



Waste Electrical and Electronic Equipment for Recycling from Molecular Devices Customers

Under the Dutch Waste Electrical and Electronic Equipment (WEEE) Management Regulations, when customers buy new electrical and electronic equipment from Molecular Devices they are entitled to send old equipment for recycling on a one-for-one, like-for-like basis. Customers are also entitled to send the new Molecular Devices equipment for recycling when it ultimately becomes waste. Molecular Devices has established a recycling contract with the European Advanced Recycling Network (EARN) and will cover any costs of recycling the equipment in this box on arrival at an EARN recycling company. Please complete this label and attach it to a suitable box for transport of old equipment to EARN.

Please confirm that the enclosed waste (EWC code 16 02 14) is:	
Old equipment that has been replaced by a purchase of new Molecular Devices equipment after 13 August 2005 (on a one-for-one, like-for-like basis); and / or	✓
Molecular Devices Equipment purchased after 13 August 2005 which has become waste	✓

Please indicate the quantity of waste in this container: _____ kg / items

Please indicate the carrier company that you will use:

	Carrier Name	Waste Carrier Registration Number	Expiry Date
✓	Koninklijke TNT Post B.V.	ZH509208VXXX	17/05/2016

Transfer from you to licensed waste carrier	Transfer from licensed waste carrier to recycler
Date:	Date:
Time:	Time:
Your Signature:	Waste Carrier Signature:
Waste Carrier Signature:	Recycler Signature:

To be sent to:

EARN Partner
Coolrec BV, Spaarpot 6
5667 KX Geldrop
P.O Box 49
NL - 5660 AA
Geldrop 916
NETHERLANDS

From (your company name and address):



Under Dutch Environmental Management Regulations you are required to keep a signed copy of this label for five years.