

November 12, 2020

**Keywords or phrases:**

Celsius® S<sup>3</sup> Benchtop System and Celsius® FFT|FFTp

# Modelling Celsius® FFT|FFTp Freeze & Thaw Temperature Profiles in the Celsius® S<sup>3</sup> Benchtop System

Sartorius Stedim FMT S.A.S, ZI des paluds, Avenue de Jouques, 13781 Aubagne, France

Correspondence

E-Mail: [leads@sartorius.com](mailto:leads@sartorius.com)

## Abstract

Celsius® S<sup>3</sup> Benchtop System is a laboratory instrument that is designed to evaluate freeze & thaw processes of new drug candidates in 30 ml and 100 ml single-use bags to match the performances of the 100L-scale Celsius® CFT (Controlled-rate Freeze & Thaw) systems. This tool allows generation of a consistent samples library that is useful for evaluating stability, storage and shipping process steps and formulations to reproduce the exact same freeze & thaw conditions obtained at production scale.

In this study, the use of the Celsius® S<sup>3</sup> Benchtop System to match typical freezing and thawing temperature profiles of two other production-scale Celsius® platforms offered by Sartorius Stedim Biotech, Celsius® FFT and Celsius® FFTp, has been evaluated. Celsius® S<sup>3</sup> Benchtop System new freeze and thaw recipes have been developed to approximate typical performances that can be obtained using Celsius® FFT | FFTp in manufacturing. This offers to the end-users an efficient tool to investigate the behavior of the bulk drug substance while submitted to different freeze and thaw conditions and to support decision for the most suitable production-scale freeze & thaw platform.

**Find out more:** [www.sartorius.com](http://www.sartorius.com)

## Introduction

The Celsius® S<sup>3</sup> Benchtop System is a tool to execute freeze & thaw process development and stability studies using minimal amount of product. Specifically designed to be scalable to the production-scale Celsius® CFT System, freeze & thaw runs are performed with as little as 30 mL of product matching the performances obtained at 100L-scale. The scalability concept of the Celsius® CFT platform comes from the use of the same freezing path length, the same material of construction of the single-use bags and the same heat transfer technology at both lab and production-scale. Freezing and thawing rates are controlled, freezing and thawing times are equivalent and the critical freeze & thaw process parameters have been characterized to be maintained throughout process scales.

Sartorius Stedim Biotech offers to the industry two other production-scale Celsius® platforms, Celsius® FFT and Celsius® FFTp. Celsius® FFT (Flexible Freeze & Thaw) and Celsius® FFTp (Flexible Freeze & Thaw for Plate Freezer) is a unique bag-in-shell system that replaces traditional freezing and thawing methods. The single-use containers are sterile, pre-assembled and ready-to-use for freezing and thawing biopharmaceutical solutions in commercially available equipment with the intent to leverage existing freeze & thaw infrastructure. Celsius® FFT can be used with any type of conventional freezers while Celsius® FFTp is specifically designed to be frozen in horizontal type plate freezers. Material of construction of the single-use containers is identical to the one used in the Celsius® CFT platform.

In horizontal plate freezers, Celsius® FFTp takes benefits of controlled-rate freeze & thaw process with fast and reproducible thermal cycling. Performances may depend upon the equipment and related manufacturer as well as to the operating conditions. No scale-down model is today proposed by the industry to match the results obtained in horizontal plate freezers.

Performances of Celsius® FFT on the other hand can be more diversified and more challenging to predict as the freezing and thawing methods are “uncontrolled” and coming from a wide range of possibilities (upright freezer, walk-in freezer, blast freezer, thawing in incubator, thawing at room temperature, thawing with or without shaking). Process conditions (freezing and thawing set point, product load) will also influence the overall production scale freezing and thawing times. Uncontrolled freezing and thawing methods using existing infrastructure are the easiest, quickest and least costly. However, freeze and thaw rates in Celsius® FFT can become slow depending on the selected methods, especially when large volume is required to be processed.

