

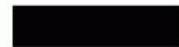


Information and Instructions on Disposal and Recycling

The packaging is made of environmentally friendly materials that can be used as secondary raw materials. If you no longer need this packaging, bring it to your local recycling and waste disposal facility according to the regulations applicable in your country. In Germany, you can dispose of this material using the VFW dual system (contract number D-59101-2009-1129).



The equipment, including accessories and empty non-rechargeable and rechargeable batteries, does not belong in your regular household waste; such equipment is manufactured from high-grade materials and can be recycled and reused. The European Directive 2002/96/EC on Waste Electrical and Electronic Equipment (WEEE) requires that electrical and electronic equipment be collected and disposed of separately from other unsorted municipal waste, with the aim of recycling it. The crossed-out waste bin symbol indicates that separate collection is required.



In Germany and many other countries, Sartorius Weighing Technology GmbH and its branch offices take care of the return and legally compliant disposal of their electrical and electronic equipment on their own. **Such equipment may not be thrown out with household waste or brought to collection centers run by local public disposal operations – not even by small commercial operators.** Please contact our on-site service staff or our central Services unit (Service Center) in Goettingen, Germany:

Sartorius Weighing Technology GmbH
Service Center
Weender Landstrasse 94-108
37075 Goettingen, Germany

Tel.: +49.551.308.3333
Fax: +49.551.308.3730
E-mail: technical-service-germany@sartorius.com

For equipment disposal in the following countries of the European Economic Area (EEA), please contact your local Sartorius affiliate, subsidiary, dealer or distributor:

Country	Person to Contact	Phone Fax	E-mail
Austria	Sartorius Mechatronics Austria GmbH	+ 43 1 796 57 60 + 43 1 796 57 60 24	info.austria@sartorius.com
Belgium	Sartorius Mechatronics Belgium N.V.	+32 2 756 06 80	info.belgium@sartorius.com
Cyprus	A.J. Vouros Ltd.	+357 22 442822	klitosth@logosnet.cy.net
Czech Republic	Sartorius Service and Consulting Center Karolina-Express	+420 475 200870	info@sartorius.cz
Denmark	VWR – Bie & Berntsen A/S	+45 43 86 87 88	info@dk.vwr.com
Estonia	Sartorius Service and Consulting Center OÜ Est-Doma	+37 27 362 716 +37 27 441 997	jaanus.vahtra@estdoma.ee
Finland	Sartorius Mechatronics UK Ltd	+44 1372 737 101	service@sartorius.com
France	Sartorius Mechatronics France SAS	+33 1 6919 2100	service.client@sartorius.com
Greece	Analytical Instruments A.E.	+30 210 67 48 973	a.nikolakis@analytical.gr

Information and Instructions on Disposal and Recycling

Country	Person to Contact	Phone Fax	E-mail
Hungary	Sartorius Mechatronics Hungária Kft.	+36 23 457 227 +36 23 457 147	mechatronika@sartorius.hu
Ireland	Sartorius Mechatronics UK Ltd	+353 1 8089050	info.ireland@sartorius.com
Italy	Sartorius Mechatronics Italy S.r.l.	+39 039 46591	info@sartorius.it
Latvia	DOMA, SIA	+371 737 62 89 +371 732 62 99	info@domagroup.lv
Lithuania	Sartorius Service & Consulting Center VILDOMA, UAB	+370 5 236 36 57 +370 5 236 36 55	dd@vildoma.lt
Luxembourg	Sartorius Mechatronics Belgium N.V.	+32 2 756 06 80	info.belgium@sartorius.com
Malta	Technoline Ltd.	+356 21 344 345	esaguna@technoline-mt.com
Netherlands	Sartorius Mechatronics Netherlands B.V.	+31 30 6052864	service.nl@sartorius.com
Norway	Sartorius Mechatronics UK Ltd	+44 1372 737 101	service@sartorius.com
Poland	Sartorius Mechatronics Poland Sp.z.o.o.	+48 61 647 3830 +48 61 647 3839	info.pl@sartorius.com
Portugal	Sartorius Mechatronics S.A.- Representação em Portugal	+351 217501000	manuel.muguiro@sartorius.com
Sweden	Sartorius Mechatronics UK Ltd	+44 1372 737 101	service@sartorius.com
Switzerland	Sartorius Mechatronics Switzerland AG	+41 44 746 50 25	roland.baumgartner@sartorius.com
Slovakia	Sartorius Service and Consulting Center Sartorius Slovensko s.r.o.	+421 31 552 6498	filtre@sartorius.sk
Slovenia	Sartorius Service and Consulting Center ALBA Trgovina in Proizvodnja, d.o.o.	+386 3 4281 800 +386 3 4281 810	alba@siol.net
Spain	Sartorius Mechatronics Spain S.A.U.	+34 91 358 60 95	spain.service@sartorius.com
UK	Sartorius Mechatronics UK Ltd	+44 1372 737 101	service@sartorius.com

