

Agilent Vibration Damper

Product Name	Part Number
VIBRATION DAMPER, ISO 63	9699340

Extremely low vibration amplitudes at the turbopump inlet flange are required for some sensitive applications, such as scanning electron microscopes, or analytical instrumentation. Agilent Turbo-V pumps have very low levels of vibration amplitude, but the vibration damper 9699340 is useful when additional dampening of the vibration amplitude is required. This vibration damper can be only used in conjunction with turbopumps with ISO 63 inlet flange. It typically reduces the main vibration amplitude by a factor of 20. It must be installed between pump and system, using the clamps provided in the kit (figure 1).

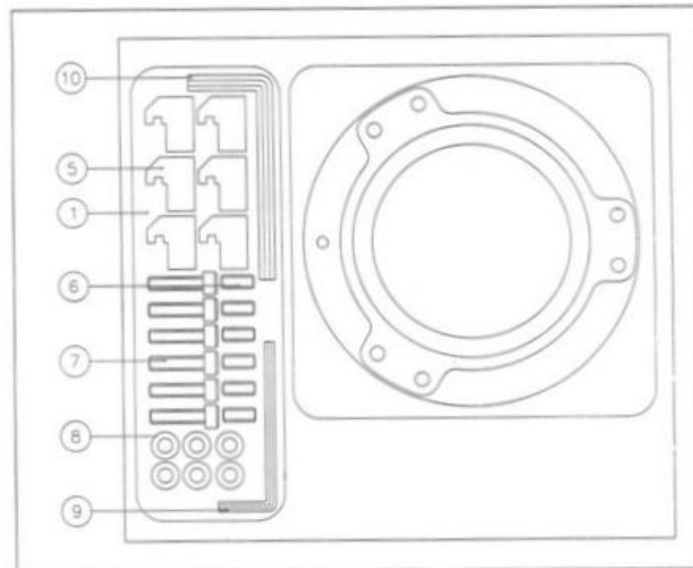


Figure 1 - Vibration damper kit

Instructions for use

Install first all six clamps (5) with the washers (8) and screws (7). Tight each screw using the Allen key (10) in turn and repeat sequential tightening until flange faces meet the damper. Then insert set screw (6) and lock it, using both Allen keys on screw and set screw.

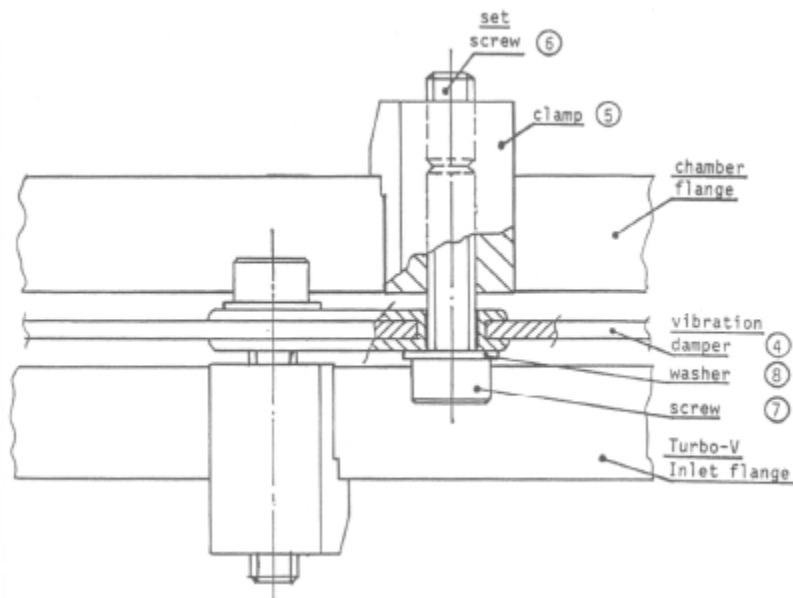


Figure 2 - Vibration damper installation

Instructions for use

Disposal

The following symbol is applied in accordance with the EC WEEE (Waste Electrical and Electronic Equipment) Directive. This symbol (valid only in countries of the European Community) indicates that the product it applies to must NOT be disposed of together with ordinary domestic or industrial waste but must be sent to a differentiated waste collection system. The end user is therefore invited to contact the supplier of the device, whether the Parent Company or a retailer, to initiate the collection and disposal process after checking the contractual terms and conditions of sale.



For more information refer to:

<http://www.agilent.com/environment/product/index.shtml>

Customer Support & Service:

NORTH AMERICA:

Toll Free: 800 882 7426
vp-l-ra@agilent.com
Lexington-service@agilent.com

EUROPE:

Toll Free: 00 800 234 234 00
vpt-customer-care@agilent.com

PACIFIC RIM:

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

Worldwide Web Site, Catalog and Order Online:

www.agilent.com/chem/vacuum
Representatives in most countries

This information is subject to change without notice.

© Agilent Technologies, Inc. 2026

Printed in Italy, 03/2026

87-900-858-01 (B.00)