This may cause degradation of the active ingredient due to cryo-concentration. Certain protein formulations are robust enough to be processed at slow rates but consideration for lab-scale studies from a process perspective requires to obtain similar time-temperature profiles throughout all process scales to confirm the impact of the freeze and thaw rates on drug substances. In this case, scalable solutions on the basis of heat and mass transfer dimensions become critical and lab-scale controlled-rate freeze and thaw processes are required to approach Celsius® FFT performances.

Even though Celsius® S<sup>3</sup> Benchtop System has not been designed to reproduce the results obtained at production-scale with Celsius® FFT|FFTp, special freezing and thawing temperature profiles can be developed to approximate those obtained with Celsius® FFT|FFTp. This would not be a truly scalable freeze & thaw process (freezing path length and heat transfer system different) but this will provide an approximation of the freeze & thaw conditions which will be experienced by the drug substances at production-scale in Celsius® FFT|FFTp.

## Material and Methods

### 1. Selection of the Celsius® FFT|FFTp Temperature profile

Celsius® FFT and Celsius® FFTp production-scale performances have been selected from a wide range of in-house and end-user data to bracket the variety of freezing and thawing times that can be obtained from the different freeze and thaw methods used in the industry.

#### a. Freezing Temperature profile

4 different process conditions for freezing have been selected and described in Table 1 to be reproduced with the Celsius® S<sup>3</sup> Benchtop System.

The following 3 process parameters from typical freezing curve (Figure 1) have been taken into account in order to develop the new temperature profiles in the Celsius® S<sup>3</sup> Benchtop System:

- time to reach the freezing point (0°C)
- time to remove the latent heat or nominal freeze time (NFT): time required for the temperature to go from +3°C to -5°C
- time to complete freeze: time required for the temperature to get stable at the vicinity of the final freezer setpoint

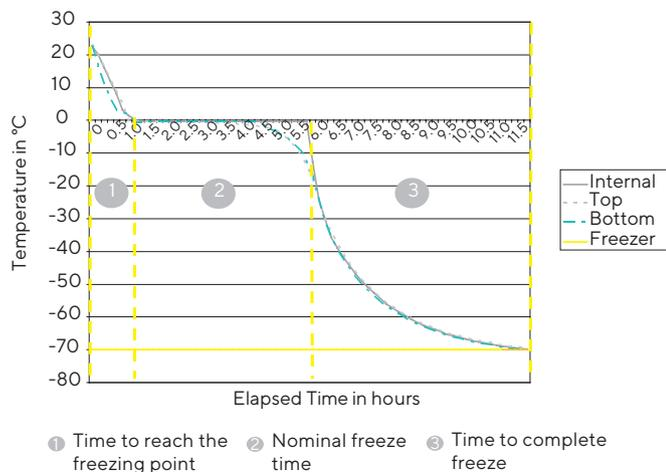


Figure 1: Celsius® FFT 2L typical freezing temperature profile

Celsius® FFT FFTp Size	Freezer Type   setpoint	Loading pattern	Freezing parameters duration	Data Source
12L Celsius® FFT	Upright freezer   -70°C	5 × containers (1 × container per shelf)	time to freezing point: ~6.5 hr nominal freeze time: ~17 hr complete freeze time: ~50 hr	end user
12L Celsius® FFT	Cold room   -25°C	20 × containers stacked on shelf (4 × 5-high containers)	time to freezing point: ~6 hr nominal freeze time: ~28 hr complete freeze time: ~60 hr	end user
12L Celsius® FFT	Blast freezer   -80°C	9-high containers stacked	time to freezing point: ~3 hr nominal freeze time: ~3 hr complete freeze time: ~11 hr	in-house SSB data
12L Celsius® FFTp	Horizontal plate freezer   -70°C	10 × containers (2 per shelf)	time to freezing point: ~1.5 hr nominal freeze time: ~2 hr complete freeze time: ~7 hr	end user

Table 1: Celsius® FFT|FFTp freezing conditions at production-scale selected for scale-down operation

### b. Thawing Temperature Profile

5 different process conditions for thawing have been selected and described in Table 2 to be reproduced with the Celsius® S<sup>3</sup> Benchtop System.

