



Agilent Technologies

Vacuum Products Division

CE

Block Valves
All Sizes

INSTRUCTION MANUAL

Manual No. 699912040
Revision L
September 2004

Block Valves: Instruction Manual for All Sizes



ConFlat Flange is a registered trademark of Agilent, Inc.

Klamp-Flange and KF Clamp are registered trademarks of Agilent, Inc.

Krytox is a registered trademark of DuPont Company

Loctite is a registered trademark of Loctite Corporation.

Scotch-Brite is a trademark of 3M.

Viton is a registered trademark of E. I du Pont de Nemours and Company.

Warranty

Products manufactured by Seller are warranted against defects in materials and workmanship for twelve (12) months from date of shipment thereof to Customer, and Seller's liability under valid warranty claims is limited, at the option of Seller, to repair, to replace, or refund of an equitable portion of the purchase price of the Product. Items expendable in normal use are not covered by this warranty. All warranty replacement or repair of parts shall be limited to equipment malfunctions which, in the sole opinion of Seller, are due or traceable to defects in original materials or workmanship. All obligations of Seller under this warranty shall cease in the event of abuse, accident, alteration, misuse, or neglect of the equipment. In-warranty repaired or replaced parts are warranted only for the remaining unexpired portion of the original warranty period applicable to the repaired or replaced parts. After expiration of the applicable warranty period, Customer shall be charged at the then current prices for parts, labor, and transportation.

Reasonable care must be used to avoid hazards. Seller expressly disclaims responsibility for loss or damage caused by use of its Products other than in accordance with proper operating procedures.

Except as stated herein, Seller makes no warranty, express or implied (either in fact or by operation of law), statutory or otherwise; and, except as stated herein, Seller shall have no liability under any warranty, express or implied (either in fact or by operation of law), statutory or otherwise. Statements made by any person, including representatives of Seller, which are inconsistent or in conflict with the terms of this warranty shall not be binding upon Seller unless reduced to writing and approved by an officer of Seller.

Warranty Replacement and Adjustment

All claims under warranty must be made promptly after occurrence of circumstances giving rise thereto, and must be received within the applicable warranty period by Seller or its authorized representative. Such claims should include the Product serial number, the date of shipment, and a full description of the circumstances giving rise to the claim. Before any Products are returned for repair and/or adjustment, written authorization from Seller or its authorized representative for the return and instructions as to how and where these Products should be returned must be obtained. Any Product returned to Seller for examination shall be prepaid via the means of transportation indicated as acceptable by Seller. Seller reserves the right to reject any warranty claim not promptly reported and any warranty claim on any item that has been altered or has been returned by non-acceptable means of transportation. When any Product is returned for examination and inspection, or for any other reason, Customer shall be responsible for all damage resulting from improper packing or handling, and for loss in transit, notwithstanding any defect or non-conformity in the Product. In all cases, Seller has the sole responsibility for determining the cause and nature of failure, and Seller's determination with regard thereto shall be final.

If it is found that Seller's Product has been returned without cause and is still serviceable, Customer will be notified and the Product returned at Customer's expense; in addition, a charge for testing and examination may be made on Products so returned.

3/1/00

This page intentionally left blank.

Contents

Preface	xi
Document Conventions	xi
Contacting Agilent	xi
Declaration of Conformity	
Introduction and Installation	1-1
Installation	1-3
Electrical Connections to Solenoid	1-4
Operation	1-5
Manually-Operated Valves.....	1-5
Pneumatically-Operated Valves	1-5
Maintenance.....	1-6
Manual Valves	1-6
Pneumatic Valves.....	1-11
Appendix A. Technical Specifications	A-1
Appendix B. Spare Parts Kits/Accessories and Replacement Parts	B-1
Request for Return Health and Safety Certification	
Sales and Service Offices	

This page intentionally left blank.

List of Figures

Figure	Caption	Page
1-1	KF Coupling.....	1-3
1-2	Solenoid Assembly.....	1-4
1-3	Solenoid Operation.....	1-5
1-4	Small and Large Manual Block Valve – Exploded View.....	1-6
1-5	Small and Large Pneumatic Block Valve – Exploded View.....	1-11
1-6	Positioning the Quad Ring in the Air Cylinder.....	1-15
1-7	Removing the Stem from the Seal Disc.....	1-15
1-8	Position of Quad Ring In its Groove.....	1-17
1-9	Position Indicator Assembly Exploded View.....	1-19
1-10	Wiring Diagram – Position Indicator.....	1-23

This page intentionally left blank.

List of Tables

Table	Title	Page
1-1	Valve Ordering Information NW16 – NW25.....	1-2
1-2	Valve Ordering Information NW50 – NW80.....	1-2
1-3	Small Block Valves: Manual Valve Parts List	1-7
1-4	Large Block Valves: Manual Valve Parts List	1-7
1-5	Small Valves: Pneumatic Valve Parts List	1-12
1-6	Large Block Valves: Pneumatic Valve Parts List.....	1-13
1-7	Position Indicator Parts List	1-20
1-8	Connector Base Wiring	1-21
A-1	Block Valve Specifications	A-1
B-1	Spare Part Kits	B-1
B-2	KF Clamp [®] and Centering Ring Ordering Information	B-2

This page intentionally left blank.

Preface

Document Conventions

This manual uses the following standard safety protocols:

WARNING



The warning messages are for attracting the attention of the operator to a particular procedure or practice which, if not followed correctly, could lead to serious injury.

CAUTION



The caution messages are displayed before procedures, which if not followed, could cause damage to the equipment.

NOTE



The notes contain important information.

This product must only be operated and maintained by trained personnel.

Before operating or servicing equipment, read and thoroughly understand all operation/maintenance manuals provided by Agilent. Be aware of the hazards associated with this equipment, know how to recognize potentially hazardous conditions, and how to avoid them. Read carefully and strictly observe all cautions and warnings. The consequences of unskilled, improper, or careless operation of the equipment can be serious.

In addition, consult local, state, and national agencies regarding specific requirements and regulations. Address any safety, operation, and/or maintenance questions to your nearest Agilent office.

Contacting Agilent

In the United States, you can contact Agilent Customer Service at 1-800-882-7426. See the back cover of this manual for a listing of our sales and service offices.

Visit our web site at: <http://www.chem.agilent.com/en-US/Products/Instruments/vacuum/pages/default.aspx>.

Declaration of Conformity
Konformitätserklärung
Déclaration de Conformité
Declaración de Conformidad
Verklaring de Overeenstemming
Dichiarazione di Conformità



Agilent Technologies

We
Wir
Nous
Nosotros
Wij
Noi

Agilent, Inc.
121 Hartwell Avenue
Lexington, MA, 02421-3133 USA

declare under our sole responsibility that the product,
erklären, in alleiniger Verantwortung, daß dieses Produkt,
déclarons sous notre seule responsabilité que le produit,
declaramos, bajo nuestra sola responsabilidad, que el producto,
verklaren onder onze verantwoordelijkheid, dat het product,
dichiariamo sotto nostra unica responsabilità, che il prodotto,

Aluminum GateKeeper Valves	NW63, NW100, NW160, NW250,
Aluminum Block Valves	NW16, NW25, NW40, NW50, NW63, NW80,
Stainless Steel Block Valves	NW16, NW25, NW40,
Stainless Steel Tube Valves	NW16, NW25, NW40

to which this declaration relates is in conformity with the following standard(s) or other normative documents.
auf das sich diese Erklärung bezieht, mit der/den folgenden Norm(en) oder Richtlinie(n) übereinstimmt.
auquel se réfère cette déclaration est conforme à la (aux) norme(s) ou au(x) document(s) normatif(s).
al que se refiere esta declaración es conforme a la(s) norma(s) u otro(s) documento(s) normativo(s).
waamaar deze verklaring verwijst, aan de volgende norm(en) of richtlijn(en) beantwoordt.
a cui se riferisce questa dichiarazione è conforme alla/e seguente/i norma/o documento/i normativo/i.

98/37/EEC, Machinery Directive

EN 60204-1 Electrical equipment of industrial machines; general requirements

John Ehmann
General Manager
Agilent, Inc.
Vacuum Products Division
Lexington, Massachusetts, USA

September 2010



Section 1. Introduction and Installation

Agilent small right-angle and in-line block valves are compact bellows-sealed valves for use in roughing and high vacuum applications. The valve body is made of aluminum or stainless steel, and incorporates Klamp-Flange[®], CFF or tube end fittings to provide reliable and convenient vacuum connections. The main seal and all vacuum seals are Viton[®].

