

Roughing Pump Service Options: New Guidelines for Certain Models Built After November 2025

Starting November 2025, the service processes for the below models have changed. Minor maintenance procedures, as described in the following documents, now require revised steps and different tools.

For more information on changes to the service processes, please contact the Agilent Contact Center.

DS Rotary Vane Pumps

Models:

9499315, 9499320, 9499325, 9499330, 9499331, 9499335, 9499336

Documents:

8790093701 | DS102/202/302/402/602 user manual

TriScroll 300 Dry Scroll Pumps

Models:

PTS03001UNIVEU, PTS03001UVPIEU, PTS03001UNIVUK, PTS03001UVPIUK, PTS03001INV

Documents:

699904280 | TriScroll 300 Tip Seal Replacement user manual

699904285 | TriScroll 300 Pump Module Replacement user manual

699904260 | TriScroll 300 Major Maintenance user manual

TriScroll 600 Dry Scroll Pumps

Models:

PTS06001UNIVEU, PTS06001UVPIEU, PTS06001UNIVUK, PTS06001UVPIUK, PTS06001INV

Documents:

699904310 | TriScroll 600 Tip Seal Replacement user manual

699904305 | TriScroll 600 Pump Module Replacement user manual

699904300 | TriScroll 600 Major Maintenance user manual

www.agilent.com

D0140973 – A.00

This information is subject to change without notice.

© Agilent Technologies, Inc. 2025

November, 2025