The following 3 process parameters from typical thawing curve have been taken into account in order to develop the new temperature profiles (Figure 2) in the Celsius® S<sup>3</sup> Benchtop System:

- time to reach the melting point (0°C)
- time at the melting point (0°C): time required for the temperature to go from -5°C to +3°C
- time to complete thaw: time required for the temperature to get stable at the vicinity of the final thawing setpoint with absence of ice



Figure 2: Celsius® FFT 2L typical thawing temperature profile

Celsius® FFT FFTp Size	Thaw method   setpoint	Loading pattern	Thawing parameters duration	Data Source	
12L Celsius® FFT	Flat table   RT	1 × container	time to melting point:	~ 6 hr	in-house SSB data
			time at melting point:	~ 28 hr	
			complete thaw time:	~ 40 hr	
12L Celsius® FFT	Flat table with forced air blowing   RT	1 × container	time to melting point:	~ 2 hr	in-house SSB data
			time at melting point:	~ 8 hr	
			complete thaw time:	~ 11 hr	
12L Celsius® FFT	Flat table with forced air blowing + 60 rpm agitation   RT	1 × container	time to melting point:	~ 1 hr	in-house SSB data
			time at melting point:	~ 3 hr	
			complete thaw time:	~ 5 hr	
12L Celsius® FFT	Blast Freezer   +22°C	9-high containers stacked	time to melting point:	~ 8 hr	in-house SSB data
			time at melting point:	~ 13 hr	
			complete thaw time:	~ 30 hr	
12L Celsius® FFTp	Horizontal plate freezer   +25°C	10 × containers (2 per shelf)	time to melting point:	~ 1.5 hr	end-user
			time at melting point:	~ 3 hr	
			complete thaw time:	~ 10 hr	

Table 2: Celsius® FFT|FFTp thawing conditions at production-scale selected for scale-down operation

## 2. Development of the new temperature profiles in Celsius® S<sup>3</sup> Benchtop System

### a. Loading samples for a freeze & thaw step

10x100ml Celsius®-Paks were filled to nominal fill volume via a syringe. The sample solution was deionized water for all experiments. Celsius®-Paks were placed in the heat exchange slots of the module of the Celsius® S<sup>3</sup> Benchtop System. Thermo couples were placed into the thermowells of the bags to be monitored and secured.

The thermocouple tip-end was located at the last point to freeze meaning  $1 \pm 0.5$  cm below the liquid surface. The module was closed.

### b. Creating new temperature profile

New temperature profiles need to be developed so that the product temperature curves obtained by using the Celsius® S<sup>3</sup> Benchtop System match the ones obtained by using the previously mentioned Celsius® FFT|FFTp production-scale performances described in table 1 and 2. Based on freezing and thawing curves from large-scale Celsius® FFT|FFTp product temperature curves, series of steps of X°C were drawn and translated into a set of numerical instructions in the text format readable by the CryoPilot software. CryoPilot software controls the chiller and the mixer by following a profile. A profile is a recipe with instructions to change the chiller temperature setpoint and the mixer mode setpoint (on| off) as a function of time.

Once the freeze & thaw runs were completed, the data files were placed into a spreadsheet application along with data from the large scale conditions. The data from both sources were plotted against each other and the parts of the set point profile that need modifications were determined in a graphical fashion. This procedure was repeated until both profiles from small and large-scale match each other to the desired degree.

### c. Running a freeze & thaw cycle

After having created and selected the desired profile, the CryoPilot software was started and used as is.

