

MYCAP® CCX is a reliable aseptic cell culture expansion technology. This closed system features a revolutionary design to ensure the media feed, inoculation, sampling and transfers are done aseptically, outside a biosafety cabinet. The MYCAP® CCX cell expansion QApp software controls peristaltic pumps for automated transfer of media, inoculum or media & inoculum. The software automatically calculates the volume(s) to be transferred, the final volume and effective cell density.

## MYCAP® CCX Processing



### Automated Transfer of Media or Inoculum to Cell Cultures

The Mycap® CCX system allows for expansion of suspension cell culture from cell banks to seed bioreactor. The expansion process is performed through multiple passages of successively larger Erlenmeyer flasks and involves the aseptic transfer of media and | or inoculum between donor and recipient flasks. Using the Mycap® CCX system for cell culture expansion eliminates the need to work in a biosafety cabinet for media or inoculum transfer since the flask's cap need not be removed for fluid transfer during passaging or sampling.

The Cubis® II software application for Mycap® CCX processing facilitates the transfer of media or inoculum between Mycap® Erlenmeyer flasks using peristaltic pumps. During the pumping process media or inoculum is transferred from a donor flask into a recipient flask and the transferred amount is gravimetrically measured. The Mycap® CCX processing application controls Watson-Marlow pump models 323Du, 530Du and 630Du via serial communication and offers a dedicated pumping process for the liquid transfer.

In general, the parameters for the Mycap® CCX expansion of cell cultures are organized in the software application in so-called “experiments”. General experiment settings are determined by users with the “Create Task” function, and process-specific experiment settings are determined by the user before the samples are processed. General settings include pump selection, definition of the experiment name, target flask volume, media density, target volume in the target flask, target cell density in the target flask, cell line, and passage number.

The process specific settings include the lot number of the medium, media volume in the target flask, cell density in the donor flask, passage date, lot number of the target flask, and number of flasks to be processed. All settings are stored in an experiment database. From the concentration and volume data, the application automatically calculates the target weight of the inoculum or media to be transferred during the transfer process and controls the peristaltic pump accordingly.

Experiment name	HeLa 12CUB
Receiving flask size	250.000 ml
Media density	0.99 g/ml
Target volume in receiving flask	100.000 ml
Target cell density (receiving flask)	1.00 10 <sup>6</sup> cls/ml
Cell line	HeLa
<span>✕</span> Parameters for experiment <span>✓</span>	

Target cell density (receiving flask)	1.00 10 <sup>6</sup> cls/ml
Actual cell density (prior flask)	10.00 10 <sup>6</sup> cls/ml
Passage date	2020-01-27
Cell line	HeLa
Passage number	2
Receiving flask lot	1234567
<span>✕</span> Experimental details 2 <span>&lt;</span> <span>&gt;</span>	

Start speed	25 rpm
Media filling speed	200 rpm
Media slow down at	90 %
Media to speed	10 rpm
Reverse run	10 rpm
Print label	On
<span>✕</span> Experimental details 3 <span>&lt;</span> <span>✓</span>	

In addition to the experimental settings, the speed for all steps during the pumping procedure is defined in rpm. The pumping procedure consists of an optional priming step to get rid of air bubbles remaining in the tube, transferring medium or inoculum to a defined percentage of the target weight, running the pump at slow speed until the final target weight is reached, and an optional reverse run to remove liquid from the tube above the weighing pan that would falsify the acquired weighing result.

The software displays the recommended default values, depending upon the type of pump used, for all steps of the pumping procedure but these can be modified by the user as needed. Additionally, depending on the selected method (add media, add inoculum or add media and inoculum) the software displays instructive text for setup of the flasks and tubes the pumping process. The pumping procedure is repeated according to the number of flasks to be filled until all samples are processed.

The screenshot shows the MYCAP® CCX cell passaging software interface. At the top, it displays "MYCAP® CCX cell passaging" and "Administrator 2020-01-27 15:36". Below this, there are fields for "isoCAL" and "Max 220 g d = 0.0001 g". The main display area shows "Gross" weight as "+ 49.9983 g" and "Target weight" as "99.000000 g". A message says "Pump is running. Wait until pump stops automatically." and there is an icon of a flask on a scale. At the bottom, there are navigation buttons: "0", "T", and "✕".

During sample processing the receiving flask is fixed on the balance weighing pan, the balance is tared and then the peristaltic pump transfers liquid from the donor flasks to the receiving flask. The volume to be transferred is automatically calculated by the software application and transformed into a target weight value using the media density.

The transferred liquid weight is measured by the balance and the pump speed is reduced at a set percentage of the target weight. The speed reduction prevents overshooting the target weight. The final weight is gravimetrically measured and the final results are calculated based on the measured weight value.

The final volume and the effective final cell density for each sample are calculated and can be documented by means of a printer connected to the balance. Labels with cell line information, lot number, passage number and date, effectively cell density, and the final media volume can optionally be printed for the receiving flasks.

Preview 1 of 1  
Printer YDP30 (only weight values) Administrator 2020-01-27 15:37

**Initial**  
 DatPas 2020-01-27  
 CellLine HeLa

**Passage** 2  
 RecLot ABC  
 Size 250.000 ml  
 TarVol 100.000 ml  
 MedVol 0.000 ml  
 Media 0.99 g/ml  
 TarMed 99.00000000 g

**Components**  
 n 1  
 Actual media weight

Result	
Media volume in receiving flask	50.000 ml
Target media weight	40.0000000 g
Target inoculum weight	10.0000000 g
Target total weight	50.0000000 g
Sample number	1
Actual media weight	49.9978 g
Actual inoculum weight	16.1915 g
Actual total weight	66.1893 g
Effective cell density	1.39 10 <sup>6</sup> cls/ml
Total volume	116.189 ml

# Sales and Service Contacts

For further contacts, visit  
[www.sartorius.com](http://www.sartorius.com)

## **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