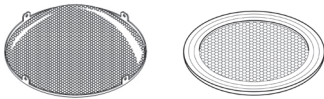


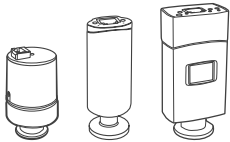
TwisTorr 305-IC Options

Optional Accessories

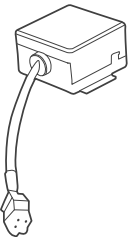
Inlet screen



Active Gauges



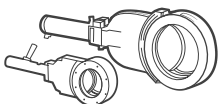
Vent Valve



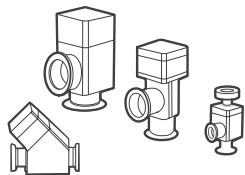
Purge Valve



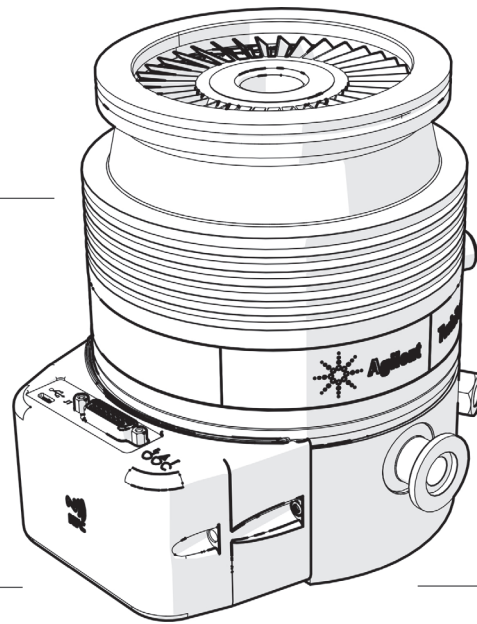
High Vacuum Valves



Aluminium Block Valves



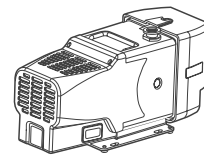
Mobile App



Suggested Agilent Foreline Pumps

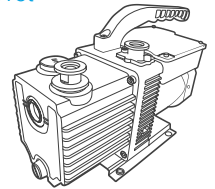
Low Gas Flow

Dry



IDP-3

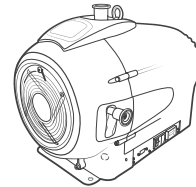
Wet



DS40M

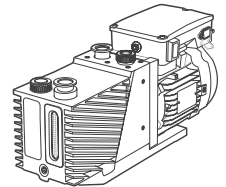
High Gas Flow

Dry



IDP-15

Wet

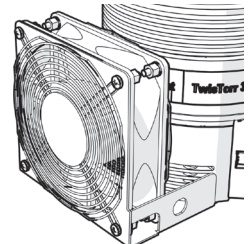
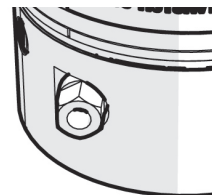


DS302

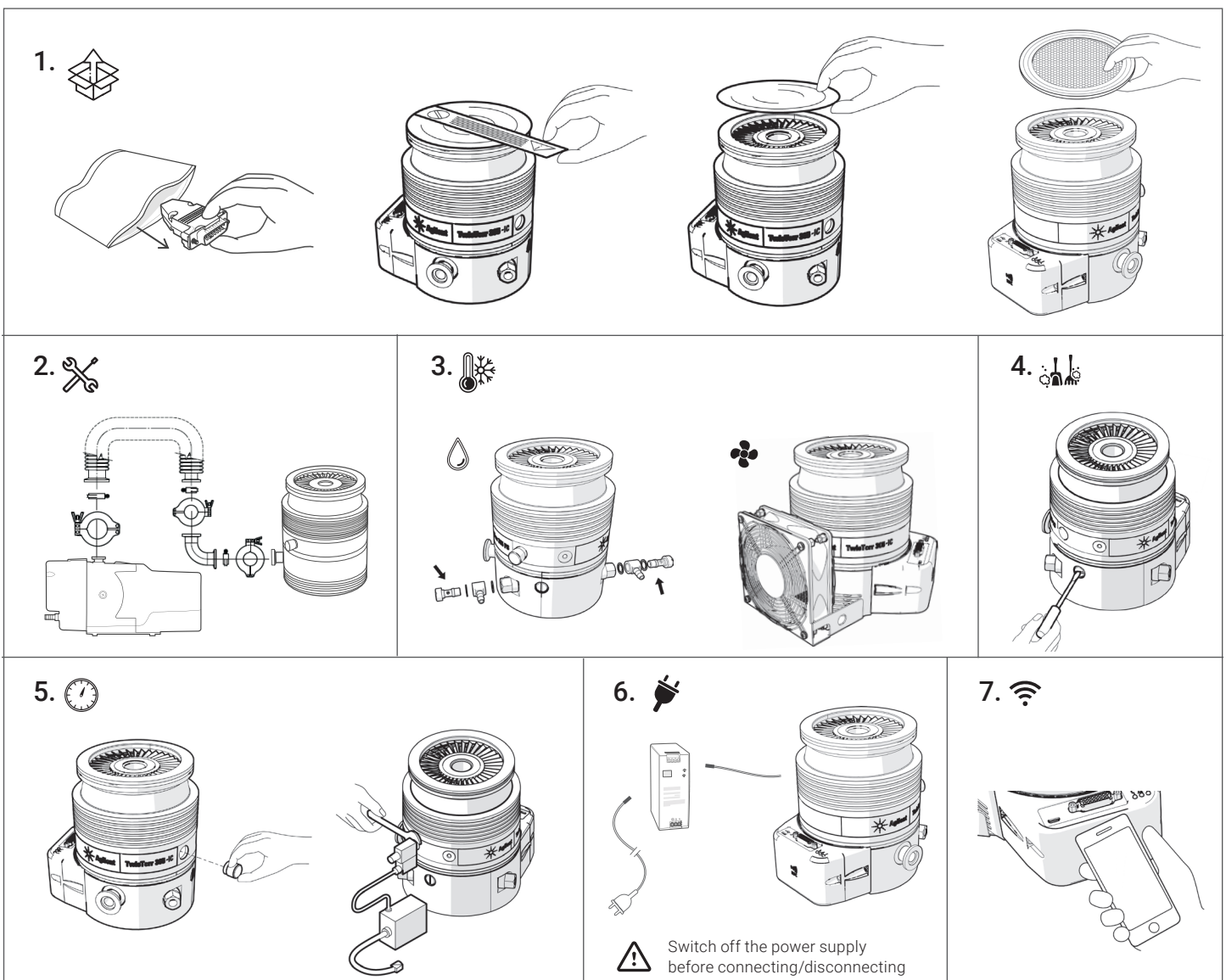
Cooling

 Water Cooling

 Air Cooling



TwisTorr 305-IC Installation



Visit the product web page:
www.agilent.com/chem/TwisTorr305-IC

This information is subject to change without notice.

© Agilent Technologies, Inc. 2020
Printed in Italy, August 3, 2020
X3513-90000 rev. B.00