# Results and Discussion

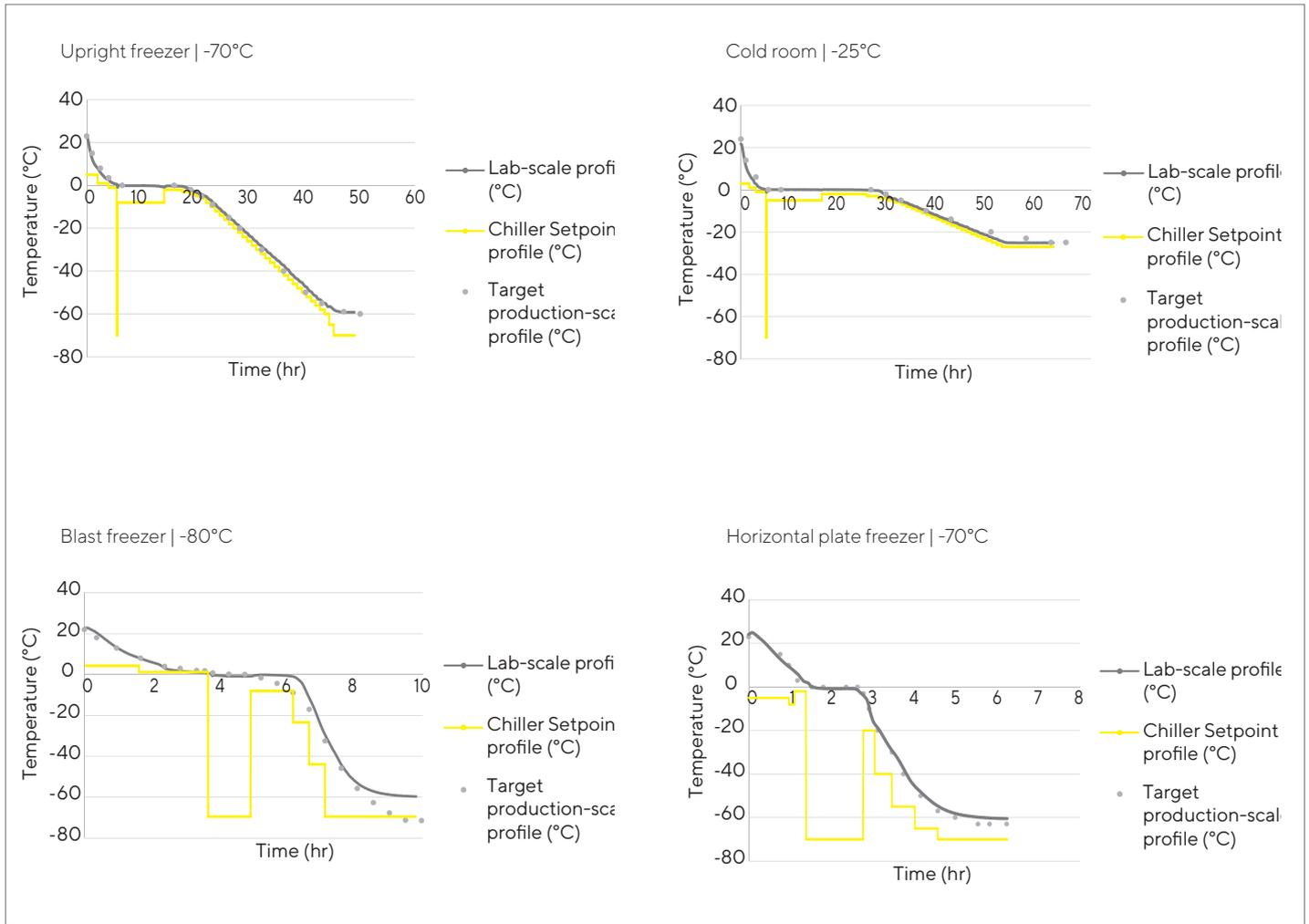


Figure 3: Superposition of production-scale and corresponding lab-scale freezing profiles generated with Celsius® S<sup>3</sup> Benchtop System. The lab-scale profile represents the average of the temperature measurement made on the 10 × 100 ml Celsius®-Pak by the 10 thermocouples. The target production-scale profile has been obtained from the average (when applicable) of the temperature measurement made from temperature sensors located inside and/or at the surface of the Celsius® FFT and FFTp containers.

