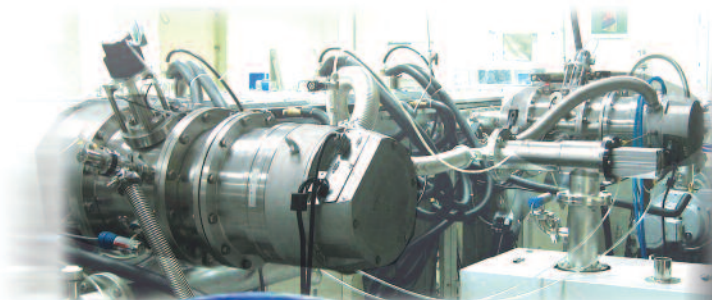


Varian, Inc.
Vacuum Technologies



Vacuum Tools for Coating System Productivity

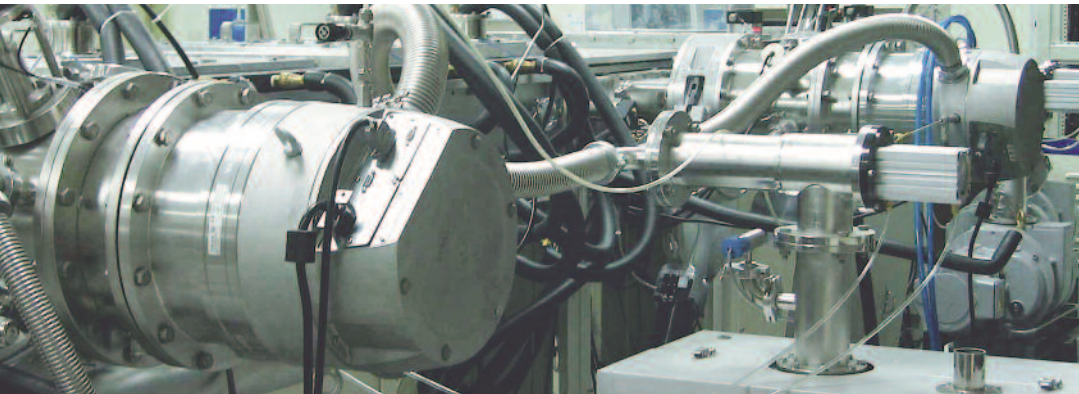


Inspiring Excellence

NOTICE: This document contains references to Varian. Please note that Varian, Inc. is now part of Agilent Technologies. For more information, go to www.agilent.com/chem.

 Agilent Technologies

Vacuum Tools for Coating System Productivity



We at Varian understand the demands of large area vacuum coating systems. We provide the vacuum equipment you need for reliable, energy efficient operations. Our innovative solutions include:

- **Turbo-V K-G Pump Systems** – high-performance, highly-reliable, and designed specifically for large area coating applications
- **MoniTorr/Remote Assist Software** – connectivity to Varian Support Engineers to monitor *and* keep you apprised of the operation of your Turbo-V K-G pumps – no matter where you are in the world
- **VS Series Leak Detectors** – intelligent and easy to use, and if you need to leak check over a distance, the portability of wireless remote control to help you keep your system up and running and leak free

Turbo-V K-G Series Pump Systems

The only truly integrated high vacuum pump system available, the Turbo-V K-G Series high performance turbo system delivers enhanced performance and efficiency for high throughput applications and features a robust mechanical design suitable for heavy duty industrial environments.

- Highest pumping speed in its class: 2200 l/s for Nitrogen
- Throughput: over 1450 sccm for Nitrogen
- Ramp up time < 6 mins
- Maintenance-free with no oil refill
- Unlimited cycle life
- MTF > 200,000 hrs
- On-board power supply
- Diagnostics and electronic interface
- Optimized power consumption and management system with active stop
- Very compact, high performance rotor design for PVD applications
- Profibus®, serial interface or remote I/O provide communication between the pump and system controls enabling operation via a PC using T-Plus Software.