The valves are available in manually-operated and pneumatically-operated versions. The pneumatic valve is also available with a position indicator. The position indicator delivers an electrical output signal when the valve cycles, and is user-set in either the normally-open or normally-closed position.

The NW 16 and NW 25 valves are identical except for their port sizes; therefore, all internal spare parts are interchangeable. The NW 40 valve, however, is a larger-bodied valve requiring different internal components.

The NW 50, NW 63 and NW 80 are available only as stainless steel, right-angle valves.

This manual contains the necessary information to install, operate and maintain the Agilent block valves listed in Table 1-1 and Table 1-2 on page 1-2.

Refer to the following sections:

- ❑ “Installation” on page 1-3
- ❑ “Operation” on page 1-5
- ❑ “Maintenance” on page 1-6

Block Valves: Instruction Manual for All Sizes

Table 1-1 Valve Ordering Information NW16 – NW25

Valve	Material	Right-Angle, All Models Part No.	In-Line, All Models Part No.
Manually-operated	Aluminum	L6280	L9180
	Stainless Steel	L9480	L9580
Pneumatically-operated (with or without solenoid, without position indicator)	Aluminum	L6281	L9181
	Stainless Steel	L9481	L9581
Pneumatically-operated (with or without solenoid, with position indicator)	Aluminum	L6282	L9182
	Stainless Steel	L9482	L9582

Table 1-2 Valve Ordering Information NW50 – NW80

Valve	Material	Right-Angle, All Models Part No.
Manually-operated	Aluminum	L72803
Pneumatically-operated (with or without solenoid, with position indicator)	Aluminum	L72823

1.1 Installation

All valves are installed in a vacuum system by brazing welding, Conflat[®] gaskets, or Klamp-Flange[®] or KF couplings.

KF coupling Consists of two symmetrical flange fittings (one of which is on the valve body), a centering ring, an O-ring and a flange clamp (Figure 1-1).

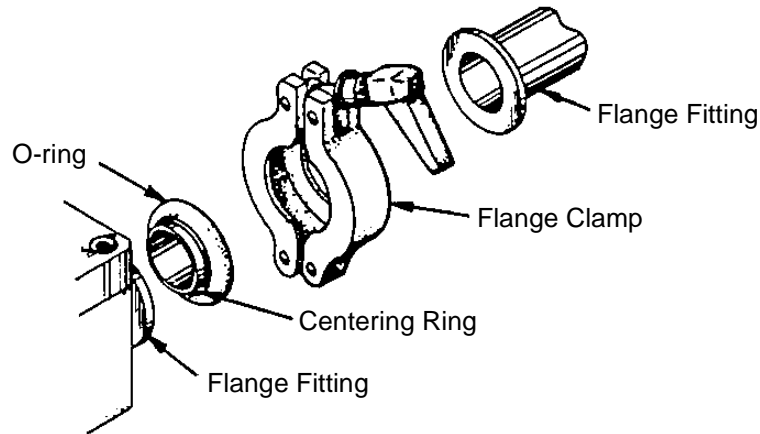


Figure 1-1 KF Coupling

NW 50 valves Installed in a vacuum system by means of a Klamp[®] (Quick Clamp) flange or ISO coupling.

ISO 63 and ISO 80 valves Installed using an ISO coupling consisting of two symmetrical flanges, a centering ring, and claw clamps.

To install the valve:

1. Wipe the O-ring clean and grease it very lightly with Dupont Krytox[®] GPL2O7 vacuum lubricant (Part No. 695400010) before installation.
2. Assemble the KF coupling and tighten the wing nut or ratchet firmly to ensure proper compression of the O-ring seal.



The pneumatic valve uses a 1/8 NPT inlet fitting and requires a minimum of 80 psig air pressure to operate.

3. Install a filter and lubricator in the air inlet line (pneumatic valve only).

1.1.1 Electrical Connections to Solenoid

To supply power to the solenoid:

1. Remove the connector from the solenoid by removing the holding screw and pulling the connector away from the solenoid.
2. Remove the terminal block from its housing by inserting a small, flat-head screwdriver in the slot provided at the terminal base and prying the housing off (Figure 1-2).

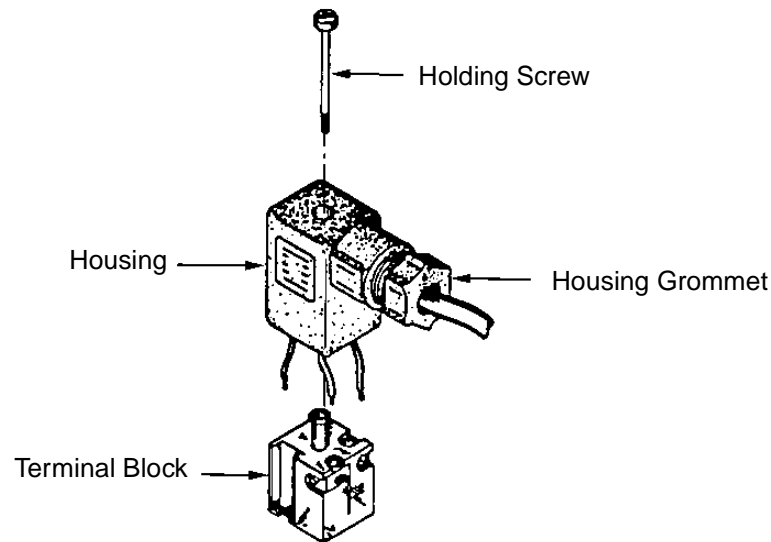


Figure 1-2 Solenoid Assembly

3. Feed the power line through the housing grommet and connect the three leads to the terminal contacts which are clearly labeled 1, 2, and $\underline{\underline{1}}$ (ground).
4. Push the terminal block back into the housing and snap it in place.



The housing can be rotated 360° in 90° increments so that the power line exits the connector in the most convenient direction.

5. Plug the connector back into the solenoid and tighten the holding screw.

1.2 Operation

This section discusses the operation of:

- ❑ “Manually-Operated Valves”
- ❑ “Pneumatically-Operated Valves”

1.2.2 Manually-Operated Valves

The manual valve operates by means of a hand-actuated knob.

To close the valve:

- ❑ Turn the knob in a clockwise direction. To properly compress the main O-ring seal and obtain a good vacuum seal, turn the knob clockwise until a slight resistance is felt, then firmly torque the knob approximately $\frac{1}{4}$ turn.

To open the valve:

- ❑ Turn the knob counterclockwise until the valve is fully open to ensure maximum conductance.

1.2.3 Pneumatically-Operated Valves

Pneumatic valves are spring-closed air-opened valves.

When electric power is:

- ❑ *Supplied* to the solenoid valve, air is supplied to the cylinder and the valve opens.
- ❑ *Removed*, the valve closes (Figure 1-3).

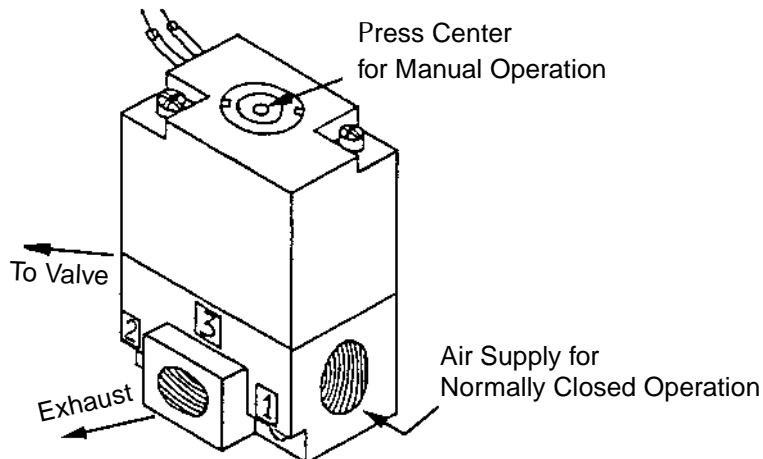


Figure 1-3 Solenoid Operation

1.3 Maintenance

This section discusses the maintenance of:

- “Manual Valves”
- “Pneumatic Valves” on page 1-11

1.3.4 Manual Valves

Manual valve maintenance consists of:

- “Disassembly and Cleaning” on page 1-8
- “Reassembly” on page 1-10

Figure 1-4 shows an exploded view of the small manual block valve and the large block valve. The numbers of the small block valve correspond to items in Table 1-3. The numbers of the large block valve correspond to items in 1-4.

