

## Agilent Fan installation

Product Name	P/N
Medium-TMP air cooling kit board	9699297

New fan mod. 9699297 is suitable for Agilent pumping systems TV 551 Navigator (models 9698832, 9698833, 9698834 and 9698835), TV 701 Navigator (models 9698836 and 9698837) and TV 1001 Navigator (models 9698838, 9698839, 9698840 and 9698841).

### 1.1 Fan outline

The following figure shows the outline of the new fan.

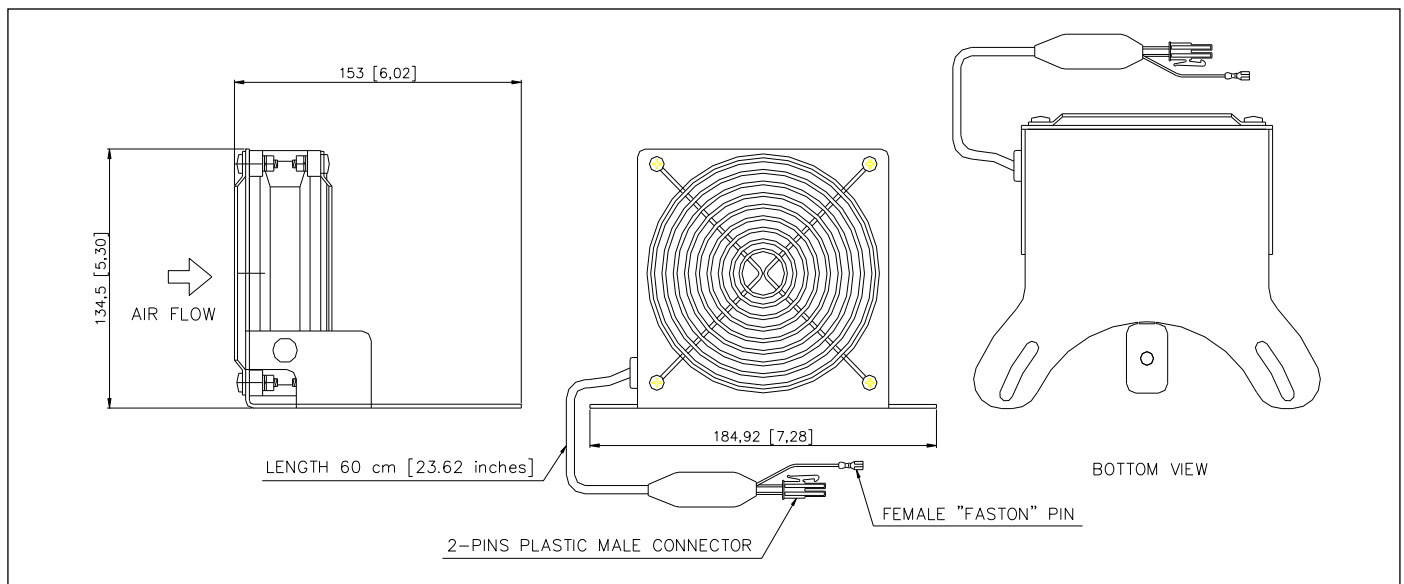


Figure 1.1 Outline drawing

## Instructions for use

### 1.2 Fan preparation

To prepare the fan for the installation it is necessary to detach the spacer from the fan installation flange (see the following figure).

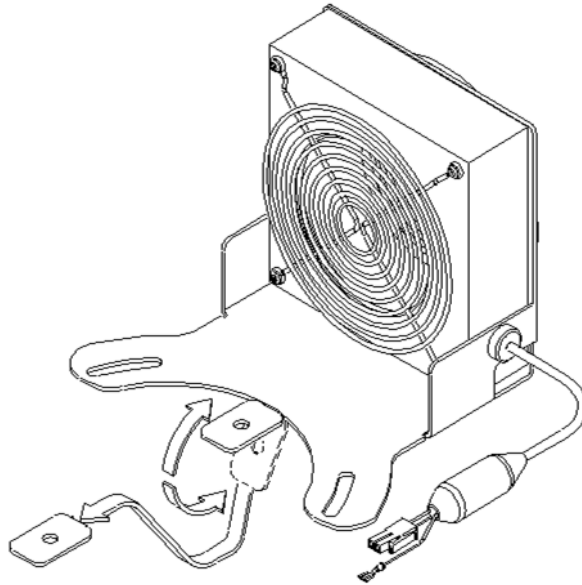


Figure 1.2 Spacer removal

#### NOTE

Do not lose the spacer: it must be used during the fan installation.

