

**Intended Use of the Equipment**

This heater is intended only for use in combination with Agilent 45 l/s SEM I.B. Ion Pumps. Do not use these heaters for other applications or in combination with other pumps.

Technical Specifications and Equipment Ratings**919-2836 and 03.704491-01**

Input Voltage: 120 Vac, 50-60Hz

Input Power: 160 W

919-2837 and 03.704491-02

Input Voltage: 230 Vac, 50-60Hz

Input Power: 160 W

Certification and Norms:

CE mark

Norm EN-60204

cCSAus mark (available only for the 9192837 and 03.704491-02 230Vac version)

Norm CAN/CSA C22.2 No.61010-1 and UL 61010-1

installation category: 2

pollution degree: II

Installation

For the installation on the ion pump, refer to the “**87-900-101-01**” “**Ion SEM Pumps**” user manual at “**HEATER INSTALLATION**” chapter.

The heater has to be connected to mains by technical people.

Connection table of the main power cord:

wire	connection
blue	neutral
brown	phase
yellow/green	earth ground
metal braid	earth ground

The heater has to be installed by providing 10A protection fuses on both lines (phase and neutral).

For different protection levels, additional evaluation shall be conducted on the end-use equipment.

Instructions for Use

The heater's supply cord is intended for internal wiring use, in the final installation it shall be fully contained into the end-use equipment.

A disconnection mean shall be foreseen into the end-use equipment where the heater is installed.



WARNING!

The temperature reached by the heater and the ion pump during the heating (bake-out) are dangerous. Do not touch the heater/ion-pump during the bake out.
These heaters must be used with a maximum ambient temperature of 40°C.

Maintenance

No maintenance is needed for the heater.

DISPOSAL

Meaning of the "WEEE" logo found in labels

The following symbol is applied in accordance with the EC WEEE (Waste Electrical and Electronic Equipment) Directive.

This symbol (**valid only in countries of the European Community**) indicates that the product it applies to must NOT be disposed of together with ordinary domestic or industrial waste but must be sent to a differentiated waste collection system.

The end user is therefore invited to contact the supplier of the device, whether the Parent Company or a retailer, to initiate the collection and disposal process after checking the contractual terms and conditions of sale.