Profiles		Parameters	Celsius® FFT   FFTp freezing profile	Celsius® S <sup>3</sup> Benchtop System freezing profile
Freezing	Upright freezer   -70°C	Time to freezing point	~6.5 hr	~ 5.5 hr
		Nominal freeze time	~17 hr	~ 18 hr
		Complete freeze time	~50 hr	~ 47 hr
	Cold room   -25°C	Time to freezing point	~ 6 hr	~ 5.5 hr
		Nominal freeze time	~28 hr	~ 29 hr
		Complete freeze time	~ 60 hr	~ 53 hr
	Blast freezer   -80°C	Time to freezing point	~ 3 hr	~ 4 hr
		Nominal freeze time	~ 3 hr	~ 4 hr
		Complete freeze time	~ 11 hr	~ 10 hr
Horizontal plate freezer   -70°C	Time to freezing point	~1.5 hr	~ 2 hr	
	Nominal freeze time	~ 2 hr	~ 2 hr	
	Complete freeze time	~ 7 hr	~ 6.5 hr	

Table 3a: Comparison of expected results and results obtained with scaled-down freezing models

Profiles		Parameters	Celsius® FFT   FFTp thawing profile	Celsius® S <sup>3</sup> Benchtop System thawing profile
Thawing	Flat table   RT	Time to melting point	~ 6 hr	~ 5.5 hr
		Time at melting point	~28 hr	~ 34 hr
		Complete thawing time	~40 hr	~ 42 hr
	Flat table with forced air blowing   RT	Time to melting point	~ 2 hr	~ 2 hr
		Time at melting point	~ 8 hr	~ 9 hr
		Complete thawing time	~ 11 hr	~ 12 hr
	Flat table with forced air blowing + 60 rpm agitation   RT	Time to melting point	~ 1 hr	~ 1 hr
		Time at melting point	~ 3 hr	~ 4 hr
		Complete thawing time	~ 5 hr	~ 6 hr
	Blast Freezer   +22°C	Time to melting point	~ 8 hr	~ 7 hr
		Time at melting point	~ 13 hr	~ 13 hr
		Complete thawing time	~ 30 hr	~ 25 hr
	Horizontal plate freezer   +25°C	Time to melting point	~ 1.5 hr	~ 1.5 hr
		Time at melting point	~ 3 hr	~ 4 hr
		Complete thawing time	~ 10 hr	~ 8 hr

Table 3b: Comparison of expected results and results obtained with scaled-down thawing models

N.B.Celsius® S<sup>3</sup> capacity limited to -60°C, even if set-point can be programmed at lower temperature.

