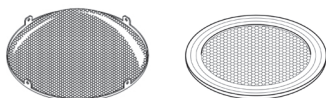


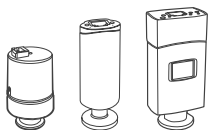
TwisTorr 305 FS Options

Optional Accessories

Inlet screen



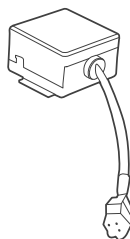
Active Gauges



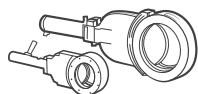
Purge Valve



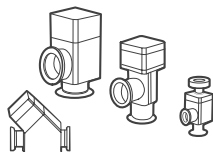
Vent Valve



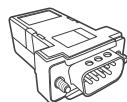
High Vacuum Valves



Aluminium Block Valves



Dongle

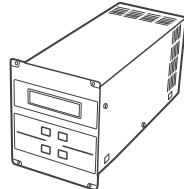


Mobile App



Controller

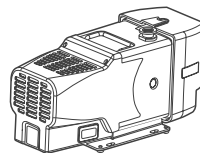
Remote Controller



Suggested Agilent Foreline Pumps

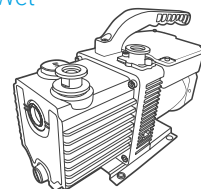
Low Gas Flow

Dry



IDP-3

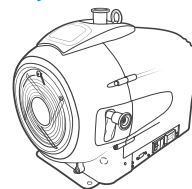
Wet



DS40M

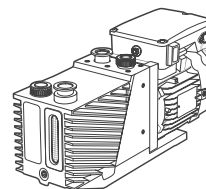
High Gas Flow

Dry



IDP-15

Wet

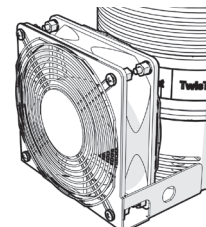


DS302

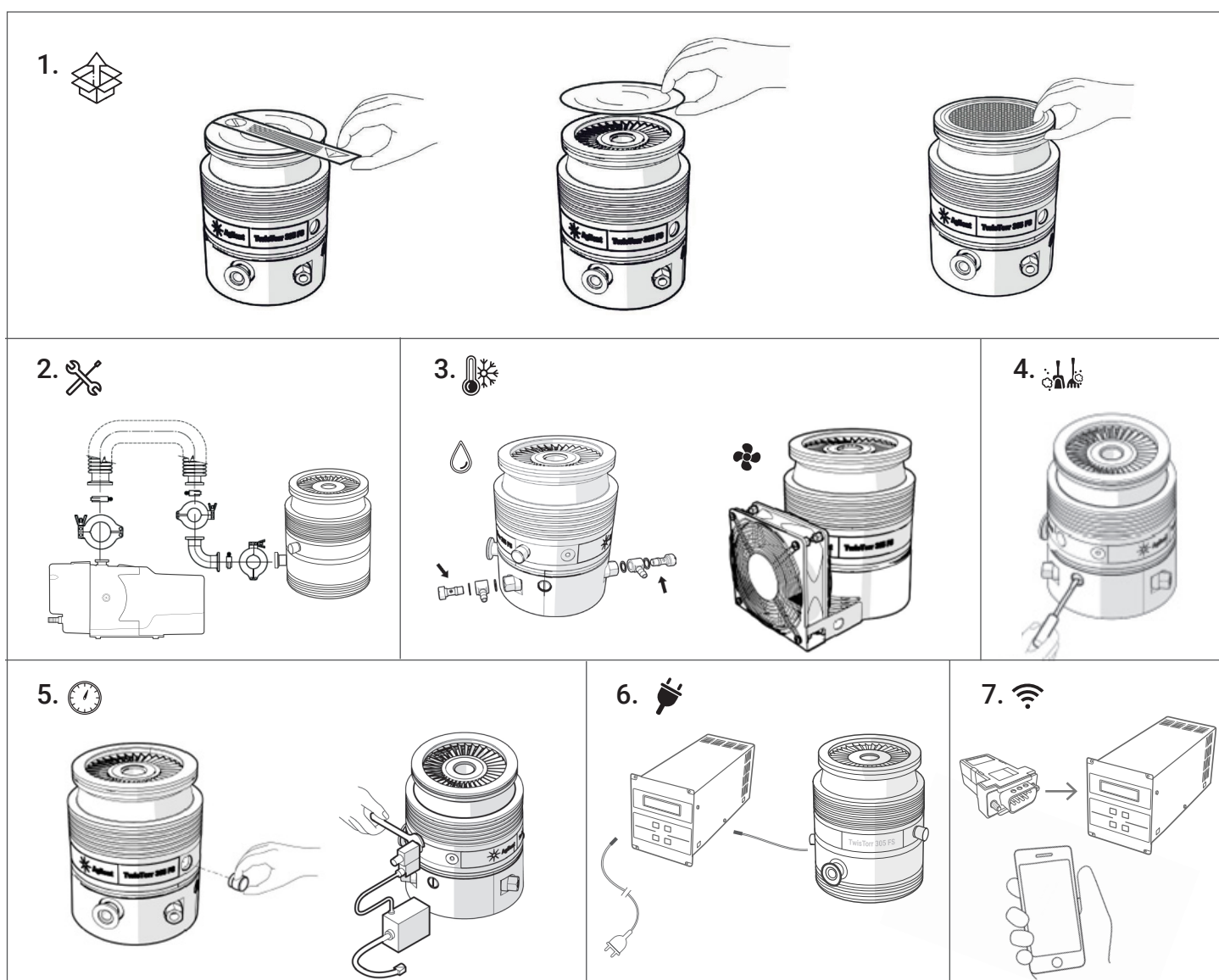
Cooling

Water Cooling

Air Cooling



TwisTorr 305 FS Installation



Visit the product web page:
www.agilent.com/chem/TwisTorr305FS