

## IDP-15 Quick Reference Card

### Start Up/Shutdown

**Warning** Read all Warnings, Cautions and Notes in the IDP-15 Dry Scroll Pump Instruction manual (X3815-90000) before pump operation. Obey all pump labeling.

**Warning** The pump is designed for operation with neutral or noncorrosive gases. It is absolutely forbidden to use it with potentially explosive or inflammable substances.

### Start Up

#### Pre-Installation Minimums

Do not install or use the pump in an environment exposed to atmospheric agents (rain, snow, ice), dust, aggressive gases, or in explosive environments or those with a high fire risk.

If placing the IDP-15 pump inside an enclosure, provide ample room to supply ambient air to the pump front and rear air intakes.

During operation, the following environmental conditions must be respected:

- Temperature: +5 °C to +40 °C (41 °F to 104 °F)
- Relative humidity: 0 to 95% (non-condensing)

- Connect the pump to the power supply using an IEC-320 style power cord of at least 10 A capacity.
- In order to reach maximum vacuum, the pump must be left running for about an hour with the inlet sealed.

#### Unpacking and Inspection

1. Orient the shipping container with *This End Up* on top.
2. Open the box and carefully lift the IDP-15 and foam blocks out of the box. Remove the foam blocks and the plywood skid.
3. Save the carton and all packing materials.
4. Inspect the pump for damage. If there is shipping damage, contact the freight carrier and your local Agilent sales office immediately.
5. Loosen the lower two front and rear cowlings screws and slide the L-brackets up so that the pump rests on its rubber feet.
6. Disengage the Shipping Locks (x 4). Beginning with the rear side, disengage the shipping locks (x 2) by inserting the 4 mm Allen Wrench provided into the holes. Turn the screw clockwise until significant resistance is met. Repeat for the front (fan side) shipping locks (x 2). When the locks are disengaged, the pump body will float on its isolation mounts.

#### Setting Mains Voltage

For IDP-15 pumps with fixed speed motors (part numbers X3815-64000 and X3815-64010) only:

- Slide the voltage selector switch to 115 V for 115-120 V mains or to 230 V for 200-230 V mains

#### Startup

1. Check that the inlet screen is installed before beginning operation.

**Warning** Do not insert a finger or any foreign object in the path of the fan; serious personal injury may result or the pump may be damaged.

2. Operate the pump within ambient temperature constraints, otherwise damage to the pump or shortened operating life may result.

#### Caution Do not block the fan ducts.

Blocking these ducts can cause pump overheating. A pump surface temperature in excess of 65 °C (150 °F) is potentially damaging. If such conditions are observed, turn the pump off and allow it to cool. Disassemble, inspect for damage, and repair, if necessary.

## IDP-15 Quick Reference Card

### Start Up/Shutdown

**Warning** Read all Warnings, Cautions and Notes in the IDP-15 Dry Scroll Pump Instruction manual (X3815-90000) before pump operation. Obey all pump labeling.

**Warning** The pump is designed for operation with neutral or noncorrosive gases. It is absolutely forbidden to use it with potentially explosive or inflammable substances.

### Start Up

#### Pre-Installation Minimums

Do not install or use the pump in an environment exposed to atmospheric agents (rain, snow, ice), dust, aggressive gases, or in explosive environments or those with a high fire risk.

If placing the IDP-15 pump inside an enclosure, provide ample room to supply ambient air to the pump front and rear air intakes.

During operation, the following environmental conditions must be respected:

- Temperature: +5 °C to +40 °C (41 °F to 104 °F)
- Relative humidity: 0 to 95% (non-condensing)

- Connect the pump to the power supply using an IEC-320 style power cord of at least 10 A capacity.
- In order to reach maximum vacuum, the pump must be left running for about an hour with the inlet sealed.

#### Unpacking and Inspection

1. Orient the shipping container with *This End Up* on top.
2. Open the box and carefully lift the IDP-15 and foam blocks out of the box. Remove the foam blocks and the plywood skid.
3. Save the carton and all packing materials.
4. Inspect the pump for damage. If there is shipping damage, contact the freight carrier and your local Agilent sales office immediately.
5. Loosen the lower two front and rear cowlings screws and slide the L-brackets up so that the pump rests on its rubber feet.
6. Disengage the Shipping Locks (x 4). Beginning with the rear side, disengage the shipping locks (x 2) by inserting the 4 mm Allen Wrench provided into the holes. Turn the screw clockwise until significant resistance is met. Repeat for the front (fan side) shipping locks (x 2). When the locks are disengaged, the pump body will float on its isolation mounts.

#### Setting Mains Voltage

For IDP-15 pumps with fixed speed motors (part numbers X3815-64000 and X3815-64010) only:

- Slide the voltage selector switch to 115 V for 115-120 V mains or to 230 V for 200-230 V mains

#### Startup

1. Check that the inlet screen is installed before beginning operation.

**Warning** Do not insert a finger or any foreign object in the path of the fan; serious personal injury may result or the pump may be damaged.

