

WEEE RoHS Compliance

English

This product is required to comply with the European Union's Waste Electrical & Electronic Equipment (WEEE) Directive 2002/96/EC. It is marked with the following symbol:



Thermo Fisher Scientific has contracted with one or more recycling or disposal companies in each European Union (EU) Member State, and these companies should dispose of or recycle this product. See www.thermoscientific.com/en/about-us/promotions/epm-weee-compliance.html?ca=WEEERoHS for further information on Thermo Fisher Scientific's compliance with these Directives and the recyclers in your country.