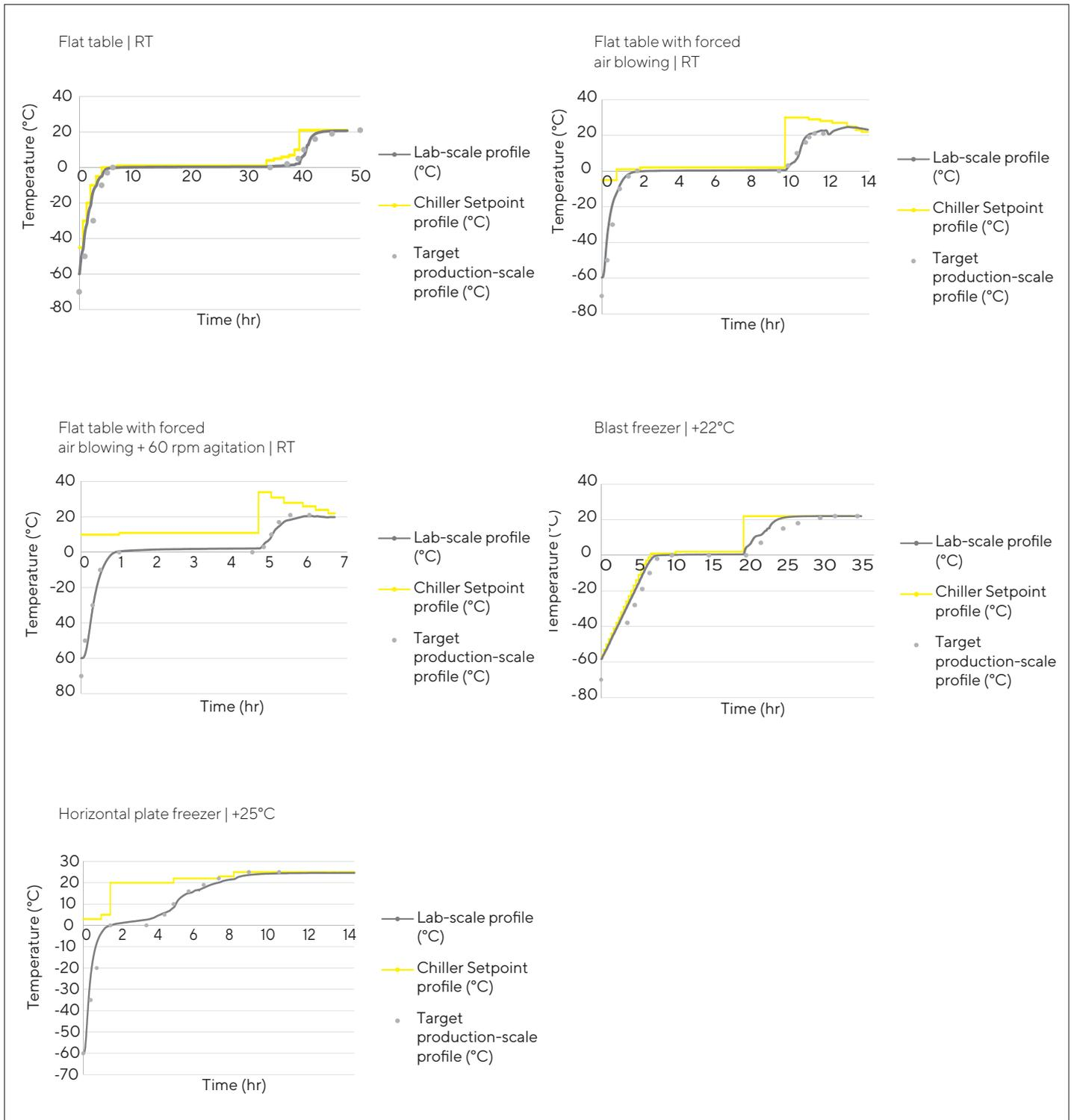


Figure 4: Superposition of production-scale and corresponding lab-scale thawing profiles generated with Celsius® S<sup>3</sup> Benchtop System. The lab-scale profile represents the average of the temperature measurement made on the 10 × 100 ml Celsius®-Pak by the 10 thermocouples. The target production-scale profile has been obtained from the average (when applicable) of the temperature measurement made from temperature sensors located inside and/or at the surface of the Celsius® FFT and FFTp containers.

Celsius® FFT|FFTp production-scale temperature profiles have been screened and selected to bracket typical conditions as well as freeze & thaw methods generally used in manufacturing environment. These selected Celsius® FFT|FFTp temperature profiles represent an example only of performances that could be expected when using the same or close process-scale conditions.

As outlined in Table 3, the newly created lab-scale temperature profiles are approaching the critical freezing and thawing process parameters of the selected Celsius® FFT|FFTp temperature profiles. Figure 1 and 2 show the temperature curves obtained with the lab-scale temperature profiles and their respective targeted production-scale Celsius® FFT|FFTp temperature profiles.

