

## TPS-compact Gas Ballast

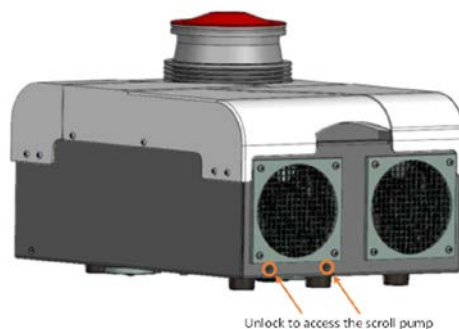
The TPS-compact packaging includes a standard gas ballast kit (P/N IDP3GBKIT) that can be used to prevent water and other condensates from accumulating within the foreline pump. The standard configuration has a seal screw installed in the 1/4-20 thread in the top of the exhaust manifold. This configuration can be used for dry applications.

When appreciable amounts of water, solvents, etc., are being pumped, remove the seal screw and replace it with the provided adapter and 1/8" NPT sintered filter plug.

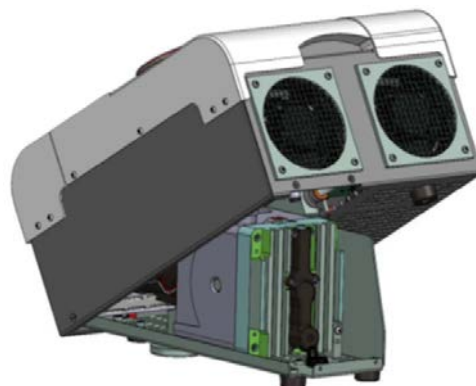
### WARNING

**Before install gas ballast adapter remove the power supply cable.**

1. Disconnect the TPS-compact from electrical power.
2. Unlock the scroll pump support by unscrewing the screws.

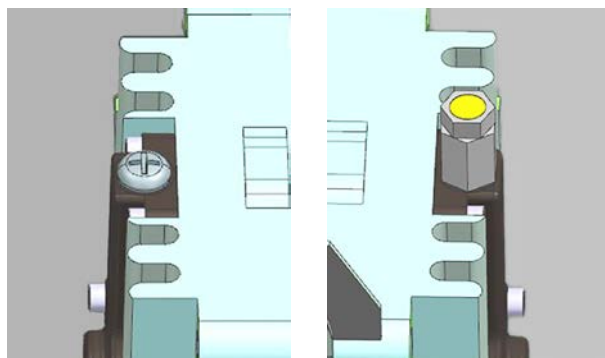


3. Lift the TPS-compact in order to access the Scroll pump site.



## Quick Reference Guide

4. Unscrew seal screw and install provided gas ballast adapter and 1/8" NPT.



5. Lock the scroll pump support by screwing the screws.

### CAUTION

**Be careful during pump support closing to prevent damage to system components (vacuum pipes and connectors).**

### NOTE

1. During the operation of the gas ballast feature the noise may increase.
2. There is not a predefined amount of time to let the gas ballast open: it depends on the amount of moisture trapped in the system.

### CAUTION

**Do not exceed 2.8 Nm of torque when installing the adapter into the manifold.**

For more details about gas ballast refer to the manuals of the single TPS-compact components.

### CAUTION

**Use the provided adapter and 1/8" NPT sintered filter plug in order to avoid dust entrance.**

### CAUTION

**The gas ballast must be sealed whenever pumping any gas not intended to be vented to atmosphere.**

#### Contacts:

U.S. and Canada (toll free):  
+1 800 882 7426  
[vpl-customer care@agilent.com](mailto:vpl-customer care@agilent.com)

Europe & other countries (toll free):  
+00 800 234 234 00  
[vpl-customer care@agilent.com](mailto:vpl-customer care@agilent.com)

China (toll free):  
400 8206778 (mobile)  
800 8206778 (landline)  
[vpd-support-china@agilent.com](mailto:vpd-support-china@agilent.com)

Visit the product web page: [www.agilent.com/chem/tpscompact](http://www.agilent.com/chem/tpscompact)

This information is subject to change without notice.  
© Agilent Technologies, Inc. 2022  
Printed in Italy, January 14th, 2022  
X3582-90000-9-PUI rev. A.00