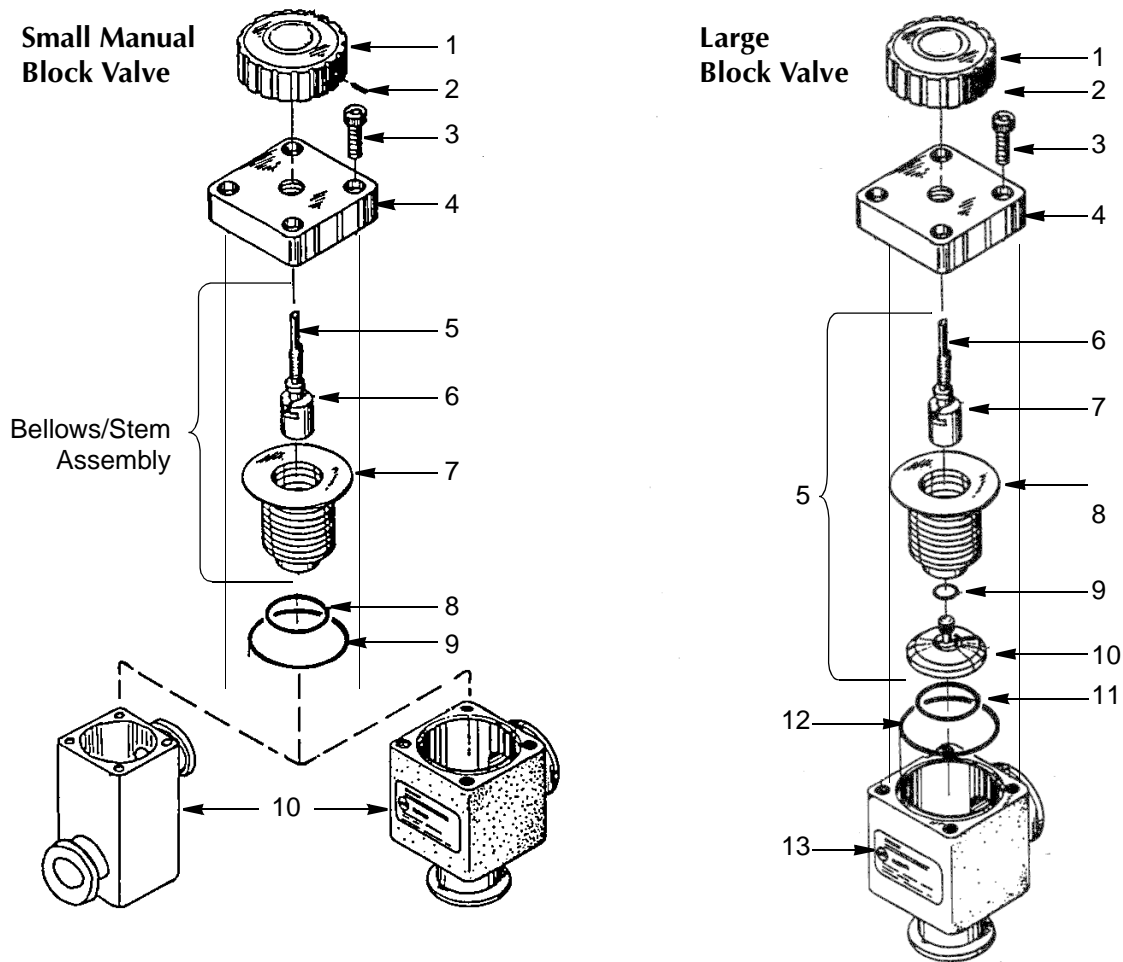


Figure 1-4 Small and Large Manual Block Valve – Exploded View

Block Valves: Instruction Manual for All Sizes

Table 1-3 Small Block Valves: Manual Valve Parts List

Item	Description	Agilent Part No. or Equivalent	
		NW16 or NW25	NW40
1	Knob	L5967301 (aluminum)	L5698301 (aluminum)
		L8819001 (plastic)	L8818001 (plastic)
2	Set Screw	#10-32 x 5/8 lg	#10-32 x 1/2 lg
3	Screw, Allen head cap	#10-32 x 5/8 lg	#10-32 x 3/4 lg
4	Top Plate	L5491001	L5686001
5	Valve Stem	L5490001	L9684001
6	Bellows Nut	85073001	85073001
7	Bellows and Seal Disc	L8739301	L8738301
8	O-ring, Viton	Parker No.2-213	Parker No.2-223
9	O-ring Viton	Parker No.2-031	Parker No 2-039
10	Valve Body	–	–

Table 1-4 Large Block Valves: Manual Valve Parts List

Item	Description	Agilent Part No. or Equivalent		
		NW50	ISO63	ISO80
1	Knob	L8819001 (plastic)	L8818001 (plastic)	L8818001 (plastic)
2	Set Screw	#10-32 x 1/2 lg	#10-32 x 1/2 lg	#10-32 x 1/2 lg
3	Screw Socket, Head Cap	#10-32 x 3/4 lg	#10-32 x 3/4 lg	#10-32 x 3/4 lg
4	Top Plate	L6606001	L6606001	L7283001
5	Bellow/Valve Stem Assembly	Ref	Ref	Ref
6	Valve Stem	L6608001	L6608001	L6606001
7	Bellows Nut	L6609001	L6609001	L6609001
8	Bellows	L6603001	L6603001	L6603001
9	O-ring, Bellows, Viton	66892014	66892014	66892014
10	Seal Disc	L6602001	L6985001	L9972001
11	O-ring, Seal Disc, Viton	660892227	660892231	660892234
12	O-ring, Bonnet, Viton	66892043	660892043	660892043

Block Valves: Instruction Manual for All Sizes

Table 1-4 Large Block Valves: Manual Valve Parts List (Continued)

Item	Description	Agilent Part No. or Equivalent		
		NW50	ISO63	ISO80
13	Valve Body	L6600001	L9970401	L9971401
	Valve Body, Nickel-plated	L6600003	L9970003	L9971003

1.3.4.1 Disassembly and Cleaning

Refer to Figure 1-4 on page 1-6 for this procedure.



In this procedure, the components (shown in Figure 1-4, identified in Table 1-3 and Table 1-4) are enclosed in parentheses using the following convention:

- A single number refers to both the small and large valve.*
- In a set of numbers separated by a semi-colon:
 - The first item number identifies an item in the small block valve.*
 - The second item number identifies an item in the large block valve.**

To disassemble and clean the valve:

1. Loosen the set screw (2) with a 3/32 Allen wrench and pull the knob (1) off the valve stem (5; 6).
2. Remove the four Allen head cap screws (3) and pull the top plate (4) and bellows/valve stem assembly (7, 8, 9, and 10; 5) from the valve body as one unit. If only the main seal O-ring (8) is to be replaced, do so at this time and then reassemble the valve ("Reassembly" on page 1-10).
3. Holding the valve stem (5; 6) of the bellows/valve stem assembly ((8, 9, and 10; 5) between the fingers of one hand, unscrew the top plate (4) with the other hand and remove the plate from the valve stem (5; 6).

The bellows/valve stem assembly (7, 8, 9, and 10; 5) is treated as a single unit; its separate components, however, are listed on the parts list.

4. Remove the O-ring (8; 11) from the seal disc and the O-ring (9; 12) from the valve body (10; 13). Use plastic tweezers or a similar non-destructive instrument to remove the O-rings so as not to mar O-ring groove surfaces.

Block Valves: Instruction Manual for All Sizes

5. Clean the internal metal components using an alcohol-dampened cloth or a light abrasive cloth such as Scotch-Brite™.

CAUTION



When cleaning the parts of the bellows/valve stem assembly (7, 8, 9, and 10; 5) , wipe only the surface area of the bellows and seal disk exposed to the vacuum system. Soaking the bellows/valve stem assembly in an alcohol bath damages the O-ring (8) and causes it to outgas under vacuum.

1.3.4.2 Reassembly

Refer to Figure 1-4 on page 1-6 for this procedure.



In this procedure, the components (shown in Figure 1-4, identified in Table 1-3 and Table 1-4) are enclosed in parentheses using the following convention:

- ❑ *A series of numbers not separated by a semi-colon refer to both the small and large valve.*
- ❑ *In a series of numbers separated by a semi-colon:*
 - ❑ *The first item number identifies an item in the small block valve.*
 - ❑ *The second item number identifies an item in the large block valve.*

To reassemble the manual valve:



Use new O-rings (8 and 9; 9, 11 and 12) for this procedure.