Turbo-V 3 K-G with MoniTorr Controller

Turbo-V 2 K-G with MoniTorr Controller

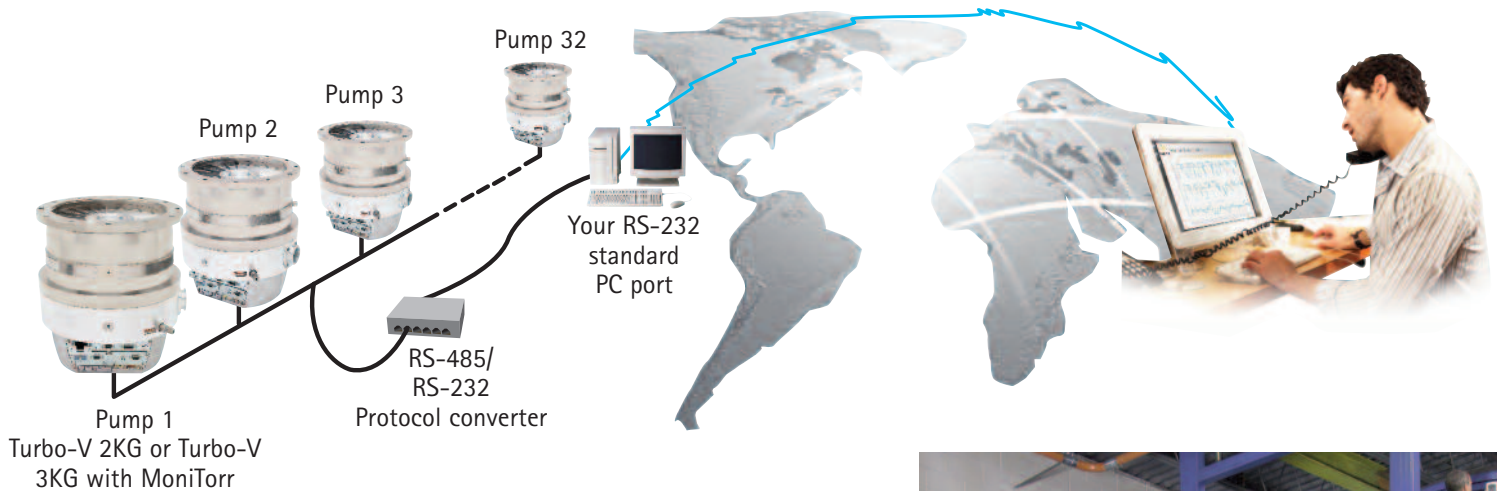


MoniTorr with Remote Assist

MoniTorr is a unique, remote diagnostic system that makes it possible for our technicians to monitor your vacuum system from our location so that you don't need to. MoniTorr is integrated into the pump's controller and continuously monitors all its key functions.

In conjunction with MoniTorr, we offer Remote Assist which is dedicated priority technical support and extended warranty featuring Varian's Advanced Exchange Program, the most comprehensive service offering for turbo pumps in the vacuum industry.

- Integrated MoniTorr card (optional) allows 24/7 monitoring of all pump parameters
- MoniTorr Software Package collects all data
- Allows remote virtual presence of Varian Support Engineers at the user site



VS Series Leak Detectors

VS Series Leak Detectors maximize your productivity and system up-time with the simplicity of two-button operation combined with advanced system intelligence. Robust design allows the VS Series to conform to rigorous industrial standards and operate dependably in the most challenging environments.

VS Wireless Remote Control/Display

Leak testing hard to reach locations is now possible up to 100 meters away. Applications in large systems in which two technicians were previously required now are performed by one, thereby reducing labor costs.

Because the wireless remote has a display and all the key buttons necessary to leak check, all major functions of the leak detector are at your fingertips – even when you are a distance from the instrument. The VS Series Wireless Remote provides truly portable leak detection solution.



VS Series leak detector and Wireless Remote in use at Carbon Composites, Inc.

