# **Gas Purge Valve**



# 1.1 General

The gas purge valve is a protection device to protect the pump bearings against particulate and corrosive gases that could move into the pump.

This device is available with:

- Small leak
- Large leak.

Both versions are available with six different fitting combination for a total of 12 different models (see Fig. 1).

The following table summarize the models part number:

Version	Small Leak	Large leak
Fittings		
NW16	9699231	9699235
M5		
NW16	9699239	9699241
M12		
NW16	9699233	9699237
Swagelock 1/4"		
Swagelock 7/16"	9699232	9699236
Swagelock 1/4"		
Swagelock 1/4"	9699234	9699238
M5		
Swagelock 1/4"	9699240	9699242
M12		

Every valve is made with a thin calibrate sapphire orifice mounted into a cartridge, with two special PTFE membranes on each tube side.

This is to avoid any particulate, powder and any liquid to come into and occlude the thin hole.

The pieces are mounted altogether into a SST leak tight body.

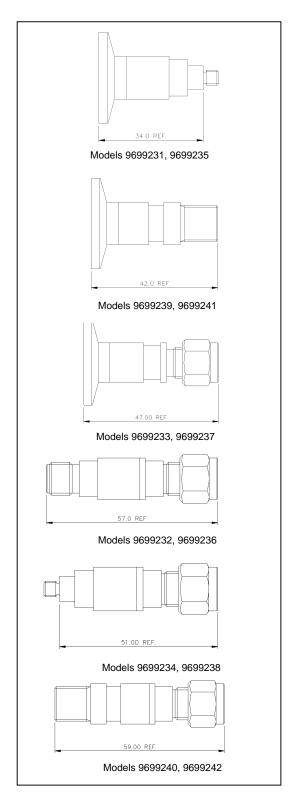


Fig. 1 - Gas Purge Valve Models

# 1.2 Gas Purge Valve Characteristics

Each version of the valve is suited for a particular application:

- the Small leak (~ 10 sccm at 1000 mbar) for mild corrosive operation when the partial pressure of the corrosive gas is maximum 10<sup>-5</sup> mbar (analytical instruments, ion implanters);
- the Large leak ( $^{\sim}$  20 sccm at 1000 mbar) for heavy corrosive operations (etching, CVD), when the partial pressure of the corrosive gas is >10<sup>-5</sup> mbar.

If you wish to use a different flowrate, you might change the supply pressure. For a given supply pressure value, the gas purge valve can give you a fixed flowrate value.

Please note that the flow amount does not depend on the vacuum side pressure of the valve, if the pressure is lower than  $\sim$  0.5 of the supply pressure.

Refers to the following diagrams to find the precise correlation. These diagrams refer to dry nitrogen and argon, for other gases you may correct the flowrate according to the calibration factor:

$$fx = \sqrt{\frac{M_x}{M_{N2}}}$$

where:

- M<sub>N2</sub> is the molecular weight of Nitrogen
- $M_x$  is the molecular weight of the other gas.

For example the gas purge flow rate for 1000 mbar of Ar for the single tube valve is:

$$Q \approx Q_{N2} \bullet fAr = Q_{N2} \bullet \sqrt{\frac{28}{40}}$$

where  $Q_{N2}$  is the corresponding nitrogen flowrate at 1000 mbar.

## 1.3 Operation Notes

Always use no less than Agilent minimum suggested purge flow rates to properly purge the pump (see the relevant pump manual). Please note that you can use all the purging flows between the minimum specified into the instruction manual and 3500 sccm, if your system is able to pump and withstand higher flow-rates.

Always operate the pump with gas purge on: during pump running, during stops, even if corrosive gases were not flowing. This provides protection against particulate that could move into the pump.

Non-observation of this basic, simple rules could affect pump warranty.

# **NOTE**

The Soft start is recommended for the installation of the pump and may be disabled when the pump is continuously operated and the interval between different run is not more than 1 week.

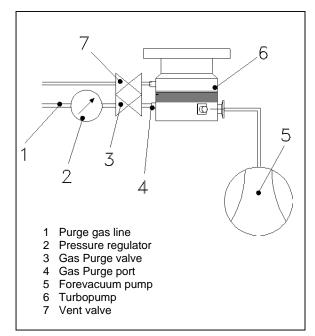


Fig. 2 - Installation diagram

87-900-890-01(B) Gas Purge Valve Page 2/4

### 1.4 Vent Procedure

Venting may be done in three different ways:

**Through the vent port** (using vent valve or vent device).

In this case the only precaution required is to keep the gas purge port on even during vent procedure.

