

ICP-OES Oil-free Air Compressor



PerkinElmer is expanding its product portfolio in response to customer requests for an industrial quiet oil-free compressor, a must for dependable and clean air supply. Each compressor comes assembled on an internally and externally coated air tank and includes a self-purging regenerative dryer with aftercooler. One version includes a sound enclosure and has a sound level of less than 60 dB. Models in 115 V - 60 Hz and 230 V - 60 Hz are equipped with ASME® rated pressure tanks. Line Pressure Regulator with gauge and Line Ball Valve are standard on all units.



Specifications

Horse Power	1.5 Hp
Output CFM	6.0 CFM
Output L.	170 L. Min
Max Pressure PSI	120 PSI
Max Pressure Bar	8 bar
Operating Pressure PSI	90-120 PSI
Operating Pressure Bar	6-8 bar
Noise Level	75 dB/A, with Enclosure 55 dB/A
Tank Size Gal.	13 Gallons
Tank Size L	50 Liters
Dimensions	16 x 26 x 35 / with Enclosure 29 x 23 x 30
Weight	119 Lbs. / with Enclosure 207 Lbs.
Packed Dimensions	33 x 28 x 43 / with Enclosure 33 x 28 x 39
Packed Weight	135 Lbs. / with Enclosure 264 Lbs.

Compressors

Description	115V/60Hz	220V/50Hz	220V/60Hz
Compressor	N0777602	N0777603	N0777604
Compressor with Enclosure	N0777605	N0777606	N0777607

Product meets both U.S. and Canadian CSA standards. ASME® Certified tank.

Please note: Enclosure cannot be sold separately.

Tested and approved.

Replacement parts

Description	Part No.
Replacement Air Intake Filter	N0777608
Micron Dryer Filter Element	N0777609
Replacement Piston Assembly	N0777610
Replacement Reed Valve	N0777611
Replacement Head Gasket	N0777612

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright ©2010, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.