

Accelerating Cell Line Development

Increasing Efficiency by Combining Ambr[®] 15
Cell Culture with Octet[®] Titer Measurements

A Powerful Duo for Process Optimization

- Screen multiple cell lines or clones in parallel
- Reduce experimental costs for media and feeds
- Perform process analytics on crude samples, enabling faster turnaround of titer analysis
- Together, these instruments speed your time to answers so you can select the best clone faster



New

Ambr® 15 Cell Culture

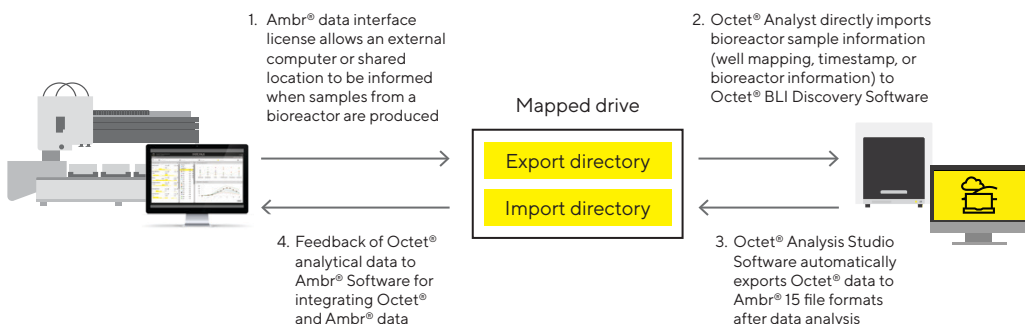
The Ambr® 15 Cell Culture is an automated, high throughput single-use bioreactor system that provides parallel operation of up to 48 single-use microbioreactors at a time and can be managed by a single operator.

Octet® BLI

Octet® systems operate on the label free, non-fluidic based BLI technology with Dip and Read assay format, that utilizes micro-titer plates and propriety biosensors coated with protein ligands to enable specific binding between the ligand and the relevant binding partner. This can be done with up to 96 samples in parallel.

The Cell Line Development workflow benefits from utilizing the Octet® BLI platform alongside the Ambr® 15 Cell Culture system with the Ambr-Octet software bridge, allowing for easy and rapid cellular product quantification. This combination allows rapid identification of the best combinations of media and clones based on product titers and cell specific productivity.

Figure: *Ambr® 15 and Octet® Workflow Utilizing the New Data Interface for Automated Data Exchange Between the Two Platforms.*



Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Corporation
565 Johnson Avenue
Bohemia, NY 11716
Phone +1 631 254 4249
Toll-free +1 800 635 2906

For further contacts, visit
www.sartorius.com/cld