# Through the gas purge port

The maximum air flow allowed for venting through the gas purge port is 3500 sccm (58.3 mbarl/sec). If the required flow for venting the turbopump is higher than the flow allowed by the gas purge valve, a "T" connector can be put on the purge port. On one side is connected the vent valve, and on the other side is connected the gas purge valve.

# **CAUTION**

When venting is performed through the gas purge port, be sure that the foreline pump is on, in order to avoid overpressure.

# 1.5 Through the chamber

The vent flow rate into the chamber must be less than three times the chamber volume (in litres) multiplied by the flow rate through the gas purge port. This is done to maintain the pressure into the pump body higher than into the chamber in order to avoid any corrosive or powder flowback.

E.g. chamber volume: 250 litres; gas purge flow: 20 sccm (0.33 mbarl/sec)

chamber venting flow rate = 3x250x20 = 15000 sccm

and the venting time will be:

chamber venting =

- = (chamber volumex1000)/chamber flow rate=
- $= (250 \times 1000)/15000 = 17 \text{ min.}$

# 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.



87-900-890-01(B) Gas Purge Valve Page 3/4

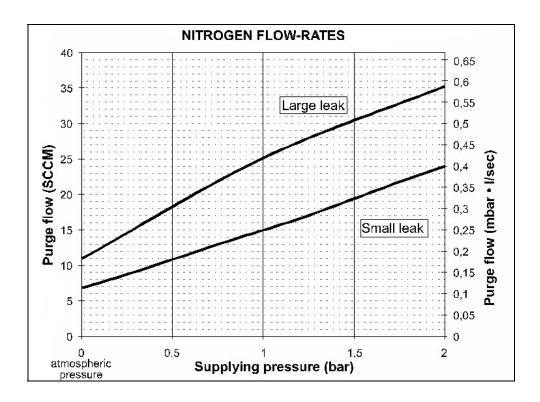


Fig. 4 - Gas Purge Valve flow rate vs supplying pressure for Nitrogen

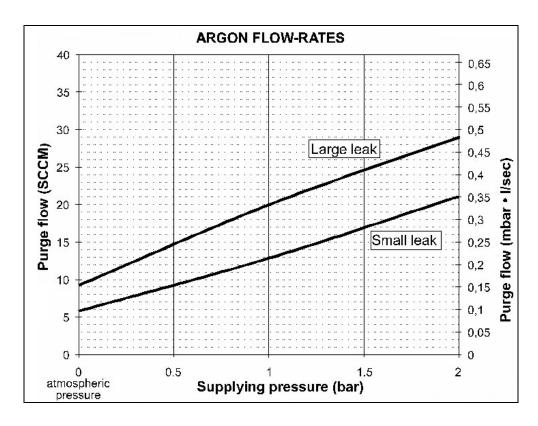


Fig. 3 - Gas purge Valve flow rate vs supplying pressure for Argon

87-900-890-01(B) Gas Purge Valve Page 4/4

#### **United States**

# **Agilent Technologies**

121 Hartwell Avenue

Lexington, MA 02421 - USA

Ph. +1 781 861 7200

Fax: +1 781 860 5437

Toll-Free: +1 800 882 7426

vpl-customercare@agilent.com

#### Netherlands

## Agilent Technologies Netherlands B.V.

**Customer Contact Center** 

Laan van Langerhuize 1, toren A-8

1186 DS Amstelveen

Tel. +31 020 547 2600

Fax +31 020 654 5748

customercare netherlands@agilent.com

#### Belaium

#### Agilent Technologies Belgium S.A./N.V.

**Customer Contact Center** 

Pegasus Park

De Kleetlaan 12A bus 12

B-1831 Diegem

Tel. +32 2 404 92 22

Fax +32 2 626 46 30

customercare belgium@agilent.com

#### Agilent Technologies Brasil

Avenida Marcos Penteado de Ulhoa

Rodrigues, 939 - 6° andar

Castelo Branco Office Park

Torre Jacarandá - Tamboré

Barueri, Sao Paulo CEP: 06460-040

Toll free: 0800 728 1405

#### China

#### Agilent Technologies (China) Co. Ltd

No.3, Wang Jing Bei Lu, Chao Yang District

Beijing, 100102, China

Tel: +86 (0)10 64397888

Fax: +86 (0)10 64392765

Toll free: 400 8206778 (mobile)

Toll free: 800 8206778 (landline)

vacuum.cnmarketing@agilent.com

vpc-customerservice@agilent.com

#### France

Agilent Technologies

Parc Technopolis - Z.A. de Courtaboeuf

3. avenue du Canada - CS 90263

91978 Les Ulis cedex, France

Tel: +33 (0) 1 64 53 61 15 Fax: +33 (0) 1 64 53 50 01

vpf.sales@agilent.com

## Southeast Asia

#### Agilent Technologies Sales Sdn Bhd

Unit 201. Level 2 uptown 2.