1. Wipe new O-rings clean and very lightly grease them with Dupont Krytox[®] GPL2O7 vacuum lubricant.
2. Place O-rings (8 and 9) into their respective grooves.
3. Take the bellows/valve stem assembly (7, 8, 9, and 10; 5) and hold it securely in one hand. Using the other hand, thread the top plate (4) onto the valve stem (5; 6) until the top plate is flush with the bellows.
4. Place the bellows/valve stem assembly (7, 8, 9, and 10; 5) with the top plate into the valve body and secure with cap screws (3). Be sure the seal disc is not in contact with the seal area of the body before tightening the cap screws.
5. Push the knob onto the valve stem and secure it by tightening the set screw on the flat of the valve stem(5; 6).

1.3.5 Pneumatic Valves

Pneumatic valve maintenance consists of:

- ❑ “Disassembly and Cleaning” on page 1-14
- ❑ “Reassembly” on page 1-16
- ❑ “Position Indicator” on page 1-18 – discusses installation and servicing of the position indicator and microswitches

Figure 1-5 shows an exploded view of the small pneumatic block valve and the large pneumatic block valve. The numbers of the small block valve correspond to the items in Table 1-5. The numbers of the large block valve correspond to the items in Table 1-6 on page 1-13.

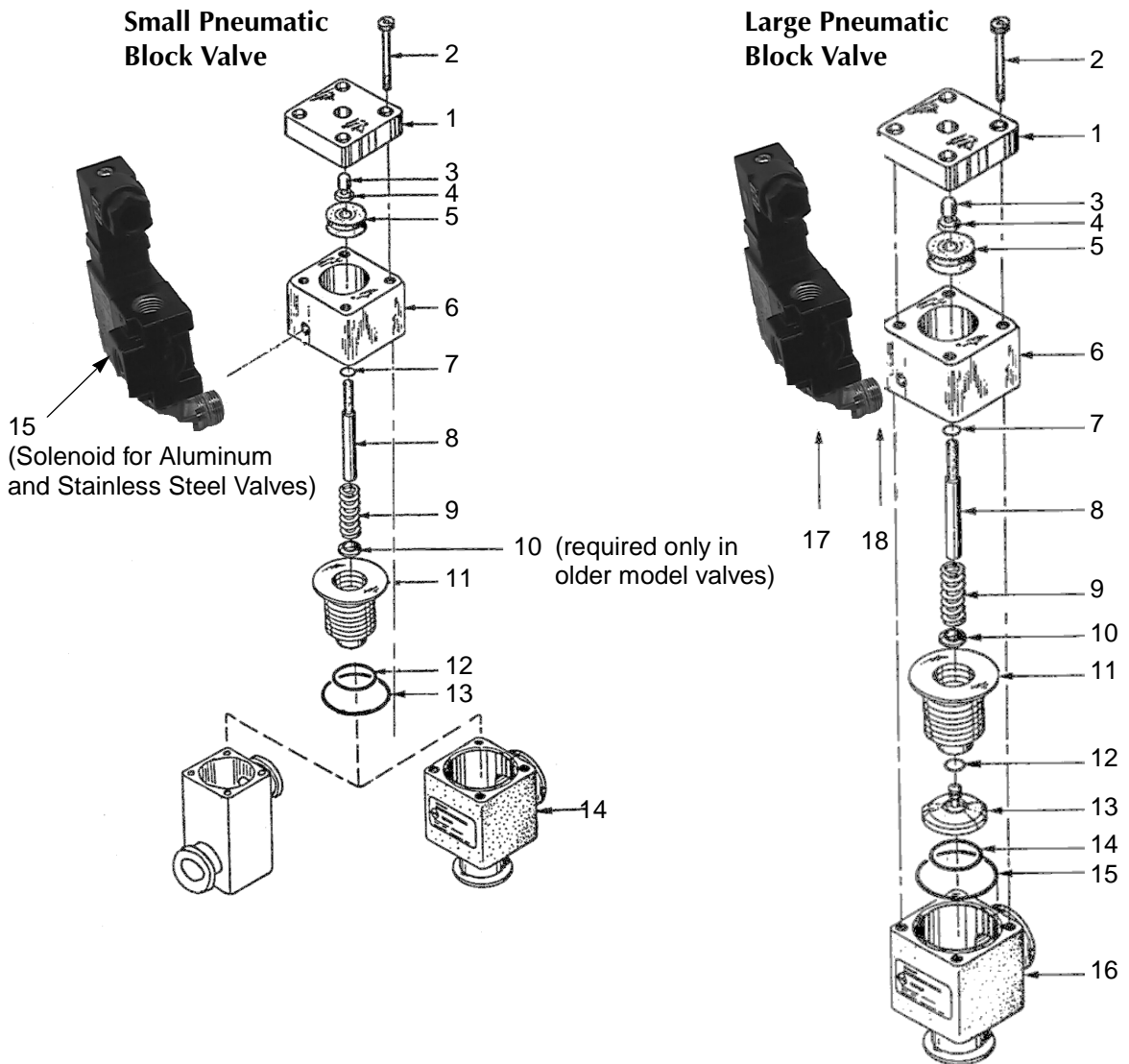


Figure 1-5 Small and Large Pneumatic Block Valve – Exploded View

Block Valves: Instruction Manual for All Sizes

Table 1-5 Small Valves: Pneumatic Valve Parts List

Item	Description	Agilent Part No. or Equivalent	
		NW16 or NW25	NW40
1	Air cylinder Cover	L5484001	L5674001
2	Screw, Allen head cap	#10-32x 2 lg	#10-32 x 2¼ lg
3	Visual Indicator Cap	L5673-001	L5673001
4	Stop Nut	ESNA 29NTE 040	ESNA 29NTE 040
5	Piston Cup	660202114	660203125
6	Air Cylinder	L5480001	L5677001
7	Quad Ring	699000025	699000025
8	Valve Stem	L5483001	L5679001
9	Spring	660285312	660265073
10	Special Washer, Aluminum (not used on new models)	L5482001	L5678001
11	Bellows & Seal Disc	L8739301	L8738301
12	O-ring	Parker No. 2-213, Viton	Parker No. 2-223, Viton
13	O-ring	Parker No. 2-031, Viton	Parker No. 2-039, Viton
14	Valve Body	–	–
		Voltage/Hertz	Part No.
15	Solenoid kit for aluminum (solenoid valve and brass elbow) and stainless steel valves (solenoid valve and straight brass fitting)	110/115 V, 50/60 Hz	BVSOLKIT110
		220/240 V, 50/60 Hz	BVSOLKIT220
		24 VDC	BVSOLKIT24D

Block Valves: Instruction Manual for All Sizes

Table 1-6 Large Block Valves: Pneumatic Valve Parts List

Item	Description		Agilent Part No. or Equivalent		
			NW50	ISO63	ISO80
1	Air Cylinder Cover		L6612001	L6612001	L9974001
2	Screw Allen Head Cap		1/4-20 x 21/4 lg	1/4-20 x 21/4 lg	1/4-20 x 21/4 lg
3	Visual Indicator Cap		L6611001	L6611001	L6611001
4	Stop Nut		617935016	617935016	617935016
5	Piston Cup		660203140	660203140	660203140
6	Air Cylinder		L6604001	L660400	L6604001
7	Quad Ring		699000026	699000026	699000026
8	Valve Stem		L6607001	L6607001	L6607001
9	Spring		660285698	660285698	660285698
10	Special Washer		L9973001	L9973001	L9973001
11	Bellows		L6603001	L6603001	L6603001
12	O-ring, Bellows, Viton		66892014	66892014	66892014
13	Seal Disc		L6602001	L6985001	L9972001
14	O-ring, Seal Disc, Viton		660892227	660892231	660892234
15	O-ring, Bonnet, Viton		660892043	660892043	660892043
16	Valve Body		L6600401	L9970401	L9971401
	Valve Body, Nickel-plated		L6600003	L9970003	L9971003
17	Solenoid	110/115 V, 50/50 Hz	626771150		
		220/240 V, 50/50 Hz	626771151		
		24 VDC	626771152		
18	Elbow 1/8 NPT	623434001	623434001	623434001	623434001

1.3.5.3 Disassembly and Cleaning

Refer to Figure 1-5 on page 1-11 for this procedure.