In countries that are not members of the European Economic Area (EEA) or where no Sartorius affiliates, subsidiaries, dealers or distributors are located, please contact your local authorities or a commercial disposal operator.

Sartorius Weighing Technology GmbH, its affiliates, subsidiaries, dealers and distributors will not take back equipment contaminated with hazardous materials (ABC contamination); please refer to our General Terms and Conditions for Services. Before returning any equipment to us, the sender must fill out the attached Non-hazardous Declaration completely and return it to us in advance. No repair or return is possible unless a completed Non-hazardous Declaration form has been submitted.

Information and Instructions on Disposal and Recycling

Prior to disposal or scrapping of the equipment, remove any rechargeable or non-rechargeable batteries and, if they are drained, dispose of them in local collection boxes. Only for customers in Germany: On request, Sartorius will have the Foundation of Manufacturers' Mutual Return System for Batteries (Stiftung Gemeinsame Rücknahmesystem Batterien) provide customers in Germany with battery collection boxes.

Note: In Germany, Sartorius Weighing Technology GmbH is entered in the EAR Registry of the Foundation for Used Electrical and Electronic Equipment (Stiftung Elektro-Altgeräte Register). The WEEE registration number for Germany is DE 49923090.

Information and Instructions on Disposal and Recycling

✂ ----- ✂

Non-hazardous Declaration

Declaration of Safety and Acceptability as Non-Hazardous to Health

The safety and health of our employees, the legal regulations governing the handling of materials that pose hazards to health and the environment and the regulations governing occupational safety mandate and require that this Non-hazardous Declaration be filled out completely for all equipment returned to the respective Sartorius affiliate, subsidiary, dealer or distributor listed for your country. No repair or return is possible unless a completed Non-hazardous Declaration form has been submitted.

Fax (at +49 551 308 3737) or mail a copy of the completed Non-hazardous Declaration form in advance to the respective Sartorius affiliate, subsidiary, dealer or distributor listed for your country. We need to have received this before the equipment/component arrives. An additional copy must accompany the equipment/component. If necessary, the shipping company must be notified.

Incomplete entries or non-compliance with this procedure will automatically lead to considerable delays in processing.

In the event of return, we do not differentiate between

- still utilizable and/or repairable used electrical and electronic equipment and no longer utilizable and/or repairable used electrical and electronic equipment or
- between waste that is subject to monitoring by the supervisory authorities and waste that is not subject to any special monitoring by supervisory authorities (used electrical and electronic equipment).

Model:	Serial No.:
I/we guarantee that the following measures have been taken: <ul style="list-style-type: none"> <input type="checkbox"/> The equipment/component has been freed of ABC contamination to ensure that persons involved with the handling/repair are not exposed to any hazard or danger whatsoever. <input type="checkbox"/> The equipment has been packaged for safety and the complete labeling has been affixed to it. <input type="checkbox"/> The shipping agent has been notified (as prescribed) about the hazardous nature of the shipment. 	
The person sending in the equipment/component is hereby notified that said person shall be liable to Sartorius Weighing Technology GmbH, its affiliates, subsidiaries, dealers and distributors as well as to any third parties – in particular to any of the employees of Sartorius Weighing Technology GmbH, its affiliates, its subsidiaries, dealers or distributors involved with the handling/repair of the equipment/component – for any damage caused by entries that are incomplete or incorrect on account of negligence, gross negligence or willful intent.	
Name: Position: Date: Signature: Company stamp:	

✂ ----- ✂