The Celsius® S<sup>3</sup> Benchtop System enables to control the chiller temperature between +45°C and -70°C with minimal 1°C possible temperature variation per step. This made possible to simulate long production-scale freezing or thawing processes that could eventually be obtained using Celsius® FFT|FFTp.

The lab-scale temperature profiles (Figure 3 and 4) obtained are close enough to their respective production-scale performances to satisfy the goal of this study.

Indeed, since the Celsius® FFT|FFTp production-scale performances may vary upon process conditions, equipment and methods, similar results rather than the exact same freezing and thawing time would be obtained by end-users when reproducing test parameters described in Table 1 and 2.

Although the scalability remains to be validated using proteins as a model, this work provides a mean to the end-user however to investigate the drug substance behavior at lab-scale as if it would be frozen and thawed at production-scale in Celsius® FFT|FFTp using different scenarios.

Coupled to the already available Celsius® CFT scalability feature, the Celsius® S<sup>3</sup> Benchtop System allows to the end-user to explore the impact of freeze & thaw on proteins through the entire portfolio of Celsius® Scalable Solutions to ultimately select the most appropriate one at production-scale to preserve the product quality.

## Conclusion

Celsius® S<sup>3</sup> Benchtop System was demonstrated to be suitable for the reproduction of temperature profiles that can be obtained when using Celsius® FFT|FFTp containers at production-scale.



Approaching Celsius® FFT|FFTp Temperature Profiles for Smart and Early Process Decision with Celsius® S<sup>3</sup> Benchtop System



Celsius® FFT



Celsius® S<sup>3</sup> Benchtop System



Celsius® FFTp

## Annex (Lab-scale freeze temperature profile)

### Profile Upright freezer | -70°C

	Time (min)	Chiller Set Point (°C)	Mixer State
START	0.00	5	OFF
	120.00	1	OFF
	240.00	-1	OFF
	330.00	-70	OFF
	340.00	-8	OFF
	850.00	-2	OFF
	1050.00	-3	OFF
	1140.00	-4	OFF
	1200.00	-5	OFF
	1260.00	-6	OFF
	1310.00	-8	OFF
	1360.00	-10	OFF
	1410.00	-12	OFF
	1460.00	-14	OFF
	1510.00	-16	OFF
	1560.00	-18	OFF
	1610.00	-20	OFF
	1660.00	-22	OFF
	1710.00	-24	OFF
	1760.00	-26	OFF
	1810.00	-28	OFF
	1860.00	-30	OFF
	1910.00	-32	OFF
	1960.00	-34	OFF
	2010.00	-36	OFF
	2060.00	-38	OFF
	2110.00	-40	OFF
	2160.00	-42	OFF
	2210.00	-44	OFF
	2260.00	-46	OFF
	2310.00	-48	OFF
	2360.00	-50	OFF
	2410.00	-52	OFF
	2460.00	-54	OFF
	2510.00	-56	OFF
	2560.00	-58	OFF
	2610.00	-60	OFF
	2660.00	-65	OFF
	2710.00	-70	OFF
			STOP

### Profile Cold room | -25°C

	Time (min)	Chiller Set Point (°C)	Mixer State
START	0.00	3	OFF
	100.00	1	OFF
	180.00	-1	OFF
	300.00	-70	OFF
	310.00	-5	OFF
	970.00	-2	OFF
	1510.00	-3	OFF
	1650.00	-4	OFF
	1770.00	-5	OFF
	1850.00	-6	OFF
	1930.00	-7	OFF
	1990.00	-8	OFF
	2050.00	-9	OFF
	2110.00	-10	OFF
	2170.00	-11	OFF
	2230.00	-12	OFF
	2290.00	-13	OFF
	2350.00	-14	OFF
	2410.00	-15	OFF
	2470.00	-16	OFF
	2530.00	-17	OFF
	2590.00	-18	OFF
	2650.00	-19	OFF
	2710.00	-20	OFF
	2770.00	-21	OFF
	2830.00	-22	OFF
	2890.00	-23	OFF
	2950.00	-24	OFF
	3010.00	-25	OFF
	3070.00	-26	OFF
	3130.00	-27	OFF
	3600.00	-27	OFF
			STOP

