

Disposal of Used ECDs

- ECD as Radioactive Waste

Environmental and Safety Regulations and Requirements for Working with ECDs

The ECD (Electron Capture Detector) is a chromatographic detector that may be used in a GC (Gas Chromatograph) instrument. ECDs from Agilent Technologies contain Ni-63 as a radioactive electrodeposited sealed source with a maximum activity of 555MBq.

Users and holders of ECDs must adhere to all applicable relevant environmental, health, safety, and nuclear regulatory requirements. The information here is for guidance only and legal responsibility remains with ECD holders and users.

This document describes how to dispose of a used ECD (an ECD as radioactive waste) when the ECD is no longer needed and/or is at the end of its operating life (EOL). Return of an unused ECD is not governed by this process.

Warning: ECDs or instruments including them must never be disposed of in WEEE or other mixed or general waste streams.

Disposal of Used ECDs

Regulatory requirements vary from country to country; some countries require the disposal of a used ECD to be done within the country and some others require the disposal to be done outside of the country. To ensure regulatory compliance and promote environment protection, Agilent strongly encourages our customers to dispose of used ECD(s) via a licensed local service provider; however, under circumstances where ECD disposal must be done by Agilent, Agilent provides two ECD disposal services - Local Disposal and Return to US Factory. Please contact the local sales representative or customer service center to obtain a quote on the relevant ECD disposal service provided by Agilent.

In countries where the removal of the ECD(s) from a GC must be done by licensed personnel, customers should have a licensed service provider remove the ECD(s) from their GC or call Agilent for ECD removal service before ECD disposal arrangement.

Agilent strives to serve our customers and comply with local regulatory requirements; as a result, country-specific service coverage is offered - please refer to the table in Appendix for the list of ECD disposal service coverage by country. The coverage is accurate as of the date of publishing and is subject to change to meet US NRC and local regulatory requirements worldwide; please always check with your Radiation Safety Officer and local Agilent customer service center to get the latest information about ECD disposal.

Packaging, Labeling, and Marking of ECD Packages

Refer to [“Agilent Electron Capture Detector, Information for General Licensees” \(5961-5664E\)](#) for Procedures on Packaging, Labeling, and Marking of ECD Packages for Return to Agilent Technologies

Appendix

Region	Country	Local Disposal by Agilent	Agilent US Factory Takeback	Comments
EMEA	Austria	No	No	ECD disposal managed by customer
	Belgium	No	No	ECD disposal managed by customer
	Denmark	No	No	ECD disposal managed by customer
	Finland	Yes	No	Agilent provides in-country ECD disposal service through 3 rd party service provider
	France	No	Yes	Customer ships ECD to the Agilent US factory
	Germany	No	No	ECD disposal managed by customer
	IDO (EMEA)	No	No	ECD disposal managed by customer or distributor
	India	No	Yes	Agilent collects ECD and ship to US factory for disposal
	Ireland	Yes	No	Agilent provides in-country ECD disposal service through 3 rd party service provider
	Italy	Yes	No	Agilent provides in-country ECD disposal service through 3 rd party service provider
	Luxemburg	No	No	ECD disposal managed by customer
	Netherlands	No	No	ECD disposal managed by customer
	Portugal	No	Yes	Customer ships ECD to the Agilent US factory
	Spain	No	Yes	Customer ships ECD to the Agilent US factory
	Sweden	No	No	ECD disposal managed by customer
	Switzerland	Yes	No	Agilent provides in-country ECD disposal service through 3 rd party service provider
	UK	Yes	No	Agilent provides in-country ECD disposal service through 3 rd party service provider

Region	Country	Local Disposal by Agilent	Agilent US Factory Takeback	Comments
North / Central / South America	Canada	No	Yes	Customer ships ECD to the Agilent US factory
	US	No	Yes	Customer ships ECD to the Agilent US factory
	Mexico	No	Yes	Customer ships ECD to the Agilent US factory
	Brazil	No	Yes	Customer ships ECD to the Agilent US factory
	Others	No	No	ECD disposal managed by customer
Asia Pacific	China	No	Yes	Agilent collects ECD and ship to US factory for disposal
	Japan	No	Yes	Agilent collects ECD and ship to US factory for disposal
	Others	No	No	ECD disposal managed by customer
Rest of World		No	No	ECD disposal managed by customer

www.agilent.com/chem

[This information is subject to change without notice.]