In this procedure, the components (shown in Figure 1-5, identified in Table 1-5 and Table 1-6) are enclosed in parentheses, using the following convention:

- ❑ *A single number refers to both the small and large valve.*
- ❑ *In a set of numbers separated by a semi-colon:*
 - ❑ *The first item number identifies an item in the small block valve.*
 - ❑ *The second item number identifies an item in the large block valve.*

To disassemble and clean the valve:

1. Using a 5/8 wrench, turn the solenoid valve fitting (15; 17) counterclockwise and remove the complete solenoid valve assembly as one unit.
2. Place the solenoid valve assembly (15; 17) aside.
3. Unscrew the four cap screws (2) and remove the air cylinder cover (1).
4. Pull the air cylinder (6) away from the valve body (14; 16). All components except for the O-ring (13; 15) remain with the air cylinder.

This group of components is collectively referred to as the trim assembly. If only the main seal is to be replaced, do so at this time, then reassemble the valve. For more information, refer to "Reassembly" on page 1-16.

5. To disassemble the trim assembly, pull the red visual indicator cap (3) off the valve stem (8).
6. Using a 7/16" deep socket wrench, turn the stop nut (4) counterclockwise while gripping the bellows seal disc (11) in the other hand.
7. Remove the stop nut (4) and pull the air cylinder (6) away from the valve stem (8).
8. Proceed to:
 - ❑ Step 9 if disassembly of the air cylinder is necessary (e.g. an air leak is suspected).
 - ❑ Step 12 on page 1-15, if disassembly of the air cylinder is not required.
9. Place the air cylinder upside down on a flat surface with the large bore and piston cup insert (5) facing down.
10. Remove the piston cup insert from the air cylinder (6) by inserting the eraser end of a pencil (or similar non-destructive object) into the center bore of the air cylinder and pushing the piston cup out the other side.

11. Ensure that the air cylinder (6) is still in the upside down position, using plastic tweezers or other nondestructive instrument, pull the quad ring (7) out of the center bore of the air cylinder (Figure 1-6).



When removing the quad ring, care must be taken not to mar any metal surfaces, especially the groove in which the quad ring is positioned.

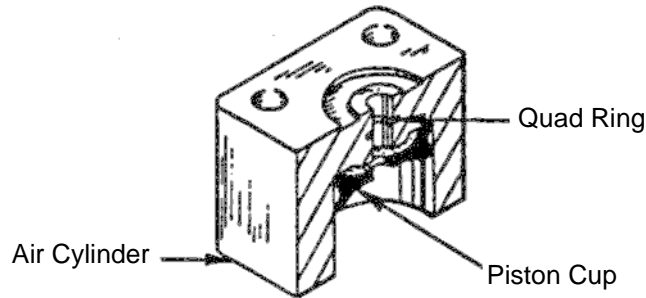


Figure 1-6 Positioning the Quad Ring in the Air Cylinder



Further disassembly requires a soft-jaw vise, two 1/4-20 nuts, and two 7/16 open end wrenches.

12. Remove the spring (9) from the valve stem (8).
13. Position the bellows/valve stem assembly in the vise with the valve stem pointed up and the seal disc between the jaws. Tighten the vise jaws on the seal disc periphery firmly but not so tight as to deform the seal disk (Figure 1-7).

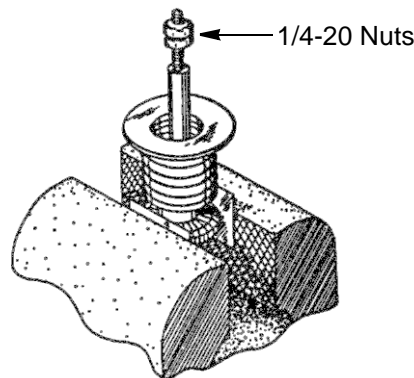


Figure 1-7 Removing the Stem from the Seal Disc

14. Thread both nuts to the midpoint of the threaded section of the stem. Using two wrenches, tighten the two nuts into each other by turning the bottom nut counterclockwise while simultaneously turning the top nut clockwise (Figure 1-7).

Block Valves: Instruction Manual for All Sizes

15. Turn the bottom nut counterclockwise using the wrench. The nuts should not rotate with respect to the stem, instead the stem rotates with respect to the seal disc. Continue turning the stem until it separates from the seal disc.
16. Remove the nuts from the stem (8).
17. Separate the remaining component: bellows and seal disc (11) by hand.
18. Using plastic tweezers or other non-destructive instrument, remove O-ring (12; 14) from the seal disc and the O-ring (13; 15) from the valve body.
19. Proceed to:
 - Step 20, if an air leak is suspected in the solenoid.
 - Otherwise, disassembly is complete.

CAUTION



When cleaning the parts of the bellows/stem assembly, wipe only the surface area of the bellows and seal disk exposed to the vacuum system. Soaking the bellows/stem assembly in an alcohol bath damages the O-ring (8) and causes it to outgas under vacuum.

20. Clean the metal components using an alcohol-dampened cloth or a light abrasive cloth such as Scotch-Brite™.
21. Wipe all surfaces with an alcohol-dampened cloth and allow components to dry thoroughly.

1.3.5.4 Reassembly

Refer to Figure 1-5 on page 1-11 for this procedure.

NOTE



In this procedure, the components (shown in Figure 1-5, identified in Table 1-5 and Table 1-6) are enclosed in parentheses, using the following convention:

- A single number or a series of numbers not separated by a semi-colon refer to both the small and large valve.*
- In a set of numbers separated by a semi-colon:*
 - The first item number identifies an item in the small block valve.*
 - The second item number identifies an item in the large block valve.*

To reassemble the valve:



When completely rebuilding the valve, be sure to reassemble with new O-rings, quad ring, and cylinder cup.

1. Wipe all O-rings (12 and 13; 14 and 15) clean and very lightly grease them with Dupont Krytox[®] GPL207 vacuum lubricant.
2. Secure the bellows and seal disc (11; 11, 12 and 13) in a vise as described in disassembly procedures, step 13.
3. Place the O-ring (12 – large valve only) into its groove on the bellows and seal disc (11).
4. Place the washer (10) into the bellow and over the seal disc threaded stud (11; 13) with the stepped side facing up. This washer is used to center the spring on bellows assemblies without a boss machined into the seal disc.
5. Take the valve stem (8) in hand, apply Loctite[®] 242 adhesive to the first three female threads.
6. Thread the valve stem (8) onto the threaded rod of the seal disc (11) by turning the stem clockwise until finger-tight.
7. Place the spring (9) over the stem so it is seated on the special washer (10). This washer is required only in older model valves.
8. Proceed to:
 - Step 9 if resassembly of the air cylinder is necessary.
 - Step 15, if resassembly of the air cylinder is not required.
9. Lubricate the quad ring (7) liberally with Sta-lube 3120 grease. The piston cup (5) has a rubber-coated bottom surface and a metal upper surface (Figure 1-8). Apply Sta-lube liberally to the side walls and the rubber-coated bottom surface, leaving the metal top surface dry.

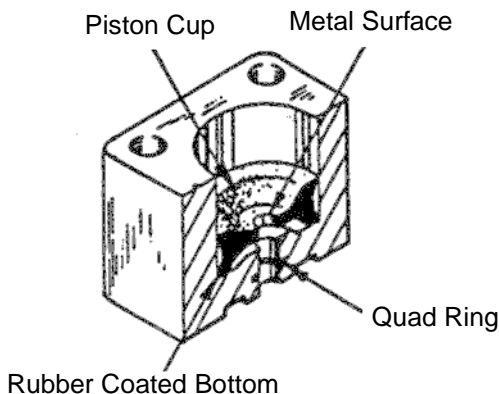


Figure 1-8 Position of Quad Ring In its Groove

10. Place the quad ring (7) in its groove.
11. Push the piston cup (5) into the air cylinder (6) with the rubber-coated side down.
12. Place the air cylinder (6) onto the valve stem (8) with the rubber side of the piston cup (5) facing up.
13. Thread the stop nut (4) onto the valve stem (8).
14. Using a torque wrench, torque the stop nut to 45 inch-pounds.
15. Push the red visual indicator cap (3) onto the valve stem (8).
16. Place the O-ring (13; 15) into its groove on the valve body.
17. Place the air cylinder/bellows assembly into the valve body.
18. Put the air cylinder cover (1) into position and secure the assembly with cap screws (2).
19. Mount the solenoid assembly to the air cylinder and tighten with a 5/8" open end wrench.

