

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

Instructions for Installation of the LD Silencer Kits: LDTS1PSILKIT LDTS2PSILKIT

FIELD INSTRUCTION SHEET

Manual No. 699909989
Revision B
Date April 2003

This page intentionally left blank.

Field Instruction Sheet

Varian Vacuum Technologies

Instructions for LD TS Silencer Kits

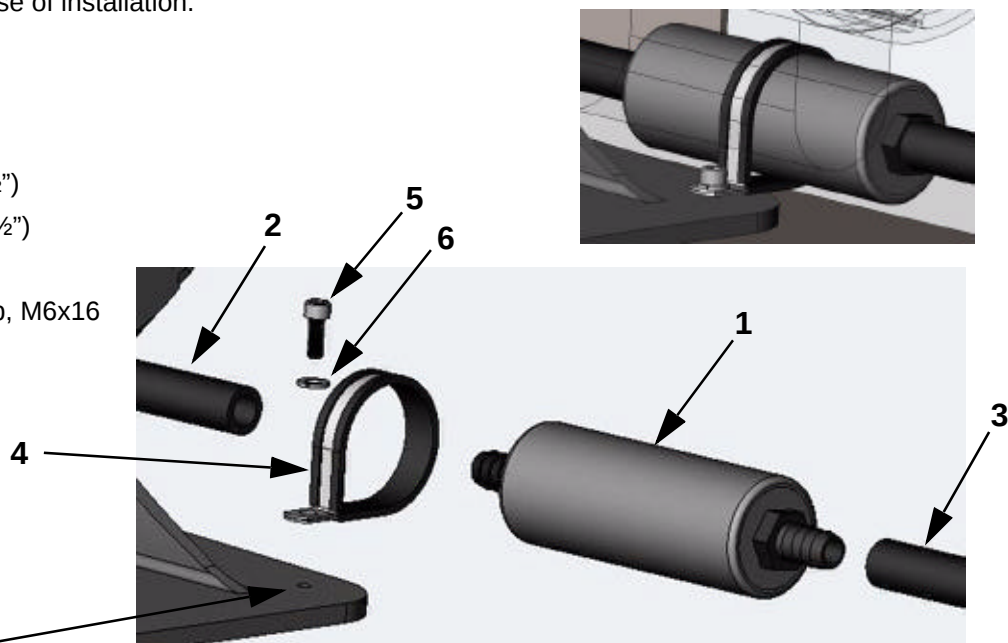
Introduction:

The **LD TS Silencer Kits** (LDTS1PSILKIT and LDTS2PSILKIT) are designed to dampen the ambient noise from the Dry pump during use on both Single- and Dual-Pump Cart units. The installation is universal, but Single-Pump systems shall be used in the description for clarity. **Note:** On Single-Pump, 4-Wheel Cart systems mount the Silencer on the Left side of the Pump for ease of installation.

Components:

- 1 - Silencer Assembly
- 2 - Short Tubing Length (7½")
- 3 - Long Tubing Length (22½")
- 4 - Clamp
- 5 - Screw, Socket Head Cap, M6x16
- 6- Washer, Split Lock, 1/2"

Mounting Hole
Located at the
Right Rear of
the Pump Base



USAGE: (For each Pump as appropriate.)

- 1- Remove Tubing between the Exhaust of Pump and the Exhaust Filter on the Cart and discard.
- 2- Wrap the Clamp (4) around the Body of the Silencer Assembly (1).
- 3- Place the Lock Washer (6) on the Screw (5) and feed the Screw through the eyehole in the Clamp.
- 4- Insert the Screw into the M6 Hole located at the right rear of the Pump Base. Hand tighten.
- 5- Connect the Short Tubing Length (2) between the Exhaust of Pump and the Silencer Assembly.
- 6- Connect the Long Tubing Length (3) between the Silencer Assembly and the Exhaust Filter. Trim if necessary to desired length.

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 Vacuum Technologies Customer Service at 1-800-8Varian or on-line at vpl.customer.support@varianinc.com.

Title

Instructions for Installation of an LD Silencer Kit

Page 1 of 1

Document No.

699909989

Revision

B

VARIAN



Sales and Service Offices

Argentina

Varian Argentina Ltd.

Sucursal Argentina
Av. Ricardo Balbin 2316
1428 Buenos Aires
Argentina
Tel: (54) 1 783 5306
Fax: (54) 1 786 5172

Australia

Varian Australia Pty Ltd.

679-701 Springvale Road
Mulgrave, Victoria ZZ 3170
Australia
Tel: (61) 395607133
Fax: (61) 395607950

Benelux

Varian Vacuum Technologies

Rijksstraatweg 269 H,
3956 CP Leersum
The Netherlands
Tel: (31) 343 469910
Fax: (31) 343 469961

Brazil

Varian Industria e Comercio Ltda.

Avenida Dr. Cardoso de Mello 1644
Vila Olimpia
Sao Paulo 04548 005
Brazil
Tel: (55) 11 3845 0444
Fax: (55) 11 3845 9350

Canada

Central coordination through:

Varian Vacuum Technologies
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 6641 1530
Fax: (86) 10 6641 1534

France and Wallonie

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 Vacuum Technologies

Via F.lli Varian, 54
10040 Leini, (Torino)
Italy
Tel: (39) 011 997 9111
Fax: (39) 011 997 9350

Japan

Varian Vacuum Technologies

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 Bldg. 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

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 Vacuum Technologies

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

Other Countries

Varian Vacuum Technologies

Via F.lli Varian, 54
10040 Leini, (Torino)
Italy
Tel: (39) 011 997 9111
Fax: (39) 011 997 9350

Internet Users:

Customer Service & Technical Support:

vpl.customer.support@varianinc.com

Worldwide Web Site:

www.varianinc.com/vacuum

Order On-line:

www.evarian.com

Representatives in most countries



VARIAN