

Instructions for use

Agilent Heater for Medium Vaclon Pumps

Intended Use of the Equipment

This heater is intended only for use in combination with Agilent 40-55-75 l/s VIPs.
Do not use these heaters for other applications or in combination with other pumps.

Product Name	Part Number
HEATER FOR VIP PLUS 40-55-75,120V CE/CSA	9192812
HEATER FOR VIP PLUS 40-55-75,230V CE/CSA	9192813

Technical Specifications and Equipment Ratings

9192812

Voltage	120 Vac, 50-60Hz
Power	180W

9192813

Voltage	230 Vac, 50-60Hz
Power	180W

Certification and Norms

CE mark

Norm EN-60204

cCSAus mark (available only for the 9192813 and 03.704492-02 230Vac version)

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

installation category: 2

pollution degree: II

Instructions for use

Installation

For the installation on the Ion pump, refer to the “87-900-105-01” “*Vac-Ion Plus 40/55/75 pumps*” user manual at “Heater installation” chapter. The heater must 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 must 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.

Maintenance

No maintenance is needed for the heater.

Instructions for use

Disposal

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.



For more information refer to:

<http://www.agilent.com/environment/product/index.shtml>

Customer Support & Service:

NORTH AMERICA:

Toll Free: 800 882 7426
vpl-ra@agilent.com
Lexington-service@agilent.com

EUROPE:

Toll Free: 00 800 234 234 00
vpt-customer-care@agilent.com

PACIFIC RIM:

please visit our website for individual office information
<http://www.agilent.com/chem/vacuum>

Worldwide Web Site, Catalog and Order Online:

www.agilent.com/chem/vacuum
Representatives in most countries

This information is subject to change without notice.

© Agilent Technologies, Inc. 2024

Printed in Italy, 11/2024

87-900-134-01 (B.00)