1.3.5.5 Position Indicator

This discussion details how to disassemble the position indicator so that:

- A defective microswitch can be fixed
(see "Replacing the Defective Microswitch" on page 1-21)

or

- A new position indicator can be installed using a kit
(see "Installing the Position Indicator Kit" on page 1-22).

In either case:

1. Disassemble the valve as explained in "Disassembly/Reassembly" on page 1-21.
2. Perform the procedure required.
3. Reassemble the valve as explained in "Disassembly/Reassembly" on page 1-21.
4. Once either procedure is complete, the switches can be tested (see "Test Procedure" on page 1-23).

Block Valves: Instruction Manual for All Sizes

Figure 1-9 shows an exploded view of the position indicator and Table 1-7 lists the associated callouts.

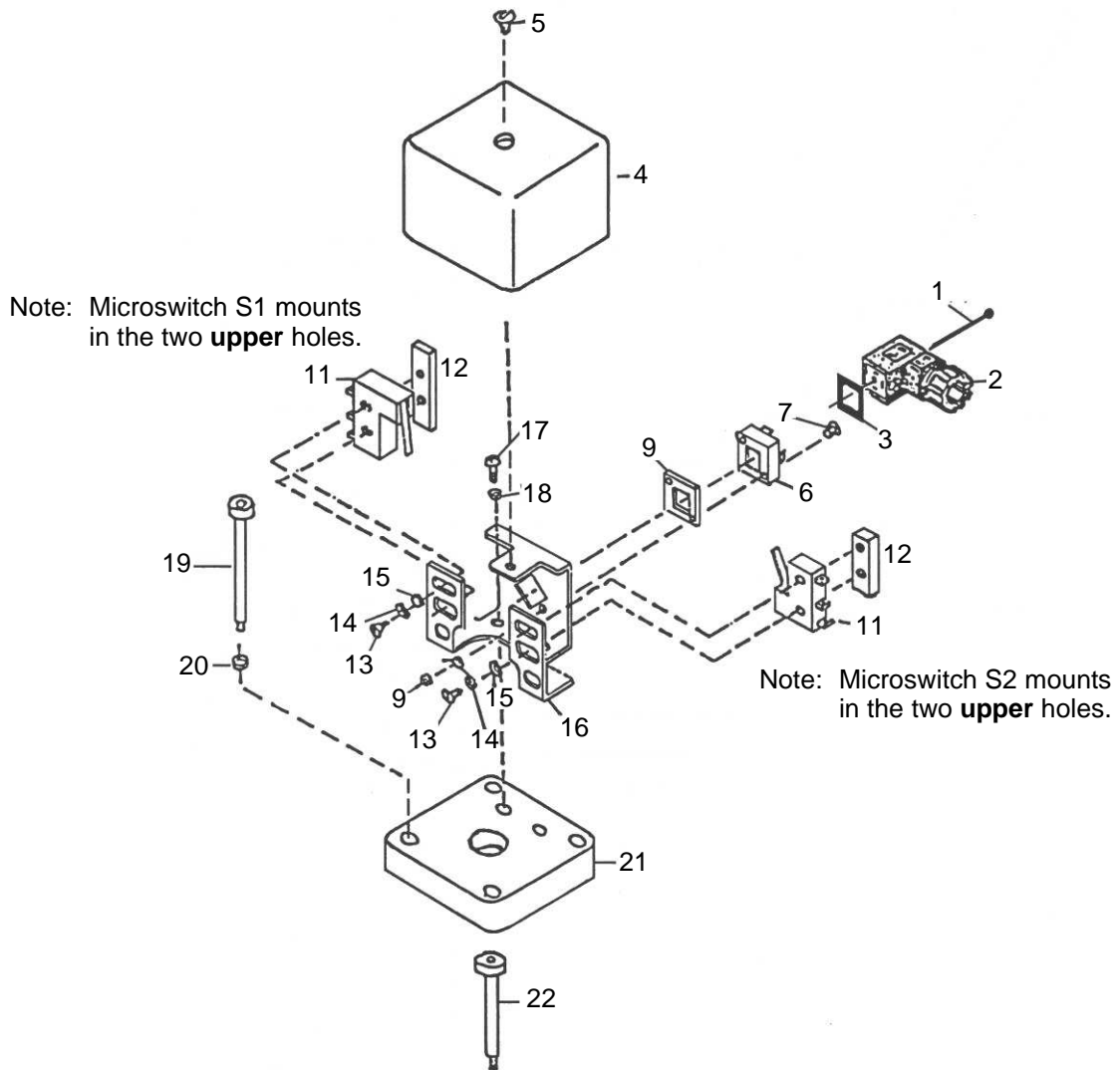


Figure 1-9 Position Indicator Assembly Exploded View

Block Valves: Instruction Manual for All Sizes

Table 1-7 Position Indicator Parts List

Item	Description	Agilent Part No. or Equivalent
1	Holding Screw	Included with item 1
2	Connector	MPM 193-07N
3	Gasket	Included with item 1
4	Position Indicator Cover	L5978001
5	Screw, Phillips head	#6-32 x 1/4 lg
6	Connector Base	MPM 193
7	Screw, Phillips head	#2-56 x 3/8 lg
8	Gasket	Included with item 6
9	Nut	#2-56
10	Lockwasher	#2
11	Microswitch	Cherry E61-10H
12	Threaded Holder	Q3033001
13	Screw, Phillips head	#2-56 x 1/2 lg
14	Lockwasher	#2
15	Washer, Flat	#2
16	Microswitch Mounting Bracket	L8673301
17	Screw, Philips head	#6-32 x 1/4 lg
18	Lockwasher	#6
19	Screw, Allen head cap	#10-32 x 2 lg (NW16, NW25, NW50, ISO63) 1/4-20 x 2 lg (NW40, ISO80)
20	Lockwasher	No.10 (NW16, NW25, NW50, ISO63) 1/4" (NW40, ISO80)
21	Air Cylinder Cover, Aluminum	L5484001 (NW16, NW25)
		L5674001 (NW40)
		L6612001 (NW50, ISO63)
		L4474001 (ISO80)
22	Screw, Shoulder, 5/16 by 1/2", Stainless Steel	L8844001

1.3.5.5.1 Disassembly/Reassembly

Refer to Figure 1-9 on page 1-19 for this procedure:

1. Remove the holding screw (1) from the connector (2) and pull the connector off the connector base (6).
2. Remove the screw (5) from the position indicator cover (4) and lift the cover off the air cylinder cover (21).


The components of the position indicator are now accessible. To reassemble the valve, reverse the steps in the procedure above.

1.3.5.5.2 Replacing the Defective Microswitch

To replace the microswitch:

1. Disconnect the soldered leads from the Microswitch pins.
2. Unscrew the mounting screws (13) and remove the Microswitch (11).
3. Reassemble a new switch on the bracket (16) as shown in Figure 1-9 on page 1-19. Note the position of the switches on the bracket and wire the microswitches as shown in Figure 1-10 on page 1-23.
4. Solder the leads to the Microswitch. Leads are wired to the connector base as indicated in Table 1-8.

Table 1-8 Connector Base Wiring

Base Pin	Wire	Dual Microswitch
1	Orange	Switch 2, NO
2	Blue	Switch 1, NO
3	White	Switch 1 and 2, Common
	Green	Ground

1.3.5.5.3 Installing the Position Indicator Kit

To install a position indicator kit:

1. Remove the cover (4) from the position indicator assembly by removing the electrical connector and the screw and lockwasher from the top cover.
2. Apply air pressure to the valve and allow the valve to remain in the open position.
3. Remove the red visual indicator cap (not shown) from the valve stem.
4. Apply a small amount of Loctite #242 to the threads of the special shoulder screw (22) before threading it into the valve stem. Install the 5/16 by 1/2" long stainless steel shoulder screw on the valve stem. Tighten the shoulder screw until it is fully threaded into the valve stem then torque it to 30 inch-pounds.
5. Remove the air pressure from the valve to close it.
6. Carefully mount the position indicator assembly on the valve. It may be necessary to move the switches back to clear the shaft. Use two #6-32 x 3/4" pan head screws and two #6 lockwashers to fasten the position indicator on the top of the valve.

1.3.5.5.3.1 Wiring the Position Indicator

Refer to Figure 1-9 on page 1-19 for this procedure.

To wire the position indicator:

1. Remove the holding screw (1) and pull the connector (2) away from the base (6).
2. Using the holding screw, push the terminal block out of the connector housing.
3. Feed the No.22 AWG wire through the housing grommet and connect the leads to the appropriate terminal block contacts for the desired operation.



