- Door Left hinges (DL)
- Copper inner chamber (Cu)
- H₂O₂ disinfection ports (GDF)
- Dual Access Point (DAP) for Cytomat 10 Hotels
- Linear Shaking (Agitation) 1-100 mm amplitude, 12-300 rpm for Cytomat 10 C450 only*

* Condensate formation is possible

One product series for every application

Developed by the inventors of microbiological incubators and taking advantage of over two decades of unequalled experience on automation, the Cytomat 10 series offers the broadest set of capabilities for incubation and storage. It delivers robust, consistent environmental conditions, with minimal effort and the fastest access times on the market.

The Cytomat 10 series is the solution for all applications - from random ambient, through low temperature and room incubation, to 37 °C incubation and hybridization, and beyond.

Model	Controlled temperature range	Low temperature storage	Ambient random access storage	Room temperature incubation	Biochemical assays	Cell Assays	Mammalian and microbial cell incubation	HTS	HCS	Short term incubation
Cytomat 10 Hotel	Ambient		●	●	●			●	●	◐
Cytomat 10 C	(Ambient + 10 °C) - 50 °C			◐	●	●	●	●	●	●
Cytomat 10 C425	4 - 25 °C	●		●	●			●	●	●
Cytomat 10 C450	4 - 50 °C	●		●	●	●	●	●	●	●

◐ Capability
● Good capability

Stacker Options

Plate Type	Stacker Version	Max. Plate Height	Capacity
1536 well-plates	Stacker P17	12 mm	280
96 or 384 well-plates	Stacker P23	18 mm	210
Transwell or insert plates	Stacker P33	28 mm	150
Deep well-plates	Stacker P50	45 mm	100

Thermo Scientific Cytomat 10 series

Discover your advanced incubation and storage solution today

Mean access time below 10 seconds

Automated decontamination routine with ContraCon

Controlled humidity through Hydra Smart technology

Wide temperature ranges available



Find out more at thermofisher.com/Cytomat

ThermoFisher
SCIENTIFIC