### Profile Blast freezer | -80°C

	Time (min)	Chiller Set Point (°C)	Mixer State
START	0.00	2	OFF
	100.00	-1	OFF
	230.00	-70	OFF
	320.00	-10	OFF
	370.00	-25	OFF
	410.00	-45	OFF
	450.00	-70	OFF
	600.00	-70	OFF STOP

### Horizontal plate freezer | -70°C

	Time (min)	Chiller Set Point (°C)	Mixer State
START	0.00	-5	OFF
	70.00	-8	OFF
	80.00	-2	OFF
	100.00	-70	OFF
	200.00	-20	OFF
	220.00	-40	OFF
	250.00	-55	OFF
	290.00	-65	OFF
	330.00	-70	OFF
	500.00	-70	OFF STOP

### Annex (Lab-scale thaw temperature profile)

#### Profile Flat Table | RT

	Time (min)	Chiller Set Point (°C)	Mixer State
START	0.00	-45	OFF
	40.00	-30	OFF
	80.00	-20	OFF
	120.00	-10	OFF
	180.00	-5	OFF
	240.00	0	OFF
	400.00	1	OFF
	2000.00	4	OFF
	2080.00	5	OFF
	2160.00	6	OFF
	2240.00	7	OFF
	2300.00	10	OFF
	2350.00	21	OFF STOP

#### Profile Blast freezer | +22°C

	Time (min)	Chiller Set Point (°C)	Mixer State
START	0.00	-56	OFF
	20.00	-53	OFF
	40.00	-50	OFF
	60.00	-47	OFF
	80.00	-44	OFF
	100.00	-41	OFF
	120.00	-38	OFF
	140.00	-35	OFF
	160.00	-32	OFF
	180.00	-29	OFF
	200.00	-26	OFF
	220.00	-23	OFF
	240.00	-20	OFF
	260.00	-17	OFF
	280.00	-14	OFF
	300.00	-11	OFF
	320.00	-8	OFF
	340.00	-5	OFF
	360.00	-3	OFF
	380.00	-1	OFF
	400.00	1	OFF
	600.00	2	OFF
	1150.00	22	OFF
	1500.00	22	OFF STOP

### Profile Flat Table with forced air blowing | RT

	Time (min)	Chiller Set Point (°C)	Mixer State	
START	0.00	-5	OFF	
	50.00	1	OFF	
	130.00	2	OFF	
	620.00	30	OFF	
	700.00	29	OFF	
	740.00	28	OFF	
	780.00	27	OFF	
	830.00	25	OFF	
	860.00	23	OFF	
	880.00	22	OFF	STOP

### Profile Horizontal plate freezer | +25°C

	Time (min)	Chiller Set Point (°C)	Mixer State	
START	0.00	3	OFF	
	60.00	5	OFF	
	90.00	20	OFF	
	300.00	22	OFF	
	450.00	23	OFF	
	500.00	25	OFF	
	600.00	25	OFF	STOP

### Profile Flat Table with forced air blowing +60 rpm agitation | RT

	Time (min)	Chiller Set Point (°C)	Mixer State	
START	0.00	10	ON	
	60.00	11	ON	
	280.00	34	ON	
	300.00	31	ON	
	320.00	28	ON	
	350.00	26	ON	
	370.00	24	ON	
	390.00	22	ON	
	400.00	22	OFF	STOP

**Germany**

Sartorius Stedim Biotech GmbH  
August-Spindler-Straße 11  
37079 Göttingen  
Phone +49 551 308 0

**USA**

Sartorius Stedim North America Inc.  
565 Johnson Avenue  
Bohemia, NY 11716  
Toll-Free +1 800 368 7178

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