The position indicator can be wired for normally-open or normally-closed operation, or both:

Normally-open The position indicator supplies an electrical output signal when the valve is open.

Normally-closed The position indicator supplies a signal when the valve is closed.

Both A signal is supplied when the valve is open or closed.

- Wire terminal contacts as detailed in Figure 1-10.

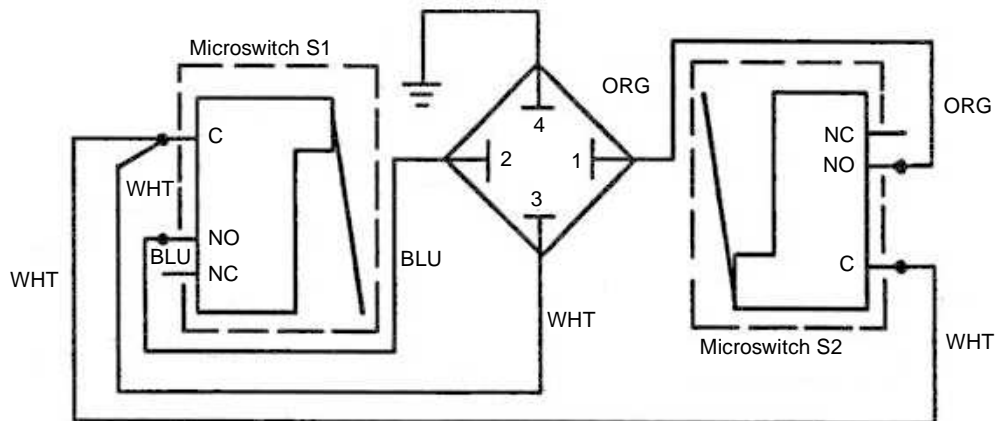


Figure 1-10 Wiring Diagram – Position Indicator

- Push the terminal block back into the connector housing and snap it into place.
- Plug the connector back into the base and secure it with the holding screw.

1.3.5.5.4 Test Procedure

Once a microswitch is replaced or a position indicator is installed, perform the test procedure for dual switches:

- Connect an Ohmmeter to pin 1 (orange wire, Figure 1-10) and pin S (white wire) of switch S2 connector, then set the meter to measure resistance (Ohms).
- With the valve in the open position (air pressure applied), adjust switch S2 (right side) until the Microswitch closes.
- Tighten the screws holding switch S2.
- Actuate the valve five times and ensure that the switch S2 opens and closes and is in the correct position.
- Close the valve by removing the air pressure from the valve.
- Connect the Ohmmeter to pin 2 (blue wire) and pin S (white wire) of the switch S1 connector. The meter should still be set to read resistance (Ohms).
- With the valve in the closed position (no air pressure), adjust switch S1 (left side) until the Microswitch closes.
- Apply air pressure to open the valve. Switch S1 should open. Readjust it if necessary by tightening the screws.
- Actuate the valve five times to ensure that the switch opens and closes and is in the correct position.
- Place the cover on the position indicator and replace the connector using the appropriate hardware.

This page intentionally left blank.

Appendix A. Technical Specifications

Table A-1 lists the block valve specifications.

Table A-1 Block Valve Specifications

	NW 16	NW 25	NW 40	NW 50	ISO 63	ISO 80
Conductance	3 liters/sec	12 liters/sec	32 liters/sec	61 liters/sec	100 liters/sec	155 liters/sec
Main Seal Leak Rate	$\leq 1 \times 10^{-9}$ std cc/sec Helium					
Body Leak Rate	$\leq 1 \times 10^{-10}$ std cc/sec Helium					
Materials						
Valve Body	Aluminum (6061-T6) or stainless steel					
Bellows	Stainless Steel T300 Series					
O-rings, vacuum	Viton V747-75					
Air Piston	Viton V747-75					
Baking Temperature						
Manual valve	80°C (176°F)			150°C (302°F)		
Pneumatic valve	80°C (176°F) without solenoid			150° C (302°F) without solenoid		
	60°C (140°F) with solenoid			60°C (140°F) with sol enoid		
Miscellaneous						
Service Life Bellows	1,000,000 cycles					
Manual valve, number of turns to open (seated to fully open)	6			9		
Pneumatic Valve						
Air connection	1/8 NPT					
Air pressure	80 psig (minimum)					
	100 psig (minimum)					
Open/close time	<1 second					
Actuation	Spring closed, air opened					
Power loss	Valve closes					

Block Valves: Instruction Manual for All Sizes

Table A-1 Block Valve Specifications (Continued)

	NW 16	NW 25	NW 40	NW 50	ISO 63	ISO 80
Position Indicator						
Actuator	Microswitch					
Electrical rating	125/250 VAC, 5 A maximum					
Signal hookup	NO or NC or both					
Microswitch life	250,000 cycles with 1 A load					

Appendix B. Spare Parts Kits/Accessories and Replacement Parts

Table B-1 lists the spare part kits by valve type.

Table B-1 Spare Part Kits

Valve Type	Valve	Part No.
Manual	NW 16 or NW 25*	L6125301
	NW 40*	L6126301
	NW 50**	L6625301
	ISO 63**	L7276301
	ISO 80**	L7278301
*Includes: Bellows, Stem, Stem nut, and O-rings **Includes: Bellows assembly, and O-rings		
Pneumatic	NW 16 or NW 25	L6127301
	NW 40	L6128301
	NW50	L6615301
	ISO 63	L7277301
	ISO 80	L7279301
Includes: Quad ring, Bellows, all O-rings, and Lubricant		

Table B-2 on page B-2 lists the accessories and replacement parts by valve type.

Block Valves: Instruction Manual for All Sizes

Table B-2 KF Clamp® and Centering Ring Ordering Information

KF Fitting	KF Clamp with Ratchet Closure Part No.	Centering Ring with Viton O-ring Part No.
NW16	KQ16AR	KC16SV
NW25	KQ25AR	KC25SV
NW40	KQ40AR	KC40SV
NW50	KQ50AR	KC50SV
ISO63	IC063100A	IC063AV
ISO80	IC063100A	IC80AV
Dupont Krytox® GPL207 Vacuum Grease		695400010
Sta-lube bearing grease		3120

Vacuum Products Division Instructions for returning products

Dear Customer:

Please follow these instructions whenever one of our products needs to be returned.

- 1) Complete the attached Request for Return form and send it to Agilent Technologies (see below), taking particular care to identify all products that have pumped or been exposed to any toxic or hazardous materials.
- 2) After evaluating the information, Agilent Technologies will provide you with a Return Authorization (RA) number via email or fax, as requested.
Note: Depending on the type of return, a Purchase Order may be required at the time the Request for Return is submitted. We will quote any necessary services (evaluation, repair, special cleaning, eg).
- 3) **Important steps for the shipment of returning product:**
 - Remove all accessories from the core product (e.g. inlet screens, vent valves).
 - Prior to shipment, drain any oils or other liquids, purge or flush all gasses, and wipe off any excess residue.
 - If ordering an Advance Exchange product, **please use the packaging from the Advance Exchange to return the defective product.**
 - Seal the product in a plastic bag, and package product carefully to avoid damage in transit. You are responsible for loss or damage in transit.
 - Agilent Technologies is not responsible for returning customer provided packaging or containers.
 - **Clearly label package with RA number.** Using the shipping label provided will ensure the proper address and RA number are on the package. Packages shipped to Agilent without a RA clearly written on the outside cannot be accepted and will be returned.
- 4) Return only products for which the RA was issued.
- 5) **Product being returned under a RA must be received within 15 business days.**
- 6) **Ship to the location specified on the printable label, which will be sent, along with the RA number, as soon as we have received all of the required information.** Customer is responsible for freight charges on returning product.
- 7) Return shipments must comply with all applicable **Shipping Regulations** (IATA, DOT, etc.) and carrier requirements.

RETURN THE COMPLETED **REQUEST FOR RETURN** FORM TO YOUR NEAREST LOCATION:

EUROPE:
Fax: 00 39 011 9979 330
Fax Free: 00 800 345 345 00
Toll Free: 00 800 234 234 00
vpt-customer@agilent.com

NORTH AMERICA:
Fax: 1 781 860 9252
Toll Free: 800 882 7426, Option 3
vpl-ra@agilent.com

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



Please read important policy information on Page 3 that applies to all returns.

