

Instructions for use

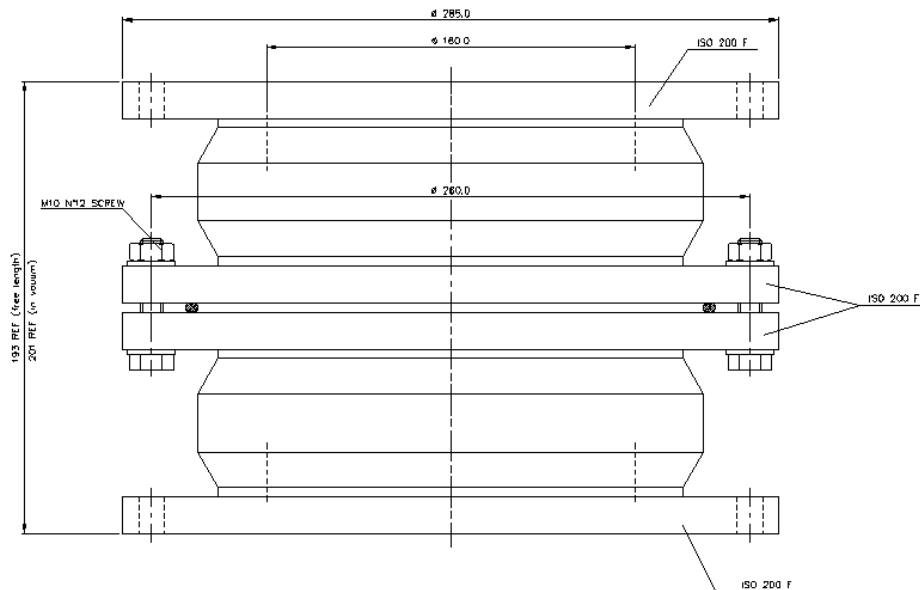
Agilent Double Vibration Damper ISO 200 F flange

Product Name	Part Number
Double Vibration Damper ISO 200 F flange	9699378



Extremely low vibration amplitudes are required for some sensitive applications, even lower than already low levels normally present in the Turbo SEM pumps. Typically scanning electron microscopes or other analytical instruments require additional dampening of vibration amplitude. To further reduce the residual amplitude (of the SEM pumps), a new double vibration damper is provided. To effectively reduce the total vibration level transmitted to the system, the Turbo pump should be suspended from the double damper. Although the actual dampening effect depends on the specific complete system, using a Agilent new double damper, the residual vibration amplitude can be typically reduced by a factor 100 (for ISO 200 F flange).

The damper's specifications are available for the V551 SEM pumps.



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
vpl-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

87-900-961-01 (B.00)