

Filter Housing Heaters

Removable Heaters
for Biopharmaceutical
Applications



Introduction

The housing heaters offer tight temperature control and advanced functionality, while complying with strict safety regulations. The ELTC-14 control unit is a stand alone solution, which features a user friendly digital display and a high | low temperature alarm relay. It is CE tested and marked for both electrical and thermal safety.

Product Information

Sartorius Stedim housing heaters are a cost-effective and low maintenance solution for biopharmaceutical processes requiring superior thermal control. The housing heaters prevent vapor condensation in sterile vent and process filters and piping systems and help maintain preferred material viscosities.

Applications

Sartorius Stedim housing heaters are used in conjunction with sterile vent and process filter housings and are ideally suited for processes such as:

- Bioprocessing and Pharmaceutical Process Fluids
- Fermentation
- Product Recovery
- Water for Injection
- Filtration and Purification Processes
- Skidded Systems
- Tanks and Vessels
- Clean room versions on request



Standard Controller ELTC/H3-14



Controller ELT-GP1 4-20mA
With an additional analogue output for an external control center or SCADA

Features

- Digital temperature control
- Digital temperature display
- Adjustable temperature limit value (80°C adjusted)
- Adjustable hysteresis ($\pm 3^\circ\text{C}$ adjusted)
- Alarm relay
- Temperature monitoring by PT 100 – 2 wire
- Silicone insulation inside-Safe to touch
- Corrosion resistant
- CE approved and marked



Heating jacket front view



Heating jacket with strip fasteners at the back

Technical Data

Specifications of the controller

Supply Voltage	230 Volt $\pm 10\%$, 50 Hz
Switching Capacity	16 Ampere 230 Volt
Measuring Input	PT 100 DIN 2-Wire
Adjustment Range	0...+100°C
Ambient Range	-30...+60°C
Outside material	Glass fabric (PTFE coated)
CE Conformity	2004/108/EC 2006/95/EC 2011/65/EC
Lengths of Cables:	
Power supply cable (black)	3 m
Data cable (red)	2 m

Ordering Information

Removable heater incl.
ELTC/H3-14 controller

Order Code	Description
Mini-housing: 7ZE---0029	5" (size 9), height 120 mm, 40 W max.
1-round housing: 7ZE---0030	5", height 155 mm, 50 W max.
7ZE---0031	10", height 270 mm, 100 W max.
7ZE---0032	20", height 540 mm, 200 W max.
7ZE---0033	30", height 810 mm, 300 W max.
7ZE---0034	40", height 1,080 mm, 400 W max.
3-round housing: 7ZE---0035	10", height 270 mm, 150 W max.
7ZE---0036	20", height 540 mm, 200 W max.
7ZE---0037	30", height 810 mm, 300 W max.
7ZE---0038	40", height 1,080 mm, 400 W max.
5-round housing: 7ZE---0039	10", height 270 mm, 200 W max.
7ZE---0040	20", height 540 mm, 300 W max.
7ZE---0041	30", height 810 mm, 400 W max.
7ZE---0042	40", height 1,080 mm, 500 W max.

Removable heater incl.
ELT-GP1 4-20mA

Order Code	Description
Mini-housing: 7ZE---0049	5" (size 9), height 120 mm, 40 W max.
1-round housing: 7ZE---0050	5", height 155 mm, 50 W max.
7ZE---0051	10", height 270 mm, 100 W max.
7ZE---0052	20", height 540 mm, 200 W max.
7ZE---0053	30", height 810 mm, 300 W max.
7ZE---0054	40", height 1,080 mm, 400 W max.
3-round housing: 7ZE---0055	10", height 270 mm, 150 W max.
7ZE---0056	20", height 540 mm, 200 W max.
7ZE---0057	30", height 810 mm, 300 W max.
7ZE---0058	40", height 1,080 mm, 400 W max.
5-round housing: 7ZE---0059	10", height 270 mm, 200 W max.
7ZE---0060	20", height 540 mm, 300 W max.
7ZE---0061	30", height 810 mm, 400 W max.
7ZE---0062	40", height 1,080 mm, 500 W max.

Spare Parts

Just the jacket without controller

Order Code	Description
Mini-housing: 7ZE---0029M	5" (size 9), height 120 mm
1-round housing: 7ZE---0030M	5", height 155 mm
7ZE---0031M	10", height 270 mm
7ZE---0032M	20", height 540 mm
7ZE---0033M	30", height 810 mm
7ZE---0034M	40", height 1,080 mm
3-round housing: 7ZE---0035M	10", height 270 mm
7ZE---0036M	20", height 540 mm
7ZE---0037M	30", height 810 mm
7ZE---0038M	40", height 1,080 mm
5-round housing: 7ZE---0039M	10", height 270 mm
7ZE---0040M	20", height 540 mm
7ZE---0041M	30", height 810 mm
7ZE---0042M	40", height 1,080 mm

Just the controller without jacket

Order Code	Description
7ZE---0022	ELTC/H3-14
7ZE---0064	ELT-GP1 4-20 mA

Germany

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen
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
[sartorius.com](https://www.sartorius.com)