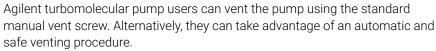


# How to Select Vent Valves for Agilent Turbo Pumps

#### **Abstract**

This technical overview is a guide to help users when selecting the proper Agilent vent valve according to the controller in use. In addition, it provides useful information about the compatibility between existing vent valve models and controller types.





This provides: (a) automatic handling of the pump's ramp down time, avoiding inrush (from wrong and excessive manual venting), and (b) protection against gas back streaming from the fore line (where impurities and hydrocarbon residuals can be present).

Automatic vent valves are normally used as accessories for safe and proper venting of the Agilent turbopumps, to maximize the pump's reliability and improve user experience.

Vent valves can be classified in different ways:

- External or Integrated
- With normally open (NO) or normally closed (NC) logic
- With 0.5 mm or 1.2 mm orifice

Different valve models are available in the Agilent portfolio, according to the turbopump controller in use. Extension cables to connect the vent valve to a specific controller are also available.

This technical overview presents the available choices, starting with the valve type, and details compatibility with controllers.

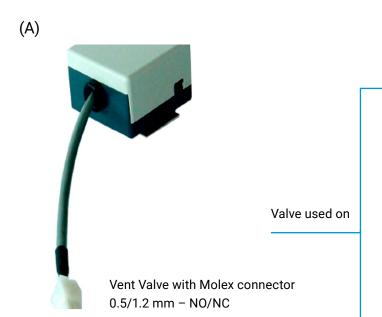


Agilent TwisTorr Turbo Pumps

### **Vent Valves for Onboard Controllers**

The appropriate vent valve type to connect to the different Agilent TwisTorr onboard controllers is displayed in Figure 1. A list of vent valves and their compatibility with the TwisTorr onboard controllers is outlined in Table 1.

Note: extension cables are not available to connect these vent valves and controllers. Instead, an Agilent TwisTorr 305-IC Fan/Vent Adapter kit is required (see Figure 2).





TwisTorr 305 FS onboard controller with TwisTorr 305-IC Fan/Vent Adapter kit\*



TwisTorr Medium-TMP onboard controller for TwisTorr 404 FS, 704 FS, and 804 FS turbo pumps



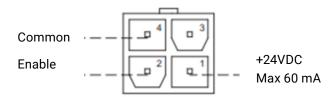


Figure 1. (A) Vent valve for Agilent TwisTorr onboard controllers. (B) A dedicated 24-VDC connector to control the optional vent valve.

# Table 1. Compatibility of Agilent vent valves with Agilent TwisTorr onboard controllers

Vent Valve Description Connector T			TwisTorr 305-IC	TwisTorr 305 FS Onboard Controller	TwisTorr Medium-TMP Onboard Controller (for 404 FS, 704 FS, 804 FS)
	Description	Connector Type	Fan/Vent Adapter Kit* Required	Part Number Part Number	Part Number
			X3507-64130	X3512-64006	
			X3507-64131	X3512-64016	
9699834	N.O. orifice 1.2 mm	Molex	Yes	Yes	Yes
9699834M006	N.O. orifice 0.5 mm	Molex	Yes	Yes	Yes
9699834M010	N.C. orifice 1.2 mm	Molex	Yes	Yes	Yes
9699834M011	N.C. orifice 0.5 mm	Molex	Yes	Yes	Yes

Note: Extension cable not available

<sup>\*</sup>Part number X3514-68001 (Figure 2).



Figure 2. Agilent TwisTorr 305-IC Fan/Vent Adapter kit required for vent valves.

### **Vent Valves for TwisTorr AG Rack Controllers**

The appropriate vent valve type to connect to the different Agilent TwisTorr AG rack controllers is displayed in Figure 3.

A list of vent valves and their compatibility with the TwisTorr AG controllers is outlined in Table 2. A list of extension cables available is outlined in Table 3.

(A)



Valve used on

Vent Valve with USB connector 0.5/1.2 mm - NO/NC

(B)

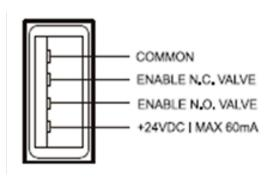


Figure 3.

(A) Vent valve for Agilent TwisTorr active gauge (AG) rack controllers.

(B) A dedicated 24-VDC connector to control the optional vent valve (see orderable parts). The vent valve can be driven automatically by controller, or by serial line.