1) CUSTOMER INFORMATION

Company Name:		Contact Name:	
Tel:	Email:	Fax:	
Customer Ship To:		Customer Bill To:	
Europe only: VAT reg. Number:		USA/Canada only: <input type="checkbox"/> Taxable <input type="checkbox"/> Non-taxable	

2) PRODUCT IDENTIFICATION

Product Description	Agilent P/N	Agilent S/N	Original Purchasing Reference

3) TYPE OF RETURN (Choose one from each row and supply Purchase Order if requesting a billable service)

- 3A.** Non-Billable Billable **➔** New PO # (hard copy must be submitted with this form):
- 3B.** Exchange Repair Upgrade Consignment/Demo Calibration Evaluation Return for Credit

4) HEALTH and SAFETY CERTIFICATION

AGILENT TECHNOLOGIES CANNOT ACCEPT ANY PRODUCTS CONTAMINATED WITH BIOLOGICAL OR EXPLOSIVE HAZARDS, RADIOACTIVE MATERIAL, OR MERCURY AT ITS FACILITY.

Call Agilent Technologies to discuss alternatives if this requirement presents a problem.

The equipment listed above (check one):

- HAS NOT** pumped or been exposed to any toxic or hazardous materials. OR
- HAS** pumped or been exposed to the following toxic or hazardous materials. If this box is checked, the following information must also be filled out. Check boxes for all materials to which product(s) pumped or was exposed:

- Toxic Corrosive Reactive Flammable Explosive Biological Radioactive

List all toxic/hazardous materials. Include product name, chemical name, and chemical symbol or formula:

NOTE: If a product is received at Agilent which is contaminated with a toxic or hazardous material that was not disclosed, **the customer will be held responsible** for all costs incurred to ensure the safe handling of the product, and **is liable** for any harm or injury to Agilent employees as well as to any third party occurring as a result of exposure to toxic or hazardous materials present in the product.

Print Name: _____ **Authorized Signature:** **Date:** _____

5) FAILURE INFORMATION:

Failure Mode (REQUIRED FIELD. See next page for suggestions of failure terms):
Detailed Description of Malfunction: (Please provide the error message)
Application (system and model):

I understand and agree to the terms of Section 6, Page 3/3.

Print Name: _____ **Authorized Signature:** **Date:** _____



Vacuum Products Division
Request for Return Form
(Health and Safety Certification)

Please use these Failure Mode to describe the concern about the product on Page 2.

TURBO PUMPS and TURBO CONTROLLERS

Table with 3 columns: APPARENT DEFECT/MALFUNCTION, POSITION, and PARAMETERS. Lists various failure modes like 'Does not start', 'Noise', 'Vertical', 'Horizontal', etc.

ION PUMPS/CONTROLLERS

Table listing failure modes for Ion Pumps/Controllers: Bad feedthrough, Vacuum leak, Error code on display, Poor vacuum, High voltage problem, Other.

VALVES/COMPONENTS

Table listing failure modes for Valves/Components: Main seal leak, Solenoid failure, Damaged sealing area, Bellows leak, Damaged flange, Other.

LEAK DETECTORS

Table listing failure modes for Leak Detectors: Cannot calibrate, Vacuum system unstable, Failed to start, No zero/high background, Cannot reach test mode, Other.

INSTRUMENTS

Table listing failure modes for Instruments: Gauge tube not working, Communication failure, Error code on display, Display problem, Degas not working, Other.

SCROLL AND ROTARY VANE PUMPS

Table listing failure modes for Scroll and Rotary Vane Pumps: Pump doesn't start, Doesn't reach vacuum, Pump seized, Noisy pump (describe), Over temperature, Other.

DIFFUSION PUMPS

Table listing failure modes for Diffusion Pumps: Heater failure, Doesn't reach vacuum, Vacuum leak, Electrical problem, Cooling coil damage, Other.

Section 6) ADDITIONAL TERMS

Please read the terms and conditions below as they apply to all returns and are in addition to the Agilent Technologies Vacuum Product Division – Products and Services Terms of Sale.

- Customer is responsible for the freight charges for the returning product. Return shipments must comply with all applicable Shipping Regulations (IATA, DOT, etc.) and carrier requirements.
Customers receiving an Advance Exchange product agree to return the defective, rebuildable part to Agilent Technologies within 15 business days. Failure to do so, or returning a non-rebuildable part (crashed), will result in an invoice for the non-returned/non-rebuildable part.
Returns for credit toward the purchase of new or refurbished Products are subject to prior Agilent approval and may incur a restocking fee. Please reference the original purchase order number.
Units returned for evaluation will be evaluated, and a quote for repair will be issued. If you choose to have the unit repaired, the cost of the evaluation will be deducted from the final repair pricing. A Purchase Order for the final repair price should be issued within 3 weeks of quotation date. Units without a Purchase Order for repair will be returned to the customer, and the evaluation fee will be invoiced.
A Special Cleaning fee will apply to all exposed products per Section 4 of this document.
If requesting a calibration service, units must be functionally capable of being calibrated.

This page intentionally left blank.

Agilent Technologies

Vacuum Product Division

United States & Canada

Agilent Technologies
Vacuum Products Division
121 Hartwell Avenue
Lexington, MA 02421 USA
Tel: +1 781 861 7200
Toll-Free: +1 800 882 7426
Fax: +1 871 860 5437

Benelux

Agilent Technologies
Vacuum Products Division
Herculesweg 8
4338 PL Middelburg
THE NETHERLANDS
Tel: +31 118 671570
Fax: +31 118 671569

China

Agilent Technologies
Vacuum Products Division
Room 1648
Central Tower South Wing
Beijing Junefield Plaza
No. 10 XuanWuMenWai Street
Beijing 100052 P.R. CHINA
Tel.: +86 (10) 6310 8550
Toll-Free: 800 820 6556
Fax: +86 (10) 6310 0141

France

Agilent Technologies
Vacuum Products Division
7 avenue des Tropiques
Z.A. de Courtaboeuf - B.P. 12
91941 Les Ulis cedex FRANCE
Tel.: +33 (0) 1 69 86 38 84
Fax: +33 (0) 1 69 86 29 88

Germany & Austria

Agilent Technologies
Vacuum Products Division
Alsfelder Strasse 6
Postfach 11 14 35
64289 Darmstadt GERMANY
Tel.: +49 (0) 6151 703 353
Fax: +49 (0) 6151 703 302

India

Agilent Technologies
Vacuum Product Division
205-A, "A" wing of Galleria,
2nd floor, Hiranandani Gardens,
Powai, Mumbai-400 076, India
Tel.: +91 22-2570 8595 / 8597
Fax: +91 22- 2570 8599

Italy

Agilent Technologies
Vacuum Products Division
via F.lli Varian 54
10040 Leini, (Torino) ITALY
Tel.: +39 011 997 9111
Toll-Free: 00 800 234 234 00
Fax: +39 011 997 9350

Japan

Agilent Technologies
Vacuum Products Division
Sumitomo Shibaura Building 4-16-36
8th Floor
4-16-36 Shibaura Minato-ku
Tokyo 108 JAPAN
Tel.: +81 3 5232 1253
Toll-Free: 0120 655 040
Fax: +81 3 5232 1710

Korea

Agilent Technologies
Vacuum Products Division
Shinsa 2nd Bldg. 2F 966-5
Daechi-dong
Kangnam-gu, Seoul
KOREA 135-280
Tel.: +82 2 3452 2452
Toll-Free: 080 222 2452
Fax: +82 2 3452 2451

Mexico

Agilent Technologies
Vacuum Products Division
Concepcion Beistegui No 109
Col Del Valle C.P. 03100
MEXICO, D.F.
Tel.: +52 5 523 9465
Fax: +52 5 523 9472

Southeast Asia

Agilent Technologies
Vacuum Products Division
South East Asia (SEA) - Alex Ho
H/P: +601 2213 1253
Fax: +603 6733 8121

Singapore

Agilent Technologies
Vacuum Products Division
Singapore
Unit 10-04 Helios
Biopolis @ one-north
11 Biopolis Way, 138667
Singapore
H/P: +65 92364988
Fax: +65 64789603

Taiwan

Agilent Technologies
Vacuum Products Division
14F-6, No. 77, Hsin Tai Wu Road,
Sec. 1
Hsi chih, Taipei Hsien, Taiwan,
R.O.C.
Tel.: +886 2 2698 9555
Toll Free: 0800 051 342
Fax: +886 2 2698 9678

UK & Ireland

Agilent Technologies
Vacuum Products Division
6 Mead Road
Oxford Industrial Park
Tel.: +44 (0) 1865 291570
Fax: +44 (0) 1865 291571



Agilent Technologies

