

Field Instructions

ConvecTorr Simulator

Product Name	P/N
ConvecTorr Simulator	L7382301

Description

The ConvecTorr Simulator (L7382301) is intended to be a convenient tool for verifying proper operation of the ConvecTorr control unit. For those cases when the target gauge tube cannot be brought to proper pressure levels for true in-situ calibration, the simulator can be used to at least ensure the control unit is calibrated to the average tube.

Operation

Normally, the simulator would be plugged into the target gauge cable at the sensor end. When the switch is set to the ATM, the controller should read approximately 1×10^5 Pa (760 Torr, 1000 mbar). If desired, the controller may be calibrated by pressing the **<CAL>** button on XGS. Sliding the switch to **VAC** should cause the controller to read approximately 1.33×10^{-1} Pa (1 mT, 1.3×10^{-3} mbar). Again, the controller may be calibrated by pressing the appropriate button on the controller keypad.

NOTE

The Simulator is not intended to take the place on NIST-traceable calibration. The product is intended to verify the functionality of control electronics.

Field Instructions

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

699908311 (B.00)