

## Instructions for use

# Thermocouple Simulator

Product Name	P/N
Thermocouple Simulator	L7383301

### Description

The Simulator (L7383301) is intended to be a convenient tool for verifying proper operation of the TC control unit. For those cases when the target gauge tube cannot be brought to proper pressure levels for true in-situ calibration, use the Simulator to at least ensure the control unit is calibrated to the average tube.

### Operation

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

## 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  
[vp1-ra@agilent.com](mailto:vp1-ra@agilent.com)  
[Lexington-service@agilent.com](mailto:Lexington-service@agilent.com)

##### **EUROPE:**

Toll Free: 00 800 234 234 00  
[vpt-customer care@agilent.com](mailto: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](http://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

699908312 (C.00)