### 1.3 Controller removal

Before the fan installation, the Navigator controller must be detached from the pump as detailed in the following procedure (see figures 1.3.1 and 1.3.2):

1. Remove the power cable from the controller.
2. Disconnect the High Voltage connector from the pump.
3. Turn the pump upside-down.
4. Unscrew the three socket head screws M5 that fix the controller to the three fixing blocks.
5. Remove the controller.
6. Unscrew the three M6 screws that fix the fixing blocks to the pump.

## Instructions for use

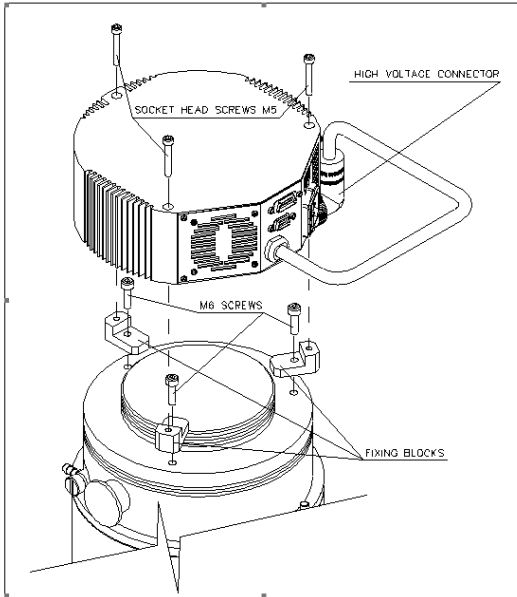


Fig. 1.3.1 TV 551/701 Navigator controller removal

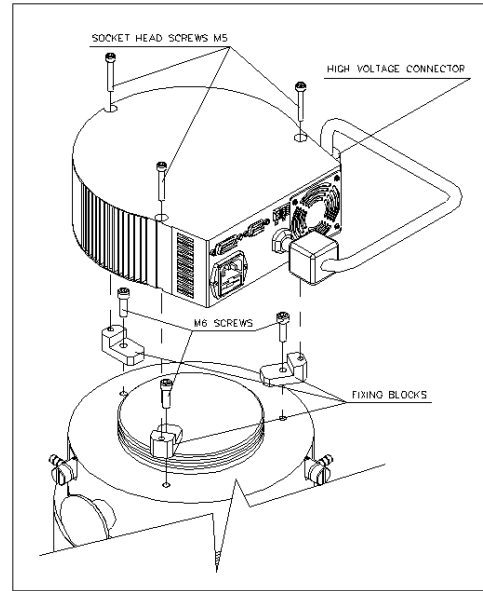


Fig. 1.3.2 TV 1001 Navigator controller removal

### 1.4 Fan installation

To install the fan the following procedure must be executed (see figures 1.4.1 and 1.4.2):

1. Position the fan and the spacer on the pump bottom.
2. Position the three fixing blocks.
3. Fix the three fixing blocks, the fan and the spacer with the three M6 screws removed during the controller removal.

