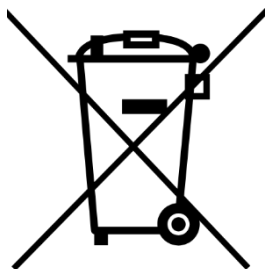


EU Directive on Waste Electrical and Electronic Equipment (WEEE) - Product Recycling Information



This product is subject to **Directive 2012/19/EU of the European Parliament and of the Council on Waste Electrical and Electronic Equipment (WEEE)**.

The purpose of the WEEE Directive is to minimise the environmental impact of electrical and electronic equipment at end-of-life by promoting separate collection and environmentally sound treatment, recycling and recovery of such equipment. Under the principle of extended producer responsibility (EPR), producers are required to ensure appropriate arrangements for the management of WEEE arising from products they place on the European market.

All products covered by WEEE Directive are marked with the **crossed-out wheeled bin** symbol, indicating that the product must not be disposed of as unsorted municipal waste and must be collected separately at end-of-life.

WEEE collection and recycling arrangements for Sartorius products

For Sartorius products marked with the crossed-out wheeled bin symbol and therefore covered by Directive 2012/19/EU, the following arrangements apply for end-of-life recycling.

Sartorius has established a recycling agreement with the **European Advanced Recycling Network (EARN)** to ensure appropriate treatment, recycling and recovery of WEEE for which Sartorius is responsible.

Under this agreement:

- Sartorius bears the costs of treatment, recycling and recovery of WEEE for which Sartorius is responsible.
- The business end-user is responsible for organising and covering the costs of transporting the WEEE to designated EARN recycling facility in the relevant country.

To arrange transport of WEEE for which Sartorius is responsible, please complete the online return form using the following link:

<https://b2btool.earn-service.com/sartorius-parcel>

The generated document will include the address of the designated EARN recycling facility, your address as the consignor, and other necessary information required for processing. Please print the document twice, attach on copy visibly to the outside of the package and place the second copy inside the package with the equipment.

Important note – Luxembourg

For customers located in Luxembourg, please do not use the online return form. Instead, contact EARN directly to organise WEEE collection and treatment. In Luxembourg, Sartoris will also cover the costs of transport from the business end-users premises to the EARN recycling facility.

[contact EARN directly](#)