TwisTorr 74 FS, 84 FS AG rack controllers



TwisTorr 74 FS, 84 FS, and 304 FS AG rack controllers and Agilent TwisTorr 305 remote controller

# Table 2. Compatibility of Agilent vent valves with Agilent AG rack controllers

		TwisTorr 74 FS AG Rack Controller	TwisTorr 84 FS AG Rack Controller	TwisTorr 304 FS AG Rack Controller	TwisTorr 305 Remote Controller
Vent Valve	Description	Part Number	Part Number	Part Number	Part Number
Part Number	Description	X3508-64301	X3508-64001	X3506-64002	X3506-64130
		X3508-64022	X3508-64002	X3506-64003	X3506-64131
9699844	N.O. orifice 0.5 mm	Yes	Yes	Yes	Yes
9699844M001	N.C. orifice 0.5 mm	Yes	Yes	Yes	Yes

Nomenclature: starting with the TwisTorr 305, the rack-mountable controllers are referred to as "remote". Connector type: USB

# Table 3. Extension cables available

Part Number	Length (m)
9699941	5
9699941M003	10
9699941M001	15
9699941M005	20
9699941M007	30
9699941M004	50

# Vent Valves for TwisTorr 74 FS and 84 FS Onboard AG Controllers

The valves mentioned in Table 2 are also compatible with the Agilent TwisTorr 74 FS and 84 FS onboard controllers, as shown in Table 4. Extension cables available for these connections are displayed in Table 5.

Table 4. Compatibility of Agilent vent valves with Agilent TwisTorr 74 FS and 84 FS onboard controllers

		TwisTorr 74 FS Onboard Controller	TwisTorr 84 FS Onboard Controller
Vent Valve Part Number	Description	Part Number	Part Number
		X3508-64030	X3509-64000
		X3508-64021	X3509-64001
9699844	N.O. orifice 0.5 mm	Yes	Yes
9699844M001	N.C. orifice 0.5 mm	Yes	Yes
Connector type: USB			

### Table 5. Extension cables available

Part Number	Length (m)
9699941	5
9699941M003	10
9699941M001	15
9699941M005	20
9699941M007	30
9699941M004	50

# **Vent Valves for Medium Rack Controllers**

The appropriate vent valve type to connect to the different Agilent TwisTorr Medium-TMP rack controllers is displayed in Figure 4.

A list of vent valves and their compatibility with the TwisTorr Medium-TMP rack controllers is outlined in Table 6. A list of extension cables available for these connections is outlined in Table 7.



Valve used on



TwisTorr Medium-TMP rack controller for TwisTorr 404 FS, 704 FS, and 804 FS turbo pumps.

Vent Valve Circular connector 1.2 mm ID - N.O/N.C

Figure 4. Vent valve for Agilent TwisTorr Medium-TMP rack controllers.

# **Vent Valves for TwisTorr Medium-TMP Rack Controllers**

# Table 6. Compatibility of Agilent vent valves with Agilent TwisTorr Medium-TMP rack controllers

	Description	Voltage	TwisTorr Medium-TMP Rack Controller (for 404 FS, 704 FS, 804 FS)	Agilent Turbo-V1K-G Rack Controller	Agilent Turbo-V2300 Rack Controller
Vent Valve Part Number			Part Number	Part Number	Part Number
			X3501-64006	X3501-64005	X3501-64004
			X3501-64016		
X3501-68002	N.O. orifice 1.2 mm	24 V	Yes	-	-
X3501-68022	N.C. Orifice 1.2 mm	24 V	Yes	-	-
Note: Ventualis has a Fire languable					

Note: Vent valve has a 5 m long cable.

# Table 7. Vent valve rack extension cables available

Part Number	Length (m)
X3501-68054	10
X3501-68064	15
X3501-68074	20
X3501-68084	30
X3501-68094	75

# Turbo-V 2K-G and 3K-G Pumps with Integrated Controller

Agilent Turbo-V 2K-G and 3K-G pump and controllers are also available (Table 8). Features of these systems include:

- The controller is integrated in the pump body.
- The purge-vent valves are integrated, factory installed, and fully controlled through the electronics (serial I/O-Profibus if available).
- Complete setting is possible for both vent and purge:
  - a. Automatic: completely operated by the microprocessor according to pump status
  - b. On command: delay time/opening delay/opening time fully configurable

Full list of commands is available in the user manual.

# Table 8. Agilent Turbo-V 2K-G and 3K-G pump and controller part numbers

Turbo-V 2K-G Pump and Controller Part Number	Turbo-V 3K-G Pump and Controller Part Number
9698871	9698888
9698873	9698889

### **Conclusions:**

Agilent can provide an automatic and safe solution for the proper venting of Agilent turbopumps.

An array of Agilent vent valves is available according to the turbopump controller in use, as well as extension cables to connect the vent valves to the controller.

These accessories can help maximize a pump's reliability and improve user experience in order to optimize your laboratory's workflow.

### **Contact information**

#### **Americas**

Agilent Technologies 121 Hartwell Avenue, Lexington, MA 02421 USA Toll free: +1 800 882 7426 vpl-customercare@agilent.com

#### Europe, Middle East, Africa, India

Agilent Technologies Italia SpA Via F.Ili Varian 54, 10040 Leini (Torino), Italy Tel: +39 011 9979 111 Toll free: 00 800 234 234 00

vpt-customercare@agilent.com

#### China

Beijing Office Agilent Technologies (China) Co. Ltd. No.3, Wang Jing Bei Lu, Chao Yang District, Beijing, 100102, China Toll free: 800 820 6778 Contacts.vacuum@agilent.com

For more information, please contact your Agilent representative or visit www.agilent.com/chem/vacuum where you can chat live with a vacuum expert.



### www.agilent.com/chem/vacuum

DE92780894

This information is subject to change without notice.

