


# On your iPad Only Go to the App store and Download the ambr 250 high throughput iBrochure App

<https://itunes.apple.com/gb/app/ambr-250-high-throughput-ibrochure/id952362252?mt=8>




**ambr 250**  
high  
throughput

**ambr 250 high throughput iBrochure** 17+

Invention Interactive Limited


Free

## iPad Screenshots



**sartorius stedim**  
BIOTECH


**ambr® 250 high throughput**  
Fully automated single-use bioreactor system



turning science into solutions

AWARDED WINNER  
Best Collaborator Award  
March & Oct 16 & 16th Innovations

AWARDED WINNER  
Best Technology Application - (Sartorius Award)  
ambr 250 automated micro bioreactor system




**Introduction**  
ambr 250 high throughput is a bioreactor system for parallel fermentation or cell culture.


The system, developed with two major industrial partners, builds on our successful micro bioreactor technology pioneered in the original ambr microscale system.

ambr 250 high throughput takes ambr technology to the next level with a platform that adds increased working fermentation volumes as well as individual bioreactor temperature and impeller control. This enhanced environment supports the culture requirements for both high demand microbial fermentation and applications where larger sample volumes are required.

ambr 250 high throughput provides an efficient and rapid high throughput scale-down system at 100-250ml volumes to explore a wide range of conditions and strains with scalability to 30L bench top stirred tank bioreactors.



**ambr 250 high throughput video gallery**  
The video clips demonstrate 'easy connect' set up, run mode and automated processing capabilities



Tap the thumbnail images to learn more about ambr 250 high throughput's key features

1. ambr 250H overview
2. Easy set up
3. Sending new vessel
4. Run mode
5. Inlet addition
6. Sampling from a vessel
7. pH calibration
8. Scale-down model