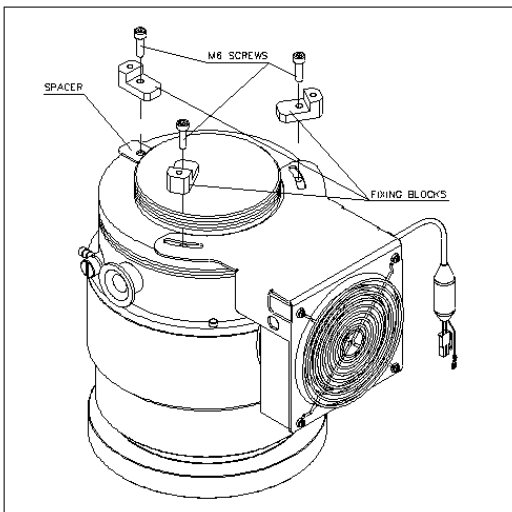


Fig. 1.4.1 TV 551/701 Navigator fan installation

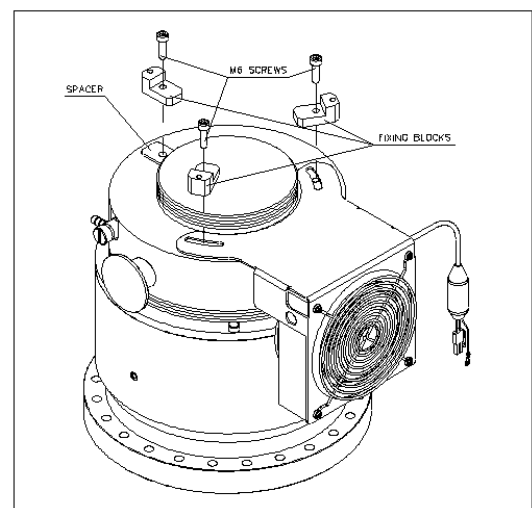


Fig. 1.4.2 TV 1001 Navigator fan installation

## Instructions for use

### 1.5 Controller installation

To complete the fan installation, the Navigator controller must be reinstalled as detailed in the following procedure (see figures 1.5.1 and 1.5.2):

1. Position the controller on the fan flange.
2. Fix the controller by means of the three socket head screws M5 removed during the controller removal.
3. Turn the pump.
4. Plug the line cord and the pump cable to the pump.
5. Connect the fan supply cable to the suitable connector of the controller as detailed in the following paragraph.
6. Connect the mating connector with the jumper on the interlock signal to start the pump.

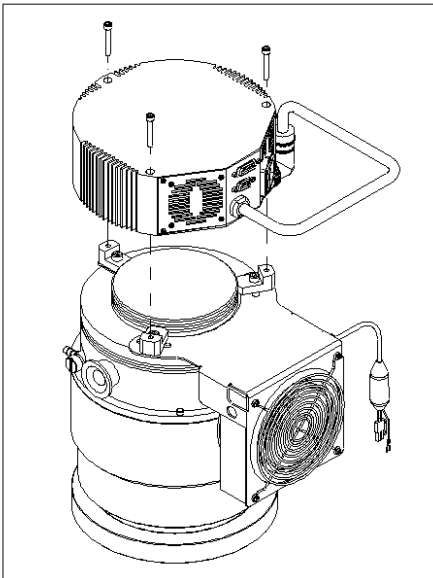


Fig. 1.5.1 TV 551/701 Navigator controller installation

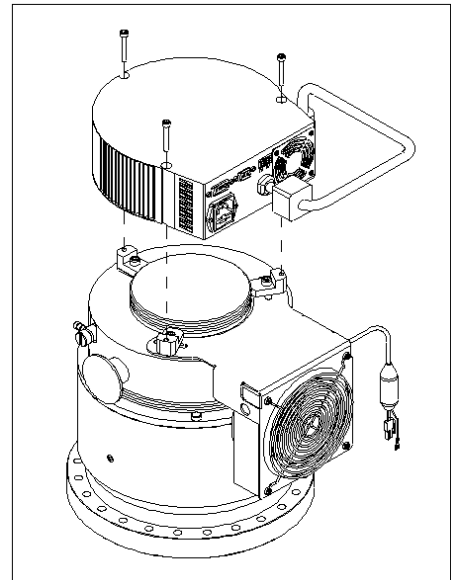


Fig. 1.5.2 TV 1001 Navigator controller installation

### 1.6 Fan to Controller Connection

The fan has one cable terminating with one 2-pins plastic male connector for the 24Vdc power supply plus a female “faston” pin for the shield.

Plug the 2-pins plastic male connector into the “FAN” named 2-pins female connector available on the controller.

If a small metal blade is present on the controller panel (shield) near the “FAN” connector, plug here the “faston” female to shield the fan’s cable, otherwise, if not present, leave it unconnected.

## 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](mailto:vpl-ra@agilent.com)  
[Lexington-service@agilent.com](mailto:Lexington-service@agilent.com)

##### EUROPE:

Toll Free: 00 800 234 234 00  
[vpt-customer-care@agilent.com](mailto: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](http://www.agilent.com/chem/vacuum)  
Representatives in most countries

This information is subject to change without notice.

© Agilent Technologies, Inc. 2026

Printed in Italy, 03/2026

87-900-966-01 (C.00)