

IDP-35 & IDP-45 UNBOXING AND INSTALLATION



UNBOXING



Quick Reference Guide

INSTALLATION (WITH OPTIONAL FILTERS)

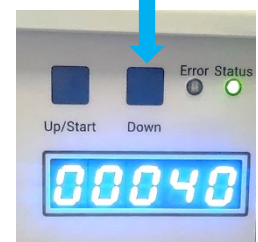
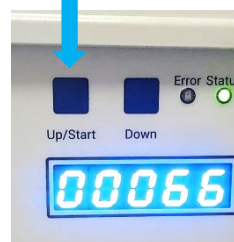


DEFAULT OPERATING MODE

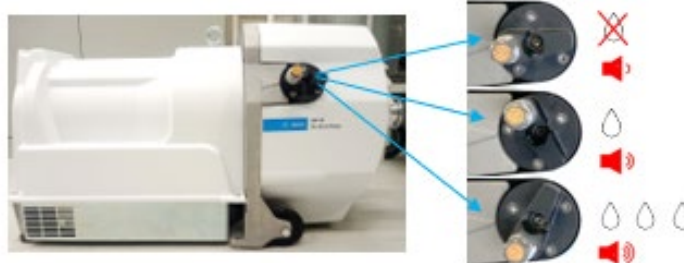
STARTING THE PUMP



SPEED TUNING



GAS BALLAST



Visit the product web page:

[IDP-35 Dry Scroll Pump](#)
[IDP-45 Dry Scroll Pump](#)

This information is subject to change without notice.

© Agilent Technologies, Inc. 2024
Printed in Italy, December 2024
X3835-90011 (A.00)



Get the IDP-35, IDP-45 Quick Reference Guide