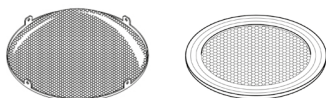


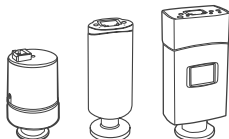
TwisTorr 305-IC, 305-ICQ & 305-IC SF Options

Optional Accessories

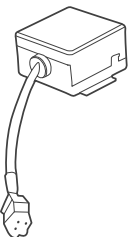
Inlet screen



Active Gauges



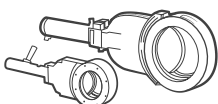
Vent Valve



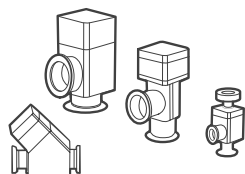
Purge Valve



High Vacuum Valves



Aluminium Block Valves



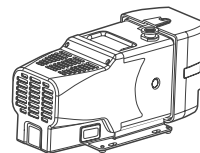
Mobile App



Suggested Agilent Foreline Pumps

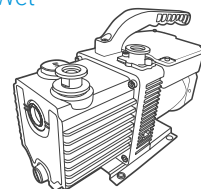
Low Gas Flow

Dry



IDP-3

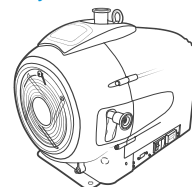
Wet



DS40M

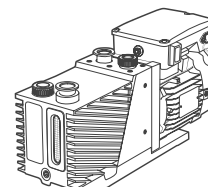
High Gas Flow

Dry



IDP-15

Wet

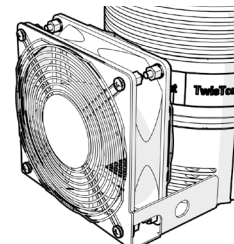
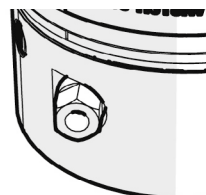


DS302

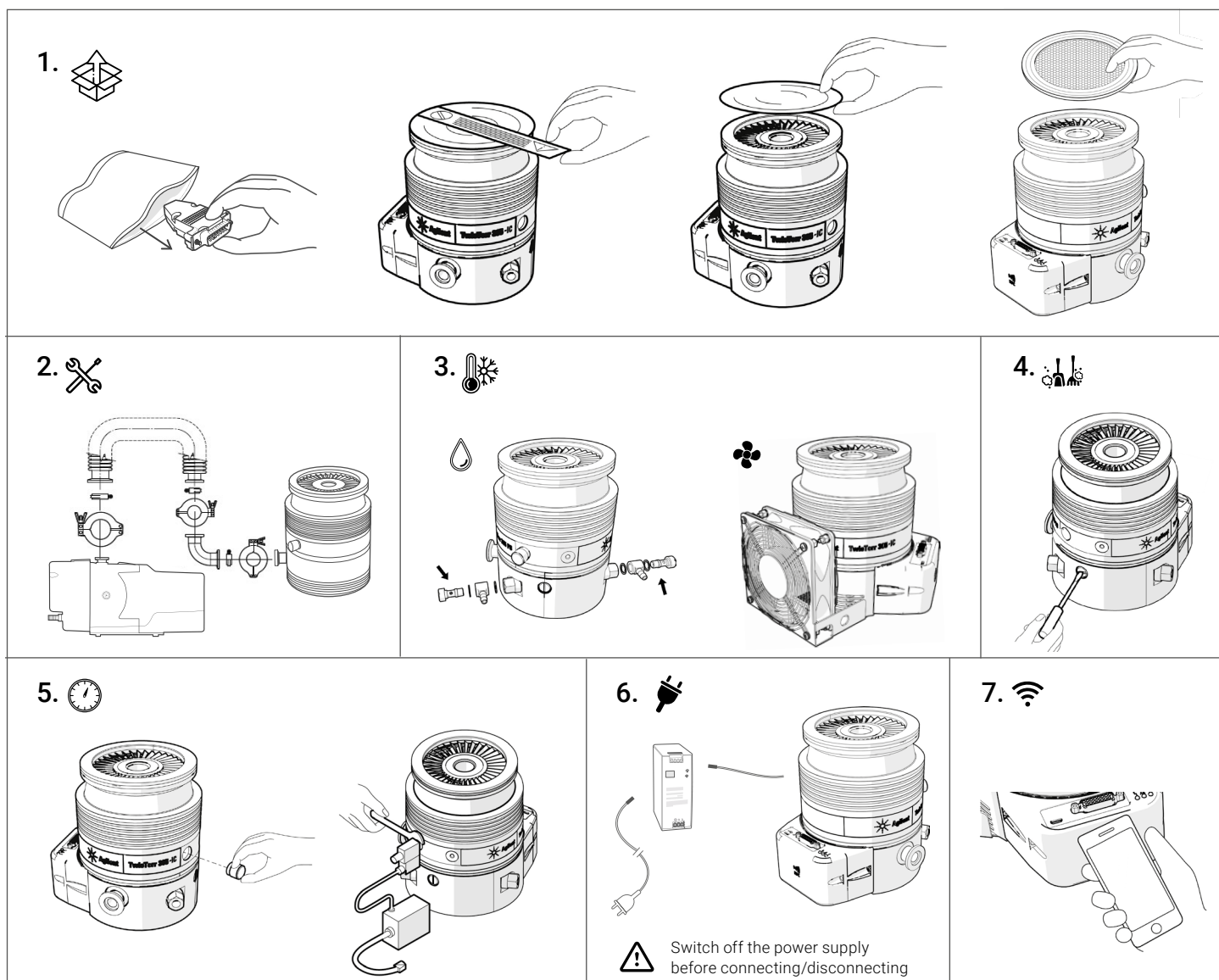
Cooling

 Water Cooling

 Air Cooling



TwisTorr 305-IC, 305-ICQ & 305-IC SF Installation



Visit the product web page:

www.agilent.com/chem/twistorr305-ic

www.agilent.com/chem/twistorr305icq

www.agilent.com/chem/twistorr305icsf

This information is subject to change without notice.

© Agilent Technologies, Inc. 2021

Printed in Italy, January 14th, 2021

X3513-90000 rev. C.00