

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

Features

- Digital temperature control
- Digital temperature display
- Adjustable temperature limit value (80°C adjusted)
- Adjustable hysteresis ($\pm 3^{\circ}\text{C}$ adjusted)
- Alarm relay
- Water proof classification IP65 (controller), IP54 (heating jacket)
- Temperature monitoring by PT100 – 2 wire
- Silicone insulation inside-Safe to touch
- Corrosion resistant
- CE approved and marked

Technical Data

Specifications

Supply Voltage	230 Volt +/- 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
Protection Type Controller	IP65 Protection Class (I)
Protection Type heating jacket	IP54 Protection Class (II)
Outside material	Glass fabric (PTFE coated)
CE Conformity	2004/108/EC 2006/95/EC 2011/65/EC
Lengths of Cables:	
Power supply cable (black)	3.0 m
Data cable (red)	2.0 m



|1



|2

Control Unit ELTC-14 (1) | Strip fastener on the back side (2)

Ordering Information


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 1080 mm, 400 W max.
3-round housing:	
7ZE---0035	10", Height 270mm, 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 1080 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 1080 mm, 500 W max.

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
www.sartorius.com