



Potter S Homogenizer
For gentle
cell disruption

Potter S Homogenizer For gentle cell disruption

The Potter S homogenizer of Sartorius is a well-established tool in laboratories worldwide, where gentle disruption of soft tissues or cultured cells is required. Due to its well-controlled action, this homogenizer can even be used for isolation of nuclei from such materials.

In cell biology, biochemistry and plant physiology, the Potter S serves to rapidly disrupt cells for the isolation of DNA, RNA and proteins, or to prepare cell organelles. In food analysis, the Potter S is used to homogenize tissues for subsequent analysis of various drugs, for instance in the retina of slaughtered animals. For testing of surface disinfectants, a suspension of single *Mycobacterium terrae* cells is generated from cell aggregates.

Homogenization with the Potter S is fast and reproducible. An integrated cooling vessel, filled with ice water or connected to an external cooling system, helps to control temperature of the sample during homogenization in order to slow down degradation of temperature-sensitive components.

If you have other homogenization tasks in your laboratory, the Micro-Dismembrators U and S, LABSONIC® M and P ultrasonic disintegrators and Dounce hand homogenizers may be an option. Please refer to our Laboratory Product Catalogue or to our website, www.sartorius.com

Benefits

Gentle action

Homogenization of cells from tissue cultures

Disruption of soft tissues

Isolation of cell nuclei

Disaggregation of bacteria

Sample can be cooled during homogenization

Controlled speed

Clamping device for cylinders or vessels protects the user

Large selection of vessels or cylinders

Plungers made of glass or PTFE

Technical Data

Ordering information

Potter S, including cooling vessel and one set of clamping rings	
853 3024	230 V/50–60 Hz
853 3032	115 V/50–60 Hz

Accessories:

For homogenization complete vessels made of borosilicate glass or homogenizer cylinders and the appropriate PTFE plungers are needed. Complete vessels consist of a glass cylinder and a glass plunger ground to match the cylinder. For this reason, these components are labelled with an individual number so that they can be properly matched by the user. All cylinders have a volume scale. The choice between glass vessels and glass cylinders with PTFE plungers depends on the intended application.

Cylinders and vessels from 2 to 15 ml have a gap of approx. 0.045–0.065 mm, the larger ones have a larger gap of approx. 0.095–0.115 mm.

Technical data

Speed:	150–1500 min ⁻¹
Accuracy:	max. –3%
Maximum orbit:	170 mm
Drive:	DC motor
Power consumption:	70 W
Adjustment of speed:	stepless, potentiometer
Admissible temperature range:	+10 °C to +60 °C
Admissible humidity:	10 to 90%, non-condensing
Power supply:	230 V/50–60 Hz 115 V/50–60 Hz
Dimensions (W×H×D) in mm:	300×850×300
Weight (net):	approx. 12.5 kg

All instruments are delivered excluding further accessories.



Accessories



Vessels made of borosilicate glass, with latticed notches, complete, cylinder with ground-in glass plunger

Cat.-No.	Description
854 0756	Homogenizer vessel 2 ml, with latticed notches
854 0705	Homogenizer vessel 5 ml, with latticed notches
854 0802	Homogenizer vessel 15 ml, with latticed notches
854 0900	Homogenizer vessel 30 ml, with latticed notches
854 1000	Homogenizer vessel 60 ml, with latticed notches



Vessels made of borosilicate glass, without latticed notches, complete, cylinder with ground-in glass plunger

Cat.-No.	Description
854 1957	Homogenizer vessel 2 ml, without latticed notches
854 1906	Homogenizer vessel 5 ml, without latticed notches
854 2007	Homogenizer vessel 15 ml, without latticed notches
854 2104	Homogenizer vessel 30 ml, without latticed notches
854 2201	Homogenizer vessel 60 ml, without latticed notches



Cylinders made of borosilicate glass, for plungers made of PTFE

Cat.-No.	Description
854 2252	Homogenizer cylinder 2 ml, for plungers made of PTFE
854 2309	Homogenizer cylinder 5 ml, for plungers made of PTFE
854 2406	Homogenizer cylinder 15 ml, for plungers made of PTFE
854 2503	Homogenizer cylinder 30 ml, for plungers made of PTFE
854 2600	Homogenizer cylinder 60 ml, for plungers made of PTFE

Plungers made of PTFE, including shaft made of stainless steel

Cat.-No.	Description
854 2651	Plunger made of PTFE 2 ml, for cylinder 854 2252
854 2708	Plunger made of PTFE 5 ml, for cylinder 854 2309
854 2805	Plunger made of PTFE 15 ml, for cylinder 854 2406
854 2902	Plunger made of PTFE 30 ml, for cylinder 854 2503
854 3003	Plunger made of PTFE 60 ml, for cylinder 854 2600

Special accessories for POTTER S

Cat.-No.	Description
853 2206	Storage rack for 10 homogenizer cylinders and plungers or complete vessels
853 3130	Clamping ring for vessels or cylinders 2 ml
853 3148	Clamping ring for vessels or cylinders 5 ml
853 3156	Clamping ring for vessels or cylinders 15 ml
853 3164	Clamping ring for vessels or cylinders 30 ml
853 3172	Clamping ring for vessels or cylinders 60 ml
853 3180	Exchange cooling vessel, glass

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-1022-e05022

Order No.: 85030-515-86