



CERTOMAT[®] H | HK
Converts your shaker
to an incubator

CERTOMAT® H | HK

Converts your shaker to an incubator

The incubation hoods CERTOMAT® H and HK of Sartorius have proven their value in laboratories worldwide. They allow control of sample temperature in an airstream with a minimum of lab space. In combination with our range of CERTOMAT® benchtop shakers, bacteria and yeasts can be cultured under optimum conditions.

When combined with an external cooler or a coolant line, the CERTOMAT® HK will even enable culturing at temperatures near or below ambient. Both hoods feature a built-in additional memory function for re-establishing the temperature set point in the event of power failure.

Hood volumes are sufficient for the mounting of a wide range of commonly used incubation flasks.

The hide door of the plexiglass hood allows easy access to shaking vessels.

Benefits

Accomodate all Sartorius benchtop shakers with EU trays

Memory function

Temperature range:
+8 °C above ambient to +60 °C
(without cooling)

+10 °C to +60 °C (with cooling)



Technical Data

Ordering Information

Incubation Hood with heater: 115/230 V/50–60 Hz	886 3202	CERTOMAT® H
Incubation Hood with heater and heat exchanger: 115/230 V/50–60 Hz	886 3245	CERTOMAT® HK

Technical data

Temperature control	Incubation temperature:	CERTOMAT® H: +8 °C above ambient to +60 °C CERTOMAT® HK: +10 °C to +60 °C at +20 °C ambient temperature, external cooling required (not included with CERTOMAT® HK)
	Air circulation	180 m ³ /h
	Heater:	550 W
	Setting of temperatures:	keys + and –
Safety	Protection class:	1
Electrical	Max. current:	2.5 A /230 V, 5.0 A /115 V
Operation conditions	Temperature:	from +10 °C to +60 °C
	Humidity:	from 10% to 90%, not condensing
Dimensions and weight	External dimensions (W × H × D):	670 × 425 × 615 mm CERTOMAT® H 670 × 515 × 615 mm CERTOMAT® HK
	Incubation chamber (W × H × D):	520 × 420 × 600 mm
	Weight:	20 kg

All units are delivered without further accessories.

Accessories

For cultivating cells or for mixing liquids a benchtop shaker incl. tray is required. Maximum tray size is 420 × 420 mm, e.g. EU-tray 885 3002. In addition further accessories are needed to mount shaking flasks, separation funnels or tubes on the tray.

Sartorius AG
Weender Landstrasse 94-108
37075 Goettingen, Germany

Phone +49.551.308.0
Fax +49.551.308.3289

www.sartorius.com

Sartorius BBI Systems GmbH
Schwarzenberger Weg 73-79
34212 Melsungen, Germany

Phone +49.5661.71.3400
Fax +49.5661.71.3702

www.sartorius-bbi-systems.com

Sartorius North America Inc.
131 Heartland Blvd.
Edgewood, New York 11717, USA

Phone +1.631.254.4249
Toll-free +1.800.3687178
Fax +1.631.254.4253

Sartorius BBI Systems, Inc.
2800 Baglyos Circle
Bethlehem, PA 18020, USA

Phone +1.610.866.4800
Fax +1.610.866.4890

Sartorius Ltd.
Longmead Business Park
Blenheim Road, Epsom
Surrey, KT19 9 QQ, U.K.

Phone +44.1372.737159
Fax +44.1372.726171

Sartorius S.A.
4, rue Emile Baudot
91127 Palaiseau Cedex, France

Phone +33.1.6919.2100
Fax +33.1.6920.0922

Sartorius S.p.A.
Via dell'Antella, 76/A
50011 Antella (FI), Italy

Phone +39.055.63.40.41
Fax +39.055.63.40.526

Sartorius, S.A.
C/Isabel Colbrand 10-12
Planta 4, Oficina 121
Polígono Industrial de Fuencarral
28050 Madrid, Spain

Phone +34.91.3586091
Fax +34.91.3588804

Sartorius Technologies N.V.
Luchthavenlaan 1-3
1800 Vilvoorde, Belgium

Phone +32.2.756.0670
Fax +32.2.756.0681

Sartorius K.K.
KY Building, 8-17
Kitashinagawa 1-chome
Shinagawa-ku
Tokyo 140-0001, Japan

Phone +81.3.3740.5407
Fax +81.3.3740.5406

Specifications subject to change
without notice.

Printed in Germany on paper that has been
bleached without any use of chlorine.

W/sart-000 · G

Publication No.: SB-1024-e05022

Order No.: 85030-517-51