2 Jalan SS21/37, Damansara Uptown

47400 Petaling Jaya, Selangor, Malaysia

Ph. +603 7712 6181 Fax: +603 7727 1239

Toll free: 1 800 880 805

vps-customerservice@agilent.com

#### Agilent Vacuum Products Division/Sales and Service Offices

#### India (Sales)

## Agilent Technologies India Pvt. Ltd.

Unit Nos 110-116, & Part of 101 & 109

First Floor, Splendor Forum,

Plot No.-3, District Centre, Jasola

New Delhi-110025

Ph. +91 11 4623 7100

Fax: +91 4623 7105 Toll Free: 1 800 180 1517

## Agilent Technologies Italia S.p.A.

Via F.Ili Varian, 54

10040 Leini, (Torino) - Italy

Tel: +39 011 9979 111

Fax: +39 011 9979 350 Toll free: 00 800 234 234 00

vpt-customerservice@agilent.com

#### .lanan

# Agilent Technologies Japan, Ltd.

9-1 Takakura-cho Hachioji-city,

Tokyo, Japan

Tel.: +81- 3-5232-1253

Fax: +81-120-565-154

Toll-Free: +81-120-477-111

jp-vvt-sales.pdl-ext@agilent.com

#### Singapore

#### Agilent Technologies Singapore Pte. Ltd,

1 Yishun Avenue 7,

Singapore 768923

Tel: (65) 6215 8045

Fax: (65) 6754 0574

Toll free: 1 800 2762622 vps-customerservice@agilent.com

#### Korea

## Agilent Technologies Korea, Ltd.

**Ilshin Builiding 4F** 

Yongsan-gu Hannam-daero

Seoul Korea 04418

Tel: +82 (0)2 2194 9449

Fax: +82 (0)2 2194 9853

Toll free: 080 222 2452

vpk-customerservice@agilent.com

## **UK and Ireland**

# **Agilent Technologies LDA UK Limited**

Lakeside Cheadle Royal Business Park

Cheadle, Cheshire SK8 3GR.

United Kingdom

Ph. +44 01865291570

Fax +44 01865291571

Toll free: 00 800 234 234 00

Toll free fax: 00 800 345 345 00 vpt-customercare@agilent.com

© Agilent Technologies, Inc. 2018 07/2018



#### India (Service)

Agilent Technologies India Pvt. Ltd.

C-Block, RM7 Centennial Plot Number- 8A, 8B, 8C,

Doddanakundi Industrial Area, ITPL Road, Mahadevapura Post, Bangalore- 560048

Ph. +91 80 40614000

Fax: +91 80 40148991

#### Taiwan

## **Agilent Technologies Taiwan Limited**

No. 20 Gao-shuang Road, Ping-zhen Dist

Tao-Yuan City

32450 Taiwan, R.O.C.

Tel: +886 3 4959004

Toll free: 0800 018 768

vpw-customerservice@agilent.com

## **Germany and Austria**

**Agilent Technologies** 

Sales & Services GmbH & Co. KG

Lyoner Str. 20

60 528 Frankfurt am Main

**GERMANY** 

Tel: +49 69 6773 43 2230

Fax: +49 69 6773 43 2250

#### Mexico

## **Agilent Technologies**

Concepcion Beistegui No 109

Col Del Valle

C.P. 03100 - Mexico, D.F.

Tel.: +52 5 523 9465 Fax: +52 5 523 9472

# Other Countries

# Agilent Technologies Italia S.p.A.

Via F.Ili Varian, 54

10040 Leini, (Torino) - Italy

Tel.: +39 011 997 9111

Fax: +39 011 997 9350

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

# **Customer Support & Service**

NORTH AMERICA:

Toll Free: 800 882 7426

vpl-ra@agilent.com Lexington-service@agilent.com

# FUROPF:

Toll Free: 00 800 234 234 00

vpt-customercare@agilent.com

# PACIFIC RIM:

please visit our website for individual office information

http://www.agilent.com/chem/vacuum

# Worldwide Web Site, Catalog and Order On-

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