Varian, Inc. Vacuum Technologies

Benelux

Varian Vacuum Technologies
Herculesweg 8, 4338 PL Middelburg
PO Box 8033, 4330 EA Middelburg
The Netherlands
Tel: (31) 118 671570
Fax: (31) 118 671569

Canada

Central coordination through:
Varian, Inc.
121 Hartwell Avenue
Lexington, MA 02421
USA
Tel: (781) 861 7200
Fax: (781) 860 5437
Toll Free # 1 (800) 882 7426

China

Varian Technologies China, Ltd.
Room 1648, Central Tower South Wing
Beijing Junefield Plaza
No. 10 XuanWuMenWai Street
Beijing 100052
P.R. China
Tel: (86) 10 63108550
Fax: (86) 10 63100141
Toll Free: 800 820 6556

Varian Technologies China, Ltd.
Unit 1801, Central Plaza
No. 227 North Huanpi Road
Shanghai 200003
P. R. China
Tel.: (86) 21 3318-4918
Fax: (86) 21 6375-6848
Toll-Free: 800 820 6556

Varian Technologies China, Ltd.
Room 2508-11 West-Tower
YangCheng
Commercial Building TiYuDong Road
Guangzhou, 510620
P. R. China
Tel.: (86) 20 38871 469
Fax: (86) 20 38871 470
Toll-Free: 800 820 8266

France

Varian s.a.
7 avenue des Tropiques
Z.A. de Courtaboeuf - B.P. 12
91941 Les Ulis cedex
France
Tel: (33) 1 69 86 38 84
Fax: (33) 1 69 86 29 88

Germany and Austria

Varian Deutschland GmbH
Alsfelder Strasse 6
Postfach 11 14 35
64289 Darmstadt
Germany
Tel: (49) 6151 703 353
Fax: (49) 6151 703 302

India

Varian India Pvt. Ltd.
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

Varian, Inc.
via F.lli Varian 54
10040 Leini, (Torino)
Italy
Tel: (39) 011 997 9 111
Fax: (39) 011 997 9 350

Japan

Varian Technologies Japan, Ltd.
Sumitomo Shibaura Building, 8th Floor
4-16-36 Shibaura Minato-ku
Tokyo 108
Japan
Tel: (81) 3 5232 1253
Fax: (81) 3 5232 1710
Toll Free: 0120 655 040

Varian Technologies Japan, Ltd.
Osaka Rep. Office, Shin-Osaka Doi Bldg.
7-5-25 Nishinakayima Yodogawa- ku
532-0011 Osaka
Japan
Tel.: (81) 6 6305 6607
Fax: (81) 6 6305 6556

Korea

Varian Technologies Korea, Ltd.
Shinsa 2nd Bldg. 2F
966-5 Daechi-dong
Kangnam-gu, Seoul
Korea 135-280
Tel: (82) 2 3452 2452
Fax: (82) 2 3452 2451
Toll Free: 080 222 2452

Mexico

Varian, S. de R.L. de C.V.
Concepcion Beistegui No 109
Col Del Valle
C.P. 03100
Mexico, D.F.
Tel: (52) 5 523 9465
Fax: (52) 5 523 9472

South East Asia

Varian Vacuum Technologies
South East Asia (SEA)
Alex Ho
H/P.: (601) 2213 1253
Fax: (603) 6733 8121
alex.ho@varianinc.com

Singapore

Varian Vacuum Technologies Singapore
Unit 10-04 Helios
Biopolis @ one-north
11 Biopolis Way
138667 Singapore
H/P.: (65) 92364988
Fax: (65) 64789603

Taiwan

Varian Technologies Asia Ltd.
14F-6, No.77, Hsin Tai Wu Rd., Sec. 1
Hsi chih, Taipei Hsien
Taiwan, R.O.C.
Tel: (886) 2 2698 9555
Fax: (886) 2 2698 9678
Toll Free: 0800 051342

Varian Technologies Asia, Ltd.
3F-11, No. 26, Tai-Yuan St.
Tai-Yuan Li, Hsinchu Hsien
Taiwan, R.O.C.
Tel.: (886) 3 552 6770
Fax: (886) 3 552 6771

UK and Ireland

Varian Ltd.
6 Mead Road
Oxford Industrial Park
Yarnton
Oxford OX5 1QU
England
Tel: (44) 1865 291570
Fax: (44) 1865 291571

United States

Varian, Inc.
121 Hartwell Avenue
Lexington, MA 02421
USA
Tel: (781) 861 7200
Fax: (781) 860 5437

Other Countries

Varian, Inc.
via F.lli Varian 54
10040 Leini, (Torino)
Italy
Tel: (39) 011 997 9 111
Fax: (39) 011 997 9 350

Toll Free Number for United States
1 (800) 882 7426

Toll Free Number for Europe
00 800 234 234 00

