

DS Pumps Quick Reference Guide

CAUTION!

Before starting the pump, fill up with lubricating oil as the pump is delivered empty.

WARNING!

The pump is designed for operation with neutral or non-corrosive fluids. It is absolutely forbidden to use potentially explosive or flammable substances.

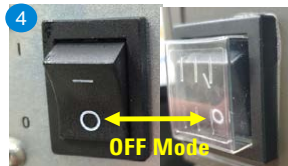
Standard for All Pumps



Step 1: Connect Flange Inlet NW25 to Dust Trap/Filter.

Step 2: Remove Safety Valve & Fill Oil to Max Level.

Step 3: Remove Cap Seal & Connect Flange Outlet NW25 to Exhaust System.



Step 4: Ensure Switch is OFF mode.

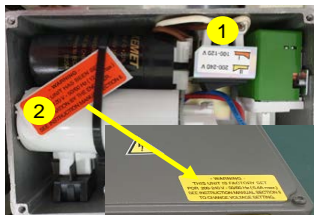
Step 5: Remove Yellow Caution Label.

Step 6: Connect the Power Cord & Lock with Retention Spring.

Step 7: Power On.

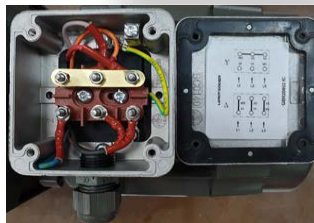


Additional Info for certain Motor Type



Single Phase Motor

1. Motor Default Voltage is 220-240Volt.
2. User may switch to 100-120V & replace Yellow Label with Orange Label on Switch Box Cover.



Three Phase Motor

1. User may connect the cable in "Y" or "Δ".

CAUTION!

Pay special attention to the three phase motor, where an inversion of polarity causes inversion of the direction of rotation of the pump with consequent possibilities of mechanical damage.



Inverter Motor with I/O Male Connector

1. User needs to plug in the I/O Female Connector (provided with Pump).



DS Pumps Maintenance

WARNING!

Disconnect the power supply to the pump before starting maintenance work.
Place a special warning sign over the power supply breaker switch: **MACHINE UNDERGOING MAINTENANCE - DO NOT POWER ON.**
When finished, remove the safety warning. If you have to perform maintenance on the pump after a considerable time in operation, leave it to cool as temperature of the outer surface may be in excess of 60 °C.

Always refer to Product User Manual for more Product Maintenance & other information. <https://www.agilent.com/en/products/vacuum-technologies>

Oil Change Procedure

CAUTION!

Do not change the oil immediately after stopping the machine as the oil temperature may still be high.

MAINTENANCE

It is recommended that the oil change is done periodically every 6 months.

Step 1: Manually remove Safety Valve.

Step 2: Remove Magnetic Plug by using M10 Hex Torque Wrench/Hex Key & drain oil from Pump.

Step 3: Check the DS FKM Gasket Condition. If damaged, replace with the new Gaskets (Note: Additional Gaskets are attached with Pumps).

Step 4: Re-install Magnetic Plug with Gasket by using M10 Hex Torque Wrench/Hex Key with 10Nm.

Step 5: Pour the oil into pump until Oil Max Level & then re-install the Safety Valve.

Step 6: Oil leak check from Magnetic Plug after operate pump for 30min, 1 hour & 2 hours.



Magnetic Plug with DS FKM Gasket



Change the DS FKM Gasket if found damaged

