

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.



979- NW25 Throttle Valve

FIELD INSTRUCTION SHEET

Manual No. 699909986
Revision A
Date October 2002

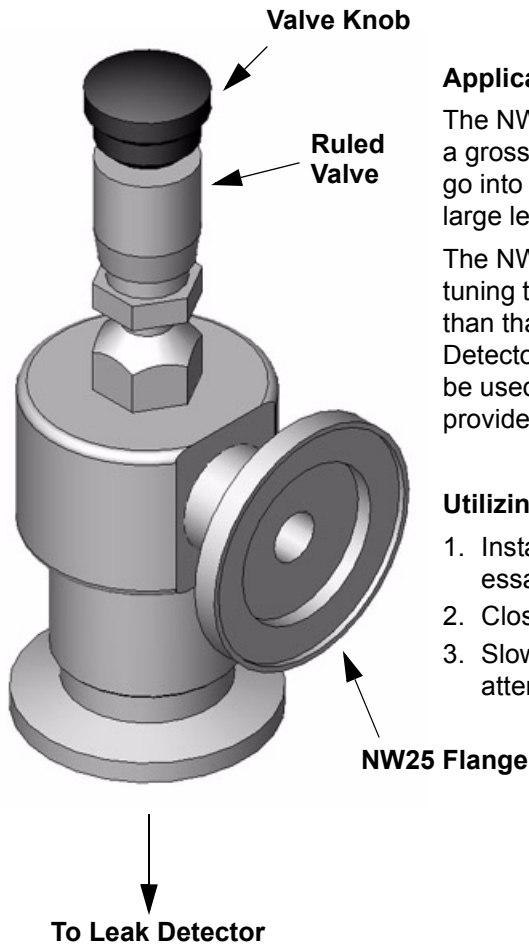
This page intentionally left blank.

Field Instruction Sheet

Varian, Inc

LD 979- NW25 Throttle Valve

The Varian NW25 Throttle Valve combines the functions of a Throttle Valve and a Tuning Leak into one effective tool. It provides an uncalibrated source of helium designed to furnish an ample leak rate signal for tuning Varian Leak Detectors. The helium used by the Tuning Leak is the helium which occurs naturally in the air. (Normally present at a concentration of about 5 parts per million. (5ppm)) It can also be attached to a system via the NW25 Flange. The NW25 Throttle Valve admits a small amount of air into the Leak Detector, which is sensitive to the trace of helium. No separate supply of helium is necessary for use with this tool. Helium flow is controlled by a ruled valve at the top of the NW25 Throttle Valve.



Application

The NW25 Throttle Valve may be used in Leak Detector applications in which a grossly leaking part or vacuum chamber will not permit the Leak Detector to go into the Test Mode. It acts as a non-quantitative approach to locating very large leaks in a vacuum system.

The NW25 Throttle Valve is not a calibrated source of helium. Its value in tuning the leak detector lies in the fact that it provides a signal a decade larger than that from a typical calibrated leak, hence it makes tuning the Leak Detector easier. Likewise, it has a ten-segment ruled needle valve which can be used as an aid to gross leak testing or for gross leak calibration. Or even provide an indication of background helium in the testing room air.

Utilizing the NW25 Throttle Valve

1. Install the NW25 Throttle Valve on the Test Port of the Leak Detector. If necessary, attach any appropriate connectors to the NW25 Flange.
2. Close the Ruled Valve by twisting the Valve Knob to the left.
3. Slowly open the Ruled Valve by twisting the Valve Knob to the right. Paying attention to the response from the Leak Detector.

Technical Assistance

The final page of this document lists all the Varian offices from which you can obtain technical assistance. In the United States, contact Varian Customer Service at 1-800-8Varian or on-line at vpl.customer.support@varianinc.com.

Title

LD979- Throttle Valve

Page 1 of 1

Document No.

699909986

Revision

A

VARIAN



Sales and Service Offices

Canada

Central coordination through:

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

China

Varian Technologies - Beijing

Room 1201, Jinyu Mansion
No. 129A, Xuanwumen Xidajie
Xicheng District
Beijing 1000031
P.R. China
Tel: (86) 10 6608 1031
Fax: (86) 10 6608 1541

France and Benelux

Varian s.a.

7 avenue des Tropiques
Z.A. de Courtaboeuf – B.P. 12
Les Ulis cedex (Orsay) 91941
France
Tel: (33) 1 69 86 38 13
Fax: (33) 1 69 28 23 08

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

101-108, 1st Floor
1010 Competent House
7, Nangal Raya Business Centre
New Delhi 110 046
India
Tel: (91) 11 5548444
Fax: (91) 11 5548445

Italy

Varian, Inc.

Via F.Ili Varian, 54
10040 Leini, (Torino)
Italy
Tel (39) 011 997 9 111
Fax (39) 011 997 9 350

Japan

Varian, Inc.

Sumitomo Shibaura Building, 8th Floor
4-16-36 Shibaura
Minato-ku, Tokyo 108
Japan
Tel: (81) 3 5232 1253
Fax: (81) 3 5232 1263

Korea

Varian Technologies Korea, Ltd.

Shinsa 2nd Building 2F
966-5 Daechi-dong
Kangnam-gu, Seoul
Korea 135-280
Tel: (82) 2 3452 2452
Fax: (82) 2 3452 2451

Mexico

Varian S.A.

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

Russia

Central coordination through:

Varian, Inc.
via F.Ili Varian 54
10040 Leini, (Torino)
Italy
Tel: (39) 011 997 9 252
Fax: (39) 011 997 9 316

Taiwan

Varian Technologies Asia Ltd.

18F-13 No.79, Hsin Tai Wu Road
Sec. 1, Hsi Chih, Taipei Hsien
Taiwan, R.O.C.
Tel: (886) 2 2698 9555
Fax: (886) 2 2698 9678

UK and Ireland

Varian Ltd.

28 Manor Road
Walton-On-Thames
Surrey KT 12 2QF
England
Tel: (44) 1932 89 8000
Fax: (44) 1932 22 8769

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.Ili Varian 54
10040 Leini, (Torino)
Italy
Tel: (39) 011 997 9 111
Fax: (39) 011 997 9 350

Customer Support and Service:

North America

Tel: 1 (800) 882-7426 (toll-free)
vtl.technical.support@varianinc.com

Europe

Tel: 00 (800) 234 234 00 (toll-free)
vtl.technical.support@varianinc.com

Japan

Tel: (81) 3 5232 1253 (dedicated line)
vtj.technical.support@varianinc.com

Korea

Tel (82) 2 3452 2452 (dedicated line)
vtk.technical.support@varianinc.com

Taiwan

Tel: 0 (800) 051 342 (toll-free)
vtw.technical.support@varianinc.com

Worldwide Web Site, Catalog and On-line Orders:

www.varianinc.com

Representatives in most countries



VARIAN