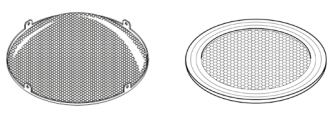


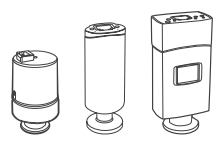
TwisTorr 305 FS, 305 FSQ & 305 SF Options

Optional Accessories

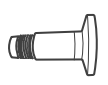
Inlet screen



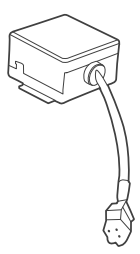
Active Gauges



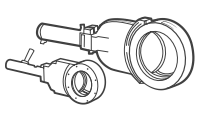
Purge Valve



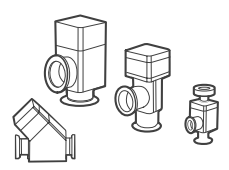
Vent Valve



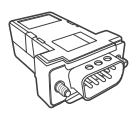
High Vacuum Valves



Aluminium Block Valves



Dongle

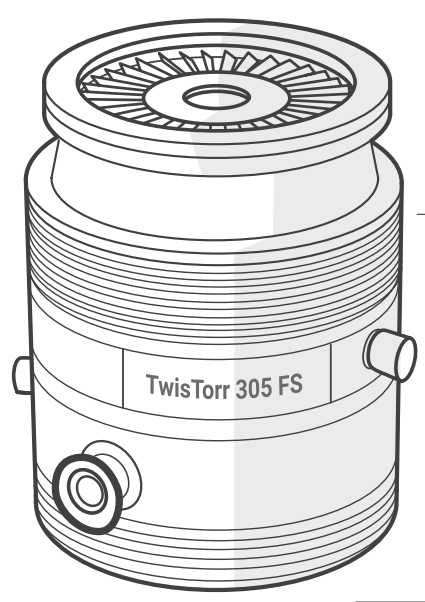
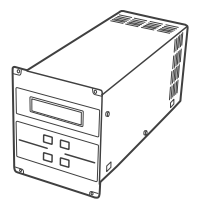


Mobile App



Controller

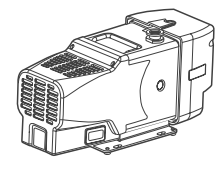
Remote Controller



Suggested Agilent Foreline Pumps

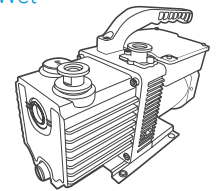
Low Gas Flow

Dry



IDP-3

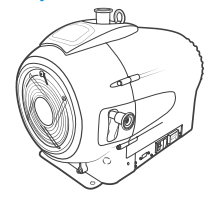
Wet



DS40M

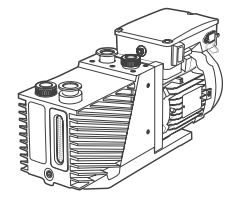
High Gas Flow

Dry



IDP-15

Wet

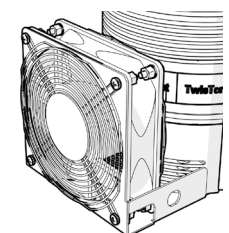


DS302

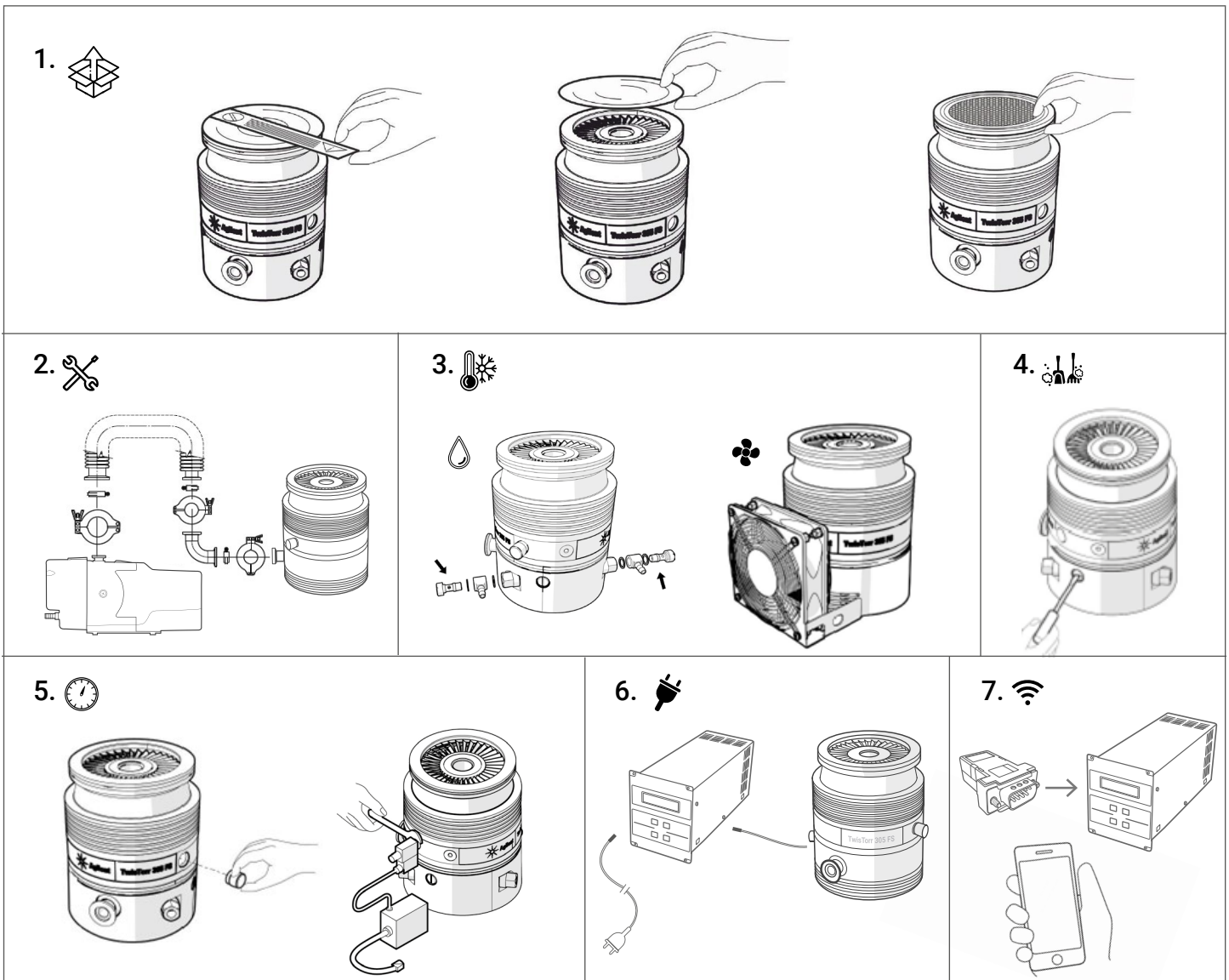
Cooling

Water Cooling

Air Cooling



Twistorr 305 FS, 305 FSQ & 305 SF Installation



Visit the product web page:

www.agilent.com/chem/twistorr305fs
www.agilent.com/chem/twistorr305fsq
www.agilent.com/chem/twistorr305sf

This information is subject to change without notice.

© Agilent Technologies, Inc. 2021
 Printed in Italy, January 14th, 2021
 X3513-90001 rev. B.00