2. Operate the pump within ambient temperature constraints, otherwise damage to the pump or shortened operating life may result.

#### Caution Do not block the fan ducts.

Blocking these ducts can cause pump overheating. A pump surface temperature in excess of 65 °C (150 °F) is potentially damaging. If such conditions are observed, turn the pump off and allow it to cool. Disassemble, inspect for damage, and repair, if necessary.

### Grounding Instructions

This pump is supplied with a power cord that has a grounding wire with an appropriate grounding plug. The plug must be inserted into an outlet that is properly installed and grounded in accordance with all local codes and ordinances.

**Warning** Improper use or modification of the grounding plug can result in a risk of electrical shock.

- Do not modify the plug provided; if it does not fit the outlet, obtain an IEC320 power cord that properly fits the electrical outlet.
- Connect the product only to an outlet that has the same configuration as the plug.
- Do not use an adapter with this product.

### Extension Cords

Do not use extension cords.

**NOTE** For all configurations, refer to the Run Currents section of the IDP-15 Dry Scroll Pump Instruction manual (699904350).

### Shutdown

To stop the pump:

- Toggle the On/Off switch.

### Mechanical Connections

#### Pump Inlet

- Use NW25, or larger, clean vacuum hardware with as short a length as practical between the pump and vacuum chamber.
- Insert a bellows between the pump and vacuum chamber to provide both vibration isolation and strain relief.

#### Pump Exhaust

- An NW16 exhaust fitting is located at the side of the pump. To avoid overheating the pump, do not restrict the exhaust flow with long lengths of small diameter tubing. Use as short as practical lengths of 16 mm ID, or larger, diameter hardware.

#### Gas Ballast

- When pumping gas loads containing water vapor or condensable gases, use the gas ballast. To activate the gas ballast, remove the solid 1/4 NPT plug from either port 1 or port 2 and install the sintered filter plug provided with the pump.

**Caution** Pumping high water vapor loads can cause a temporary increase in ultimate pressure due to adsorption and absorption of water vapor by the internal surfaces of the pump. Pumping water vapor loads in excess of the water vapor handling capability can cause reduced time between tip seal replacements.

### Tip Seal Replacement

Refer to the IDP-15 Dry Scroll Pump Instruction manual (X3815-90000).

### Troubleshooting

Refer to the IDP-15 Dry Scroll Pump Instruction manual (X3815-90000). Visit the Agilent Literature Library at <http://www.chem.agilent.com/en-US/Search/Library/Pages/default.aspx>. Keyword *IDP-15*, choose Publication Type *User Manual*.



Agilent Technologies

Manual No. X3815-90001 Rev. B  
July 2015

### Grounding Instructions

This pump is supplied with a power cord that has a grounding wire with an appropriate grounding plug. The plug must be inserted into an outlet that is properly installed and grounded in accordance with all local codes and ordinances.

**Warning** Improper use or modification of the grounding plug can result in a risk of electrical shock.

- Do not modify the plug provided; if it does not fit the outlet, obtain an IEC320 power cord that properly fits the electrical outlet.
- Connect the product only to an outlet that has the same configuration as the plug.
- Do not use an adapter with this product.

### Extension Cords

Do not use extension cords.

**NOTE** For all configurations, refer to the Run Currents section of the IDP-15 Dry Scroll Pump Instruction manual (699904350).

### Shutdown

To stop the pump:

- Toggle the On/Off switch.

### Mechanical Connections

#### Pump Inlet

- Use NW25, or larger, clean vacuum hardware with as short a length as practical between the pump and vacuum chamber.
- Insert a bellows between the pump and vacuum chamber to provide both vibration isolation and strain relief.

#### Pump Exhaust

- An NW16 exhaust fitting is located at the side of the pump. To avoid overheating the pump, do not restrict the exhaust flow with long lengths of small diameter tubing. Use as short as practical lengths of 16 mm ID, or larger, diameter hardware.

#### Gas Ballast

When pumping gas loads containing water vapor or condensable gases, use the gas ballast. To activate the gas ballast, remove the solid 1/4 NPT plug from either port 1 or port 2 and install the sintered filter plug provided with the pump.

**Caution** Pumping high water vapor loads can cause a temporary increase in ultimate pressure due to adsorption and absorption of water vapor by the internal surfaces of the pump. Pumping water vapor loads in excess of the gas ballast water vapor handling capability can cause reduced time between tip seal replacements.

### Tip Seal Replacement

Refer to the IDP-15 Dry Scroll Pump Instruction manual (X3815-90000).

### Troubleshooting

Refer to the IDP-15 Dry Scroll Pump Instruction manual (X3815-90000). Visit the Agilent Literature Library at <http://www.chem.agilent.com/en-US/Search/Library/Pages/default.aspx>. Keyword *IDP-15*, choose Publication Type *User Manual*.



Agilent Technologies

Manual No. X3815-90001 Rev. B  
July 2015